

**Listing of antibodies printed on the Kinex™ KAM-2025 antibody microarray.**



[www.kinexus.ca](http://www.kinexus.ca)

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
14-3-3 (KCIP-1)	Pan	P31946											●					Kinexus	NN441-3	RpAb
14-3-3 (KCIP-1)	Pan	P31946											●					Kinexus	NN441-2	RpAb
14-3-3 (KCIP-1)	Pan	P31946											●					No	sc-1657	MmAb
14-3-3-F (YWHAH)	Pan	Q04917											●					Kinexus	NN450-1	RpAb
14-3-3-G (YWHAG)	Pan	P61981											●					Kinexus	NN453-2	RpAb
14-3-3-S (YWHAS; SFN)	Pan	P31947											●					Kinexus	NN454-2	RpAb
14-3-3-S (YWHAS; SFN)	Pan	P31947											●					Kinexus	NN454-1	RpAb
14-3-3-T (YWHAQ)	Pan	P27348											●					Kinexus	NN455-2	RpAb
14-3-3-T (YWHAQ)	Pan	P27348											●					Kinexus	NN455-1	RpAb
14-3-3-Z (YWHAZ)	Pan	P63104											●					Kinexus	NN001-3	RpAb
14-3-3-Z (YWHAZ)	Pan	P63104											●					Kinexus	NN001-2	RpAb
14-3-3b (YWHAB)	Pan	P31946											●					Kinexus	NN452-2	RpAb
14-3-3b (YWHAB)	Pan	P31946											●					Kinexus	NN452-1	RpAb
14-3-3e (YWHAE)	Pan	P62258											●					Kinexus	NN451-3	RpAb
14-3-3e (YWHAE)	Pan	P62258											●					Kinexus	NN451-1	RpAb
4E-BP1 (PHAS1)	Pan	Q13541						●										No	Y021215	RpAb
4E-BP1 (PHAS1)	Pan	Q13541						●										No	sc-6024	GpAb
4E-BP1 (PHAS1)	T37+T46	Q13541						●										Kinexus	PN550	RpAb
A6 (Twinfilin-1)	Y309	Q12792	• DSK										●					Kinexus	PK501	RpAb
A6r (Twinfilin-2)	Y309	Q61BS0	• DSK										●					Kinexus	PK502	RpAb
AAK1	S637+Pan	Q2M218	• PSTK															Kinexus	PK503	RpAb
Abl (Abl1)	Pan	P00519	• PYK															Kinexus	NK001-4	RpAb
Abl (Abl1)	Pan	P00519	• PYK															Kinexus	NK001-6	RpAb
Abl (Abl1)	Pan	P00519	• PYK															No	sc-887	RpAb
Abl (Abl1)	Pan	P00519	• PYK															No	sc-131	RpAb
Abl (Abl1)	Y139	P00519	• PYK															Kinexus	PK504	RpAb
Abl (Abl1)	Y226	P00519	• PYK															Kinexus	PK505	RpAb
Abl (Abl1)	Y257	P00519	• PYK															Kinexus	PK506	RpAb
Abl (Abl1)	Y264	P00519	• PYK															Kinexus	PK507	RpAb
Abl (Abl1)	Y393+T394	P00519	• PYK															Kinexus	PK873	RpAb
Abl (Abl1)	Y413	P00519	• PYK															Kinexus	PK896	RpAb
Abl (Abl1)	Y469	P00519	• PYK															Kinexus	PK508	RpAb
ACACA (ACC1; ACCA)	S29	Q13085								●								Kinexus	PN733	RpAb
ACACA (ACC1; ACCA)	S80	Q13085								●								Kinexus	PN863	RpAb
ACACA (ACC1; ACCA)	S80	Q13085								●								No	07-303	RpAb
ACACB (ACC2)	T1342	O00763								●								Kinexus	PN734	RpAb
AChE	Y164	P22303								●								Kinexus	PN931	RpAb
ACK1 (TNK2)	Pan	Q07912	• PYK															No	sc-323	RpAb
ACK1 (TNK2)	Y284	Q07912	• PYK															Kinexus	PK511	RpAb
ACK1 (TNK2)	Y518	Q07912	• PYK															Kinexus	PK512	RpAb
ACK1 (TNK2)	Y859+Y860	Q07912	• PYK															Kinexus	PK513	RpAb
ACLY	S455	P53396								●								Kinexus	PN735	RpAb
ACLY	Y682	P53396								●								Kinexus	PN686	RpAb
ACO1 IREB1)	S711+Pan	P21399								●								Kinexus	PN736	RpAb
ACO1 IREB1)	S806	P21399								●								Kinexus	PN737	RpAb
ACP1	Y132+Y133	P24666	• PSTP															Kinexus	PN687	RpAb
ACS1 (ACSL1)	Y567	P33121								●								Kinexus	PN688	RpAb
ACTA1 (alpha-actin)/ACTB	Pan	P68133												●				No	sc-8432	MmAb
ACTB (beta-actin)	Y53	P60709												●				Kinexus	PN501	RpAb
ACTB (beta-actin)	Y294	P60709												●				Kinexus	PN500	RpAb
ACTN1	Y246	P12814												●				Kinexus	PN502	RpAb
ADAM22	Pan	Q9P0K1				●												No	NN189-2	MmAb
ADD1 (Adducin-a)	Y35	Q9UEY8												●				Kinexus	PN867	RpAb
ADD1 (Adducin-a)	Y550	Q9UEY8												●				Kinexus	PN868	RpAb
ADD1 (Adducin-a)	T724+S726	Q9UEY8												●				Kinexus	PN866	RpAb
ADD1 (Adducin-a)	S726	Q9UEY8												●				Kinexus	PN807	RpAb
ADK	S86	P55263								●								Kinexus	PN689	RpAb
ADRA2C (ADRA2L2;	Pan	P18825				●												No	NN190-2	MmAb
ADRA2C (ADRA2L2;	Pan	P18825				●												No	NN190-1	MmAb
ADRB2	S355+S356	P07550				●												Kinexus	PN784	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
AGTR1 (AT1)	S335+T336+S338	P30556				●												Kinexus	PN785	RpAb
AHSA1 (AHA1)	Pan	O95433									●							No	NN191-2	RatmAb
AIF	Pan	O95831										●						No	sc-9416	GpAb
AKAP12	S627+S629	Q02952											●					Kinexus	PN1009	RpAb
AKAP12	S696	Q02952											●					Kinexus	PN933	RpAb
Akt1 (PKBa)	Pan	P31749	• PSTK															Kinexus	NK129-4	RpAb
Akt1 (PKBa)	Pan	P31749	• PSTK															Kinexus	NK129-5	RpAb
Akt1 (PKBa)	Pan	P31749	• PSTK															Kinexus	NK129-3	RpAb
Akt1 (PKBa)	Pan	P31749	• PSTK															No	sc-5298	MmAb
Akt1 (PKBa)	T308	P31749	• PSTK															Kinexus	PK515	RpAb
Akt1 (PKBa)	Y315	P31749	• PSTK															Kinexus	PK516	RpAb
Akt1 (PKBa)	Y326	P31749	• PSTK															Kinexus	PK517	RpAb
Akt1 (PKBa)	S473	P31749	• PSTK															Kinexus	PK869	RpAb
Akt1 (PKBa)	S473	P31749	• PSTK															Kinexus	PK958	RpAb
Akt1 (PKBa)	Y474	P31749	• PSTK															No	06-0020	RpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															Kinexus	NK130-4	RpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															Kinexus	NK130-8	RpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															Kinexus	NK130-9	RpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															Kinexus	NK130-5	RpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															No	sc-7125	GpAb
Akt2 (PKBb)	Pan	P31751	• PSTK															No	sc-7127	GpAb
Akt3 (PKBg)	Pan	Q9Y243	• PSTK															Kinexus	NK131-3	RpAb
ALK	Pan	Q9UM73	• PYK			●												Kinexus	NK003-2	RpAb
ALK	Pan	Q9UM73	• PYK			●												Kinexus	NK003-3	RpAb
ALK	Pan	Q9UM73	• PYK			●												No	18-0266	RpAb
ALK	Y1092	Q9UM73	• PYK			●												Kinexus	PK518	RpAb
ALK	Y1096	Q9UM73	• PYK			●												Kinexus	PK519	RpAb
ALK	Y1507	Q9UM73	• PYK			●												Kinexus	PK520	RpAb
ALOX5 (5-LO)	S272	P09917								●								Kinexus	PN738	RpAb
ALS2 (ALS2CR6)	S483	Q96Q42																Kinexus	PN869	RpAb
ALS2 (ALS2CR6)	S492	Q96Q42																Kinexus	PN870	RpAb
AML1 (RUNX1)	T273	Q01196			●													Kinexus	PN568	RpAb
AML2 (RUNX3)	T231	Q13761			●													Kinexus	PN569	RpAb
AMPKa1 (PRKAA1)	Pan	Q13131	• PSTK															Kinexus	NK259-2	RpAb
AMPKa1 (PRKAA1)	Pan	Q13131	• PSTK															Kinexus	NK259-1	RpAb
AMPKa1 (PRKAA1)	T172	Q13131	• PSTK															Kinexus	PK897	RpAb
AMPKa1 (PRKAA1)	T183+S184	Q13131	• PSTK															Kinexus	PK521	RpAb
AMPKa1 (PRKAA1)	T490	Q13131	• PSTK															Kinexus	PK918	RpAb
AMPKa2 (PRKAA2)	Pan	P54646	• PSTK															Kinexus	NK260-1	RpAb
AMPKa2 (PRKAA2)	Pan	P54646	• PSTK															Kinexus	NK260-2	RpAb
AMPKa2 (PRKAA2)	S377	P54646	• PSTK															Kinexus	PK522	RpAb
AMPKa2 (PRKAA2)	Y436	P54646	• PSTK															Kinexus	PK919	RpAb
ANAX11	Y482	P50995														●		Kinexus	PN871	RpAb
ANKRD3 (RIPK4)	Pan	P57078	• PSTK															No	AP7808b	RpAb
ANKRD3 (RIPK4)	S438+Pan	P57078	• PSTK															Kinexus	PK523	RpAb
Ankyrin (ANK1)	Pan	P16157											●					No	NN195-2	MmAb
Ankyrin B (ANK2)	Pan	Q01484											●					No	NN196-1	MmAb
ANXA1	Y207	P04083														●		Kinexus	PN551	RpAb
ANXA1	Y207	P04083														●		Kinexus	PN503	RpAb
ANXA2	Y238	P07355														●		Kinexus	PN504	RpAb
AP2B1 (ADTB2)	Y276	P63010															●	Kinexus	PN934	RpAb
AP2M1 (AP-2 mu 2)	T156	Q96CW1											●				●	Kinexus	PN935	RpAb
AP3B1	S276	O00203															●	Kinexus	PN936	RpAb
AP3B2	S272	Q13367															●	Kinexus	PN937	RpAb
APC	S2129	P25054											●					Kinexus	PN739	RpAb
APG1	Pan	O95757									●							No	sc-6242	RpAb
APG2	Pan	P34932									●							No	sc-6240	RpAb
APP (A4; ABPP; Amyloid)	Pan	P05067			●	●					●							Kinexus	NN192-2	RpAb
APP (A4; ABPP; Amyloid)	Pan	P05067			●	●					●							Kinexus	NN192-3	RpAb
APP (A4; ABPP; Amyloid)	Pan	P05067			●	●					●							Kinexus	NN192-4	RpAb
APP (A4; ABPP; Amyloid)	Pan	P05067			●	●					●							No	NN194-1	RpAb
APP (A4; ABPP; Amyloid)	T743	P05067			●	●					●							Kinexus	PN808	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
APP (A4; ABPP; Amyloid)	Y757	P05067			●	●					●							Kinexus	PN809	RpAb
Aquaporin 3 (AQP3)	Pan	Q92482					●											No	NN201-1	RpAb
AR	S96	P10275			●	●												Kinexus	PN570	RpAb
AR	S310	P10275			●	●												Kinexus	PN571	RpAb
Arg (Abl2)	Pan	P42684	• PYK															Kinexus	NK238-1	RpAb
Arg (Abl2)	Y439	P42684	• PYK															Kinexus	PK509	RpAb
Arg (Abl2)	Y439+T440	P42684	• PYK															Kinexus	PK510	RpAb
ARID1A	S363	O14497			●													Kinexus	PN740	RpAb
ARRB1	Pan	P49407											●					No	610551	MmAb
ARRB1	S412	P49407											●					No	44-200	RpAb
ASK1 (MAP3K5)	Pan	Q99683	• PSTK															No	sc-6368	GpAb
ASK1 (MAP3K5)	Pan	Q99683	• PSTK															No	sc-7931	RpAb
ASK1 (MAP3K5)	T838	Q99683	• PSTK															Kinexus	PK525	RpAb
ASK1 (MAP3K5)	S1033	Q99683	• PSTK															Kinexus	PK524	RpAb
ASK1 (MAP3K5)	T1326	Q99683	• PSTK															Kinexus	PK920	RpAb
ASS1	T174+S180	P00966								●								Kinexus	PN741	RpAb
ATAD1 ATPase	Pan	Q8NBU5																No	NN204-1	MmAb
ATASE (PPAT)	T356	Q06203								●								Kinexus	PN742	RpAb
ATF2 (CRE-BP1)	T69+T71	P15336			●													Kinexus	PN572	RpAb
ATF2 (CRE-BP1)	T69+T71	P15336			●													Kinexus	PN552	RpAb
ATF2 (CRE-BP1)	T71	P15336			●													No	sc-8398	MmAb
ATG2A	Pan	Q2TAZ0							●									No	NN209-1	RpAb
ATG4B	Pan	Q9Y4P1							●									No	NN374-1	RpAb
ATG4C (APG4C; AURL1;	Pan	Q96DT6							●									No	NN210-2	RpAb
ATG4C (APG4C; AURL1;	Pan	Q96DT7							●									No	NN210-1	RpAb
ATG4D (PG4D; AURL4)	Pan	Q86TL0							●									No	NN211-2	RpAb
ATG5 (APG5L; ASP)	Pan	Q9H1Y0							●									No	NN212-1	RpAb
ATG12 (APG12; APG12L)	Pan	O94817							●									No	NN213-1	RpAb
ATG13	Pan	O75143							●									No	NN214-1	RpAb
ATM	Pan	Q13315	• PSTK															Kinexus	NK230-2	RpAb
ATM	Pan	Q13315	• PSTK															Kinexus	NK230-1	RpAb
ATM	Pan	Q13315	• PSTK															Kinexus	NK230-3	RpAb
ATM	Pan	Q13315	• PSTK															Kinexus	NK230	RpAb
ATM	Pan	Q13315	• PSTK															No	sc-7230	RpAb
ATM	Pan	Q13315	• PSTK															No	sc-1214	GpAb
ATM	S1981	Q13315	• PSTK															Kinexus	PK526	RpAb
ATM	Y2969	Q13315	• PSTK															Kinexus	PK527	RpAb
ATP7B	Pan	P35670													●			No	NN215-2	MmAb
ATR	Pan	Q13535	• PSTK															Kinexus	NK237-2	RpAb
ATR	Pan	Q13535	• PSTK															Kinexus	NK237-3	RpAb
ATR	Pan	Q13535	• PSTK															Kinexus	NK237-1	RpAb
ATR	S435+S436	Q13535	• PSTK															Kinexus	PK528	RpAb
AurKA (Aurora A; ARK-1; AIK;	Pan	O14965	• PSTK															Kinexus	NK008-4	RpAb
AurKA (Aurora A; ARK-1; AIK;	Pan	O14965	• PSTK															Kinexus	NK008-5	RpAb
AurKA (Aurora A; ARK-1; AIK;	Pan	O14965	• PSTK															Kinexus	NK008-3	RpAb
AurKA (Aurora A; ARK-1; AIK;	T287+T288	O14965	• PSTK															Kinexus	PK529	RpAb
AurKB (AurB; AIK2)	Pan	Q96GD4	• PSTK															Kinexus	NK193-3	RpAb
AurKB (AurB; AIK2)	Pan	Q96GD4	• PSTK															Kinexus	NK193-4	RpAb
AurKB (AurB; AIK2)	Pan	Q96GD4	• PSTK															Kinexus	NK193-2	RpAb
AurKB (AurB; AIK2)	S227	Q96GD4	• PSTK															Kinexus	PK530	RpAb
AurKB (AurB; AIK2)	T232	Q96GD4	• PSTK															Kinexus	PK531	RpAb
AurKC (Aurora C; AurC; AIK3)	Pan	Q9UQB9	• PSTK															Kinexus	NK009-3	RpAb
AurKC (Aurora C; AurC; AIK3)	Pan	Q9UQB9	• PSTK															Kinexus	NK009-4	RpAb
AurKC (Aurora C; AurC; AIK3)	Pan	Q9UQB9	• PSTK															Kinexus	NK009-2	RpAb
AurKC (Aurora C; AurC; AIK3)	S193	Q9UQB9	• PSTK															Kinexus	PK532	RpAb
AVPR1A	T378+S380	P37288				●												Kinexus	PN787	RpAb
Axl (UFO)	Pan	P30530	• PYK			●												Kinexus	NK010-3	RpAb
Axl (UFO)	Pan	P30530	• PYK			●												Kinexus	NK010-2	RpAb
Axl (UFO)	Y702+Y703	P30530	• PYK			●												Kinexus	PK533	RpAb
Axl (UFO)	Y779	P30530	• PYK			●												Kinexus	PK898	RpAb
B-Myb (MYBL2)	T487	P10244			●													Kinexus	PN573	RpAb
B23 (Nucleophosmin; NPM)	Pan	P06748																No	sc-6013-R	RpAb
Bad	S99	Q92934										●						No	07-007	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Bad	S99	Q92934										●						No	AB10424	RpAb
Bak	Pan	Q16611										●						No	sc-832	RpAb
BCKD (BCKDK)	Pan	O14874	• PSTK															Kinexus	NK257-2	RpAb
BCKD (BCKDK)	Pan	O14874	• PSTK															Kinexus	NK257-3	RpAb
BCKD (BCKDK)	Pan	O14874	• PSTK															Kinexus	NK257-1	RpAb
BCKDHA	S337+Pan	P12694								●								Kinexus	PN810	RpAb
Bcl-xL	Pan	Q07817										●						No	DB002	RpAb
Bcl2	S70	P10415										●						No	sc-293128	MmAb
Bcl2	S70+Pan	P10415										●						Kinexus	PN811	RpAb
BCLAF1	S512	Q9NYF8			●													Kinexus	PN574	RpAb
Bcr	Y177	P11274	• PSTK															Kinexus	PK538	RpAb
Bcr	Y591	P11274	• PSTK															Kinexus	PK539	RpAb
Bcr	Y644	P11274	• PSTK															Kinexus	PK540	RpAb
BDNF (Abrineurin)	Pan	P23560																No	NN216-1	RpAb
Beclin 1 (BECN1; GT197)	Pan	Q14457											●					No	NN217-1	RpAb
Beclin 2 (BECN1L1)	Pan	A8MW95											●					No	NN372-1	RpAb
Bid	Pan	P55957										●						No	sc-6539	GpAb
BK Beta2 (KCNMB2)	Pan	Q9Y691					●											No	NN219-1	MmAb
BKR2	S366+T369	P30411				●												Kinexus	PN788	RpAb
BLK	Pan	P51451	• PYK															No	AP7697a	RpAb
BLK	Y187	P51451	• PYK															Kinexus	PK541	RpAb
BLK	Y188	P51451	• PYK															Kinexus	PK542	RpAb
BLK	Y389	P51451	• PYK															Kinexus	PK543	RpAb
BLVRA (Billiverdin-IX alpha-	Pan	P53004								●								No	NN220-1	RpAb
BMPR2 (BMPR-II)	S375	Q13873				●												Kinexus	PK544	RpAb
Bmx (Etk)	Pan	P51813	• PYK															No	B63620	RpAb
Bmx (Etk)	Y40	P51813	• PYK															Kinexus	PK545	RpAb
BRCA1	T509	P38398			●													Kinexus	PN690	RpAb
BRCA2	S70	P51587											●					Kinexus	PN789	RpAb
BRD2	S37	P25440	• PSTK															Kinexus	PK546	RpAb
BRK	Y342	Q13882	• PYK															Kinexus	PK548	RpAb
BRK	S446+Y447	Q13882	• PYK															Kinexus	PK547	RpAb
BRSK1	T205	Q8TDC3	• PSTK															Kinexus	PK549	RpAb
Btk	Pan	Q06187	• PYK															No	sc-1696	RpAb
Btk	Pan	Q06187	• PYK															No	sc-1107	GpAb
Btk	Y223+Y225	Q06187	• PYK															Kinexus	PK550	RpAb
Btk	Y551	Q06187	• PYK															Kinexus	PK551	RpAb
BUB1 (BUB1A)	Pan	Q43683	• PSTK															No	AP8058b	RpAb
BUBR1 (BUB1B)	S670+Pan	O60566	• PSTK															Kinexus	PK552	RpAb
C-EBP-alpha	T226+T230	P49715			●													Kinexus	PN575	RpAb
CACNB1 (CACB1)	T499	Q02641					●											Kinexus	PN938	RpAb
CACNG2 (Stargazin; TARP	Pan	Q9Y698					●											No	NN224-2	MmAb
CACNG2 (Stargazin; TARP	Pan	Q9Y699					●											No	NN224-1	MmAb
CACNG3 (TARP3)	Pan	O60359					●											Kinexus	NN456	RpAb
CAD	S1859	P27708								●								Kinexus	PN691	RpAb
Calnexin (CANX)	Pan	P27824														●		No	NN136-2	RpAb
Calnexin (CANX)	Pan	P27824														●		No	NN136-5	RpAb
CaMK1a (CAMK1)	T177	Q14012	• PSTK															Kinexus	PK553	RpAb
CaMK1a (CAMK1)	Y235	Q14012	• PSTK															Kinexus	PK874	RpAb
CaMK1a/d (CaMKI)	Pan	Q8IU85	• PSTK															No	sc-1544	GpAb
CaMK1d	Pan	Q8IU85	• PSTK															No	sc-1543	GpAb
CaMK1d	T180	Q8IU85	• PSTK															Kinexus	PK554	RpAb
CaMK2a (CaMKII)	Pan	Q9UQM7	• PSTK															No	sc-9035	RpAb
CaMK2a (CaMKII)	Pan	Q9UQM7	• PSTK															No	sc-5390	GpAb
CaMK2a (CaMKII)	Pan	Q9UQM7	• PSTK															No	NK302-2	MmAb
CaMK2a (CaMKII)	Pan	Q9UQM7	• PSTK															No	NK302-1	MmAb
CaMK2a (CaMKII)	Y230	Q9UQM7	• PSTK															Kinexus	PK875	RpAb
CaMK2a (CaMKII)	T286	Q9UQM7	• PSTK															Kinexus	PK555	RpAb
CAMK2b	Pan	Q13554	• PSTK															Kinexus	NK018-3	RpAb
CAMK2d	Pan	Q13557	• PSTK															Kinexus	NK019-3	RpAb
CaMK2g	Pan	Q13555	• PSTK															Kinexus	NK020-1	RpAb
CaMK4 (CaMPK4)	Pan	Q16566	• PSTK															Kinexus	NK021-1	RpAb
CaMK4 (CaMPK4)	Pan	Q16566	• PSTK															Kinexus	NK021-2	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
CaMK4 (CaMPK4)	T200	Q16566	• PSTK															Kinexus	PK556	RpAb
CaMKK1 (CaMKK)	Pan	Q8N5S9	• PSTK															Kinexus	NK022-2	RpAb
CaMKK1 (CaMKK)	Pan	Q8N5S9	• PSTK															Kinexus	NK022	RpAb
CaMKK1 (CaMKK)	S74	Q8N5S9	• PSTK															Kinexus	PK557	RpAb
Cas-L	Y166+Pan	Q14511											●					Kinexus	PN505	RpAb
CASP1 (Caspase-1)	Pan	P29466										●						No	sc-622	RpAb
CASP3 (Caspase-3; Yama)	Pan	P42574										●						No	160745	RpAb
CAV1	Pan	Q03135											●					No	Y021112	RpAb
Cav1.2 (CACNA1C)	Pan	Q13936					●											No	NN225-1	MmAb
Cav1.3 (CACNA1D; CACH3)	Pan	Q01668					●											No	NN226-1	MmAb
CavBeta2 (CACNB2; CAB2)	Pan	Q08289					●											No	NN228-1	MmAb
Cbl	Y674	P22681											●					Kinexus	PN743	RpAb
CBS	S227+Pan	P35520								●								Kinexus	PN744	RpAb
CD45 (PTPRC)	Y1216	P08575		• PYP		●												Kinexus	PP527	RpAb
CD74	Pan	P04233																No	NN229-1	MmAb
CDC25A	Pan	P30304		• DSP														Kinexus	NP038-2	RpAb
CDC25A	Pan	P30304		• DSP														Kinexus	NP038-3	RpAb
CDC25A	Pan	P30304		• DSP														Kinexus	NP038-1	RpAb
CDC25B	Pan	P30305		• DSP														Kinexus	NP002-2	RpAb
CDC25B	Pan	P30305		• DSP														Kinexus	NP002-3	RpAb
CDC25C	Pan	P30307		• DSP														Kinexus	NP003-2	RpAb
CDC25C	Pan	P30307		• DSP														Kinexus	NP003-3	RpAb
CDC25C	Pan	P30307		• DSP														No	sc-5620	RpAb
CDC2L5 (CHED; CDK13)	Pan	Q14004	• PSTK															No	AP7514a	RpAb
CDC2L5 (CHED; CDK13)	Pan	Q14004	• PSTK															Kinexus	NK024-2	RpAb
CDC37	Pan	Q16543											●					No	NN230-1	RpAb
CDC5L	T385	Q99459			●													Kinexus	PN576	RpAb
CDC7	Pan	O00311	• PSTK															No	sc-7519	GpAb
CDC7	Pan	O00311	• PSTK															No	sc-7518	GpAb
CDC7	T376	O00311	• PSTK															Kinexus	PK558	RpAb
CDK1 (CDC2)	Pan	P06493	• PSTK															No	AHZ0122	MmAb
CDK1 (CDC2)	Pan	P06493	• PSTK															Kinexus	NK025-4	RpAb
CDK1 (CDC2)	Pan	P06493	• PSTK															Kinexus	NK025-7	RpAb
CDK1 (CDC2)	Pan	P06493	• PSTK															No	sc-954	RpAb
CDK1 (CDC2)	Pan	P06493	• PSTK															No	sc-54	MmAb
CDK1 (CDC2)	T14	P06493	• PSTK															No	2543	RpAb
CDK1 (CDC2)	T14	P06493	• PSTK															Kinexus	PK559	RpAb
CDK1 (CDC2)	T14+Y15	P06493	• PSTK															Kinexus	PK560	RpAb
CDK1 (CDC2)	Y15	P06493	• PSTK															Kinexus	PK562	RpAb
CDK1 (CDC2)	Y19	P06493	• PSTK															Kinexus	PK563	RpAb
CDK1 (CDC2)	T161	P06493	• PSTK															Kinexus	PK561	RpAb
CDK1 (CDC2)	T161	P06493	• PSTK															No	11134	RpAb
CDK2	Pan	P24941	• PSTK															No	AP7518b	RpAb
CDK2	Pan	P24941	• PSTK															No	sc-163	RpAb
CDK2	Pan	P24941	• PSTK															No	sc-748	RpAb
CDK2	Pan	P24941	• PSTK															No	sc-6248	MmAb
CDK2	Pan	P24941	• PSTK															No	610145	MmAb
CDK2	T160	P24941	• PSTK															Kinexus	PK568	RpAb
CDK4	Pan	P11802	• PSTK															No	sc-260	RpAb
CDK4	Pan	P11802	• PSTK															No	sc-601	RpAb
CDK4	T172	P11802	• PSTK															Kinexus	PK569	RpAb
CDK5	Pan	Q00535	• PSTK															Kinexus	NK028-2	RpAb
CDK5	Pan	Q00535	• PSTK															No	sc-6247	MmAb
CDK5	Y15	Q00535	• PSTK															Kinexus	PK570	RpAb
CDK5	S159+Pan	Q00535	• PSTK															Kinexus	PK899	RpAb
CDK6	Pan	Q00534	• PSTK															No	C150M	MmAb
CDK6	Pan	Q00534	• PSTK															No	sc-7961	MmAb
CDK6	Pan	Q00534	• PSTK															No	sc-177	RpAb
CDK6	Pan	Q00534	• PSTK															No	AP-CC006	RpAb
CDK6	Y13	Q00534	• PSTK															Kinexus	PK571	RpAb
CDK6	Y24	Q00534	• PSTK															Kinexus	PK572	RpAb
CDK7 (MO15)	Pan	P50613	• PSTK															Kinexus	NK030-1	RpAb
CDK7 (MO15)	Pan	P50613	• PSTK															No	sc-7344	MmAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
CDK7 (MO15)	Pan	P50613	• PSTK															No	13-8700	RpAb
CDK7 (MO15)	T170	P50613	• PSTK															Kinexus	PK573	RpAb
CDK8	Pan	P49336	• PSTK															No	AP7524a	RpAb
CDK8	Pan	P49336	• PSTK															Kinexus	NK031-4	RpAb
CDK8	Pan	P49336	• PSTK															No	sc-1521	GpAb
CDK9	Pan	P50750	• PSTK															No	sc-7331	GpAb
CDK9	Pan	P50750	• PSTK															No	sc-484	RpAb
CDK9	T186	P50750	• PSTK															Kinexus	PK575	RpAb
CDK9	S347	P50750	• PSTK															Kinexus	PK574	RpAb
CDK10 (PISSLRE)	Pan	Q15131	• PSTK															Kinexus	NK033-3	RpAb
CDK10 (PISSLRE)	Pan	Q15131	• PSTK															Kinexus	NK033-2	RpAb
CDK10 (PISSLRE)	T196	Q15131	• PSTK															Kinexus	PK564	RpAb
CDK11A (Cdc2L2; PITSLRE)	T583	Q9UQ88	• PSTK															Kinexus	PK565	RpAb
CDK11B (PITSLRE; p58/GTA;)	Pan	P21127	• PSTK															No	sc-928	RpAb
CDK12 (Cdc2L7; CRK7)	S383+S385	Q9NYV4	• PSTK															Kinexus	PK566	RpAb
CDK12 (Cdc2L7; CRK7)	T893	Q9NYV4	• PSTK															Kinexus	PK567	RpAb
CDK14 (PFTAIRE1)	Pan	O94921	• PSTK															Kinexus	NK287-1	RpAb
CDK15 (PFTAIRE2;	Pan	Q96Q40	• PSTK															No	AP7512a	RpAb
CDK15 (PFTAIRE2;	Pan	Q96Q40	• PSTK															Kinexus	NK004-2	RpAb
CDK15 (PFTAIRE2;	Pan	Q96Q40	• PSTK															Kinexus	NK004-3	RpAb
CDK15 (PFTAIRE2;	Y114	Q96Q40	• PSTK															Kinexus	PK957	RpAb
CDKL1 (KKIALRE)	Pan	Q00532	• PSTK															Kinexus	NK199	RpAb
CDKL1 (KKIALRE)	Pan	Q00532	• PSTK															No	sc-7588	GpAb
CDKL1 (KKIALRE)	Pan	Q00532	• PSTK															No	sc-7589	GpAb
CDKL2 (KKIAMRE)	Pan	Q92772	• PSTK															Kinexus	NK261-1	RpAb
CDKL2 (KKIAMRE)	Pan	Q92772	• PSTK															No	sc-6287	GpAb
CDKL2 (KKIAMRE)	Pan	Q92772	• PSTK															No	sc-6286	GpAb
CDKL3	Pan	Q8IVW4	• PSTK															Kinexus	NK262-1	RpAb
CDKL5	Y171	O76039	• PSTK															Kinexus	PK576	RpAb
CDKN1A (p21; WAF1; Cip1;	Pan	P38936	• PSTK-															No	sc-397	RpAb
CDKN1B (p27; Kip1)	Pan	P46527	• PSTK-															No	sc-528	RpAb
CDKN2A (p14ARF; p19ARF)	Pan	P42771	• PSTK-															Kinexus	PN693	RpAb
CDKN2A (p16; p16-INK4a)	Pan	P42771	• PSTK-															Kinexus	PN694	RpAb
CDKN2A (p16; p16-INK4a)	Pan	P42771	• PSTK-															No	sc-9968	MmAb
CDKN2C (p18; p18 INK4c;	Pan	P42773	• PSTK-															No	sc-865	RpAb
CDKN2C (p18; p18 INK4c)	Pan	P42773	• PSTK-															No	AHZ0352	RpAb
CDKN2D (p19; p19 INK4d)	Pan	P55273	• PSTK-															No	sc-1063-G	GpAb
CENPA	Pan	P49450																No	NN231-1	MmAb
CHAK1 (TRPM7)	Pan	Q96QT4				●	●											No	AP8052b	RpAb
CHAK1 (TRPM7)	Pan	Q96QT4				●	●											No	AP8052a	RpAb
CHAK1 (TRPM7)	S1504	Q96QT4				●	●											Kinexus	PK921	RpAb
CHAK2 (TRPM6)	Pan	Q9BX84				●	●											No	AP8053c	RpAb
CHAK2 (TRPM6)	Pan	Q9BX84				●	●											No	AP8053b	RpAb
Chapsyn-110 (DLG2; PSD-93)	Pan	Q15700											●					No	NN232-1	MmAb
CHAT (Choline acetylase)	Pan	P28329								●								No	NN233-1	RpAb
Chk1 (CHEK1)	Pan	O14757	• PSTK															No	sc-7898	RpAb
Chk1 (CHEK1)	Pan	O14757	• PSTK															No	sc-8408	MmAb
Chk1 (CHEK1)	Pan	O14757	• PSTK															No	sc-7235	GpAb
Chk1 (CHEK1)	S280	O14757	• PSTK															Kinexus	PK577	RpAb
Chk1 (CHEK1)	S317	O14757	• PSTK															Kinexus	PK578	RpAb
Chk1 (CHEK1)	S345	O14757	• PSTK															Kinexus	PK579	RpAb
Chk2 (CHEK2)	Pan	O96017	• PSTK															No	sc-8812	GpAb
Chk2 (CHEK2)	Pan	O96017	• PSTK															No	sc-9064	RpAb
Chk2 (CHEK2)	T68	O96017	• PSTK															Kinexus	PK581	RpAb
Chk2 (CHEK2)	T383	O96017	• PSTK															Kinexus	PK580	RpAb
CHMP2B	S199	Q9UQN3																Kinexus	PN872	RpAb
CHRM1	S451+T455+	P11229				●												Kinexus	PN790	RpAb
CHRNA9	Y374	Q9UGM1				●												Kinexus	PN939	RpAb
CHRNA9	Y430	Q9UGM1				●												Kinexus	PN940	RpAb
CK1a (CSNK1A1)	Pan	P48729	• PSTK															No	AP7400a	RpAb
CK1a (CSNK1A1)	Pan	P48729	• PSTK															No	sc-6478	GpAb
CK1d (CSNK1D)	Pan	P48730	• PSTK															No	AP7402a	RpAb
CK1d (CSNK1D)	Pan	P48730	• PSTK															No	sc-6474	GpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
CK1d (CSNK1D)	Pan	P48730	• PSTK															No	sc-6475	GpAb
CK1d (CSNK1D)	T176	P48730	• PSTK															Kinexus	PK923	RpAb
CK1d (CSNK1D)	T347	P48730	• PSTK															Kinexus	PK924	RpAb
CK1d (CSNK1D)	S370	P48730	• PSTK															Kinexus	PK922	RpAb
CK1e (CSNK1E)	Pan	P49674	• PSTK															No	AP7403a	RpAb
CK1e (CSNK1E)	Pan	P49674	• PSTK															No	sc-6472	GpAb
CK1g3 (CSNK1G3)	Pan	Q9Y6M4	• PSTK															No	AP7406a	RpAb
CK2a1 (CSNK2A1)	Pan	P68400	• PSTK															Kinexus	NK041-2	RpAb
CK2a1 (CSNK2A1)	Pan	P68400	• PSTK															No	sc-373894	MmAb
CK2a1 (CSNK2A1)	Y255	P68400	• PSTK															Kinexus	PK582	RpAb
CK2a1 (CSNK2A1)	T360+S362	P68400	• PSTK															No	11572	RpAb
CK2b1 (Casein Kinase IIb)	Pan	P67870	• PSTK-															No	sc-6483	GpAb
CICN3 ( CIC-3; Clcn3)	Pan	P51790					●											No	NN234-1	MmAb
CLK1	Pan	P49759	• PSTK															No	AP7529a	RpAb
CLK1	S337	P49759	• PSTK															Kinexus	PK583	RpAb
CLK1	S337+T338	P49759	• PSTK															Kinexus	PK584	RpAb
CLK2	Pan	P49760	• PSTK															No	AP7530a	RpAb
CLK3	Pan	P49761	• PSTK															No	AP7531a	RpAb
CLK4	Pan	Q9HAZ1	• PSTK															No	AP7532a	RpAb
CLTC	Y634	Q00610															●	Kinexus	PN941	RpAb
CNR2	S335+S336+	P34972				●												Kinexus	PN792	RpAb
CNTN1	Y742	Q12860																Kinexus	PN942	RpAb
Cofilin 1 (CFL1)	S3	P23528												●				Kinexus	PN812	RpAb
Cofilin 1 (CFL1)	Y140	P23529												●				Kinexus	PN813	RpAb
Cofilin 1 (CFL1)	Y140	P23529												●				Kinexus	PN506	RpAb
COT (MAP3K8; TPL2)	Pan	P41279	• PSTK															Kinexus	NK042-2	RpAb
COT (MAP3K8; TPL2)	Pan	P41279	• PSTK															Kinexus	NK042-1	RpAb
COT (MAP3K8; TPL2)	T290	P41279	• PSTK															Kinexus	PK956	RpAb
COT (MAP3K8; TPL2)	S334	P41279	• PSTK															Kinexus	PK585	RpAb
COX1 (PTGS1)	Pan	P23219								●								No	NN381-1	RpAb
COX2	Pan	P35354								●								No	610204	MmAb
COX2	Y446	P35354								●								Kinexus	PN695	RpAb
COX4/COX4I1	Pan	P13073								●								No	NN379-1	RpAb
CPD	Pan	O75976								●								Kinexus	NN462-3	RpAb
CPD	Pan	O75976								●								Kinexus	NN462-2	RpAb
Cpn10 (HSPE1)	Pan	P61604									●							No	NN235-1	RpAb
CPR (POR)	Pan	P16435																No	NN236-1	RpAb
CREB1	S133	P16220			●													Kinexus	PN553	RpAb
CREB1	S133	P16220			●													No	06-519	RpAb
CREB1	S129+S133	P16220			●													Kinexus	PN577	RpAb
CREBBP	S121+S124	Q92793			●													Kinexus	PN793	RpAb
CRMP2 (DPYSL2)	Y499	Q16555												●				Kinexus	PN944	RpAb
CRMP2 (DPYSL2)	T509	Q16555												●				Kinexus	PN814	RpAb
CRMP2 (DPYSL2)	T514	Q16555												●				Kinexus	PN943	RpAb
CRMP2 (DPYSL2)	S522	Q16555												●				Kinexus	PN815	RpAb
CSF1R (Fms)	Pan	P07333	• PYK			●												Kinexus	NK234-3	RpAb
CSF1R (Fms)	Pan	P07333	• PYK			●												Kinexus	NK234-4	RpAb
CSF1R (Fms)	Pan	P07333	• PYK			●												Kinexus	NK234-2	RpAb
CSF1R (Fms)	Y699	P07333	• PYK			●												Kinexus	PK587	RpAb
CSF1R (Fms)	S807+Y809	P07333	• PYK			●												Kinexus	PK586	RpAb
CSF1R (Fms)	Y809	P07333	• PYK			●												Kinexus	PK588	RpAb
Csk	Pan	P41240	• PYK															No	sc-9898	GpAb
Csk	Pan	P41240	• PYK															No	sc-286	RpAb
Csk	Y184	P41240	• PYK															Kinexus	PK589	RpAb
CTNNA1	S641	P35221												●				No	11330	RpAb
CTNNB1 (Catenin-b)	Pan	P35222			●									●				No	DB077	RpAb
CTNNB1 (Catenin-b)	Pan	P35222			●									●				No	KalCTNB5	RpAb
CTNNB1 (Catenin-b)	S33	P35222			●									●				No	11218	RpAb
CTNNB1 (Catenin-b)	S33+S37	P35222			●									●				Kinexus	PN816	RpAb
CTNNB1 (Catenin-b)	S33+S37+T4	P35222			●									●				Kinexus	PN578	RpAb
CTNNB1 (Catenin-b)	Y333	P35222			●									●				No	11574	RpAb
CTNNB1 (Catenin-b)	Y489	P35222			●									●				Kinexus	PN745	RpAb
CTNNB1 (Catenin-b)	S552	P35222			●									●				Kinexus	PN945	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
CTNNB1 (Catenin-b)	Y654	P35222			●									●				Kinexus	PN579	RpAb
CTNNB1 (Catenin-b)	S675	P35222			●									●				Kinexus	PN946	RpAb
CTSS	Pan	P25774							●									Kinexus	NN461-2	RpAb
CTSS	Pan	P25774							●									Kinexus	NN461-3	RpAb
CTSS	Pan	P25774							●									Kinexus	NN461-1	RpAb
CUX1 (CUTL1)	S1270	P39880			●													Kinexus	PN580	RpAb
Cyclin A1 (CCNA; CCN1)	Pan	P20248	• PSTK-															No	100-401-15	RpAb
Cyclin B1 (CCNB1; CLNB1)	Pan	P14635	• PSTK-															No	sc-245	MmAb
Cyclin C (CCNC)	Pan	P24863	• PSTK-															No	sc-1061	RpAb
Cyclin D1 (CCND1; CLND1)	Pan	P24385	• PSTK-															No	sc-753	RpAb
Cyclin D2 (CCND2; CLND2)	Pan	P30279	• PSTK-															No	sc-754	RpAb
Cyclin G1 (CCNG1; CLNG1)	Pan	P51959	• PSTK-															No	sc-7865	RpAb
DAG1	T790	Q14118												●				Kinexus	PN947	RpAb
DAPK2	Pan	Q9UIK4	• PSTK															No	AP7218a	RpAb
DARPP32 (PPP1R1B)	Y74	Q9UD71		• PSTP-														Kinexus	PN794	RpAb
DARPP32 (PPP1R1B)	T75	Q9UD71		• PSTP-														Kinexus	PN818	RpAb
DCAMLK1 (DCLK1)	Pan	O15075	• PSTK															No	AP7219a	RpAb
DCK (dCK)	S74+Pan	P27707							●									Kinexus	PN696	RpAb
DCTN1	S105+T108	Q14203																Kinexus	PN873	RpAb
DCTN1	S541	Q14203																Kinexus	PN874	RpAb
DCTN1	T1152+S115	Q14203																Kinexus	PN875	RpAb
DDR1	Pan	Q08345	• PYK			●												Kinexus	NK263-1	RpAb
DDR1	Y796+Y797	Q08345	• PYK			●												Kinexus	PK591	RpAb
DDR1	Y797	Q08345	• PYK			●												Kinexus	PK592	RpAb
DDR2 (Tyro10)	Y736	Q16832	• PYK			●												Kinexus	PK593	RpAb
DDR2 (Tyro10)	Y740	Q16832	• PYK			●												Kinexus	PK594	RpAb
DFF45	Pan	O00273																No	M037-3	MmAb
DGKd (DGK-D)	Pan	Q16760							●									No	AP8126a	RpAb
DGKd (DGK-D)	Pan	Q16760							●									No	AP8126b	RpAb
DGKz (DGK-Z)	Pan	Q13574							●									No	AP8060b	RpAb
DICER1 (Dicer; HERNA;	Pan	Q9UPY3																No	NN237-1	MmAb
DLG1 (SAP97)	Y760	Q12959											●					Kinexus	PN948	RpAb
DLG1 (SAP97)	Y784	Q12959											●					Kinexus	PN949	RpAb
DLG3 (SAP102; Synapse-	Pan	Q92796											●					No	NN238-1	MmAb
DLG4 (PSD95)	Pan	P78352											●					No	NN239-1	MmAb
DLG4 (PSD95)	Pan	P78352											●					No	NN142	MmAb
DLG4 (PSD95)	Pan	P78352											●					No	NN239-2	MmAb
DLG4 (PSD95)	S73	P78352											●					Kinexus	PN900	RpAb
DLG4 (PSD95)	Y240	P78352											●					Kinexus	PN901	RpAb
DLG4 (PSD95)	S295	P78352											●					Kinexus	PN846	RpAb
DLG4 (PSD95)	S561	P78352											●					Kinexus	PN899	RpAb
DLK (DAK; ZPK; MAP3K12;	Pan	Q12852	• PSTK															No	AP7217b	RpAb
DLK (DAK; ZPK; MAP3K12;	Pan	Q12852	• PSTK															No	AP7214a	RpAb
DLK (DAK; ZPK; MAP3K12;	Pan	Q12852	• PSTK															No	sc-8124	GpAb
DLK (DAK; ZPK; MAP3K12;	Pan	Q12852	• PSTK															No	sc-8125	GpAb
DLK (DAK; ZPK; MAP3K12;	Pan	Q12852	• PSTK															No	sc-10805	RpAb
DLK (DAK; ZPK; MAP3K12;	S269	Q12852	• PSTK															Kinexus	PK590	RpAb
DLK (DAK; ZPK; MAP3K12;	S308+Pan	Q12852	• PSTK															Kinexus	PK900	RpAb
DMPK1 (DMPK)	Pan	Q09013	• PSTK															No	AP7034b	RpAb
DMPK1 (DMPK)	Pan	Q09013	• PSTK															Kinexus	NK264-1	RpAb
DMPK2 (MRCK-g)	Pan	Q6DT37	• PSTK															Kinexus	NK265-1	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															Kinexus	NK048-4	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															Kinexus	NK048-4	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															Kinexus	NK048-6	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															Kinexus	NK048-7	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															Kinexus	NK048-5	RpAb
DNAPK (PRKDC)	Pan	P78527	• PSTK															No	sc-9051	RpAb
DNAPK (PRKDC)	Y883	P78527	• PSTK															Kinexus	PK596	RpAb
DNAPK (PRKDC)	S2056	P78527	• PSTK															Kinexus	PK925	RpAb
DNAPK (PRKDC)	T2609	P78527	• PSTK															Kinexus	PK595	RpAb
DNM2	Y597	P50570														●		Kinexus	PN951	RpAb
DNM2	S764	P50570														●		Kinexus	PN950	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
DNMT1 (AIM; CXXC9; DNMT;	Pan	P26358								●								No	NN241-2	MmAb
DNMT1 (AIM; CXXC9; DNMT;	Pan	P26358								●								No	NN241-3	MmAb
DNMT3A	S105	Q9Y6K1																Kinexus	PN746	RpAb
DOK1	Pan	Q99704											●					No	AP2001a	RpAb
DOK3	Y398	Q7L591											●					Kinexus	PN508	RpAb
DOK7	Y395	Q18PE1											●					Kinexus	PN952	RpAb
DOK7	Y405	Q18PE1											●					Kinexus	PN953	RpAb
DOR1	T358+T361+	P41143				●												Kinexus	PN795	RpAb
DPYD	S766+T768	Q12882								●								Kinexus	PN747	RpAb
DTNA	T504	Q9Y4J8																Kinexus	PN954	RpAb
DUSP1 (MKP1; CL100; VH1)	Pan	P28562		• DSP														Kinexus	NP006-3	RpAb
DUSP1 (MKP1; CL100; VH1)	Pan	P28562		• DSP														Kinexus	NP006-4	RpAb
DUSP1 (MKP1; CL100; VH1)	Pan	P28562		• DSP														No	sc-1199	RpAb
DUSP1 (MKP1)	Pan	P28562		• DSP														Kinexus	NP006-2	RpAb
DUSP2 (PAC1)	Pan	Q05923		• DSP														Kinexus	NP008-3	RpAb
DUSP2 (PAC1)	Pan	Q05923		• DSP														Kinexus	NP008-4	RpAb
DUSP2 (PAC1)	Pan	Q05923		• DSP														Kinexus	NP008-2	RpAb
DUSP2 (PAC1)	Pan	Q05923		• DSP														No	sc-1620	GpAb
DUSP2 (PAC1)	Pan	Q05923		• DSP														No	sc-1621	GpAb
DUSP3 (VHR)	Pan	P51452		• DSP														Kinexus	NP030-3	RpAb
DUSP3 (VHR)	Pan	P51452		• DSP														Kinexus	NP030-4	RpAb
DUSP3 (VHR)	Pan	P51452		• DSP														No	sc-8890	GpAb
DUSP3 (VHR)	Pan	P51452		• DSP														No	sc-8889	GpAb
DUSP3 (VHR)	Pan	P51452		• DSP														No	V47020	MmAb
DUSP4 (MKP2; VH2)	Pan	Q13115		• DSP														Kinexus	NP007-3	RpAb
DUSP4 (MKP2; VH2)	Pan	Q13115		• DSP														Kinexus	NP007-4	RpAb
DUSP4 (MKP2; VH2)	Pan	Q13115		• DSP														Kinexus	NP007-2	RpAb
DUSP5	Pan	Q16690		• DSP														Kinexus	NP039-2	RpAb
DUSP5	Pan	Q16690		• DSP														Kinexus	NP039-3	RpAb
DUSP5	Pan	Q16690		• DSP														Kinexus	NP039-1	RpAb
DUSP6	Pan	Q16828		• DSP														Kinexus	NP040-2	RpAb
DUSP6	Pan	Q16828		• DSP														Kinexus	NP040-1	RpAb
DUSP7	Pan	Q16829		• DSP														Kinexus	NP041-2	RpAb
DUSP7	Pan	Q16829		• DSP														Kinexus	NP041-3	RpAb
DUSP7	Pan	Q16829		• DSP														Kinexus	NP041-1	RpAb
DUSP8	Pan	Q13202		• DSP														Kinexus	NP042-3	RpAb
DUSP8	Pan	Q13202		• DSP														Kinexus	NP042-1	RpAb
DUSP9	Pan	Q99956		• DSP														Kinexus	NP043-3	RpAb
DUSP9	Pan	Q99956		• DSP														Kinexus	NP043-2	RpAb
DUSP10 (MKP5)	Pan	Q9Y6W6		• DSP														Kinexus	NP047-2	RpAb
DUSP10 (MKP5)	Pan	Q9Y6W6		• DSP														Kinexus	NP047-1	RpAb
DUSP11	Pan	O75319		• DSP														Kinexus	NP048-2	RpAb
DUSP11	Pan	O75319		• DSP														Kinexus	NP048-3	RpAb
DUSP11	Pan	O75319		• DSP														Kinexus	NP048-1	RpAb
DUSP12	Pan	Q9UNI6		• DSP														Kinexus	NP046-2	RpAb
DUSP12	Pan	Q9UNI6		• DSP														Kinexus	NP046-3	RpAb
DUSP12	Pan	Q9UNI6		• DSP														Kinexus	NP046-1	RpAb
DYRK1A	Y321	Q13627	• DSK															Kinexus	PK597	RpAb
DYRK2	Pan	Q92630	• DSK															No	AP7534a	RpAb
DYRK2	Pan	Q92630	• DSK															Kinexus	NK266-1	RpAb
DYRK2	Y382	Q92630	• DSK															Kinexus	PK598	RpAb
DYRK3	Pan	O43781	• DSK															No	AP7535a	RpAb
DYRK4	Pan	Q9NR20	• DSK															No	AP7536b	RpAb
EEF1A1	Y141	P68104						●										Kinexus	PN509	RpAb
EEF2	T57	P13639						●										Kinexus	PN555	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												No	AP7628b	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												No	AP7628a	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												Kinexus	NK052-5	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												Kinexus	NK052-6	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												Kinexus	NK052-4	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												Kinexus	NK052-7	RpAb
EGFR (ErbB1)	Pan	P00533	• PYK			●												Kinexus	NK052-8	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
EGFR (ErbB1)	Pan	P00533	• PYK			●												No	100-401-145	RpAb
EGFR (ErbB1)	Y869	P00533	• PYK			●												Kinexus	PK602	RpAb
EGFR (ErbB1)	Y998	P00533	• PYK			●												Kinexus	PK603	RpAb
EGFR (ErbB1)	Y1069	P00533	• PYK			●												Kinexus	PK599	RpAb
EGFR (ErbB1)	Y1110	P00533	• PYK			●												Kinexus	PK600	RpAb
EGFR (ErbB1)	Y1172	P00533	• PYK			●												Kinexus	PK601	RpAb
eIF2a	S52+Pan	Q9BY44						●										Kinexus	PN820	RpAb
EIF2AK3 (PERK; PEK; PKR)	S555	Q9NZJ5	• PSTK															Kinexus	PK926	RpAb
EIF2AK3 (PERK; PEK; PKR)	S715	Q9NZJ5	• PSTK															Kinexus	PK927	RpAb
EIF2AK3 (PERK; PEK; PKR)	T982	Q9NZJ5	• PSTK															Kinexus	PK604	RpAb
eIF4B	S422	P23588						●										No	11513	RpAb
eIF4B	S504	P23588						●										Kinexus	PN819	RpAb
eIF4E	Pan	P06730						●										No	610270	MmAb
eIF4E	S209	P06730						●										Kinexus	PN956	RpAb
eIF4G1	T205+T207	Q04637						●										Kinexus	PN1008	RpAb
eIF4G1	S1185+S118	Q04637						●										Kinexus	PN1004	RpAb
eIF4G1	S1231	Q04637						●										No	11514	RpAb
Elk1	S324+Pan	P19419			●													Kinexus	PN581	RpAb
EML4	Y226	Q9HC35												●				Kinexus	PN510	RpAb
ENO1 (NNE)	Y44	P06733								●								Kinexus	PN511	RpAb
ENO2 (ENOG)	Y25	P09104								●								Kinexus	PN512	RpAb
EP300 (p300)	S2279	Q09472																Kinexus	PN748	RpAb
EphA1	Pan	P21709	• PYK			●												No	AP7606b	RpAb
EphA1	Y599	P21709	• PYK			●												Kinexus	PK928	RpAb
EphA1	Y781	P21709	• PYK			●												Kinexus	PK605	RpAb
EphA2	Pan	P29317	• PYK			●												No	AP7607b	RpAb
EphA2	Pan	P29317	• PYK			●												No	AP7701a	RpAb
EphA2	Y588	P29317	• PYK			●												Kinexus	PK606	RpAb
EphA2	Y772	P29317	• PYK			●												Kinexus	PK607	RpAb
EphA3	Pan	P29320	• PYK			●												No	AP7608a	RpAb
EphA3	Pan	P29320	• PYK			●												No	AP7608b	RpAb
EphA3	Y779	P29320	• PYK			●												Kinexus	PK608	RpAb
EphA4	Pan	P54764	• PYK			●												No	AP7609a	RpAb
EphA4	Pan	P54764	• PYK			●												No	AP7609b	RpAb
EphA4	Y602	P54764	• PYK			●												Kinexus	PK930	RpAb
EphA4	Y779	P54764	• PYK			●												Kinexus	PK931	RpAb
EphA5	Pan	P54756	• PYK			●												No	AP7610a	RpAb
EphA5	Pan	P54756	• PYK			●												No	AP7610b	RpAb
EphA6	Pan	Q9UF33	• PYK			●												No	AP7611a	RpAb
EphA6	Pan	Q9UF33	• PYK			●												No	AP7611b	RpAb
EphA7	Pan	Q15375	• PYK			●												No	AP7612a	RpAb
EphA8	Pan	P29322	• PYK			●												No	AP7613a	RpAb
EphA8	Pan	P29322	• PYK			●												No	AP7613b	RpAb
EphB1	Pan	P54762	• PYK			●												No	AP7622a	RpAb
EphB1	Y594	P54762	• PYK			●												Kinexus	PK609	RpAb
EphB2	Pan	P29323	• PYK			●												No	AP7623a	RpAb
EphB2	Pan	P29323	• PYK			●												No	AP7623b	RpAb
EphB2	Pan	P29323	• PYK			●												Kinexus	NK267-1	RpAb
EphB2	Pan	P29323	• PYK			●												Kinexus	NK267-2	RpAb
EphB2	Y780	P29323	• PYK			●												Kinexus	PK610	RpAb
EphB3	Pan	P54753	• PYK			●												No	AP7624a	RpAb
EphB3	Y600	P54753	• PYK			●												Kinexus	PK611	RpAb
EphB4	Pan	P54760	• PYK			●												No	AP7625a	RpAb
EphB4	Y596	P54760	• PYK			●												Kinexus	PK612	RpAb
EphB6	Pan	Q15197	• PYK			●												No	AP7627a	RpAb
EphB6	Pan	Q15197	• PYK			●												No	AP7627b	RpAb
EPM2A (LAFPTase)	Pan	O95278		• DSP														No	NN245-1	MmAb
ERa (ESR1; ER-alpha)	S104	P03372			●	●												No	11070	RpAb
ERa (ESR1; ER-alpha)	S167	P03372			●	●												Kinexus	PN583	RpAb
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●												No	AP7629b	RpAb
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●												No	AP7629a	RpAb
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●												No	DB049	RpAb
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●												Kinexus	NK054-4	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type	
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●												Kinexus	NK054-6	RpAb	
ErbB2 (Neu; HER2)	Pan	P04626	• PYK			●													Kinexus	NK054-7	RpAb
ErbB2 (Neu; HER2)	Y735	P04626	• PYK			●													Kinexus	PK614	RpAb
ErbB2 (Neu; HER2)	Y877	P04626	• PYK			●													Kinexus	PK615	RpAb
ErbB2 (Neu; HER2)	Y1248	P04626	• PYK			●													Kinexus	PK613	RpAb
ERBB2IP (Erbin)	Y1104	Q96RT1											●						Kinexus	PN513	RpAb
ErbB3 (HER3)	Pan	P21860	• PYK			●													Kinexus	NK231-2	RpAb
ErbB3 (HER3)	Pan	P21860	• PYK			●													Kinexus	NK231-3	RpAb
ErbB3 (HER3)	Pan	P21860	• PYK			●													Kinexus	NK231-4	RpAb
ErbB3 (HER3)	Pan	P21860	• PYK			●													Kinexus	NK231-2	RpAb
ErbB3 (HER3)	Y1289	P21860	• PYK			●													Kinexus	PK616	RpAb
ErbB3 (HER3)	Y1307	P21860	• PYK			●													Kinexus	PK617	RpAb
ErbB3 (HER3)	Y1328	P21860	• PYK			●													Kinexus	PK618	RpAb
ErbB4 (HER4)	Pan	Q15303	• PYK			●													No	AP7631a	RpAb
ErbB4 (HER4)	Pan	Q15303	• PYK			●													Kinexus	NK235-1	RpAb
ErbB4 (HER4)	Pan	Q15303	• PYK			●													Kinexus	NK235-3	RpAb
ErbB4 (HER4)	Y733	Q15303	• PYK			●													Kinexus	PK619	RpAb
ErbB4 (HER4)	Y875	Q15303	• PYK			●													Kinexus	PK620	RpAb
ERF	T526	P50548			●														Kinexus	PN584	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																No	AP7500a	RpAb
ERK1 (MAPK3)	Pan	P28482	• PSTK																No	AP7501a	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-1	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-2	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-3	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-5	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-4	RpAb
ERK1 (MAPK3)	Pan	P27361	• PSTK																Kinexus	NK055-6	RpAb
ERK1 (MAPK3)	Pan	P28482	• PSTK																Kinexus	NK056-2	RpAb
ERK1 (MAPK3)	Pan	P28482	• PSTK																Kinexus	NK056-3	RpAb
ERK1 (MAPK3)	Pan	P28482	• PSTK																Kinexus	NK056-4	RpAb
ERK1 (MAPK3)	Y204	P27361	• PSTK																Kinexus	PK864	RpAb
ERK1 (MAPK3)	Y204	P27361	• PSTK																No	sc-7383	MmAb
ERK1 (MAPK3)	T207	P27361	• PSTK																Kinexus	PK950	RpAb
ERK1 (MAPK3)	T207	P27361	• PSTK																Kinexus	PK865	RpAb
ERK1 (MAPK3)	S265	P27361	• PSTK																Kinexus	PK878	RpAb
ERK1 (MAPK3)	S283	P27361	• PSTK																Kinexus	PK879	RpAb
ERK1 (MAPK3)	T202+Y204	P27361	• PSTK																Kinexus	PK621	RpAb
ERK1 (MAPK3)	Y204+T207	P27361	• PSTK																Kinexus	PK866	RpAb
ERK1 (MAPK3)	Y263+S266	P28482	• PSTK																Kinexus	PK880	RpAb
ERK2 (MAPK1)	T185+Y187	P28482	• PSTK																Kinexus	PK622	RpAb
ERK3 (MAPK6)	Pan	Q16659	• PSTK																No	AP7502a	RpAb
ERK3 (MAPK6)	S189	Q16659	• PSTK																Kinexus	PK623	RpAb
ERK4 (MAPK4)	Pan	P31152	• PSTK																No	AP7503a	RpAb
ERK4 (MAPK4)	S186	P31152	• PSTK																Kinexus	PK624	RpAb
ERK5 (MAPK7; BMK)	Pan	Q13164	• PSTK																Kinexus	NK206-1	RpAb
ERK5 (MAPK7; BMK)	Pan	Q13164	• PSTK																Kinexus	NK206-4	RpAb
ERK5 (MAPK7; BMK)	Pan	Q13164	• PSTK																Kinexus	NK206-5	RpAb
ERK5 (MAPK7; BMK)	Pan	Q13164	• PSTK																No	sc-1285	GpAb
ERK5 (MAPK7; BMK)	Pan	Q13164	• PSTK																No	sc-1284	GpAb
ERK5 (MAPK7; BMK)	T219+Y221	Q13164	• PSTK																Kinexus	PK625	RpAb
ERK5 (MAPK7; BMK)	Y221	Q13164	• PSTK																Kinexus	PK626	RpAb
ESRRA (ERR1)	S19+S22	P11474			●	●													Kinexus	PN585	RpAb
ESYT1	Y822	Q9BSJ8														●			Kinexus	PN514	RpAb
Ets-1	S282	P14921			●														Kinexus	PN586	RpAb
ETV6 (Tel)	S22	P41212			●														Kinexus	PN587	RpAb
EZH2	T487	Q15910																	Kinexus	PN697	RpAb
Ezrin (EZR; Cytovillin)	Y354	P15311												●					No	11063	RpAb
FADD	Pan	Q13158											●						No	sc-6035	GpAb
FAK (PTK2)	Pan	Q05397	• PYK																No	610087	MmAb
FAK (PTK2)	Pan	Q05397	• PYK																No	sc-558	RpAb
FAK (PTK2)	Pan	Q05397	• PYK																No	sc-932	RpAb
FAK (PTK2)	Y397	Q05397	• PYK																Kinexus	PK627	RpAb
FAK (PTK2)	Y576+Y577	Q05397	• PYK																Kinexus	PK628	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
FAK (PTK2)	Y577	Q05397	• PYK															Kinexus	PK629	RpAb
FAK (PTK2)	S722	Q05397	• PYK															No	sc-16662-R	RpAb
FAS (Apo1; CD95)	Pan	P25445				●												No	sc-7886	RpAb
FasL (FasLG)	Pan	P48023										●						No	F37720	MmAb
FASN (FAS)	Y45	P49327								●								Kinexus	PN749	RpAb
FASN (FAS)	S207	P49327								●								Kinexus	PN698	RpAb
FBPase (FBP1)	S88	P09467								●								Kinexus	PN750	RpAb
FBPase (FBP1)	Y265	P09467								●								Kinexus	PN699	RpAb
FBPase 2 (FBP2)	Y216	O00757								●								Kinexus	PN700	RpAb
FBPase 2 (FBP2)	Y259	O00757								●								Kinexus	PN751	RpAb
Fer (TYK3)	Pan	P16591	• PYK															No	AP7704b	RpAb
Fer (TYK3)	Y402	P16591	• PYK															Kinexus	PK630	RpAb
β	Y714	P16591	• PYK															Kinexus	PK631	RpAb
Fes	Pan	P07332	• PYK															No	sc-7670	GpAb
Fes	Y713	P07332	• PYK															Kinexus	PK632	RpAb
Fes	Y713+S716	P07332	• PYK															Kinexus	PK633	RpAb
FGF13 (FHF2)	Pan	Q92913																No	NN246-1	MmAb
FGFR1 (FLT2)	Pan	P11362	• PYK			●												No	AP7643a	RpAb
FGFR1 (FLT2)	Pan	P11362	• PYK			●												No	AP7636a	RpAb
FGFR1 (FLT2)	Pan	P11362	• PYK			●												Kinexus	NK062-2	RpAb
FGFR1 (FLT2)	Pan	P11362	• PYK			●												Kinexus	NK062-3	RpAb
FGFR1 (FLT2)	Y653+Y654	P11362	• PYK			●												Kinexus	PK634	RpAb
FGFR2 (BEK)	Pan	P21802	• PYK			●												No	AP7637b	RpAb
FGFR2 (BEK)	Pan	P21802	• PYK			●												Kinexus	NK063-3	RpAb
FGFR2 (BEK)	Pan	P21802	• PYK			●												Kinexus	NK063-4	RpAb
FGFR2 (BEK)	Pan	P21802	• PYK			●												Kinexus	NK063-2	RpAb
FGFR2 (BEK)	Y656+Y657	P21802	• PYK			●												Kinexus	PK635	RpAb
FGFR3	Pan	P22607	• PYK			●												No	AP7638a	RpAb
FGFR3	Pan	P22607	• PYK			●												Kinexus	NK236-1	RpAb
FGFR3	Pan	P22607	• PYK			●												Kinexus	NK236-2	RpAb
FGFR3	Pan	P22607	• PYK			●												Kinexus	NK236-3	RpAb
FGFR3	Y647+Y648	P22607	• PYK			●												Kinexus	PK637	RpAb
FGFR3	Y647+Y648	P22607	• PYK			●												Kinexus	PK636	RpAb
FGFR4	Pan	P22455	• PYK			●												No	AP7639a	RpAb
FGFR4	Pan	P22455	• PYK			●												Kinexus	NK239-1	RpAb
FGFR4	Pan	P22455	• PYK			●												Kinexus	NK239-3	RpAb
Fgr	Pan	P09769	• PYK															Kinexus	NK268-1	RpAb
Fgr	Y208+Y209	P09769	• PYK															Kinexus	PK638	RpAb
Fgr	Y412	P09769	• PYK															Kinexus	PK639	RpAb
FIH (HIF1; HIF1AN)	Pan	Q9NWT6									●							No	NN248-1	MmAb
FKBP52 (FKBP4; FKBP51;	Pan	Q02790																No	NN127-2	MmAb
FLT3 (STK1; CD135)	Pan	P36888	• PYK			●												No	AP7644b	RpAb
FLT3 (STK1; CD135)	Pan	P36888	• PYK			●												No	AP7644a	RpAb
FLT3 (STK1; CD135)	Pan	P36888	• PYK			●												Kinexus	NK240-1	RpAb
FLT3 (STK1; CD135)	Pan	P36888	• PYK			●												Kinexus	NK240-2	RpAb
FLT3 (STK1; CD135)	Y842	P36888	• PYK			●												Kinexus	PK640	RpAb
Fos	Pan	P01100			●													No	DB025	RpAb
Fos	T232	P01100			●													No	44-280G	RpAb
Fos	T232	P01100			●													Kinexus	PN588	RpAb
Fos	S362+S363	P01100			●													Kinexus	PN589	RpAb
FOXK1	S441+S445	P85037			●													Kinexus	PN590	RpAb
FOXK2	S424+S428	Q01167			●													Kinexus	PN591	RpAb
FOXM1	T611	Q08050			●													Kinexus	PN592	RpAb
FOXO1A (FKHR; FKHL1)	S256	Q12778			●													Kinexus	PN593	RpAb
FOXO1A (FKHR; FKHL1)	S256	Q12778			●													No	11115	RpAb
FOXO1A (FKHR; FKHL1)	S319	Q12778			●													No	11136	RpAb
FOXO1A (FKHR; FKHL1)	S329	Q12778			●													Kinexus	PN595	RpAb
FOXO1A (FKHR; FKHL1)	S319+Pan	Q12778			●													Kinexus	PN594	RpAb
FOXO3 (FKHL1)	T32	O43524			●													Kinexus	PN596	RpAb
FOXO3 (FKHL1)	S253	O43524			●													Kinexus	PN821	RpAb
FOXO3 (FKHL1)	S294	O43524			●													Kinexus	PN597	RpAb
FRA1	S265	P15407			●													Kinexus	PN599	RpAb
Frk	Pan	P42685	• PYK															No	AP7708a	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Frk	Pan	P42685	• PYK															Kinexus	NK269-1	RpAb
Frk	Pan	P42685	• PYK															Kinexus	NK269-2	RpAb
Frk	Y387	P42685	• PYK															Kinexus	PK641	RpAb
Frk	Y497	P42685	• PYK															Kinexus	PK642	RpAb
FUS	S277	P35637																Kinexus	PN876	RpAb
FUS	S340	P35637																Kinexus	PN877	RpAb
FUS	Y468	P35637																Kinexus	PN878	RpAb
Fused (STK36)	S159	Q9NRP7	• PSTK															Kinexus	PK643	RpAb
FXR (NR1H4)	Pan	Q96R11				●												No	sc-13063	RpAb
Fyn	Pan	P06241	• PYK															No	AP7709a	RpAb
Fyn	Pan	P06241	• PYK															Kinexus	NK065-2	RpAb
Fyn	Pan	P06241	• PYK															No	sc-161	RpAb
Fyn	T12	P06241	• PYK															Kinexus	PK901	RpAb
Fyn	Y213+Y214	P06241	• PYK															Kinexus	PK644	RpAb
Fyn	Y420	P06241	• PYK															Kinexus	PK881	RpAb
Fyn	Y531	P06241	• PYK															Kinexus	PK645	RpAb
G6PD	Y401	P11413								●								Kinexus	PN515	RpAb
G6PD	Y503+Y507	P11413								●								Kinexus	PN701	RpAb
Gab1	Y406	Q13480											●					Kinexus	PN516	RpAb
GABA A Receptor (GABRA2)	Pan	Q9UBS5				●												No	NN250-1	MmAb
GABA A Receptor (GABRB3)	Pan	Q9UBS5				●												No	NN254-1	MmAb
GABA B Receptor 2	Pan	O75899				●												No	NN256-2	MmAb
GABARAP (FLC3B)	Pan	O95166				●												No	NN257-1	RpAb
GABBR1 (GABA B Receptor	Pan	Q9UBS5				●												No	NN259-1	MmAb
GABBR1 (GABA B Receptor	T873	Q9UBS5				●												Kinexus	PN796	RpAb
GABPA	S303	Q06546			●													Kinexus	PN957	RpAb
GABPA	Y380+Y381	Q06546			●													Kinexus	PN958	RpAb
GABPB1	Y364	Q06547			●													Kinexus	PN959	RpAb
GADD153 (DDIT3; CHOP)	Pan	P35638									●							No	sc-7351	MmAb
GALNT7	Pan	Q86SF2								●								Kinexus	NN463-1	RpAb
GATA1	S142	P15976			●													No	11041	RpAb
GATA3	S369	P23771			●													Kinexus	PN702	RpAb
GATAD2B	T120+S122	Q8WXI9			●													Kinexus	PN600	RpAb
GCK (Glucokinase)	S411+Pan	P35557								●								Kinexus	PK893	RpAb
GCK (MAP4K2)	Pan	Q12851	• PSTK															No	AP7901a	RpAb
GCK (MAP4K2)	Pan	Q12851	• PSTK															No	sc-6161	GpAb
GCK (MAP4K2)	S170	Q12851	• PSTK															Kinexus	PK646	RpAb
GCLC	S5+S8	P48506																Kinexus	PN752	RpAb
GCN2 (EIF2AK4)	Pan	Q9P2K8	• PSTK															No	AP8062a	RpAb
GCN2 (EIF2AK4)	T667	Q9P2K8	• PSTK															Kinexus	PK877	RpAb
GFAP	Pan	P14136												●				Kinexus	NN260-3	RpAb
GFAP	Pan	P14136												●				No	NN260-2	MmAb
GFAP	Pan	P14136												●				No	NN260-1	MmAb
GFAP	S133	P14136												●				Kinexus	PN932	RpAb
GIT1	Y545	Q9Y2X7																Kinexus	PN517	RpAb
GluR1	S849	P42261				●												Kinexus	PN822	RpAb
GluR2	Y876	P42262				●												Kinexus	PN823	RpAb
GLUT2 (SLC2A2)	Pan	P11168														●		No	NN262-1	RpAb
GPAT3 (LPAAT-THETA)	S68	Q53EU6								●								Kinexus	PN753	RoAb
GR (NR3C1)	S226+Pan	P04150			●	●												Kinexus	PN601	RpAb
GRK1 (Rhodopsin kinase)	Pan	Q15835	• PSTK															No	AP7003a	RpAb
GRK1 (Rhodopsin kinase)	Pan	Q15835	• PSTK															No	sc-561	MmAb
GRK1 (Rhodopsin kinase)	Pan	Q15835	• PSTK															No	sc-8004	MmAb
GRK2 (BARK1; ADRBK1)	Y356	P25098	• PSTK															Kinexus	PK537	RpAb
GRK2 (BARK1; ADRBK1)	S670	P25098	• PSTK															Kinexus	PK536	RpAb
GRK5 (GPRK5)	Pan	P34947	• PSTK															No	AP7007a	RpAb
Grp75 (HspA9)	Pan	P38646									●							No	NN047-2	MmAb
Grp78 (Hspa5; BiP)	Pan	P11021									●							No	NN048-4	MmAb
Grp78 (Hspa5; BiP)	Pan	P11021									●							No	NN048-5	MmAb
Grp94 (HSP90B1)	Pan	P14625									●							No	NN049-3	RatmAb
Grp94 (HSP90B1)	Pan	P14625									●							No	NN049-4	RpAb
Grp170 (HYOU1; ORP-150)	Pan	Q9Y4L1									●							Kinexus	NN265-3	RpAb
Grp170 (HYOU1; ORP-150)	Pan	Q9Y4L1									●							Kinexus	NN265-4	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Grp170 (HYOU1; ORP-150)	Pan	Q9Y4L1									●							Kinexus	NN265-5	RpAb
Grp170 (HYOU1; ORP-150)	Pan	Q9Y4L1									●							No	NN265-1	MmAb
GSK3a	Pan	P49840	• PSTK															Kinexus	NK069-3	RpAb
GSK3a	T19+S21	P49840	• PSTK															Kinexus	PK648	RpAb
GSK3a	S278+Y279	P49840	• PSTK															Kinexus	PK647	RpAb
GSK3a	Y279	P49840	• PSTK															Kinexus	PK649	RpAb
GSK3a	Y279	P49840	• PSTK															Kinexus	PK882	RpAb
GSK3a	Y284+Y285	P49840	• PSTK															Kinexus	PK650	RpAb
GSK3b	Pan	P49841	• PSTK															Kinexus	NK270-3	RpAb
GSK3b	Pan	P49841	• PSTK															Kinexus	NK070	RpAb
GSK3b	Pan	P49841	• PSTK															No	sc-7291	MmAb
GSK3b	Pan	P49841	• PSTK															No	NK070-4	RpAb
GSK3b	T43	P49841	• PSTK															Kinexus	PK933	RpAb
GSK3b	T275+T277	P49841	• PSTK															Kinexus	PK883	RpAb
GSK3b	T390	P49841	• PSTK															Kinexus	PK932	RpAb
GTF2F1	S385+T389	P35269			●													Kinexus	PK651	RpAb
GTF2I	S412	P78347			●													Kinexus	PN602	RpAb
GUK1	Y53	Q16774								●								Kinexus	PK652	RpAb
GYS1	S641+S645	P13807								●								Kinexus	PN703	RpAb
GYS2	Y45	P54840								●								Kinexus	PN704	RpAb
H2AFX (H2AX; Histone	S140	P16104																Kinexus	PN824	RpAb
H3F3A (HIST1H3A; Histone	T4	P84243																No	05-746	RpAb
H3F3A (HIST1H3A; Histone	S10	P84243																Kinexus	PN826	RpAb
H3F3A (HIST1H3A; Histone	S10	P84243																No	06-570	RpAb
Haspin (GSG2)	Pan	Q8TF76	• PSTK															No	AP8064c	RpAb
HBS1L	Y56+Y58	Q9Y450			●													Kinexus	PN603	RpAb
HCA59	Y147	Q9NZ63																Kinexus	PN518	RpAb
HCFC1	S1507	P51610			●													Kinexus	PN604	RpAb
Hck	Pan	P08631	• PYK															Kinexus	NK271-1	RpAb
HDAC4	S632	P56524																Kinexus	PN825	RpAb
HDAC4/5/9	S246/259/220	P56524																No	11517	RpAb
HePTP (PTPN7)	S44	P35236	• PYP															Kinexus	PP500	RpAb
HePTP (PTPN7)	T66	P35236	• PYP															Kinexus	PP501	RpAb
HePTP (PTPN7)	S143	P35236	• PYP															Kinexus	PP528	RpAb
HGK (MAP4K4; NIK; ZC1)	Pan	O95819	• PSTK															Kinexus	NK300-1	RpAb
HGK (MAP4K4; NIK; ZC1)	T187	O95819	• PSTK															Kinexus	PK653	RpAb
HGS (Hrs)	Y216	O14964	• PYK															Kinexus	PN519	RpAb
HIF1A (HIF1 alpha)	Pan	Q16665									●							No	sc-13515	MmAb
HIF1A (HIF1 alpha)	Pan	Q16665									●							No	NN267-1	MmAb
HIPK1	Y352	Q86Z02	• PSTK															Kinexus	PK654	RpAb
HIPK2	Pan	Q9H2X6	• PSTK															Kinexus	NK272-1	RpAb
Histone H2B (HIST1H2BN)	S14	Q99877																No	07-191	RpAb
HMGA1	S36+T39	P17096			●													Kinexus	PN605	RpAb
HMGA1	T53	P17096			●													Kinexus	PN606	RpAb
HMGB1	S35+S39	P09429			●													Kinexus	PN607	RpAb
HMGCR	S872	P04035								●								Kinexus	PN705	RpAb
HMGCS1	S495	Q01581								●								Kinexus	PN754	RpAb
HNF4A	S167	P41235			●													Kinexus	PN608	RpAb
HNRNPA1	S4+S6	P09651																Kinexus	PN879	RpAb
HNRNPA1	Y347	P09651																Kinexus	PN880	RpAb
HO1 (HO; HMOX1)	Pan	P09601									●							No	NN052-6	RpAb
HO2 (HMOX2)	Pan	P30519								●								No	OSA-200	RpAb
Hpk1	Pan	Q92918	• PSTK															No	sc-6230	GpAb
HRAS (H-Ras)	Pan	P01112																Kinexus	NN281-3	RpAb
HRAS (H-Ras)	Y157	P01112																Kinexus	PN755	RpAb
HRI (HCR; EIF2AK1)	S258	Q9BQI3	• PSTK															Kinexus	PK876	RpAb
Hsc70 (HSPA8; Hsc70;	Pan	P11142									●							No	sc-7298	MmAb
Hsc70 (HSPA8; Hsc70;	Pan	P11142									●							No	NN060-12	MmAb
Hsc70 (HSPA8; Hsc70;	Pan	P11142									●							No	NN054-4	MmAb
HSF1	Pan	Q00613			●													No	NN268-2	RatmAb
HSF1	Pan	Q00613			●													No	NN268-1	RatmAb
HSF1	S303+S307	Q00613			●													Kinexus	PN609	RpAb
HSF4 (HSTF4)	Pan	Q9ULV5			●						●							No	H65520	MmAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
HSP27 (HSP28; HSPB1)	Pan	P04792									●							No	NN152-2	RpAb
HSP27 (HSP28; HSPB1)	S15	P04792									●							No	SPA-525	RpAb
HSP40 (DNAJB1; DNAJ1;	Pan	P25685									●							No	NN057-2	MmAb
HSP60 (HSPD1; CPN60;	Pan	P10809									●							No	sc-13115	MmAb
HSP60 (HSPD1; CPN60;	Pan	P10809									●							No	SPA-875	RpAb
HSP60 (HSPD1; CPN60;	Pan	P10809									●							No	NN059-4	MmAb
HSP60 (HSPD1; CPN60;	Pan	P10809									●							No	NN059-3	MmAb
HSP70 (Hsc70; Hsp71;	Pan	P0DMV8									●							No	SPA-810	MmAb
HSP70 (Hsc70; Hsp71;	Pan	P0DMV8									●							No	NN060-11	MmAb
HSP70 (Hsc70; Hsp71;	Pan	P0DMV8									●							No	NN060-7	MmAb
HSP70 (Hsc70; Hsp71;	Pan	P0DMV8									●							No	NN060-9	MmAb
HSP90a (HSP90AA1; HSP90;	Pan	P07900									●							No	sc-13119	MmAb
HSP90a (HSP90AA1; HSP90;	Pan	P07900									●							No	NN061-4	MmAb
HSP90a (HSP90AA1; HSP90;	Pan	P07900									●							No	NN061-2	MmAb
HSP86; HSPC; HSPCA;											●							No	NN061-2	MmAb
HSP90AB1 (HSP90; HSP84;	Pan	P08238									●							No	NN061-3	MmAb
HSP90AB1 (HSP90; HSP84;	Pan	P08238									●							No	NN061-16	RpAb
HSP90AB1 (HSP90; HSP84;	Y484	P08238									●							Kinexus	PN520	RpAb
HSP105 (HSPH1; HSP110)	Pan	Q92598									●							No	sc-6241	RpAb
HSP105 (HSPH1; HSP110)	Pan	Q92598									●							No	NN269-1	RpAb
HspBP1	Pan	Q9NZL4									●							No	H98620	MmAb
HSPC121	S114	Q9P035								●								Kinexus	PN756	RpAb
Huntingtin (HTT)	S13+S16+Pa	P42858												●				Kinexus	PN827	RpAb
Huntingtin (HTT)	S417+S419+	P42858												●				Kinexus	PN828	RpAb
Huntingtin (HTT)	S421	P42858												●				Kinexus	PN829	RpAb
IAP1 (BIRC2; API2)	Pan	Q13490										●						No	sc-7943	RpAb
ICK	Pan	Q9UPZ9	• PSTK															Kinexus	NK073-5	RpAb
ICK	Pan	Q9UPZ9	• PSTK															Kinexus	NK073-3	RpAb
ICK	Pan	Q9UPZ9	• PSTK															Kinexus	NK073-4	RpAb
ICK	Y159	Q9UPZ9	• PSTK															Kinexus	PK656	RpAb
ICK	Y156+T157	Q9UPZ9	• PSTK															Kinexus	PK656	RpAb
IDH1	T75+T77	O75874								●								Kinexus	PN706	RpAb
IDH1	Y391	O75874								●								Kinexus	PN757	RpAb
IGF1R	Pan	P08069	• PYK			●												No	AP7649a	RpAb
IGF1R	Pan	P08069	• PYK			●												Kinexus	NK074-2	RpAb
IGF1R	Pan	P08069	• PYK			●												Kinexus	NK074-3	RpAb
IGF1R	Y1161+T116	P08069	• PYK			●												Kinexus	PK657	RpAb
IGF1R	Y1165+Y116	P08069	• PYK			●												No	11088	RpAb
IGF1R	Y1280	P08069	• PYK			●												No	11302	RpAb
IGF1R	Y1346	P08069	• PYK			●												Kinexus	PK658	RpAb
IkBα (MAD3; IκBa)	Pan	P25963	• PSTK-															No	DB075	RpAb
IkBα (MAD3; IκBa)	Pan	P25963	• PSTK-															No	sc-371	RpAb
IkBα (MAD3; IκBa)	Pan	P25963	• PSTK-															No	sc-8404	MmAb
IκBβ (NFKBIB)	Pan	Q15653																No	DB076	RpAb
IκBε (NFKBIE)	S161	H3BNC2																No	11213	RpAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	AHO0352	MmAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	sc-7606	MmAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	sc-7190	GpAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	sc-7218	RpAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	KAP-TF115	RpAb
IKKα (IκBKA; CHUK)	Pan	O15111	• PSTK															No	71-2300	RpAb
IKKα (IκBKA; CHUK)	T179+S180	O15111	• PSTK															Kinexus	PK659	RpAb
IKKβ (IκBKB; IKKB)	Pan	O14920	• PSTK															Kinexus	NK076-6	RpAb
IKKβ (IκBKB; IKKB)	Pan	O14920	• PSTK															Kinexus	NK076-5	RpAb
IKKβ (IκBKB; IKKB)	Pan	O14920	• PSTK															No	KAP-TF118	RpAb
IKKe (IκBKE)	S172	Q14164	• PSTK															Kinexus	PK660	RpAb
IKKg (NEMO)	S377	Q9Y6K9	• PSTK															Kinexus	PN758	RpAb
IKZF1	S361+S364	Q13422			●													Kinexus	PN797	RpAb
IKZF1	Y413	Q13422			●													Kinexus	PN707	RpAb
ILK1 (ILK)	Pan	Q13418	• PSTK															No	AP7651b	RpAb
ILK1 (ILK)	Pan	Q13418	• PSTK															No	sc-7516	GpAb
ILK1 (ILK)	Pan	Q13418	• PSTK															No	KAP-ST203	RpAb
ILK1 (ILK)	S343	Q13418	• PSTK															Kinexus	PK661	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
ILK1 (ILK)	Y351	Q13418	• PSTK															Kinexus	PK662	RpAb
InsR (IR)	Pan	P06213	• PYK			●												Kinexus	NK079-2	RpAb
InsR (IR)	Y1185	P06213	• PYK			●												Kinexus	PK951	RpAb
InsR (IR)	Y1189	P06213	• PYK			●												Kinexus	PK663	RpAb
InsR (IR)	Y1361	P06213	• PYK			●												Kinexus	PK952	RpAb
Integrin b1 (ITGB1; CD29)	Pan	P05556				●												No	sc-9970	MmAb
IRAK1	Pan	P51617	• PSTK															No	sc-7883	RpAb
IRAK1	T387	P51617	• PSTK															Kinexus	PK664	RpAb
IRAK2	Pan	O43187	• PSTK															No	KAP-ST20E	RpAb
IRAK3 (IRAK-M)	Pan	Q9Y616	• PSTK															No	AP7804b	RpAb
IRAK4	T345+S346	Q9NWZ3	• PSTK															Kinexus	PK665	RpAb
IRF3	T135	Q14653			●													Kinexus	PN610	RpAb
IRR (INSRR)	Pan	P14616	• PYK			●												Kinexus	NK273-1	RpAb
IRS1	Y612	P35568											●					Kinexus	PN960	RpAb
IRS1	S636+S639	P35568											●					Kinexus	PN831	RpAb
ITGA4 (Integrin a4)	S1021	P13612												●				No	44-864	RpAb
ITK	Pan	Q08881	• PYK															Kinexus	NK274-1	RpAb
ITK	Pan	Q08881	• PYK															Kinexus	NK274-3	RpAb
ITK	Y512	Q08881	• PYK															Kinexus	PK666	RpAb
ITSN2	Y968	Q9NZM3											●					Kinexus	PN521	RpAb
IWS1	S720+T721+	Q96ST2			●													Kinexus	PN611	RpAb
JAK1	Pan	P23458	• PYK															Kinexus	NK084-3	RpAb
JAK1	Pan	P23458	• PYK															Kinexus	NK084-5	RpAb
JAK1	Pan	P23458	• PYK															No	sc-1677	MmAb
JAK1	Pan	P23458	• PYK															No	sc-295	RpAb
JAK1	Y1034+Y103	P23458	• PYK															Kinexus	PK884	RpAb
JAK1	T1107	P23458	• PYK															Kinexus	PK895	RpAb
JAK2	Pan	O60674	• PYK															Kinexus	NK085-3	RpAb
JAK2	Pan	O60674	• PYK															Kinexus	NK085-4	RpAb
JAK2	Pan	O60674	• PYK															Kinexus	NK085-2	RpAb
JAK2	Y570	O60674	• PYK															Kinexus	PK668	RpAb
JAK2	Y1007+Y100	O60674	• PYK															No	ab68268	RpAb
JAK2	Y1007+Y100	O60674	• PYK															Kinexus	PK667	RpAb
JAK2	Pan	O60674	• PYK															No	sc-7229	RpAb
JAK2	Pan	O60674	• PYK															No	sc-294	RpAb
JAK3	Pan	P52333	• PYK															Kinexus	NK086-3	RpAb
JAK3	Pan	P52333	• PYK															Kinexus	NK086-4	RpAb
JAK3	Pan	P52333	• PYK															Kinexus	NK086-2	RpAb
JAK3	Pan	P52333	• PYK															No	sc-1078	GpAb
JAK3	Pan	P52333	• PYK															No	sc-6932	MmAb
JAK3	Y785	P52333	• PYK															Kinexus	PK934	RpAb
JAK3	Y980+Y981	P52333	• PYK															Kinexus	PK669	RpAb
JIP2 (Mapk8ip2)	Pan	Q13387											●					No	NN272-1	MmAb
JNK1 (MAPK8; SAPK1)	Pan	P45983	• PSTK															Kinexus	NK217-2	RpAb
JNK1 (MAPK8; SAPK1)	Pan	P45983	• PSTK															Kinexus	NK217-3	RpAb
JNK1 (MAPK8; SAPK1)	Pan	P45983	• PSTK															No	sc-571	RpAb
JNK1 (MAPK8; SAPK1)	Pan	P45983	• PSTK															No	sc-474-G	GpAb
JNK1 (MAPK8; SAPK1)	Y185	P45983	• PSTK															Kinexus	PK670	RpAb
JNK2 (MAPK9)	Pan	P45984	• PSTK															Kinexus	NK189-2	RpAb
JNK2 (MAPK9)	Pan	P45984	• PSTK															Kinexus	NK189-3	RpAb
JNK2 (MAPK9)	Pan	P45984	• PSTK															Kinexus	NK189-4	RpAb
JNK2 (MAPK9)	Pan	P45984	• PSTK															Kinexus	NK196	RpAb
JNK2 (MAPK9)	Pan	P45984	• PSTK															No	sc-7345	MmAb
JNK2 (MAPK9)	T183+Y185	P45984	• PSTK															Kinexus	PK902	RpAb
JNK3 (MAPK10)	Pan	P53779	• PSTK															Kinexus	NK197-4	RpAb
JNK3 (MAPK10)	Pan	P53779	• PSTK															Kinexus	NK197	RpAb
JNK3 (MAPK10)	Pan	P53779	• PSTK															Kinexus	NK197-2	RpAb
Jun (c-Jun)	Pan	P05412			●													No	J31920	MmAb
Jun (c-Jun)	S63	P05412			●													No	9261L	RpAb
Jun (c-Jun)	S63	P05412			●													Kinexus	PN557	RpAb
Jun (c-Jun)	S63	P05412			●													No	sc-822	MmAb
Jun (c-Jun)	S73	P05412			●													Kinexus	PN612	RpAb
Jun (c-Jun)	T239	P05412			●													Kinexus	PN613	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Jun (c-Jun)	S243	P05412			●													Kinexus	PN614	RpAb
KAP (KAP1; CDKN3; Cdi1)	Pan	Q16667		• DSP														No	sc-475	RpAb
KCC2 (SLC12A5)	S963	Q9H2X9														●		Kinexus	PN832	RpAb
KCNQ1 (KCNA8; KCNA9;	Pan	P51787					●											No	NN274-1	MmAb
KCNQ2	Pan	O43526					●											No	NN275-1	MmAb
KDEL Receptor (Kdelr1;	Pan	P24390				●												No	NN153	MmAb
KHS1 (MAP4K5; KHS)	Pan	Q9Y4K4	• PSTK															Kinexus	NK089-2	RpAb
KHS1 (MAP4K5; KHS)	Pan	Q9Y4K4	• PSTK															No	sc-6429	GpAb
KHS1 (MAP4K5; KHS)	Y31	Q9Y4K4	• PSTK															Kinexus	PK672	RpAb
KHS1 (MAP4K5; KHS)	S174	Q9Y4K4	• PSTK															Kinexus	PK671	RpAb
KIF5A	S155	Q12840																Kinexus	PN881	RpAb
KIF5A	S176	Q12840																Kinexus	PN882	RpAb
Kir2.2 (KCNJ12; IRK2;	Pan	Q14500					●											No	NN278-1	MmAb
Kir6.1 (KCNJ8; uKATP-1)	Pan	Q15842					●											No	NN280-1	MmAb
Kit	Pan	P10721	• PYK			●												Kinexus	NK241-1	RpAb
Kit	Pan	P10721	• PYK			●												Kinexus	NK241-2	RpAb
Kit	Pan	P10721	• PYK			●												Kinexus	NK241-3	RpAb
Kit	Y721	P10721	• PYK			●												Kinexus	PK885	RpAb
Kit	S821+Y823	P10721	• PYK			●												Kinexus	PK674	RpAb
Kit	Y936	P10721	• PYK			●												Kinexus	PK673	RpAb
KMT2C	S4267	Q8NEZ4																Kinexus	PN798	RpAb
KOR-1 (MSL-1)	Y369	P41145				●												Kinexus	PN799	RpAb
KRAS (K-Ras; KRAS2;	Pan	P01116																Kinexus	NN281-2	RpAb
KRAS (K-Ras; KRAS2;	S181	P01116																Kinexus	PN862	RpAb
Ksr1	Pan	Q8IVT5	• PSTK															Kinexus	NK090-1	10320
Ksr1	Pan	Q8IVT5	• PSTK															No	sc-6212	GpAb
Ksr1	S406	Q8IVT5	• PSTK															Kinexus	PK675	RpAb
Ksr2	S490	Q6VAB6	• PSTK															Kinexus	PK676	RpAb
Kv3.1 (KCNC1; Kv4)	Pan	P48547					●											No	NN282-1	RatmAb
Kv4.2	T602+T607	Q9NZV8					●											Kinexus	PN833	RpAb
L1CAM (CD171)	T1172	P32004													●			Kinexus	PN961	RpAb
L1CAM (CD171)	Y1176	P32004													●			Kinexus	PN962	RpAb
LASS2	T346+S348	Q96G23																Kinexus	PN1005	RpAb
LATS1	Pan	O95835	• PSTK															No	AP7035a	RpAb
LATS1	Pan	O95835	• PSTK															Kinexus	NK091-2	RpAb
LATS1	S464	O95835	• PSTK															Kinexus	PK677	RpAb
LATS1	S909	O95835	• PSTK															Kinexus	PK678	RpAb
LATS2 (KPM)	Pan	Q9NRM7	• PSTK															Kinexus	NK092-2	RpAb
LATS2 (KPM)	Pan	Q9NRM7	• PSTK															Kinexus	NK092-1	RpAb
Lck	Pan	P06239	• PYK															No	sc-433	MmAb
Lck	Y192	P06239	• PYK															Kinexus	PK679	RpAb
Lck	Y263+Y264	P06239	• PYK															Kinexus	PK680	RpAb
Lck	Y394	P06239	• PYK															Kinexus	PK903	RpAb
LEDGF	S273+S275	O75475			●													Kinexus	PN615	RpAb
LEF1	T155+Pan	Q9UJU2			●													Kinexus	PN616	RpAb
LG1 (Epitempin-1)	Pan	O95970																No	NN285-1	MmAb
LIMK1	T508	P53667	• PSTK															Kinexus	PK681	RpAb
LKB1 (STK11; STRAD)	Pan	Q15831	• PSTK															Kinexus	NK227-3	RpAb
LKB1 (STK11; STRAD)	Pan	Q15831	• PSTK															Kinexus	NK227-4	RpAb
LKB1 (STK11; STRAD)	Pan	Q15831	• PSTK															Kinexus	NK227-2	RpAb
LKB1 (STK11; STRAD)	S31	Q15831	• PSTK															Kinexus	PK682	RpAb
LKB1 (STK11; STRAD)	S428	Q15831	• PSTK															Kinexus	PK683	RpAb
LMNA	T19+S22	P02545												●				Kinexus	PN966	RpAb
LMNA	S301	P02545												●				Kinexus	PN963	RpAb
LMNA	S390+S392	P02545												●				Kinexus	PN964	RpAb
LMNA	S458	P02545												●				Kinexus	PN965	RpAb
LMR2 (LMTK2; KPI-2)	S1450	Q8IWU2	• PSTK															Kinexus	PK684	RpAb
LOK	S191	O94804	• PSTK															Kinexus	PK685-2	RpAb
LOK	S191	O94804	• PSTK															Kinexus	PK685	RpAb
LOK	T952	O94804	• PSTK															Kinexus	PK686	RpAb
LRP4 (LRP10; MEGF7)	Pan	O75096				●												No	NN287-2	MmAb
LRP4 (LRP10; MEGF7)	Pan	O75096				●												No	NN286-1	MmAb
LRP4 (LRP10; MEGF7)	S1887	O75096				●												Kinexus	PN967	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
LRRK2 (PARK8)	Pan	Q5S007	• PSTK															No	NK303-1	MmAb
LRRK2 (PARK8)	S910	Q5S007	• PSTK															Kinexus	PK904	RpAb
LRRK2 (PARK8)	S935	Q5S007	• PSTK															Kinexus	PK905	RpAb
LRRK2 (PARK8)	S935	Q5S007	• PSTK															Kinexus	PK935	RpAb
LRRK2 (PARK8)	S973	Q5S007	• PSTK															Kinexus	PK936	RpAb
LRRK2 (PARK8)	T1410	Q5S007	• PSTK															Kinexus	PK937	RpAb
LRRK2 (PARK8)	T1503	Q5S007	• PSTK															Kinexus	PK938	RpAb
LRRK2 (PARK8)	T2031+S203	Q5S007	• PSTK															Kinexus	PK939	RpAb
LTB4R	T308+S310	Q15722				●												Kinexus	PN800	RpAb
LTK	Y672	P29376				●												Kinexus	PK687	RpAb
Lyn	Pan	P07948	• PYK															Kinexus	NK095-2	RpAb
Lyn	Pan	P07948	• PYK															No	sc-7274	MmAb
Lyn	Y508	P07948	• PYK															Kinexus	PK688	RpAb
MAFG	S124+Pan	O15525			●													Kinexus	PN617	RpAb
MAK	Pan	P20794	• PSTK															No	AP7542b	RpAb
MAK	T157	P20794	• PSTK															Kinexus	PK689	RpAb
Malin (NHLRC1; EPM2B)	Pan	Q6VVV1							●									No	NN288-1	MmAb
MAP2	T1613+T161	P11137												●				Kinexus	PN834	RpAb
MAP2	S1782	P11137												●				Kinexus	PN835	RpAb
MAPKAPK2 (RPS6KC1)	Pan	P49137	• PSTK															Kinexus	NK097-2	RpAb
MAPKAPK2 (RPS6KC1)	Pan	P49137	• PSTK															No	sc-6221	GpAb
MAPKAPK2 (RPS6KC1)	Pan	P49137	• PSTK															No	sc-7871	RpAb
MAPKAPK2 (RPS6KC1)	T222	P49137	• PSTK															Kinexus	PK690	RpAb
MAPKAPK2 (RPS6KC1)	Y225+T226	P49137	• PSTK															Kinexus	PK691	RpAb
MAPKAPK3	Y76	Q16644	• PSTK															Kinexus	PK692	RpAb
MAPKAPK5 (PRAK)	Pan	Q8IW41	• PSTK															No	sc-8251	GpAb
MAPKAPK5 (PRAK)	T186	Q8IW41	• PSTK															Kinexus	PK693	RpAb
MARK1	Pan	Q9P0L2	• PSTK															Kinexus	NK098-2	RpAb
MARK1	T215	Q9P0L2	• PSTK															Kinexus	PK694	RpAb
MARK1	T215	Q9P0L2	• PSTK															Kinexus	PK695	RpAb
MARK2	Pan	Q7KZ17	• PSTK															Kinexus	NK275-2	RpAb
MARK2	Pan	Q7KZ17	• PSTK															Kinexus	NK275-1	RpAb
MARK3	Pan	P27448	• PSTK															Kinexus	NK276-2	RpAb
MARK3	Pan	P27448	• PSTK															Kinexus	NK276-1	RpAb
MARK3	T507	P27448	• PSTK															Kinexus	PK697	RpAb
MARK4	Pan	Q96L34	• PSTK															Kinexus	NK277-2	RpAb
MARK4	Pan	Q96L34	• PSTK															Kinexus	NK277-1	RpAb
MAT1A	T341	Q00266								●								Kinexus	PN759	RpAb
MATR3	T130	P43243																Kinexus	PN887	RpAb
MATR3	S188	P43243																Kinexus	PN883	RpAb
MATR3	Y219	P43243																Kinexus	PN888	RpAb
MATR3	S248+Y250	P43243																Kinexus	PN884	RpAb
MATR3	S533	P43243																Kinexus	PN885	RpAb
MATR3	S596+S598	P43243																Kinexus	PN886	RpAb
MBP	Y203	P02686												●				Kinexus	PN968	RpAb
MBP	T232	P02686												●				Kinexus	PN558	RpAb
Mcl1	Pan	Q07820										●						No	sc-819	RpAb
MCM2	S40+S41	P49736			●													Kinexus	PN618	RpAb
MCM2	Y137+S139	P49736			●													Kinexus	PN620	RpAb
MDC1 (NFB1)	Pan	Q14676												●				No	NN289-1	MmAb
MDM2	S166	Q00987							●									No	11550	RpAb
MEF2A	Pan	Q02078			●													No	sc-313	RpAb
MEF2A	T108	Q02078			●													Kinexus	PN622	RpAb
MEF2C	S396	Q06413			●													Kinexus	PN623	RpAb
MEF2D	S121	Q14814			●													Kinexus	PN624	RpAb
MEF2D	S180	Q14814			●													Kinexus	PN625	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															Kinexus	NK099-8	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															Kinexus	NK099-9	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															Kinexus	NK099-7	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															Kinexus	NK099-3	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															Kinexus	NK099-4	RpAb
MEK1 (MAP2K1; MKK1)	Pan	Q02750	• DSK															No	610122	MmAb
MEK1 (MAP2K1; MKK1)	S222+Pan	Q02750	• DSK															Kinexus	PK698	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
MEK1 (MAP2K1; MKK1)	T286	Q02750	• DSK															Kinexus	PK886	RpAb
MEK1 (MAP2K1; MKK1)	S298	Q02750	• DSK															No	KAP-MA020	RpAb
MEK2 (MAP2K2; MKK2)	Pan	P36507	• DSK															No	AP7905a	RpAb
MEK2 (MAP2K2; MKK2)	Pan	P36507	• DSK															Kinexus	NK100-5	RpAb
MEK2 (MAP2K2; MKK2)	Pan	P36507	• DSK															Kinexus	NK100-4	RpAb
MEK2 (MAP2K2; MKK2)	Pan	P36507	• DSK															Kinexus	NK100-6	RpAb
MEK2 (MAP2K2; MKK2)	Pan	P36507	• DSK															No	610236	MmAb
MEK5 (MAP2K5; MKK5)	Pan	Q13163	• DSK															Kinexus	NK104-4	RpAb
MEK5 (MAP2K5; MKK5)	Pan	Q13163	• DSK															Kinexus	NK104-5	RpAb
MEK5 (MAP2K5; MKK5)	Pan	Q13163	• DSK															Kinexus	NK104-2	RpAb
MEK5 (MAP2K5; MKK5)	Pan	Q13163	• DSK															Kinexus	NK104-3	RpAb
MEK5 (MAP2K5; MKK5)	Pan	Q13163	• DSK															No	sc-1287	GpAb
MEK5 (MAP2K5; MKK5)	S311	Q13163	• DSK															Kinexus	PK699	RpAb
MEKK1 (MAP3K1)	Pan	Q13233	• PSTK															No	AP7907a	RpAb
MEKK1 (MAP3K1)	Pan	Q13233	• PSTK															Kinexus	NK107-3	RpAb
MEKK1 (MAP3K1)	Pan	Q13233	• PSTK															No	sc-252	RpAb
MEKK2 (MAP3K2)	Pan	Q9Y2U5	• PSTK															No	AP7908a	RpAb
MEKK2 (MAP3K2)	Pan	Q9Y2U5	• PSTK															Kinexus	NK108-3	RpAb
MEKK2 (MAP3K2)	Pan	Q9Y2U5	• PSTK															Kinexus	NK108-4	RpAb
MEKK2 (MAP3K2)	Pan	Q9Y2U5	• PSTK															Kinexus	NK108-5	RpAb
MEKK2 (MAP3K2)	Pan	Q9Y2U5	• PSTK															No	sc-1089	RpAb
MEKK2 (MAP3K2)	S239+Pan	Q9Y2U5	• PSTK															Kinexus	PK700	RpAb
MEKK4 (MAP3K4)	Pan	Q9Y6R4	• PSTK															No	AP7910a	RpAb
MEKK6 (MAP3K6; ASK2)	Pan	O95382	• PSTK															Kinexus	NK225-2	RpAb
MEKK6 (MAP3K6; ASK2)	Pan	O95382	• PSTK															Kinexus	NK225-3	RpAb
MEKK6 (MAP3K6; ASK2)	Pan	O95382	• PSTK															Kinexus	NK225-4	RpAb
MELK	Pan	Q14680	• PSTK															Kinexus	NK229-3	RpAb
MELK	Pan	Q14680	• PSTK															Kinexus	NK229-4	RpAb
MELK	Pan	Q14680	• PSTK															Kinexus	NK229-2	RpAb
MELK	Y438	Q14680	• PSTK															Kinexus	PK701	RpAb
MERTK (MER)	Y749	Q12866	• PYK			●												Kinexus	PK702	RpAb
MERTK (MER)	Y749+Y753	Q12866	• PYK			●												Kinexus	PK703	RpAb
MERTK (MER)	Y753	Q12866	• PYK			●												Kinexus	PK704	RpAb
Met (HGF receptor)	Pan	P08581	• PYK			●												Kinexus	NK110-3	RpAb
Met (HGF receptor)	Pan	P08581	• PYK			●												Kinexus	NK110-4	RpAb
Met (HGF receptor)	Pan	P08581	• PYK			●												Kinexus	NK110-2	RpAb
Met (HGF receptor)	Y1003	P08581	• PYK			●												Kinexus	PK708	RpAb
Met (HGF receptor)	Y1230+Pan	P08581	• PYK			●												Kinexus	PK709	RpAb
Met (HGF receptor)	Y1234	P08581	• PYK			●												Kinexus	PK710	RpAb
Met (HGF receptor)	S1236	P08581	• PYK			●												Kinexus	PK705	RpAb
Met (HGF receptor)	Y1234+Y123	P08581	• PYK			●												Kinexus	PK711	RpAb
Met (HGF receptor)	Y1234+Y123	P08581	• PYK			●												Kinexus	PK712	RpAb
Met (HGF receptor)	T1355+Y135	P08581	• PYK			●												Kinexus	PK707	RpAb
Met (HGF receptor)	T1241	P08581	• PYK			●												Kinexus	PK706	RpAb
MITF	S414	O75030			●													Kinexus	PN626	RpAb
MKK3 (MAP2K3; MEK3)	Pan	P46734	• DSK															Kinexus	NK101-4	RpAb
MKK3 (MAP2K3; MEK3)	Pan	P46734	• DSK															Kinexus	NK101-5	RpAb
MKK3 (MAP2K3; MEK3)	Pan	P46734	• DSK															Kinexus	NK101-6	RpAb
MKK3 (MAP2K3; MEK3)	Pan	P46734	• DSK															No	sc-959	RpAb
MKK3 (MAP2K3; MEK3)	Pan	P46734	• DSK															No	sc-960	RpAb
MKK3 (MAP2K3; MEK3)	S218	P46734	• DSK															Kinexus	PK713	RpAb
MKK3 (MAP2K3; MEK3)	Y230	P46734	• DSK															Kinexus	PK714	RpAb
MKK3 (MAP2K3; MEK3);	Pan	P46734	• DSK															No	sc-8407	MmAb
MKK4 (MAP2K4; MEK4)	Pan	P45985	• DSK															Kinexus	NK103-4	RpAb
MKK4 (MAP2K4; MEK4)	Pan	P45985	• DSK															Kinexus	NK103-5	RpAb
MKK4 (MAP2K4; MEK4)	Pan	P45985	• DSK															Kinexus	NK103-6	RpAb
MKK4 (MAP2K4; MEK4)	Pan	P45985	• DSK															No	sc-964	RpAb
MKK4 (MAP2K4; MEK4)	S257	P45985	• DSK															Kinexus	PK715	RpAb
MKK4 (MAP2K4; MEK4)	S80+Pan	P45985	• DSK															Kinexus	PK716	RpAb
MKK6 (MAP2K6; MEK6)	Pan	P52564	• DSK															Kinexus	NK105-4	RpAb
MKK6 (MAP2K6; MEK6)	Pan	P52564	• DSK															Kinexus	NK105-5	RpAb
MKK6 (MAP2K6; MEK6)	Pan	P52564	• DSK															Kinexus	NK105-3	RpAb
MKK7 (MAP2K7; MEK7)	Pan	O14733	• DSK															Kinexus	NK106-4	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
MKK7 (MAP2K7; MEK7)	Pan	O14733	• DSK															Kinexus	NK106-5	RpAb
MKK7 (MAP2K7; MEK7)	Pan	O14733	• DSK															Kinexus	NK106-3	RpAb
MKK7 (MAP2K7; MEK7)	T275	O14733	• DSK															Kinexus	PK717	RpAb
MLK1 (MAP3K9)	Pan	P80192	• PSTK															Kinexus	NK278-1	RpAb
MLK2 (MAP3K10)	Pan	Q02779	• PSTK															Kinexus	NK279-1	RpAb
MLK3 (MAP3K11)	S281	Q16584	• PSTK															Kinexus	PK718	RpAb
MLK3 (MAP3K11)	T277+S281	Q16584	• PSTK															Kinexus	PK870	RpAb
MLK4 (MAP3K21)	Pan	Q5TCX8	• PSTK															Kinexus	NK280-1	RpAb
MLL	S2196	Q03164																Kinexus	PN801	RpAb
MLTK	T161+T162	Q9NYL2	• PSTK															Kinexus	PK719	RpAb
Mnk1	T250+T255	Q9BUB5	• PSTK															No	2111L	RpAb
Mnk2	Pan	Q9HBH9	• PSTK															No	sc-7445	GpAb
MOK (RAGE)	Pan	Q9UQ07	• PSTK															Kinexus	NK281-2	RpAb
MOK (RAGE)	Pan	Q9UQ07	• PSTK															Kinexus	NK281-1	RpAb
MOK (RAGE)	Y167	Q9UQ07	• PSTK															Kinexus	PK721	RpAb
MOK (RAGE)	T159+Y161	Q9UQ07	• PSTK															Kinexus	PK720	RpAb
MOR1 (mu opioid receptor)	T372+S377+	P35372				●												Kinexus	PN802	RpAb
Mos	Pan	P00540	• PSTK															Kinexus	NK112	RpAb
Mos	Y263	P00540	• PSTK															Kinexus	PK722	RpAb
MPL	Y591	P40238				●												Kinexus	PN760	RpAb
MRCKa (PK428)	Pan	Q5VT25	• PSTK															Kinexus	NK282-1	RpAb
MRCKb (CDC42BPB)	Pan	Q9Y5S2	• PSTK															Kinexus	NK283-1	RpAb
MRLC1	T19+S20	P24844												●				Kinexus	PN836	RpAb
MSH6	S14	P52701									●							Kinexus	PN708	RpAb
MSK1 (RPS6KA5)	S212	O75582	• PSTK															Kinexus	PK723	RpAb
MSK1 (RPS6KA5)	S376	O75582	• PSTK															No	9595L	RpAb
MSK2 (RPS6KA4)	T194+S196	O75676	• PSTK															Kinexus	PK868	RpAb
MSK2 (RPS6KA4)	T687	O75676	• PSTK															Kinexus	PK725	RpAb
MST1 (STK4; Krs2)	Pan	Q13043	• PSTK															No	sc-6213	GpAb
MST1 (STK4; Krs2)	Pan	Q13043	• PSTK															No	M77120	MmAb
MST1 (STK4; Krs2)	T183	Q13043	• PSTK															Kinexus	PK871	RpAb
MST1 (STK4; Krs2)	T187	Q13043	• PSTK															Kinexus	PK726	RpAb
MST3 (STK24)	T184+Pan	Q9Y6E0	• PSTK															Kinexus	PK727	RpAb
MST3 (STK24)	T190	Q9Y6E0	• PSTK															Kinexus	PK728	RpAb
mTOR (FRAP)	Pan	P42345	• PSTK															Kinexus	NK116-4	RpAb
mTOR (FRAP)	Pan	P42345	• PSTK															Kinexus	NK116-3	RpAb
mTOR (FRAP)	Pan	P42345	• PSTK															Kinexus	NK116-5	RpAb
mTOR (FRAP)	S2448	P42345	• PSTK															Kinexus	PK729	RpAb
mTOR (FRAP)	S2478+S248	P42345	• PSTK															Kinexus	PK730	RpAb
MuSK	Pan	O15146	• PYK			●												No	sc-6009	GpAb
MuSK	Y554	O15146	• PYK			●												Kinexus	PK953	RpAb
MuSK	Y756	O15146	• PYK			●												Kinexus	PK872	RpAb
Myb	S532+Pan	P10242			●													Kinexus	PN627	RpAb
Myc (c-Myc)	T58	P01106			●													No	11034	RpAb
Myc (c-Myc)	T58+S62	P01106			●													Kinexus	PN628	RpAb
Myf-5	Pan	P13349			●													No	sc-302	RpAb
MYT1 (PLPB1)	S143	Q99640	• PSTK															Kinexus	PK887	RpAb
NALCN (VGCNL1)	Pan	Q8IZF0					●											No	NN293-1	MmAb
Nav1.2	S568	Q99250					●											Kinexus	PN837	RpAb
NCAM1 (CD56)	S784	P13591													●			Kinexus	PN969	RpAb
NCOA3 (SRC-3)	S867	Q9Y6Q9			●													Kinexus	PN629	RpAb
NDR1 (NDR; STK38)	S281+T282	Q15208	• PSTK															Kinexus	PK731	RpAb
Nek1	S1052	Q96PY6	• PSTK															Kinexus	PK940	RpAb
Nek2	Pan	P51955	• PSTK															No	NK117-5	RpAb
Nek2	Pan	P51955	• PSTK															No	sc-7440	GpAb
Nek2	Pan	P51955	• PSTK															No	sc-7439	GpAb
Nek2	T170+S171	P51955	• PSTK															Kinexus	PK733	RpAb
Nek2	S171	P51955	• PSTK															Kinexus	PK732	RpAb
Nek4	Pan	P51957	• PSTK															No	AP8076b	RpAb
Nek6	S206	Q9HC98	• PSTK															Kinexus	PK734	RpAb
Nek7	T191+S195	Q8TDX7	• PSTK															Kinexus	PK735	RpAb
NF-H (NEFH)	Pan	P12036												●				Kinexus	NN583-1	RpAb
NF-H (NEFH)	S532	P12036												●				Kinexus	PN970	RpAb



Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
p38a MAPK (MAPK14; CSBP;	T180+Y182	Q16539	• PSTK															Kinexus	PK739	RpAb
p38a MAPK (MAPK14; CSBP;	T180+Y182	Q15759	• PSTK															Kinexus	PK740	RpAb
p38a MAPK (MAPK14; CSBP;	Y182	Q15759	• PSTK															Kinexus	PK959	RpAb
p38b MAPK (MAPK11)	Pan	Q15759	• PSTK															Kinexus	NK248-2	RpAb
p38b MAPK (MAPK11)	Pan	Q15759	• PSTK															Kinexus	NK248-1	RpAb
p38b MAPK (MAPK11)	T180+Y182	Q15759	• PSTK															Kinexus	PK741-2	RpAb
p38b MAPK (MAPK11)	T180+Y182	Q15759	• PSTK															Kinexus	PK741	RpAb
p38d MAPK (MAPK13)	Pan	O15264	• PSTK															Kinexus	NK121-3	RpAb
p38d MAPK (MAPK13)	Pan	O15264	• PSTK															Kinexus	NK121-4	RpAb
p38d MAPK (MAPK13)	Pan	O15264	• PSTK															Kinexus	NK121-2	RpAb
p38d MAPK (MAPK13)	T180+Y182	O15264	• PSTK															Kinexus	PK960	RpAb
p38d MAPK (MAPK13)	Y182	O15264	• PSTK															Kinexus	PK743	RpAb
p38d MAPK (MAPK13)	S261+T265	O15264	• PSTK															Kinexus	PK888	RpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															No	NK059-2	RpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															Kinexus	NK059-3	RpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															Kinexus	NK059-4	RpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															Kinexus	NK059-5	RpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															No	sc-6023	GpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															No	sc-6022	GpAb
p38g MAPK (MAPK12; ERK6;	Pan	P53778	• PSTK															No	07-139	RpAb
p70S6K (S6Ka; RPS6KB1)	Pan	P23443	• PSTK															Kinexus	NK223-2	RpAb
p70S6K (S6Ka; RPS6KB1)	Pan	P23443	• PSTK															Kinexus	NK223-4	RpAb
p70S6K (S6Ka; RPS6KB1)	Pan	P23443	• PSTK															No	sc-230	RpAb
p70S6K (S6Ka; RPS6KB1)	T252	P23443	• PSTK															Kinexus	PK744	RpAb
p70S6K (S6Ka; RPS6KB1)	T412	P23443	• PSTK															No	9205L	RpAb
p70S6K (S6Ka; RPS6KB1)	T412	P23443	• PSTK															Kinexus	PK745	RpAb
p70S6K (S6Ka; RPS6KB1)	S434	P23443	• PSTK															No	11269	RpAb
p70S6K (S6Ka; RPS6KB1)	T444+S447	P23443	• PSTK															Kinexus	PK746	RpAb
p70S6K (S6Ka; RPS6KB1)	S447	P23443	• PSTK															No	11284	RpAb
p70S6Kb (S6Kb2; RPS6KB2)	S423	Q9UBS0	• PSTK															Kinexus	PK747	RpAb
p73 (TRP73)	Pan	O15350			●													No	KAM-CC24	MmAb
p73 (TRP73)	Y99	O15350			●													Kinexus	PN861	RpAb
PAH	S16	P00439								●								Kinexus	PN713	RpAb
PAK1 (PAKa)	Pan	Q13153	• PSTK															Kinexus	NK122-2	RpAb
PAK1 (PAKa)	Pan	Q13153	• PSTK															No	sc-881	RpAb
PAK1 (PAKa)	S144	Q13153	• PSTK															Kinexus	PK748	RpAb
PAK1 (PAKa)	T212	Q13153	• PSTK															No	Y011154	RpAb
PAK1 (PAKa)	T423	Q13153	• PSTK															Kinexus	PK749	RpAb
PAK2 (PAKg)	Pan	Q13177	• PSTK															Kinexus	NK200	RpAb
PAK2 (PAKg)	Pan	Q13177	• PSTK															No	sc-1872	GpAb
PAK2 (PAKg)	Y130	Q13177	• PSTK															Kinexus	PK751	RpAb
PAK2 (PAKg)	S141	Q13177	• PSTK															Kinexus	PK750	RpAb
PAK4	S474	O96013	• PSTK															Kinexus	PK752	RpAb
PAK5 (PAK7)	Pan	Q9P286	• PSTK															Kinexus	NK190-3	RpAb
PAK5 (PAK7)	Pan	Q9P286	• PSTK															Kinexus	NK190-2	RpAb
PAK5 (PAK7)	S602	Q9P286	• PSTK															Kinexus	PK753	RpAb
PAK6	Pan	Q9NQU5	• PSTK															Kinexus	NK124-2	RpAb
Paxillin 1 (PXN; PAX1)	Y118	P49023												●				No	44-722G	RpAb
PBK (TOPK)	Pan	Q96KB5	• PSTK															Kinexus	NK284-1	RpAb
PBK (TOPK)	Y74	Q96KB5	• PSTK															Kinexus	PK754	RpAb
PBK (TOPK)	Y272	Q96KB5	• PSTK															Kinexus	PK889	RpAb
PBRM1	S303+Pan	Q86U86			●													Kinexus	PN767	RpAb
PBRM1	S948	Q86U86			●													Kinexus	PN714	RpAb
PCDH-gamma-A3	Pan	Q9Y5H0												●				No	NN308-2	MmAb
PCDH-gamma-C3 (PCDH2)	Pan	Q9UN70												●				No	NN304-1	MmAb
PCK1	S118	P35558								●								Kinexus	PN768	RpAb
PCTAIRE1 (CDK16; PCTK1)	Pan	Q00536	• PSTK															No	AP7547b	RpAb
PCTAIRE1 (CDK16; PCTK1)	Y176	Q00536	• PSTK															Kinexus	PK755	RpAb
PCTK2 (PCTAIRE2; CDK17)	Pan	Q00537	• PSTK															Kinexus	NK285-1	RpAb
PCTK2 (PCTAIRE2; CDK17))	S180	Q00537	• PSTK															Kinexus	PK756	RpAb
PCTK3 (PCTAIRE3; CDK18)	Pan	Q07002	• PSTK															Kinexus	NK286-1	RpAb
PCYT1A (CTPCT; CCTA)	Pan	P49585								●								Kinexus	NN456-2	RpAb
PCYT1A (CTPCT; CCTA)	Pan	P49585								●								Kinexus	NN606-1	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
PCYT1A (CTPCT; CCTA)	S329+S331	P49585								●								Kinexus	PN561	RpAb
PCYT1A (CTPCT; CCTA)	T342+S343	P49585								●								Kinexus	PN547	RpAb
PCYT1A (CTPCT; CCTA)	Y359+S362	P49585								●								Kinexus	PN548	RpAb
PCYT1B (CCTB)	Pan	Q9Y5K3								●								Kinexus	NN605-1	RpAb
PCYT1B (CCTB)	S315+S319	Q9Y5K3								●								Kinexus	PN546	RpAb
PDGFRA	Pan	P16234	• PYK			●												Kinexus	NK242-2	RpAb
PDGFRA	Pan	P16234	• PYK			●												Kinexus	NK242-1	RpAb
PDGFRA	Y762	P16234	• PYK			●												Kinexus	PK758	RpAb
PDGFRA	Y768	P16234	• PYK			●												Kinexus	PK759	RpAb
PDGFRA	S847+Y849	P16234	• PYK			●												Kinexus	PK757	RpAb
PDGFRB	Pan	P09619	• PYK			●												Kinexus	NK243-3	RpAb
PDGFRB	Pan	P09619	• PYK			●												Kinexus	NK243-1	RpAb
PDGFRB	Y716	P09619	• PYK			●												No	07-021	RpAb
PDK1 (PDHK1)	Pan	Q15118	• PSTK															Kinexus	NN179-2	RpAb
PDK1 (PDHK1)	Pan	Q15118	• PSTK															Kinexus	NN179-1	RpAb
PDK1 (PDPK1)	Pan	O15530	• PSTK															Kinexus	NK126-3	RpAb
PDK1 (PDPK1)	Pan	O15530	• PSTK															Kinexus	NK126-4	RpAb
PDK1 (PDPK1)	S241	O15530	• PSTK															Kinexus	PK760	RpAb
PDK2 (PDHK2; PyDK2)	Pan	Q15119	• PSTK															Kinexus	NN180-2	RpAb
PDK2 (PDHK2; PyDK2)	Pan	Q15119	• PSTK															Kinexus	NN180-1	RpAb
PDK3 (PDHK3)	Pan	Q15120	• PSTK															Kinexus	NN181-2	RpAb
PDK4 (PDHK4)	Pan	Q16654	• PSTK															Kinexus	NN178-3	RpAb
PDK4 (PDHK4)	Pan	Q16654	• PSTK															Kinexus	NN178-2	RpAb
PDLIM5 (LIM)	Y251	Q96HC4											●					Kinexus	PN522	RpAb
PECAM-1	Y713	P16284													●			Kinexus	PN523	RpAb
PEX6 (Peroxin-6; PXAAA1)	Pan	Q13608																No	NN305-1	RpAb
PFKFB3	S461	Q16875								●								Kinexus	PN715	RpAb
PFKFB3	S467	Q16875								●								Kinexus	PN769	RpAb
PFKL	S775	P17858								●								Kinexus	PN770	RpAb
PFKP	S386	Q01813								●								Kinexus	PN716	RpAb
PFKP	Y651+Pan	Q01813								●								Kinexus	PN771	RpAb
PFN1	Y60	P07737												●				Kinexus	PN898	RpAb
PFN1	Y129	P07737												●				Kinexus	PN524	RpAb
PGK1	Y196	P00558								●								Kinexus	PN525	RpAb
PHLPP1 (PHLPP)	Pan	O60346	• PSTP															Kinexus	NP049-1	RpAb
PHLPP1 (PHLPP)	Y1712+Y171	O60346	• PSTP															Kinexus	PP529	RpAb
PHLPP2	Pan	Q6ZVD8	• PSTP															Kinexus	NP050-1	RpAb
PHYH (Phytanic acid oxidase;)	Pan	O14832								●								No	NN306-1	MmAb
PI3KR4	Pan	Q99570	• PSTK															No	AP8026a	RpAb
PI4KCB	Pan	Q9UBF8								●								No	AP8030b	RpAb
Piccolo (Pclo)	Pan	Q9Y6V0												●				No	NN307-2	RpAb
PIK3CA	Y317	P42336	• PI-K															Kinexus	PK894	RpAb
PIK3R1 (PI3K p85)	Y467	P27986	• PI-K															Kinexus	PN526	RpAb
PIK3R1 (PI3K p85)	Y580	P27986	• PI-K															Kinexus	PN527	RpAb
PIK3R2	Y464	O00459	• PI-K															Kinexus	PN528	RpAb
Pim1	Pan	P11309	• PSTK															Kinexus	NK258-2	RpAb
Pim2	Pan	Q9P1W9	• PSTK															Kinexus	NK288-2	RpAb
Pim2	Pan	Q9P1W9	• PSTK															Kinexus	NK288-1	RpAb
Pim2	T195	Q9P1W9	• PSTK															Kinexus	PK761	RpAb
Pim3	Pan	Q86V86	• PSTK															Kinexus	NK289-1	RpAb
PINK1 (BRPK)	Pan	Q9BXM7	• PSTK															No	NK304-1	MmAb
PINK1 (BRPK)	S228	Q9BXM7	• PSTK															Kinexus	PK906	RpAb
PINK1 (BRPK)	T257	Q9BXM7	• PSTK															Kinexus	PK942	RpAb
PINK1 (BRPK)	T313	Q9BXM7	• PSTK															Kinexus	PK907	RpAb
PINK1 (BRPK)	S402	Q9BXM7	• PSTK															Kinexus	PK908	RpAb
PIP5K	S307	Q9Y217	• PSTK															Kinexus	PK762	RpAb
PIP5K	Y1772+Y177	Q9Y217	• PSTK															Kinexus	PK943	RpAb
PIP5K2a (PIP5K2A)	Pan	P48426	• PI-K															No	NK209	RpAb
PKCa (PRKCA)	Pan	P17252	• PSTK															Kinexus	NK201	RpAb
PKCa (PRKCA)	Pan	P17252	• PSTK															No	P16520	MmAb
PKCa (PRKCA)	Y195	P17252	• PSTK															Kinexus	PK764	RpAb
PKCa (PRKCA)	S226+T228	P17252	• PSTK															Kinexus	PK944	RpAb
PKCa (PRKCA)	T497	P17252	• PSTK															Kinexus	PK763	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
PKCa (PRKCA)	T638	P17252	• PSTK															Kinexus	PK909	RpAb
PKCa (PRKCA)	T638/T641	P17252	• PSTK															No	9375L	RpAb
PKCb (PRKCB1)	Pan	P05771	• PSTK															Kinexus	NK133-2	RpAb
PKCb (PRKCB1)	Pan	P05771	• PSTK															No	sc-209	RpAb
PKCb (PRKCB1)	T500	P05771	• PSTK															Kinexus	PK766	RpAb
PKCb (PRKCB1)	S661	P05771	• PSTK															Kinexus	PK765	RpAb
PKCb2 (PRKCB2)	Pan	P05771	• PSTK															No	sc-210	RpAb
PKCd (PRKCD)	Pan	Q05655	• PSTK															No	sc-213	RpAb
PKCd (PRKCD)	Y313	Q05655	• PSTK															Kinexus	PK768	RpAb
PKCd (PRKCD)	Y334	Q05655	• PSTK															Kinexus	PK769	RpAb
PKCd (PRKCD)	T507	Q05655	• PSTK															Kinexus	PK767	RpAb
PKCd (PRKCD)	S664	Q05655	• PSTK															No	44-976G	RpAb
PKCe (PRKCE)	Pan	Q02156	• PSTK															No	sc-214	RpAb
PKCe (PRKCE)	Pan	Q02156	• PSTK															No	sc-214-G	GpAb
PKCg (PRKCG)	Pan	P05129	• PSTK															No	sc-211	RpAb
PKCg (PRKCG)	T514	P05129	• PSTK															No	KAP-PK011	RpAb
PKCg (PRKCG)	T655	P05129	• PSTK															No	44-965	RpAb
PKCh (PRKCH; PRKCeta)	Pan	P24723	• PSTK															No	sc-215	RpAb
PKCi (PRKCiota)/PKCI	Pan	P41743	• PSTK															No	sc-32020	GpAb
PKCt (PRKCQ; PKC-q; PKC-	Y545	Q04759	• PSTK															Kinexus	PK773	RpAb
PKCt (PRKCQ; PKC-q; PKC-	S676	Q04759	• PSTK															No	11297	RpAb
PKCt (PRKCQ; PKC-q; PKC-	S685	Q04759	• PSTK															Kinexus	PK961	RpAb
PKCt (PRKCQ; PKC-q; PKC-	S695	Q04759	• PSTK															Kinexus	PK772	RpAb
PKCt (PRKCQ; PKC-q; PKC-	S695	Q04759	• PSTK															No	11173	RpAb
PKCz (PRKCZ)	Pan	Q05513	• PSTK															No	sc-216	RpAb
PKCz (PRKCZ)	S262+Y263	Q05513	• PSTK															Kinexus	PK774	RpAb
PKCz (PRKCZ)	T410	Q05513	• PSTK															Kinexus	PK775	RpAb
PKD1 (PRKCM; PKCm;	Pan	Q15139	• PSTK															No	sc-639	RpAb
PKD1 (PRKCM; PKCm;	S205	Q15139	• PSTK															Kinexus	PK770	RpAb
PKD1 (PRKCM; PKCm;	S738+S742	Q15139	• PSTK															Kinexus	PK771	RpAb
PKD1 (PRKCM; PKCm;	S910	Q15139	• PSTK															Kinexus	PK912	RpAb
PKD2 (PRKD2)	S197+S198	Q9BZL6	• PSTK															Kinexus	PK784	RpAb
PKD3 (PRKCN)	Pan	O94806	• PSTK															Kinexus	NK139-2	RpAb
PKG1a (PRKG1A)	Pan	Q13976	• PSTK															Kinexus	NK202-2	RpAb
PKG1a (PRKG1A)	Pan	Q13976	• PSTK															Kinexus	NK202	RpAb
PKG1a (PRKG1A)	Pan	Q13976	• PSTK															No	sc-10344	GpAb
PKG1a (PRKG1A)	T515+T517	Q13976	• PSTK															Kinexus	PK776	RpAb
PKG1a (PRKG1A)	T517	Q13976	• PSTK															Kinexus	PK776-2	RpAb
PKG2 (PRKG2)	Pan	Q13237	• PSTK															Kinexus	NK290-2	RpAb
PKG2 (PRKG2)	Pan	Q13237	• PSTK															Kinexus	NK290-1	RpAb
PKM2	S37	P14618								●								Kinexus	PN718	RpAb
PKM2	Y105	P14618								●								Kinexus	PN772	RpAb
PKM2	Y148	P14618								●								Kinexus	PN717	RpAb
PKM2	Y370	P14618								●								Kinexus	PN773	RpAb
PKM2	Y390	P14618								●								Kinexus	PN529	RpAb
PKN1 (PRK1)	Pan	Q16512	• PSTK															No	sc-1842	GpAb
PKN1 (PRK1)	T774	Q16512	• PSTK															Kinexus	PK781	RpAb
PKN2 (PRK2)	Pan	Q16513	• PSTK															Kinexus	NK149-3	RpAb
PKR1 (PRKR; EIF2AK2)	Pan	P19525	• PSTK															No	sc-6282	MmAb
PKR1 (PRKR; EIF2AK2)	T446+Pan	P19525	• PSTK															Kinexus	PK777	RpAb
PKR1 (PRKR; EIF2AK2)	T451	P19525	• PSTK															Kinexus	PK910	RpAb
PLCB3	S1105	Q01970								●								Kinexus	PN719	RpAb
PLCD1	S460	P51178								●								Kinexus	PN720	RpAb
PLCE1	S1096+T110	Q9P212								●								Kinexus	PN721	RpAb
PLCG1	Y771	P19174								●								No	11523	RpAb
PLCG1	Y771+Y775	P19174								●								Kinexus	PN774	RpAb
PLCG1	Y783	P19174								●								Kinexus	PN530	RpAb
PLCG1	Y977	P19174								●								Kinexus	PN722	RpAb
PLCG2 (PLC R)	Pan	P16885								●								No	sc-407	RpAb
PLCG2 (PLC R)	Y753	P16885								●								Kinexus	PN723	RpAb
PLCG2 (PLC R)	Y759	P16885								●								Kinexus	PN775	RpAb
PLCG2 (PLC R)	Y759	P16885								●								Kinexus	PN531	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
PLCG2 (PLC R)	Y1217	P16885								●								Kinexus	PN776	RpAb
Plk1 (PLK)	Pan	P53350	• PSTK															Kinexus	NK145-2	RpAb
Plk1 (PLK)	Pan	P53350	• PSTK															No	33-1700	RpAb
Plk1 (PLK)	T210	P53350	• PSTK															Kinexus	PK778	RpAb
Plk1 (PLK)	T210	P53350	• PSTK															Kinexus	PK778-2	RpAb
Plk1 (PLK)	Y217	P53350	• PSTK															Kinexus	PK779	RpAb
Plk2 (SNK)	Pan	Q9NYY3	• PSTK															No	NK146	RpAb
Plk2 (SNK)	Pan	Q9NYY3	• PSTK															No	sc-9577	GpAb
Plk3 (CNK; FNK)	Pan	Q9H4B4	• PSTK															No	AP3213a	RpAb
Plk3 (CNK; FNK)	Pan	Q9H4B4	• PSTK															Kinexus	NK147-2	RpAb
Plk3 (CNK; FNK)	Pan	Q9H4B4	• PSTK															No	sc-9578	GpAb
Plk3 (CNK; FNK)	Pan	Q9H4B4	• PSTK															No	sc-9580	GpAb
Plk4 (SAK; STK18)	Pan	O00444	• PSTK															Kinexus	NK291-1	RpAb
Plk4 (SAK; STK18)	T170	O00444	• PSTK															Kinexus	PK780	RpAb
PML	S518+Pan	P29590			●													Kinexus	PN641	RpAb
POU2F1	S385	P14859			●													Kinexus	PN643	RpAb
PPARg-1	S112	P37231			●	●												Kinexus	PN644	RpAb
PPFIBP1	S599+S601	Q86W92											●					Kinexus	PN777	RpAb
PPP1CA (PP1/Ca)	Pan	P62136	• PSTP															No	sc-7482	MmAb
PPP1CA (PP1/Ca)	Y306	P62136	• PSTP															Kinexus	PP542	RpAb
PPP1CA (PP1/Ca)	Y134+Y137	P62136	• PSTP															Kinexus	PP541	RpAb
PPP1CB (PP1/Cb)	Y306	P62140	• PSTP															Kinexus	PP530	RpAb
PPP1CB (PP1/Cb)	T316	P62140	• PSTP															Kinexus	PP502	RpAb
PPP1R11 (HCG V)	Y64	O60927	• PSTP-															Kinexus	PN532	RpAb
PPP1R12A (MYPT1; MBS)	Y496	O14974	• PSTP-															Kinexus	PP543	RpAb
PPP1R12A (MYPT1; MBS)	Y766	O14974	• PSTP-															Kinexus	PP544	RpAb
PPP1R12B (MYPT2)	T646	O60237	• PSTP-															Kinexus	PP503	RpAb
PPP1R14B (PHI-1)	Y29	Q96C90	• PSTP-															Kinexus	PP546	RpAb
PPP1R14B (PHI-1)	T57	Q96C90	• PSTP-															Kinexus	PP545	RpAb
PPP1R16A (MYPT3)	Y434	Q96I34	• PSTP-															Kinexus	PP547	RpAb
PPP1R1A (I-1; Inhibitor-1)	T35	Q13522	• PSTP-															Kinexus	PP548	RpAb
PPP2CA (PP2A/Ca)	Pan	P67775	• PSTP															No	P47720	MmAb
PPP2CA (PP2A/Ca)	Y284	P67775	• PSTP															Kinexus	PP549	RpAb
PPP2CA (PP2A/Ca)	Y307	P67775	• PSTP															Kinexus	PP504	RpAb
PPP2CB	T304	P62714	• PSTP															Kinexus	PP505	RpAb
PPP2R4 (PP2A subunit B';	Y223	Q15257	• PSTP-															Kinexus	PP550	RpAb
PPP2R5D (2A5D; B56D)	Y580	Q14738	• PSTP-															Kinexus	PP551	RpAb
PPP2R5E	Y99	Q16537	• PSTP-															Kinexus	PP552	RpAb
PPP3CA (PP2B-Aa;	Pan	Q08209	• PSTP															No	sc-17808	MmAb
PPP3CC (Calcinerin Ag)	S463+Pan	P48454	• PSTP															Kinexus	PP506	RpAb
PPP4R1 (PP4/A'2 (PPP4/A'2)	Pan	Q8TF05	• PSTP-															No	P215P	RpAb
PPP5C (PP5C; PP5; PPT)	Y119	P53041	• PSTP															Kinexus	PP507	RpAb
PPP6C	Y261	O00743	• PSTP															Kinexus	PP553	RpAb
PPPM1A (PP2Ca;	Y362	P35813	• PSTP															Kinexus	PP508	RpAb
PPPM1B (PP2Cb; PPM1B)	Y367	O75688	• PSTP															Kinexus	PP540	RpAb
PRKACA (PKA; PKACA)	Pan	P17612	• PSTK															No	610981	MmAb
PRKACA/B (PKA; PKACA/B)	T196+T198	P17612	• PSTK															Kinexus	PK782	RpAb
PRKACB (PKA)	Y69	P22694	• PSTK															Kinexus	PK783	RpAb
PRKACB (PKA)	S339	P22694	• PSTK															No	44-992G	RpAb
PRKAR2B)	Pan	P31323	• PSTK															No	AP8096a	RpAb
PRKX	Pan	P51817	• PSTK															Kinexus	NK292-1	RpAb
PRKX	T201+T203	P51817	• PSTK															Kinexus	PK785	RpAb
PRKY	Pan	O43930	• PSTK															Kinexus	NK293-1	RpAb
PRMT5	T634	O14744																Kinexus	PN549	RpAb
PRP4K (PRP4; PRPF4B)	Y849	Q13523	• PSTK															Kinexus	PK786	RpAb
PSEN1	T354	P49768							●									Kinexus	PN977	RpAb
PTEN	Pan	P60484	• PI-P															Kinexus	NP023-3	RpAb
PTEN	Pan	P60484	• PI-P															Kinexus	NP023-4	RpAb
PTEN	Pan	P60484	• PI-P															Kinexus	NP023-5	RpAb
PTEN	Pan	P60484	• PI-P															No	sc-7974	MmAb
PTEN	Y315	P60484	• PI-P															Kinexus	PP531	RpAb
PTEN	Y336	P60484	• PI-P															Kinexus	PP556	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
PTPN1 (PTP1B)	Pan	P18031		• PYP														No	P18020	MmAb
PTPN1 (PTP1B)	Y46	P18031		• PYP														Kinexus	PP533	RpAb
PTPN1 (PTP1B)	S50	P18031		• PYP														Kinexus	PP509	RpAb
PTPN1 (PTP1B)	Y66	P18031		• PYP														Kinexus	PP510	RpAb
PTPN1 (PTP1B)	Y20+Pan	P18031		• PYP														Kinexus	PP532	RpAb
PTPN2 (TC-PTP)	S304	P17706		• PYP														Kinexus	PP515	RpAb
PTPN3 (PTP-H1)	S357+S359	P26045		• PYP														Kinexus	PP519	RpAb
PTPN6 (PTP1C; SH-PTP1;	Y536	P29350		• PYP														Kinexus	PP520	RpAb
PTPN6 (PTP1C; SH-PTP1;	Y564	P29350		• PYP														Kinexus	PP538	RpAb
PTPN6 (PTP1C; SH-PTP1)	Pan	P29350		• PYP														No	sc-7289	MmAb
PTPN11 (PTP1D; PTP2C;	Pan	Q06124		• PYP														No	B50279-51	RpAb
PTPN11 (PTP1D; PTP2C;	Pan	Q06124		• PYP														No	B50270-51	RpAb
PTPN11 (PTP1D; PTP2C;	Y62	Q06124		• PYP														Kinexus	PP512	RpAb
PTPN11 (PTP1D; PTP2C;	Y546	Q06124		• PYP														Kinexus	PP534	RpAb
PTPN11 (PTP1D; PTP2C;	Y584	Q06124		• PYP														Kinexus	PP535	RpAb
PTPN12 (PTP-PEST; PTPG1)	S39	Q05209		• PYP														Kinexus	PP513	RpAb
PTPN12 (PTP-PEST; PTPG1)	S435	Q05209		• PYP														Kinexus	PP536	RpAb
PTPN14	S486+Pan	Q15678		• PYP														Kinexus	PP514	RpAb
PTPN21	S637	Q16825		• PYP														Kinexus	PP516	RpAb
PTPN22	Y499	Q9Y2R2		• PYP														Kinexus	PP517	RpAb
PTPN23 (HD-PTP)	Y1165	Q9H3S7		• PYP														Kinexus	PP518	RpAb
PTPN23 (HD-PTP)	Y1229	Q9H3S7		• PYP														Kinexus	PP537	RpAb
PTPRA (PTP-alpha)	Y798	P18433		• PYP		●												Kinexus	PP521	RpAb
PTPRB (PTP-beta)	Y1981	P23467		• PYP		●												Kinexus	PP522	RpAb
PTPRE (PTP-epsilon)	Y696	P23469		• PYP		●												Kinexus	PP554	RpAb
PTPRF (LAR)	Y1621	P10586		• PYP		●												Kinexus	PP523	RpAb
PTPRK (PTP-kappa)	Y858	Q15262		• PYP		●												Kinexus	PP539	RpAb
PTPRK (PTP-kappa)	Y916	Q15262		• PYP		●												Kinexus	PP524	RpAb
PTPRM (PTPRL1)	Y929	P28827		• PYP		●												Kinexus	PP526	RpAb
PTRF	Y308	Q6NZI2			●													Kinexus	PN646	RpAb
PTRF	Y308	Q6NZI2			●													Kinexus	PN533	RpAb
PU.1	S146	P17947			●													Kinexus	PN647	RpAb
PYK2 (PTK2B; FAK2)	Pan	Q14289	• PYK															No	AP7703a	RpAb
PYK2 (PTK2B; FAK2)	Pan	Q14289	• PYK															No	sc-1514	GpAb
PYK2 (PTK2B; FAK2)	Y402	Q14289	• PYK															Kinexus	PK788	RpAb
PYK2 (PTK2B; FAK2)	Y402	Q14289	• PYK															Kinexus	PK787	RpAb
PYK2 (PTK2B; FAK2)	Y579+Y580	Q14289	• PYK															Kinexus	PK789	RpAb
QKI (HKQ)	Pan	Q96PU8																No	NN311-1	MmAb
RAB1B	Pan	Q9H0U4																No	NN378-1	RpAb
RAB5A (Rab5)	Pan	P20339																No	NN150	RpAb
RACK1 (GNB2L1)	Pan	P63244											●					No	NN045	RpAb
Raf-A (ARaf)	Pan	P10398	• PSTK															Kinexus	NK205-3	RpAb
Raf-A (ARaf)	Pan	P10398	• PSTK															Kinexus	NK205-4	RpAb
Raf-A (ARaf)	Pan	P10398	• PSTK															Kinexus	NK205-5	RpAb
Raf-A (ARaf)	Pan	P10398	• PSTK															No	sc-408	RpAb
Raf-A (ARaf)	Y302	P10398	• PSTK															Kinexus	PK500	RpAb
Raf-B (B-Raf; BRaf; RafB)	Pan	P15056	• PSTK															Kinexus	NK156-4	RpAb
Raf-B (B-Raf; BRaf; RafB)	Pan	P15056	• PSTK															Kinexus	NK156-6	RpAb
Raf-B (B-Raf; BRaf; RafB)	Pan	P15056	• PSTK															No	sc-9002	RpAb
Raf-B (B-Raf; BRaf; RafB)	Pan	P15056	• PSTK															No	sc-5284	MmAb
Raf-B (B-Raf; BRaf; RafB)	Pan	P15056	• PSTK															No	/H000673M	MmAb
Raf-B (B-Raf; BRaf; RafB)	S446+S447	P15056	• PSTK															Kinexus	PK534	RpAb
Raf-B (B-Raf; BRaf; RafB)	S729	P15056	• PSTK															Kinexus	PK535	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															No	610151	MmAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															Kinexus	NK155-5	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															Kinexus	NK155-7	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															Kinexus	NK155-8	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															Kinexus	NK155-2	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															Kinexus	NK155-6	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															No	sc-376142	MmAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															No	sc-7198	RpAb
Raf1 (c-Raf; RafC)	Pan	P04049	• PSTK															No	sc-227	RpAb
Raf1 (c-Raf; RafC)	S259	P04049	• PSTK															Kinexus	PK790	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Raf1 (c-Raf; RafC)	S296	P04049	• PSTK															Kinexus	PK791	RpAb
Raf1 (c-Raf; RafC)	S301+T303	P04049	• PSTK															Kinexus	PK792	RpAb
Raf1 (c-Raf; RafC)	S338	P04049	• PSTK															Kinexus	PK913	RpAb
Raf1 (c-Raf; RafC)	Y340+Y341	P04049	• PSTK															Kinexus	PK914	RpAb
RARA	S77	P10276			●	●												Kinexus	PN648	RpAb
Rb (RB1)	S249+T252	P06400											●					Kinexus	PN724	RpAb
Rb (RB1)	T373	P06400											●					Kinexus	PN978	RpAb
Rb (RB1)	S612	P06400											●					No	44-572	RpAb
Rb (RB1)	S807+S811	P06400											●					Kinexus	PN847	RpAb
RBM9iso6	T7	O43251			●								●					Kinexus	PN649	RpAb
RDBP	S89+T91	P18615			●													Kinexus	PN650	RpAb
REEP1	Pan	Q9H902																No	NN312-1	MmAb
REEP2 (SGC32445)	Pan	Q9BRK0																No	NN312-2	MmAb
RelB (RELB)	S573	Q01201			●													No	Y011255	RpAb
Ret (GDNF receptor)	Pan	P07949	• PYK			●												Kinexus	NK244-2	RpAb
Ret (GDNF receptor)	Pan	P07949	• PYK			●												Kinexus	NK244-3	RpAb
Ret (GDNF receptor)	Pan	P07949	• PYK			●												Kinexus	NK244-1	RpAb
Ret (GDNF receptor)	Y905	P07949	• PYK			●												Kinexus	PK793	RpAb
Ret (GDNF receptor)	Y1062	P07949	• PYK			●												Kinexus	PK915	RpAb
Rhodopsin (OPN2; RHO;	Pan	P08100				●												No	NN313-2	MmAb
Rhodopsin (OPN2; RHO;	Pan	P08100				●												No	NN313-1	MmAb
Rictor	S21	Q6R327											●					Kinexus	PN979	RpAb
Rictor	T1135	Q6R327											●					Kinexus	PN980	RpAb
RIOK1	Y466	Q9BRS2	• PSTK															Kinexus	PK794	RpAb
RIOK2	S332+S335+	Q9BVS4	• PSTK															Kinexus	PK890	RpAb
RIPK1 (RIP; RIPK)	Pan	Q13546	• PSTK															No	sc-133102	MmAb
RIPK1 (RIP; RIPK)	S166	Q13546	• PSTK															Kinexus	PK945	RpAb
RIPK1 (RIP; RIPK)	S320	Q13546	• PSTK															Kinexus	PK946	RpAb
RIPK1 (RIP; RIPK)	Y384	Q13546	• PSTK															Kinexus	PK795	RpAb
RIPK2 (RICK; RIP2; CARD3)	S176	O43353	• PSTK															Kinexus	PK796	RpAb
RIPK2 (RICK; RIP2; CARD3)	Y381	O43353	• PSTK															Kinexus	PK797	RpAb
ROCK1 (ROKb)	T233	Q13464	• PSTK															Kinexus	PK947	RpAb
ROCK1 (ROKb)	Y913	Q13464	• PSTK															Kinexus	PK798	RpAb
ROCK2 (ROKa)	Pan	O75116	• PSTK															No	sc-1851	GpAb
ROCK2 (ROKa)	Pan	O75116	• PSTK															No	07-443	RpAb
ROCK2 (ROKa)	Y722	O75116	• PSTK															Kinexus	PK799	RpAb
Ron (RONa; MST1R)	Pan	Q04912	• PYK			●												Kinexus	NK161-3	RpAb
Ron (RONa; MST1R)	Pan	Q04912	• PYK			●												Kinexus	NK161-4	RpAb
Ron (RONa; MST1R)	Pan	Q04912	• PYK			●												Kinexus	NK161-2	RpAb
Ron (RONa; MST1R)	Y1238	Q04912	• PYK			●												Kinexus	PK800	RpAb
Ron (RONa; MST1R)	Y1238+Y123	Q04912	• PYK			●												Kinexus	PK801	RpAb
ROR2 (RON2)	Y645+Y646	Q01974	• PYK			●												Kinexus	PK802	RpAb
RORA	S35	P35398			●	●												Kinexus	PN651	RpAb
Ros (ROS1)	Pan	P08922	• PYK			●												Kinexus	NK163-3	RpAb
Ros (ROS1)	Pan	P08922	• PYK			●												Kinexus	NK163-4	RpAb
Ros (ROS1)	Pan	P08922	• PYK			●												Kinexus	NK163-2	RpAb
Ros (ROS1)	Y2114+Y211	P08922	• PYK			●												Kinexus	PK803	RpAb
RPS6	Pan	P62753						●										No	sc-74576	MmAb
RPS6	Pan	P62753						●										No	sc-74459	MmAb
RPS6	S235	P62753						●										No	07-433	RpAb
RPS6	S235+S236	P62753						●										Kinexus	PN685	RpAb
RPS6	S235+S236+	P62753						●										Kinexus	PN559	RpAb
RPTOR (RAPTOR)	T706	Q8N122											●					Kinexus	PN982	RpAb
RPTOR (RAPTOR)	S859+S863	Q8N122											●					Kinexus	PN981	RpAb
RSK1 (RPS6KA1; p90RSK)	Pan	Q15418	• PSTK															Kinexus	NK164-3	RpAb
RSK1 (RPS6KA1; p90RSK)	Pan	Q15418	• PSTK															No	sc-231	RpAb
RSK1 (RPS6KA1; p90RSK)	Y220+S221	Q15418	• PSTK															Kinexus	PK807	RpAb
RSK1 (RPS6KA1; p90RSK)	S221	Q15418	• PSTK															Kinexus	PK804	RpAb
RSK1 (RPS6KA1; p90RSK)	S380	Q15418	• PSTK															Kinexus	PK805	RpAb
RSK1 (RPS6KA1; p90RSK)	T573	Q15418	• PSTK															Kinexus	PK806	RpAb
RSK2 (RPS6KA3)	Pan	P51812	• PSTK															Kinexus	NK165-2	RpAb
RSK2 (RPS6KA3)	Pan	P51812	• PSTK															Kinexus	NK165-3	RpAb
RSK2 (RPS6KA3)	Pan	P51812	• PSTK															No	AF1518	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
RSK3 (RPS6KA2)	Y217+S218	Q15349	• PSTK															Kinexus	PK808	RpAb
RSK4 (RPS6KA6)	Pan	Q9UK32	• PSTK															Kinexus	NK166-2	RpAb
RTN4 (NOGO)	S107	Q9NQC3																Kinexus	PN983	RpAb
RUNX1 (AML1)	S276+Pan	Q01196			●													Kinexus	PN778	RpAb
RXRa	S260	P19793			●	●												Kinexus	PN652	RpAb
RYK	Pan	P34925	• PYK															No	AP7677b	RpAb
S100A9 (CAGB; MRP14)	Pan	P06702														●		Kinexus	NN459-3	RpAb
SAA2	Pan	P0DJ19									●							Kinexus	NN460-3	RpAb
SAC3 (FIG4)	Y631	Q92562								●								Kinexus	PN984	RpAb
SATB1	S38	Q01826			●													Kinexus	PN653	RpAb
SCN2B (NaVbeta2; Gm183)	Pan	O60939					●											No	NN314-1	MmAb
SCN3B (NaVbeta3)	Pan	Q9NY72					●											No	NN315-1	MmAb
SCNN1A (ENaCA; Alpha-	Pan	P37088					●											No	NN316-1	RpAb
SCYL1	S754	Q96KG9	• PSTK															Kinexus	PK809	RpAb
SETD2	S2080+S208	Q9BYW2																Kinexus	PN779	RpAb
SFC (KITLG)	Pan	P21583																No	10004254-M	MmAb
SGK1	Pan	O00141	• PSTK															Kinexus	NK294-1	RpAb
Sgk223	Y413	Q86YV5	• PSTK															Kinexus	PK810	RpAb
Sgk269 (PEAK1)	Y635	Q9H792	• PSTK															Kinexus	PK811	RpAb
Sgk288 (ANKK1)	Pan	Q8NFD2	• PSTK															Kinexus	NK295-1	RpAb
SGK3 (CISK; SGK1)	Pan	Q96BR1	• PSTK															Kinexus	NK170-4	RpAb
SGK3 (CISK; SGK1)	Pan	Q96BR1	• PSTK															Kinexus	NK170-3	RpAb
SH2BP1	T925	Q6PD62			●													Kinexus	PN654	RpAb
SHANK2 (CortBP1; Cortactin-	Pan	Q9UPX8											●					No	NN318-1	MmAb
SHANK3 (PROSAP2; PSAP2)	Pan	Q9BYB0											●					No	NN319-1	MmAb
Shc1 (Shc)	Y427	P29353											●					Kinexus	PN986	RpAb
Shc1 (Shc)	Y349+Y350	P29353											●					Kinexus	PN985	RpAb
Shc1 (Shc)	Y349+Y350	P29353											●					No	sc-23764	GpAb
Shc3	Y406	Q92529											●					Kinexus	PN987	RpAb
Shc4	Y424	Q6S5L8											●					Kinexus	PN988	RpAb
SHIP1 (INPP5D)	Pan	Q92835	• PI-P															Kinexus	NP044-2	RpAb
SHIP1 (INPP5D)	Pan	Q92835	• PI-P															Kinexus	NP044-3	RpAb
SHIP1 (INPP5D)	Pan	Q92835	• PI-P															Kinexus	NP044-1	RpAb
SHIP1 (INPP5D)	Y187	Q92835	• PI-P															Kinexus	PN560	RpAb
SHIP2 (INPPL1)	Pan	O15357	• PI-P															Kinexus	NP045-2	RpAb
SHIP2 (INPPL1)	Pan	O15357	• PI-P															Kinexus	NP045-3	RpAb
SHIP2 (INPPL1)	Pan	O15357	• PI-P															Kinexus	NP045-1	RpAb
SHIP2 (INPPL1)	Y886	O15357	• PI-P															Kinexus	PN534	RpAb
SIK (SNF1LK)	Pan	P57059	• PSTK															Kinexus	NK251-2	RpAb
SIK (SNF1LK)	Pan	P57059	• PSTK															Kinexus	NK251-3	RpAb
SIK (SNF1LK)	Pan	P57059	• PSTK															Kinexus	NK251-1	RpAb
SIK (SNF1LK)	T182	P57059	• PSTK															Kinexus	PK812	RpAb
SIK2 (QIK)	Pan	Q9H0K1	• PSTK															Kinexus	NK249-2	RpAb
SIK2 (QIK)	Pan	Q9H0K1	• PSTK															Kinexus	NK249-3	RpAb
SIK2 (QIK)	Pan	Q9H0K1	• PSTK															Kinexus	NK249-1	RpAb
SIK2 (QIK)	S358	Q9H0K1	• PSTK															Kinexus	PK813	RpAb
SIK3 (QSK)	Pan	Q9Y2K2	• PSTK															Kinexus	NK250-2	RpAb
SIK3 (QSK)	Pan	Q9Y2K2	• PSTK															Kinexus	NK250-3	RpAb
SIK3 (QSK)	Pan	Q9Y2K2	• PSTK															Kinexus	NK250-1	RpAb
SIK3 (QSK)	T163	Q9Y2K2	• PSTK															Kinexus	PK814	RpAb
SIK3 (QSK)	T411	Q9Y2K2	• PSTK															Kinexus	PK815	RpAb
SIN3A	S832+Pan	Q96ST3			●													Kinexus	PN655	RpAb
SIT	Y90	Q9Y3P8											●					Kinexus	PN535	RpAb
SIT	Y95	Q9Y3P8											●					Kinexus	PN536	RpAb
SLC12A1 (NKCC2;	Pan	Q13621														●		No	NN320-1	RpAb
SLC12A3 (NCC; Thiazide-	Pan	P55017														●		No	NN321-1	RpAb
SLC12A5 (KCC2; K-Cl	Pan	Q9H2X9														●		No	NN322-1	MmAb
SLC2A1 (GLUT-1)	T478	P11166														●		Kinexus	PN989	RpAb
SLC38A1 (ATA1; NAT2)	Pan	Q9H2H9														●		No	NN324-1	MmAb
SLC9A3 (NHE3)	Pan	P48764														●		No	NN326-1	RpAb
SLK	S189	Q9H2G2	• PSTK															Kinexus	PK816	RpAb
Smad1	S465	Q15797			●													No	11321	RpAb
Smad2	T220	Q15796			●													No	11323	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
Smad3	T8+Pan	P84022			●													Kinexus	PN658	RpAb
Smad4	Y513	Q13485			●													Kinexus	PN725	RpAb
SMARCA4	S610+S613	P51532			●													Kinexus	PN726	RpAb
SMARCB1	T134	Q12824			●													Kinexus	PN727	RpAb
SMG1	Pan	Q96Q15	• PYK															Kinexus	NK233-1	RpAb
SMG1	Pan	Q96Q15	• PYK															Kinexus	NK233-3	RpAb
SMG1	Pan	Q96Q15	• PYK															Kinexus	NK233-2	RpAb
SMG1	T3550	Q96Q15	• PYK															Kinexus	PK817	RpAb
SNAP25	S28+T29	P60880															●	Kinexus	PN1003	RpAb
SNAP25	T138	P60880															●	Kinexus	PN991	RpAb
SNAP25	S187	P60880															●	Kinexus	PN990	RpAb
SNCA (a-Synuclein)	Y39	P37840												●				Kinexus	PN850	RpAb
SNCA (a-Synuclein)	S129+Pan	P37840												●				Kinexus	PN849	RpAb
SNCA (a-Synuclein)	Y136	P37840												●				No	11286	RpAb
SND1	Y908	Q7KZF4			●													Kinexus	PN661	RpAb
snRNP70	Y126	P08621																Kinexus	PN537	RpAb
SOCS7	Y561	O14512			●													Kinexus	PN728	RpAb
SOD1 (hSod1)	Pan	P00441									●							No	NN098-4	RpAb
SOD1 (hSod1)	T40	P00441									●							Kinexus	PN905	RpAb
SOD1 (hSod1)	S99	P00441									●							Kinexus	PN904	RpAb
SOD1 (hSod1)	S108	P00441									●							Kinexus	PN903	RpAb
SOD2 (Mn-SOD)	Pan	P04179									●							No	NN068-1	RpAb
SOD2 (Mn-SOD)	Pan	P04179									●							No	NN068-2	RpAb
SOD3 (Cu/Zn; ECSOD)	Pan	P08294									●							No	NN328-1	MmAb
SOX9	S181+Pan	P48436			●													Kinexus	PN662	RpAb
SP4	S770	Q02446			●													Kinexus	PN851	RpAb
SPG11	S1955	Q96JI7																Kinexus	PN906	RpAb
SPG11	S1961	Q96JI7																Kinexus	PN907	RpAb
SPRED1	S105	Q7Z699																Kinexus	PN566	RpAb
SPT5	T791	O00267			●													Kinexus	PN663	RpAb
SQSTM1 (Sequestosome-1)	Pan	Q13501											●					No	NN375-1	RpAb
SQSTM1 (Sequestosome-1)	T269+S272	Q13501											●					Kinexus	PN909	RpAb
SQSTM1 (Sequestosome-1)	T339+S342	Q13501											●					Kinexus	PN910	RpAb
SQSTM1 (Sequestosome-1)	S403	Q13501											●					Kinexus	PN908	RpAb
Src	Pan	P12931	• PYK															No	sc-19	RpAb
Src	Pan	P12931	• PYK															No	KAP-TK121	RpAb
Src	Pan	P12931	• PYK															No	05-184	MmAb
Src	Y419	P12931	• PYK															Kinexus	PK818	RpAb
Src	Y419	P12931	• PYK															Kinexus	PK818-2	RpAb
Src	Y530	P12931	• PYK															Kinexus	PK948	RpAb
SRF	S224	P11831			●													Kinexus	PN664	RpAb
SRPK1	S222	Q96SB4	• PSTK															Kinexus	PK819	RpAb
SRPK1	S587	Q96SB4	• PSTK															Kinexus	PK891	RpAb
SRPK2	Pan	P78362	• PSTK															Kinexus	NK296-1	RpAb
SRPK2	Y319	P78362	• PSTK															Kinexus	PK820	RpAb
SSTR2	S341+S343	P30874				●												Kinexus	PN804	RpAb
STAG2	Y433	Q8N3U4																Kinexus	PN729	RpAb
STAM2	Y374	O75886											●					Kinexus	PN538	RpAb
STAT1	Pan	P42224			●													No	DB037	RpAb
STAT1	Y701	P42224			●													Kinexus	PN666	RpAb
STAT1	Y701	P42224			●													No	sc-8394	MmAb
STAT1	S727+Pan	P42224			●													Kinexus	PN667	RpAb
STAT2	Y690	P52630			●													Kinexus	PN668	RpAb
STAT3	Y705+T708	P40763			●													Kinexus	PN539	RpAb
STAT3	S727	P40763			●													Kinexus	PN669	RpAb
STAT4	Pan	Q14765			●													No	DB030	RpAb
STAT4	Y693	Q14765			●													Kinexus	PN670	RpAb
STAT5A	Y694	P42229			●													Kinexus	PN671	RpAb
STAT5A	S780	P42229			●													Kinexus	PN672	RpAb
STAT5B	Pan	P51692			●													No	DB032	RpAb
STAT6	Pan	P42226			●													No	DB038	RpAb
STAT6	Y641	P42226			●													Kinexus	PN673	RpAb
STI1 (STIP1)	Pan	P31948									●							No	S65720	MmAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
STLK3 (DCHT)	T231	Q9UEW8	• PSTK															Kinexus	PK916	RpAb
STLK6 (ALS2CR2)	S304+S306	Q9C0K7	• PSTK															Kinexus	PN911	RpAb
STX1B	S14	P61266															●	Kinexus	PN852	RpAb
STXBP1 (HUNC18)	S313	P61764															●	Kinexus	PN992	RpAb
STXBP1 (HUNC18)	Y473	P61764															●	Kinexus	PN994	RpAb
STXBP1 (HUNC18)	T574	P61764															●	Kinexus	PN993	RpAb
SUR1 (Abcc8)	Pan	Q09428					●											No	NN340-1	MmAb
Syk	Pan	P43405	• PYK															No	sc-1240	MmAb
Syk	Y323	P43405	• PYK															Kinexus	PK821	RpAb
Syk	Y352	P43405	• PYK															Kinexus	PK822	RpAb
Syk	Y525+Y526	P43405	• PYK															Kinexus	PK823	RpAb
SYN1 (Synapsin 1)	Pan	P17600															●	No	Y021259	RpAb
SYN1 (Synapsin 1)	S605	P17600															●	Kinexus	PN854	RpAb
SYN1 (Synapsin 1)	S551+S553	P17600															●	Kinexus	PN853	RpAb
Synaptophysin (SYP)	Pan	P08247															●	No	NN370-1	MmAb
SYNJ1 (INPP5G)	Y784+Y786	O43426															●	Kinexus	PN996	RpAb
SYNJ1 (INPP5G)	S1084	O43426															●	Kinexus	PN995	RpAb
SYT1 (SVP65)	T129	P21579															●	Kinexus	PN997	RpAb
SYT1 (SVP65)	Y312	P21579															●	Kinexus	PN998	RpAb
SYT1 (SVP65)	Y381	P21579															●	Kinexus	PN999	RpAb
SYT3	Pan	Q9BQG1															●	No	NN366-1	MmAb
SYT6	Pan	Q5T7P8															●	No	NN367-1	MmAb
SYT9 (Synaptotagmin 9)	Pan	Q86SS6															●	No	NN343-1	MmAb
SYT10 (Synaptotagmin 10)	Pan	Q6XYQ8															●	No	NN344-1	MmAb
SYT12 (Synaptotagmin 12)	Pan	Q8IV01															●	No	NN345-1	MmAb
TAK1 (MAP3K7)	Pan	O43318	• PSTK															Kinexus	NK175-3	RpAb
TAK1 (MAP3K7)	Pan	O43318	• PSTK															No	sc-7967	MmAb
TAK1 (MAP3K7)	Pan	O43318	• PSTK															No	AP-ST0091	RpAb
TAK1 (MAP3K7)	T184+T187	O43318	• PSTK															Kinexus	PK825	RpAb
TAK1 (MAP3K7)	S439	O43318	• PSTK															Kinexus	PK824	RpAb
TAO1 (TAOK1)	S181	Q7L7X3	• PSTK															Kinexus	PK826	RpAb
TAO1 (TAOK1)	Y309	Q7L7X3	• PSTK															Kinexus	PK827	RpAb
TAO2 (TAOK2)	Pan	Q9UL54	• PSTK															No	sc-514268	MmAb
TAO3 (JIK, TAOK3)	Pan	Q9H2K8	• PSTK															No	AP7902a	RpAb
TAO3 (JIK, TAOK3)	Pan	Q9H2K8	• PSTK															Kinexus	NK087-2	RpAb
TARDBP	S409+S410	Q13148			●													Kinexus	PN674	RpAb
Tau (MAPT)	S202	P10636												●				No	28017	RpAb
Tau (MAPT)	T498	P10636												●				Kinexus	PN855	RpAb
Tau (MAPT)	S515+S516+	P10636												●				Kinexus	PN856	RpAb
Tau (MAPT)	T522 (T205)	P10636												●				No	Y021108	RpAb
Tau (MAPT)	T529+T534	P10636												●				Kinexus	PN857	RpAb
Tau (MAPT)	S713+Pan	P10636												●				Kinexus	PN858	RpAb
TBC1D7	Y14	Q9P0N9																Kinexus	PN540	RpAb
TBK1 (IKKd)	Pan	Q9UHD2	• PSTK															No	sc-9085	RpAb
TBK1 (IKKd)	S172	Q9UHD2	• PSTK															Kinexus	PK828	RpAb
TBK1 (IKKd)	S716	Q9UHD2	• PSTK															Kinexus	PK949	RpAb
TBXA2R	S329+S331	P21731				●												Kinexus	PN805	RpAb
TDP-43 (TADBP)	S48	Q13148																Kinexus	PN914	RpAb
TDP-43 (TADBP)	S183	Q13148																Kinexus	PN912	RpAb
TDP-43 (TADBP)	S403+S404	Q13148																Kinexus	PN913	RpAb
Tec	Y519	P42680	• PYK															Kinexus	PK829	RpAb
TERF1	T371	P54274			●													Kinexus	PN675	RpAb
TGFβ1	Pan	P01137																No	sc-146	RpAb
TGM2	Y369	P21980																Kinexus	PN541	RpAb
TH (TY3H)	S19	P07101					●											Kinexus	PN731	RpAb
TH (TY3H)	S71	P07101					●											Kinexus	PN730	RpAb
THRAP3	S253	Q9Y2W1			●													Kinexus	PN676	RpAb
TIE2 (TEK)	Pan	Q02763	• PYK			●												No	AP7684a	RpAb
TIE2 (TEK)	Y897	Q02763	• PYK			●												Kinexus	PK830	RpAb
TIE2 (TEK)	Y992	Q02763	• PYK			●												Kinexus	PK831	RpAb
TLK1	Pan	Q9UKI8	• PSTK															No	AP8101b	RpAb
TLN1	Y70	Q9Y490												●				Kinexus	PN542	RpAb
TLR4 (CD284)	Pan	O00206				●												No	NN368-1	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
TMEM205	Pan	Q6UW68																No	sc-514568	MmAb
TNFR1 (CD120a; TNFRSF1A;	Pan	P19438				●												No	NN347-1	RpAb
TNK1	Y277	Q13470	• PYK															Kinexus	PK832	RpAb
TORC2 (CRTC2)	S433	Q53ET0			●													Kinexus	PN677	RpAb
TP53 (p53)	Pan	P04637			●													No	sc-126	MmAb
TP53 (p53)	Pan	P04637			●													No	NN082-2	MmAb
TP53 (p53)	Pan	P04637			●													No	NN082-3	RpAb
TP53 (p53)	S6	P04637			●													No	11092	RpAb
TP53 (p53)	S6+S9	P04637			●													Kinexus	PN637	RpAb
TP53 (p53)	S15	P04637			●													Kinexus	PN859	RpAb
TP53 (p53)	T18+S20	P04637			●													Kinexus	PN638	RpAb
TP53 (p53)	S392	P04637			●													Kinexus	PN640	RpAb
TP53BP1	T1056	Q12888			●													Kinexus	PN678	RpAb
TPH1	T225+S228	P17752								●								Kinexus	PN780	RpAb
TPPb	Pan	NA								●								No	B50263-51	RpAb
TRADD	Pan	Q15628											●					No	T50320	MmAb
Trail (TNSF10)	Pan	P50591																No	DB040	RpAb
TRAK2 (ALS2CR3)	Pan	O60296																No	SMC-483D	MmAb
TRAP1 (HSP75)	Pan	Q12931									●							No	NN349-2	RpAb
TRAP1 (HSP75)	Pan	Q12931									●							No	NN349-1	MmAb
TRHR	S364+T365	P34981				●												Kinexus	PN806	RpAb
TRIM28 (TIF1B)	Y458	Q13263	• PSTK		●													Kinexus	PK834	RpAb
TRIM28 (TIF1B)	S473	Q13263	• PSTK		●													Kinexus	PK833	RpAb
TRIM28 (TIF1B)	Y517	Q13263	• PSTK		●													Kinexus	PK835	RpAb
TRIM33 (TIF1G)	S1119	Q9UPN9	• PSTK		●													Kinexus	PK836	RpAb
TrkA (NGFR; NTRK1)	Y496	P04629	• PYK			●												No	sc-8058	MmAb
TrkA (NGFR; NTRK1)	Y680+Y681	P04629	• PYK			●												Kinexus	PK837	RpAb
TrkB (NTRK2)	Y516	Q16620	• PYK			●												Kinexus	PK838	RpAb
TrkB (NTRK2)	Y702	Q16620	• PYK			●												Kinexus	PK839	RpAb
TrkB (NTRK2)	Y706+Y707	Q16620	• PYK			●												Kinexus	PK917	RpAb
TrkC (NTRK3)	Y709+Y710	Q16288	• PYK			●												Kinexus	PK840	RpAb
TRPV4	Pan	Q9HBA0					●											Kinexus	NN464-2	RpAb
TRRAP	Pan	Q9Y4A5	• PSTK															Kinexus	NK232-2	RpAb
TRRAP	Pan	Q9Y4A5	• PSTK															Kinexus	NK232	RpAb
TRRAP	Pan	Q9Y4A5	• PSTK															Kinexus	NK232-3	RpAb
TSC1	S505	Q92574											●					Kinexus	PN781	RpAb
TSSK3	T168	Q96PN8	• PSTK															Kinexus	PK841	RpAb
TTBK1	Pan	Q5TCY1	• PSTK															Kinexus	NK306-1	RpAb
TTBK1	Pan	Q5TCY1	• PSTK															Kinexus	NK306-2	RpAb
TTBK1	T186	Q5TCY1	• PSTK															Kinexus	PK955	RpAb
TTBK2	Pan	Q6IQ55	• PSTK															Kinexus	NK307-2	RpAb
TTK (MPS1)	Pan	P33981	• DSK															No	AP8103a	RpAb
TTK (MPS1)	S677	P33981	• DSK															Kinexus	PK842	RpAb
TTK (MPS1)	Y833+Y836	P33981	• DSK															Kinexus	PK843	RpAb
TUBA4A	Pan	P68366												●				No	NN380-1	RpAb
TUBA4A	Y103	P68366												●				Kinexus	PN915	RpAb
TUBA4A	Y210	P68366												●				Kinexus	PN916	RpAb
TXK	Y420	P42681	• PYK															Kinexus	PK844	RpAb
TYK2	Pan	P29597	• PYK															Kinexus	NK181-3	RpAb
TYK2	Pan	P29597	• PYK															Kinexus	NK181-4	RpAb
TYK2	Pan	P29597	• PYK															Kinexus	NK181-5	RpAb
TYK2	Pan	P29597	• PYK															No	sc-169	RpAb
TYK2	Pan	P29597	• PYK															No	sc-7205	RpAb
TYK2	Y292	P29597	• PYK															Kinexus	PK846	RpAb
TYK2	Y1054+Y105	P29597	• PYK															Kinexus	PK845	RpAb
Tyro3	Y681	Q06418	• PYK			●												Kinexus	PK847	RpAb
Tyro3	Y685+Y686	Q06418	• PYK			●												Kinexus	PK848	RpAb
UBF	T201	P17480			●													Kinexus	PN680	RpAb
UBF	S484	P17480			●													Kinexus	PN679	RpAb
Ubiquitin (Ub; UBB)	Pan	NA							●									No	sc-8017	MmAb
UBQLN2	T128	Q9UHD9							●									Kinexus	PN920	RpAb
UBQLN2	S222	Q9UHD9							●									Kinexus	PN918	RpAb
UBQLN2	S256	Q9UHD9							●									Kinexus	PN919	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
UGDH	Y352	O60701								●								Kinexus	PN782	RpAb
ULK1	Pan	O75385	• PSTK															Kinexus	NK298-1	RpAb
ULK1	Pan	O75385	• PSTK															No	NK298-2	RpAb
ULK2	Pan	Q8IYT8	• PSTK															No	NK354-1	RpAb
ULK3	Pan	Q6PHR2	• PSTK															No	NK355-1	RpAb
ULK4	Pan	Q96C45	• PSTK															Kinexus	NK299-1	RpAb
UTRN	Y3197	P46939												●				Kinexus	PN1000	RpAb
UVRAG	Pan	Q9P2Y5															●	No	NN356-1	RpAb
VACAMKL	Y245	Q8NCB2	• PSTK															Kinexus	PK892	RpAb
VAMP1	S63	P23763															●	Kinexus	PN1001	RpAb
VAMP4	S30	O75379															●	Kinexus	PN1002	RpAb
VAPB	T46	O95292															●	Kinexus	PN924	RpAb
VAPB	T150	O95292															●	Kinexus	PN922	RpAb
VAPB	T164	O95292															●	Kinexus	PN923	RpAb
VAPB	S206	O95292															●	Kinexus	PN921	RpAb
VAV1 (Vav)	Y826	P15498																Kinexus	PN543	RpAb
VCP	S37	P55072																Kinexus	PN925	RpAb
VCP	T56	P55072																Kinexus	PN929	RpAb
VCP	T316	P55072																Kinexus	PN927	RpAb
VCP	T509	P55072																Kinexus	PN928	RpAb
VCP	S702	P55072																Kinexus	PN926	RpAb
VCP	T761	P55072																Kinexus	PN930	RpAb
VEGFR1 (Flt1)	Pan	P17948	• PYK															No	AP7642b	RpAb
VEGFR1 (Flt1)	Pan	P17948	• PYK															Kinexus	NK226-2	RpAb
VEGFR1 (Flt1)	Pan	P17948	• PYK															Kinexus	NK226-1	RpAb
VEGFR1 (Flt1)	Pan	P17948	• PYK															No	AP7642a	RpAb
VEGFR1 (Flt1)	Y1048	P17948	• PYK															Kinexus	PK850	RpAb
VEGFR1 (Flt1)	Y1053	P17948	• PYK															Kinexus	PK851	RpAb
VEGFR2 (KDR)	Pan	P35968	• PYK															Kinexus	NK245-2	RpAb
VEGFR2 (KDR)	Pan	P35968	• PYK															Kinexus	NK245-3	RpAb
VEGFR2 (KDR)	Pan	P35968	• PYK															Kinexus	NK245-1	RpAb
VEGFR2 (KDR)	Y1054	P35968	• PYK															Kinexus	PK852	RpAb
VEGFR3 (Flt4)	Pan	P35916	• PYK															Kinexus	NK064-3	RpAb
VEGFR3 (Flt4)	Pan	P35916	• PYK															Kinexus	NK064-2	RpAb
VEGFR3 (Flt4)	Y1063	P35916	• PYK															No	11531	RpAb
VEGFR3 (Flt4)	Y1068	P35916	• PYK															Kinexus	PK853	RpAb
Versican (GHAP; CSPG2)	Pan	P13611														●		No	NN360-1	MmAb
VGLUT1 (SLC17A7; BNPI)	Pan	Q9P2U7														●		No	NN361-1	MmAb
VGLUT3 (SLC17A8; Solute)	Pan	Q8NDX2														●		No	NN362-1	MmAb
VHL	S111	P40337							●									Kinexus	PN783	RpAb
VIM (Vimentin)	Y117	P08670												●				Kinexus	PN544	RpAb
Vrk1	Pan	Q99986	• PSTK															No	AP7408c	RpAb
WASP	Y291	P42768											●					Kinexus	PN545	RpAb
Wee1	Pan	P30291	• DSK															Kinexus	NK185	RpAb
Wee1	S642	P30291	• DSK															Kinexus	PK854	RpAb
Wip1 (PPM1D)	Pan	O15297	• PSTP															No	sc-7263	GpAb
Wip1 (PPM1D)	Pan	O15297	• PSTP															No	sc-7264	GpAb
WNK1	Pan	Q9H4A3	• PSTK															Kinexus	NK252-2	RpAb
WNK1	Pan	Q9H4A3	• PSTK															Kinexus	NK252-3	RpAb
WNK1	Pan	Q9H4A3	• PSTK															Kinexus	NK252-1	RpAb
WNK1	T60	Q9H4A3	• PSTK															Kinexus	PK856	RpAb
WNK1	S382+Pan	Q9H4A3	• PSTK															Kinexus	PK855	RpAb
WNK1	T2245	Q9H4A3	• PSTK															Kinexus	PK857	RpAb
WNK2	Pan	Q9Y3S1	• PSTK															Kinexus	NK253-2	RpAb
WNK2	Pan	Q9Y3S1	• PSTK															Kinexus	NK253-3	RpAb
WNK2	Pan	Q9Y3S1	• PSTK															Kinexus	NK253-1	RpAb
WNK3 (PRKWNK3)	Pan	Q9BYP7	• PSTK															Kinexus	NK254-2	RpAb
WNK3 (PRKWNK3)	Pan	Q9BYP7	• PSTK															Kinexus	NK254-3	RpAb
WNK3 (PRKWNK3)	Pan	Q9BYP7	• PSTK															Kinexus	NK254-1	RpAb
WNK4 (PRKWNK4)	Pan	Q96J92	• PSTK															Kinexus	NK255-3	RpAb
WNK4 (PRKWNK4)	Pan	Q96J92	• PSTK															Kinexus	NK255-1	RpAb
WT1	S365	P19544			●													Kinexus	PN732	RpAb
YAP1	S109	P46937			●													Kinexus	PN681	RpAb

Target Protein Name	Pan-specific or P-Site	Human UniProt ID	Protein Kinase	Protein Phosphatase	Transcription	Receptor	Ion Channel	Protein Synthesis	Protein Degradation	Metabolic Enzyme	Stress Protein	Apoptosis	Adapter/ Scaffold	Cytoskeletal	Cell Adhesion	Ca/ Lipid / Solute	Vesicle Protein	Kinexus Anti-body	Kinexus Antibody Code	Anti-body Type
YAP1	T119	P46937			●													Kinexus	PN682	RpAb
YAP1	S127	P46937			●													Kinexus	PN683	RpAb
YDJ1 (MAS5; Yeast dnaJ	Pan	P25491																No	NN057-6	MmAb
Yes	Pan	P07947	• PYK															No	sc-8403	MmAb
Yes	Pan	P07947	• PYK															No	sc-46674	MmAb
Yes	Pan	P07947	• PYK															No	sc-14	RpAb
Yes	Y222+Y223	P07947	• PYK															Kinexus	PK858	RpAb
YLPM1	S561	P49750			●													Kinexus	PN684	RpAb
YSK1 (STK25; SOK1)	Pan	O00506	• PSTK															No	sc-6865	GpAb
YSK1 (STK25; SOK1)	Pan	O00506	• PSTK															No	sc-6884	GpAb
YSK1 (STK25; SOK1)	T174	O00506	• PSTK															Kinexus	PK859	RpAb
YSK4 (MAP3K19)	Pan	Q56UN5	• PSTK															Kinexus	NK256-2	RpAb
YSK4 (MAP3K19)	Pan	Q56UN5	• PSTK															Kinexus	NK256-3	RpAb
YSK4 (MAP3K19)	Pan	Q56UN5	• PSTK															Kinexus	NK256-1	RpAb
ZAP70	Pan	P43403	• PYK															No	sc-1526	GpAb
ZAP70	Pan	P43403	• PYK															No	sc-574	RpAb
ZAP70	Y248	P43403	• PYK															Kinexus	PK860	RpAb
ZAP70	Y292	P43403	• PYK															Kinexus	PK861	RpAb
ZAP70	Y319	P43403	• PYK															Kinexus	PK862	RpAb
ZAP70	Y492+Y493	P43403	• PYK															Kinexus	PK863	RpAb
ZC2 (TNIK; GK)	Pan	Q9UKE5	• PSTK															Kinexus	NK301-1	RpAb