

Sequence Alignment of Activating Phosphorylation Sites in Human MAPK

Erk1 DFGLARVADPDHDHTGFLTE[•]Y[•]VATRWRAPPE
 Erk2 DFGLARVADPDHDHTGFLTE[•]Y[•]VATRWRAPPE
 Erk3 DFGLARIMDPHYSHKGHLSEGLVTKWYRSPE
 Jnk1 DFGLARTACTNFM-----MTPYVVTRYRRAPE
 p38 DFGLARHTDDE-----MTGYVATRWRAPPE

