



Superagonist TCR/MHC complex [pdb 1G6R]: T-cell receptor alpha (orange) and beta (yellow) extracellular chains; Major histocompatibility complex Class 1 molecule (green); Beta-2 microglobulin chain (red); & antigen peptide Siyr (blue).