

KinSight™ Phosphatase Selectivity Profiling – Target List

Powered by the PhosphoSens® Platform



HGNC (Common) Name	HGNC Approved Full Name	Ser (S) Thr (T) Tyr (Y)	Sensor Peptide Substrate
CDC25A	cell division cycle 25A	Dual (Y)	AQT0523
CDC25B	cell division cycle 25B	Dual (Y)	AQT0523
CDC25C	cell division cycle 25C	Dual (Y)	AQT0523
DUSP22 (JSP-1, LMW-DSP2, MKPX)	dual specificity phosphatase 22	Dual	AQT0737
DUSP3 (VHR)	dual specificity phosphatase 3	Dual (Y)	AQT0264
PPM1A (PP2Ca, PP2CA)	protein phosphatase, Mg2+/Mn2+ dependent 1A	S/T	AQT0267
PPM1D (WIP1, PP2C-delta)	protein phosphatase, Mg2+/Mn2+ dependent 1D	S/T	AQT0525
PPM1G (PP2Cg)	protein phosphatase, Mg2+/Mn2+ dependent 1G	S/T	AQT0267
PPM1F (CamKPase)	protein phosphatase, Mg2+/Mn2+ dependent 1F	S/T	AQT0523
PPP1CA (PP1A/PPP1A)	protein phosphatase 1 catalytic subunit alpha	S/T	AQT0480
PPP1CB (PP1B, PPP1B)	protein phosphatase 1 catalytic subunit beta	S/T	AQT0523
PPP2CA (PP2A α, RP-C)	protein phosphatase 2 catalytic subunit alpha	S/T	AQT0523
PPP2CB (PP2Abeta)	protein phosphatase 2 catalytic subunit beta	S/T	AQT0523
PPP3CA/B (Calcineurin A/B, CNA alpha/beta, CALNA/B)	protein phosphatase 3 catalytic subunit alpha/beta	S/T	AQT0670
PTPN1 (PTP1B, PTB1B)	protein tyrosine phosphatase non-receptor type 1	Y	AQT0266
PTPN2 (TCPTP, PTPT)	protein tyrosine phosphatase non-receptor type 2	Y	AQT0264
PTPN4 (MEG1)	protein tyrosine phosphatase non-receptor type 4	Y	AQT0737
PTPN5 (STEP)	protein tyrosine phosphatase non-receptor type 5	Y	AQT0664
PTPN6 (SHP 1, PTP-1C)	protein tyrosine phosphatase non-receptor type 6	Y	AQT0737
PTPN7 (HEPTP, LC-PTP)	protein tyrosine phosphatase non-receptor type 7	Y	AQT0772
PTPN9 (MEG2)	protein tyrosine phosphatase non-receptor type 9	Y	AQT0737
PTPN11 (SHP 2, PTP2C, PTP1D)	protein tyrosine phosphatase non-receptor type 11	Y	AQT0737
PTPN12 (PTP-PEST, PTPG1)	protein tyrosine phosphatase non-receptor type 12	Y	AQT0737
PTPN13 (FAP-1, PTP-BAS, PTP-E1, PNP1)	protein tyrosine phosphatase non-receptor type 13	Y	AQT0737
PTPN22 (PTPN8, Lyp)	protein tyrosine phosphatase non-receptor type 22	Y	AQT0737
PTPRA (R-PTP-alpha, PTPA, LRP)	protein tyrosine phosphatase receptor type A	Y	AQT0667
PTPRB (R-PTP-beta, VE-PTP)	protein tyrosine phosphatase receptor type B	Y	AQT0667
PTPRC (L-CA, CD45, T200)	protein tyrosine phosphatase receptor type C	Y	AQT0630
PTPRE (R-PTP-epsilon)	protein tyrosine phosphatase receptor type E	Y	AQT0667
PTPRF (LAR)	protein tyrosine phosphatase receptor type F	Y	AQT0737
PTPRG (R-PTP-gamma)	protein tyrosine phosphatase receptor type G	Y	AQT0667
PTPRJ (R-PTP-eta, DEP1, CD148)	protein tyrosine phosphatase receptor type J	Y	AQT0737
PTPRM (PTPμ, PTPRL1, R-PTP-mu)	protein tyrosine phosphatase receptor type M	Y	AQT0737
PTPRS (R-PTP-sigma)	protein tyrosine phosphatase receptor type S	Y	AQT0737