

PhosphoSens Sensor Peptides for Continuous Phosphatase Assays

A Powerful Approach for Understanding Phosphatase and Phosphatase Inhibitor Activity

Continuous • Direct • Catalytic

Search below to match your PHOSPHATASE with our SENSOR PEPTIDE substrate!

Bulk Sensor Peptide: Designed for higher throughput applications. 1mg is sufficient for ~1,200 assays in a low-volume 384-well plate (10µM sensor peptide in 25µL reaction volume). Available from 1mg to +200mg aliquots. Optimized ancillary reagents sold separately.

Assay Kits: Designed for streamlined technology evaluation. Each kit enables 96 reactions in a low-volume 384-well plate (10µM sensor peptide in 25µL reaction volume). Includes all assay components (excluding your kinase).

See your target of interest? Contact us here. Or email us at hello@assayquant.com!

Don't see your target of interest? Contact us. Our PhosphoSens Technology enables rapid assay development.

HGNC Common Name	HGNC Approved Full Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Class or Family	Enzyme used in Sensor Peptide Development	Assay Strength*
PHLPP1 (PHLPP, SCOP)	PH domain Leucine-Rich Repeat-Containing Protein Phosphatase 1	CSKS-AQT0481B CSKS-AQT0481K	S/T	Soluble	Academic Lab	M
PHLPP2 (PHLPL)	PH domain Leucine-Rich Repeat-Containing Protein Phosphatase 2	CSKS-AQT0481B CSKS-AQT0481K	S/T	Soluble	Academic Lab	M
PPP1CA (PP1A, PPP1A)	Serine/threonine-protein phosphatase PP1-alpha catalytic subunit	CSKS-AQT0480B CSKS-AQT0480K	S/T	Soluble	BPS (41106)	M
PPM1A (PP2Calpha, PP2CA)	Protein Phosphatase 1A	CSKS-AQT0267B CSKS-AQT0267K	S/T	Soluble	SignalChem (P02-20G)	VS
PPM1G (PP2Cgamma)	Protein Phosphatase 1G	CSKS-AQT0267B CSKS-AQT0267K	S/T	Soluble	SignalChem (P07-20G)	S
PPP1CB (PP1B, PPP1B)	Serine/threonine-protein phosphatase PP1-alpha catalytic subunit	CSKS-AQT0523B CSKS-AQT0523K	S/T	Soluble	SignalChem (P14-20G)	VS
PPP2CA (PP2Aalpha/PP2AR, RP-C)	Serine/threonine-protein phosphatase 2A catalytic subunit alpha isoform	CSKS-AQT0523B CSKS-AQT0523K	S/T	Soluble	BPS (30056)	VS
PPP3CA/B (Calcineurin A/B, CNA alpha/beta, CALNA/B)	Protein Phosphatase 3 catalytic subunit alpha/beta	CSKS-AQT0670B CSKS-AQT0670K	S/T	Soluble	R&D Systems (3160-CA20)	S
PPM1D (WIP1, PP2C-delta)	Protein phosphatase 1D	CSKS-AQT0525B CSKS-AQT0525K	S/T	Soluble	SignalChem (P04-20G)	S
CDC25A	M-phase inducer phosphatase 1	CSKS-AQT0523B CSKS-AQT0523K	Dual (Y)	Soluble	SignalChem (C04-20G)	W
CDC25B	M-phase inducer phosphatase 2	CSKS-AQT0523B CSKS-AQT0523K	Dual (Y)	Soluble	SignalChem (C04-20BG)	W
CDC25C	M-phase inducer phosphatase 3	CSKS-AQT0523B CSKS-AQT0523K	Dual (Y)	Soluble	SignalChem (C04-20CG)	W
DUSP3 (VHR)	Dual specificity protein phosphatase 3	CSKS-AQT0264B CSKS-AQT0264K	Dual (Y)	Soluble	R&D Systems (1654-VH-050)	M
DUSP22 (JSP-1, LMW-DSP2, MKPX)	Dual specificity protein phosphatase 22	CSKS-AQT0737B CSKS-AQT0737K	Dual	Soluble	SignalChem (D22-20G)	S
PTPN1 (PTP1B, PTB1B)	Tyrosine-protein phosphatase non-receptor type 1	CSKS-AQT0266B CSKS-AQT0266K	Y	Soluble	BPS (30010)	VS
PTPN2 (TCPTP, PTPT)	Tyrosine-protein phosphatase non-receptor type 2	CSKS-AQT0264B CSKS-AQT0264K	Y	Soluble	BPS (30013)	VS
PTPN4 (MEG1)	Tyrosine-protein phosphatase non-receptor type 4	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	BPS (30015)	VS
PTPN5 (STEP)	Tyrosine-protein phosphatase non-receptor type 5	CSKS-AQT0664B CSKS-AQT0664K	Y	Soluble	Academic Lab	M
PTPN6 (SHP-1, PTP-1C)	Tyrosine-protein phosphatase non-receptor type 6	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	BPS (100342)	VS
PTPN7 (HEPTP, LC-PTP)	Tyrosine-protein phosphatase non-receptor type 7	CSKS-AQT0772B CSKS-AQT0772K	Y	Soluble	SignalChem (P34-20G)	VS

HGNC Common Name	HGNC Approved Full Name	Product Number (Bulk, Kit)	Ser (S) Thr (T) Tyr (Y)	Class or Family	Enzyme used in Sensor Peptide Development	Assay Strength*
PTPN9 (MEG2)	Tyrosine-protein phosphatase non-receptor type 9	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	BPS (30020)	VS
PTPN11 (SHP-2, PTP2C, PTP1D)	Tyrosine-protein phosphatase non-receptor type 11	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	BPS (30022)	VS
PTPN12 (PTP-PEST, PTPG1)	Tyrosine-protein phosphatase non-receptor type 12	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	BPS (30023)	VS
PTPN13 (FAP-1, PTP-BAS, PTP-E1, PNP1)	Tyrosine-protein phosphatase non-receptor type 13	CSKS-AQT0737B CSKS-AQT0737K	Y	Soluble	SignalChem (P40-21G)	VS
PTPRA (R-PTP-alpha, PTPA, LRP)	Receptor-type tyrosine-protein phosphatase alpha	CSKS-AQT0667B CSKS-AQT0667K	Y	Receptor	SignalChem (P48-21G)	VS
PTPRB (R-PTP-beta, VE-PTP)	Receptor-type tyrosine-protein phosphatase beta	CSKS-AQT0667B CSKS-AQT0667K	Y	Receptor	SignalChem (P49-21G)	M
PTPRC (L-CA, CD45, T200)	Receptor-type tyrosine-protein phosphatase C	CSKS-AQT0630B CSKS-AQT0630K	Y	Receptor	BPS (30044)	VS
PTPRE (R-PTP-epsilon)	Receptor-type tyrosine-protein phosphatase epsilon	CSKS-AQT0667B CSKS-AQT0667K	Y	Receptor	SignalChem (P52-21G)	S
PTPRF (LAR)	Receptor-type tyrosine-protein phosphatase F	CSKS-AQT0737B CSKS-AQT0737K	Y	Receptor	SignalChem (P53-21G)	VS
PTPRJ (R-PTP-eta, DEP1, CD148)	Receptor-type tyrosine-protein phosphatase eta	CSKS-AQT0737B CSKS-AQT0737K	Y	Receptor	SignalChem (P56-21G)	VS
PTPRS (R-PTP-sigma)	Receptor-type tyrosine-protein phosphatase S	CSKS-AQT0737B CSKS-AQT0737K	Y	Receptor	SignalChem (P64-21G)	VS

Summary of PhosphoSens Kinase Target Assay Strength Based on Reaction Rates*

Categories	Total	Very Strong (VS)	Strong (S)	Medium (M)	Weak (W)
Initial Rate (RFU/pmole/min)		>1000	300-999	100-299	30-99
# Assays by Category	31	17	5	6	3
% Assays by Category		55%	16%	19%	10%
% Assays Very Strong or Strong		71%		Being Improved	