

Technical Data Sheet

Product Name	<i>Kinase Substrates Library, Group II, biotinylated, 18 distinct peptide mixtures</i>
Size	1 Set
Catalog #	AS-62335
Purity	Unpurified
Description	<p>This positional scanning synthetic peptide combinatorial library (PS-SPCL) has been used in rapid identification of serine/threonine kinase phosphorylation motif. The use of this Group II of distinct 18 peptide mixtures, in combination with Group I of 180 distinct peptide mixtures (cat# AS-62017-1), can provide sequence preference information of different kinases. Amount provided is 1 mg of peptide mixture x 18 vials. Sequences: Y-A-pT-X-X-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-pT-X-X-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-pT-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-pT-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-pT-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-pT-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-X-pT-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-X-X-X-pT-A-G-K-K(LC-biotin)-NH₂ Y-A-pY-X-X-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-pY-X-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-pY-X-X-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-pY-S/T-X-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-pY-X-X-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-X-pY-X-A-G-K-K(LC-biotin)-NH₂ Y-A-X-X-X-X-X-S/T-X-X-X-pY-A-G-K-K(LC-biotin)-NH₂ X = degenerate mixture of the 17 natural amino acids excluding cysteine, serine, and threonine. LC = 'long chain' version with an additional aminohexanoic acid spacer between the biotin and lysine side chain. S/T = equimolar mixture of serine and threonine. pT is phosphothreonine, and pY is phosphotyrosine.</p>
Storage	-20°C
References	Turk, BB. et al. <i>Nature Protocols</i> 1 , 375 (2006); Hutti, JE. et al. <i>Nature Methods</i> 1 , 27 (2004); Pinilla, C. et al. <i>Biotechniques</i> 13 , 901 (1992).