

PP1/PP2A Substrate

Synthetic peptide substrate derived from human GSK3 β

Catalog # P50-58

Lot # F480-5

Product Description

The PP1/PP2A Substrate peptide [GRPRTS(p)SFAEG] is derived from human GSK3 β (glycogen synthase kinase-3 beta) (amino acids 3-13) and is suitable as PP1 and PP2A substrate.

Molecular Weight

The theoretical molecular weight is 1,242.6.

Purity

The purity was determined to be 95.42% by HPLC analysis.

Formulation

1mg of peptide supplied as a lyophilized powder.

Storage and Stability

Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Reconstitution Protocol

Dilute peptide in 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

PP1/PP2A Substrate

Synthetic peptide substrate derived from human GSK3 β

Catalog #	P50-58
Quantity	1 mg
Lot #	F480-5
Purity	95.42%
Format	1 mg lyophilized powder
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C . For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.