

S6K Substrate

Synthetic peptide substrate derived from human 40S ribosomal protein S6

Catalog # S05-58

Lot # K175-2

Product Description

The S6K Substrate peptide sequence (KRRRLASLR) is based on human 40S ribosomal protein S6 (amino acid 230-238).

Molecular Weight

The theoretical molecular weight is 1155.4

Purity

The purity was determined to be 99.1% 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 distilled H₂O to a final concentration of 1mg/ml.

Related Kinases

S6K Substrate can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
p70S6K, Active	R21-10H
RSK1, Active	R15-10G
RSK3, Active	R16-10G
PIM1, Active	P35-10G
PIM2, Active	P36-10G

S6K Substrate

Synthetic peptide substrate derived from human 40S ribosomal protein S6

Catalog Number	S05-58
Quantity	1 mg
Specific Lot Number	K175-2

Purity	99.1%
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.