

Modified EIF2S Peptide

Synthetic peptide substrate derived from human EIF2S1

Catalog # E23-58B

Lot # I339-6

Product Description

The Modified EIF2S Peptide sequence (Modified-CILLSELSRRRIR) is derived from human EIF2S1 (46-57) and is suitable for use as the substrate for EIF2AK kinase family and MNK1.

Molecular Weight

The theoretical molecular weight is 1614.99

Purity

The purity was determined to be 95.88% by HPLC analysis.

Formulation

1mg of peptide supplied as a 1ml vial.

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.

Related Kinases

Modified EIF2S Peptide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
MNK1, Active	M54-10G
EIF2AK1, Active	H07-10G
EIF2AK2, Active	P80-11G
EIF2AK3, Active	E11-11G
EIF2AK4, Active	E12-11G

Modified EIF2S Peptide

Synthetic peptide substrate derived from human EIF2S1

Catalog Number	E23-58B
Quantity	1 mg in 1 ml volume
Specific Lot Number	I339-6
Purity	95.88%
Concentration	1 mg/ml
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.