

EIF2S Peptide

Synthetic peptide substrate derived from human EIF2S1

Catalog # E23-58

Lot # I267-2

Product Description

The 13 amino acids of EIF2S peptide sequence (CILLSELSRRRIR) is derived from human protein EIF2S1 between residues 46-57. This peptide is suitable for use as a substrate for MNK1 and EIF2AK family kinases.

Molecular Weight

The theoretical molecular weight is 1614.99

Purity

The purity was determined to be more than 95.88% 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.

Related Kinases

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

Product Name	Catalog Number
EIF2K 2, Active	P80-11G
EIF2K 3, Active	E11-11G
EIF2K4, Active	E12-11G
MNK1, Active	M54-10G

EIF2S Peptide

Synthetic peptide substrate derived from human EIF2S1

Catalog Number E23-58

Quantity 1 mg

Specific Lot Number I267-2

Purity >95.88%

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.