

IGF1Rtide

Synthetic peptide substrate derived from human IRS-1

Catalog # I15-58

Lot # N140-6

Product Description

The 14 amino acids of IGF1Rtide peptide sequence (KKKSPGEYVNIEFG) is derived from human IRS-1 protein residues 891-902.

Molecular Weight

The theoretical molecular weight is 1595.83

Purity

The purity was determined to be 97.2% 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

IGF1Rtide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
FLT1, active	F11-11G
IGF1R, active	I02-11H
JAK2, active	J02-11H
RET, active	R02-11G

IGF1Rtide

Synthetic peptide substrate derived from human IRS-1

Catalog # I15-58

Quantity 1 mg

Lot # N140-6

Purity 97.2%

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.