

IGF1Rtide

Synthetic peptide substrate derived from human IRS-1

Catalog # I15-58

Lot # G069-1

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 more than 95% 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
JAK2, active	J02-11H
RET, active	R02-11G

IGF1Rtide

Synthetic peptide substrate derived from human IRS-1

Catalog Number I15-58

Quantity 1 mg

Specific Lot Number G069-1

Purity	>95%
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.