

IRS1 (Y608) Peptide

Synthetic peptide substrate derived from mouse IRS1

Catalog # I40-58

Lot # G1309-5

Product Description

The synthetic IRS1 (Y608) Peptide [KKHTDDGYMPMSPGVA] is derived from mouse insulin receptor substrate 1 (amino acids 603-616) and is suitable as substrate for JAK1 and JAK2.

Molecular Weight

The theoretical molecular weight is 1,813.05.

Purity

The purity was determined to be 95.78% 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.

IRS1 (Y608) Peptide

Synthetic peptide substrate derived from mouse IRS1

Catalog #	I40-58
Quantity	1 mg
Specific Lot #	G1309-5
Purity	95.78%
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.