

GSK3 Substrate

Peptide derived from human muscle glycogen synthase 1

Catalog # G50-58

Lot # O1059-3

Product Description

The GSK3 Substrate peptide sequence (YRRAAVPPSPSLR-RHSSPHQ(pS)EDEEE) is based on human muscle glycogen synthase 1 (amino acid 636-661).

Molecular Weight

The theoretical molecular weight is 3029.13

Purity

The purity was determined to be 95.85% 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 25mM Tris-HCl buffer (pH 7.5) to a final concentration of 1mg/ml.

Related Kinases

GSK3 Substrate can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
GSK3 β , Active	G09-10G
GSK3 α , Active	G08-10G

GSK3 Substrate

Synthetic peptide substrate derived from human muscle glycogen synthase 1

Catalog # G50-58

Quantity 1 mg

Lot # O1059-3

Purity 95.11%

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.