

## GSK3 Substrate

Peptide derived from human muscle glycogen synthase 1

**Catalog # G50-58**

Lot # M071-3

### 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  $-20^{\circ}\text{C}$ . 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 $\beta$ , Active	G09-10G
GSK3 $\alpha$ , Active	G08-10G

## GSK3 Substrate

Synthetic peptide substrate derived from human muscle glycogen synthase 1

Catalog Number	G50-58
Quantity	1 mg
Specific Lot Number	M071-3
Purity	95.85%
Format	1 mg lyophilized powder
Stability	1yr at $-20^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-20^{\circ}\text{C}$ . 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**