

Crosstide

Synthetic peptide substrate derived from GSK3

Catalog # C51-58

Lot # K151-1

Product Description

The Crosstide peptide sequence (GRPRTSSFAEG) is based on the GSK3.

Molecular Weight

The theoretical molecular weight is 1164.2

Purity

The purity was determined to be 90.2% 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 distilled H₂O to a final concentration of 1mg/ml.

Related Kinases

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

| Product Name | Catalog Number |
|----------------|----------------|
| AKT1, Active | A16-10G |
| AKT2, Active | A17-10G |
| AKT3, Active | A18-10G |
| SGK1, Active | S06-11G |
| SGK2, Active | S07-10G |
| p70S6K, Active | R21-10H |
| RSK2, Active | R17-10G |

Crosstide

Synthetic peptide substrate derived from GSK3

| | |
|---------------------|---|
| Catalog Number | C51-58 |
| Quantity | 1 mg |
| Specific Lot Number | K151-1 |
| Purity | 90.2% |
| 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. |

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.