

## Crosstide

Synthetic peptide substrate derived from GSK3

**Catalog # C51-58**

Lot # Z1456-7

### 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 96.39% 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 H<sub>2</sub>O 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 #	C51-58
Quantity	1 mg
Lot #	Z1456-7
Purity	96.39%
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](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.**