

## SRC Substrate

Synthetic peptide substrate derived from the p34cdc2

**Catalog # S30-58**

Lot # K199-1

### Product Description

The SRC Substrate peptide sequence (KVEKIGEGTYGVVYK-amide) is based on the p34cdc2 (amino acid 6-20).

### Molecular Weight

The theoretical molecular weight is 1668.9

### Purity

The purity was determined to be 95.9% 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

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

Product Name	Catalog Number
SRC, Active	S19-10G
LYN B, Active	L13-10G

## SRC Substrate

Synthetic peptide substrate derived from the p34cdc2

Catalog Number S30-58

Quantity 1 mg

Specific Lot Number K199-1

Purity 95.9%

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.**