

SRC Substrate

Synthetic peptide substrate derived from the p34cdc2

Catalog # S30-58

Lot # N140-7

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 1668.97

Purity

The purity was determined to be 98.15% 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 20mM Tris-HCl, pH 7.5 solution 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	S19-18G
LYN B, Active	L13-10G

SRC Substrate

Synthetic peptide substrate derived from the p34cdc2

Catalog #	S30-58
Quantity	1 mg
Lot #	N140-7
Purity	98.15%
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.