

PKA Substrate

Synthetic peptide substrate derived from cAMP-dependent protein kinase inhibitor alpha

Catalog # C01-58

Lot # K115-2

Product Description

The PKA Substrate peptide sequence (CGRTGRRNSI-amide) is based on the cAMP-dependent protein kinase inhibitor alpha (amino acid 15-23).

Molecular Weight

The theoretical molecular weight is 1119.3

Purity

The purity was determined to be 98% 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

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

Product Name	Catalog Number
AKT2, Active	A17-10G

PKA Substrate

Synthetic peptide substrate derived from cAMP-dependent protein kinase inhibitor alpha

Catalog Number C01-58

Quantity 1 mg

Specific Lot Number K115-2

Purity 98%

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.