

AMARA Peptide

Synthetic peptide substrate evaluated as a substrate for SIK and AMPK

Catalog # A11-58

Lot # D196-6

Product Description

The AMARA substrate peptide sequence (AMARAASAAALARRR) is routinely evaluated as a substrate for SIK and AMPK.

Molecular Weight

The theoretical molecular weight is 1542.84

Purity

The purity was determined to be 98.4% 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 25mM Tris-HCl buffer (pH 7.5) to a final concentration of 1mg/ml.

Related Kinases

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

Product Name	Catalog Number
SIK	S14-11H
AMPK (A1/B1/G1)	P47-10H
AMPK (A1/B1/G2)	P55-10H
AMPK (A1/B1/G3)	P56-10H
AMPK (A1/B2/G1)	P50-10H
AMPK (A2/B1/G1)	P48-10H

AMARA Peptide

Synthetic peptide substrate evaluated as a substrate for SIK and AMPK

Catalog Number	A11-58
Quantity	1 mg
Specific Lot Number	D196-6
Purity	98.4%
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.