

Abltide

Synthetic peptide substrate derived from Abl

Catalog # A02-58

Lot # K055-2

Product Description

The Abltide peptide sequence (EAIYAAPFAKKK) is based on the C-terminus of Abl.

Molecular Weight

The theoretical molecular weight is 1336.6

Purity

The purity was determined to be 97.7% 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

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

Product Name	Catalog Number
ABL1, Active	A03-11H

Abltide

Synthetic peptide substrate derived from Abl

Catalog Number	A02-58
Quantity	1 mg
Specific Lot Number	K055-2
Purity	97.7%
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.