

## Abltide

Synthetic peptide substrate derived from Abl

**Catalog # A02-58**

Lot # M360-3

### Product Description

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

### Molecular Weight

The theoretical molecular weight is 1336.61

### 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 20mM Tris-HCl, pH 7.5 solution 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
ABL1, Active	A03-18H
ABL2, Active	A04-11H

## Abltide

Synthetic peptide substrate derived from Abl

Catalog Number      A02-58

Quantity      1 mg

Specific Lot Number      M360-3

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.