

MLC Peptide

Synthetic peptide substrate derived from human MYL9

Catalog # M09-58

Lot # D196-8

Product Description

The MLC Peptide sequence (AKRPQRATSNVFS) represents the phosphate-accepting domain of human MYL9 (amino acids 12-23) and can be used as a substrate for PHKG1, PHKG2 and DAPK2.

Molecular Weight

The theoretical molecular weight is 1461.66

Purity

The purity was determined to be 96.9% 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

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

Product Name	Catalog Number
DAPK2, Active	D01-11G
PHKG2, Active	P20-10G

MLC Peptide

Synthetic peptide substrate derived from human MYL9

Catalog Number	M09-58
Quantity	1 mg
Specific Lot Number	D196-8
Purity	96.9%
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.