

MRCL3 Peptide

Synthetic peptide substrate derived from human MRCL3

Catalog # M56-58

Lot # F528-5

Product Description

The synthetic substrate MRCL3 Peptide (KKRPQRATSN-VFAM-NH₂) is derived from human myosin regulatory light chain MRCL3 (amino acid 11-24) and can be used for MLCK, MYLK2 and MYLK3 kinase assay.

Molecular Weight

The theoretical molecular weight is 1631.90

Purity

The purity was determined to be 96.43% 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

MRCL3 Peptide substrate can be utilized as a substrate for MLCK, MYLK2 and MYLK3 kinase assay.

Product Name	Catalog Number
MLCK, Active	M62-11G
MYLK2, Active	M63-10H
MYLK3, Active	M64-10G

MRCL3 Peptide

Synthetic peptide substrate derived from human MRCL3

Catalog Number	M56-58
Quantity	1 mg
Specific Lot Number	F528-5
Purity	96.43%
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.