

PKCtide

Synthetic peptide substrate derived from protein kinase C epsilon

Catalog # P15-58

Lot # M360-4

Product Description

The PKCtide peptide sequence (ERM₁PRKRQGSVRRRV) is based on protein kinase C epsilon (amino acid 149-164).

Molecular Weight

The theoretical molecular weight is 2067.5

Purity

The purity was determined to be 95.1% 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

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

Product Name	Catalog Number
PKCβI, Active	P62-10G
PKCε, Active	P65-10G
PKCη, Active	P67-10G
PKCγ, Active	P66-10G
PKCα, Active	P61-10G
PKCβII, Active	P63-10G
PKCδ, Active	P64-10G
PKCι, Active	P68-10G
PKCμ, Active	P72-10G
PKCν, Active	P73-10G
PKCζ, Active	P75-10G

PKCtide

Synthetic peptide substrate derived from protein kinase C epsilon

Catalog Number	P15-58
Quantity	1 mg
Specific Lot Number	M360-4

Purity	95.1%
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.