

CSKtide

Synthetic peptide substrate evaluated for CSK and TNK1 assay

Catalog # C63-58

Lot # F528-4

Product Description

The synthetic peptide CSKtide (KKKEEYFFFG-NH₂) contains a tyrosine protein kinase phosphorylation site and can be used for CSK and TNK1 assay.

Molecular Weight

The theoretical molecular weight is 1435.50

Purity

The purity was determined to be 97.41% 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

CSKtide peptide substrate can be utilized as a substrate for CSK and TNK1.

CSKtide

Synthetic peptide substrate evaluated for CSK and TNK1 assay

Catalog Number C63-58

Quantity 1 mg

Specific Lot Number F528-4

Purity 97.41%

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.