

DYRKtide

Synthetic peptide substrate evaluated for DYRK family assay

Catalog # D96-58

Lot # W235-4

Product Description

The synthetic peptide (RRRFRPASPLRGPPK) contains a serine/threonine protein kinase phosphorylation site and is routinely evaluated as a substrate for DYRK family kinases.

Molecular Weight

The theoretical molecular weight is 1791.15

Purity

The purity was determined to be 95.31% 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.

DYRKtide

Synthetic peptide substrate evaluated for DYRK family assay

Catalog Number D96-58

Quantity 1 mg

Specific Lot Number W235-4

Purity 95.31%

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.