

Tyr-phosphopeptide-2

Synthetic peptide substrate derived from epidermal growth factor receptor isoform 1

Catalog # T70-58

Lot # K211-1

Product Description

The Tyr-phosphopeptide-2 sub peptide sequence (DADEY(p)LIPDQGG) is based on mouse epidermal growth factor receptor isoform 1 (amino acid 1014-1024).

Molecular Weight

The theoretical molecular weight is 1315.3

Purity

The purity was determined to be 95.3% 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

Tyr-phosphopeptide-2 can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
PTPN6, Active	P33-20G
PTPN11, Active	P38-21G
PTPRF, Active	P53-21G

Tyr-phosphopeptide-2

Synthetic peptide substrate derived from epidermal growth factor receptor isoform 1

Catalog Number T70-58

Quantity 1 mg

Specific Lot Number K211-1

Purity 95.30%

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.