

Poly (6:2:5:1 Ala,Glu,Lys,Tyr) Peptide

Synthetic peptide used as tyrosine kinase substrate

Catalog # P60-58

Lot # F724-3

Product Description

The Poly (6:2:5:1 Ala,Glu,Lys,Tyr) substrate peptide is used as universal substrate for protein tyrosine kinases.

Molecular Weight

The peptide molecular weight is around 20,000 to 30,000.

Purity

The purity was determined to be 95%.

Concentration

1mg of peptide supplied in 1 ml of 25mM Tris-HCl buffer (pH 7.5) to a final concentration of as 1mg/ml.

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.

Poly (6:2:5:1 Ala,Glu,Lys,Tyr) Peptide

Synthetic peptide used as tyrosine kinase substrate

Catalog Number P60-58

Quantity 1 mg

Specific Lot Number F724-3

Purity 95%

Concentration 1 mg/ml

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.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.