

Poly (4:1 Glu, Tyr) Peptide

Synthetic peptide used as tyrosine kinase substrate

Catalog # P61-58

Lot # A1316-5

Product Description

The Poly (4:1 Glu, Tyr) substrate peptide is used as universal substrate for protein tyrosine kinases.

Molecular Weight

The peptide molecular weight is around 20,000 Da to 50,000 Da.

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 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 (4:1 Glu, Tyr) Peptide

Synthetic peptide used as tyrosine kinase substrate

Catalog #	P61-58
Quantity	1 mg
Lot #	A1316-5
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 frozen.

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.