

RS Repeat Peptide

Synthetic peptide substrate evaluated for CLK family assay

Catalog # R55-58

Lot # F722-5

Product Description

The synthetic peptide (GRSRSRSRSRSRSR) contains serine/threonine protein kinase phosphorylation sites and is routinely evaluated as a substrate for CLK, EIF2AK and SRPK family kinases.

Molecular Weight

The theoretical molecular weight is 1934.13

Purity

The purity was determined to be 97.52% 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

RS Repeat Peptide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
CLK4, Active	C61-10G
EIF2AK1, Active	H07-10G
EIF2AK2, Active	P80-11G
EIF2AK3, Active	E11-11G
EIF2AK4, Active	E12-11G
MNK1, Active	M54-10G
SRPK1, Active	S21-10G
SRPK2, Active	S22-10G

RS Repeat Peptide

Synthetic peptide substrate evaluated for CLK family assay

Catalog #	R55-58
Quantity	1 mg
Lot #	F722-5
Purity	97.52%
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.