

## GRKtide

Synthetic peptide substrate evaluated for GRK family assay

**Catalog # G46-58**

Lot # R339-6

### Product Description

The synthetic peptide sequence (CRRREEEESAAA) is routinely evaluated as a substrate for GRK family kinases, including GRK2, GRK3, GRK4 and GRK7.

### Molecular Weight

The theoretical molecular weight is 1536.2

### Purity

The purity was determined to be 96.1% 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

GRKtide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
GRK2, Active	A14-10G
GRK7, Active	G05-10G

## GRKtide

Synthetic peptide substrate evaluated for GRK family assay

Catalog Number	G46-58
Quantity	1 mg
Specific Lot Number	R339-6
Purity	96.1%
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.