

## GRK4 Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # G02-31G**

Lot # P1749-5

### Product Description

Recombinant human GRK4 (140-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number for GRK4 is [BC117320](#).

### Gene Aliases

GPRK2L, GPRK4, GRK4a, IT11

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

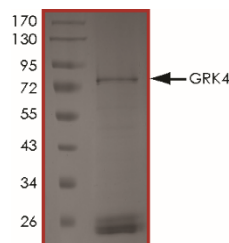
### Scientific Background

GRK4 or G protein-coupled receptor kinase 4 is a member of the guanine nucleotide-binding protein (G protein)-coupled receptor kinase subfamily of the Ser/Thr protein kinase family which phosphorylates the activated forms of G protein-coupled receptors thus initiating its deactivation and it also play an important role in receptor desensitization. GRK4 protein has kinase activity and that it interacts with and is inhibitable by calmodulin (1). GRK4 has been linked to both genetic and acquired hypertension. GRK4 is highly expressed in testis (2).

### References

1. Salles, M. et al: G protein-coupled receptor kinase GRK4: molecular analysis of the four isoforms and ultrastructural localization in spermatozoa and germinal cells. J. Biol. Chem. 272: 10188-10195, 1997.
2. Premont, R. T. et al: Characterization of the G protein-coupled receptor kinase GRK4: identification of four splice variants. J. Biol. Chem. 271: 6403-6410, 1996.

### Purity



The purity of GRK4 was determined to be **>70%** by densitometry, GRK4 approx. MW **75kDa**.

## GRK4 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	G02-31G
Lot #	P1749-5
Purity	>70%
Concentration	0.05 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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

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