

Data Sheet

GRK6

Human, recombinant, N-terminal GST tag

Catalog #: 40066

Lot #: 110215 **Conc.:** 0.1 mg/ml

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

Stability: >12 months at -80°C

References:

1. Gaudreau, R., *et al. J. Biol Chem.* **277**: 31567-31576 (2002).
2. Gainetdinov, R.R., *et al. Neuron* **38**: 291-303 (2003).

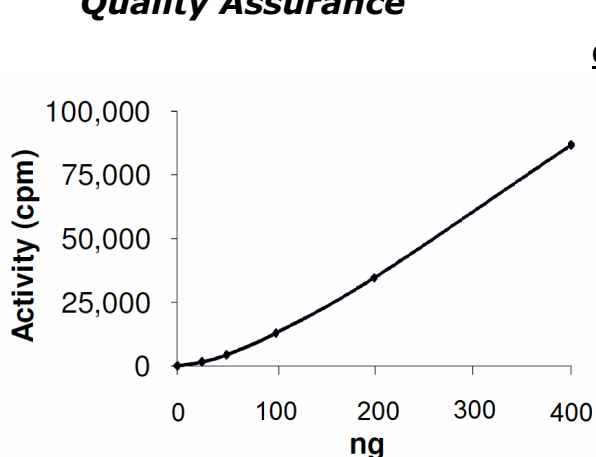
Description: Human GRK6, also known as GPRK6 or FLJ3213 (GenBank Accession No. BC_017272), full length, with N-terminal GST tag, MW ~93 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 11 pmol/min/μg

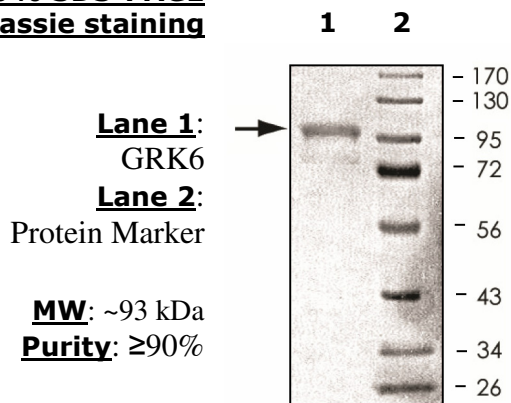
Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate GRK6 with PLKtide substrate and [³³P]-ATP at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



10% SDS-PAGE
Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com