

## Data Sheet

### GRK5

Human, recombinant, N-terminal GST tag

**Catalog #:** 40063

**Lot #:** 110419      **Conc.:** 0.1 mg/ml

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

**Stability:** >12 months at -80°C

**References:**

1. Haribabu, B., *et al. Proc. Natl. Acad. Sci. USA* **90**: 9398-9402 (1993).
2. Fan, J., *et al. Nature Med.* **9**: 315-321 (2003).

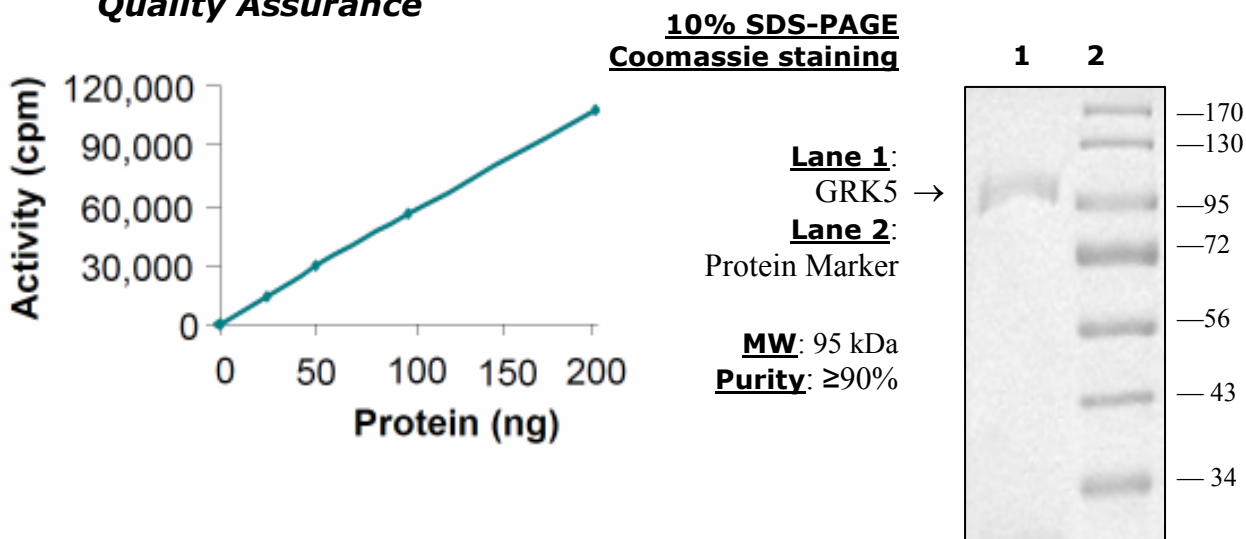
**Description:** Human GRK5, also known as GPRK5 (GenBank Accession No. NM\_005308), full length, with N-terminal GST tag, MW ~95 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

**Specific Activity:** 40 pmol/min/μg

Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate GRK5 with 0.5 mg/ml casein substrate, 12.5 mg/ml phosphatidylcholine and [<sup>33</sup>P]-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

### Quality Assurance



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)