

Data Sheet

GCK

Human, recombinant, N-terminal GST tag

Catalog #: 40107

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. Pombo, C.M., *et al. Nature.* **377:** 750-754 (1995).
2. Ren, M., *et al. Proc. Natl. Acad. Sci.* **93:** 5151-5155 (1996).

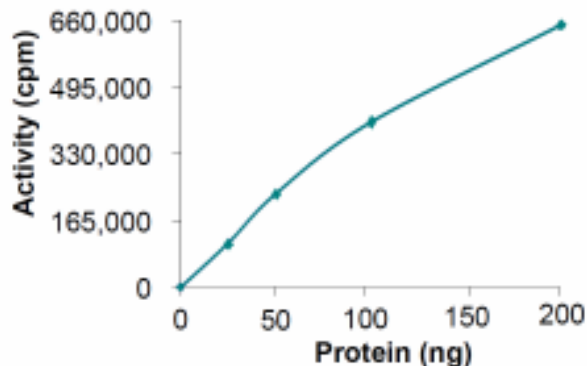
Description: Human GCK, also known as MAP4K2, GenBank Accession No. NM_004579, full length, with N-terminal GST tag, MW ~125 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 290 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 GCK with 0.25 mg/ml Myelin Basic Protein (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

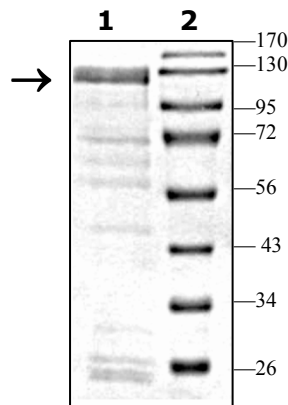


10% SDS-PAGE Coomassie staining

Lane 1:
GCK

Lane 2:
Protein Marker

MW: 125 kDa
Purity: ≥80%



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