

Data Sheet

ROCK2, Active

Human, recombinant, N-terminal GST tag

Catalog #: 40086

Lot#: 110506

Conc.: 0.1 mg/ml

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

Stability: >6 months at -80°C

References:

1. Zhao, Y M. *et al. Dev Biol.* 2004 Nov 1;**275(1)**:183-91.
2. Noma, K. *et al. Am J Physiol Cell Physiol.* 2006 Mar;**290(3)**:C661-8.

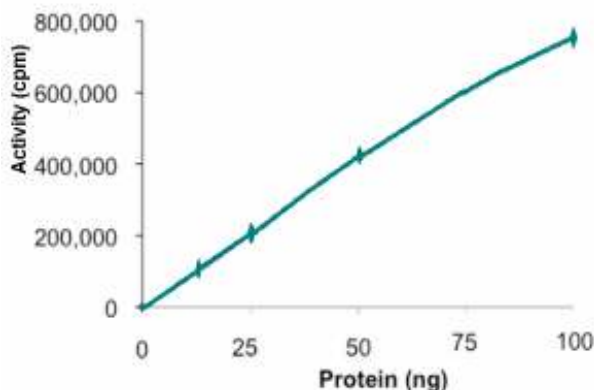
Description: Recombinant human ROCK2 (5-554) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is NM_004850.

Specific Activity: 51 pmol/min/μg

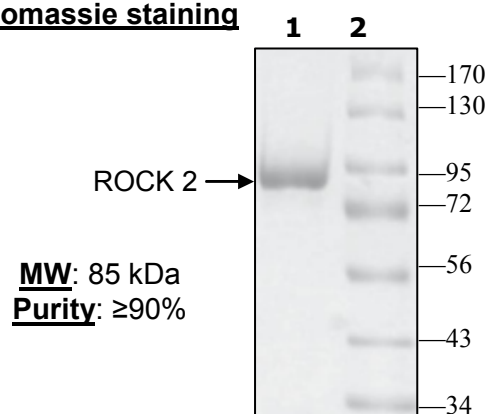
Assay Buffer: 5 mM MOPS, 2.5 mM β-glycerophosphate, 5 mM MgCl₂, 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 2 mM ATP. Incubate ROCK1 with 0.25 mg/ml S6K peptide substrate and [³³P]-ATP at 30°C for 15 min., then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation 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 **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com