

Data Sheet

ACK1

Human, recombinant, N-terminal GST-tag

Catalog #: 40400

Lot#: 110420-20 **Conc.:** 0.67 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 16 mM
glutathione, 3 mM DTT, and 20% glycerol.

Stability: >6 months at -80°C

References:

1. Eley, L. *et al.*, *Biochem. Biophys. Res. Commun.* **371** (4), 877-882 (2008).
2. Grovdal, L.M. *et al.*, *Exp. Cell Res.* **314** (6), 1292-1300 (2008).

Description:

Human Ack1 (GenBank Accession No. NM_005781), (a-a 110-476) with N-terminal GST tag, MW=68 kDa, expressed in Sf9 insect cells via a baculovirus expression system.

Assay conditions: Add 50 µl reaction mix (ACK1 diluted in 1x kinase buffer supplemented with 0.2 mM ATP) to wells coated with substrate. Incubate for 1h. Add Europium-labeled anti-p-Tyr antibody and incubate for 1h. Time resolved fluorescence is measured at room temperature at $\lambda_{exc}=330\pm 40$ nm and $\lambda_{em}=620\pm 10$ nm.

Application:

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

Quality Assurance

10% SDS-PAGE Coomassie staining

Lane 1:

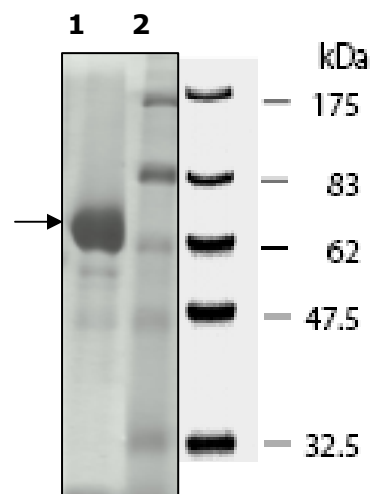
ACK1

Lane 2:

Protein Marker
BioLabs (#P7708L)

MW: 68 kDa

Purity: ≥62%



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