

Data Sheet

ERK3, GST-Tag

Human, recombinant, N-terminal GST tag

Catalog #: 100366

Lot #: 190423

Conc.: 0.05 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: At least 12 months at -80°C.
 Avoid multiple freeze/thaw cycles.

References:

1. Boulton, T.G., *et al.*, *Cell*. **65**: 663-675, (1991).
2. Meloche, S., *et al.*, *Oncogene*. **13**: 1575-1579, (1996).

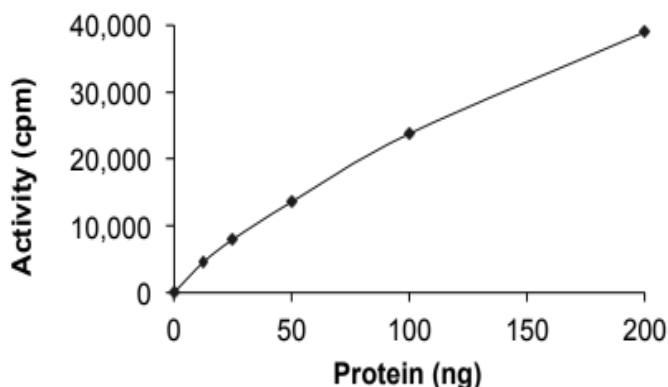
Description: Human ERK3, also known as MAPK6, HsT17250, p97MAPK, and PRKM6, GenBank Accession No. NM_002748, a.a. 1-721(end), with N-terminal GST tag, expressed in a Baculovirus infected Sf9 cell expression system. This protein runs at a higher M.W. by SDS-PAGE. Calculated MW =109 kDa.

Specific Activity: 22 pmol/min/μg.

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

Quality Assurance

Specific Activity



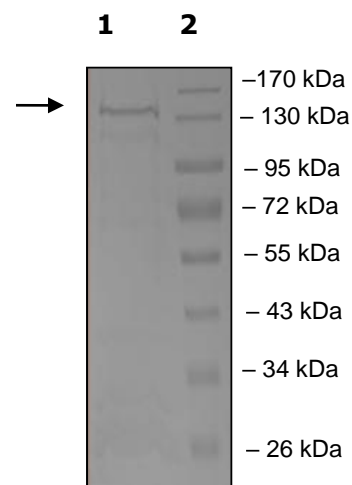
4-20% SDS-PAGE Coomassie staining

Lane 1:
ERK3

Lane 2:
Protein Marker

MW: 109 kDa

Purity: ≥75%



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** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

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** 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)