

## Data Sheet

### **ERK7, GST-Tag (Mouse)**

Mouse, recombinant, N-terminal GST tag

**Catalog #:** 100365

**Lot #:** 190422

**Conc.:** 0.10 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:**

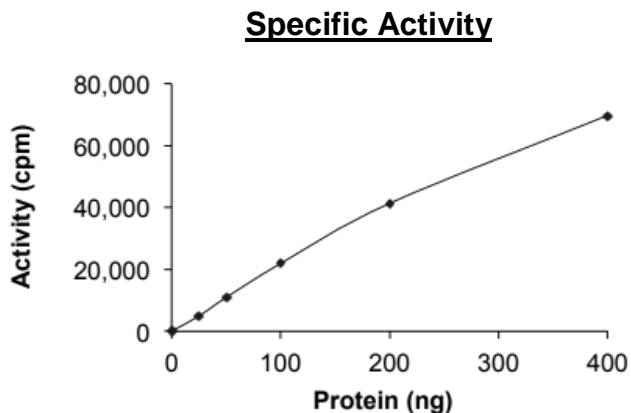
1. Rossi, M., *et al.*, *J. Biol Chem.* **286(10)**: 8507-8522, (2011).
2. Abe, M.K., *et al.*, *J. Biol Chem.* **(276)24**: 21272-21279, (2001).

**Description:** Mouse ERK7, also known as MAPK15, GenBank Accession No. BC048082, a.a. 1-549 (full length), with N-terminal GST tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=86 kDa.

**Specific Activity:** 7 pmol/min/μg

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

## **Quality Assurance**



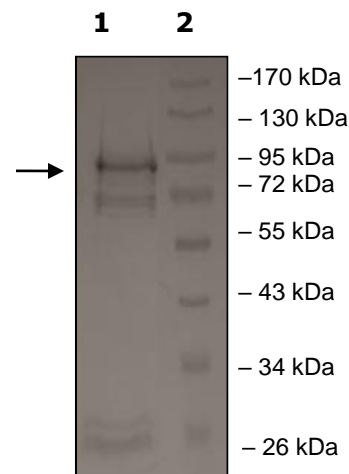
**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
ERK7

**Lane 2:**  
Protein Marker

**MW:** 86 kDa

**Purity:** ≥70%



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)