

## Data Sheet

### MEG-2, GST-tag

Human, recombinant, N-terminal GST-tag  
**Catalog #:** 30020  
**Lot#:** 171205      **Conc.:** 0.20 mg/ml

**Formulated in:** 25 mM Tris-HCl, pH 8.0, 67 mM NaCl, 1.4 mM KCl, 50% glycerol, 10 mM glutathione, and 3 mM DTT.

**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles.

**References:**

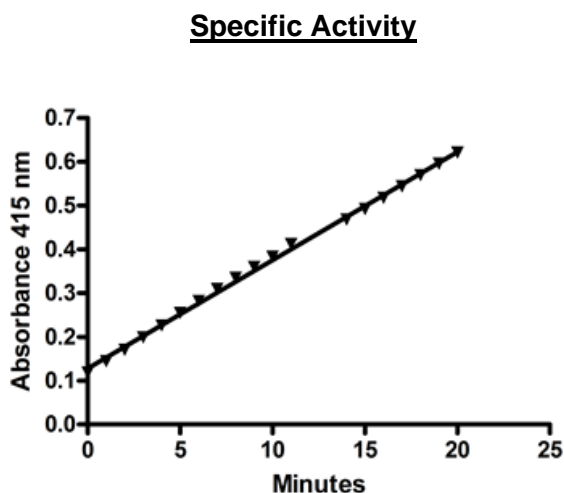
1. Q. Ying, *et al.*, *J. Cell Biochem.* 2002, **86**: 79
2. H. Huynh, *et al.* *J. Immunol.* 2003, **171**: 6661

**Description:** Human MEG-2, also known as, Protein Tyrosine Phosphatase Non-Receptor Type 9 (PTP-9) and PTPase-MEG2, Genbank Accession No.: M83738, a.a. 285-593(end) with N-terminal GST-tag, expressed in an *E. coli* expression system. MW=60 kDa.

**Specific Activity:** 1000 pmole/min/μg. Assay conditions: Enzyme reaction 400 ng/rxn (20 μl) containing 20 mM HEPES, pH 7.2, 100 mM NaCl, 2 mM EDTA, 2 mM DTT, 1 mM pNPP and PTPase was monitored at 415 nm continuously with readings at 1-minute intervals for 20 minutes.

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

## Quality Assurance

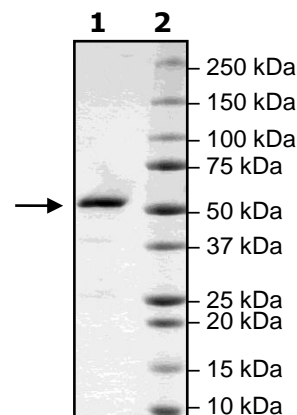


**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2 μg MEG2  
**Lane 2:**  
Protein Marker

**MW:** 60 kDa

**Purity:** ≥90%



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)



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto: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)