

## Data Sheet

### DRAK2, GST-tag (Mouse)

Mouse, Recombinant N-terminal GST-tag

**Catalog #:** 100444

**Lot #:** 190717-1      **Conc:** 0.17 mg/ml

**Formulated in:** 50 mM HEPES, pH 7.5, 0.001% Brij-35, 1 mM DTT, 10 mM MgCl<sub>2</sub>, 150 mM KCl, 5% Glycerol, 0.1 mM EDTA, and 1mM Glutathione.

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

**References:**

1. McGargill, M. *et al.*, *Immunity* **21** (6): 781-791 (2004).
2. Mao, J. *et al.*, *J. Biol. Chem.* **281**: 12587-12595 (2006).

**Description:** Mouse DRAK2, also known as DAP kinase-related apoptosis-inducing protein kinase 2, and death-associated protein kinase-related 2, GenBank Accession No. NM\_133810, a.a. 2-372 (end), expressed in a baculovirus-infected Sf9 cell expression system, MW = 68 kDa.

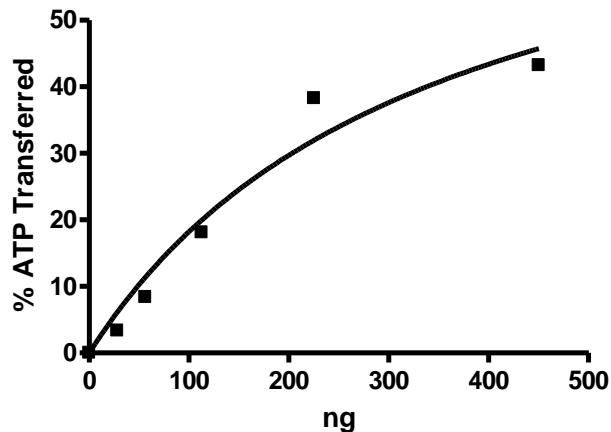
**Specific Activity:** 5 pmole/min/μg

**Assay Conditions:** Reaction was performed in 50 mM Tris pH 7.4, 5 mM MgCl<sub>2</sub>, 1 mM DTT, 20 μM ATP, 0.2 mg/ml MRCL3 Peptide. Reaction was initiated with the addition of Mouse DRAK2 and was incubated at 30°C for 45 min. The luminescence was read with the Kinase Glo Max reagent (Promega).

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

### Quality Assurance

**Specific Activity**



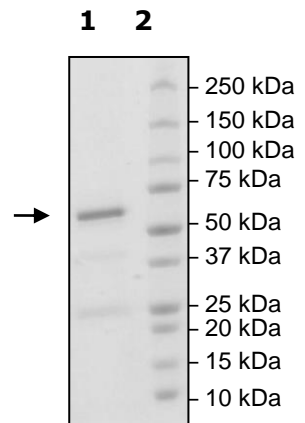
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
DRAK2

**Lane 2:**  
Protein Marker

**Purity:** ≥75%

**MW:** 68 kDa



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)

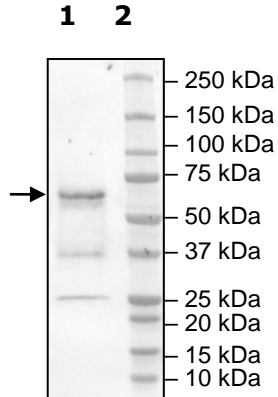
**Western Blot**

**4-20% SDS-PAGE**  
**Anti-GST**

**Lane 1:**  
**Mouse DRAK2**

**Lane 2:**  
**Protein Marker**

**MW:** 68 kDa



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)