

## Data Sheet

### mCot

Mouse, recombinant, N-terminal His-GST-tag

**Catalog #:** 41050

Lot#: 110901

Conc.: 0.1 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.4 mM KCl, 5 mM glutathione, 134 mM imidazole, 3 mM DTT and 20% glycerol.

**Stability:** >6 months at -80 °C

**References:**

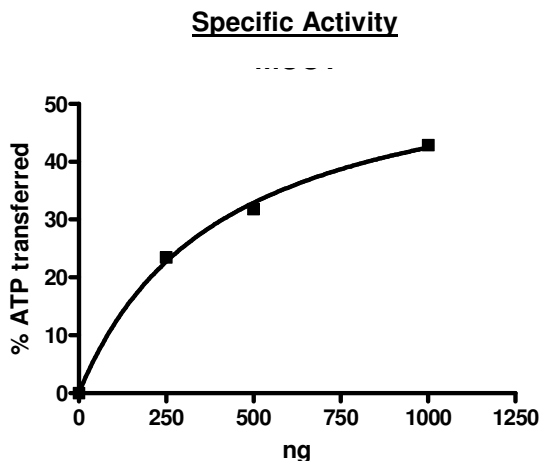
1. Jeong, J.H. *et al.*, *PLoS One*. 2011 Jan 18; **6(1)**:e16205.
2. Soria-Castro, I. *et al.* *J Biol Chem*. 2010 Oct 29; **285(44)**:33805-15.

**Description:** Mouse Cot kinase (MAP3K8) GenBank Accession No. NM\_007746, a.a. 30-397, with N-terminal His-GST-tag, expressed in a Baculovirus infected Sf9 cell expression system. The theoretical MW is 70 kDa.

**Specific Activity:** 8.5 pmole/min/mg  
**Assay Conditions:** Reaction was done in a kinase buffer (40 mM Tris, 10 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA) containing 5 µg inactive MEK1 (BPS #40075), 20 µM ATP, and mCOT for 60 min at 30°C. The amount of ATP transferred was measured by Kinase-Glo plus (Promega).

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

## Quality Assurance

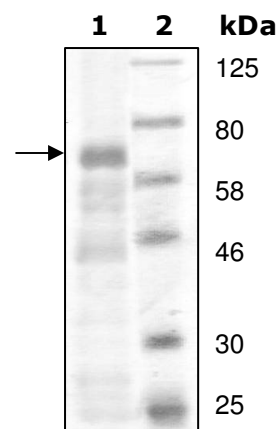


**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 2.5 µg mCot

**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 70 kDa  
**Purity:** ≥38%



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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)