

Data Sheet

Mouse Cot (MAP3K8), His-GST-Tags Recombinant

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

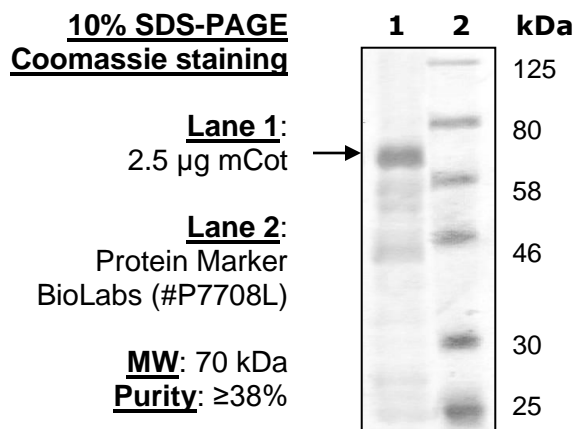
References:

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.

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

Quality Assurance



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: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com