

Data Sheet

Cot (MAP3K8)

Human, recombinant, N-terminal GST-tag

Catalog #: 40050

Lot#: 130311G **Conc.:** 0.25 mg/ml

Formulated in: 25 mM Tris-HCl, pH 8.0,
 100 mM NaCl, 50% glycerol, and 3 mM
 DTT.

Stability: >6 months at -80 °C

References:

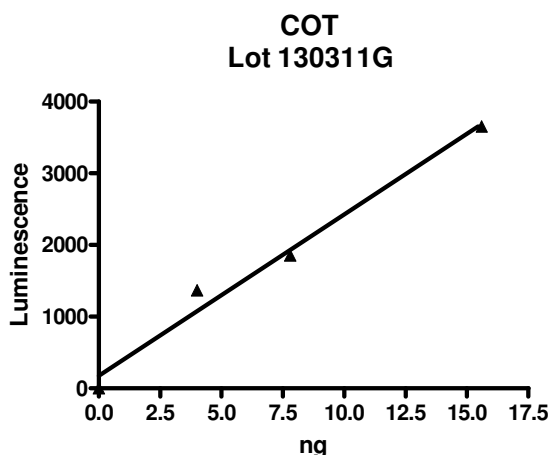
1. Stafford, M.J. *et al.*, *FEBS Lett.* **580**,
 4010-4014 (2006)
2. Babu, G. *et al.*, *J. Biol. Chem.* **281**,
 14041-14047 (2006)

Description: Human Cot (GenBank
 Accession No. NM_005204), (a.a. 30-
 397), with N-terminal GST-HIS₆-tag,
 expressed in a Baculovirus infected Sf9 cell
 expression system. The theoretical MW is
 72kDa.

Specific Activity: 47 pmol/min/μg
 Assay conditions: 100 μl of 240 nM
 biotinylated MEK1 peptide was coated onto
 a neutravidin plate for 1 h at 37 °C. After
 overnight blocking with TBS superblock, the
 reaction was done with 10 μM ATP for 60
 min at RT, and phospho-MEK1 was detected
 using an anti-phospho-MEK1 antibody.
 Specific activity was determined relatively by
 comparing with a standard kinase under the
 same condition.

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

Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:

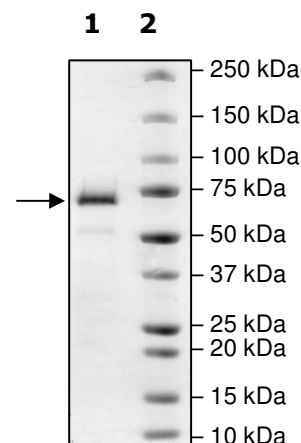
2 μg Cot

Lane 2:

Protein Marker

MW: 72 kDa

Purity: ≥95%



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

Please visit our website at: www.bpsbioscience.com