

Data Sheet

MAPK14 (p38 α), His-tag, active

Human, Recombinant, N-terminal His-tag
Catalog #: 40243
Lot #: 140124-P **Conc.:** 0.65 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
 20, 3 mM DTT, and 20% glycerol

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

References:

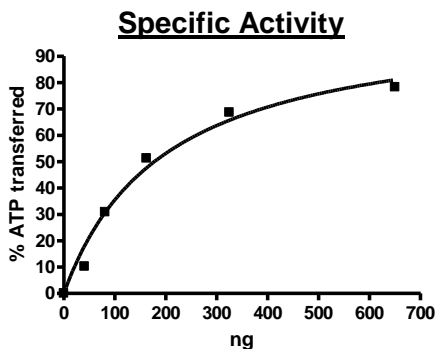
1. Hatherley, D., *et al.*, *Mol. Cell*, 2008,
 31:266-77
2. Baba, N., *et al.*, *J. Exp. Med.*, 2013,
 210(6):1251-63

Description: Mitogen-activated protein
 kinase 14 (MAPK14) GenBank Accession
 No. NM_001315, a.a. 2-360 (end), with
 N-terminal His-tag, MW 42.3 kDa,
 expressed in a baculovirus-infected Sf9
 cell expression system. This protein is
 activated by *in vitro* phosphorylation and
 SEC.

Specific Activity: ≥ 100 pmol/min/ μg
 Assay conditions: Assay was performed
 in kinase buffer containing 0.2 mM DTT
 using P38 peptide substrate (0.1 mg/ml)
 and 20 μM ATP. Reaction was done at
 30 $^{\circ}\text{C}$ for 40 min. Amount of ATP
 transferred was calculated using Kinase-
 Glo reagent (Promega).

Application: Useful for studying enzyme
 kinetics, screening inhibitors, and
 selectivity profiling.

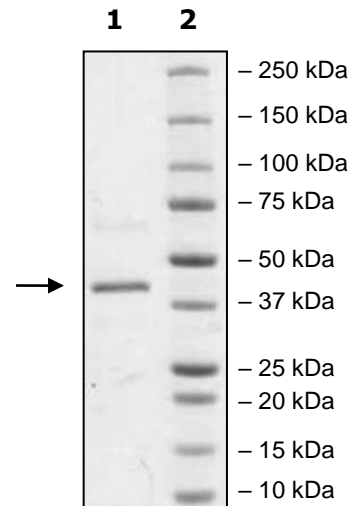
Quality Assurance



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

Lane 1:
 1 μg MAPK14
Lane 2:
 Protein Marker

MW: 42.3 kDa
Purity: $\geq 80\%$



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 **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Sequence

MHHHHHSQERPTFYRQELNKTIVEVPERYQNLSPVGS GAYGSVCAAFDTKTGLRVAVKKLSRPFQ
SIIHAKRTYRELRLLLKMKHENVIGLLDVFTPARSLEEFNDVYLVTHLMGADLNNIVKCQKLTDDH
VQFLIYQILRGLKYIHSADIIHRDLKPSNLAVNEDCELKILDFGLARHTDDEMTGYVATRWRAP
IMLNWMHYNQTVDIWSVGCIMAELLTGRTLFPGTDHINQLQQIMRLTGTPPAYLINRMP SHEARNY
IQSLTQMPKMN FANVFIGANPLAVDLLEKMLVLDSDKRITAAQALAHAYFAQYHDPDDEPVADPYD
QSFESRDL LIDEWKSLTYDEVISFVPPPLDQEEMES

Red – His-tag

Black – MAPK14 (a.a. 2-360(end))

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 **1.858.259.5734**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com