

## Data Sheet

### MAPK14/P38 alpha inactive

Human, recombinant, N-terminal GST tag  
**Catalog #:** 40091  
**Lot#:** 111012      **Conc.:** 0.2 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5,  
150mM NaCl, 0.1mM EGTA, 0.1mM EDTA,  
0.1mM PMSF, 25% glycerol, and 0.25 mM  
DTT.

**Stability:** 12 months at  $-70^{\circ}\text{C}$

**References:**

1. Bell C.E. et al., *Hum Reprod.* 2009 Jun;24(6):1373-86.
2. Carraro G. et al., *Dev Biol.* 2009 Sep 15;333(2):238-50.

**Description:**

Human MAPK14 (GenBank Accession No. NM\_139012), full length, with N-terminal GST tag, MW=65 kDa, expressed in an E.coli expression system.

**Application:**

MAPK14 is a substrate for MKK6 kinase assay.

### Quality Assurance

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

**Lane 1:**  
MAPK14  
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 65 kDa  
**Purity:**  $\geq 90\%$

