

## Product Information

### MAP2K4

Product Number : **07-144**

#### Product description

Full-length human MAP2K4 [1-399(end) amino acids of accession number NP\_003001.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP\_002392.2] using baculovirus expression system. GST-MAP2K4 was purified by using glutathione sepharose chromatography.

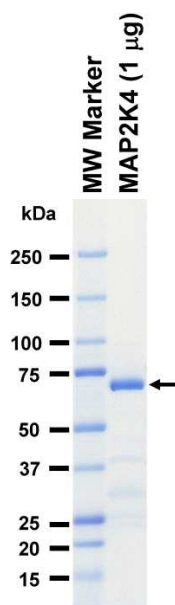
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.  
Avoid repeating freeze-thaws.

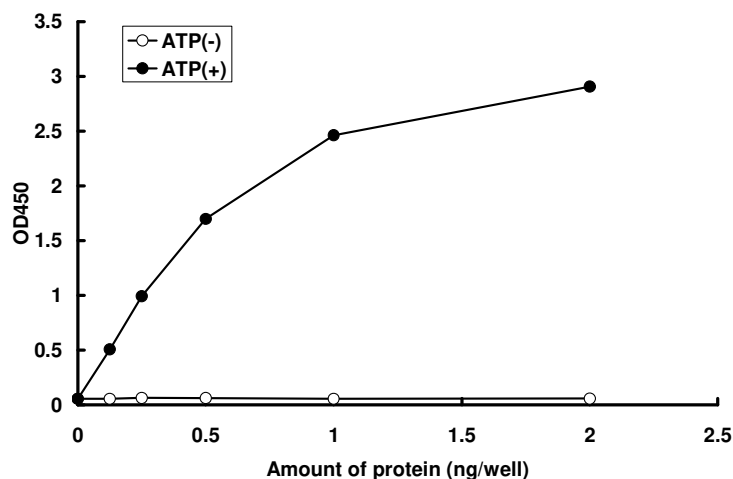
#### SDS-PAGE



Purity: 81 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with GST-fused substrate protein, and after stopping kinase reaction by EDTA, the reaction solution was transferred into glutathione-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG) .

Substrate : JNK1 [inactive mutant]

ATP : 100  $\mu$ M