

## Product Information

### MAP3K4

 Product Number : **07-106**

#### Product description

Human MAP3K4, catalytic domain [1312-1608(end) amino acids of accession number NP\_005913.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system. GST-MAP3K4 was purified by using glutathione sepharose chromatography.

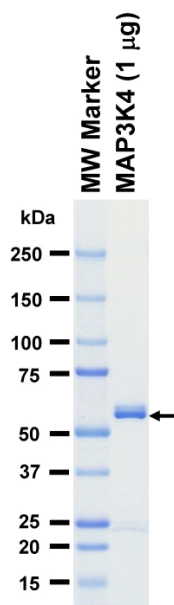
#### Storage buffer:

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.

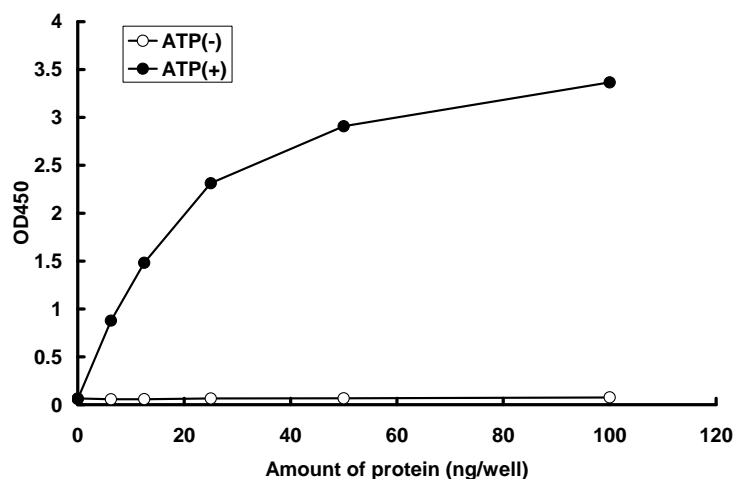
#### SDS-PAGE



Purity: 91 %

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 : MAP2K7 [inactive mutant]

ATP : 100  $\mu$ M