

## Product Information

### MAP3K14

 Product Number : **07-102**

#### Product description

Human MAP3K14, catalytic domain [319-947(end) amino acids of accession number NP\_003945.2] was expressed as N-terminal GST-fusion protein (96 kDa) using baculovirus expression system. GST-MAP3K14 was purified by using glutathione sepharose chromatography.

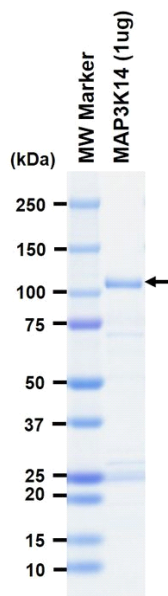
#### Storage buffer:

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.

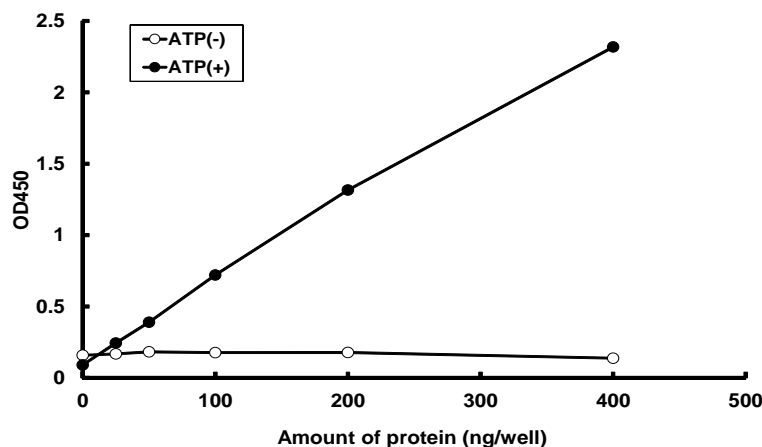
#### SDS-PAGE



Purity: 52 %

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

#### Activity data



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

Substrate : BTN-IKKa[inactive mutant]

 ATP : 100  $\mu$ M