

## Product Information

### MAP3K2

Product Number : **07-004**

#### Product description

Human MAP3K2, catalytic domain [337-620(end) amino acids of accession number NP\_006600.2] was expressed as N-terminal His-tagged protein (35 kDa) using baculovirus expression system. His-tagged MAP3K2 was purified by using Ni-NTA affinity chromatography.

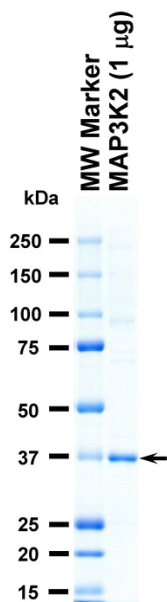
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

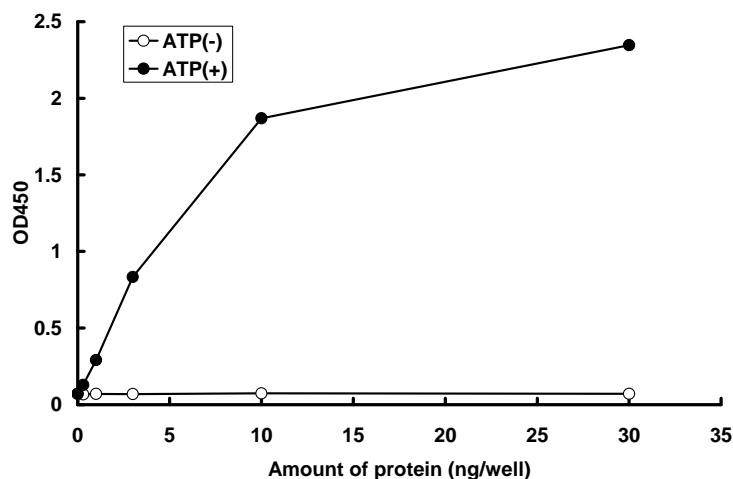
#### SDS-PAGE



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 : MAP2K7

ATP : 2  $\mu$ M