

Product Information

MAP3K2(FL)

Product Number : **07-154**

Product description

Full-length human MAP3K2 [1-619(end) amino acids of accession number NP_006600.3] was expressed as N-terminal GST-fusion protein (97 kDa) using baculovirus expression system. GST-MAP3K2 was purified by using glutathione sepharose chromatography.

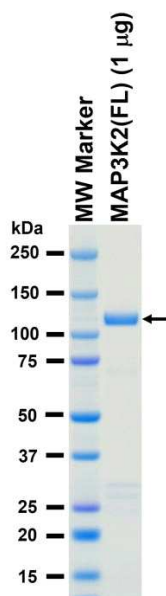
Storage buffer:

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.

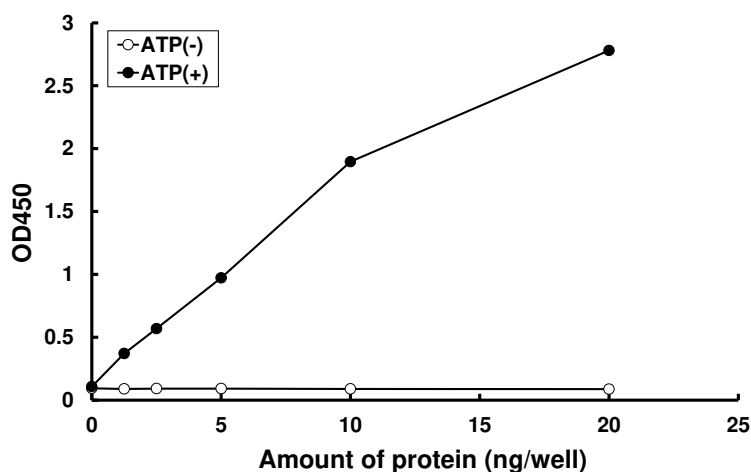
SDS-PAGE



Purity: 88 %

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-MAP2K7[inactive]

ATP : 100 µM