

Product Information

MAP2K7

 Product Number : **07-148**

Product description

Full-length human MAP2K7 [1-419(end) amino acids of accession number NP_660186.1] was co-expressed as N-terminal GST-fusion protein (75 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] using baculovirus expression system. GST-MAP2K7 was purified by using glutathione sepharose chromatography.

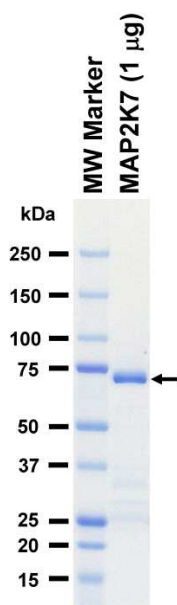
Storage buffer:

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.

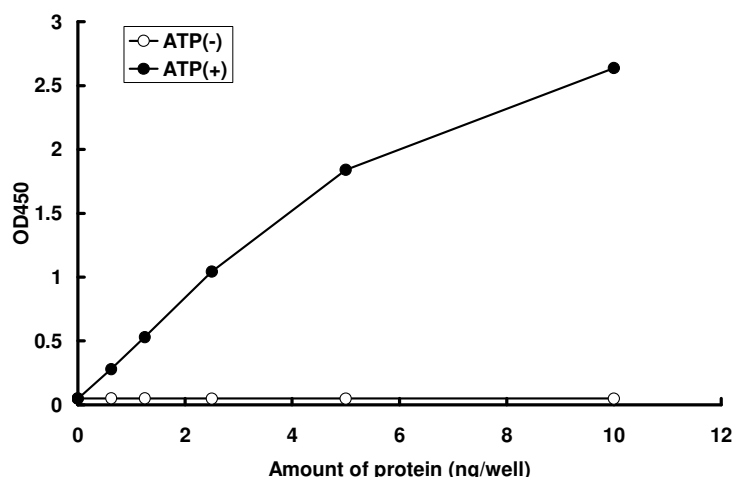
SDS-PAGE



Purity: 87 %

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 μ M