

Product Information

MAP2K6

 Product Number : **07-146**

Product description

Full-length human MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human His-tagged MLK3 [99-398 amino acids of accession number NP_002410.1] using baculovirus expression system. GST-MAP2K6 was purified by using glutathione sepharose chromatography.

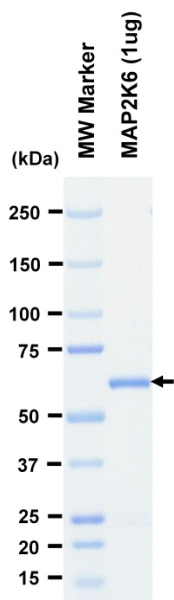
Storage buffer:

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.

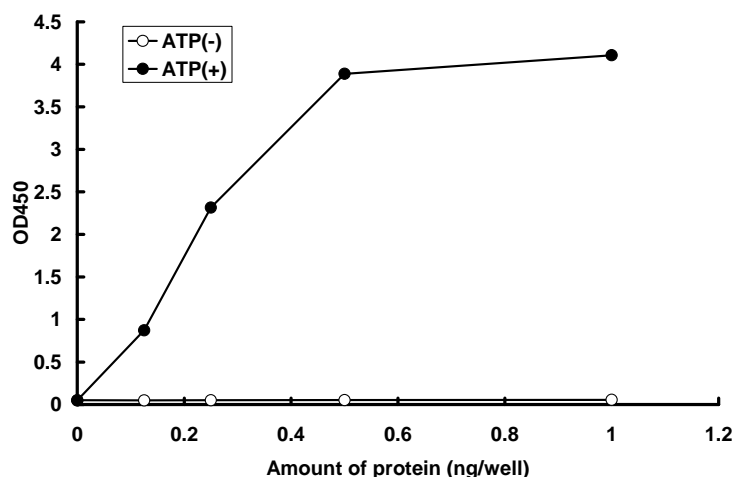
SDS-PAGE



Purity: 94 %

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 : p38 α (9-352) [inactive mutant]

 ATP : 100 μ M