

Product Information

MAP2K6[inactive]

Product Number : **07-146-10**

Product description

Full-length human MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] was expressed as N-terminal GST-fusion protein (65kDa) using E. coli expression system. GST-MAP2K6 was purified by using glutathione sepharose chromatography.

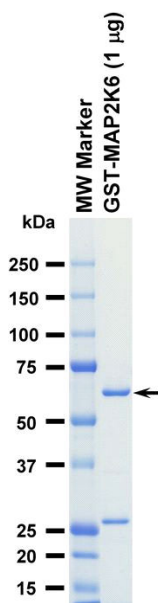
Storage buffer:

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.

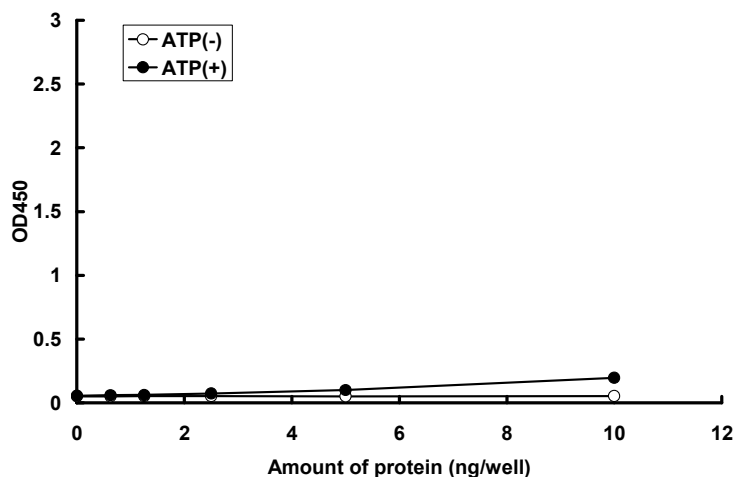
SDS-PAGE



Purity: 59 %

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.

Substrate : p38α(9-352)

ATP : 100 µM