

Product Information

MAP2K1[inactive]

 Product Number : **07-141-10**

Product description

Full-length human MAP2K1 [1-393(end) amino acids of accession number NP_002746.1] was expressed as N-terminal GST-fusion protein (70 kDa) using E.coli expression system. GST-MAP2K1 was purified by using glutathione sepharose chromatography.

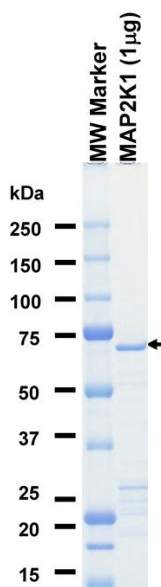
Storage buffer:

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

Storage and Handling:

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

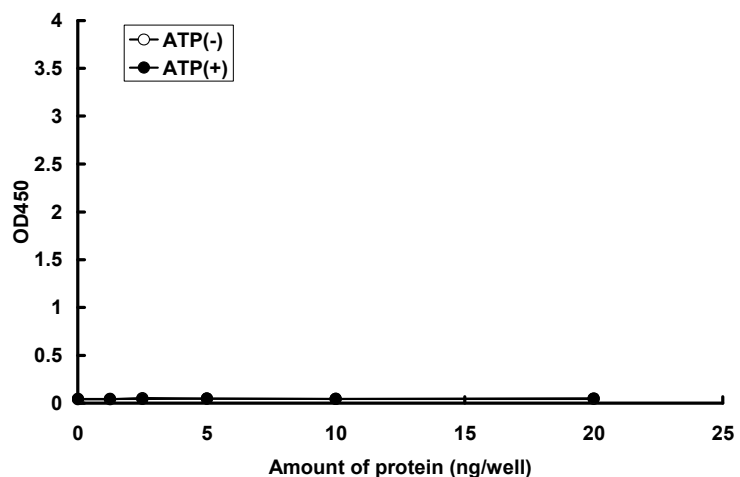
SDS-PAGE



Purity: 56 %

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 : Erk2

 ATP : 100 μ M