

Product Information

MAP2K1

 Product Number : **07-041**

Product description

Full-length human MAP2K1 [1-393(end) amino acids of accession number NP_002746.1] was coexpressed as N-terminal His-tagged protein (47 kDa) with human GST-RAF1 [306-648(end) amino acids and Y340D, Y341D of accession number NP_002871.1] using baculovirus expression system. His-tagged MAP2K1 was purified by using Ni-NTA affinity chromatography and glutathione sepharose chromatography.

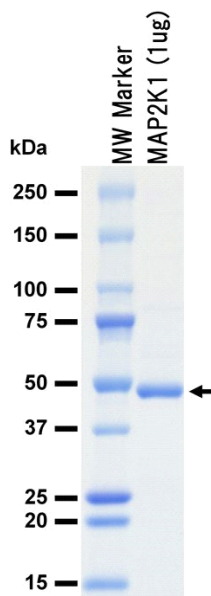
Storage buffer:

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.

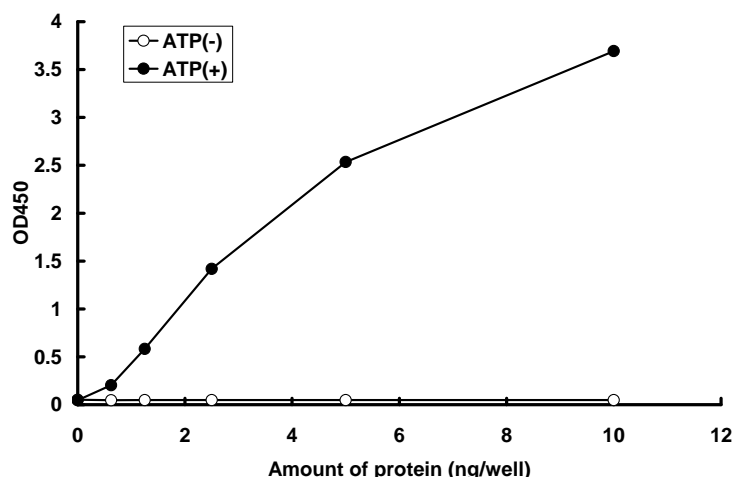
SDS-PAGE



Purity: 97 %

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 : Erk2 [inactive mutant]

 ATP : 20 μ M