

Product Information

MAP2K2

Product Number : **07-142**

Product description

Full-length human MAP2K2 [1-400(end) amino acids of accession number NP_109587.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged RAF1 [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] using baculovirus expression system. GST-MAP2K2 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

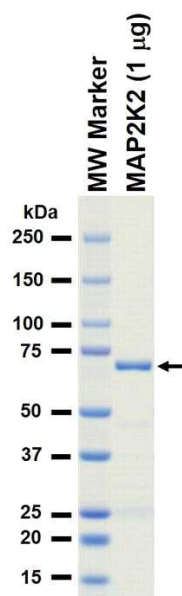
Storage buffer:

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.

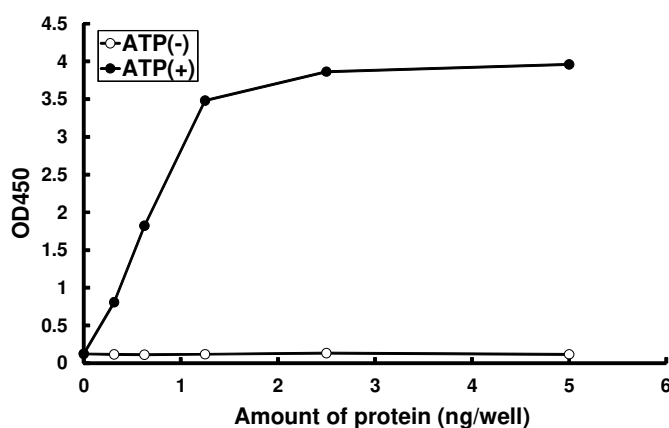
SDS-PAGE



Purity: 84 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated substrate protein/peptide, and after stopping kinase reaction by EDTA, the reaction solution was transferred into streptavidin-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG).

Substrate : BTN-Erk2[inactive]

ATP : 100 µM