

Product Information

MAP2K5

Product Number : **07-145**

Product description

Full-length human MAP2K5 [1-448(end) amino acids of accession number NP_660143.1] was co-expressed as N-terminal GST-fusion protein (77 kDa) with human His-tagged MAP3K3[1-626(end) amino acids of accession number NP_002392.2], CDC37 and HSP90 using baculovirus expression system. GST-MAP2K5 was purified by using glutathione sepharose chromatography.

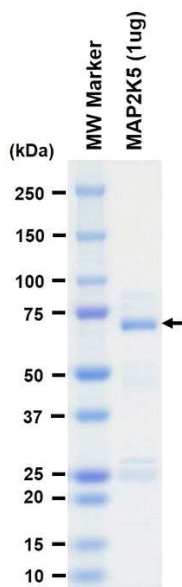
Storage buffer:

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.

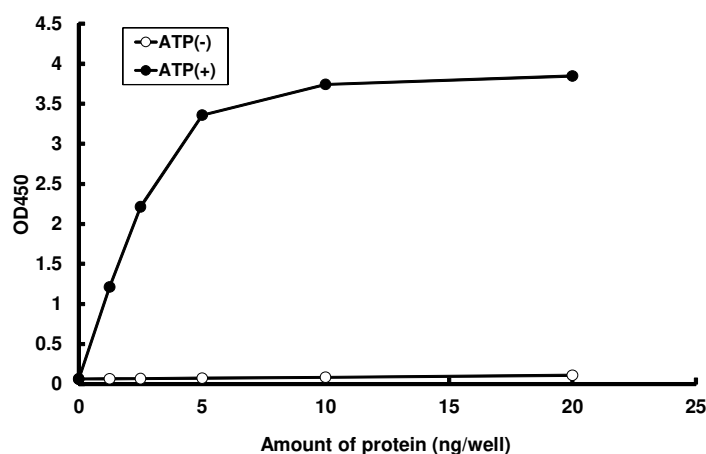
SDS-PAGE



Purity: 58 %

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 : Erk5 [inactive mutant]

ATP : 100 μ M