

## Product Information

### MAP2K1[P124L]

 Product Number : **07-150**

#### Product description

Full-length human MAP2K1 [1-393(end) amino acids and P124L of accession number NP\_002746.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with human His-tagged RAF1 [306-648 amino acids and Y340D and Y341D of accession number NP\_002871.1] using baculovirus expression system. GST-MAP2K1[P124L] was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

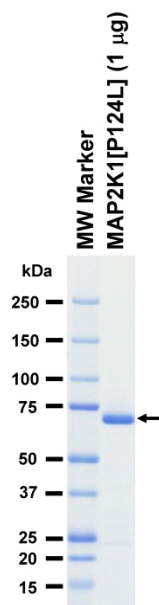
#### Storage buffer:

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.

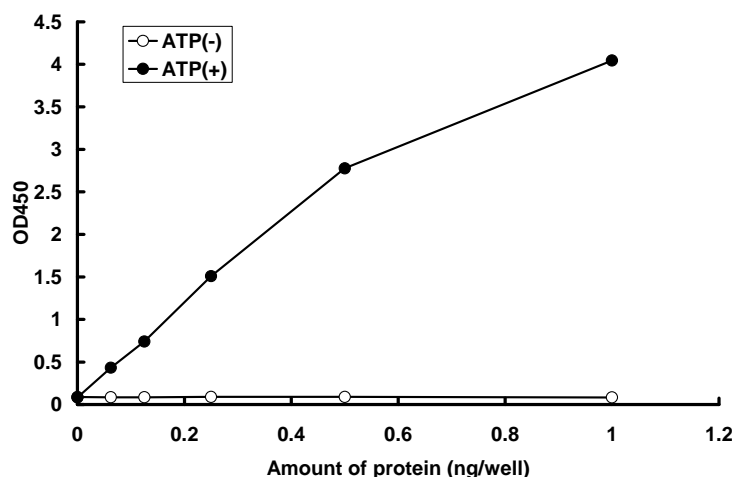
#### SDS-PAGE



Purity: 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 : 100 µM