

## Product Information

### p38 $\alpha$ (MAPK14)

Product Number : **04-152**

#### Product description

Truncated human p38 $\alpha$  [9-352 amino acids of accession number NP\_620581.1] was expressed as N-terminal GST-fusion protein (66 kDa) using E. coli expression system. GST-p38 $\alpha$  was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38 $\alpha$  was purified using glutathione sepharose chromatography.

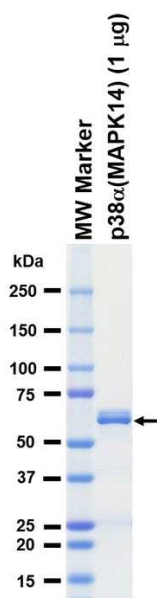
#### Storage buffer:

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.

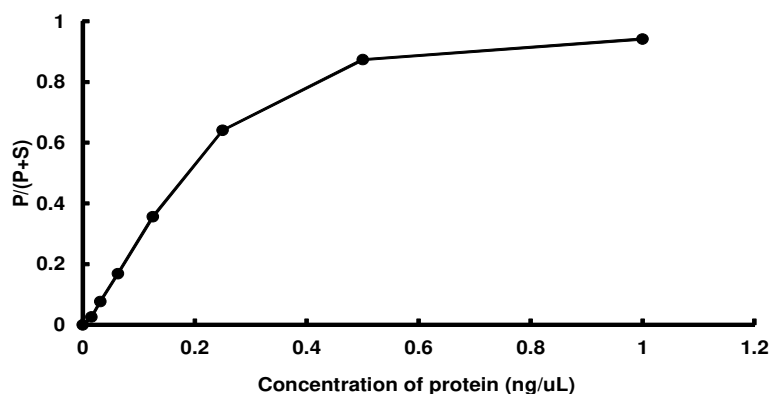
#### SDS-PAGE



Purity: 89 %

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

#### Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Modified Erk14

ATP : 100  $\mu$ M