

Product Information

p38 γ (MAPK12)

Product Number : **04-155**

Product description

Full-length human p38 γ [1-367(end) amino acids of accession number NP_002960.2] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38 γ was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38 γ was purified using glutathione sepharose chromatography.

Storage buffer:

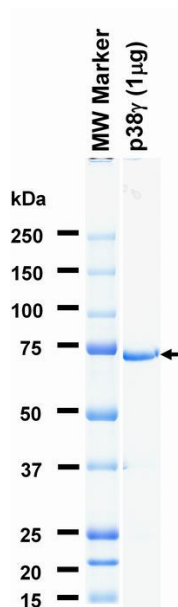
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

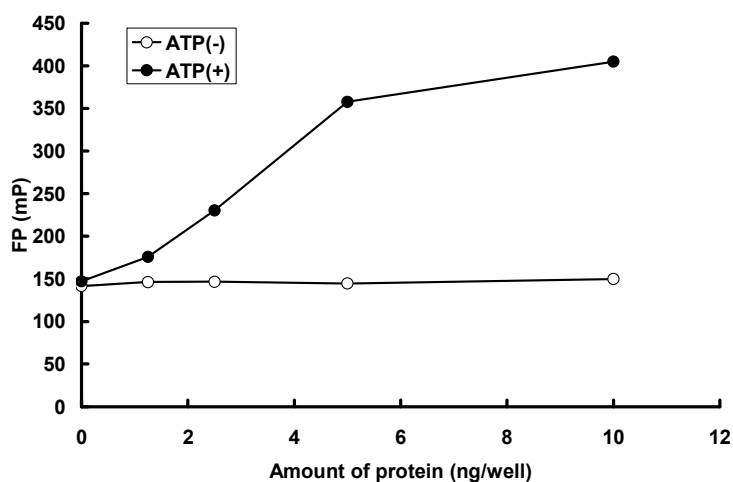
SDS-PAGE



Purity: 97 %

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

Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : Erktide

ATP : 3 μ M