

Product Information

BTN-p38 α (MAPK14)

 Product Number : **04-452-20N**

Product description

Truncated human p38 α [9-352 amino acids of accession number NP_620581.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (43 kDa) with human His-MAP2K6[1-334(end) amino acids of accession number NP_002749.2] and His-MLK3[99-398 amino acids of accession number NP_002410.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

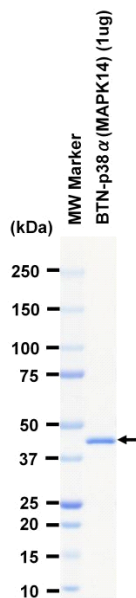
Storage buffer:

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.

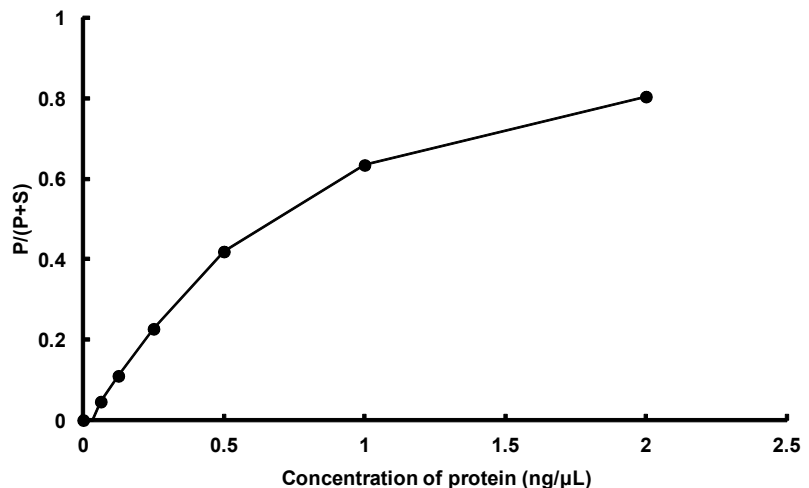
SDS-PAGE



Purity: 100 %

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

Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : Modified Erktide

 ATP : 100 μ M