

Product Information

TAK1-TAB1(MAP3K7)

 Product Number : **09-019**

Product description

Fused gene of human TAK1 [1-303 amino acids of accession number NP_663304.1] and human TAB1 [437-504 amino acids of accession number NP_006107.1] was expressed as N-terminal His-tagged protein (45kDa) using baculovirus expression system. His-tagged TAK1-TAB1 was purified by using Ni-NTA affinity chromatography.

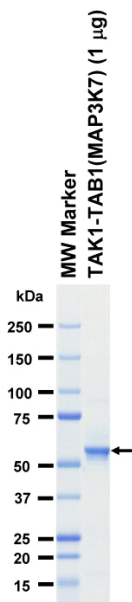
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
 1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

Store at -80C.
 Avoid repeating freeze-thaws.

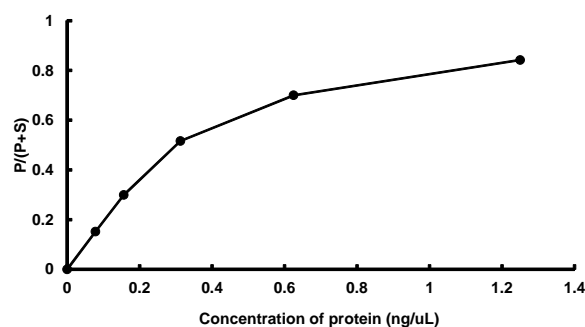
SDS-PAGE



Purity: 96 %

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 : LRRKtide

ATP : 100 μM