

Product Information

IKK β (IKBKB)

Product Number : **05-084**

Product description

Truncated human IKK β [1-662 amino acids of accession number NP_001547.1] was expressed as N-terminal His-tagged protein (77 kDa) using baculovirus expression system. His-tagged IKK β was purified by using Ni-NTA affinity chromatography and anion exchange chromatography.

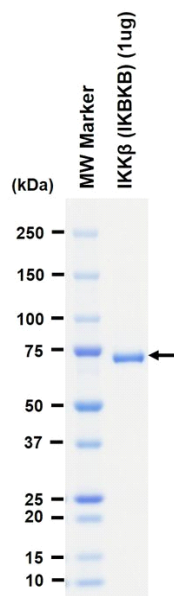
Storage buffer:

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.

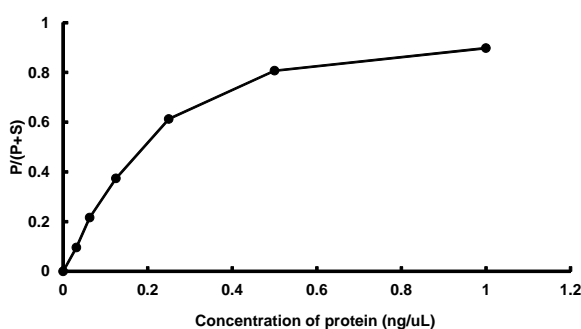
SDS-PAGE



Purity: 98 %

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 I κ B α -derived peptide

ATP : 100 μ M