

Product Information

ZAP70

Product Number : **08-377**

Product description

Full-length human ZAP70 [1-619(end) amino acids of accession number NP_001070.2] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-ZAP70 was purified by using glutathione sepharose chromatography. GST-ZAP70 was cleaved by PreScission protease and GST-free ZAP70 (71 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

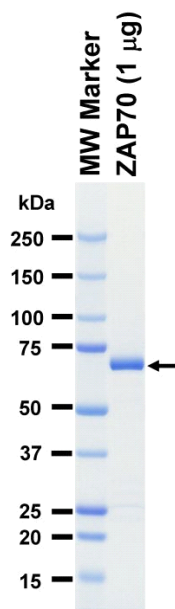
Storage buffer:

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.

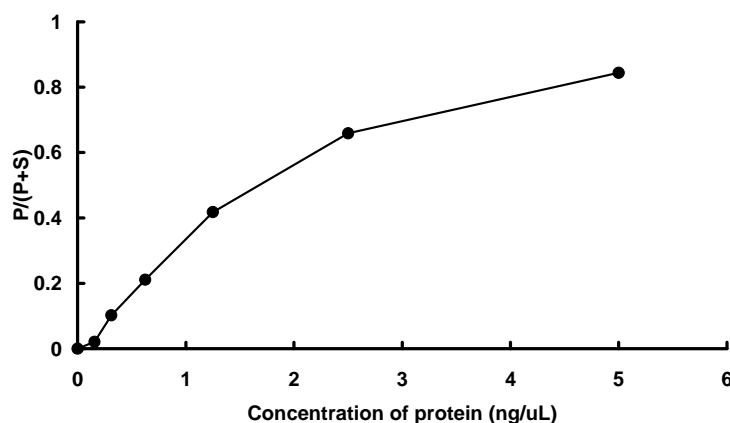
SDS-PAGE



Purity: 96 %

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

Activity data



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

Substrate : Blk/Lyntide

ATP : 100 µM