

Product Information

BUB1/BUB3

Product Number : **05-187**

Product description

Full-length human BUB1 [1-1085 (end) amino acids of accession number NP_004327] was co-expressed as N-terminal GST-fusion protein (149 kDa) with DYKDDDDK tagged BUB3 [1-328 (end) amino acids of accession number NP_004716] using baculovirus expression system. GST-BUB1 was purified by using glutathione sepharose chromatography.

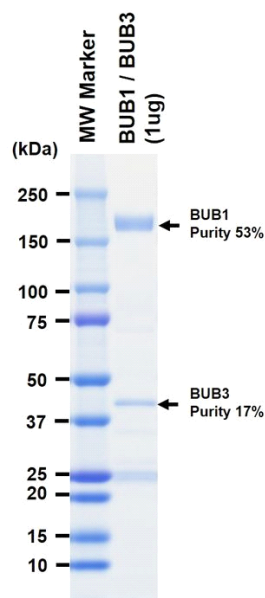
Storage buffer:

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.

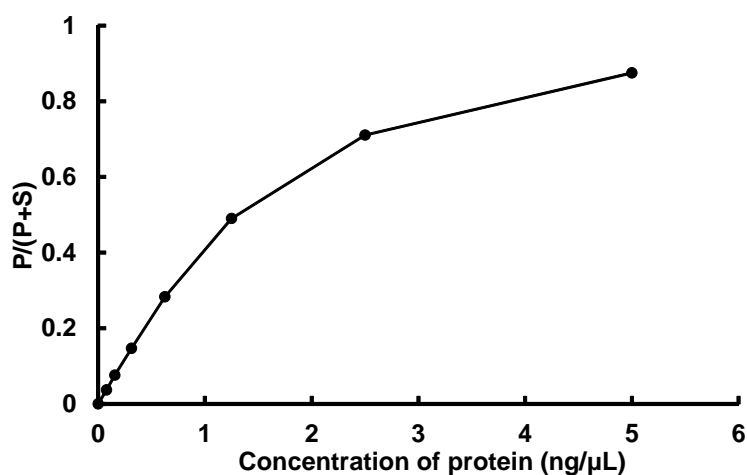
SDS-PAGE



Purity: 70 %

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 : H2A peptide

ATP : 100 μM