

Product Information

ALK[C1156Y]

Product Number : **08-530**

Product description

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and C1156Y of accession number BAG10812.1] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[C1156Y] was purified by using glutathione sepharose chromatography.

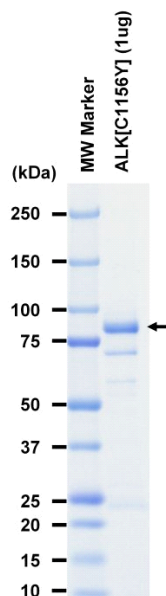
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

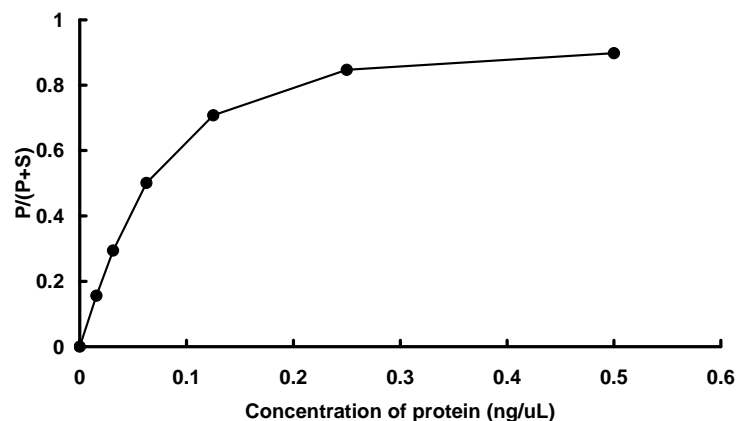
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 71 %
 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 : Srctide

ATP : 100 μM