

## Product Information

### NPM1-ALK

Product Number : **08-517**

#### Product description

Fused gene of human fusion NPM1-ALK [1-680 amino acids of accession number BAA08343.1] was expressed as N-terminal GST-fusion protein (103kDa) using baculovirus expression system. GST-NPM1-ALK 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

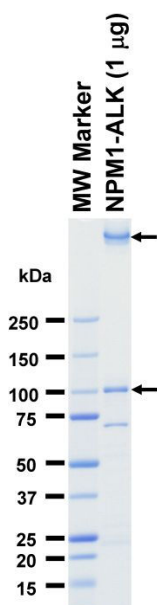
#### Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

Use within 12 months from the date of shipment.

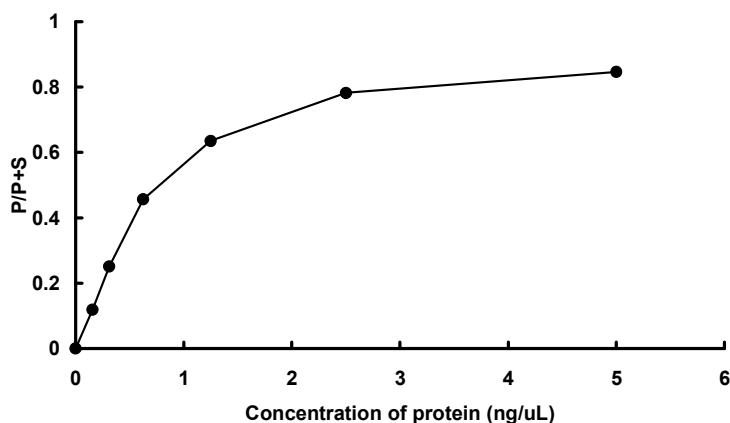
#### SDS-PAGE



Purity: 81 %

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 : Src tide

ATP : 100 µM