

## Product Information

### AKT2

Product Number : **01-102**

#### Product description

Human AKT2, catalytic domain [120-481(end) amino acids of accession number NP\_001617.1] was co-expressed as N-terminal GST-fusion protein (69 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP\_002604.1] using baculovirus expression system. GST-AKT2 was purified by using glutathione sepharose chromatography.

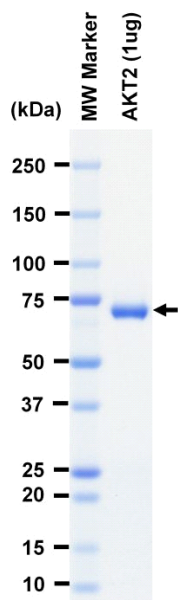
#### Storage buffer:

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.

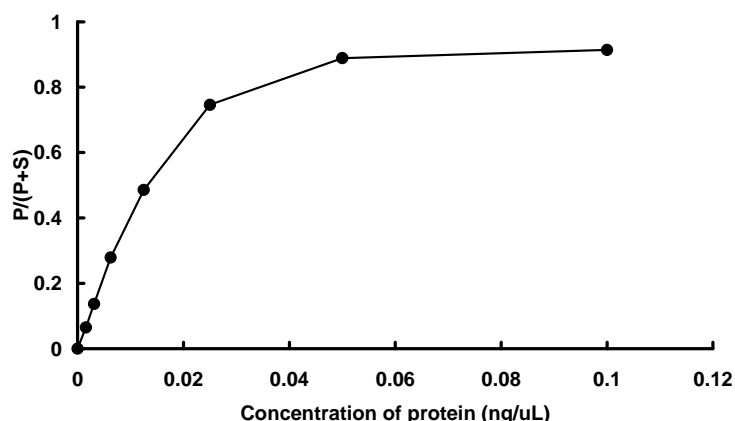
#### SDS-PAGE



Purity: 99 %

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 : Crosstide

ATP : 100  $\mu$ M