

Product Information

AKT3

Product Number : **01-103**

Product description

Human AKT3, catalytic domain [108-479(end) amino acids of accession number NP_005456.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-AKT3 was purified by using glutathione sepharose chromatography.

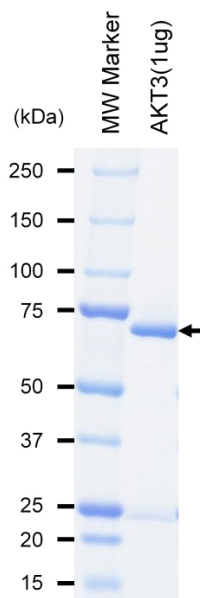
Storage buffer:

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.

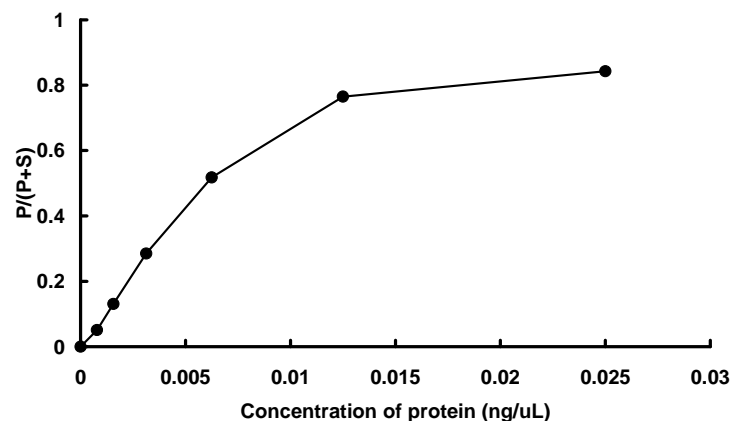
SDS-PAGE



Purity: 87 %

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 μ M