

Product Information

DAPK3

Product Number : **02-136**

Product description

Full-length human DAPK3 [1-454(end) amino acids of accession number NP_001339.1] was expressed as N-terminal GST-fusion protein (80 kDa) using baculovirus expression system. GST-DAPK3 was purified by using glutathione sepharose chromatography.

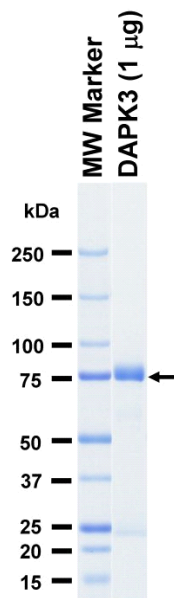
Storage buffer:

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.

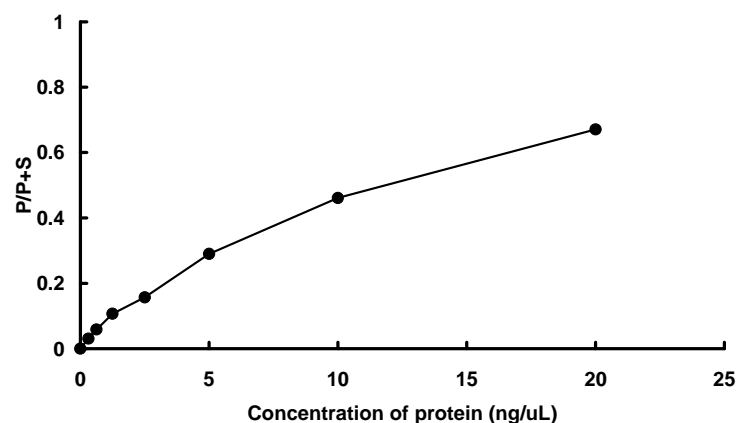
SDS-PAGE



Purity: 91 %

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 : ZIPTide

ATP : 1000 μ M

Product Information

DAPK3

 Product Number : **02-136**

Product description

Full-length human DAPK3 [1-454(end) amino acids of accession number NP_001339.1] was expressed as N-terminal GST-fusion protein (80 kDa) using baculovirus expression system. GST-DAPK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

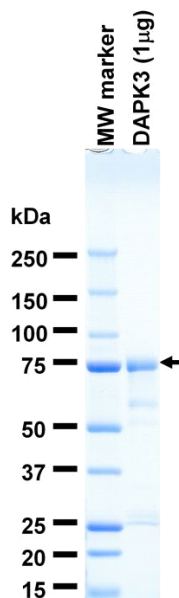
Storage buffer:

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.

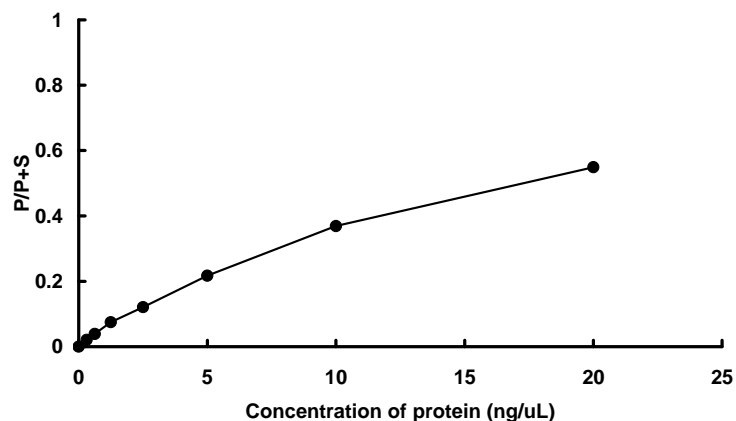
SDS-PAGE



Purity: 67 %

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 : ZIPtide

 ATP : 1000 μ M