

Product Information

TAOK2

Product Number : **07-133**

Product description

Human TAOK2, catalytic domain [1-319 amino acid of accession number NP_004774.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-TAOK2 was purified by using glutathione sepharose chromatography.

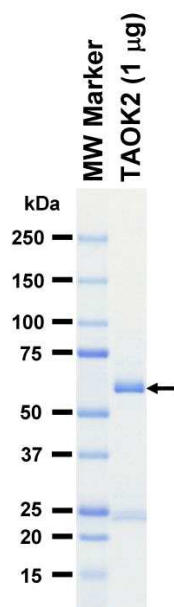
Storage buffer:

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.

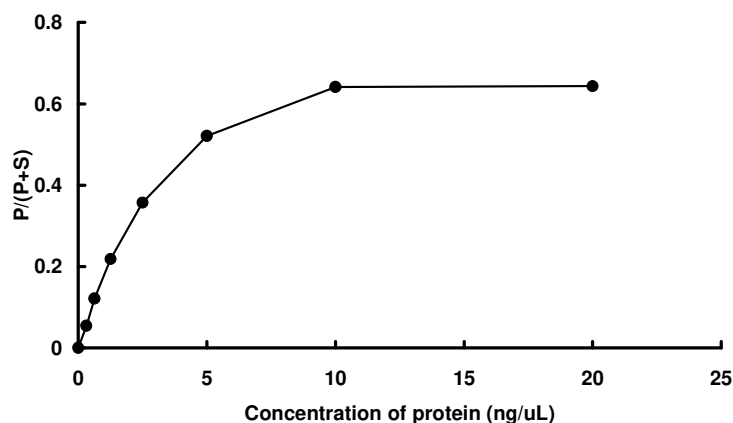
SDS-PAGE



Purity: 73 %

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 : TAOKtide

ATP : 100 µM