

Product Information

TAOK3

 Product Number : **07-135**

Product description

Full-length human TAOK3 [1-898(end) amino acids of accession number NP_057365.3] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system. GST-TAOK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

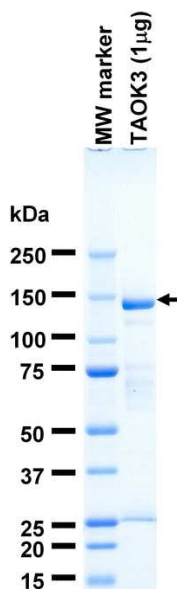
Storage buffer:

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.

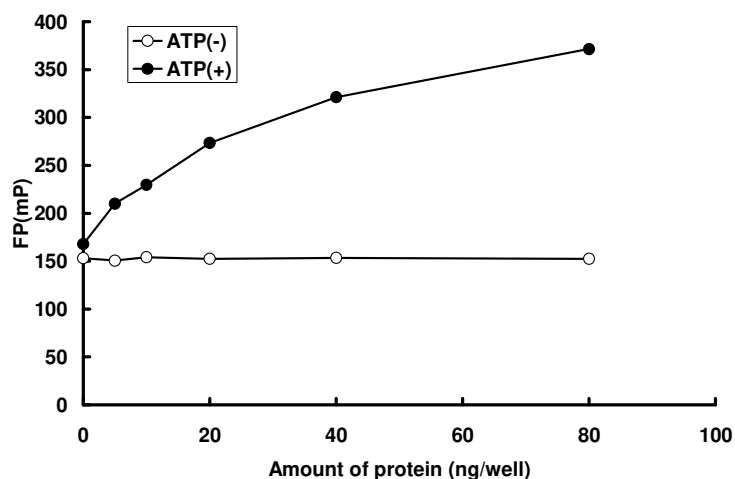
SDS-PAGE



Purity: 73 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : PKCε peptide

ATP : 100 μM