

Product Information

ZAK

Product Number : **09-120**

Product description

Full-length human ZAK [1-455(end) amino acids of accession number NP_598407.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-ZAK was purified by using glutathione sepharose chromatography and gel filtration chromatography.

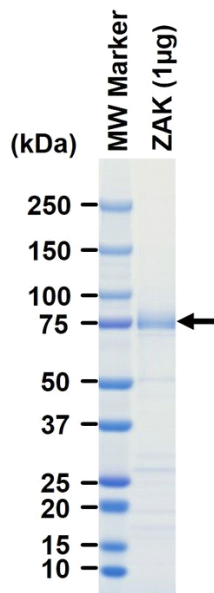
Storage buffer:

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.

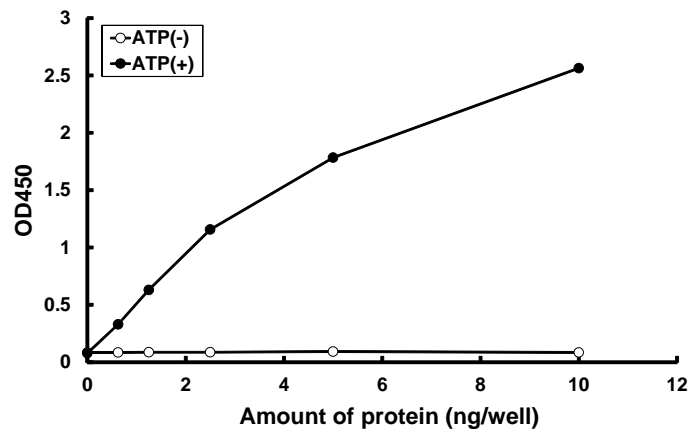
SDS-PAGE



Purity: 63 %

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

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated substrate protein/peptide, and after stopping kinase reaction by EDTA, the reaction solution was transferred into streptavidin-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG (or HRP-labeled anti-mouse IgG).

Substrate : BTN-MAP2K7[inactive]

ATP : 100 μ M