

Product Information

ULK3

Product Number : **05-173**

Product description

Full-length human ULK3 [1-472(end) amino acids of accession number NP_001092906.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-ULK3 was purified by using glutathione sepharose chromatography.

Storage buffer:

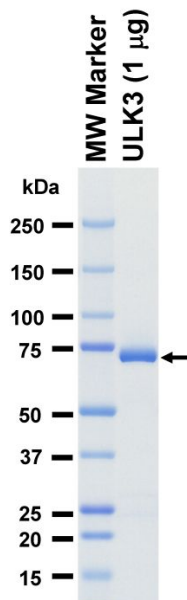
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

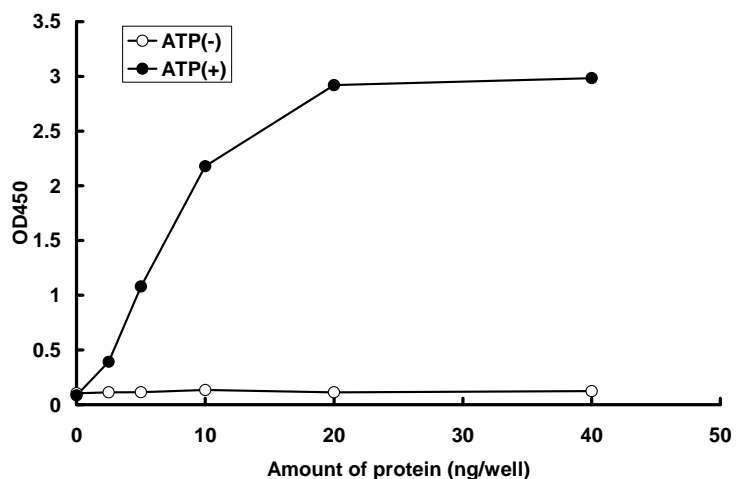
SDS-PAGE



Purity: 96 %

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, 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.

Substrate : MBP

ATP : 100 µM