

## Product Information

### ULK1(ATG1)(1-331aa)

Product Number : 05-136

#### Product description

Human ULK1, catalytic domain [1-331 amino acids of accession number NP\_003556.2] was expressed as N-terminal GST-fusion protein (64 kDa) using baculovirus expression system. GST-ULK1 was purified by using glutathione sepharose chromatography.

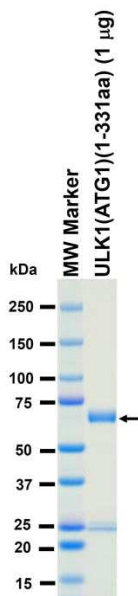
#### Storage buffer:

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.

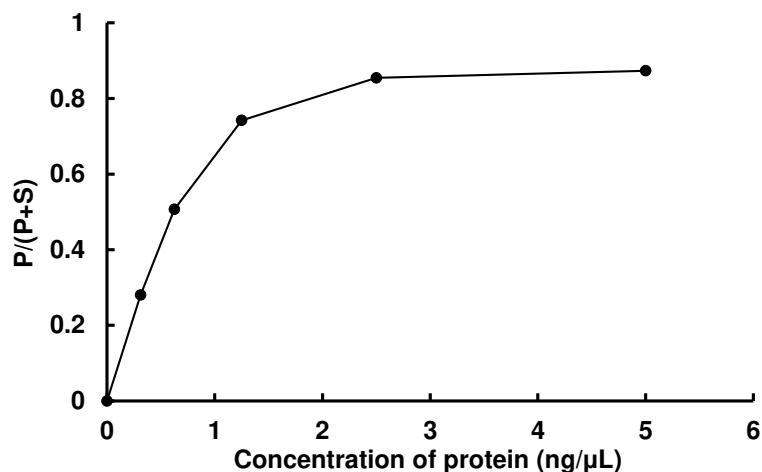
#### SDS-PAGE



Purity: 75 %

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 : CDC25c peptide

ATP : 100 μM