

## Product Information

### LZK(MAP3K13)

 Product Number : **09-114**

#### Product description

Full-length human LZK(MAP3K13) [1-966(end) amino acids of accession number NP\_004712.1] was expressed as N-terminal GST-fusion protein (136 kDa) using baculovirus expression system. GST-LZK(MAP3K13) was purified by using glutathione sepharose chromatography

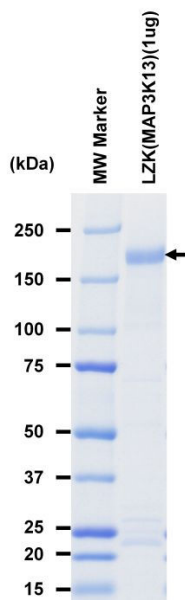
#### Storage buffer:

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.  
 Use within 12 months from the date of shipment.

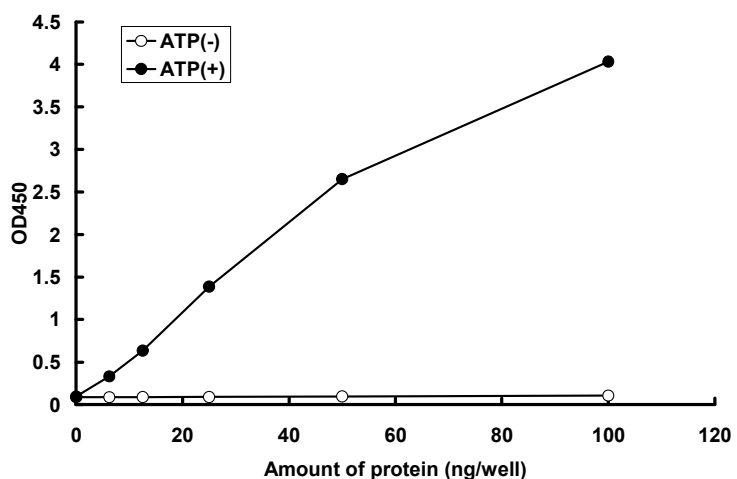
#### SDS-PAGE



Purity: 72 %

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

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with GST-fused substrate protein, and after stopping kinase reaction by EDTA, the reaction solution was transferred into glutathione-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG.

Substrate : MAP2K7 [inactive mutant]

 ATP : 100  $\mu$ M