

Product Information

LIMK2

 Product Number : **09-106**

Product description

Full-length human LIMK2 [1-638(end) amino acids of accession number NP_005560.1] was co-expressed as N-terminal GST-fusion protein (99 kDa) with human His-tagged ROCK2 [1-553 amino acids of accession number NP_004841.2] using baculovirus expression system. GST-LIMK2 was purified by using glutathione sepharose chromatography.

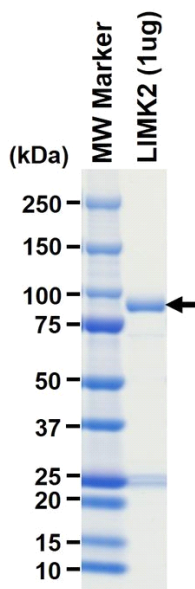
Storage buffer:

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.

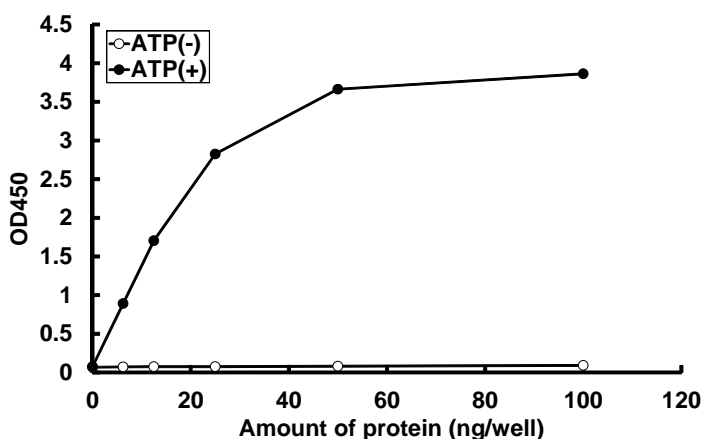
SDS-PAGE



Purity: 70 %

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(or HRP-labeled anti-mouse IgG) .

Substrate : Cofilin2

 ATP : 100 μ M