

Product Information

LIMK1

Product Number : **09-105**

Product description

Truncated human LIMK1 [285-638 amino acids of accession number NP_002305.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with human His-tagged ROCK2 [1-553 amino acids of accession number NP_004841.2] using baculovirus expression system. GST-LIMK1 was purified by using glutathione sepharose chromatography.

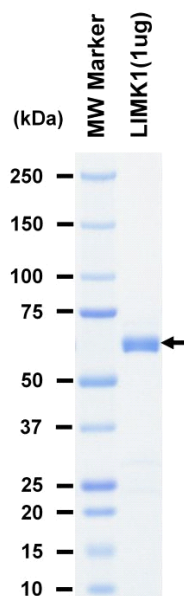
Storage buffer:

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.

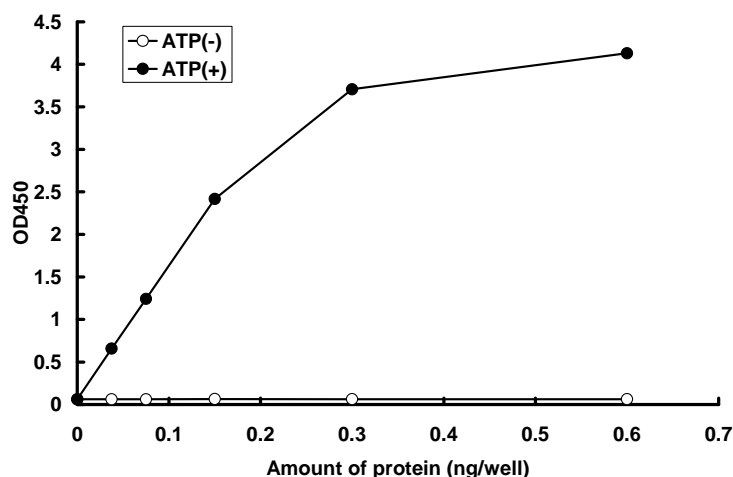
SDS-PAGE



Purity: 94 %

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