

Product Information

LKB1(STK11)/MO25α/STRADα

 Product Number : **02-119**

Product description

Full-length human LKB1 [1-433(end) amino acids of accession number NP_000446.1] , MO25α [1-341(end) amino acids of accession number NP_057373.1] and STRADα [1-431(end) amino acids of accession number NP_001003787.1] were co-expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-LKB1 was purified by using glutathione sepharose chromatography.

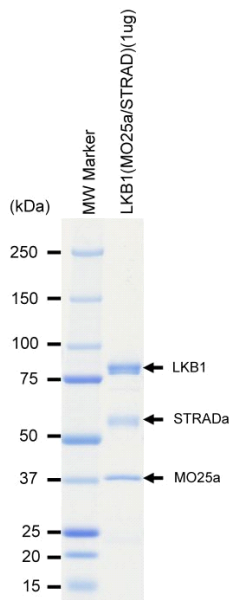
Storage buffer:

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.

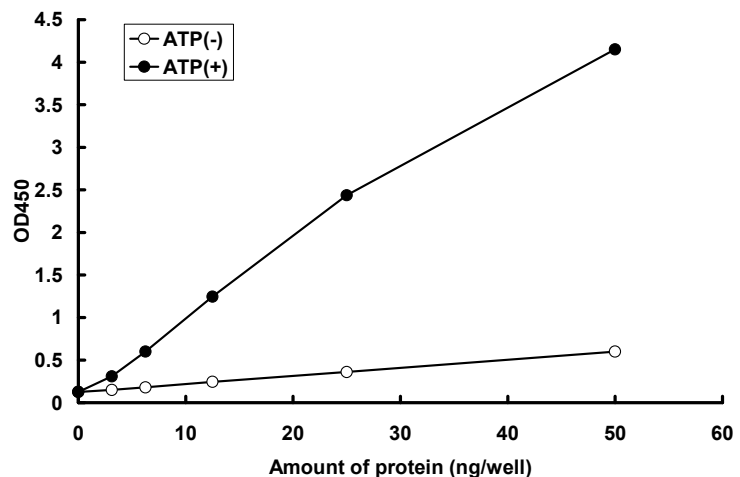
SDS-PAGE



Purity: 97 %

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/peptide, 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 : LKBtide

ATP : 100 μM