

Product Information

MYLK4

 Product Number : **02-161**

Product description

Full-length human MYLK4 [1-388(end) amino acids of accession number NP_001012418.2] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-MYLK4 was purified by using glutathione sepharose chromatography.

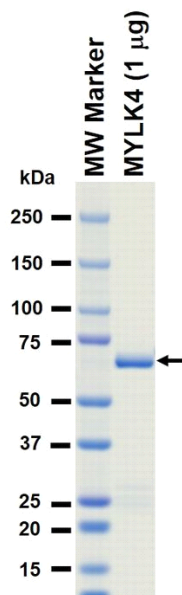
Storage buffer:

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.

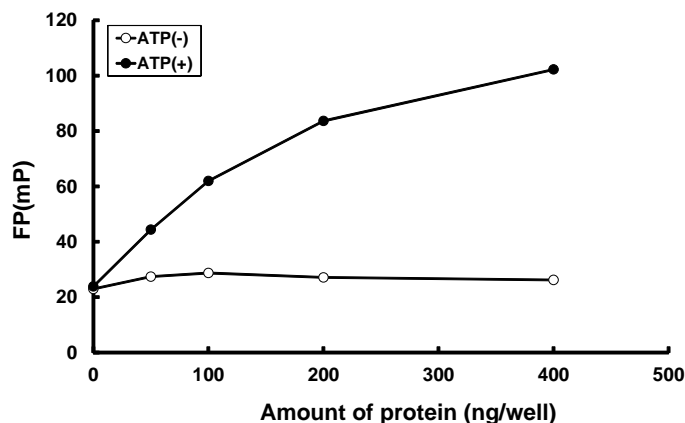
SDS-PAGE



Purity: 91 %

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

Activity data



The activity was determined by IMAP™ assay. The enzyme was incubated with fluorescein labeled peptide and phosphorylation was detected by IMAP™ technology (fluorescence polarization).

Substrate : MLCTide

 ATP : 100 μ M