

Product Information

BTN-MYLK4

 Product Number : **02-461-20N**

Product description

Full-length human MYLK4 [1-388(end) amino acids of accession number NP_001012418.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (49 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

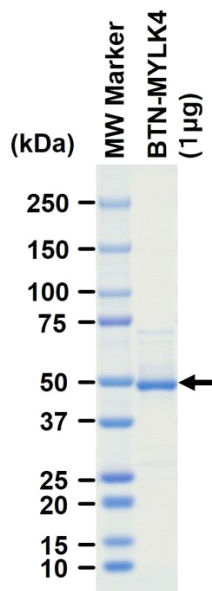
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

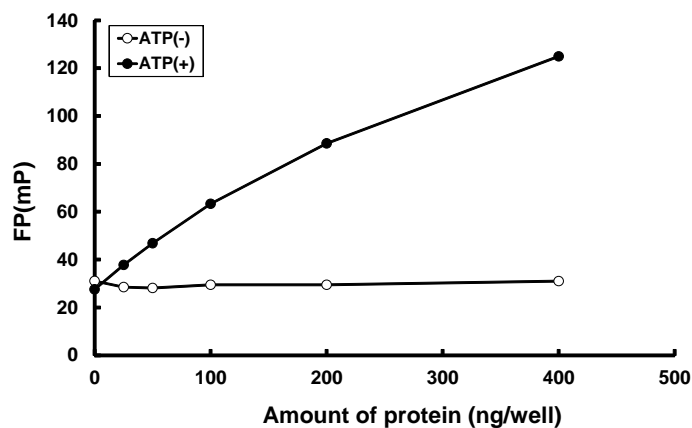
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 88 %
 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 µM