

## Product Information

### smMLCK(MYLK)

 Product Number : **02-151**

#### Product description

Human smMLCK, catalytic domain [1428-1771 amino acids of accession number NP\_444253.2] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-smMLCK was purified by using glutathione sepharose chromatography.

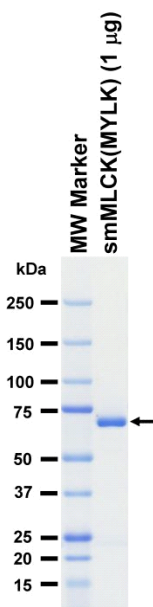
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
 1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

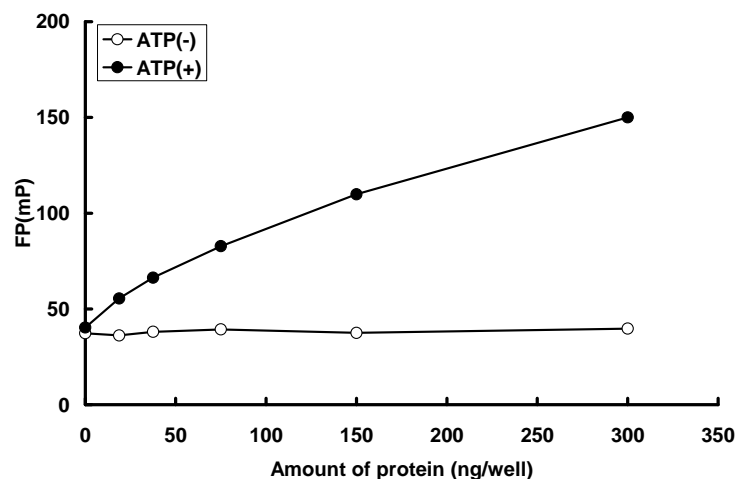
#### SDS-PAGE



Purity: 97 %

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

#### Activity data



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

Substrate : GS peptide

 ATP : 100  $\mu$ M