

## Product Information

### BTN-LATS2/MOBKL1A

 Product Number : **01-424-20N**

#### Product description

Human LATS2, catalytic domain [553-1088(end) amino acids of accession number NP\_055387.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (65 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP\_775739.1] and human GST-tagged MST2 [1-491(end) amino acids of accession number NP\_006272.2] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and glutathione sepharose chromatography.

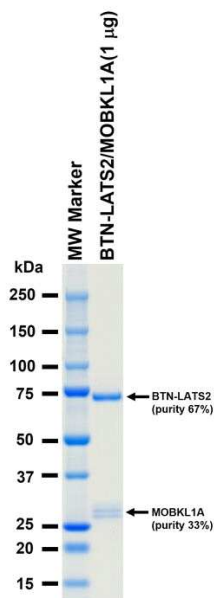
#### Storage buffer:

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.

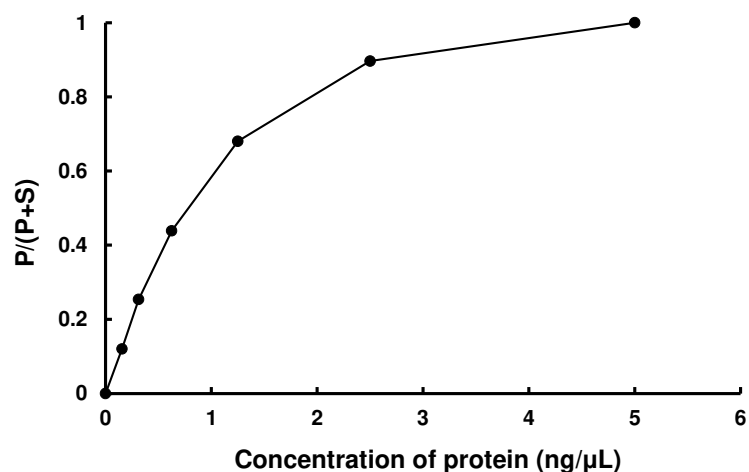
#### SDS-PAGE



Purity: 100 %

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

#### Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : SGKtide

ATP : 1000 µM