

Product Information

LATS1/MOBKL1A

 Product Number : **01-523**

Product description

Human LATS1, catalytic domain [589-1130(end) amino acids of accession number NP_004681.1] was co-expressed as N-terminal GST-fusion protein (90 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] and human BTN-MST2 [1-491(end) amino acids of accession number NP_006272.2] using baculovirus expression system. GST-LATS1/MOBKL1A was purified by using glutathione sepharose chromatography and DYKDDDDK tag antibody agarose.

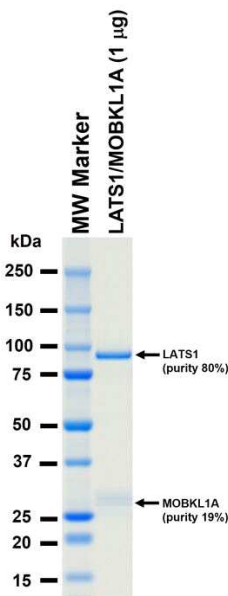
Storage buffer:

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

Storage and Handling:

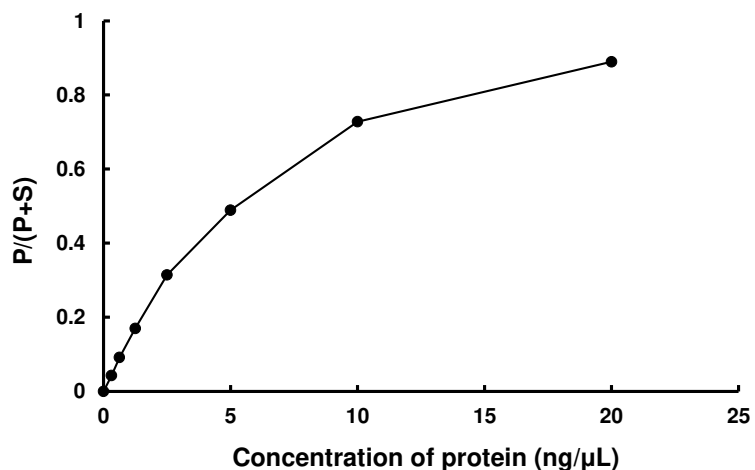
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 99 %
 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