
LATS1/MOBKL1A

Product code: 01-523

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.

Assay platform : MSA
Substrate : SGKtide

ATP (μ M) Km app / Bin : 23 / 25
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 4.2
IC50 at 1 mM ATP (nM) : 11
