
LATS2/MOBKL1A

Product code: 01-524

Human LATS2, catalytic domain [553-1088(end) amino acids of accession number NP_055387.2] was co-expressed as N-terminal GST-fusion protein (89 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 : 38 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.7
IC50 at 1 mM ATP (nM) : 4.0
