

---

# 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. GST-LATS2/MOBKL1A was purified by using glutathione sepharose chromatography and DYKDDDDK tag antibody agarose.

Assay platform	: MSA
Substrate	: SGKtide
ATP ( $\mu$ 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

