
LATS2

Product code 01-124

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] using baculovirus expression system. GST-LATS2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : SGKtide

ATP (μ M) Km app / Bin : 380 / 400

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 1.6

IC50 at 1 mM ATP (nM) : 2.2
