
MSK1(RPS6KA5)

Product code 01-147

Full-length human MSK1 [1-802(end) amino acids of accession number NP_004746.2] was co-expressed as N-terminal GST-fusion protein (117 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP_002736.3] using baculovirus expression system. GST-MSK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : Crosstide

ATP (μ M) Km app / Bin : 13 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 1.6

IC50 at 1 mM ATP (nM) : 30
