
MSK2(RPS6KA4)

Product code: 01-148

Full-length human MSK2 [1-772(end) amino acids of accession number NP_003933.1] was co-expressed as N-terminal GST-fusion protein (114 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP_002736.3] using baculovirus expression system. GST-MSK2 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Crosstide
ATP (μ M) Km app / Bin	: 40 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.3
IC50 at 1 mM ATP (nM)	: 20
