
MAP2K4

Product code: 07-144

Full-length human MAP2K4 [1-399(end) amino acids of accession number NP_003001.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] using baculovirus expression system. GST-MAP2K4 was purified by using glutathione sepharose chromatography.

Assay platform : MSA
Substrate : JNK2
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 4600

* JNK2/Modified Erktide
