
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
