
MAP2K7

Product code: 07-148

Full-length human MAP2K7 [1-419(end) amino acids of accession number NP_660186.1] was co-expressed as N-terminal GST-fusion protein (75 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] using baculovirus expression system. GST-MAP2K7 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) : 1100

* JNK2/Modified Erktide
