
MAP2K1

Product code: 07-141

Full-length human MAP2K1 [1-393(end) amino acids of accession number NP_002746.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged RAF1 [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] using baculovirus expression system. GST-MAP2K1 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

Assay platform : MSA
Substrate : Erk2
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 58

* Erk2/Modified Erktide
