
MAP2K6

Product code: 07-146

Full-length human MAP2K6 [2-334(end) amino acids of accession number NP_002749.2] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human His-tagged MLK3 [99-398 amino acids of accession number NP_002410.1] using baculovirus expression system.

Assay platform : MSA
Substrate : p38 α
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 140

* p38 α /Modified Erktide
