
MAP3K4

Product code: 07-106

Human MAP3K4, catalytic domain [1312-1608(end) amino acids of accession number NP_005913.3] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : MAP2K6
Cascade Assay *
ATP (μ M) : 1000
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 100

* MAP2K6/p38 α /Modified Erktide
