
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
