
MAP3K3

Product code: 07-105

Full-length human MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] was expressed as N-terminal GST-fusion protein (98 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)	: 72

* MAP2K6/p38 α /Modified Erktide
