
MAP3K5

Product code: 07-107

Human MAP3K5, catalytic domain [654-971 amino acids of accession number NP_005914.1] was expressed as N-terminal GST-tagged protein (62 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) : 14

* MAP2K6/p38 α /Modified Erktide
