
MAP4K2

Product code: 07-111

Full-length human MAP4K2 [1-820(end) amino acid of accession number NP_004570.2] was expressed as N-terminal GST-fusion protein (119 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : S6K2 peptide

ATP (μ M) Km app / Bin : 93 / 100
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 1.3
IC50 at 1 mM ATP (nM) : 0.87
