
MAP4K5(KHS1)

Product code: 07-113

Full-length human MAP4K5 [1-846(end) amino acids of accession number NP_006566.2] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP (μ M) Km app / Bin	: 28 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 3.1
