
HPK1(MAP4K1)

Product code: 07-410

Human HPK1, catalytic domain [1-346 amino acids of accession number NP_009112.1] was expressed as N-terminal DYKDDDDK tagged protein (41 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: S6K2 peptide
ATP (μ M) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 6.9
IC50 at 1 mM ATP (nM)	: 32

