
CHK1(CHEK1)

Product code: 02-117

Full-length human CHK1 [1-476(end) amino acids and I471V of accession number NP_001265.2] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 50 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.24
IC50 at 1 mM ATP (nM)	: 1.1

