
CHK2(CHEK2)

Product code: 02-162

Full-length human CHK2 [1-543(end) amino acids of accession number NP_009125.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	CHKtide
ATP (μ M) Km app / Bin	:	51 / 50
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at ATP Bin (nM)	:	11
IC50 at 1 mM ATP (nM)	:	25

