
CDC7/ASK

Product code: 05-109

Full-length human CDC7 [1-574(end) amino acids of accession number NP_003494.1] was co-expressed as N-terminal GST-fusion protein (92 kDa) with Dbf4(ASK) [1-674(end) amino acids of accession number NP_006707.1] using baculovirus expression system. GST-CDC7 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: MCM2 peptide
ATP (μ M) Km app / Bin	: 2.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: 1600
