
CDK7/CycH/MAT1

Product code: 04-108

Full-length human CDK7 [1-346(end) amino acids of accession number NP_001790.1] was co-expressed as N-terminal GST-fusion protein (66 kDa) with CyclinH [1-323(end) amino acids of accession number NP_001230.1] and MAT1 [1-309(end) amino acids of accession number NP_002422.1] using baculovirus expression system. GST-CDK7 was purified by using glutathione sepharose chromatography.

Assay platform : MSA
Substrate : CTD3 peptide
ATP (μ M) Km app / Bin : 32 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 17
IC50 at 1 mM ATP (nM) : 120
