
CDK2 / CycA2

Product code 04-103

Full-length human CDK2 [1-298(end) amino acids of accession number NP_001789.2] was co-expressed as N-terminal GST-tagged protein (61 kDa) with GST-CyclinA2 [1-432(end) amino acids of accession number NP_001228.1] using baculovirus expression system. GST-CDK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Modified Histone H1

ATP (μ M) K_{mapp} / B_{in} : 27 / 25
Metal : Mg
Reference compound : Staurosporine
IC₅₀ at ATP B_{in} (nM) : 1.0
IC₅₀ at 1 mM ATP (nM) : 7.1
