
CDK9/CycT1

Product code: 04-110

Full-length human CDK9 [1-372(end) amino acids of accession number NP_001252.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-CyclinT1 [1-726(end) amino acids of accession number NP_001231.2] using baculovirus expression system.

Assay platform : MSA
Substrate : CDK9 substrate
ATP (μ M) K_m app / B_{in} : 9.4 / 10
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP B_{in} (nM) : 5.2
IC50 at 1 mM ATP (nM) : 130
