
PLK1

Product code: 05-157

Full-length human PLK1 [1-603(end) amino acids of accession number NP_005021.2] was expressed as N-terminal GST-fusion protein (95 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : CDC25ctide

ATP (μ M) Km app / Bin : 5.6 / 5
Metal : Mg
Reference compound : GW843682X
IC50 at ATP Bin (nM) : 3.6
IC50 at 1 mM ATP (nM) : 47
