
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. GST-PLK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CDC25ctide

ATP (μ M) K_{mapp} / B_{in} : 5.6 / 5

Metal : Mg

Reference compound : GW843682X

IC50 at ATP B_{in} (nM) : 3.6

IC50 at 1 mM ATP (nM) : 47
