

---

# 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 ( $\mu$ 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

---