

---

# 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

---