

---

## ***PLK2***

Product code 05-158

Full-length human PLK2 [1-685(end) amino acids of accession number NP\_006613.2] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PLK2 was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: CHK2 peptide
ATP ( $\mu$ M) Kmapp / Bin	: 30 / 30
Metal	: Mg
Reference compound	: GW843682X
IC50 at ATP Bin (nM)	: 4.8
IC50 at 1 mM ATP (nM)	: n.a.

---