

---

---

# PLK3

Product code: 05-159

Human PLK3, catalytic domain [58-340 amino acids of accession number NP\_004064.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CDC25ctide

ATP ( $\mu$ M)  $K_m$  app /  $B_{in}$  : 6.8 / 5

Metal : Mg

Reference compound : GW843682X

IC50 at ATP  $B_{in}$  (nM) : 33

IC50 at 1 mM ATP (nM) : 450

---

---