

---

---

# CHK1(CHEK1)

Product code: 02-117

Full-length human CHK1 [1-476(end) amino acids and I471V of accession number NP\_001265.2] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CHKtide  
  
ATP ( $\mu$ M) Km app / Bin : 50 / 50  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.24  
IC50 at 1 mM ATP (nM) : 1.1

---

---