

---

## CK1γ2(CSNK1G2)

Product code: 03-106

Full-length human CK1γ2 [1-415(end) amino acids of accession number NP\_001310.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : CKtide

ATP (μM) Km app / Bin : 10 / 10

Metal : Mg

Reference compound : 5-Iodotubercidin

IC50 at ATP Bin (nM) : 510

IC50 at 1 mM ATP (nM) : >10000

---