

---

# CK1γ1(CSNK1G1)

Product code 03-105

Full-length human CK1γ1 [1-422(end) amino acids of accession number NP\_071331.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-CK1γ1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : CKtide

ATP (μM) Km app / Bin : 6.3 / 5  
Metal : Mg  
Reference compound : 5-Iodotubercidin  
IC50 at ATP Bin (nM) : 1300  
IC50 at 1 mM ATP (nM) : n.a.

---