

---

## CK1γ3(CSNK1G3)

Product code: 03-107

Full-length human CK1γ3 [1-447(end) amino acids of accession number NP\_004375.2] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : CKtide

ATP (μM) Km app / Bin : 3.2 / 5  
Metal : Mg  
Reference compound : 5-Iodotubercidin  
IC50 at ATP Bin (nM) : 920  
IC50 at 1 mM ATP (nM) : n.a.

---