
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. GST-CK1γ2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
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) : n.a.
