
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. GST-CK1γ3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
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.
