
CK1 γ 2(CSNK1G2)

Product code: 03-106

Full-length human CK1 γ 2 [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 (μ M) Km app / Bin	: 10 / 10
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 510
IC50 at 1 mM ATP (nM)	: >10000
