
CK1 γ 1(CSNK1G1)

Product code: 03-105

Full-length human CK1 γ 1 [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.

Assay platform	: MSA
Substrate	: CKtide
ATP (μ M) 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.
