
CK1 ϵ (CSNK1E)

Product code: 03-104

Human CK1 ϵ , catalytic domain [1-348 amino acids of accession number NP_001885.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CKtide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 300
IC50 at 1 mM ATP (nM)	: 5800
