
CK1δ(CSNK1D)

Product code: 03-103

Human CK1δ, catalytic domain [1-294 amino acids of accession number NP_001884.2] was expressed as N-terminal GST-fusion protein (61 kDa) using E. coli expression system.

Assay platform : MSA
Substrate : CKtide

ATP (μM) Km app / Bin : 7.7 / 10
Metal : Mg
Reference compound : 5-Iodotubercidin
IC50 at ATP Bin (nM) : 25
IC50 at 1 mM ATP (nM) : 570
