
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. GST-CK1δ was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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
