

---

## ***CK1ε(CSNK1E)***

Product code 03-104

Human CK1ε, 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. GST-CK1ε was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CKtide

ATP (μM) Kmapp / Bin : 16 / 25

Metal : Mg

Reference compound : 5-Iodotubercidin

IC50 at ATP Bin (nM) : 300

IC50 at 1 mM ATP (nM) : 5800

---