

---

---

## CK1 $\alpha$ (CSNK1A1)

Product code: 03-101

Full-length human CK1 $\alpha$  [1-337(end) amino acids of accession number NP\_001883.4] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CKtide
ATP ( $\mu$ M) Km app / Bin	: 4.1 / 5
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 150
IC50 at 1 mM ATP (nM)	: >10000

---

---