
CK2 α 1/ β (CSNK2A1/B)

Product code: 05-184

Full-length human CK2 α 1 [1-391(end) amino acids of accession number NP_001886.1] was co-expressed as N-terminal GST-fusion protein (72 kDa) with human His-tagged CK2 β [1-215 amino acids of accession number NP_001311.3] using baculovirus expression system.

Assay platform	: MSA
Substrate	: CK2tide
ATP (μ M) Km app / Bin	: 2.9 / 5
Metal	: Mg
Reference compound	: TBB
IC50 at ATP Bin (nM)	: 60
IC50 at 1 mM ATP (nM)	: 4800

