
CK2 α 2/ β (CSNK2A2/B)

Product code: 05-185

Full-length human CK2 α 2 [1-350(end) amino acids of accession number NP_001887.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with human His-tagged CK2 β [1-215 amino acids of accession number NP_001311.3] using baculovirus expression system. GST-CK2 α 2 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: CK2tide
ATP (μ M) Km app / Bin	: 2.1 / 5
Metal	: Mg
Reference compound	: TBB
IC50 at ATP Bin (nM)	: 50
IC50 at 1 mM ATP (nM)	: 5800
