

---

## CK2 $\alpha$ 2/ $\beta$ (CSNK2A2/B)

Product code 05-185

Full-length human CK2 $\alpha$ 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 $\beta$  [1-215 amino acids of accession number NP\_001311.3] using baculovirus expression system. GST-CK2 $\alpha$ 2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : CK2tide

ATP ( $\mu$ M) Km app / Bin : 2.1 / 5  
Metal : Mg  
Reference compound : TBB  
IC50 at ATP Bin (nM) : 50  
IC50 at 1 mM ATP (nM) : n.a.

---