

---

---

## ***CK2 $\alpha$ 1* / *$\beta$* (CSNK2A1 / B)**

Product code 05-184

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

Assay platform : Mobility Shift Assay

Substrate : CK2tide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 2.9 / 5

Metal : Mg

Reference compound : TBB

IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 60

IC<sub>50</sub> at 1 mM ATP (nM) : 4800

---

---