

---

---

## 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              | : MSA     |
| Substrate                   | : CK2tide |
| ATP ( $\mu$ 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    |

