
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.

| | |
|-----------------------------|-----------|
| 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) | : n.a. |
