

---

## CaMK2α(CAMK2A)

Product code 02-109

Full-length human CaMK2α [1-478(end) amino acids of accession number NP\_741960.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CaMK2α was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : GS peptide

ATP (μM) Km app / Bin : 33 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.75

IC50 at 1 mM ATP (nM) : 3.2

---