

---

---

## CaMK2 $\beta$ (CAMK2B)

Product code: 02-110

Full-length human CaMK2 $\beta$  [1-503 amino acids of accession number NP\_742078.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP ( $\mu$ M) Km app / Bin	: 19 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.54
IC50 at 1 mM ATP (nM)	: 5.8

---

---