
CaMK2β(CAMK2B)

Product code: 02-110

Full-length human CaMK2β [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 (μM) 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
