
CaMK2 β (CAMK2B)

Product code 02-110

Full-length human CaMK2 β [1-503 amino acids of accession number NP_742078.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-CaMK2 β was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : GS peptide

ATP (μ 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
