
CaMK2γ(CAMK2G)

Product code 02-112

Full-length human CaMK2γ[1-518(end) amino acids of accession number NP_751910.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-CaMK2γ was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : GS peptide

ATP (μM) Km app / Bin : 23 / 25

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.39

IC50 at 1 mM ATP (nM) : 1.1
