
CaMK2δ(CAMK2D)

Product code 02-111

Full-length human CaMK2δ [1-478 amino acids of accession number NP_742113.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CaMK2δ was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : GS peptide

ATP (μM) Km app / Bin : 6.3 / 5

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.26

IC50 at 1 mM ATP (nM) : 0.64
