
CaMK2 δ (CAMK2D)

Product code: 02-111

Full-length human CaMK2 δ [1-478 amino acids of accession number NP_742113.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 6.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.26
IC50 at 1 mM ATP (nM)	: 3.7
