
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.

Assay platform	: MSA
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
