
CaMK2 α (CAMK2A)

Product code: 02-109

Full-length human CaMK2 α [1-478(end) amino acids of accession number NP_741960.1] 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	: 33 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.75
IC50 at 1 mM ATP (nM)	: 3.2
