
CaMK4

Product code: 02-108

Full-length human CaMK4 [1-473(end) amino acids of accession number NP_001735.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 140
IC50 at 1 mM ATP (nM)	: 1000
