
CaMK1 α (CAMK1)

Product code: 02-104

Full-length human CaMK1 α [1-370(end) amino acids of accession number NP_003647.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform : MSA

Substrate : GS peptide

ATP (μ M) Km app / Bin : 750 / 1000

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 16

IC50 at 1 mM ATP (nM) : 16
