
PKC α (PRKCA)

Product code: 01-133

Full-length human PKC α [1-672(end) amino acids of accession number NP_002728.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system.

Assay platform : MSA
Substrate : PKC peptide

ATP (μ M) Km app / Bin : 36 / 50
Metal : Mg+Ca
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 0.33
IC50 at 1 mM ATP (nM) : 3.6
