
PKC α (PRKCA)

Product code: 01-133

Full-length human PKC α [1-672(end) amino acids and V568I of accession number NP_002728.2] 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
