
PKC γ (PRKCG)

Product code: 01-137

Full-length human PKC γ [1-697(end) amino acids of accession number NP_002730.1] was expressed as N-terminal GST-fusion protein (106 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 74 / 75
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 11
