
PKC θ (PRKCQ)

Product code: 01-140

Full-length human PKC θ [1-706(end) amino acids and P330L of accession number NP_006248.1] was expressed as N-terminal GST-fusion protein (109 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 18 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.72
IC50 at 1 mM ATP (nM)	: n.a.

