
PKCζ(PRKCZ)

Product code: 01-141

Full-length human PKCζ [1-592(end) amino acids of accession number NP_002735.3] was expressed as N-terminal GST-fusion protein (94kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP (μM) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 55
IC50 at 1 mM ATP (nM)	: n.a.
