
PKD3(PRKD3)

Product code: 02-159

Full-length human PKD3 [1-890(end) amino acids of accession number NP_005804.1] was expressed as N-terminal GST-fusion protein (127 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.80
IC50 at 1 mM ATP (nM)	: n.a.

