
PKD1(PRKD1)

Product code: 02-157

Full-length human PKD1 [1-912(end) amino acids of accession number NP_002733.2] was expressed as N-terminal GST-fusion protein (129 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: n.a.

