
PKD1(PRKD1)

Product code 02-157

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

Assay platform : Mobility Shift Assay

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.
