

---

## PKD2(PRKD2)

Product code: 02-158

Full-length human PKD2 [1-878(end) amino acids of accession number NP\_057541.2] was expressed as N-terminal GST-fusion protein (124 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: GS peptide
ATP ( $\mu$ M) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 16

---