

---

---

# PKN1

Product code: 01-144

Full-length human PKN1 [1-942(end) amino acids of accession number NP\_002732.3] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system.

Assay platform	: IMAP
Substrate	: S6K peptide
ATP ( $\mu$ M) Km app / Bin	: 19 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.15
IC50 at 1 mM ATP (nM)	: n.a.

