

---

---

# PKC $\beta$ 1(PRKCB1)

Product code: 01-134

Full-length human PKC $\beta$ 1 [1-671(end) amino acids of accession number NP\_997700.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP ( $\mu$ M) Km app / Bin	: 79 / 75
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.71
IC50 at 1 mM ATP (nM)	: n.a.

