

---

---

## PKC $\beta$ 2(PRKCB2)

Product code: 01-165

Full-length human PKC $\beta$ 2 [1-673(end) amino acids of accession number NP\_002729.2] 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	: 41 / 50
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.43
IC50 at 1 mM ATP (nM)	: n.a.

