

---

---

# PKAC $\beta$ (PRKACB)

Product code: 01-128

Full-length human PKAC $\beta$  [1-351(end) amino acids of accession number NP\_002722.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	Kemptide
ATP ( $\mu$ M) Km app / Bin	:	4.7 / 5
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at ATP Bin (nM)	:	1.0
IC50 at 1 mM ATP (nM)	:	37

