

---

---

# 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

---

---