

---

## PKC $\alpha$ (PRKCA)

Product code 01-133

Full-length human PKC $\alpha$  [1-672(end) amino acids of accession number NP\_002728.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC $\alpha$  was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : PKC peptide

ATP ( $\mu$ M) Km app / Bin : 36 / 50

Metal : Mg+Ca

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.33

IC50 at 1 mM ATP (nM) : 3.6

---