

---

## ***PKC $\epsilon$ (PRKCE)***

Product code 01-136

Full-length human PKC $\epsilon$  [1-737(end) amino acids of accession number NP\_005391.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-PKC $\epsilon$  was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay  
Substrate : PKC peptide

ATP ( $\mu$ M)  $K_{mapp}$  /  $B_{in}$  : 16 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC<sub>50</sub> at ATP  $B_{in}$  (nM) : 0.56  
IC<sub>50</sub> at 1 mM ATP (nM) : 5.6

---