

---

---

## 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.

Assay platform	: MSA
Substrate	: PKC peptide
ATP ( $\mu$ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.37
IC50 at 1 mM ATP (nM)	: 1.8

