

---

---

## 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           |

---

---