

---

---

## PKAC $\alpha$ (PRKACA)

Product code: 01-127

Full-length human PKAC $\alpha$  [1-351(end) amino acids of accession number NP\_002721.1] 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 | : 2.6 / 5       |
| Metal                       | : Mg            |
| Reference compound          | : Staurosporine |
| IC50 at ATP Bin (nM)        | : 0.80          |
| IC50 at 1 mM ATP (nM)       | : 86            |

---

---