

---

---

## PKAC $\gamma$ (PRKACG)

Product code: 01-129

Full-length human PKAC $\gamma$  [1-351(end) amino acids of accession number NP\_002723.2] 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	: 4.5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: n.a.

---

---