

---

---

## PKC $\eta$ (PRKCH)

Product code: 01-138

Full-length human PKC $\eta$  [1-683(end) amino acids of accession number NP\_006246.2] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: PKC peptide
ATP ( $\mu$ M) Km app / Bin	: 42 / 50
Metal	: Mg+Zn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.78
IC50 at 1 mM ATP (nM)	: n.a.

---

---