

---

---

## CaMK1 $\delta$ (CAMK1D)

Product code: 02-106

Full-length human CaMK1 $\delta$ [1-357(end) amino acid of accession number NP\_065130.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: Synapsin peptide
ATP ( $\mu$ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.5
IC50 at 1 mM ATP (nM)	: n.a.

---

---