

---

## CaMK2δ(CAMK2D)

Product code: 02-111

Full-length human CaMK2δ [1-478 amino acids of accession number NP\_742113.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : GS peptide

ATP (μM) Km app / Bin : 6.3 / 5  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 0.26  
IC50 at 1 mM ATP (nM) : 0.64

---