

---

# CaMK4

Product code: 02-108

Full-length human CaMK4 [1-473(end) amino acids of accession number NP\_001735.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : GS peptide

ATP ( $\mu$ M) Km app / Bin : 20 / 25  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 140  
IC50 at 1 mM ATP (nM) : 1000

---