

---

# 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. GST-CaMK4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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

---