

## Data Sheet

### CAMK2δ

Human recombinant, with N-terminal GST-tag

**Catalog #:** 41109

**Lot #:** 141211

**Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

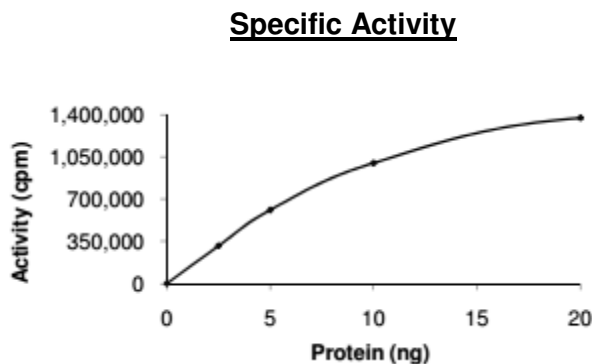
1. Hoch, B., *et al.*, *Circ. Res.* 1999; **84**: 713-721.
2. Zhu, W.-Z., *et al.*, *J. Clin. Invest.* 2003; **111**: 617-625.

**Description:** Recombinant human calcium/calmodulin-dependent protein kinase 2 delta (CAMK2δ), also known as CAMKIIδ, CAMK2d and CAMKIIδ, GenBank Accession No. NM\_172115, full-length with an N-terminal GST tag and expressed in a baculovirus-infected Sf9 insect cell expression system. MW = 77 kDa.

**Specific Activity:** 5,860 pmol/min/μg.  
 Assay conditions: Varying amounts of enzyme were incubated in 5 mM MOPS, pH 7.2, 2.5 mM β-glycerophosphate, 4 mM MgCl<sub>2</sub>, 2.5 mM MnCl<sub>2</sub>, 1 mM EGTA, 0.4 mM EDTA, 0.01 mM DTT, 0.05 mM ATP, 7.5 μg of Autocamide 2 synthetic peptide substrate, Ca<sup>2+</sup>/Calmodulin solution, and 32 μCi of [<sup>33</sup>P]-ATP was incubated at 30°C for 15 minutes. Then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

**Application:** Useful for studying enzyme kinetics, screening inhibitors, and selectivity profiling.

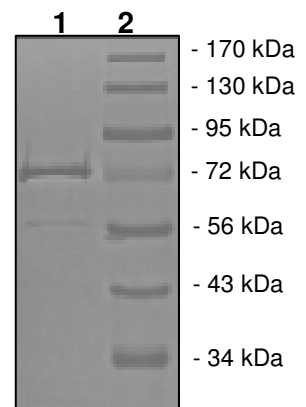
## Quality Assurance



**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:** CAMK2δ  
**Lane 2:** Protein Marker

**MW:** 77 kDa  
**Purity:** ≥85%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)