

Data Sheet

CK1 δ GST -Tag

Human, Recombinant, with N terminal
GST-tag

Catalog #: 79347

Lot #: 171023 **Conc.:** 0.1 mg/ml

Description: Full length Recombinant CK1 δ also known as CSNK1D, HCKID, and CK1D. Genbank Accession NM_001893 with N-terminal GST-tag, expressed in Sf9 cells via a baculovirus expression system. MW = 72kDa.

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

Applications: Useful for the kinase assay kits and western blotting.

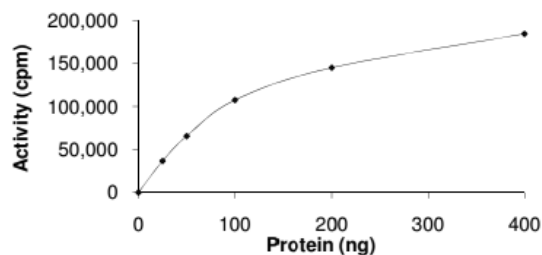
Stability: At least 6 months at -80°C .
Avoid freeze/thaw cycles.

References: (add references)

1. von Blume, J., *et al.*, *EMBO J.* 2007;**26(22)**:4619-33.
2. Smal, C., *et al.*, *Nature.* 2010;**502(1)**:44-52.

Quality Assurance

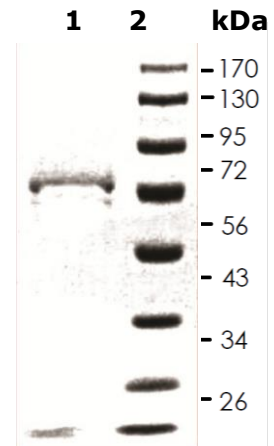
Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
CK1
Lane 2:
Protein Marker

MW: 72 kDa
Purity: $\geq 80\%$



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
Please visit our website at: www.bpsbioscience.com