

Data Sheet

CK1 α 1, GST-Tag

Human, Recombinant, with N-terminal
GST-tag

Catalog #: 79346

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

Description: Full length recombinant human CK1 α 1, also known as CSNK1A1, CK1, HLCDGP1, PRO2975, and CK1A1. Genbank Accession NM_001892 with an N-terminal GST-tag, expressed in Sf9 cells via a baculovirus expression system. MW = 62kDa.

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

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

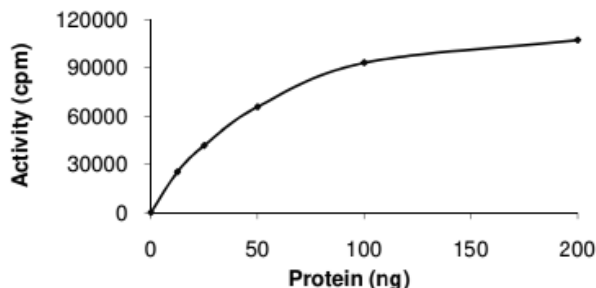
Stability: At least 6 months at -80°C .
Avoid freeze/thaw cycles.

References: (add references)

1. Jia, J., *et al.*, *Nature* 2004;**432**:1045-1050.
2. Huart, A.S., *et al.*, *J Biol Chem.* 2009;**475**:284(47):32384-94.

Quality Assurance

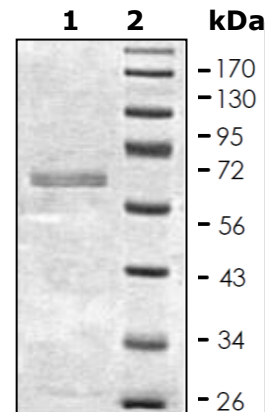
Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
CK1 α 1
Lane 2:
Protein Marker

MW: 62 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