

Data Sheet

CDK4, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100052

Lot #: 180523

Conc.: 0.56 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 20% glycerol,
3mM DTT, 100ug/ml flag peptide

Stability: At least 6 months at -80°C .
Avoid freeze/thaw cycles. Storing diluted
enzyme is not recommended, if necessary,
use carrier protein (BSA 0.1 – 0.5%).

References:

1. Bullrich, F. *et al.*, *Cancer Res.* (1995);
55:1199-1205.

2. Ye, X. *et al.*, *Proc. Nat. Acad. Sci.*
(2001); **98**:1682-1686.

Description:

Human CDK4, also known as cyclin-dependent kinase 4, PSK-J3, and cell division protein kinase 4, with N-terminal FLAG-tag, a.a 2-303(end), GenBank Accession No. NM_000075, expressed in a Baculovirus-infected Sf9 cell expression system.

Quality Assurance

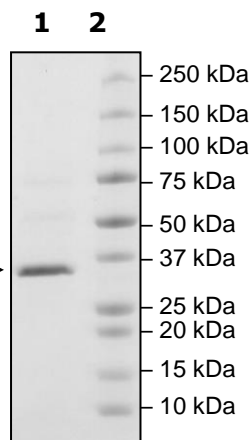
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
Cdk4

Lane 2:
Protein Marker

Purity: $\geq 90\%$ →

MW:
CDK4: 36 kDa



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