

## Data Sheet

### CDK19/Cyclin C, GST-tags

Human, Recombinant, N-terminal GST-tags

**Catalog #:** 100593

**Lot #:** 191216

**Conc.:** 0.1 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Lahti, J. M. *et al*, *Nature Genet.* **7**: 370-375, 1994.
2. Bunnell, B. *et al*, *Proc. Nat. Acad. Sci.* **87**: 7467-7471, 199

**Description:** CDK19/Cyclin C complex of human CDK19, also known as CDK11, CDC2L6, CDK8-like, and KIAA1028, GenBank Accession No. NM\_015076, a.a. 1-502(end) with N-terminal GST-tag, MW= 92 kDa, and human Cyclin C, also known as CCNC and CycC, GenBank Accession No. NM\_005190, a.a. 1-238(end) with N-terminal GST-tag. MW= 55 kDa. Co-expressed in a Baculovirus infected Sf9 cell expression system.

## Quality Assurance

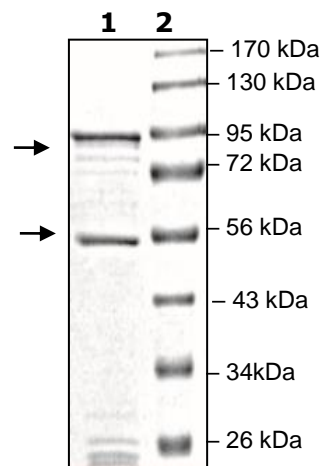
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
CDK19/Cyclin C

**Lane 2:**  
Protein Marker

**Purity:**  $\geq 90\%$

**MW:**  
CDK19: 92 kDa  
CyclinC: 55 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** 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)