

## Data Sheet

### CDC7/DBF4 Complex

Human, Recombinant, N-terminal GST-tags

**Catalog #:** 100538

**Lot #:** 191024

**Conc.:** 0.1 mg/ml

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

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

**Reference:**

1. Hess, G. F. *et. al. Gene* **211**: 133-140, 1998.
2. Sheu, Y.-J *et. al. Nature* **463**: 113-117, 2010

**Description:** CDC7/DBF4 complex of human **CDC7**, also known as HsCDC7 and Hsk 1, GenBank Accession No. NM\_003503 a.a. 1-574(end) with N-terminal GST-tag and MW= 94 kDa, and human **DBF4**, also known as ASK and CHIF, GenBank Accession No. NM\_006716, a.a. 1-674(end) with N-terminal GST-tag and MW= 125 kDa, co-expressed expressed in an Sf9 infected baculovirus expression system.

**Specific Activity:**  $\geq 8$  pmol/min/ $\mu\text{g}$

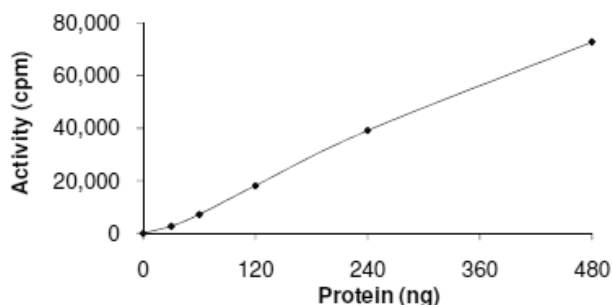
**Assay Conditions:** 5 mM MOPS, 2.5 mM

$\beta$ -glycerophosphate, 5 mM  $\text{MgCl}_2$ , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 50 ng/ml BSA, 2 mM ATP, 200  $\mu\text{g}$  of PDKtide synthetic peptide substrate. Add [ $^{32}\text{P}$ ]-ATP and incubate at  $30^{\circ}\text{C}$  for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

**Applications:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

**Specific Activity**



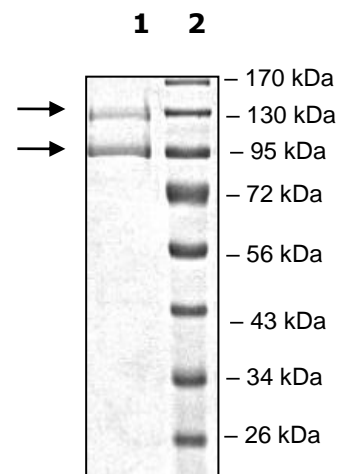
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
CDC7/DBF4

**Lane 2:**  
Protein Marker

**MW:** CDC7 94 kDa  
DBF4 125 kDa

**Purity:**  $\geq 85\%$



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

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