

## Data Sheet

### PKN1 (PRK1), GST-Tag

Human, Recombinant, N-terminal GST-Tag

**Catalog #:** 79678

**Lot #:** 181205

**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, 25% glycerol

**Stability:** At least 6 months at  $-70^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### References:

1. Amano, M. *et al.*, *Science*. 1996; **271**: 648-651.
2. Metzger, E. *et al.*, *EMBO J*. 2003; **22**: 270-280.

#### Description:

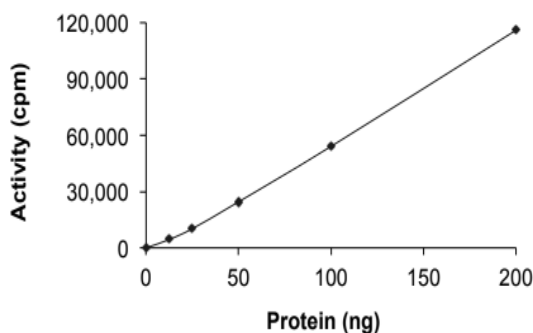
Human PKN1 protein, also known as PRK1, DBK, and PKN, GenBank Accession No. BC040061, a.a. 289-end, with N-terminal GST-Tag, expressed in an Sf9 insect cell using a baculovirus expression system. MW = 115 kDa.

**Specific Activity:** 30 pmol/min/ $\mu\text{g}$

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

### Quality Assurance

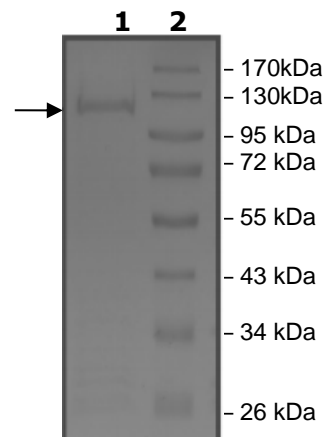
#### Specific Activity



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

**Lane 1:**  
PKN1/PRK1  
**Lane 2:**  
Protein Marker

**MW:** 115 kDa  
**Purity:**  $\geq 90\%$



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)