

## Data Sheet

### Plk3 Polo Box Domain

Human, recombinant, N-terminal MBP tag and C-terminal His tag

Catalog #: 50302

Lot #: 120403MN      Conc.: 0.72 mg/ml

**Formulated in:** 50mM Tris-HCL 8.0, 138mM NaCl, 2.7mM KCL, 10% glycerol, 250mM Imidazole, and 3mM DTT.

**Stability:** >6 months at -80°C

**References:**

1. Li,Z., et al., J. Biol. Chem. **280** (17), 16843-16850 (2005)
2. Xie,S., et al., Oncogene **23** (21), 3822-3829 (2004)

**Description:** Human N-terminal MBP tagged Plk3 polo box domain (GenBank Accession No.NM\_004073/NP\_004064), (a.a. 335-646) with C-terminal His tag, MW=80.5 kDa, expressed in an E.coli expression system.

**Application:**

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

### Quality Assurance

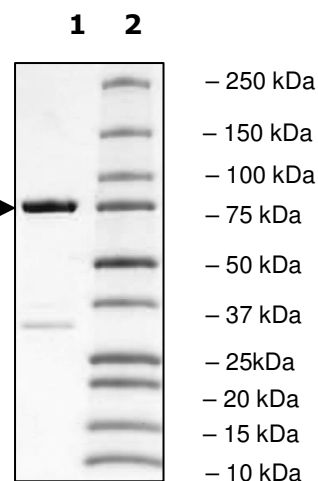
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2.0 µg PLK3

**Lane 2:** →  
Protein Marker  
(BioRad 161-0374)

**MW:** 80.5 kDa

**Purity:** ≥90%



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

To place your order, please contact us by Phone **1.858.829.3082** 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)