

Data Sheet

Aurora B Protein Kinase

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

Catalog #: 40002

Lot#: 120522-E250B **Conc.:** 0.11 mg/ml

Formulated in: 40mM Tris pH8.0, 138 mM NaCl, 2.2 mM KCl, 220 mM Imidazole 3mM DTT, and 10% Glycerol.

Stability: >6 months at -80°C

References:

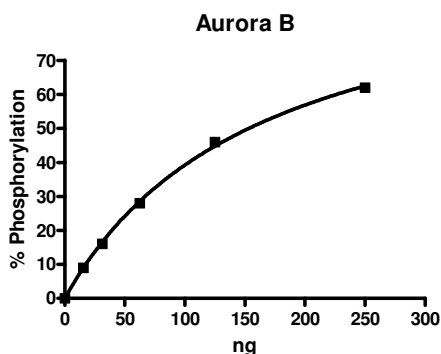
1. Delaval, B. *et al.*, *Oncogene* **23** (26), 4516-4522 (2004)
2. Yasui, Y. *et al.*, *J. Biol. Chem.* **279**, 12997-13003 (2004)

Description: Human Aurora kinase B (GenBank Accession No. NM_004217), (a.a. 2-344), with N-terminal GST tag and C-terminal His tag, MW= 66kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 6.7 pmole/min/ug.
Assay buffer: 50 mM HEPES, pH 7.4, 3 mM MgCl₂, 3 mM MnCl₂, 1 mM DTT, 3 uM Na-orthovanadate, 0.5 mM ATP, 4 uM aurora B substrate, and 4 ug/ml Aurora B kinase.

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

Quality Assurance

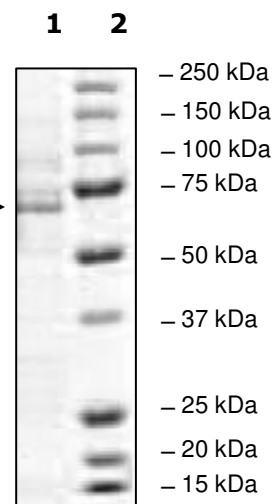


10% SDS-PAGE Coomassie staining

Lane 1:
1.5 ug Aurora B

Lane 2: →
Protein Marker

MW: 66 kDa
Purity: ≥50%



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