



Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

### For research use only

Not intended or approved for  
diagnostic or therapeutic use.

## Aurora Kinase B, active

(Recombinant human protein expressed in Sf9 cells)

Catalog # **E-K057**  
Size: 10 µg  
Specific activity: 22 U/mg

### Product Description

Aurora-B is a serine/threonine protein kinase. Recombinant human Aurora kinase B (AURKB, amino acids 2-344) with N-terminal GST-6xHis tag was expressed in a baculovirus infected Sf9 insect cell expression system, purified using Ni-NTA beads. The GenBank Accession number is NM\_004217. MW = 70 kDa.

### Specific Activity

22 U/mg. One unit of Aurora-B activity is defined as 1 nmol phosphate incorporated into Ser/Thr 1 peptide substrate per minute at pH 7.4 and 30 °C.

### Assay Condition

50 mM HEPES, pH 7.4, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 1 mM DTT, 3 µM Na-orthovanadate, 0.1 mM ATP, 4 µM Aurora B substrate, and 4 µg/ml Aurora kinase B.

### Purity

> 80% by Coomassie-stained SDS-PAGE.

### Formulation

Recombinant protein is in storage buffer (25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 3 mM DTT, and 50% glycerol).

### Storage and Stability

Store product frozen at or below -70 °C. Enzyme will be stable for at least 6 months at -70 °C as undiluted stock. Freeze in working aliquots to avoid repeated thawing and freezing.

### Application

Aurora kinase B is useful for the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

### References

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

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

TDS E-K057 Rev: 1 (03/15/06)