

## Data Sheet

### Aurora A Protein Kinase, His-Tag

Human, Recombinant, N-terminal His-tag  
**Catalog #:** 100112  
**Lot#:** 180622-3      **Conc.:** 0.12 mg/ml

**Formulated in:** 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 200mM Imidazole, 20% glycerol, and 3 mM DTT.

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

**References:**

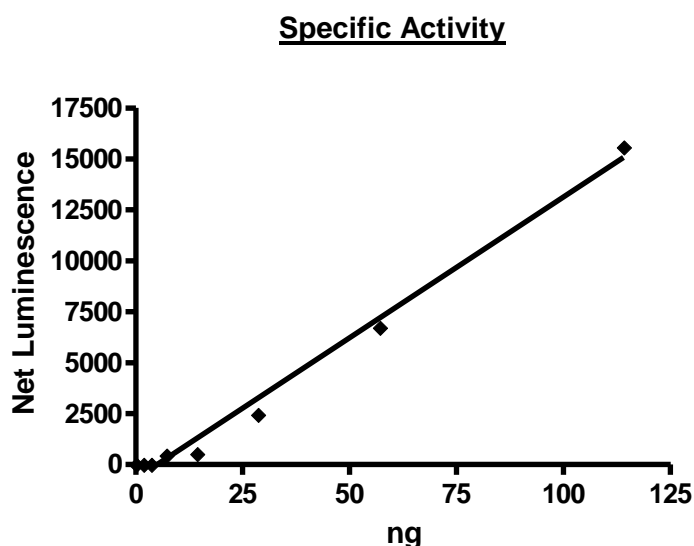
1. Ouchi, M. *et al.*, *J. Biol Chem.* **279** (19), 19643-19648 (2004).
2. Kunitoku, N. *et al.*, *Dev. Cell* **5**(6), 853-864 (2003).

**Description:** Human Aurora kinase A also known as, AURKA, Serine/Threonine-Protein Kinase Aurora-A, and Breast Tumor-Amplified Kinase, GenBank Accession No. NM\_003600, a.a. 2-403 (end) with N-terminal His-tag, expressed in a HEK 293 cell expression system. MW = 46 kDa.

**Specific Activity:** 45 pmole/min/ $\mu\text{g}$ .  
**Assay Conditions:** Assay was done in Kinase buffer containing 2 mM DTT using KEMPTide substrate (0.2 mg/ml) and 10  $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  for 45 min. Amount of ATP transferred was calculated using ADP-Glo reagent.

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

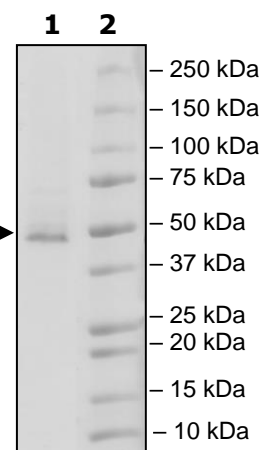
## Quality Assurance



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

**Lane 1:**  
2 $\mu\text{g}$  Aurora A  
**Lane 2:**  
Protein Marker

**MW:** 46 kDa  
**Purity:**  $\geq 75\%$



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)