

Data Sheet

Aurora A Protein Kinase, His-Tag

Human, Recombinant, N-terminal His-tag
Catalog #: 40004
Lot#: 190402 **Conc.:** 0.25 mg/ml

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

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

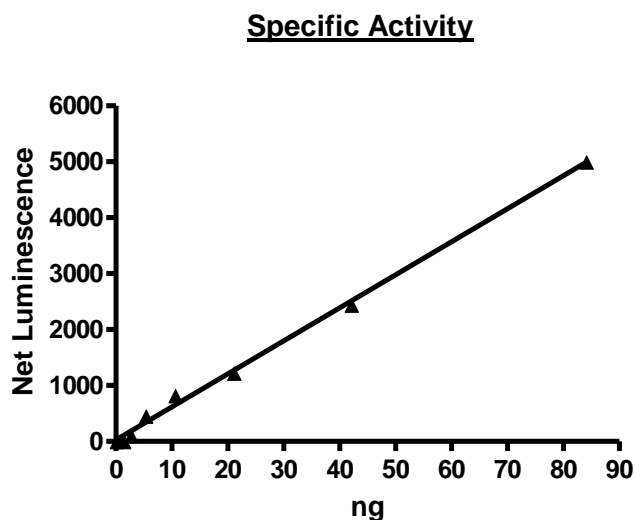
1. Ouchi, M. *et al.*, *J. Biol Chem.* **279** (19), 19643-19648 (2004).
2. Kunitoku, N. *et al.*, *Dev. Cell* **5**(6), 853-864 (2003).

Description: Human Aurora Kinase A also known as STK15, STK6, BTAK, HARK1, ARK1, and AURA, GenBank Accession No. NM_003600, a.a. 1-403 (end) with N-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system, MW=50 kDa.

Specific Activity: ≥45 pmole/min/μg. Assay Conditions: Assay was done in Kinase buffer containing 2 mM DTT using KEMPTide substrate (0.2 mg/ml) and 10 μM ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using ADP-Glo reagent (Promega).

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

Quality Assurance



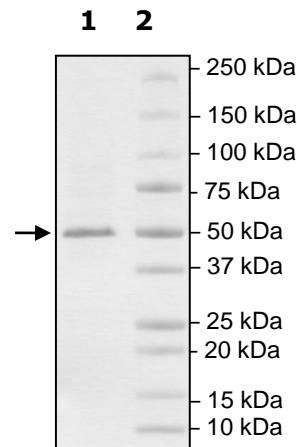
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
Aurora A

Lane 2:
Protein Marker

Purity: ≥90%

MW:
50 kDa



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
 Please visit our website at: www.bpsbioscience.com