

# PRODUCT INFORMATION



## Aurora Kinase Inhibitor II

Item No. 17541

CAS Registry No.: 331770-21-9

Formal Name: N-[4-[(6,7-dimethoxy-4-quinazolyl)amino]phenyl]-benzamide

Synonym: 4-(4'-Benzamidoanilino)-6,7-dimethoxyquinazoline

MF: C<sub>23</sub>H<sub>20</sub>N<sub>4</sub>O<sub>3</sub>

FW: 400.4

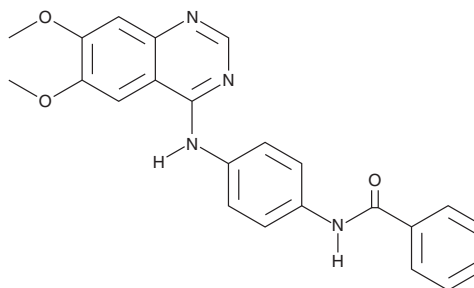
Purity: ≥98%

UV/Vis.: λ<sub>max</sub>: 223, 262, 354 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### Laboratory Procedures

Aurora kinase inhibitor II is supplied as a crystalline solid. A stock solution may be made by dissolving the Aurora kinase inhibitor II in the solvent of choice. Aurora kinase inhibitor II is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of Aurora kinase inhibitor II in these solvents is approximately 2.5 and 0.25 mg/ml, respectively.

Aurora kinase inhibitor II is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Aurora kinase inhibitor II should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Aurora kinase inhibitor II has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

The Aurora kinases are a family of serine/threonine kinases that are key regulators of mitosis and cytokinesis. Aurora kinase inhibitor II is a cell-permeable anilinoquinazoline that blocks the activity of Aurora A (IC<sub>50</sub> = 0.39 μM).<sup>1</sup> In MCF-7 cells, it blocks proliferation with an IC<sub>50</sub> value of 1.25 μM.<sup>1</sup>

### Reference

1. Heron, N.M., Anderson, M., Blowers, D.P., *et al.* SAR and inhibitor complex structure determination of a novel class of potent and specific Aurora kinase inhibitors. *Bioorg. Med. Chem. Lett.* **16(5)**, 1320-1323 (2006).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/05/2016

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM