

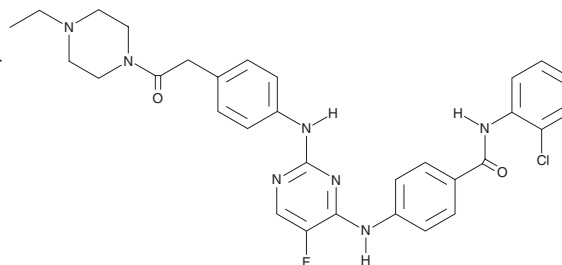
# PRODUCT INFORMATION



## Aurora A Inhibitor I

Item No. 21600

**CAS Registry No.:** 1158838-45-9  
**Formal Name:** N-(2-chlorophenyl)-4-[[2-[[4-[2-(4-ethyl-1-piperazinyl)-2-oxoethyl]phenyl]amino]-5-fluoro-4-pyrimidinyl]amino]-benzamide  
**MF:** C<sub>31</sub>H<sub>31</sub>ClFN<sub>7</sub>O<sub>2</sub>  
**FW:** 588.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 281, 338 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Aurora A inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the Aurora A inhibitor I in the solvent of choice, which should be purged with an inert gas. Aurora A inhibitor I is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Aurora A inhibitor I in these solvents is approximately 33 mg/ml.

### Description

Aurora A inhibitor I is a potent and selective inhibitor of Aurora A kinase (IC<sub>50</sub>s = 3.4 and 3,400 nM for Aurora A and Aurora B, respectively).<sup>1</sup> It prevents centrosome separation and reduces phosphorylated Aurora A, but not phosphorylated histone H3, levels at the centrosomes, indicating inhibition is specific to Aurora A kinase. Aurora A inhibitor I reduces the proliferation of HCT116 and HT-29 colorectal carcinoma cells *in vitro* (IC<sub>50</sub>s = 0.19 and 2.9 μM, respectively).

### Reference

1. Aliagas-Martin, I., Burdick, D., Corson, L., *et al.* A class of 2,4-bisanilinopyrimidine Aurora A inhibitors with unusually high selectivity against Aurora B. *J. Med. Chem.* **52(10)**, 3300-3307 (2009).

#### 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, 11/11/2022

#### 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