

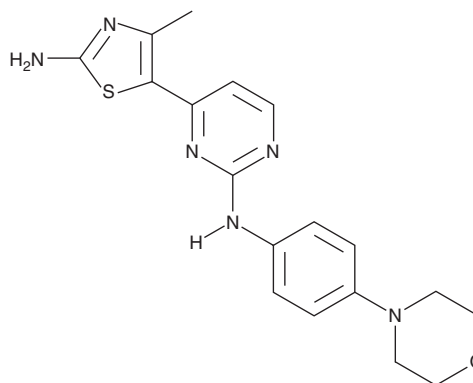
PRODUCT INFORMATION



CYC-116

Item No. 21205

CAS Registry No.: 693228-63-6
Formal Name: 4-(2-amino-4-methyl-5-thiazolyl)-
N-[4-(4-morpholinyl)phenyl]-2-
pyrimidinamine
MF: C₁₈H₂₀N₆OS
FW: 368.5
Purity: ≥98%
UV/Vis.: λ_{max}: 295, 337 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

CYC-116 is supplied as a crystalline solid. A stock solution may be made by dissolving the CYC-116 in the solvent of choice. CYC-116 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of CYC-116 in these solvents is approximately 10 and 1.25 mg/ml, respectively.

CYC-116 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CYC-116 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CYC-116 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CYC-116 is a potent Aurora kinase inhibitor exhibiting K_i values of 8.0 and 9.2 nM for Aurora kinases A and B, respectively.¹ It is significantly less potent against a panel of cyclin-dependent kinases and other related kinases, with the exception of Flt-3 (K_i = 44 nM). CYC-116 exhibits broad-spectrum antiproliferative activity against a panel of human cancer cell lines with a mean IC₅₀ value of 0.54 μM (range = 0.09 – 1.6 μM). It is orally bioavailable and inhibits the growth of human tumor xenografts in mice.

Reference

1. Wang, S., Midgley, C.A., Scaërou, F., *et al.* Discovery of N-phenyl-4-(thiazol-5-yl)pyrimidin-2-amine aurora kinase inhibitors. *J. Med. Chem.* **53**(11), 4367-4378 (2010).

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, 12/07/2022

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