

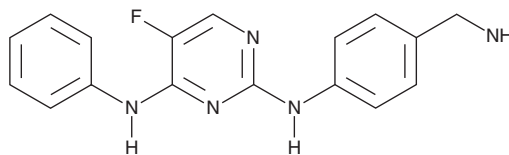
PRODUCT INFORMATION



CZC-8004

Item No. 16718

CAS Registry No.: 916603-07-1
Formal Name: N²-[4-(aminomethyl)phenyl]-5-fluoro-N⁴-phenyl-2,4-pyrimidinediamine
Synonym: Dianilinopyrimidine-01
MF: C₁₇H₁₆FN₅
FW: 309.3
Purity: ≥98%
UV/Vis.: λ_{max}: 274, 321 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

CZC-8004 is supplied as a crystalline solid. A stock solution may be made by dissolving the CZC-8004 in the solvent of choice, which should be purged with an inert gas. CZC-8004 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CZC-8004 in these solvents is approximately 0.1, 5, and 2 mg/ml, respectively.

CZC-8004 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CZC-8004 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CZC-8004 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CZC-8004 is an aminopyrimidine that binds a range of tyrosine kinases, including ABL, BTK, FAK, FER, JAK1, SRC, SYK, TEC, TNK1, TYK2, and YES.¹ It inhibits these kinases at low micromolar concentrations and can be immobilized to capture and purify these enzymes.¹

Reference

1. Klutchko, S.R., Hamby, J.M., Boschelli, D.H., *et al.* 2-Substituted aminopyrido[2,3-*d*]pyrimidin-7(8*H*)-ones. Structure-activity relationships against selected tyrosine kinases and in vitro and in vivo anticancer activity. *J. Med. Chem.* **41**(17), 3276-3292 (1998).

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, 11/09/2022

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