

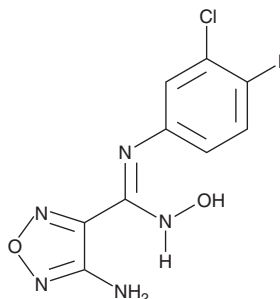
PRODUCT INFORMATION



INCB024360 analog

Item No. 18265

CAS Registry No.: 914471-09-3
Formal Name: 4-amino-N'-(3-chloro-4-fluorophenyl)-N-hydroxy-1,2,5-oxadiazole-3-carboximidamide
Synonym: IDO5L
MF: C₉H₇ClFN₅O₂
FW: 271.6
Purity: ≥98%
UV/Vis.: λ_{max}: 246, 285 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

INCB024360 is supplied as a crystalline solid. A stock solution may be made by dissolving the INCB024360 in the solvent of choice. INCB024360 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of INCB024360 in these solvents is approximately 30 mg/ml.

INCB024360 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, INCB024360 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. INCB024360 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Indoleamine 2,3-dioxygenase (IDO) metabolizes tryptophan to kynurenine, leading to the production of NAD⁺ via the kynurenine pathway. It has been implicated in mediating pathological immunosuppression associated with certain diseases, including cancer. INCB024360 analog is a potent IDO1 inhibitor that demonstrates IC₅₀ values of 67 and 19 nM for human IDO in enzymatic activity and HeLa cell assays, respectively.¹⁻³ It is inactive against tryptophan 2,3-dioxygenase (IC₅₀ > 10 μM).² At 25-75 mg/kg, subcutaneous administration of INCB024360 analog to mice significantly decreased kynurenine levels in plasma and dose-dependently reduced the growth of GM-CSF-secreting B16 tumor xenografts.² A fluorine-18 moiety has been used to label this compound for use as a probe for imaging IDO1 expression via positron emission tomography.⁴

Reference

1. Liu, X., Shin, N., Koblisch, H.K., et al. *Blood* **115**(17), 3520-3530 (2010).
2. Yue, E.W., Douty, B., Wayland, B., et al. *J. Med. Chem.* **52**(23), 7364-7367 (2009).
3. Tomek, P., Palmer, B.D., Flanagan, J.U., et al. *Anal. Bioanal. Chem.* **405**(8), 2515-2524 (2013).
4. Huang, X., Gillies, R.J., and Tian, H. *J. Labelled Comp. Radiopharm.* **58**(4), 156-162 (2015).

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, 12/02/2022

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