

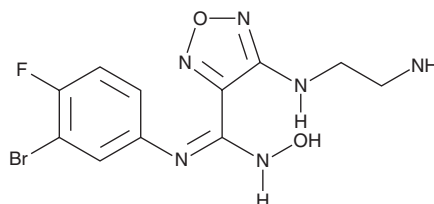
PRODUCT INFORMATION



IDO-IN-3

Item No. 28432

CAS Registry No.: 2070018-30-1
Formal Name: 4-[(2-aminoethyl)amino]-N'-(3-bromo-4-fluorophenyl)-N-hydroxy-1,2,5-oxadiazole-3-carboximidamide
MF: C₁₁H₁₂BrFN₆O₂
FW: 359.2
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 290 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

IDO-IN-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the IDO-IN-3 in the solvent of choice, which should be purged with an inert gas. IDO-IN-3 is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml.

Description

IDO-IN-3 is an inhibitor of indoleamine-2,3-dioxygenase 1 (IDO1; IC₅₀s = 290 and 98 nM for the human recombinant enzyme in a cell-free assay and in HeLa cells, respectively).¹

Reference

1. Yue, E.W., Sparks, R.B., Polam, P., *et al.* INCB24360 (epacadostat), a highly potent and selective indoleamine-2,3-dioxygenase 1 (IDO1) inhibitor for immuno-oncology. *Med. Chem. Lett.* **8**(5), 486-491 (2017).

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, 10/11/2022

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