

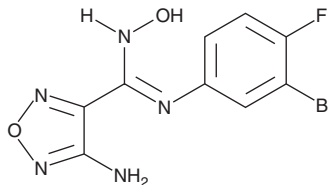
PRODUCT INFORMATION



IDO-IN-1

Item No. 19402

CAS Registry No.: 914638-30-5
Formal Name: 4-amino-N-(3-bromo-4-fluorophenyl)-N'-hydroxy-1,2,5-oxadiazole-3-carboximidamide
MF: C₉H₇BrFN₅O₂
FW: 316.1
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 282 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-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the IDO-IN-1 in the solvent of choice, which should be purged with an inert gas. IDO-IN-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of IDO-IN-1 in these solvents is approximately 10 mg/ml.

IDO-IN-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IDO-IN-1 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. IDO-IN-1 has a solubility of approximately 0.025 mg/ml in a 1:30 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

IDO-IN-1 is a potent indoleamine 2,3-dioxygenase (IDO) inhibitor that demonstrates IC₅₀ values of 59 and 12 nM for human IDO enzymatic activity and HeLa cell assays, respectively.¹ It is inactive against tryptophan 2,3-dioxygenase (TOD; IC₅₀ > 10 μM).

References

1. Yue, E.W., Douty, B., Wayland, B., *et al.* Discovery of potent competitive inhibitors of indoleamine 2,3-dioxygenase with in vivo pharmacodynamic activity and efficacy in a mouse melanoma model. *J. Med. Chem.* **52**(23), 7364-7367 (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, 10/27/2022

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