

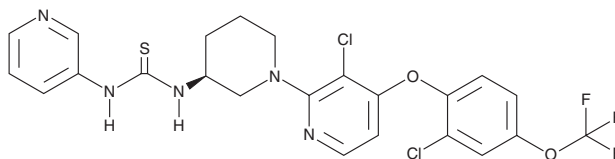
PRODUCT INFORMATION



(S)-DO271

Item No. 30038

CAS Registry No.: 2301865-01-8
Formal Name: N-[(3S)-1-[3-chloro-4-[2-chloro-4-(trifluoromethoxy)phenoxy]-2-pyridinyl]-3-piperidinyl]-N'-3-pyridinyl-thiourea
MF: C₂₃H₂₀Cl₂F₃N₅O₂S
FW: 558.4
Purity: ≥98%
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

(S)-DO271 is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-DO271 in the solvent of choice, which should be purged with an inert gas. (S)-DO271 is soluble in the organic solvent DMSO.

Description

(S)-DO271 is an inactive control for the α/β -hydrolase domain-containing protein 12 (ABHD12) inhibitor DO264 (Item No. 29402).^{1,2} Unlike DO264, (S)-DO271 does not inhibit ABHD12-dependent hydrolysis of lysophosphatidylserine (lyso-PS) in mouse brain membrane lysates when used at a concentration of 1 μ M or modify inflammatory cytokine production in M1-polarized THP-1 macrophages.

References

- Ogasawara, D., Ichu, T.-A., Vartabedian, V.F., *et al.* Selective blockade of the lyso-PS lipase ABHD12 stimulates immune responses in vivo. *Nat. Chem. Biol.* **14**(12), 1099-1108 (2018).
- Ogasawara, D., Ichu, T.-A., Jing, H., *et al.* Discovery and optimization of selective and in vivo active inhibitors of the lysophosphatidylserine lipase α/β -hydrolase domain-containing 12 (ABHD12). *J. Med. Chem.* **62**(3), 1643-1656 (2019).

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, 03/21/2024

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