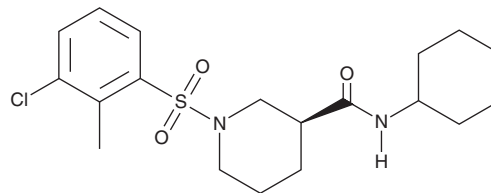


PRODUCT INFORMATION



INCB 13739
Item No. 40223

CAS Registry No.: 872506-67-7
Formal Name: (3S)-1-[(3-chloro-2-methylphenyl) sulfonyl]-N-cyclohexyl-3-piperidinecarboxamide
MF: C₁₉H₂₇ClN₂O₃S
FW: 398.9
Purity: ≥98%
Supplied as: A 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

INCB 13739 is supplied as a solid. A stock solution may be made by dissolving the INCB 13739 in the solvent of choice, which should be purged with an inert gas. INCB 13739 is soluble in methanol and DMSO.

Description

INCB 13739 is an inhibitor of 11 β -hydroxysteroid dehydrogenase 1 (11 β -HSD1; IC₅₀ = 3.2 nM).¹ It is greater than 1,000-fold selective for 11 β -HSD1 over 11 β -HSD2, the mineralocorticoid receptor, and the glucocorticoid receptor at 5 μ M. INCB 13739 inhibits 11 β -HSD1 activity in peripheral blood mononuclear cells (PBMCs; IC₅₀ = 1.1 nM).

Reference

1. Zhang, C., Xu, M., He, C., *et al.* Discovery of 1'-(1-phenylcyclopropane-carbonyl)-3H-spiro [isobenzofuran-1,3'-pyrrolidin]-3-one as a novel steroid mimetic scaffold for the potent and tissue-specific inhibition of 11 β -HSD1 using a scaffold-hopping approach. *Bioorg. Med. Chem. Lett.* **69**:128782, (2022).

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, 02/01/2024

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