

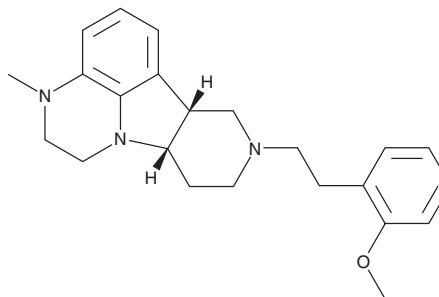
PRODUCT INFORMATION



IHCH-7079

Item No. 41988

CAS Registry No.: 2957888-63-8
Formal Name: 2,3,6bR,7,8,9,10,10aS-octahydro-8-[2-(2-methoxyphenyl)ethyl]-3-methyl-1H-pyrido[3',4':4,5]pyrrolo[1,2,3-de]quinoxaline
MF: C₂₃H₂₉N₃O
FW: 363.5
Purity: ≥98%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

IHCH-7079 is a β -arrestin-biased agonist of the serotonin (5-HT) receptor subtype 5-HT_{2A} (K_i = 16.98 nM).¹ It reduces the time spent immobile in the tail suspension and forced swim tests in mice in a model of acute restraint stress-induced depression when administered at a dose of 10 mg/kg, and this effect can be blocked by the 5-HT_{2A} receptor antagonist MDL100907 (volinanserin; Item No. 15936). IHCH-7079 (2 and 10 mg/kg) does not induce the head-twitch response in mice, indicating a lack of hallucinogenic potential.

Reference

1. Cao, D., Yu, J., Wang, H., *et al.* Structure-based discovery of nonhallucinogenic psychedelic analogs. *Science* **375**(6579), 403-411 (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, 06/26/2026

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