

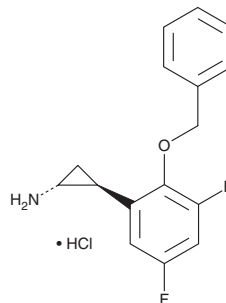
PRODUCT INFORMATION



S 2101

Item No. 19802

CAS Registry No.: 1239262-36-2
Formal Name: (1R,2S)-rel-2-[3,5-difluoro-2-(phenylmethoxy)phenyl]-cyclopropanamine, monohydrochloride
Synonyms: KDM1 Inhibitor II, LSD1 Inhibitor II, Lysine-specific Demethylase 1 Inhibitor II
MF: C₁₆H₁₅F₂NO • HCl
FW: 311.8
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

S 2101 is supplied as a solid. A stock solution may be made by dissolving the S 2101 in water. We do not recommend storing the aqueous solution for more than one day.

Description

S 2101 is an inhibitor of lysine-specific demethylase 1 (LSD1; $K_i = 0.61 \mu\text{M}$).¹ It is selective for LSD1 over monoamine oxidase A (MAO-A) and MAO-B (K_i s = 17 and 110 μM , respectively).

Reference

1. Mimasu, S., Umezawa, N., Sato, S., *et al.* Structurally designed trans-2-phenylcyclopropylamine derivatives potently inhibit histone demethylase LSD1/KDM1. *Biochemistry* **49(30)**, 6494-6503 (2010).

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, 11/06/2023

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