

PRODUCT INFORMATION

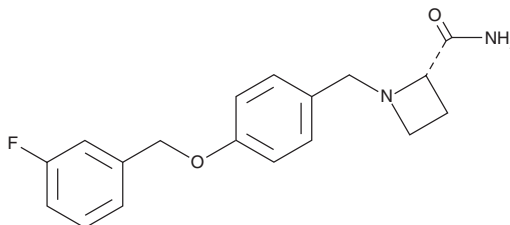


C3

Item No. 33166

CAS Registry No.: 2286257-25-6
Formal Name: 1-[[4-[(3-fluorophenyl)methoxy]phenyl]methyl]-2S-azetidinecarboxamide

MF: C₁₈H₁₉FN₂O₂
FW: 314.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

C3 is supplied as a crystalline solid. A stock solution may be made by dissolving the C3 in the solvent of choice, which should be purged with an inert gas. C3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of C3 in ethanol and DMSO is approximately 2 mg/ml and approximately 15 mg/ml in DMF.

Description

C3 is an inhibitor of monoamine oxidase B (MAO-B; IC₅₀ = 0.021 μM).¹ It is selective for MAO-B over MAO-A (IC₅₀ = 26.8 μM). C3 (0.1, 1, and 3 mg/kg) prevents MPTP-induced decreases in brain dopamine levels in mice.

Reference

1. Jin, C.-F., Wang, Z.-Z., Chen, K.-Z., *et al.* Computational fragment-based design facilitates discovery of potent and selective monoamine oxidase-B (MAO-B) inhibitor. *J. Med. Chem.* **63**(23), 15021-15036 (2020).

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, 10/08/2025

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