

PRODUCT INFORMATION



CC-90011

Item No. 32727

CAS Registry No.: 2097523-60-7

Formal Name: 4-[2-(4-amino-1-piperidiny)-5-(3-fluoro-4-methoxyphenyl)-1,6-dihydro-1-methyl-6-oxo-4-pyrimidinyl]-2-fluoro-benzonitrile, monobenzenesulfonate

Synonyms: LSD1-IN-7, LSD1 Inhibitor 7, Lysine-specific Demethylase 1-IN-7, Lysine-specific Demethylase 1 Inhibitor 7

MF: C₂₄H₂₃F₂N₅O₂ • C₆H₆O₃S

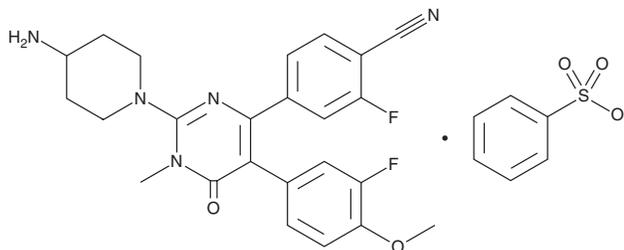
FW: 609.6

Purity: ≥95%

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

CC-90011 is supplied as a solid. A stock solution may be made by dissolving the CC-90011 in the solvent of choice, which should be purged with an inert gas. CC-90011 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

CC-90011 is an inhibitor of lysine-specific demethylase 1 (LSD1; IC₅₀ = 0.3 nM).¹ It is selective for LSD1 over LSD2, monoamine oxidase A (MAO-A), and MAO-B at 10 μM. CC-90011 induces differentiation in THP-1 cells (EC₅₀ = 7 nM) and inhibits the proliferation of NCI H1417 small cell lung cancer (SCLC) cells (EC₅₀ = 6 nM). It also inhibits lipid accumulation induced by palmitic acid (Item No. 10006627) and oleic acid (Item Nos. 90260 | 24659) in Huh7 hepatocytes when used at a concentration of 20 μM.² *In vivo*, CC-90011 (2.5 mg/kg) reduces tumor volume in an NCI H1417 mouse xenograft model.¹

References

1. Kanouni, T., Severin, C., Cho, R.W., *et al.* Discovery of CC-90011: A potent and selective reversible inhibitor of lysine specific demethylase 1 (LSD1). *J. Med. Chem.* **63**(23), 14522-14529 (2020).
2. Yang, Z., Zhang, S., Liu, X., *et al.* Histone demethylase KDM1A promotes hepatic steatosis and inflammation by increasing chromatin accessibility in NAFLD. *J. Lipid Res.* **65**(3), 100513 (2024).

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, 07/07/2025

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