

PRODUCT INFORMATION



DS9300

Item No. 43812

CAS Registry No.: 2259641-46-6
Formal Name: (2R,4R)-1-[[4,4-difluoro-1-(4-methoxyphenyl)cyclohexyl]carbonyl]-4-fluoro-N-1H-pyrazolo[4,3-b]pyridin-5-yl-2-pyrrolidinecarboxamide

MF: C₂₅H₂₆F₃N₅O₃

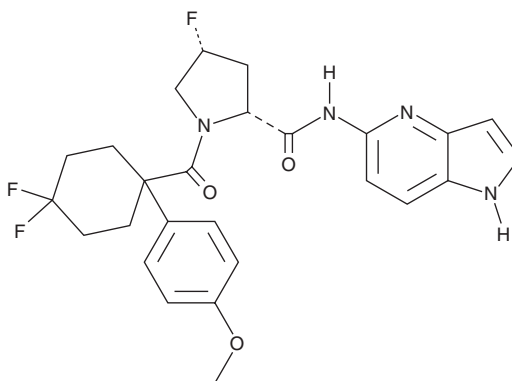
FW: 501.5

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

DS9300 is supplied as a solid. A stock solution may be made by dissolving the DS9300 in the solvent of choice, which should be purged with an inert gas. DS9300 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

DS9300 is an inhibitor of histone acetyltransferase p300 (IC₅₀ = 28 nM).¹ It inhibits acetylation of lysine 27 on histone H3 (H3K27) in LK-2 cells (IC₅₀ = 54 nM). DS9300 inhibits the growth of VCaP prostate cancer cells (GI₅₀ = 0.6 nM). *In vivo*, DS9300 (0.3-3 mg/kg) reduces tumor volume in a VCaP castrated mouse xenograft model.

Reference

1. Kanada, R., Kagoshima, Y., Suzuki, T., *et al.* Discovery of DS-9300: A highly potent, selective, and once-daily oral EP300/CBP histone acetyltransferase inhibitor. *J. Med. Chem.* **66**(1), 695-715 (2023).

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, 07/21/2025

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