

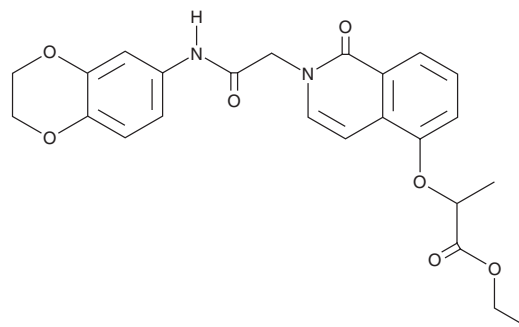
PRODUCT INFORMATION



SID 7969543

Item No. 43705

CAS Registry No.: 868224-64-0
Formal Name: 2-[[2-[2-[(2,3-dihydro-1,4-benzodioxin-6-yl)amino]2-oxoethyl]-1,2-dihydro-1-oxo-5-isoquinolinyl]oxy]-propanoic acid, ethyl ester
MF: C₂₄H₂₄N₂O₇
FW: 452.5
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

SID 7969543 is supplied as a solid. A stock solution may be made by dissolving the SID 7969543 in the solvent of choice, which should be purged with an inert gas. SID 7969543 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SID 7969543 is an inhibitor of nuclear receptor subfamily 5 group A member 1 (NR5A1), also known as steroidogenic factor 1 (SF-1; IC₅₀ = 760 nM).¹ It is selective for NR5A1/SF-1 over retinoic acid receptor-related orphan receptor α (ROR α) and the herpes simplex virus (HSV) transcriptional activator VP16 (IC₅₀s = >33.333 μ M for both). SID 7969543 is selectively cytotoxic to PER-485, THP-1, and MOLM-13 leukemia cells, which contain mixed-lineage leukemia 1 (MLL1) rearrangement mutations (IC₅₀s = 1.4, 2.6, and 1 μ M, respectively), over a panel of wild-type MLL1-containing leukemia cells and MRC-5 fibroblasts (IC₅₀s = >20 μ M for all).²

References

1. Madoux, F., Li, X., Chase, P., *et al.* Potent, selective and cell penetrant inhibitors of SF-1 by functional ultra-high-throughput screening. *Mol. Pharmacol.* **73(6)**, 1776-1784 (2008).
2. Karsa, M., Ronca, E., Bongers, A., *et al.* Systematic *in vitro* evaluation of a library of approved and pharmacologically active compounds for the identification of novel candidate drugs for KMT2A-rearranged leukemia. *Front. Oncol.* **11**, 779859 (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, 07/24/2025

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