

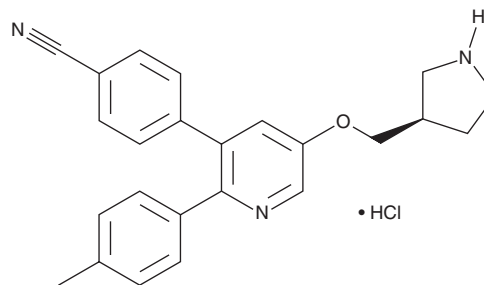
PRODUCT INFORMATION



GSK690 (hydrochloride)

Item No. 45320

CAS Registry No.: 2436760-79-9
Formal Name: 4-[2-(4-methylphenyl)-5-[(3R)-3-pyrrolidinylmethoxy]-3-pyridinyl]-benzonitrile, monohydrochloride
MF: C₂₄H₂₃N₃O • HCl
FW: 405.9
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

GSK690 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the GSK690 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GSK690 (hydrochloride) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

GSK690 is a lysine-specific demethylase 1 (LSD1) inhibitor (IC₅₀ = 0.09 μM in a cell-free assay).¹ It is selective for LSD1 over monoamine oxidase A (MAO-A; IC₅₀ = >200 μM). GSK690 inhibits LSD1 activity in THP-1 acute myeloid leukemia cells (IC₅₀ = 1.4 μM). GSK690 potentiates reductions in viability of SJCRH30 rhabdomyosarcoma cells induced by the histone deacetylase (HDAC) inhibitor JNJ-26481585 (Item No. 14088).²

References

1. Hitchin, J.R., Blagg, J., Burke, R., *et al.* Development and evaluation of selective, reversible LSD1 inhibitors derived from fragments. *MedChemComm* **4**, 1513 (2013).
2. Haydn, T., Metzger, E., Schuele, R., *et al.* Concomitant epigenetic targeting of LSD1 and HDAC synergistically induces mitochondrial apoptosis in rhabdomyosarcoma cells. *Cell Death Dis.* **8(6)**, e2879 (2017).

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, 04/24/2026

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