

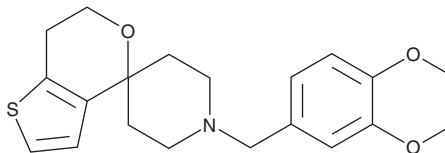
PRODUCT INFORMATION



GSK2200150A

Item No. 36620

CAS Registry No.: 1443138-53-1
Formal Name: 1-[(2,3-dihydro-1,4-benzodioxin-6-yl)methyl]-6',7'-dihydro-spiro[piperidine-4,4'-[4H]thieno[3,2-c]pyran
MF: C₂₀H₂₃NO₃S
FW: 357.5
Purity: ≥98%
UV/Vis.: λ_{max}: 229 nm
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

GSK2200150A is supplied as a solid. A stock solution may be made by dissolving the GSK2200150A in the solvent of choice, which should be purged with an inert gas. GSK2200150A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GSK2200150A in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

GSK2200150A is a spirocyclic antimycobacterial agent.¹ It is active against *M. tuberculosis* and *M. bovis* (MICs = 0.38 and 1.6 μM, respectively).

References

1. Balcells, L., Bates, R.H., Young, R.J., *et al.* Fueling open-source drug discovery: 177 small-molecule leads against tuberculosis. *ChemMedChem* **8**(2), 313-321 (2013).

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, 06/14/2023

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