

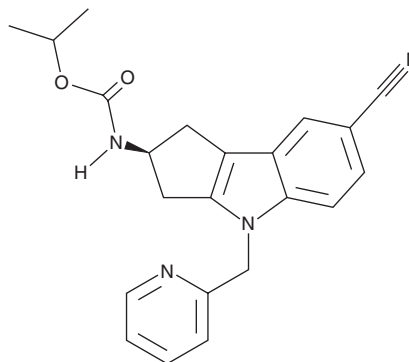
PRODUCT INFORMATION



LY2452473

Item No. 43312

CAS Registry No.: 1029692-15-6
Formal Name: N-[(2S)-7-cyano-1,2,3,4-tetrahydro-4-(2-pyridinylmethyl)cyclopent[b]indol-2-yl]-carbamic acid, 1-methylethyl ester
Synonym: TT 701
MF: C₂₂H₂₂N₄O₂
FW: 374.4
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

LY2452473 is supplied as a solid. A stock solution may be made by dissolving the LY2452473 in the solvent of choice, which should be purged with an inert gas. LY2452473 is soluble (≥10 mg/ml) in DMSO.

Description

LY2452473 is an androgen receptor antagonist (IC₅₀ = 8.565 nM).¹

Reference

1. Popović, L., Brankatschk, B., Palladino, G., *et al.* Polypharmacological profiling across protein target families and cellular pathways using the multiplexed cell-based assay platform safetyProfiler reveals efficacy, potency and side effects of drugs. *Biomed. Pharmacother.* **180**, 117523 (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, 04/29/2025

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