

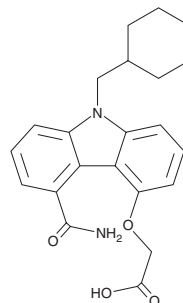
PRODUCT INFORMATION



LY433771

Item No. 35406

CAS Registry No.: 220862-64-6
Formal Name: 2-[[5-(aminocarbonyl)-9-(cyclohexylmethyl)-9H-carbazol-4-yl]oxy]-acetic acid
Synonym: LSN433771
MF: C₂₂H₂₄N₂O₄
FW: 380.4
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 250, 291, 346 nm
Supplied as: A crystalline 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

LY433771 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY433771 in the solvent of choice, which should be purged with an inert gas. LY433771 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LY433771 in these solvents is approximately 3 mg/ml.

Description

LY433771 is an inhibitor of type X secretory phospholipase A₂ (sPLA₂; IC₅₀ = 3 nM).¹

Reference

1. Ono, T., Ueno, M., and Hanasaki, K. V type and/or x type sPLA₂ inhibitors. *Nogi Pharmaceutical Co., Ltd. WO2001026653A1* (2001).

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, 10/09/2025

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