

PRODUCT INFORMATION



LY2922470

Item No. 43109

CAS Registry No.: 1423018-12-5
Formal Name: (βS)-4-[[5-[(3,4-dihydro-8-methoxy-1(2H)-quinolinyl)methyl]-2-thienyl]methoxy]-β-1-propyn-1-yl-benzenepropanoic acid

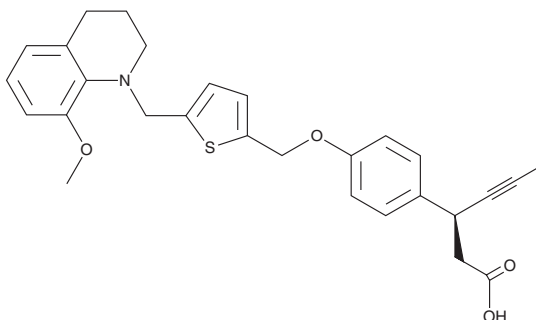
MF: C₂₈H₂₉NO₄S

Purity: ≥95%

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

LY2922470 is supplied as a solid. A stock solution may be made by dissolving the LY2922470 in the solvent of choice, which should be purged with an inert gas. LY2922470 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

LY2922470 is an agonist of free fatty acid receptor 1 (FFAR1/GPR40).¹ It induces calcium flux in HEK293 cells expressing FFAR1/GPR40 (EC₅₀s = 7, 1, and 3 nM for the human, mouse, and rat receptor, respectively). LY2922470 (30 mg/kg) increases blood levels of glucagon-like peptide 1 (GLP-1) in fasted mice. It also decreases cerebral infarct area in a mouse model of ischemia-reperfusion injury induced by middle cerebral artery occlusion (MCAO) when administered at doses of 10, 20, or 40 mg/kg.²

References

1. Hamdouchi, C., Kahl, S.D., Lewis, A.P., *et al.* The discovery, preclinical, and early clinical development of potent and selective GPR40 agonists for the treatment of type 2 diabetes mellitus (LY2881835, LY2922083, and LY2922470). *J. Med. Chem.* **59**(24), 10891-10916 (2016).
2. Lu, Y., Zhou, W., Cui, Q., *et al.* G protein-coupled receptor 40 agonist LY2922470 alleviates ischemic-stroke-induced acute brain injury and functional alterations in mice. *Int. J. Mol. Sci.* **24**(15), 12244 (2023).

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, 04/16/2025

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