

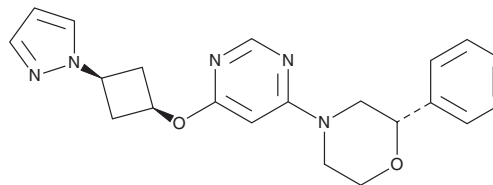
PRODUCT INFORMATION



ELOVL1-22

Item No. 36142

Formal Name: (S)-4-(6-((1S,3R)-3-(1H-pyrazol-1-yl)cyclobutoxy)pyrimidin-4-yl)-2-phenylmorpholine
MF: C₂₁H₂₃N₅O₂
FW: 377.4
Purity: ≥98%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

ELOVL1-22 is an inhibitor of ELOVL fatty acid elongase 1 (ELOVL1).¹ It reduces lysohexacosanoylphosphatidylcholine (LPC(C26:0)) synthesis in HEK293 cells (IC₅₀ = 400 pM), a marker of ELOVL1 inhibition, and is selective for ELOVL1 over ELOVL4, -6, and -7 (IC₅₀s = >50 μM for all), as well as a panel of 34 receptors at 10 μM. ELOVL1-22 (1-32 mg/kg) reduces blood and CNS levels of LPC(C26:0) in an *Abcd1*^{-/-} mouse model of adrenoleukodystrophy (ALD), a neurodegenerative disease characterized by rapid inflammatory demyelination of the CNS.

Reference

1. Boyd, M.J., Collier, P.N., Clark, M.P., *et al.* Discovery of novel, orally bioavailable pyrimidine ether-based inhibitors of ELOVL1. *J. Med. Chem.* **64**(24), 17777-17794 (2021).

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, 12/01/2025

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