

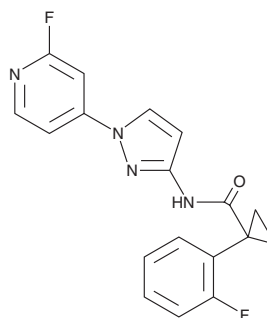
# PRODUCT INFORMATION



## ELOVL1-27

Item No. 35323

**CAS Registry No.:** 2227482-41-7  
**Formal Name:** 1-(2-fluorophenyl)-N-[1-(2-fluoro-4-pyridinyl)-1H-pyrazol-3-yl]-cyclopropanecarboxamide  
**MF:** C<sub>18</sub>H<sub>14</sub>F<sub>2</sub>N<sub>4</sub>O  
**FW:** 340.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 289 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

ELOVL1-27 is supplied as a crystalline solid. A stock solution may be made by dissolving the ELOVL1-27 in the solvent of choice, which should be purged with an inert gas. ELOVL1-27 is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

### Description

ELOVL1-27 is an inhibitor of ELOVL fatty acid elongase 1 (ELOVL1; IC<sub>50</sub> = 900 nM).<sup>1</sup> It reduces lysohexacosanoylphosphatidylcholine (LPC(C26:0)) synthesis in HEK293 cells (IC<sub>50</sub> = 13 nM), as well as in primary fibroblasts isolated from patients with cerebral adrenoleukodystrophy (CALD) or adrenomyeloneuropathy (AMN; IC<sub>50</sub>s = 15-43 nM), neurodegenerative diseases characterized by rapid inflammatory demyelination of the CNS. ELOVL1-27 (2, 16, and 32 mg/kg) reduces blood LPC(C26:0) levels in wild-type mice and rats. It also reduces CNS levels of LPC(C26:0) in an ABCD1<sup>-/-</sup> mouse model of ALD when administered at a dose of 1 or 10 mg/kg per day for at least three months.

### Reference

1. Come, J.H., Senter, T.J., Clark, M.P., *et al.* Discovery and optimization of pyrazole amides as inhibitors of ELOVL1. *J. Med. Chem.* **64**(24), 17753-17776 (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, 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