

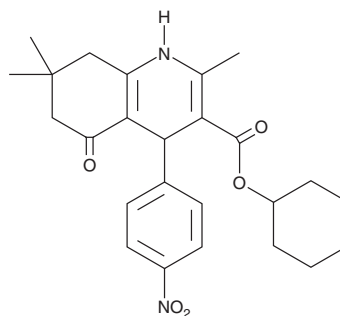
PRODUCT INFORMATION



FLI-06

Item No. 21272

CAS Registry No.: 313967-18-9
Formal Name: 1,4,5,6,7,8-hexahydro-2,7,7-trimethyl-4-(4-nitrophenyl)-5-oxo-3-quinolinecarboxylic acid, cyclohexyl ester
MF: C₂₅H₃₀N₂O₅
FW: 438.5
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 282, 374 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

FLI-06 is supplied as a crystalline solid. A stock solution may be made by dissolving the FLI-06 in the solvent of choice, which should be purged with an inert gas. FLI-06 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of FLI-06 in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

FLI-06 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FLI-06 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. FLI-06 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

FLI-06 is a first-of-its-kind inhibitor of Notch signaling (EC₅₀ = 2.3 μM) that acts upstream of α-secretase and β-secretase cleavage.¹ It disrupts the intracellular trafficking and processing of the Notch signaling pathway, inhibiting general secretion at a stage before exit from the endoplasmic reticulum (ER) and converting ER morphology from tubule to sheet.¹ FLI-06 is reported to alter the maturation pattern of amyloid precursor protein (APP) and prevent the shedding of APPs in HEK293 cells stably expressing a mutated APP that yields robust amounts of amyloid β.¹

Reference

1. Krämer, A., Mentrup, T., Kleizen, B., et al. Small molecules intercept Notch signaling and the early secretory pathway. *Nat. Chem. Biol.* **9**(11), 731-738 (2013).

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, 10/14/2022

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