

PRODUCT INFORMATION



Autogramin-2

Item No. 43556

CAS Registry No.: 2375541-45-8
Formal Name: 6,7-dihydro-2-[[4-(1-methylethoxy)benzoyl]amino]-thiazolo[5,4-c]pyridine-5(4H)-carboxylic acid, 1,1-dimethylethyl ester

MF: C₂₁H₂₇N₃O₄S

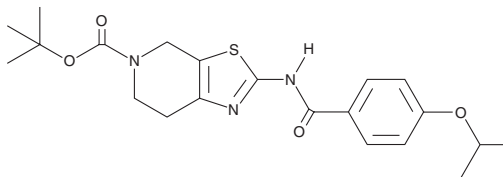
FW: 417.5

Purity: ≥98%

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

Autogramin-2 is supplied as a solid. A stock solution may be made by dissolving the autogramin-2 in the solvent of choice, which should be purged with an inert gas. Autogramin-2 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Autogramin-2 is an inhibitor of Aster-A (IC₅₀ = 1.93 μM), a protein that transports cholesterol from the plasma membrane to the endoplasmic reticulum (ER).¹ Autogramin-2 is selective for Aster-A over Aster-B and -C (IC₅₀s = >20 μM for both). It inhibits autophagy induced by amino acid starvation or rapamycin-mediated mammalian target of rapamycin 1 (mTORC1) inhibition in MCF-7 cells (IC₅₀s = 0.27 and 0.14 μM, respectively).²

References

1. Whitmarsh-Everiss, T., Olsen, A.H., and Laraia, L. Identification of inhibitors of cholesterol transport proteins through the synthesis of a diverse, sterol-inspired compound collection. *Angew. Chem. Int. Ed. Engl.* **60(51)**, 26755-26761 (2021).
2. Laraia, L., Friese, A., Corkery, D.P., et al. The cholesterol transfer protein GRAMD1A regulates autophagosome biogenesis. *Nat. Chem. Biol.* **15(7)**, 710-720 (2019).

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, 08/08/2025

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