PRODUCT INFORMATION



Liproxstatin-1 analog

Item No. 28591

CAS Registry No.: 1170643-61-4

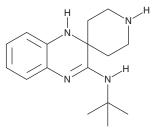
Formal Name: N-(1,1-dimethylethyl)-spiro[piperidine-

4,2'(1'H)-quinoxalin]-3'-amine

MF: $C_{16}H_{24}N_4$ FW: 272.4 **Purity:** ≥98% λ_{max} : 236 nm A crystalline solid UV/Vis.: Supplied as:

-20°C Storage: Stability: ≥2 years

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



Laboratory Procedures

Liproxstatin-1 analog is supplied as a crystalline solid. A stock solution may be made by dissolving the liproxstatin-1 analog in the solvent of choice, which should be purged with an inert gas. Liproxstatin-1 analog is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of liproxstatin-1 analog in these solvents is approximately 5, 15, and 25 mg/ml, respectively.

Description

Liproxstatin-1 analog is an analog of the ferroptosis inhibitor liproxstatin-1 (Item No. 17730).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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.

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, 01/15/2020

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