

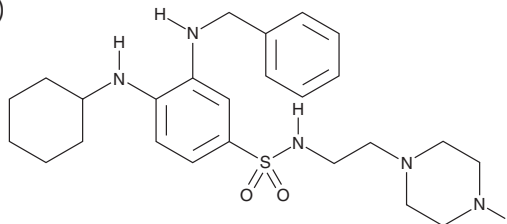
PRODUCT INFORMATION



UAMC-3203

Item No. 26525

CAS Registry No.: 2271358-64-4
Formal Name: 4-(cyclohexylamino)-3-[(phenylmethyl)amino]-N-[2-(1-piperazinyl)ethyl]benzenesulfonamide
MF: C₂₅H₃₇N₅O₂S
FW: 471.7
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 286 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

UAMC-3203 is supplied as a crystalline solid. A stock solution may be made by dissolving the UAMC-3203 in the solvent of choice, which should be purged with an inert gas. UAMC-3203 is soluble in the organic solvent ethanol.

Description

UAMC-3203 is an inhibitor of ferroptosis that has an IC₅₀ value of 10 nM for inhibition of erastin-induced ferroptosis in IMR-32 neuroblastoma cells.¹ It decreases iron-induced plasma lactate dehydrogenase (LDH) levels in a mouse model of acute iron poisoning when administered at a dose of 20 μmol/kg. It is not toxic to mice following chronic administration of a 20 μmol/kg dose for four weeks. UAMC-3203 has increased solubility and a longer half-life in mouse, rat, and human microsomes and isolated plasma than the ferroptosis inhibitor ferrostatin-1 (Item No. 17729). In an *in silico* membrane dynamics study, UAMC-3203 was incorporated into a phospholipid bilayer.

Reference

1. Devisscher, L., Van Coillie, S., Hofmans, S., *et al.* Discovery of novel, drug-like ferroptosis inhibitors with in vivo efficacy. *J. Med. Chem.* **61**(22), 10126-10140 (2018).

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, 06/04/2021

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