

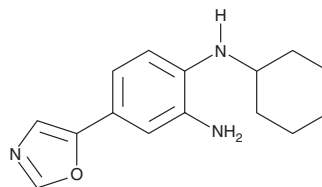
PRODUCT INFORMATION



UAMC-4821

Item No. 43146

Formal Name: N¹-cyclohexyl-4-(oxazol-5-yl)benzene-1,2-diamine
MF: C₁₅H₁₉N₃O
FW: 257.3
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

UAMC-4821 is supplied as a solid. A stock solution may be made by dissolving the UAMC-4821 in the solvent of choice, which should be purged with an inert gas. UAMC-4821 is soluble (≥10 mg/ml) in DMSO.

Description

UAMC-4821 is an inhibitor of ferroptosis.¹ It inhibits ferroptosis induced by ML162 (Item No. 20455) in HT-1080 fibrosarcoma cells (IC₅₀ = 5.2 nM). UAMC-4821 (4 μM) traps lipid radicals induced by di-*tert*-undecylhyponitrite (DTUN; Item No. 32742) in a fluorescence-enabled inhibited autoxidation (FENIX) assay.

Reference

- Scarpellini, C., Klejborowska, G., Lanthier, C., *et al.* Oxazole-based ferroptosis inhibitors with promising properties to treat central nervous system diseases. *J. Med. Chem.* **68**(4), 4908-4928 (2025).

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, 05/22/2025

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