

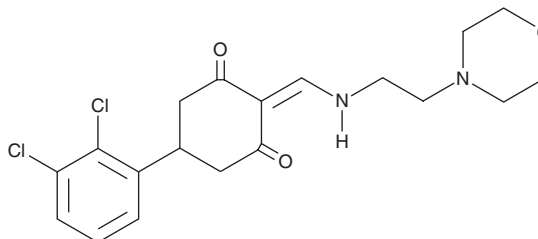
PRODUCT INFORMATION



DC-LC3in-D5

Item No. 38339

CAS Registry No.: 2868312-73-4
Formal Name: 5-(2,3-dichlorophenyl)-2-[[[2-(4-morpholinyl)ethyl]amino]methylene]-1,3-cyclohexanedione
MF: $C_{19}H_{22}Cl_2N_2O_3$
FW: 397.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 257, 305 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

DC-LC3in-D5 is supplied as a crystalline solid. A stock solution may be made by dissolving the DC-LC3in-D5 in the solvent of choice, which should be purged with an inert gas. DC-LC3in-D5 is soluble in chloroform. DC-LC3in-D5 is slightly soluble in methanol.

Description

DC-LC3in-D5 is an inhibitor of LC3 ($IC_{50} = 200$ nM).¹ It selectively binds to LC3-I and LC3-II over GABA type A receptor-associated protein (GABARAP), scavenger receptor class B member 2 (SCARB2), and paraoxonase 2 (PON2) in an activity-based protein profiling (ABPP) assay using HeLa cells. DC-LC3in-D5 (25 μ M) inhibits LC3-II lipidation, autophagosome formation, and the protein-protein interaction between LC3-II and ATG7 in HeLa cells.

Reference

1. Fan, S., Yue, L., Wan, W., *et al.* Inhibition of autophagy by a small molecule through covalent modification of the LC3 protein. *Angew. Chem. Int. Ed.* **60**(50), 26105-26114 (2021).

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, 10/09/2025

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