

PRODUCT INFORMATION



HCy-Lyso

Item No. 41179

Formal Name: (E)-4-(4-(2-(1-ethyl-3,3-dimethylindolin-2-yl)vinyl)phenyl)morpholine

MF: C₂₄H₃₀N₂O

FW: 362.5

Purity: ≥98%

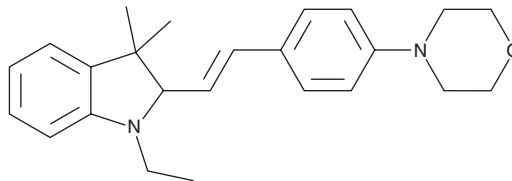
Ex./Em. Max: 510/592 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Light sensitive



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

Laboratory Procedures

HCy-Lyso is supplied as a solid. A stock solution may be made by dissolving the HCy-lyso in the solvent of choice, which should be purged with an inert gas. HCy-Lyso is sparingly soluble (1-10 mg/ml) in chloroform.

Description

HCy-Lyso is a lysosome-specific fluorescent reporter for the detection of hydroxyl radicals.¹ It selectively fluoresces in low pH environments in the presence of hydroxy radicals over hypochlorite, superoxide, nitric oxide, or peroxynitrite radicals. HCy-Lyso (10 μM) displays lysosomal fluorescence in the presence of the ferroptosis inducers erastin (Item No. 17754) or (1S,3R)-RSL3 (Item No. 19288), an effect that can be blocked by the ferroptosis inhibitor ferrostatin-1 (Item No. 17729), in 4T1 cells. HCy-Lyso displays excitation/emission maxima of 510/592 nm, respectively, in low pH environments in the presence of hydroxy radicals.

Reference

1. Zhong, L., Fu, D., Xu, J., *et al.* Rational design of a lysosome-targeted fluorescent probe for monitoring the generation of hydroxyl radicals in ferroptosis pathways. *RSC Adv.* **14**(18), 12864-12872 (2024).

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, 07/01/2024

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