

PRODUCT INFORMATION



Cy-OH-LP

Item No. 42571

Formal Name: (E)-2-((E)-2-(2-chloro-3-((E)-2-(1,3,3-trimethylindolin-2-yl)vinyl)cyclohex-2-en-1-ylidene)ethylidene)-1,3,3-trimethylindoline

Synonym: Cy-hydroxy-LP

MF: C₃₂H₃₇ClN₂

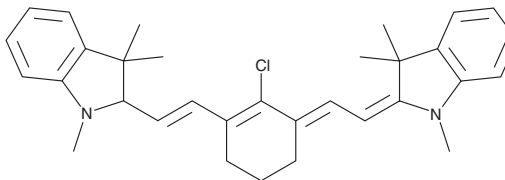
FW: 485.1

Purity: ≥95%

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

Cy-OH-LP is supplied as a solid. A stock solution may be made by dissolving the Cy-OH-LP in the solvent of choice, which should be purged with an inert gas. Cy-OH-LP is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Cy-OH-LP is a turn-on fluorescent and photoacoustic probe for the detection of lipid-associated hydroxyl radicals.¹ In the presence of hydroxyl radicals, Cy-OH-LP is oxidized to Cy7, which displays weak fluorescent and photoacoustic signals in water-soluble environments and strong fluorescent and photoacoustic signals in lipid-rich environments due to its lipid solubility. Cy-OH-LP (1 mg/kg) has been used for photoacoustic imaging of hydroxyl radicals in atherosclerotic plaques in *ApoE*^{-/-} mice fed a high-fat diet.

References

1. Yang, F.-K., Cao, J., Zhang, T., *et al.* Dual-activated photoacoustic probe for reliably detecting hydroxyl radical in ischemic cardiovascular disease in mouse and human samples. *ACS Sens.* **9(10)**, 5445-5453 (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, 02/04/2025

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