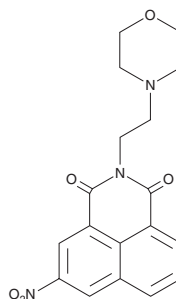


PRODUCT INFORMATION



LysoFP-NO₂
Item No. 27024

CAS Registry No.: 69408-75-9
Formal Name: 2-[2-(4-morpholinyl)ethyl]-5-nitro-1H-benz[de]isoquinoline-1,3(2H)-dione
MF: C₁₈H₁₇N₃O₅
FW: 355.3
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 275, 332 nm
EX./Em. Max: 440/528 nm
Supplied as: A crystalline 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

LysoFP-NO₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the lysoFP-NO₂ in the solvent of choice. LysoFP-NO₂ is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

Description

LysoFP-NO₂ is a turn-on fluorescent probe for carbon monoxide (CO) that localizes to the lysosome.¹ In the presence of lysosomal CO, LysoFP-NO₂ is transformed into LysoFP-NH₂, which is highly fluorescent. LysoFP-NO₂ is selective for CO over various reactive nitrogen, oxygen, and sulfur species. It displays excitation/emission maxima of 440/528 nm, respectively, and is not cytotoxic to HepG2 cells for up to five hours when used at a concentration of 30 μM.

Reference

1. Dhara, K., Lohar, S., Patra, A., *et al.* A new lysosome-targetable turn-on fluorogenic probe for carbon monoxide imaging in living cells. *Anal. Chem.* **90(4)**, 2933-2938 (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, 02/16/2024

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