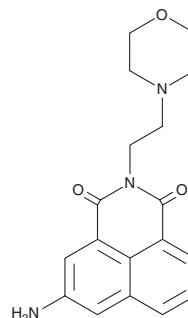


PRODUCT INFORMATION



LysoFP-NH₂ Item No. 27023

CAS Registry No.: 69408-85-1
Formal Name: 5-amino-2-[2-(4-morpholinyl)ethyl]-1H-benz[de]isoquinoline-1,3(2H)-dione
MF: C₁₈H₁₉N₃O₃
FW: 325.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 250, 343, 423 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-NH₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the lysoFP-NH₂ in the solvent of choice. LysoFP-NH₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of lysoFP-NH₂ in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

LysoFP-NH₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lysoFP-NH₂ should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LysoFP-NH₂ has a solubility of approximately 0.03 mg/ml in a 1:30 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LysoFP-NH₂ is the fluorescent form of the lysosomal turn-on fluorescent probe for carbon monoxide (CO) lysoFP-NO₂ (Item No. 27024).¹ LysoFP-NH₂ is formed when lysoFP-NO₂ reacts with CO. It displays excitation/emission maxima of 440/528 nm, respectively.

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