

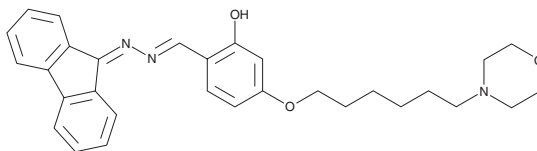
PRODUCT INFORMATION



AIE-LysoY

Item No. 40641

CAS Registry No.: 2056920-87-5
Formal Name: 2-hydroxy-4-[[6-(4-morpholinyl)hexyl]oxy]-benzaldehyde, 2-(9H-fluoren-9-ylidene)hydrazone
MF: C₃₀H₃₃N₃O₃
FW: 483.6
Purity: ≥95%
Ex./Em. Max: 390/565 nm
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.

Description

AIE-LysoY is a lysosome-specific fluorescent probe.¹ It displays excitation/emission maxima of 390/565 nm, respectively. AIE-LysoY has been used to visualize lysosome-autophagosome fusion induced by rapamycin (Item No. 13346) in HeLa cells.

Reference

1. Leung, C.W.T., Wang, Z., Zhao, E., *et al.* A lysosome-targeting AIEgen for autophagy visualization. *Adv. Healthc. Mater.* **5**(4), 427-431 (2016).

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, 10/01/2025

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