

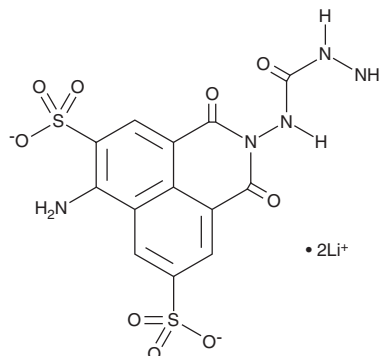
PRODUCT INFORMATION



Lucifer Yellow CH

Item No. 25573

CAS Registry No.: 67769-47-5
Formal Name: 6-amino-2-[(hydrazinylcarbonyl)amino]-2,3-dihydro-1,3-dioxo-1H-benz[de]isoquinoline-5,8-disulfonic acid, dilithium salt
Synonym: Lucifer Yellow carbohydrazide
MF: $C_{13}H_9N_5O_9S_2 \cdot 2Li$
FW: 457.2
Purity: $\geq 80\%$
Ex./Em. Max: 430/540 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Lucifer yellow CH is supplied as a solid. A stock solution may be made by dissolving the lucifer yellow CH in water. The solubility of lucifer yellow CH in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Lucifer yellow CH is a fluorescent probe that contains a free hydrazido group that reacts with aliphatic aldehydes at room temperature.¹ Lucifer yellow CH displays excitation/emission maxima of 430/540 nm, respectively. It has been used as a biological tracer to monitor the branching pattern and regeneration of neurons, gap junction detection and characterization, and selective ablation of cells following aldehyde fixation.^{1,2}

References

1. Stewart, W.W. Lucifer dyes—highly fluorescent dyes for biological tracing. *Nature* **292**(5818), 17-21 (1981).
2. Hanani, M. Lucifer yellow – an angel rather than the devil. *J. Cell. Mol. Med.* **16**(1), 22-31 (2012).

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, 11/29/2022

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