

PRODUCT INFORMATION

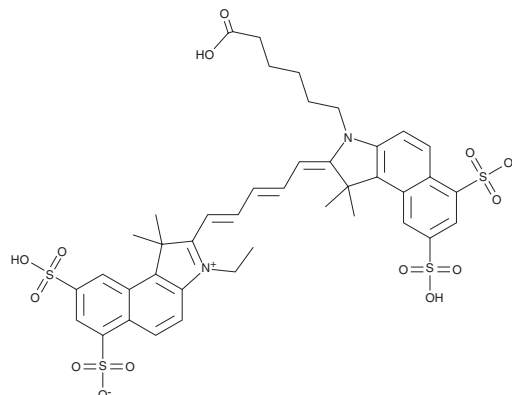


Sulfo-Cyanine 5.5

Item No. 37369

CAS Registry No.: 210892-23-2
Formal Name: 2-[5-[3-(5-carboxypentyl)-1,3-dihydro-1,1-dimethyl-6,8-disulfo-2H-benz[e]indol-2-ylidene]-1,3-pentadien-1-yl]-3-ethyl-1,1-dimethyl-6,8-disulfo-1H-benz[e]indolium, inner salt

Synonym: Cy5.5
MF: C₄₁H₄₄N₂O₁₄S₄
FW: 917.1
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Ex./Em. Max: 675/694 nm



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

Laboratory Procedures

Sulfo-cyanine 5.5 is supplied as a solid. A stock solution may be made by dissolving the sulfo-cyanine 5.5 in the solvent of choice, which should be purged with an inert gas. Sulfo-cyanine 5.5 is soluble in acetonitrile and DMSO (sonicated).

Description

Sulfo-cyanine 5.5 is an amine-reactive fluorescent probe.¹ It reacts with amine groups to form conjugates with peptides and proteins for detection by various fluorescent-based applications.^{1,2} It has been used in cancer-targeted nanovesicles encapsulating perfluorooctyl bromide (PFOB) to identify their localization in mice.² Sulfo-cyanine 5.5 displays excitation/emission maxima of 675/694 nm, respectively.³

References

1. Josephson, L., Kircher, M.F., Mahmood, U., *et al.* Near-infrared fluorescent nanoparticles as combined MR/optical imaging probes. *Bioconjug. Chem.* **13(3)**, 554-560 (2002).
2. Wang, R., Yao, Y., Gao, Y., *et al.* CD133-targeted hybrid nanovesicles for fluorescent/ultrasonic imaging-guided HIFU pancreatic cancer therapy. *Int. J. Nanomedicine* **18**, 2539-2552 (2023).
3. Kwon, Y.-D., Oh, J.-M., Chun, S., *et al.* Synthesis and evaluation of multivalent nitroimidazole-based near-infrared fluorescent agents for neuroblastoma and colon cancer imaging. *Bioorg. Chem.* **113**, 104990 (2021).

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, 11/08/2023

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