

PRODUCT INFORMATION

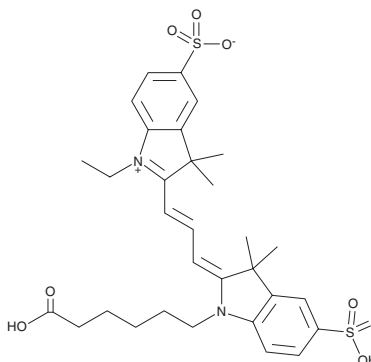


Cy3

Item No. 30392

CAS Registry No.: 146368-13-0
Formal Name: 2-[3-[1-(5-carboxypentyl)-1,3-dihydro-3,3-dimethyl-5-sulfo-2H-indol-2-ylidene]-1-propen-1-yl]-1-ethyl-3,3-dimethyl-5-sulfo-3H-indolium, inner salt

MF: C₃₁H₃₈N₂O₈S₂
FW: 630.8
Purity: ≥98%
UV/Vis.: λ_{max}: 525, 558 nm
Ex./Em. Max 550/570 nm, respectively
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

Cy3 is supplied as a crystalline solid. A stock solution may be made by dissolving the Cy3 in the solvent of choice, which should be purged with an inert gas. Cy3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Cy3 in these solvents is approximately 1, 15, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Cy3 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Cy3 in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cy3 is a cyanine-containing fluorochrome.¹ It has commonly been conjugated to secondary antibodies for use in immunocytochemistry and immunohistochemistry applications. Cy3 displays excitation/emission maxima of 550/570 nm, respectively.

Reference

1. Suzuki, T., Matsuzaki, T., Hagiwara, H., *et al.* Recent advances in fluorescent labeling techniques for fluorescence microscopy. *Acta Histochem. Cytochem.* **40(5)**, 131-137 (2007).

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, 10/14/2022

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