

PRODUCT INFORMATION



Cy3-SE

Item No. 30393

CAS Registry No.: 146368-16-3
Formal Name: 2-[3-[1-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-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

Synonyms: Cy3 NHS ester, N-hydroxysuccinimidyl Ester Cy3

MF: C₃₅H₄₁N₃O₁₀S₂

FW: 727.8

Purity: ≥95%

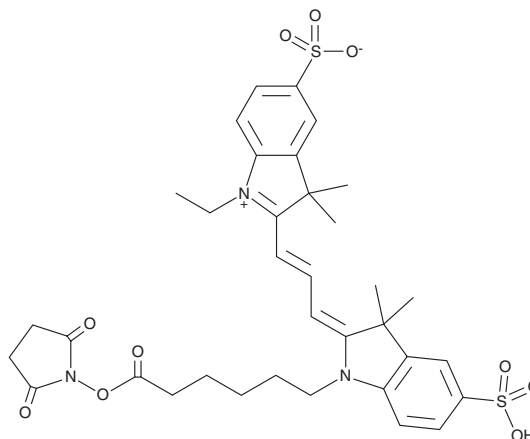
UV/Vis.: λ_{max}: 524, 557 nm

Ex./Em. Max.: 550/570 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

Cy3-SE is supplied as a crystalline solid. A stock solution may be made by dissolving the Cy3-SE in the solvent of choice, which should be purged with an inert gas. Cy3-SE is soluble in DMSO.

Description

Cy3-SE is a hydrophilic amine-reactive fluorescent probe.^{1,2} It displays excitation/emission maxima of 550/570 nm, respectively. It has commonly been used to label nucleic acids.

References

1. Randolph, J.B. and Waggoner, A.S. Stability, specificity and fluorescence brightness of multiply-labeled fluorescent DNA probes. *Nucleic Acids Res.* **25(14)**, 2923-2929 (1997).
2. Harvey, B.J. and Levitus, M. Nucleobase-specific enhancement of Cy3 fluorescence. *J. Fluoresc.* **19(3)**, 443-448 (2009).

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, 11/23/2022

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