

PRODUCT INFORMATION

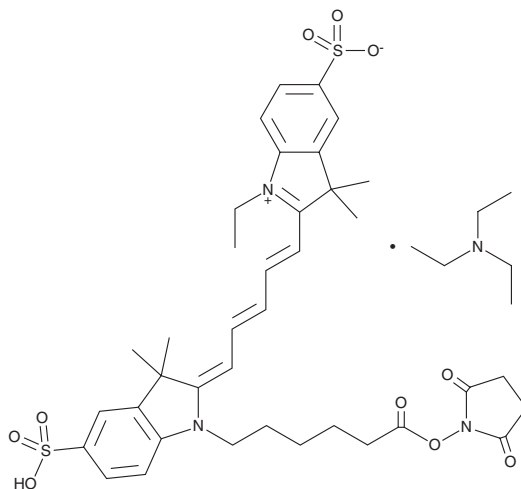


Cy5-SE (triethylamine salt)

Item No. 30387

CAS Registry No.: 1497420-70-8
Formal Name: 2-[5-[1-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-1,3-dihydro-3,3-dimethyl-5-sulfo-2H-indol-2-ylidene]-1,3-pentadien-1-yl]-1-ethyl-3,3-dimethyl-5-sulfo-3H-indolium, inner salt, compd. with N,N-diethylethanamine

Synonym: Sulfo-Cy5
MF: C₃₇H₄₃N₃O₁₀S₂ • C₆H₁₅N
FW: 855.1
Purity: ≥95%
Ex./Em. Max: 646/662 nm, respectively
UV/Vis.: λ_{max}: 649 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

Cy5-SE (triethylamine salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Cy5-SE (triethylamine salt) in the solvent of choice, which should be purged with an inert gas. Cy5-SE (triethylamine salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Cy5-SE (triethylamine salt) in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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 Cy5-SE (triethylamine salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Cy5-SE (triethylamine salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cy5-SE is a hydrophilic amine-reactive fluorescent probe.¹ It displays excitation/emission maxima of 646/662 nm, respectively. Cy5-SE-conjugated ligands have been used in the characterization of hematopoietic tumor microenvironments.

Reference

1. Soodgupta, D., Zhou, H., Beaino, W., *et al.* Ex vivo and in vivo evaluation of overexpressed VLA-4 in multiple myeloma using LLP2A imaging agents. *J. Nucl. Med.* **57(4)**, 640-645 (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, 10/11/2022

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