

PRODUCT INFORMATION



Cy5-Yne (triethylamine salt)

Item No. 40134

Formal Name: 2-[5-[1,3-dihydro-3,3-dimethyl-1-[6-oxo-6-(2-propyn-1-ylamino)hexyl]-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-alkyne

MF: $C_{36}H_{43}N_3O_7S_2 \cdot C_6H_{15}N$

FW: 795.1

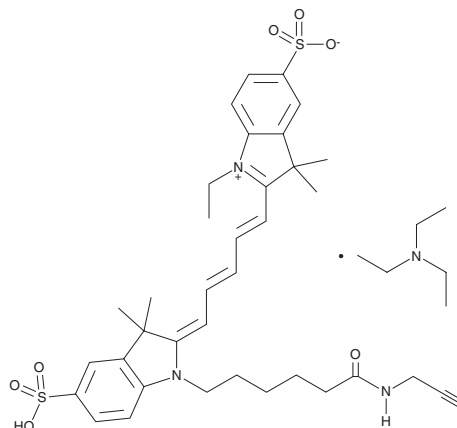
Purity: $\geq 95\%$

Ex./Em. Max: 650/680 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

Cy5-Yne (triethylamine salt) is supplied as a solid. A stock solution may be made by dissolving the Cy5-Yne (triethylamine salt) in the solvent of choice, which should be purged with an inert gas. Cy5-Yne (triethylamine salt) is soluble in DMSO. Cy5-Yne (triethylamine salt) is slightly soluble in acetonitrile.

Cy5-Yne (triethylamine salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Cy5-Yne is a clickable azido-reactive fluorescent probe.¹ It displays excitation/emission maxima of 650/680 nm, respectively. When conjugated to peptides, Cy5-Yne has been used as a pH-independent fluorescent probe for the activity of lysosomal cathepsin proteases in *S. typhimurium*-infected and bystander bone marrow-derived macrophages (BMDMs).

Reference

1. Sanman, L.E., van der Linden, W.A., Verdoes, M., *et al.* Bifunctional probes of cathepsin protease activity and pH reveal alterations in endolysosomal pH during bacterial infection. *Cell Chem. Biol.* **23(7)**, 793-804 (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, 04/02/2024

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