

# PRODUCT INFORMATION



## Cy7-SE

Item No. 30753

**CAS Registry No.:** 477908-53-5  
**Formal Name:** 2-[7-[1-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-1,3-dihydro-3,3-dimethyl-5-sulfo-2H-indol-2-ylidene]-1,3,5-heptatrien-1-yl]-1-ethyl-3,3-dimethyl-5-sulfo-3H-indolium

**Synonyms:** Cy7 NHS Ester,  
N-hydroxysuccinimidyl Ester Cy7

**MF:** C<sub>39</sub>H<sub>45</sub>N<sub>3</sub>O<sub>10</sub>S<sub>2</sub>

**FW:** 779.9

**Purity:** ≥98%

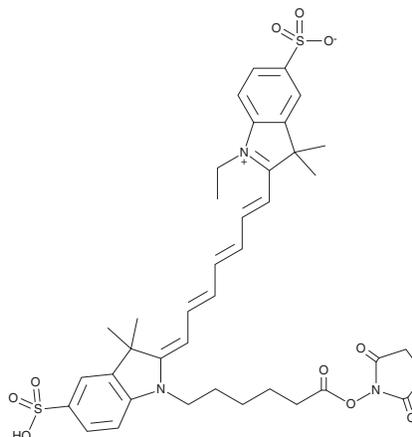
**UV/Vis.:** λ<sub>max</sub>: 755, 758 nm

**Ex./Em. Max:** 747/774 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

Cy7-SE is supplied as a crystalline solid. A stock solution may be made by dissolving the Cy7-SE in the solvent of choice, which should be purged with an inert gas. Cy7-SE is soluble in ethanol, DMSO, and dimethyl formamide.

### Description

Cy7-SE is a hydrophilic amine-reactive fluorescent probe.<sup>1</sup> It displays excitation/emission maxima of 747/774 nm, respectively.<sup>2</sup> Cy7-SE has commonly been used in the labeling and detection of nucleic acids, antibodies, and proteins.

### References

1. Berlier, J.E., Rothe, A., Buller, G., *et al.* Quantitative comparison of long-wavelength alexa fluor dyes to Cy dyes: Fluorescence of the dyes and their bioconjugates. *J. Histochem. Cytochem.* **51(12)**, 1699-1712 (2003).
2. Jiang, C., Cai, H., Peng, X., *et al.* Targeted imaging of tumor-associated macrophages by cyanine 7-labeled mannose in xenograft tumors. *Mol. Imaging* **16**, 1536012116689499 (2017).

#### 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, 12/16/2022

#### 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