

PRODUCT INFORMATION



Cy5 NHS ester

Item No. 45624

CAS Registry No.: 1263093-76-0
Formal Name: 2-[5-(1,3-dihydro-1,3,3-trimethyl-2H-indol-2-ylidene)-1,3-pentadien-1-yl]-1-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-3,3-dimethyl-3H-indolium, monotetrafluoroborate

Synonyms: Cy5 N-hydroxysuccinimide ester, Cyanine 5 NHS ester

MF: $C_{36}H_{42}N_3O_4 \cdot BF_4$

FW: 667.6

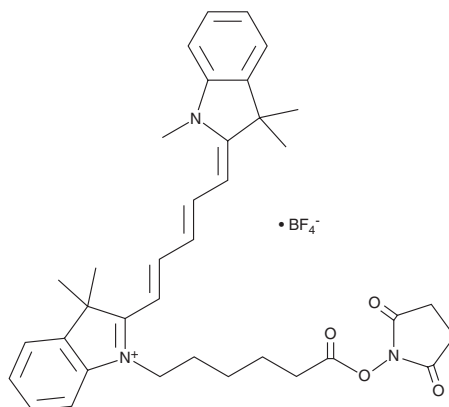
Purity: $\geq 95\%$

Ex./Em. Max: 646/662 nm

Supplied as: A 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 NHS ester is supplied as a solid. A stock solution may be made by dissolving the Cy5 NHS ester in the solvent of choice, which should be purged with an inert gas. Cy5 NHS ester is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Cy5 NHS ester is an amine-reactive fluorochrome. It displays excitation/emission maxima of 646/662 nm, respectively. Cy5 NHS ester has been conjugated to a peptide motif based on a Factor VII-activating protease (FSAP) substrate for use as an activity-based probe for the FSAP serine protease domain (SPD).¹

Reference

1. Rut, W., Nielsen, N.V., Czarna, J., *et al.* Fluorescent activity-based probe for the selective detection of Factor VII activating protease (FSAP) in human plasma. *Thromb. Res.* **182**, 124-132 (2019).

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, 05/27/2026

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