

PRODUCT INFORMATION

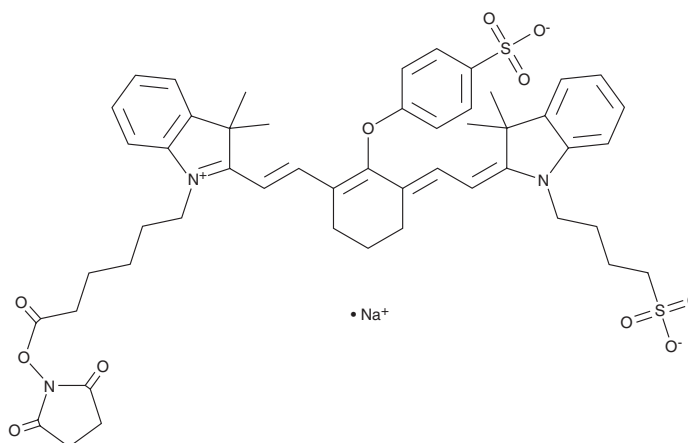


iFluor™ IRDye 800RS NHS Ester

Item No. 44614

CAS Registry No.: 918324-94-4
Formal Name: 2-[2-[3-[2-[1,3-dihydro-3,3-dimethyl-1-(4-sulfobutyl)-2H-indol-2-ylidene]ethylidene]-2-(4-sulfophenoxy)-1-cyclohexen-1-yl]ethenyl]-1-[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]-3,3-dimethyl-3H-indolium, inner salt, monosodium salt

MF: $C_{50}H_{56}N_3O_{11}S_2 \cdot Na$
FW: 962.1
Purity: $\geq 90\%$
Ex./Em. Max: 767/785 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

iFluor™ IRDye 800RS NHS ester is supplied as a solid. A stock solution may be made by dissolving the iFluor™ IRDye 800RS NHS ester in the solvent of choice, which should be purged with an inert gas. iFluor™ IRDye 800RS NHS ester is soluble (≥ 10 mg/ml) in DMSO.

Description

iFluor™ IRDye 800RS NHS ester is a near-infrared amine-reactive fluorescent dye. It displays excitation/emission maxima of 767/785 nm. iFluor™ IRDye 800RS NHS ester can be used to label nucleic acids for use in fluorescence-based applications, including hybridization assays and optical imaging.

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, 11/06/2025

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