

PRODUCT INFORMATION



iFluor™ IRD 800 CW SE

Item No. 44618

Formal Name: 2-[2-[3-[2-[1,3-dihydro-3,3-dimethyl-5-sulfo-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-5-sulfo-3H-indolium, inner salt

MF: $C_{50}H_{54}N_3O_{17}S_4 \cdot 3C_6H_{16}N$

FW: 1,403.8

Purity: $\geq 90\%$

Ex./Em. Max: 774/792 nm

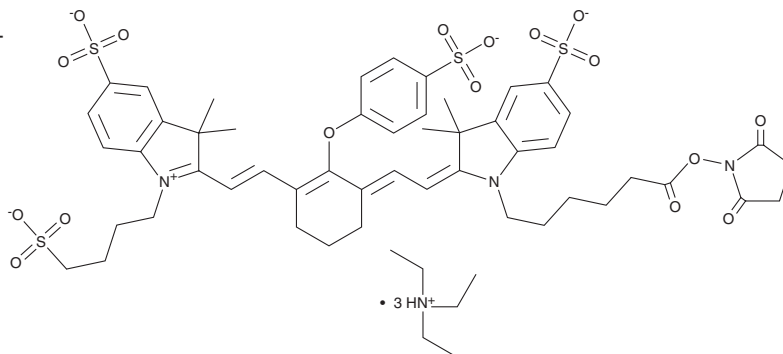
Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Light sensitive

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



Description

iFluor™ IRD 800 CW SE is an amine-reactive near-infrared fluorescent probe. It displays excitation/emission maxima of 774 and 792 nm, respectively. iFluor™ IRD 800 CW SE can be used to label peptides, antibodies, and amine-modified oligonucleotides for use in flow cytometry (FC), immunocytochemistry (ICC), microscopy, protein array, and Western blot (WB) applications. Formulations containing iFluor™ IRD 800 CW SE-conjugated antibodies have been used as imaging agents in fluorescence-guided surgery in patients with cancer.

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, 10/21/2025

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