

PRODUCT INFORMATION



Fluorescent SM-102

Item No. 43088

Formal Name: heptadecan-9-yl
8-((2-hydroxyethyl)
(6-((6-((7-nitrobenzo[c][1,2,5]
oxadiazol-4-yl)amino)hexanoyl)
oxy)hexyl)amino)octanoate

Synonym: NBD-SM-102

MF: $C_{45}H_{79}N_5O_8$

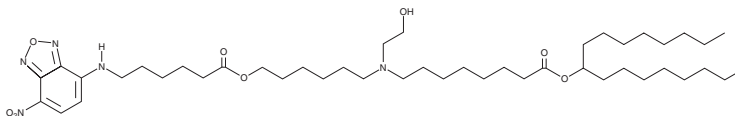
FW: 818.2

Purity: $\geq 95\%$

Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

Fluorescent SM-102 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol, chloroform, and DMSO purged with an inert gas can be used. Fluorescent SM-102 is slightly soluble (0.1-1 mg/ml) in methanol, chloroform, and DMSO.

Description

Fluorescent SM-102 is a derivative of SM-102 (Item No. 33474) containing the fluorescent probe nitrobenzofurazan (NBD).

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, 05/12/2026

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