

PRODUCT INFORMATION

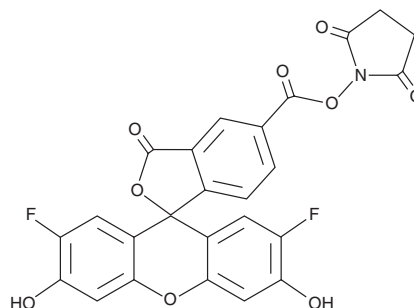


Oregon Green™ 488 succinimidyl ester

Item No. 43466

CAS Registry No.: 198139-51-4
Formal Name: 2',7'-difluoro-3',6'-dihydroxy-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-5-carboxylic acid, 2,5-dioxo-1-pyrrolidinyl ester
Synonyms: 2',7'-Difluorofluorescein succinimidyl ester, OG 488 SE, OG 488 succinimidyl ester

MF: C₂₅H₁₃F₂NO₉
FW: 509.4
Purity: ≥85%
UV/Vis.: λ_{max}: 214, 278, 458 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

Oregon Green™ 488 succinimidyl ester is supplied as a solid. A stock solution may be made by dissolving the Oregon Green™ 488 succinimidyl ester in the solvent of choice, which should be purged with an inert gas. Oregon Green™ 488 succinimidyl ester is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Oregon Green™ 488 succinimidyl ester is an amine-reactive fluorescent probe.¹ It has been used in the synthesis of fluorescent DNA probes and to fluorescently label proteins.^{1,2} A related derivative, Oregon Green™ 488 (Item No. 35372), displays absorbance/emission maxima of 490/514 nm, respectively.³

References

1. Bordello, J., Sánchez, M.I., Vázquez, M.E., *et al.* Single-molecule approach to DNA minor-groove association dynamics. *Angew. Chem. Int. Ed. Engl.* **51(30)**, 7541-7544 (2012).
2. Lehto, M.T. and Sharom, F.J. Proximity of the protein moiety of a GPI-anchored protein to the membrane surface: A FRET study. *Biochemistry* **41(26)**, 8386-8376 (2002).
3. Sun, W.-C., Gee, K.R., Klaubert, D.H., *et al.* Synthesis of fluorinated fluoresceins. *J. Org. Chem.* **62(19)**, 6469-6475 (1997).

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, 02/03/2026

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