

PRODUCT INFORMATION



N-Butylfluorescein

Item No. 20909

CAS Registry No.: 335193-91-4

Formal Name: 3'-butoxy-6'-hydroxy-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one

MF: $C_{24}H_{20}O_5$

FW: 388.4

Purity: $\geq 98\%$

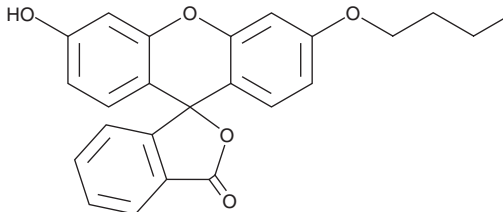
Ex./Em. Max: 467/512 nm

UV/Vis.: λ_{max} : 226, 277 nm

Supplied as: A crystalline 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

N-Butylfluorescein is supplied as a crystalline solid. A stock solution may be made by dissolving the N-butylfluorescein in the solvent of choice, which should be purged with an inert gas. N-Butylfluorescein is slightly soluble in methanol and chloroform.

Description

N-Butylfluorescein is a fluorescent compound that displays excitation/emission maxima of 467/512 nm, respectively.¹ It has been used in the synthesis of butyl fluorescein myo-inositol phosphate (butyl FLIP), a fluorogenic substrate of phosphatidylinositol-specific phospholipase C (PI-PLC).²

References

1. Li, X. and Taylor, J.S. General strategy for the preparation of membrane permeable fluorogenic peptide ester conjugates for in vivo studies of ester prodrug stability. *Bioorg. Med. Chem.* **12**(3), 545-552 (2004).
2. Zaikova, T.O., Rukavishnikov, A.V., Birrell, G.B., *et al.* Synthesis of fluorogenic substrates for continuous assay of phosphatidylinositol-specific phospholipase C. *Bioconjug. Chem.* **12**(2), 307-313 (2001).

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, 12/16/2022

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