

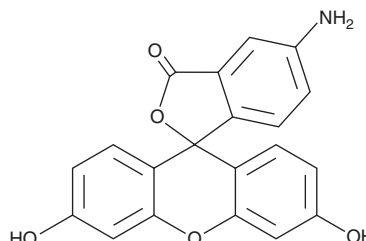
PRODUCT INFORMATION



5-Aminofluorescein

Item No. 36270

CAS Registry No.: 3326-34-9
Formal Name: 5-amino-3',6'-dihydroxy-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one
Synonyms: 5-AF, Fluoresceinamine isomer 1
MF: C₂₀H₁₃NO₅
FW: 347.3
Purity: ≥95%
UV/Vis.: λ_{max}: 226 nm
Ex./Em. Max: 490/515 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

5-Aminofluorescein is supplied as a solid. A stock solution may be made by dissolving the 5-aminofluorescein in the solvent of choice, which should be purged with an inert gas. 5-Aminofluorescein is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5-aminofluorescein in these solvents is approximately 5 mg/ml.

Description

5-Aminofluorescein is a fluorescent compound and an inhibitor of the mRNA N⁶-methyladenosine (m⁶A) demethylase fat mass and obesity-associated (FTO) protein (IC₅₀ = 6.55 μM).^{1,2} It displays excitation/emission maxima of 490 and 515 nm, respectively, and has been conjugated to human serum albumin for fluorescence imaging of glioblastomas.^{1,3}

References

1. Al-Natour, M.A., Yousif, M.D., Cavanagh, R., *et al.* Facile dye-initiated polymerization of lactide-glycolide generates highly fluorescent poly(lactic-co-glycolic acid) for enhanced characterization of cellular delivery. *ACS Macro Lett.* **9**(3), 431-437 (2020).
2. Wang, T., Hong, T., Huang, Y., *et al.* Fluorescein derivatives as bifunctional molecules for the simultaneous inhibiting and labeling of FTO protein. *J. Am. Chem. Soc.* **137**(43), 13736-13739 (2015).
3. Ding, R., Frei, E., Fardanesh, M., *et al.* Pharmacokinetics of 5-aminofluorescein-albumin, a novel fluorescence marker of brain tumors during surgery. *J. Clin. Pharmacol.* **51**(5), 672-678 (2011).

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, 11/11/2022

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