

PRODUCT INFORMATION



5-(4,6-Dichloro-s-triazin-2-ylamino)fluorescein (hydrochloride)

Item No. 34080

CAS Registry No.: 21811-74-5
Formal Name: 5-[(4,6-dichloro-1,3,5-triazin-2-yl)amino]-3',6'-dihydroxy-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one, monohydrochloride

Synonyms: Dichlorotriazinylaminofluorescein, DTAF

MF: C₂₃H₁₂Cl₂N₄O₅ • HCl

FW: 531.7

Purity: ≥98%

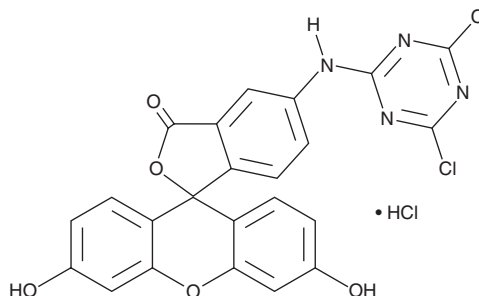
UV/Vis.: λ_{max}: 211, 225, 277 nm

Ex./Em. Max: 492/513 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-(4,6-Dichloro-s-triazin-2-ylamino)fluorescein (DTAF) (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the DTAF (hydrochloride) in the solvent of choice, which should be purged with an inert gas. DTAF (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DTAF (hydrochloride) in these solvents is approximately 1 mg/ml.

Description

DTAF is an amine-reactive fluorescent probe for labeling proteins.¹ It displays excitation/emission maxima of 492/513 nm, respectively. DTAF has been conjugated to IgG for use as fluorescent antibodies and as a pre-column or on-column reagent to label amino acids and amino acid-containing compounds.¹⁻³ It has also been used as a derivatizing agent in the analysis of ephedrine and pseudoephedrine from various sources.³

References

1. Bakeslee, D. and Baines, M.G. Immunofluorescence using dichlorotriazinylaminofluorescein (DTAF). I. Preparation and fractionation of labelled IgG. *J. Immunol. Methods* **13(3-4)**, 305-320 (1976).
2. De Boer, S.H. and Wiczorek, A. Production of monoclonal antibodies to *Corynebacterium sepedonicum*. *Phytopathology* **74(12)**, 1431-1434 (1984).
3. Wang, W., Li, C., Li, Y., et al. Rapid and ultrasensitive determination of ephedrine and pseudoephedrine derivatized with 5-(4,6-dichloro-s-triazin-2-ylamino) fluorescein by micellar electrokinetic chromatography with laser-induced fluorescence detection. *J. Chromatogr. A*. **1102(1-2)**, 273-279 (2005).

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, 11/17/2022

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