

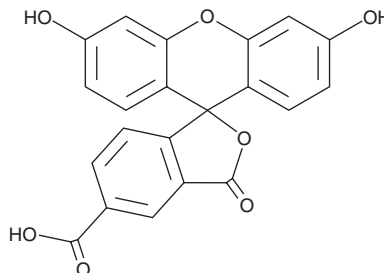
# PRODUCT INFORMATION



## 5-Carboxyfluorescein

Item No. 19581

**CAS Registry No.:** 76823-03-5  
**Formal Name:** 3',6'-dihydroxy-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-5-carboxylic acid  
**Synonym:** 5-FAM  
**MF:** C<sub>21</sub>H<sub>12</sub>O<sub>7</sub>  
**FW:** 376.3  
**Purity:** ≥98%  
**Ex./Em. Max:** 492/518 nm  
**UV/Vis.:** λ<sub>max</sub>: 209, 225, 277 nm  
**Supplied as:** A crystalline 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-Carboxyfluorescein is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-carboxyfluorescein in the solvent of choice, which should be purged with an inert gas. 5-Carboxyfluorescein is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 5-carboxyfluorescein in these solvents is approximately 5, 0.5, and 1 mg/ml, respectively.

5-Carboxyfluorescein is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-carboxyfluorescein should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 5-Carboxyfluorescein has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

5-Carboxyfluorescein is a single isomer derivative of 5(6)-carboxyfluorescein (Item No. 17172) that can be used to fluorescently label biomolecules through the interaction of carboxylic acid with primary amines.<sup>1</sup> It demonstrates excitation/emission maxima of 492 and 518 nm, respectively.

### Reference

1. Fischer, R., Mader, O., Jung, G., *et al.* Extending the applicability of carboxyfluorescein in solid-phase synthesis. *Bioconjugate Chem.* **14**(3), 653-660 (2003).

#### 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, 10/27/2022

#### 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