

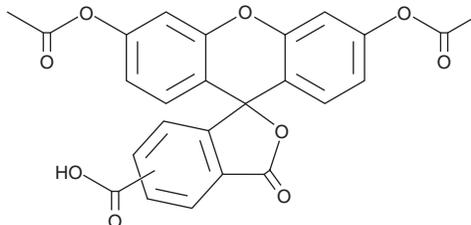
# PRODUCT INFORMATION



## 5(6)-Carboxyfluorescein diacetate

Item No. 19582

**CAS Registry No.:** 124387-19-5  
**Formal Name:** 3',6'-bis(acetyloxy)-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-ar-carboxylic acid  
**Synonym:** CFDA  
**MF:** C<sub>25</sub>H<sub>16</sub>O<sub>9</sub>  
**FW:** 460.4  
**Purity:** ≥98% (mixture of isomers)  
**UV/Vis.:** λ<sub>max</sub>: 220, 291 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(6)-Carboxyfluorescein diacetate (CFDA) is supplied as a crystalline solid. A stock solution may be made by dissolving the CFDA in the solvent of choice. CFDA is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CFDA in these solvents is approximately 30 and 50 mg/ml, respectively.

CFDA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CFDA should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CFDA has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CFDA is a fluorescent probe used to stain cells with intact membranes for analysis in applications such as flow cytometry or fluorescence microscopy. It has relatively slow cellular efflux with a half-life of ~94 minutes and displays fluorescence with excitation/emission maxima of 494/521 nm.<sup>1</sup>

### Reference

1. Dive, C., Cox, H., Watson, J. V., *et al.* Polar fluorescein derivatives as improved substrate probes for flow cytoenzymological assay of cellular esterases. *Mol. Cell Probe* **2**(2), 131-145 (1988).

#### 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, 05/02/2023

#### 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