

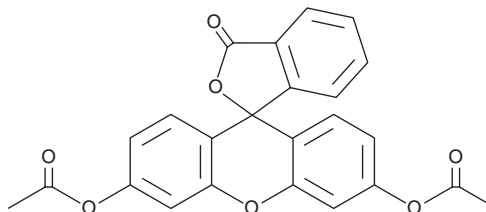
# PRODUCT INFORMATION



## Fluorescein diacetate

Item No. 35701

**CAS Registry No.:** 596-09-8  
**Formal Name:** 3',6'-bis(acetyloxy)-spiro[isobenzofuran-1(3H),9'-(9H)xanthen]-3-one  
**Synonyms:** 3',6'-Diacetylfluorescein, FDA, MFCD 5062, NSC 4726, NSC 667259  
**MF:** C<sub>24</sub>H<sub>16</sub>O<sub>7</sub>  
**FW:** 416.4  
**Purity:** ≥98%  
**Ex./Em. Max:** 490/520 nm  
**UV/Vis.:** λ<sub>max</sub>: 220 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

Fluorescein diacetate is supplied as a solid. A stock solution may be made by dissolving the fluorescein diacetate in the solvent of choice, which should be purged with an inert gas. Fluorescein diacetate is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of fluorescein diacetate in these solvents is approximately 1 and 5 mg/ml, respectively.

### Description

Fluorescein diacetate is a cell-permeable fluorogenic probe to quantify cell viability.<sup>1</sup> It is cleaved by intracellular esterases to produce fluorescein which displays excitation/emission maxima of 490/520 nm, respectively.<sup>2</sup>

### References

1. Brunius, G. Technical aspects of the use of 3',6'-diacetyl fluorescein for vital fluorescent staining of bacteria *Curr. Microbiol.* **4**, 321-323 (1980).
2. 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, 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