

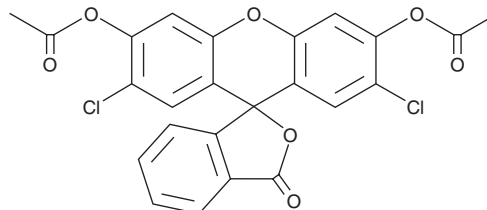
PRODUCT INFORMATION



2',7'-Dichlorofluorescein diacetate

Item No. 20656

CAS Registry No.: 2044-85-1
Formal Name: 3',6'-bis(acetyloxy)-2',7'-dichloro-spiro[isobenzofuran-1(3H),9'-[9H]xanthen]-3-one
Synonym: MFCD 37501
MF: C₂₄H₁₄Cl₂O₇
FW: 485.3
Purity: ≥98%
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

2',7'-Dichlorofluorescein diacetate is supplied as a crystalline solid. A stock solution may be made by dissolving the 2',7'-dichlorofluorescein diacetate in the solvent of choice, which should be purged with an inert gas. 2',7'-Dichlorofluorescein diacetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2',7'-dichlorofluorescein diacetate in ethanol and DMF is approximately 25 mg/ml and approximately 33 mg/ml in DMSO. The solubility of 2',7'-dichlorofluorescein diacetate in 0.1 M Na₂CO₃ is approximately 5 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2',7'-dichlorofluorescein diacetate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2',7'-dichlorofluorescein diacetate in PBS (pH 7.2) is approximately 0.02 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2',7'-Dichlorofluorescein diacetate is a cell-permeable fluorescent dye and an oxidized form of 2,7-dichlorodihydrofluorescein diacetate (Item No. 85155).

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, 09/01/2023

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