

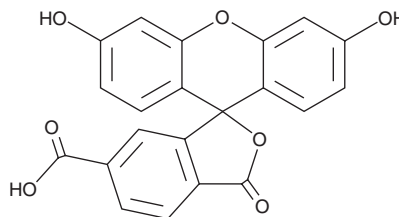
PRODUCT INFORMATION



6-Carboxyfluorescein

Item No. 21233

CAS Registry No.: 3301-79-9
Formal Name: 3',6'-dihydroxy-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-6-carboxylic acid
Synonym: 6-FAM
MF: C₂₁H₁₂O₇
FW: 376.3
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 289 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

6-Carboxyfluorescein (6-FAM) is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-FAM in the solvent of choice. 6-FAM is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 6-FAM in these solvents is approximately 5, 0.5, and 1 mg/ml, respectively.

6-FAM is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 6-FAM should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 6-FAM 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

6-FAM 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.¹ 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**, 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, 12/22/2022

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