

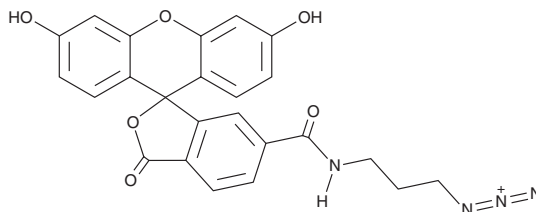
PRODUCT INFORMATION



FAM Azide 6-isomer

Item No. 42389

CAS Registry No.: 1386385-76-7
Formal Name: N-(3-azidopropyl)-3',6'-dihydroxy-3-oxo-spiro[isobenzofuran-1(3H),9'-[9H]xanthene]-6-carboxamide
Synonym: 6-FAM Azide
MF: C₂₄H₁₈N₄O₆
FW: 458.4
Purity: ≥95%
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

FAM azide 6-isomer is supplied as a solid. A stock solution may be made by dissolving the FAM azide 6-isomer in the solvent of choice, which should be purged with an inert gas. FAM azide 6-isomer is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

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 FAM azide 6-isomer can be prepared by directly dissolving the solid in aqueous buffers. FAM azide 6-isomer is soluble (≥10 mg/ml) in PBS (pH 7.2) .

Description

FAM azide 6-isomer is a clickable form of 6-carboxyfluorescein (6-FAM; Item No. 21233).¹ It has been used to label 5-ethynyl-2'-deoxyuridine (EdU; Item No. 20518) for use in cell viability, proliferation, and cell cycle progression assays in cancer cells. A related derivative, 6-FAM, displays excitation/emission maxima of 492/518 nm, respectively.²

References

1. Stengl, A., Hörl, D., Leonhaedt, H., *et al.* A simple and sensitive high-content assay for the characterization of antiproliferative therapeutic antibodies. *SLAS Discov.* **22**(3), 309-315 (2017).
2. Fischer, R., Mader, O., Jung, G., *et al.* Extending the applicability of carboxyfluorescein in solid-phase synthesis. *Bioconjug. 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, 12/10/2024

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