

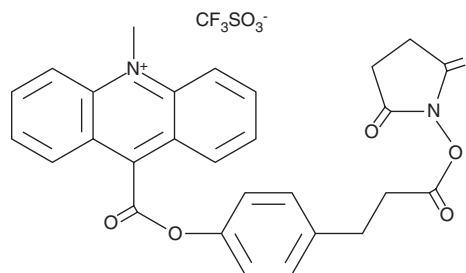
PRODUCT INFORMATION



Acridinium NHS ester

Item No. 200200

CAS Registry No.: 177332-37-5
Formal Name: 9[[4-[3-[(2,5-dioxo-1-pyrrolidinyl)oxy]-3-oxopropyl]phenoxy]carbonyl]-10-methyl-acridinium 1,1,1-trifluoromethanesulfonate
Synonym: Acridinium Compound I
MF: C₂₈H₂₃N₂O₆ • CF₃SO₃
FW: 632.6
Purity: ≥95%
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

Acridinium NHS ester is supplied as a crystalline solid. A stock solution may be made by dissolving the acridinium NHS ester in the solvent of choice, which should be purged with an inert gas. Acridinium NHS ester is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of acridinium NHS ester in these solvents is approximately 25 mg/ml.

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

Description

The acridinium NHS ester can be used to label proteins and nucleic acids. The covalently bound acridinium NHS ester will produce chemiluminescence in the presence of hydrogen peroxide. Acridinium labeled proteins can be used as a detection method in immunoassays.^{1,2}

References

1. Weeks, I., Beheshti, I., McCapra, F., *et al.* Acridinium esters as high-specific-activity labels in immunoassay. *Clin. Chem.* **29**, 1474-1479 (1983).
2. Weeks, I., Sturgess, M., Brown, R.C., *et al.* Immunoassays using acridinium esters. *Methods Enzymol.* **133**, 367-387 (1986).

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, 02/14/2024

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