

Chemiluminescent labeling kit

Chemiluminescent acridinium esters are commonly used to label antibodies, proteins, nucleic acids, and peptides containing primary amine groups. These chemiluminescent molecules are frequently for use as detection reagents in immunoassays as a safe alternative to radioactive tracers. The Chemiluminescent labeling kit provides the necessary components for quick, efficient, and easy acridinium labeling.

Citations: 1

[View Online »](#)

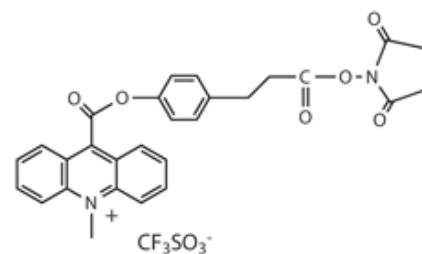
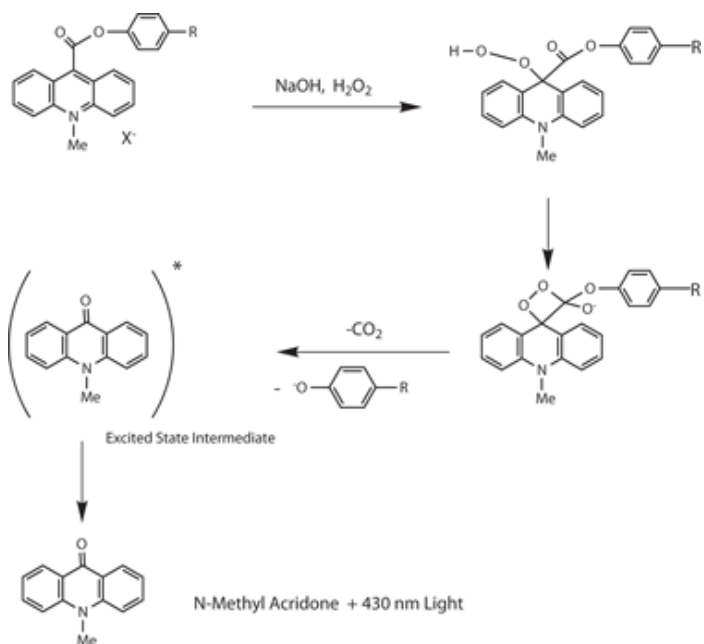
Ordering Information

[Order Online »](#)

ADI-907-001	5Reactions
-------------	------------

Manuals, SDS & CofA

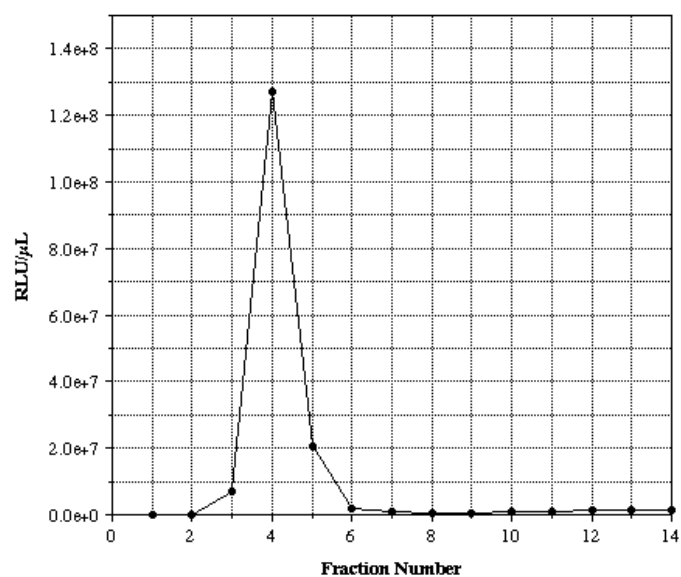
[View Online »](#)



Acridinium C₂ NHS Ester

4-(2-succinimidyl-oxycarbonyl-ethyl)-phenyl-10-acridinium-9-carboxylate trifluoromethyl sulfonate

C₂₉H₂₃F₃N₂O₉S Formula Weight= 632.55



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application

Chemiluminescence

Application Notes

A complete kit for the labeling of antibodies, proteins, peptides, and nucleic acids with Acridinium NHS Ester.

Contents

Acridinium NHS Ester, Dry DMF, Bicarbonate labeling buffer, 10% Lysine, Gel filtration column, Column buffer concentrate, Set of Trigger solutions

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-](#)
uk@enzolifesciences.com