

Acridinium Ester, (trigger solution)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-906-001	2x250ml
-------------	---------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stable for one year upon date of receipt when stored as indicated.

Long Term Storage Ambient

Shipping Ambient Temperature – Dangerous Good

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Chemiluminescent Acridinium Esters are commonly used to label antibodies, proteins, nucleic acids, and some peptides, frequently for use as a detection reagent in immunoassays; Trigger solution for Acridinium NHS Ester.

Formulation Liquid. Solution 1 contains hydrogen peroxide in nitric acid. Solution 2 contains cetyl trimethylammonium chloride in sodium hydroxide.

Source Species Independent

Technical Info / Product Notes Chemiluminescent Acridinium Esters are commonly used to label antibodies, proteins, nucleic acids, and some peptides, frequently for use as a detection reagent in immunoassays.



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](mailto:info-uk@enzolifesciences.com)