

Acridinium C2, NHS Ester

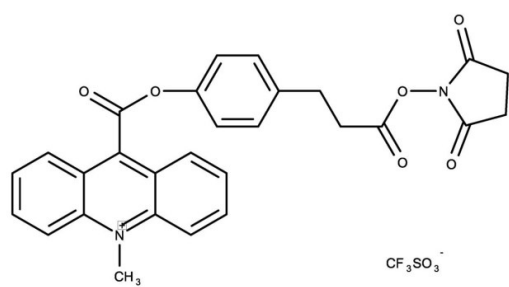
Ordering Information

[Order Online »](#)

ADI-906-000-1	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	For chemiluminescent labeling of antibodies, antigens, and nucleic acids.
CAS	177332-37-5
Formula	$C_{29}H_{23}F_3N_2O_9S$
Formulation	Solid.
MW	632.56
Solubility	Soluble in DMSO or methanol.
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.

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