

NAO (ultra pure)

Mitochondria dye

10-N-Nonyl acridine orange (NAO) is an acridine orange derivative and is generally used as a fluorescent marker of the inner mitochondrial membrane in whole cells. It is believed to bind to negatively charged phospholipids. NAO accumulation in the cell seems to be related to specific interactions with mitochondrial membrane proteins and/or lipids, such as cardiolipin, and is largely independent of mitochondrial membrane potential. With respect to apoptosis, the presence of mitochondrial membrane potential can be probed with Rhodamine123 while the structure and integrity of mitochondria can be assessed using 10-N-nonyl-acridine orange. Wavelength Maxima: Excitation 495nm, Emission 519nm

Citations: 10

[View Online »](#)

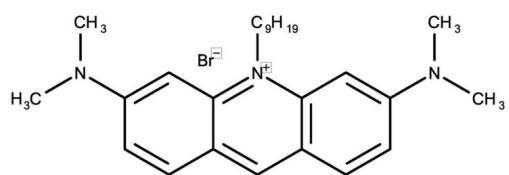
Ordering Information

[Order Online »](#)

ENZ-52306	25mg
-----------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Nonyl Acridine Orange, 3,6-bis(Dimethylamino)-10-nonyl-acridinium bromide
CAS	75168-11-5
MW	472.5
Purity	≥95% (HPLC).
Solubility	Soluble in DMSO.

Technical Info / Product Notes This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.



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)