

DAR-4M AM (cell permeable)

Fluorescent probe for NO detection. Cell permeable.

Cell permeable analog of DAR-4M (Prod. No. ALX-620-067). Can be used to detect nitric oxide (NO) directly in cell specimens prepared from various sources, such as the brain (Ex: ~560nm; Em: 575nm). Stable in a pH range of 4-12. Low decrease in fluorescence intensity over time. Detection limit 7nM. *Not for sale in Japan.*

Citations: 6

[View Online »](#)

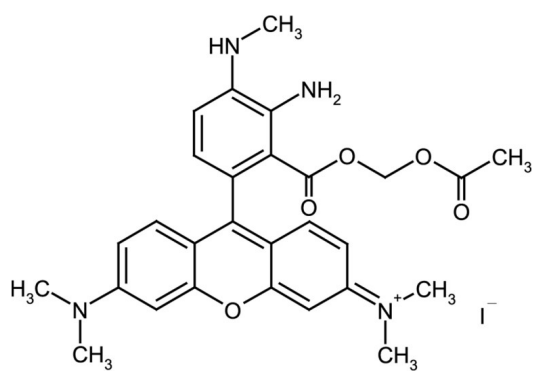
Ordering Information

[Order Online »](#)

ALX-620-069-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Prepare 500-fold dilution (~10µM) in phosphate buffer (0.1M phosphate, pH 7.4) immediately before use. BSA, phenol red and amines may affect the fluorescence and must be used with caution. Do not store diluted solutions.
Handling	Warm to room temperature before opening. Protect from light. Keep under inert gas. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Diaminorhodamine-4M AM
Appearance	Reddish-violet liquid.
Formula	$C_{28}H_{31}N_4O_5 \cdot I$
Formulation	Dissolved in 0.32ml DMSO.
MW	503.6 . 126.9
Purity	≥98% (HPLC)

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)