

DAF-2 DA (cell permeable)

Fluorescent probe for NO detection. Cell permeable.

Cell permeable analog of DAF-2 (Prod. No. ALX-620-052) that is hydrolyzed to DAF-2 by intracellular esterases. Can be used in fluorescence microscopy to measure real-time changes in nitric oxide (NO) levels *in vivo* (Ex(max): 495nm; Em(max): 515nm).

Citations: 16

[View Online »](#)

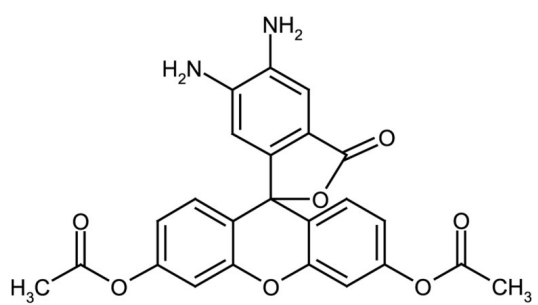
Ordering Information

[Order Online »](#)

ALX-620-056-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	Protect from light. Keep under inert gas. Hygroscopic. 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	4,5-Diaminofluorescein . diacetate
Appearance	Liquid.
CAS	205391-02-2
Formula	$C_{24}H_{18}N_2O_7$
Formulation	Dissolved in 0.45ml DMSO.
MW	446.4
Purity	≥98% (HPLC)
Technical Info / Product Notes	Replacement for ADI-906-042

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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