

DAF-2

Fluorescent probe for NO detection

Highly sensitive fluorescent probe for the real-time detection of nitric oxide (NO) *in vivo*. In the cell, the product reacts rapidly with NO in the presence of O₂ to form the highly fluorescent compound triazolofluorescein (DAF-2T, Ex 495nm; Em 515nm). *Not for sale in Japan.*

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-620-052-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

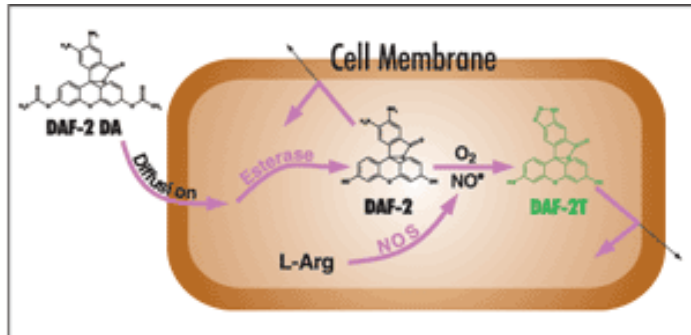
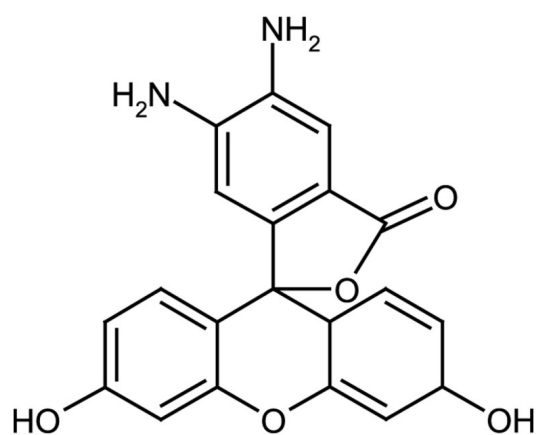


Figure: The cell-permeable diacetate derivative, DAF-2 DA (Prod. No. [ALX-620-056](#)), is used to load cells. Subsequent hydrolysis by cytosolic esterases releases DAF-2 (Prod. No. ALX-620-052)



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. 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
Appearance	Pale yellow liquid.
CAS	205391-01-1
Formula	$C_{20}H_{16}N_2O_5$
Formulation	Dissolved in 0.55ml DMSO.
MW	364.3
Purity	≥98% (HPLC)



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)