

Nitric Oxide Indicators: DAF-FM™ and DAF-FM™ Diacetate

Catalog Numbers D-23841, D-23842, and D-23844

Doc. Part No. MP23841 Pub. No. MAN0002276 Rev. A.0

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Quick facts

Storage upon receipt:

- -20°C
- Desiccate
- Protect from light

Ex/Em: 495/515

Introduction

DAF-FM™ (4-amino-5-methylamino-2',7'-difluorescein; Cat. No. D-23841) and DAF-FM™ diacetate (4-amino-5-methylamino-2',7'-difluorescein diacetate; Cat. Nos. D-23842 and D-23844) represent two important new reagents for quantitating low concentrations of nitric oxide (Figure 1). Developed by Kojima and collaborators^{1,2}, these compounds are essentially nonfluorescent until they react with NO to form a fluorescent benzotriazole. DAF-FM™ diacetate is cell-permeant and passively diffuses across cellular membranes. Once inside cells, it is deacetylated by intracellular esterases to become DAF-FM™. The fluorescence quantum yield of DAF-FM™ is ~0.005, but increases about 160-fold, to ~0.81, after reacting with nitric oxide³. With excitation/emission maxima of 495/515 nm, DAF-FM™ can be detected by any instrument that can detect fluorescein, including flow cytometers, microscopes, fluorescent microplate readers and fluorometers.

Probably the most successful indicator for nitric oxide has been 4,5-diaminofluorescein diacetate (DAF-2 diacetate)^{1,2}, which was also developed by Kojima and collaborators. DAF-2 has been used to identify individual nitric oxide-producing neurons in brain slices^{4,5}, in mitochondria⁶ and in living plant cells⁷. Simultaneous measurements of intracellular Ca²⁺ with fura-2 and nitric oxide production with DAF-2 have been reported⁸.

The DAF-FM™ reagent has some important advantages over DAF-2. The spectra of the NO adduct of DAF-FM™ are independent of pH above pH 5.5³. Also, the NO adduct of DAF-FM™ is significantly more photostable than that of DAF-2³, which means additional time for image capture. Finally, DAF-FM™ is a more sensitive reagent for NO than is DAF-2 (NO detection limit for DAF-FM™ ~3 nM³ versus ~5 nM for DAF-2²).

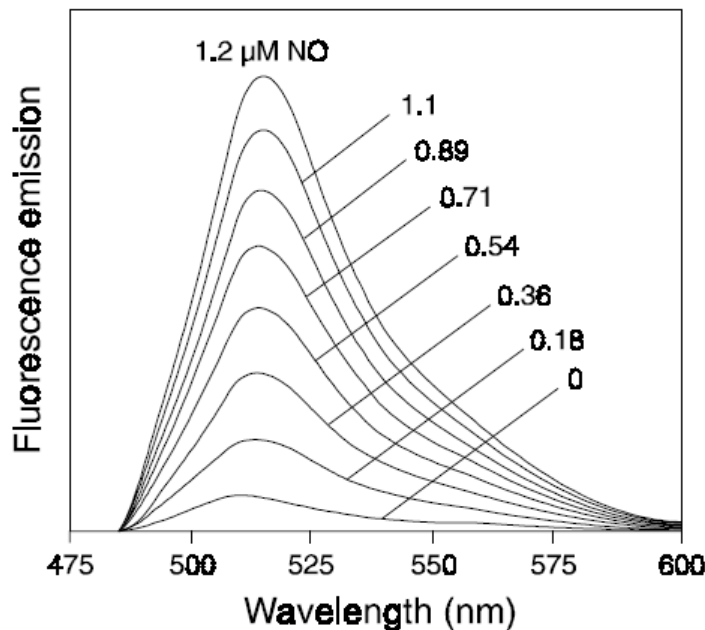


Figure 1 Fluorescence emission spectra of DAF-FM™ in solutions containing zero to 1.2 μM nitric oxide (NO) radical.

Storage and handling

Store DAF-FM™ (Cat. No. D-23841) and DAF-FM™ diacetate (Cat. Nos. D-23841 and D-23844) at -20°C, desiccated and protected from light upon receipt.

A ~7 mM stock solution of DAF-FM™ can be prepared by dissolving the entire contents of the vial in 0.35 mL of high-quality anhydrous DMSO.

A ~5 mM stock solution of DAF-FM™ diacetate can be made by dissolving the 1 mg packaging (Cat. No. D-23841) in 0.4 mL of high-quality anhydrous DMSO or the 50 μg packaging (Cat. No. D-23844) in 20 μL of high-quality anhydrous DMSO.

For long-term storage of the DMSO stock solutions, divide the solution into aliquots in order to minimize freeze-thaw cycles. These aliquoted solutions should be stable for at least six months. Please allow the solutions to warm to room temperature before opening. Working solutions of these reagents should be prepared immediately before use. The diluted reagent should not be stored for later use.

Application

Both the DAF-FM™ and the diacetate DMSO stock solutions can be diluted into aqueous buffers. Bovine serum albumin (BSA) and phenol red may affect the fluorescence and should be used with caution. DAF-FM™ and DAF-FM™ diacetate are sold for research use only. They should not be used for *in vitro* diagnostics.

DAF-FM™

The optimal dilution buffer and working concentration should be determined empirically. A suggested starting concentration range is between 1–10 μ M.

DAF-FM™ can be loaded into cells by pressure injection or perfusion from a patch-clamp pipette. With these methods it is advisable to use a dead-cell stain, such as propidium iodide, to identify cells that do not recover from the loading procedure. Fluorescence excitation and emission maxima are 495 and 515 nm, respectively. Because these wavelengths are very similar to fluorescein, detection systems designed for fluorescein or FITC can be used.

DAF-FM™ diacetate

The following loading protocol is provided as an introductory guide. Optimal loading concentration, time and temperature will need to be determined empirically. In general, it is desirable to use the minimum dye concentration required to yield fluorescence signals with adequate signal to noise ratios. Subcellular compartmentalization is usually lessened by lowering the incubation temperature.

1. Prepare viable cells in suspension or on a slide.
2. Dilute the DMSO stock solution into a suitable buffer. A suggested starting concentration range is between 1–10 μ M.
3. Incubate the cells with the diluted DAF-FM™ diacetate for 20–60 minutes at 4°C to 37°C. Adherent cultures do not need to be trypsinized for loading.
4. Wash the cells to remove excess probe. Replace with fresh buffer or medium, and then incubate for an additional 15–30 minutes to allow complete de-esterification of the intracellular diacetates.
5. Fluorescence excitation and emission maxima are 495 and 515 nm, respectively. Because these wavelengths are very similar to fluorescein, detection systems designed for fluorescein or FITC can be used.

Product list

Unless otherwise indicated, all materials are available through thermofisher.com.

Cat. No.	Product Name	Unit Size
D-23841	DAF-FM™ (4-amino-5-methylamino- 2',7'-difluorofluorescein)	1 mg
D-23842	DAF-FM™ diacetate (4-amino-5-methylamino- 2',7'-difluorofluorescein diacetate)	1 mg
D-23844	DAF-FM™ diacetate (4-amino-5-methylamino- 2',7'-difluorofluorescein diacetate) *special packaging*	10 × 50 μ g

References

1. Chem Pharm Bull 46, 373 (1998);
2. Anal Chem 70, 2446 (1998);
3. Angew Chem Int Ed Engl 21, 3209 (1999);
4. J Neurosci Methods 92, 101 (1999);
5. Brain Res 852, 239 (2000);
6. Biochem Biophys Res Commun 272, 129 (2000);
7. Plant J 23, 817 (2000);
8. Cell Calcium 27, 281 (2000).

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Manufacturer: Life Technologies Corporation | 29851 Willow Creek | Eugene, OR 97402

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

16 August 2018

ThermoFisher
SCIENTIFIC