## 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** 

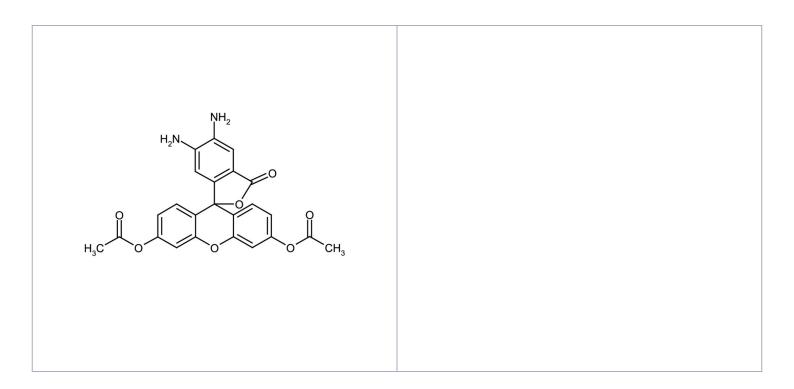
Order Online »

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 ( $\sim$ 10 $\mu$ M) 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

