DAF-FM DA (cell permeable)

Fluorescent probe for NO detection. Cell permeable.

Cell permeable photo-stable nitric oxide (NO) fluorescent indicator, that is hydrolyzed to DAF-FM by intracellular esterases (Ex: 500nm; Em: 515nm). Stable in a pH range of 5.5-12. High yield (quantum efficiency). Detection limit 3nM. *Not for sale in Japan*.

Citations: 9

View Online »

Ordering Information

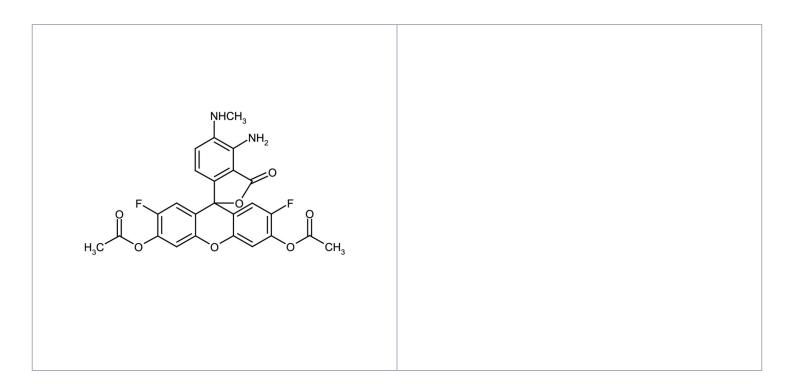
Order Online »

ALX-620-071-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 μ 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. 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 Diaminofluorescein-FM . diacetate

Appearance Liquid.

CAS 254109-22-3

Formula $C_{25}H_{18}F_2N_2O_7$

Formulation Dissolved in 0.4ml DMSO.

MW 496.4

Purity ≥98% (HPLC)

Last modified: May 29, 2024