DAR-4M

Fluorescent probe for NO detection

Fluorescent probe (Ex: ~560nm; Em: 575nm) for the detection of nitric oxide (NO). It can be used to detect NO directly in biopsy tissues, cultured cells, and tissue sections. Stable in a pH range of 4-12. Low decrease in fluorescence intensity over time. Detection limit 7nM. *Not for sale in Japan*.

Citations: 2

View Online »

Ordering Information

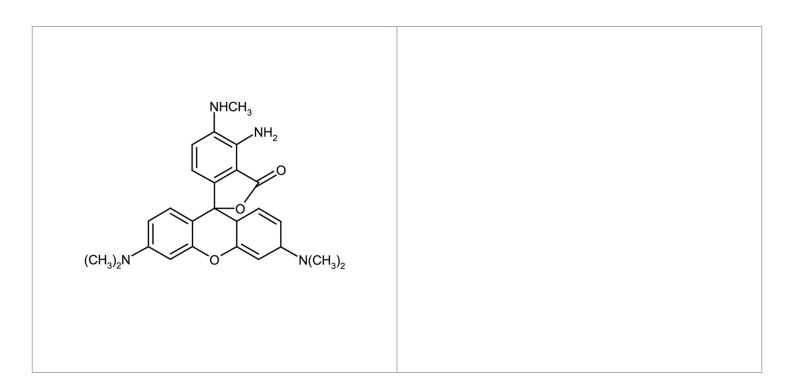
Order Online »

ALX-620-067-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 (~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 Warm to room temperature before opening. 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 Diaminorhodamine-4M

Appearance Reddish-violet liquid.

Couple Target Nitric oxide

Couple Type Ligand

Formula $C_{25}H_{26}N_4O_3$

Formulation Dissolved in 0.47ml DMSO.

MW 430.5

Purity ≥98% (HPLC)

Last modified: May 29, 2024

