

## Fluo-5N, AM \*Cell permeant\*

Catalog number: 20566  
Unit size: 10x50 ug

Component	Storage	Amount (Cat No. 20566)
Fluo-5N, AM *Cell permeant*	Freeze (< -15 °C), Minimize light exposure	10x50 ug

### OVERVIEW

Fluo-5N is an analog of Fluo-4 with lower calcium-binding affinity ( $K_d$  = ~90  $\mu$ M), making it suitable for detecting intracellular calcium levels in the range of 1  $\mu$ M to 1 mM that would saturate the response of Fluo-4. Fluo-5N AM ester may be directly loaded into live cells by adding the dissolved indicator directly to dishes containing the cultured cells. It is compatible with excitation at 488 nm by argon-ion laser sources, making Fluo-5N useful for confocal microscopy, flow cytometry, and microplate screening applications. It has excitation and emission wavelengths at 494 and 516 nm respectively. Upon calcium binding, its fluorescence intensity increases by >100 fold.

### KEY PARAMETERS

#### Flow cytometer

Emission	530/30 nm filter
Excitation	488 nm laser
Instrument specification(s)	FITC channel

#### Fluorescence microscope

Emission	FITC
Excitation	FITC
Recommended plate	Black wall/clear bottom

#### Fluorescence microplate reader

Cutoff	515
Emission	525
Excitation	490
Recommended plate	Black wall/clear bottom
Instrument specification(s)	Bottom read mode/Programmable liquid handling

### PREPARATION OF STOCK SOLUTIONS

*Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles*

#### Fluo-5N AM Stock Solution

1. Prepare a 2 to 5 mM stock solution of Fluo-5N AM in high-quality, anhydrous DMSO.

### PREPARATION OF WORKING SOLUTION

#### Fluo-5N AM Working Solution

1. On the day of the experiment, either dissolve Fluo-5N AM in DMSO or thaw an aliquot of the indicator stock solution to room temperature.
2. Prepare a 2 to 20  $\mu$ M Fluo-5N AM working solution in a buffer of your choice (e.g., Hanks and Hepes buffer) with 0.04% Pluronic® F-127. For most cell lines, Fluo-5N AM at a final concentration of 4-5

$\mu$ M is recommended. The exact concentration of indicators required for cell loading must be determined empirically.

**Note:** The nonionic detergent Pluronic® F-127 is sometimes used to increase the aqueous solubility of Fluo-5N AM. A variety of [Pluronic® F-127 solutions](#) can be purchased from AAT Bioquest.

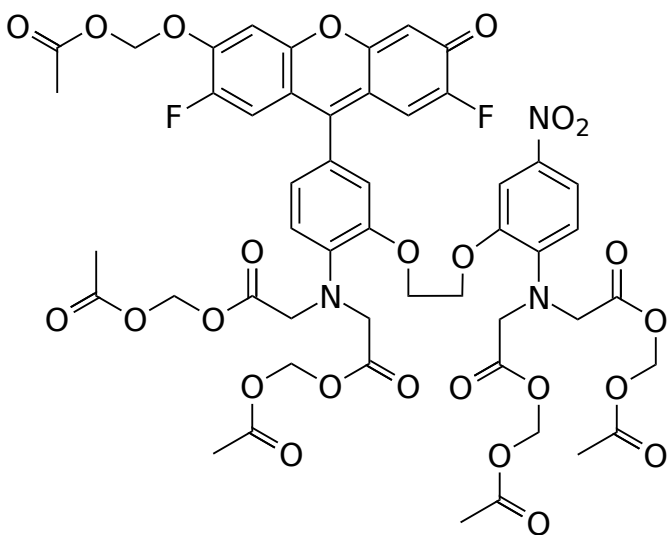
**Note:** If your cells contain organic anion-transporters, probenecid (1-2 mM) may be added to the dye working solution (final in well concentration will be 0.5-1 mM) to reduce leakage of the de-esterified indicators. A variety of [ReadiUse™ Probenecid products](#), including water-soluble, sodium salt, and stabilized solutions, can be purchased from AAT Bioquest.

### SAMPLE EXPERIMENTAL PROTOCOL

Following is our recommended protocol for loading AM esters into live cells. This protocol only provides a guideline and should be modified according to your specific needs.

1. Prepare cells in growth medium overnight.
2. On the next day, add 1X Fluo-5N AM working solution to your cell plate.  
  
**Note:** If your compound(s) interfere with the serum, replace the growth medium with fresh HHBS buffer before dye-loading.
3. Incubate the dye-loaded plate in a cell incubator at 37 °C for 30 to 60 minutes.  
  
**Note:** Incubating the dye for longer than 2 hours can improve signal intensities in certain cell lines.
4. Replace the dye working solution with HHBS or buffer of your choice (containing an anion transporter inhibitor, such as 1 mM probenecid, if applicable) to remove any excess probes.
5. Add the stimulant as desired and simultaneously measure fluorescence using either a fluorescence microscope equipped with a FITC filter set or a fluorescence plate reader containing a programmable liquid handling system such as an FDSS, FLIPR, or FlexStation, at 490/525 nm cutoff 515 nm.

## EXAMPLE DATA ANALYSIS AND FIGURES



**Figure 1.** Chemical structure for Fluo-5N, AM \*Cell permeant\*

#### DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.