

# Fluo-3 AM (ultra pure)

## Calcium dye

Fluo-3 AM is widely used in flow cytometry, confocal laserscanning microscopy and fluorescence spectroscopy. The probe has proved particularly valuable for determining GPCR function via cell-based high-throughput screening. Fluo-3 AM is a membrane permeable probe and can be passively loaded in cells by simple incubation. Once inside the cell, Fluo-3 becomes fluorescent in the presence of free Ca (II) . Fluo-3 is a long wavelength calcium indicator probe (488 nm excitable). The longer wavelength of absorption is convenient for minimizing photodamage in cells. Wavelength Maxima: Excitation 506nm, Emission 526nm

Citations: 4

[View Online »](#)

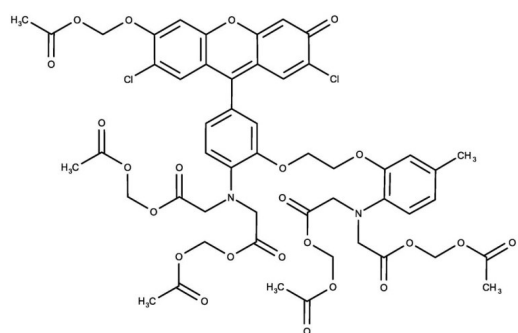
## Ordering Information

[Order Online »](#)

ENZ-52004	1mg
-----------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	1-[Amino-5-(2,7-dichloro-6-acetomethoxy-3-oxo-3H-xanthen-9-yl)phenoxy]-2-(2'-amino-5'-methylphenoxy)ethane-N,N,N',N'-tetraacetic acid, pentaacetoxymethyl ester
CAS	121714-22-5
MW	1129.8
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.
Technical Info / Product Notes	This product is a member of the CELLESTIAL <sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL <sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)