

Fluo-3 (ultra pure)

Calcium dye

Fluo-3 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 is a membrane impermeable probe and can be loaded in cells by microinjection, electroporation or scrape loading. The acetomethoxy (AM) version is necessary for passive cell loading. Once inside the cell, Fluo-3 becomes fluorescent, in the presence of free Ca (II). Fluo-3 is a long wavelength calcium probe (488 nm excitable). The longer wavelength of absorption is convenient for minimizing photodamage in cells. Wavelength Maxima: Excitation 506nm, Emission 526nm

Citations: 2

[View Online »](#)

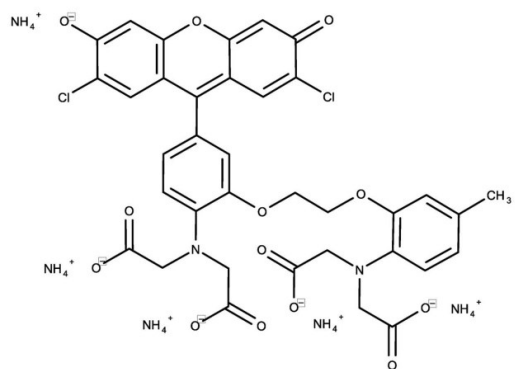
Ordering Information

[Order Online »](#)

ENZ-52005	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-[2-Amino-5-(2,7-dichloro-6-hydroxy-3-oxo-3H-xanthen-9-yl)phenoxy]-2-(2'-amino-5'-methylphenoxy)ethane-N,N,N',N'-tetraacetic acid . 5NH ₃
CAS	339221-91-9
MW	854.7
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

Technical Info / Product Notes

This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] 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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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