

FLUO 3/AM

Fluorescent probe for Ca^{2+} . Cell permeable.

Cell permeable analog of FLUO 3. Fluorescent probe for Ca^{2+} . After crossing the membrane, the product is quickly metabolized by cytoplasmic esterases to the membrane impermeant FLUO 3.

Citations: 7

[View Online »](#)

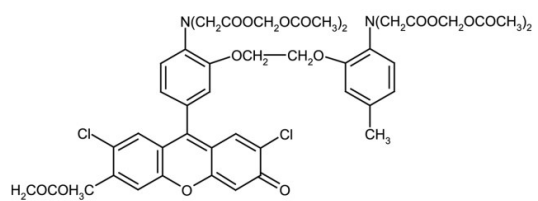
Ordering Information

[Order Online »](#)

ALX-620-003-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

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

Appearance Red solid.

CAS 121714-22-5

Formula $C_{51}H_{50}Cl_2N_2O_{23}$

Identity Determined by 1H -NMR and fluorescence.

MW 1129.9

Purity ≥85% (HPLC)

Solubility Soluble in DMSO (0.9mg/ml) and acetonitrile (2mg/ml).

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-be@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