

# INDO-1 AM (ultra pure)

## Calcium dye

Indo-1 acetomethoxy (AM) is a popular UV-excitable calcium indicator, similar to Fura-2. In contrast to Fura-2, Indo-1 has a dual emission peak. The main emission peak in calcium-free solution is 475nm while in the presence of calcium the emission is shifted to 400nm. Indo-1 AM is a membrane permeable probe and can be easily loaded into cells by passive incubation. The probe is widely used in flow cytometry. For this application, it is critical to have an instrument with an argon laser tuned to UV or a heliumcadmium laser available. Wavelength Maxima: Excitation 346nm, Emission 475nm

Citations: 2

[View Online »](#)

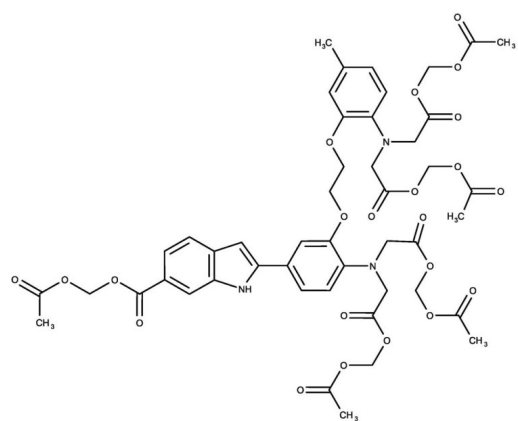
## Ordering Information

[Order Online »](#)

ENZ-52008	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-(6-carboxyindol-2-yl)phenoxy]-2-(2'-amino-5'-methylphenoxy)ethane-N,N,N',N'-tetraacetic acid, pentaacetoxymethylester
Appearance	Yellow powder.
CAS	112926-02-0
Formula	$C_{47}H_{51}N_3O_{22}$
MW	1009.9
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.

Last modified: May 29, 2024



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)