

FURA-2 AM (ultra pure)

Calcium dye

Fura-2 acetomethoxy (AM) is one of the most widely used calcium indicators for homogenous quantitative dual wavelength ratiometric cell measurements. Fura-2 AM is particularly useful for digital imaging microscopy. It is less susceptible to photobleaching than Indo-1. The probe is excited only by UV light, which results in significantly less interference by visible wavelength excitable fluorescent compounds. One application of the probe is as a useful counter screen tool for GPCR and calcium channel hits identified using calcium indicators excited at 488nm in primary screens. Fura-2 AM can be easily loaded into cells by passive incubation. Wavelength Maxima: Excitation 370nm, Emission 476nm

Citations: 8

[View Online »](#)

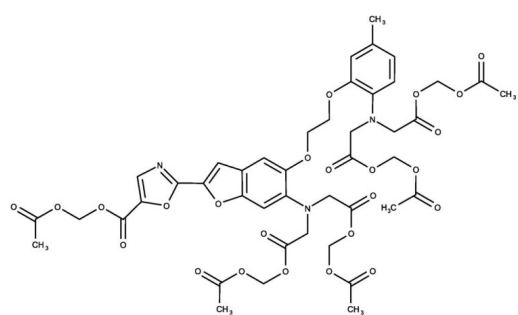
Ordering Information

[Order Online »](#)

ENZ-52006	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-(5-Carboxyoxazol-2-yl)-6-aminobenzofuran-5-oxy]-2-(2'-amino-5'-methyl-phenoxy)ethane-N,N,N',N'-tetraacetic acid, pentaacetoxymethyl ester
Appearance	Off-white to yellow solid.
CAS	108964-32-5
Formula	$C_{44}H_{47}N_3O_{24}$
MW	1001.9
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