

FURA-2 (ultra pure)

Calcium dye

Fura-2 is one of the most widely used calcium indicators for homogenous quantitative dual wavelength ratiometric cell measurements. Fura-2 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 channels identified using calcium indicators excited at 488nm in primary screens. Fura-2 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. Wavelength Maxima: Excitation 363nm, Emission 512nm

Citations: 4

[View Online »](#)

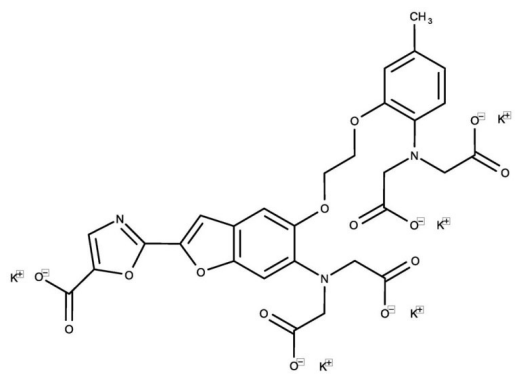
Ordering Information

[Order Online »](#)

ENZ-52007	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 . 5K
Appearance	Light yellow solid.
CAS	113694-64-7
Formula	$C_{29}H_{22}K_5N_3O_{14}$
MW	832.0
Purity	≥95% (HPLC)
Solubility	Soluble in water.

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](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)