

# FLUOFORTE®

## Reagent

### Calcium dye

Used in a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets. Fluorescent calcium indicator dye developed to improve cell loading and calcium response while maintaining the convenient Fluo-3, Fluo-4 and Calcium 4 spectral wavelength of maximum excitation at ~490 nm and maximum emission at ~520 nm. Relative to the cited dyes, FLUOFORTE® AM dye yields the brightest signal and largest assay window.

Citations: 24

[View Online »](#)

### Ordering Information

[Order Online »](#)

ENZ-52015	1mg
-----------	-----

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Red solid.
Application	Flow Cytometry, Fluorescence microscopy, Fluorescent detection, HTS
Application Notes	Provides a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.
Technical Info / Product Notes	The FLUOFORTE <sup>®</sup> Reagent 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-be@enzolifesciences.com](mailto:info-be@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)