

# GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Reagent

## Calcium dye

The GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Reagent is the brightest and most sensitive fluorescent calcium indicator relative to other commercially available red fluorescent dyes with maximum excitation at ~530nm and maximum emission at ~555nm. This product is used in a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets.

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-52016-M001	1mg
ENZ-52016-5C50	5x50µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Protect from moisture. 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	Lyophilized, orange.
Application	Flow Cytometry, Fluorescence microscopy, Fluorescent detection, HTS
MW	~1'000
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

**Technical Info / Product Notes** The GFP-CERTIFIED<sup>®</sup> 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 cell analysis applications involving confocal microscopy, flow cytometry, microplate readers and HCS/HTS, where consistency and reproducibility are required.



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)