

# GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Reagent (Non-AM)

## Calcium dye

The GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Reagent (Non-AM) is the most sensitive fluorescent calcium indicator relative to other commercially available red fluorescent dyes. It can be used for detecting intracellular calcium mobilization across a broad spectrum of biological targets. Esterase hydrolysis is not required for live cell analysis via flow cytometry or fluorescence microscopy.

## Ordering Information

[Order Online »](#)

|                |        |
|----------------|--------|
| ENZ-52555-0250 | 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.      |
| Emission Maximum   | 551nm             |
| Excitation Maximum | 524nm             |
| MW                 | ~800              |
| Purity             | ≥95% (HPLC)       |
| Solubility         | Soluble in water. |

## 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.

Based on application requirement, we also offer ENZ-52016.

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)