

GFP-CERTIFIED[®] FLUOFORTE[®] Reagent

Calcium dye

The GFP-CERTIFIED[®] FLUOFORTE[®] 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	≥90% (HPLC)
Solubility	Soluble in DMSO.

Technical Info / Product Notes

The GFP-CERTIFIED[®] FLUOFORTE[®] Reagent 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 cell analysis applications involving confocal microscopy, flow cytometry, microplate readers and HCS/HTS, where consistency and reproducibility are required.

Last modified: October 14, 2025



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-be@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