

GFP-CERTIFIED[®] FLUOFORTE[®] Reagent (Non-AM)

Calcium dye

The GFP-CERTIFIED[®] FLUOFORTE[®] 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.

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[®] 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.

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

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