

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: 21

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|-----------|--------|
| ENZ-52014 | 5x50µg |
|-----------|--------|

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 | Lyophilized, orange. |
| 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 [®] 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 imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required. |



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