## FLUOFORTE<sup>®</sup> 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<sup>®</sup> AM dye yields the brightest signal and largest assay window.

Citations: 24

View Online »

**Ordering Information** 

**Order Online** »

**ENZ-52015** 1mg

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 Red solid.

**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<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 imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency

and reproducibility are required.

Last modified: May 29, 2024

