# 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® 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 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

