# GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Reagent

#### Calcium dye

The GFP-CERTIFIED® FLUOFORTE® Reagent is the brightest and most sensitive fluorescent calcium indictor 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<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.

Last modified: October 14, 2025

