

# GFP-CERTIFIED<sup>®</sup>

# FLUOFORTE<sup>®</sup>

## Calcium assay kit

### A brighter, more robust and GFP-compatible (orange dye) fluorescent assay for calcium mobilization

Enzo Life Sciences' GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> Calcium Assay Kit (high-throughput) provides a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets. Relative to other commercially available dyes, the GFP-CERTIFIED<sup>®</sup> FLUOFORTE<sup>®</sup> dye is the brightest and most sensitive fluorescent calcium indicator. The kit provides a homogeneous mix-and-read, no-wash calcium mobilization assay. The homogeneous cell-based assay for calcium offers fewer steps, lower variability and an easier protocol for adherent and non-adherent cell lines. In addition, it requires neither a washing step, nor exogenous addition of a quencher dye, which could adversely effect receptor-ligand interaction kinetics.

Citations: 19

[View Online »](#)

### Ordering Information

[Order Online »](#)

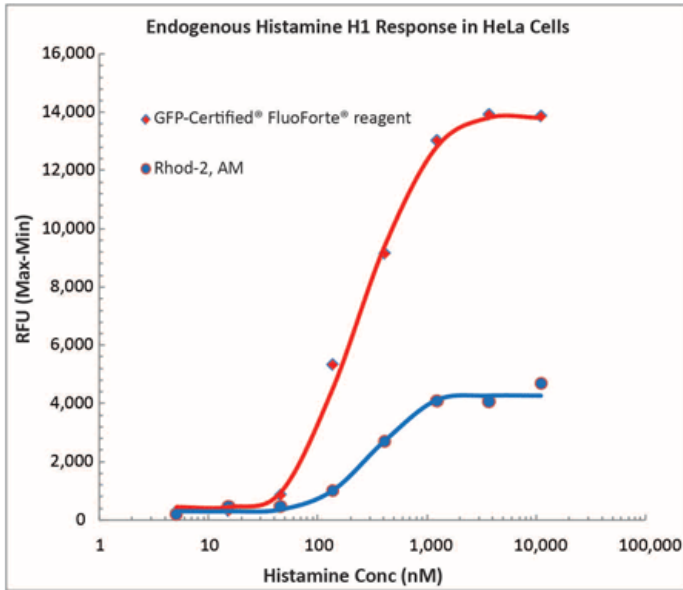
ENZ-51020-KP010

10x96 wells

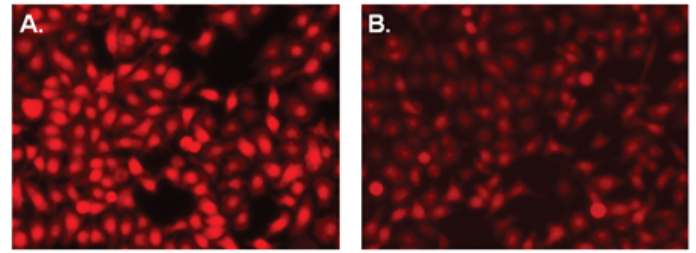
### Manuals, SDS & CofA

[View Online »](#)

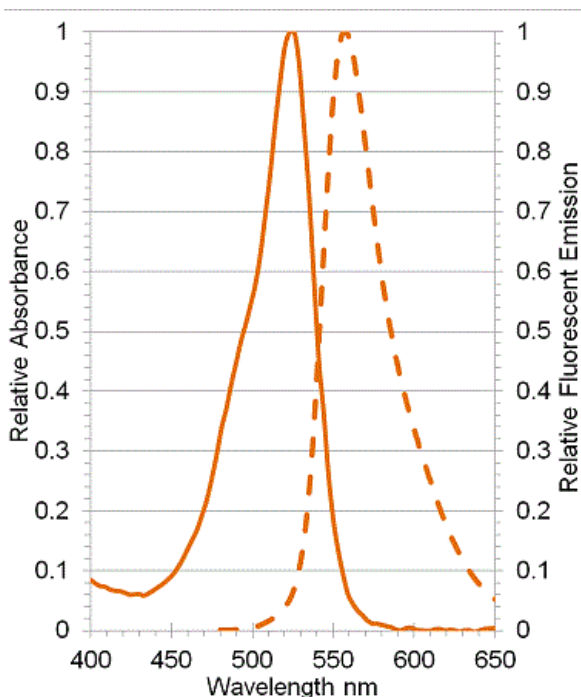
- Monitor calcium mobilization in GFP-expressing cell lines
- The brightest and most sensitive red-shifted fluorescent Ca<sup>2+</sup> indicator
- Provides comparable EC50 and IC50 values as obtained using Rhod-4 AM and Fluo-4 AM
- Larger assay window and higher Z' factor values, allowing measurements of challenging cell lines and receptors



Histamine H1 receptors: HeLa cells were seeded overnight at 40,000 cells per 100  $\mu$ L per well in a 96-well black wall/clear bottom microplate. The cells were incubated with 100  $\mu$ L of GFP-CERTIFIED® FLUOFORTE® reagent or Rhod-2 AM dye for 1 hour at 37°C. Histamine (20 $\mu$ L/well) was added using a BioTek two-syringe pump dispenser to achieve the final indicated concentrations. No significant difference in EC50 of Histamine for GFP-CERTIFIED® FLUOFORTE® and Rhod2-AM dye was observed. GFP-CERTIFIED® FLUOFORTE® reagent generated higher intensity signal and a larger assay window.



U-2 OS cells were incubated with (A) GFP-CERTIFIED® FLUOFORTE® AM dye or (B) Rhod-2 AM dye. Cells were washed and treated with a final concentration of 100nM ATP and imaged immediately. GFP-CERTIFIED® FLUOFORTE® AM dye exhibits brighter fluorescence intensity than Rhod-2 AM.



## 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

Application	Flow Cytometry, Fluorescence microscopy, Fluorescent detection, HTS
Application Notes	The GFP-CERTIFIED® FLUOFORTE® Calcium Assay kit provides a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets.
Contents	<p><b>Starter pack</b> (Prod. No. ENZ-51020-KP010):</p> <p><b>Reagent A</b> (lyophilized GFP-CERTIFIED® FLUOFORTE® dye), 1 vial</p> <p><b>Reagent B</b> (dye efflux inhibitor), 10ml</p> <p><b>Reagent C</b> (Hanks' buffer with 20mM HEPES, 100ml)</p> <p>Sufficient for 10 x 96-well plates</p> <p><b>High-throughput</b> (Prod. No. ENZ-51020-KP100):</p> <p><b>Reagent A</b> (lyophilized GFP-CERTIFIED® FLUOFORTE® dye), 10 vials</p> <p><b>Reagent B</b> (dye efflux inhibitor), 10 x 10ml</p> <p>Sufficient for 100 x 96-well plates</p>
Quality Control	% purity of Reagent A (GFP-CERTIFIED® FLUOFORTE® dye) by HPLC is ≥95%.
Quantity	ENZ-51020-KP010: for 10 x 96-well plates ENZ-51020-KP100: for 100 x 96-well plates

The GFP-CERTIFIED® FLUOFORTE® Calcium assay kit 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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)