



## **FLUOFORTE<sup>®</sup> Calcium Assay Kit**

**Green FLUOFORTE<sup>®</sup> Dye**

Catalog #: ENZ-51016

High-Throughput, 100 plates kit

Catalog #: ENZ-51017

Starter Pack, 10 plates kit

## **USE FOR RESEARCH PURPOSES ONLY**

The 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. Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Enzo, CELLestial, and FLUOFORTE are trademarks of Enzo Life Sciences, Inc. FLIPR is a trademark of Molecular Devices. CellLux is a trademark of PerkinElmer. FDSS is a trademark of Hamamatsu Photonics. Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## TABLE OF CONTENTS

Introduction .....	2
Reagents Provided and Storage .....	2
Other Materials Needed .....	3
Safety Warnings and Precautions .....	3
Methods and Procedures .....	4
A. Cell Preparation .....	4
B. Preparation of FLUOFORTE® Dye-Loading Solution Using <b>ENZ-51017</b> (10 plate kit) .....	5
C. Preparation of FLUOFORTE® Dye-Loading Solution Using <b>ENZ-51016</b> (100 plate kit) .....	5
D. Calcium Mobilization Assay .....	6
E. Calcium Measurement in Flow Cytometry .....	6
F. Calcium Mobilization Microscopy Assay .....	7
Expected Results .....	8
References .....	9
Troubleshooting Guide .....	11
Contact Information .....	18

## INTRODUCTION

The calcium ion is an important second messenger involved in many physiological and signal transduction processes within cells. Fluo-3, Fluo-4 and Calcium 4 dyes are widely used calcium ion indicators for in-cell measurement of agonist-stimulated and antagonist-inhibited calcium signaling in high-throughput screening applications.<sup>1-4</sup> However, their relatively weak fluorescence signals have limited their application in some challenging cell lines and with certain membrane receptors.

Enzo Life Sciences' FLUOFORTE<sup>®</sup> Calcium Assay Kit provides a homogeneous fluorescence-based assay for detecting intracellular calcium mobilization across a broad spectrum of biological targets. Relative to other commercially available dyes, FLUOFORTE<sup>®</sup> dye yields the brightest signal and largest assay window. The kit provides a homogeneous mix-and-read, no-wash calcium mobilization assay. The homogenous 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 affect receptor-ligand interaction kinetics.<sup>5</sup>

## REAGENTS PROVIDED AND STORAGE

All reagents are shipped on dry ice. Upon receipt, the kit should be stored at -20°C, protected from light. When stored properly, these reagents are stable for six months from date of receipt. **Avoid repeated freezing and thawing.**

Reagent	Quantity	
	ENZ-51017 (10 plate kit)	ENZ-51017 (100 plate kit)
Reagent A: FLUOFORTE <sup>®</sup> dye, lyophilized	1 vial	10 vials
Reagent B: Dye efflux inhibitor	10 x 1 mL	10 x 10 mL
Reagent C: Hanks' buffer with 20 mM HEPES (HHBS)	1 x 100 mL	Not included

## OTHER MATERIALS NEEDED

- A fluorometric imaging plate reader capable of performing quantitative optical screening for cell-based kinetic assays. (Molecular Devices FLIPR, PerkinElmer CellLux, Hamamatsu FDSS system, or similar instrumentation)
- Calibrated, adjustable precision pipettors, preferably with disposable plastic tips
- Deionized water
- Anhydrous DMSO
- Serum (optional).
- Growth medium (e.g., Dulbecco's modified Eagle medium, D-MEM)
- 10X Hanks' Balanced Salt Solution, HBSS (e.g., Invitrogen #14065-056)
- 1M HEPES Buffer (e.g., Invitrogen #15630-080)
- Assay Plates: 96- or 384-well black-wall, clear bottom plates or 1536-well low base black-wall, clear bottom plates, 1536-well lids
- Compound plates: 96-well or 384-well polypropylene plates, 1536-well polystyrene plates
- Polypropylene tubes compatible with a flow cytometer

## SAFETY WARNINGS AND PRECAUTIONS

- This product is for research use only and is not intended for diagnostic purposes.
- Some components of this kit may contain hazardous substances. They can be harmful if ingested or absorbed through the skin and may cause irritation to the eyes. The reagents of the kit should be treated as possible mutagens and should be handled with care and disposed of properly.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas. All blood components and biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.
- To avoid photobleaching, perform all manipulations in low light environments or protected from light by other means.

## METHODS AND PROCEDURES

### Brief Summary of Assay Work Flow

- *Prepare cells.*
- *Remove medium.*
- *Add FLUOFORTE<sup>®</sup> dye-loading solution.*
- *Incubate plate for 1 hour.*
- *Add test agents.*
- *Read fluorescence.*

**NOTE:** PLEASE READ THE ENTIRE PROCEDURE BEFORE STARTING. Allow all reagents to be used to warm to room temperature before proceeding. Upon thawing of solutions, gently hand-mix or vortex the reagents prior to use to ensure a homogenous solution.

### A. CELL PREPARATION

1. **Adherent Cells.** The day before the experiment, plate the cells overnight in growth medium using  $4 \times 10^4$  to  $8 \times 10^4$  cells per well at a plating volume of 100  $\mu$ L per well for 96-well plates, or using  $1 \times 10^4$  to  $2 \times 10^4$  cells per well at a plating volume of 25  $\mu$ L per well for a 384-well plates.

After overnight incubation, remove the growth medium from the cell plates. Then proceed to section D, page 6.

**NOTE:** It is important to remove the growth medium in order to minimize background fluorescence, and compound interference with serum or culture media.

2. **Non-adherent Cells.** On the day of the experiment, centrifuge the cells from the culture medium and then resuspend the cell pellets in FLUOFORTE<sup>®</sup> dye-loading solution (see section B). Plate the cells using  $1.25 \times 10^5$  to  $2.5 \times 10^5$  cells per well at a plating volume of 100  $\mu$ L per well for 96-well plates, or  $3 \times 10^4$  to  $6 \times 10^4$  cells per well at a plating volume of 25  $\mu$ L per well for 384-well plates. Centrifuge the plates at 800 rpm for 2 minutes, **with brake off**, prior to starting the experiments. Proceed to section D.3 page 6.

**NOTE:** Each cell line should be evaluated on an individual basis to determine the optimal cell density for the intracellular calcium mobilization assay.

## B. PREPARATION OF FLUOFORTE® DYE-LOADING SOLUTION USING ENZ-51017 (10 plates kit)

The following procedure is for preparation of FLUOFORTE® dye-loading solution **for use in 1 plate**. Before starting, equilibrate Reagent A (FLUOFORTE® dye), a vial of Reagent B and Reagent C (HBSS) to room temperature.

1. **Reagent A Stock Solution.** Add 100 µL DMSO to the vial containing **Reagent A**. Mix well.

***NOTE:** 10 µL of the reconstituted Reagent A is enough for 1 plate. The remaining unused, reconstituted Reagent A can be aliquoted and stored at < -20°C for at least one month if stored properly. The tubes (preferably amber vials) should be capped tightly. Avoid exposure to light and repeated freeze-thaw cycles.*

2. **1X Assay Buffer.** Mix well 9 mL of **Reagent C** (HHBS) with the contents of 1 vial (1 mL) of **Reagent B**.

***NOTE:** 10 mL of 1X Assay Buffer is sufficient for one plate. Unused buffer may be stored at < -20°C up to 1 month. Avoid exposure to light and repeated freeze-thaw cycles.*

3. **FLUOFORTE® Dye-Loading Solution.** Add 10 µL of Reagent A Stock Solution (from step B-1) to 10 mL of 1X Assay Buffer (from step B-2). Mix well. This working solution is stable for at least **2 hours** at room temperature.

## C. PREPARATION OF FLUOFORTE® DYE-LOADING SOLUTION USING ENZ-51016 (100 plates kit)

The following procedure is for preparation of FLUOFORTE® dye-loading solution **for use in 10 plates**. Before starting, equilibrate Reagent A (FLUOFORTE® dye) and a vial of Reagent B to room temperature.

1. **Hanks' Buffer with 20 mM HEPES (HHBS).** Prepare 100 mL of HHBS by mixing the following:

10 mL 10X Hanks' Balanced Salt Solution (HBSS)

2 mL HEPES Buffer

88 mL Deionized water

2. **1X Assay Buffer.** Mix well 90 mL of **HHBS** (from step C-1) with the contents of 1 vial (10 mL) of **Reagent B**.

3. **FLUOFORTE® Dye-Loading Solution.**

- a. Dissolve the contents of 1 vial of **Reagent A** in 100 µL of DMSO. Mix well.



- b. Add 100  $\mu$ L of **Reagent A** solution to 100 mL of 1X Assay Buffer (from step C-2). Mix well. This working solution is stable for at least **2 hours** at room temperature.

#### D. CALCIUM MOBILIZATION PLATE ASSAY

1. Obtain prepared cell plates (see section A).
2. Add FLUOFORTE® Dye-Loading Solution to each well (100  $\mu$ L/well for 96-well plates or 25  $\mu$ L/well for 384-well plates).
3. Incubate the cell plates for 1 hour at room temperature, or incubate about 45 minutes at 37°C then incubate for 15 minutes at room temperature prior to assay.

**NOTE:** *The incubation time should be optimized for each cell line. The incubation time should be limited to 1~2 hours. DO NOT wash the cells after dye loading.*

4. Prepare the compound plates by dissolving the compound in the buffer of choice. The FLUOFORTE® Calcium Assay is optimized for an agonist addition between 1/5-1/6 of the total final volume (e.g., Add 20-25  $\mu$ L/well of agent when using a 96-well plate that contains 100  $\mu$ L of FLUOFORTE® Dye-Loading Solution<sup>2</sup>).
5. Run the calcium flux assay by monitoring the fluorescence at Ex=490 nm/Em=525 nm with a fluorometric imaging plate reader.

**NOTE:** *Faster addition speeds can lead to better mixing of compounds and lower signal variance across the plate. Make sure to follow the recommended experimental setup parameters provided by the instrument manufacturer before reading the plate. It is also important to run the signal test before the experiment. Different instruments have their own intensity range. Adjust the signal test intensity to the level of 10% to 15% of the maximum instrument intensity counts. For example, the maximum fluorescence intensity count for FLIPR-384 is 65,000, so the instrument settings should be adjusted to have its signal test intensity around 7,000 to 10,000.*

#### E. CALCIUM MEASUREMENT IN FLOW CYTOMETRY

1. Aliquot between  $1.25 \times 10^5$  to  $2.5 \times 10^5$  cells per well at a volume of 250  $\mu$ L in each polypropylene tube.
2. Add 250  $\mu$ L FLUOFORTE® Dye-Loading Solution to each well.
3. Incubate the cell plates for 1 hour at room temperature, or incubate about 45 minutes at 37°C then incubate for 15 minutes at minutes at room temperature prior to assay.



**NOTE:** The incubation time should be optimized for each cell line. The incubation time should be limited to 1~2 hours. **DO NOT** wash the cells after dye loading.

4. Add agonist at desired concentration in small volume. Use a vehicle control.
5. Incubate the cell plates for 1 hour at room temperature, or incubate about 45 minutes at 37°C then incubate for 15 minutes at room temperature prior to assay.

**NOTE:** The incubation time should be optimized for each cell line. The incubation time should be limited to 1~2 hours. **DO NOT** wash the cells after dye loading.

6. Run samples on flow cytometer

**NOTE:** Calcium flux is a transient event. The timing of the sample analysis should be consistent and may need optimization.

## F. CALCIUM MOBILIZATION MICROSCOPY ASSAY

1. Obtain prepared cell slides (see section A).
2. Add FLUOFORTE® Dye-Loading Solution to each region on the slide.
3. Incubate the cell slides for 1 hour at room temperature, or incubate about 45 minutes at 37°C then incubate for 15 minutes at room temperature prior to assay.

**NOTE:** The incubation time should be optimized for each cell line. The incubation time should be limited to 1~2 hours. **DO NOT** wash the cells after dye loading.

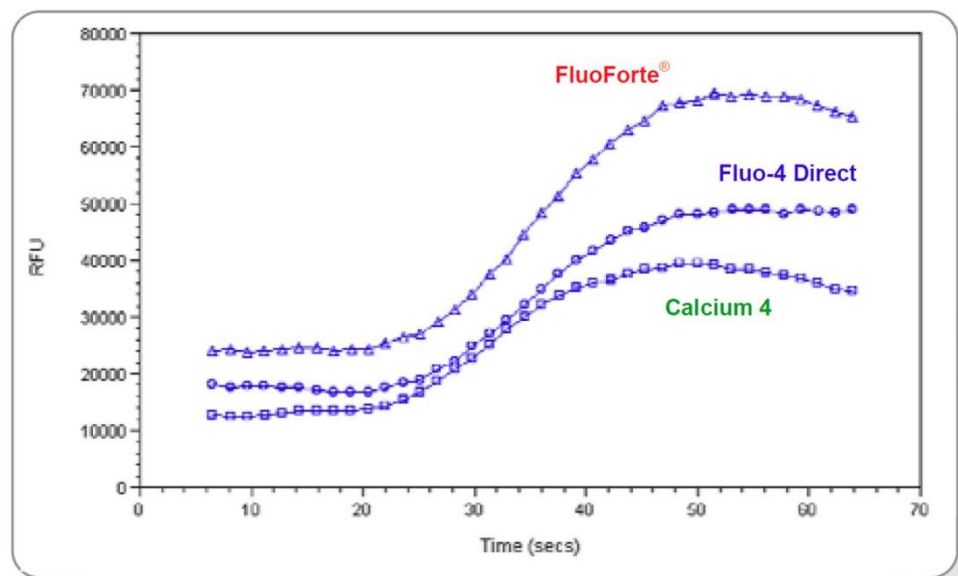
4. Prepare the compound plates by dissolving the compound in the buffer of choice. The FLUOFORTE® Calcium Assay is optimized for an agonist addition between 1/5-1/6 of the total final volume (e.g., Add 20-25 µL/well of agent when using a 96-well plate that contains 100 µL of FLUOFORTE® Dye-Loading Solution<sup>2</sup>).
5. Run the calcium flux assay by monitoring the fluorescence with a fluorescence microscope with the appropriate filter.

## EXPECTED RESULTS

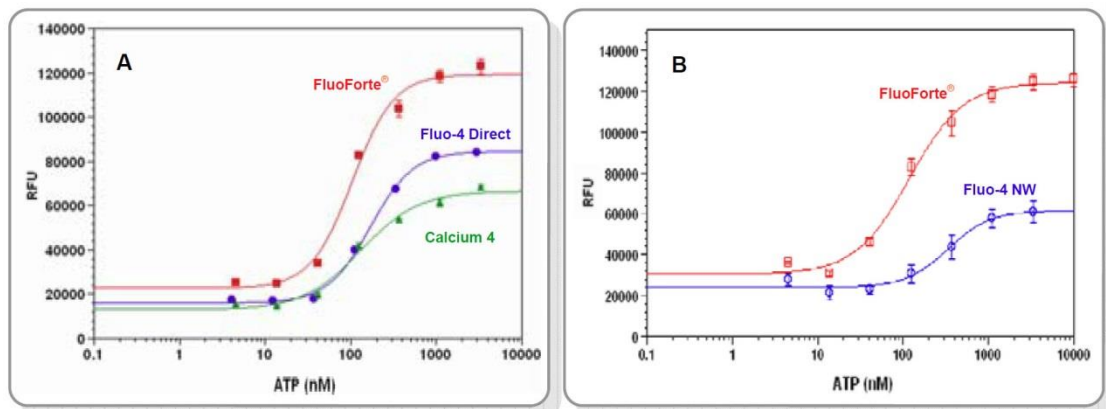
In a side-by-side comparison of FLUOFORTE<sup>®</sup>, Fluo-4 Direct assay and Calcium 4 assay, CHO M1 cells were stimulated with 200 nM of ATP. FLUOFORTE<sup>®</sup> yields the brightest signal and largest assay window. (shown in Figure 1). This facilitates measurements of challenging cell lines and receptors.

Dose responses for ATP in CHO M1 cells gave similar EC<sub>50</sub> for all the assays (shown in Figure 2). This demonstrates consistent pharmacology among the assays. However, relative fluorescence units (RFU) of FLUOFORTE<sup>®</sup> are much higher than Fluo 4 and Calcium 4.

Overall FLUOFORTE<sup>®</sup> Calcium Assay kit provides an optimized assay method for monitoring G-protein-coupled receptors (GPCRs) and calcium channels.<sup>6</sup> Its ability to generate very strong signal enables researchers to perform calcium mobilization assays with a wide range of receptor and calcium channel targets.



**Figure 1: Comparisons of FLUOFORTE<sup>®</sup> Fluo-4 Direct, and Calcium 4 detection of intracellular calcium mobilization in CHO-M1 cells.** CHO cells were seeded overnight in 40,000 cells per 100  $\mu$ L per well in a 96-well black wall, clear bottom costar plate. The cells were incubated with 100  $\mu$ L of Life Technologies' Fluo-4 Direct kit, Molecular Devices' Calcium 4 kit (both based on manufactures' protocol) or Enzo's FLUOFORTE<sup>®</sup> kit. ATP (20  $\mu$ L/well) was added by FlexStation to achieve concentrations of 200 nM.



**Figure 2: ATP Dose Response Curves in CHO-M1 cells.** CHO cells were seeded overnight at 40,000 cells per 100  $\mu$ L per well in a 96-well black wall/clear bottom microplate. **Panel A:** The cells were incubated with 100  $\mu$ L of Life Technologies' Fluo-4 Direct kit, Molecular Devices' Calcium 4 kit (both based upon manufacturer's protocol) or Enzo's FLUOFORTE<sup>®</sup> dye. **Panel B:** The cells were incubated with 100  $\mu$ L of Enzo's FLUOFORTE<sup>®</sup> Calcium assay kit, or with Life Technologies' Fluo-4 NW kit (based upon manufacturer's protocol).

ATP (20  $\mu$ L/well) was added by FlexStation to achieve the final indicated concentrations. No significant difference in EC<sub>50</sub> of ATP for FLUOFORTE<sup>®</sup>, Fluo-4 Direct and Calcium 4 was observed (shown in panel A), and also comparable EC<sub>50</sub> values were observed between FLUOFORTE<sup>®</sup> and Fluo-4 NW (shown in panel B). In all cases, FLUOFORTE<sup>®</sup> generated the highest intensity signal.

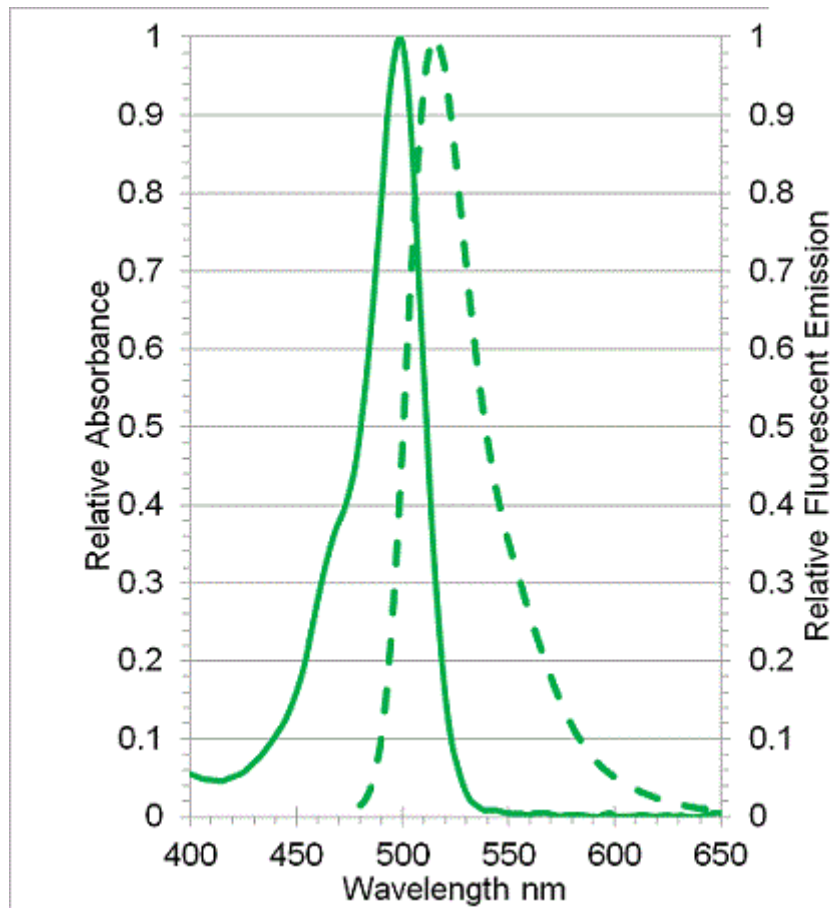
## REFERENCES

1. Martin VV, Beierlein M, Morgan JL, Rothe A, Gee KR. (2004) Novel Fluo-4 analogs for fluorescent calcium measurements. *Cell Calcium*, 36, 509.
2. do Ceu Monteiro M, Sansonetty F, Goncalves MJ, O'Connor JE. (1999) Flow cytometric kinetic assay of calcium mobilization in whole blood platelets using Fluo-3 and CD41. *Cytometry*, 35, 302.
3. Perez-Terzic C, Stehno-Bittel L, Clapham DE. (1997) Nucleoplasmic and cytoplasmic differences in the fluorescence properties of the calcium indicator Fluo-3. *Cell Calcium*, 21, 275.
4. Tretyn A, Kado RT, Kendrick RE. (1997) Loading and localization of Fluo-3 and Fluo-3/AM calcium indicators in sinapis alba root tissue. *Folia Histochem Cytobiol*, 35, 41.

5. Zheng, W. et al. (2004) High throughput assay technologies for ion channel drug discovery. *Assay Drug. Dev. Technol.* 5, 543–552.
6. Kiselyov K, Shin DM, Muallem S. (2003) Signalling specificity in GPCRdependent Ca<sup>2+</sup> signalling. *Cell Signal.* 15(3), 243-253.

## TROUBLESHOOTING GUIDE

Problem	Potential Cause	Suggestion
High baseline fluorescence	Contributions to baseline fluorescence by growth medium and organic anion transport	Remove the medium before adding the indicator dye to the wells. Use the dye solution within 2 hours at room temperature.
Untreated cells have calcium response	Inconsistent DMSO concentration	Make sure that the buffer used for the negative control wells have the same final concentration of DMSO as those in the test compounds.
Response is smaller than expected	Agonists and antagonists may stick to the pipette tips. Experimental setup parameters and dye loading time are not optimized.	Use 0.1% of BSA in all compound buffer diluents. Fast addition speeds are recommended to ensure better mixing of compounds and improved cell response. Dye loading typically takes between 30 minutes and one hour. Optimizing the conditions for each cell line is recommended.
Well to well variation observed	Incorrect dispenser and experimental setup parameter used.	Use instrument manufacturer's recommended dispenser and setup parameters (i.e., volume, height and speed of dispensing) for compound addition.
Fluorescence drop upon compound addition	Dislodging the cells during addition	Decrease the rate of addition or seed fewer cells in the wells to avoid this problem.



Excitation and emission spectra for the FLUOFORTE<sup>®</sup> dye.



# Product Manual

## NOTES





# Product Manual

## NOTES



# Product Manual

## NOTES



# Product Manual

## NOTES



# Product Manual

## NOTES



# Product Manual

**[www.enzolifesciences.com](http://www.enzolifesciences.com)**

***Enabling Discovery in Life Science®***

**GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**EUROPE/ASIA**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
[info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

For local distributors and detailed product information visit us online:

**[www.enzolifesciences.com](http://www.enzolifesciences.com)**