

LYSO-ID[®] Red Cytotoxicity Kit (GFP CERTIFIED[®])

Catalog #: ENZ-51015-KP002

2 x 96-well plates

USE FOR RESEARCH PURPOSES ONLY

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

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**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
Precautions	3
Reagents Provided and Storage	3
Additional Materials Required	3
Reagent Preparation	4
Staining Live, Adherent Cells	5
Appendices	6
References.....	8
Trouble Shooting.....	9
Contact Information.....	10

INTRODUCTION

Cationic amphiphilic drugs, toxic agents and other basic compounds often show signs of large apparent accumulation in organisms, consistent with various forms of cellular sequestration. A variety of cell types are known to respond to these agents by the formation of multiple large vacuoles, which have been variously referred to as lamellar bodies or lysosomal inclusion bodies. Drug concentration into acidic vacuoles is thought to be due to ion trapping and the vacuole enlargement that follows is considered to arise from osmotic phenomena.¹⁻⁴ The vacuoles generated by cationic amphiphilic drugs, such as procainamide and chloroquine, have been formally classified as autophagosomes.²⁻⁴ The persistence of giant autophagosome structures in cells is thought to be due to the retention of the stable cationic drugs by continuing ion trapping, as well as inhibition of autodigestion pathways. Repeated cycles of autophagy may also add layers of membrane to the vacuoles.⁴ The cytopathological sequestration of drugs within cells, leading to an excess accumulation of phospholipid-rich membranes in vacuoles, is referred to as phospholipidosis.

Enzo Life Sciences' LYSO-ID® Red Cytotoxicity Kit (GFP CERTIFIED®) is a 96-well cell-based assay that provides a rapid and quantitative approach for determining drug- or toxic agent-induced lysosome and lysosome-like organelle perturbations in live cells. Unlike conventional lysosome stains, this kit is effective for detecting phospholipidosis, which is induced by cationic amphiphilic drugs, such as chloroquine and verapamil. In addition to compounds that cause phospholipidosis, agents that cause the accumulation of autophagosomes by blocking the downstream lysosomal pathway and/or intracellular trafficking of autophagosomes also lead to increases in the accumulation of intracellular LYSO-ID® Red dye signal in the described assay. This microplate assay offers several advantages over alternative methods based upon electron microscopy, fluorescence microscopy, flow cytometry or long term incubation with fluorescent phospholipid analogs. Chief among these advantages are the ability to analyze drug response in cells without co-incubation with artificial phospholipid analogs and the ability to perform drug screening in a rapid and quantitative high-throughput manner using a conventional fluorescence microplate reader. A lysosome-perturbation agent, verapamil, is provided as a positive control for monitoring changes in vacuole number and volume. A blue nuclear counterstain is integrated into the detection reagent to identify cell death or loss.

SAFETY WARNINGS & PRECAUTIONS

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



Handle with care



Avoid freeze / thaw cycles

- This product is for research use only and is not intended for diagnostic purposes.
- Some components of this kit may contain hazardous substances. Reagents can be harmful if ingested or absorbed through the skin and may cause irritation to the eyes. They 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.

REAGENTS PROVIDED AND STORAGE

All reagents are shipped on dry ice. Upon receipt, the kit should be stored upright at $\leq -20^{\circ}\text{C}$, protected from light. When stored properly, these reagents are stable for at least twelve months. Avoid repeated freezing and thawing.

Reagents provided in the kit are sufficient for two 96-well plates using live, adherent cells.



Reagents require separate storage conditions.

Reagent	Quantity
10X Dual Color Detection Reagent	2 x 1ml
Detection Buffer	20ml
Verapamil Control	3 μ mol
10X Assay Buffer	15ml

ADDITIONAL MATERIALS REQUIRED

1. Fluorescence microplate reader
2. Calibrated, adjustable precision pipetters, preferably with disposable plastic tips
3. Glass microscope slides (optional)
4. Glass cover slips (optional)
5. Deionized water
6. Anhydrous DMSO (optional)
7. Serum (e.g., Fetal Bovine Serum)
8. Growth medium (e.g., Dulbecco's Modified Eagle Medium, D-MEM)

METHODS AND PROCEDURES

NOTE: Allow all reagents to thaw at room temperature before starting with the procedures. Upon thawing, gently hand-mix or vortex the reagents prior to use to ensure a homogenous solution. Briefly centrifuge the vials at the time of first use, as well as for all subsequent uses, to gather the contents at the bottom of the tube.

REAGENT PREPARATION

Positive Control

Verapamil is a cationic amphiphilic drug known to cause phospholipidosis, the accumulation of phospholipids in acidic organelles in the cells.⁵

Verapamil is supplied as a positive control for increasing lysosome number and volume. Reconstitute the lyophilized verapamil (3 μ moles) in 30 μ l DMSO for a 100mM stock solution. It is recommended that overnight treatment with the agent be performed using 50-120 μ M final concentration in medium in order to observe changes in lysosomal signal intensity. Unused stock solution of verapamil may be stored at -20°C for several weeks.

1X Assay Buffer

Allow the 10X Assay Buffer to warm to room temperature. Make sure that the reagent is free of any crystallization before dilution. The addition of fetal bovine serum (FBS) to 2% is highly recommended.

Prepare 1X Assay Buffer by mixing the following:

- 15ml 10X Assay Buffer
- 3ml Fetal Bovine Serum (FBS)
- 132ml Deionized Water

1X Dual Color Detection Reagent

The following procedure is for preparation of 1X Dual Color Detection Reagent for use in one 96-well plate. The addition of fetal bovine serum (FBS) to 2% is highly recommended.

In an appropriate size container, mix the following:

- 1ml 10X Dual Color Detection Reagent
- 8.8ml Detection Buffer
- 0.2ml Deionized Water or Fetal Bovine Serum (FBS)

STAINING LIVE, ADHERENT CELLS

The procedure described below was developed for U2OS, epithelial cell line (MCDK) and HeLa cells for which it is recommended that cells be seeded on plates at a density of 2.0×10^5 to 2.5×10^5 cells/mL, using 100 μ l cells/well. Any cell number and plate coating requirements should be optimized for the chosen cell model.

Positive control cells should be pretreated with verapamil (see step page 4) overnight. Response to verapamil is time and concentration dependent and may also vary significantly among cell types and cell lines. Negative control cells should be treated with a vehicle (water, DMSO, media or other solvent used to reconstitute or dilute an inducer or inhibitor) for an equal length of time under similar conditions.

1. Seed cells in 96-well microplates, using 100 μ l cells/well, the day before addition of test compound/probe. The cells should be plated such that at the end of the experiment the well should be about 90% confluent. Incubate overnight to allow the cells to grow and adhere to the plate under standard tissue culture practices.
2. After overnight incubation, treat the cells with the compound of interest under normal culture conditions for a time period sufficient for assessing the effects of the agent.

For positive control, treat the cells overnight with 50-100 μ M final concentration of verapamil in medium. The negative control (untreated cells) should be treated with the same vehicle used (e.g., DMSO) as the test cells.

3. After incubation with the compound of interest, carefully aspirate the medium and dispense 100 μ l of 1X Assay Buffer (from page 4) to each well.
4. Carefully aspirate all the buffer and dispense 100 μ l of the 1X Dual Color Detection Reagent (from page 4) to each well.
5. Protect the samples from light and incubate for 30 minutes at room temperature.
6. Wash the cells twice with 200 μ l of 1X Assay Buffer each wash. Remove all excess buffer and add 80 μ l 1X Assay Buffer to each well.
7. Analyze the plate with a fluorescence microplate reader. It is recommended to acquire data as soon after completing the assay as feasible. The red lysosome stain can be read with a Texas Red filter (Excitation ~540nm, Emission ~680) and the blue nuclear counterstain can be read with a DAPI filter

set (Excitation ~340, Emission ~480). If the blue nuclear counterstain signal decreases by more than 30%, the compound is considered generally cytotoxic. Increases in the red lysosome signal, without significant loss of blue signal, indicates the accumulation of the probe within the cells arising from an increase in lysosome or lysosome-like vesicle size and/or number.

APPENDICES

Filter Set Selection

The selection of optimal filter sets for a fluorescence microplate application requires matching the optical filter specifications to the spectral characteristics of the dyes employed in the analysis. Consult the microplate reader or filter set manufacturer for assistance in selecting optimal filter sets for your microplate reader.

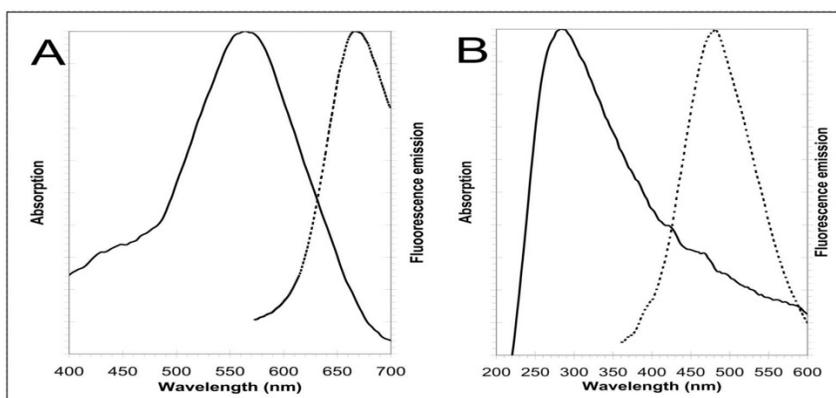


Figure 1. Absorption and fluorescence emission spectra for the red lysosome stain (A) and the blue nuclear counterstain (B). All spectra were determined in 1X Assay Buffer.

Validation of Detection Reagent and Pertubing agent

Lysosomes are membrane-bound subcellular organelles involved in the degradation of macromolecules and pathogens in diverse processes including endocytosis, phagocytosis and autophagy. The red fluorescent probe is selectively sequestered in acidic organelles by a mechanism involving protonation and retention within the organelles. Through careful selection of titratable groups on the probe, labeling has been expanded into the lysosome-like vacuoles of cells pretreated with weakly basic, cell-permeant compounds, such as the anti-arrythmic drug verapamil. The U2OS cells, pretreated overnight with 100 μ M verapamil show a dramatic increase in lysosome-like vesicle number and volume, confirming that the probe associates with this subcellular compartment. The probe can be employed for highlighting lysosome-like organelles under certain conditions, wherein cells produce lysosome-like bodies that contain most of the degradative enzymes of the lysosome, but are not as acidic as the parent organelle. Using a

conventional fluorescence microplate reader, the half maximal effective concentration (EC50) of verapamil of 36 μ M was estimated (see **Figure 2**). The high Z-factor scores obtained, using this assay, demonstrate excellent signal-to-noise and signal-to-background ratios.

This organelle profiling toolkit can provide a screening tool to help researchers in selecting a successful candidate compound with low or weak lysosome-perturbing activity for further drug development efforts, as well as providing preliminary benchmarking of dosing limits in preclinical toxicity studies. It may also be effective in screening compounds in cell-based models of various lysosomal storage diseases.

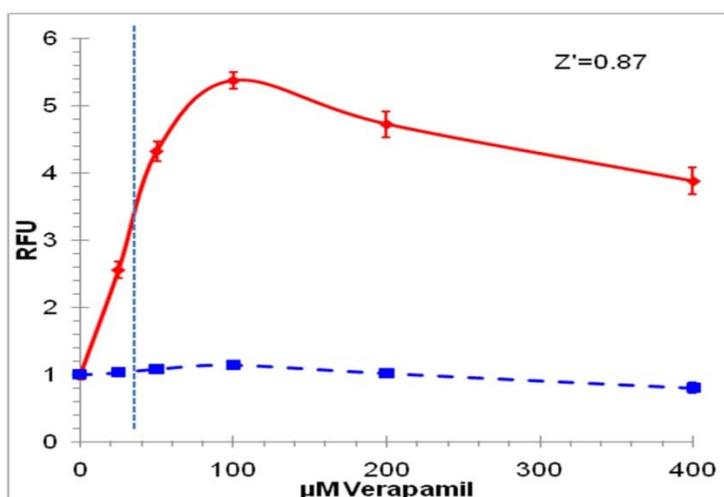


Figure 2: Using a conventional fluorescence microplate reader, the half maximal effective concentration (EC50) of verapamil in U2OS cells was estimated. The high Z-factor (0.87 for 100 μ M verapamil) obtained using the assay demonstrates excellent signal-to-noise and signal-to-background ratios. The error bars denote the standard deviation of at least six determinations. The red line with diamond markers denotes the red fluorescent lysosomal signal, and the blue line with square markers denotes the blue nuclear signal.

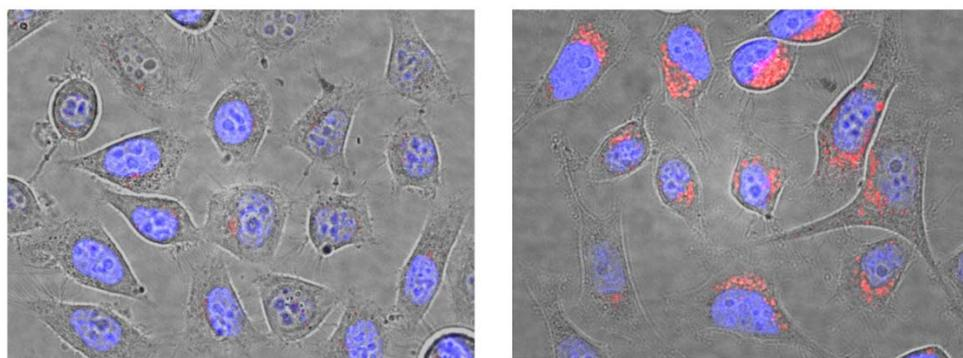


Figure 3: Composite bright-field and fluorescence microscopy images demonstrating staining of U2OS cells with Dual Color Detection Reagent. Cells pretreated overnight with a perturbing agent are shown in the right panel

REFERENCES

1. Anderson N. and Borlak J: Drug-induced phospholipidosis FEBS Lett. 2006. 580: 5533–5540.
2. Morissette, G., Moreau, E., C.-Gaudreault, R., Marceau, F., 2004. Massive cell vacuolization induced by organic amines such as procainamide. J. Pharmacol. Exp. Ther. 310, 395–406
3. Morissette, G., Moreau, E., C.-Gaudreault, R., Marceau, F., 2005. N-substituted 4-aminobenzamides (procainamide analogs): an assessment of multiple cellular effects concerning ion trapping. Mol. Pharmacol. 68, 1576–1589.
4. Morissette, G., Lodge, R. and Marceau, F. 2008 Toxicol. and Appl. Pharmacol. 228, 364–377.
5. Lemieux, B., Percival, M.D., Falguyret, J.-P., 2004. Quantitation of the lysosomotropic character of cationic amphiphilic drugs using the fluorescent basic amine Red DND-99. Anal. Biochem. 327, 247-251.

TROUBLESHOOTING GUIDE

Problem	Potential Cause	Suggestion
Acidic organelles are not sufficiently observed.	A low concentration of the Dual Color Detection Reagent was used or the reagent was incubated with the cells for an insufficient length of time.	Either increase the reagent concentration or increase the time allowed for the dye to accumulate in the lysosome.
Acidic organelle signals are too low to be readily quantified.	Insufficient cell concentration	Increase the number of cells per well. The ideal concentration is ~90% confluency.
Insignificant difference between untreated cells and induced cells observed	Insufficient removal of excess stain	Carefully remove all liquid during the wash steps by aspiration. A second aspiration is recommended to remove all remaining liquid. If the problem persists, additional wash steps may be added.
Dual Color Detection Reagent fails to stain acidic organelles in fixed and/or permeabilized cells.	The dye is only suitable for live-cell staining.	Use the dye only for live-cell analysis.
Precipitate is observed in the 10X Assay Buffer	Precipitate forms at low temperatures.	Allow solution to warm to room temperature or 37°C, then vortex to dissolve all precipitate.
Blue nuclear counterstain signal is dramatically reduced in treated cells.	The compound is generally cytotoxic.	If the blue nuclear signal decreases by more than 30%, the compound is generally cytotoxic. Optimize the concentration of the perturbing reagent.
Cells do not appear healthy by microscopic examination.	Some cells require serum to remain healthy.	Add serum to the detection reagent and wash solutions. Serum improves staining. Typical amounts of serum to add range from 2% to 10%.
Verapamil-treated cells appear dead or are no longer attached to the plate surface.	The EC ₅₀ of Verapamil may differ with different cell lines	Try lowering the dose of verapamil, or shortening the time of exposure.



Product Manual

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

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

For local distributors and detailed product information visit us online:
www.enzolifesciences.com