

EarlyTox Cell Integrity Kit

The EarlyTox™ Cell Integrity Kit from Molecular Devices is an optimized set of reagents that simplifies the measurement of live and dead cells in a single well. The assay uses two nuclear dyes that allow you to detect changes in outer cell membrane permeability resulting from cell damage or cell death via necrosis, apoptosis, or other mechanisms. Cell viability is assessed by counting the number or percent of live and damaged cells. Use the assay for studies of the effects of different treatments on cell viability, evaluation of toxicity effects of pharmaceutical compounds or other chemicals, studies of necrosis and apoptosis, and many other applications. The EarlyTox Cell Integrity Kit is designed to work with many cell types, both adherent and non-adherent.

The simple workflow and reagent performance of the protocol make it amenable to high-throughput screening using the ImageXpress® Micro XLS Widefield High Content Analysis System, ImageXpress® Micro 4 Automated Imaging System, ImageXpress® Nano Automated Imaging System, or the SpectraMax® i3x Multi-Mode Microplate Reader with the SpectraMax® MiniMax™ 300 Imaging Cytometer. See [Setup for Imaging Instruments on page 11](#) and [Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer on page 9](#).

Available Kits

Assay Kit	Explorer Kit	Bulk Kit
EarlyTox™ Cell Integrity Kit (200 Tests, two (2) 96-well or 384-well plates)	R8213	
EarlyTox™ Cell Integrity Kit (1000 Tests, ten (10) 96-well or 384-well plates)		R8214

Contents

Chapter 1: About the EarlyTox Cell Integrity Kit	3
Assay Principle	3
Chapter 2: Materials and Equipment	5
Kit Components	5
Materials Required But Not Provided	5
Storage and Handling	5
Chapter 3: Assay Protocols	7
Standard Protocol	7
Homogeneous Assay Protocol	8
Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer	9
Setup for Imaging Instruments	11
Chapter 4: Data Analysis Examples	13
Obtaining Support	15

EarlyTox Cell Integrity Kit

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/patents>

Product manufactured by Molecular Devices, LLC.
3860 N. First Street, San Jose, California, 95134, United States of America.
Molecular Devices, LLC is ISO 9001 registered.
©2023 Molecular Devices, LLC.
All rights reserved.



Chapter 1: About the EarlyTox Cell Integrity Kit

1

The EarlyTox Cell Integrity Kit provides a fast, simple, and reliable fluorescence-based assay to detect changes in cell viability based on permeability of outer cell membranes (cell integrity).

The kit is based on two nuclear dyes:

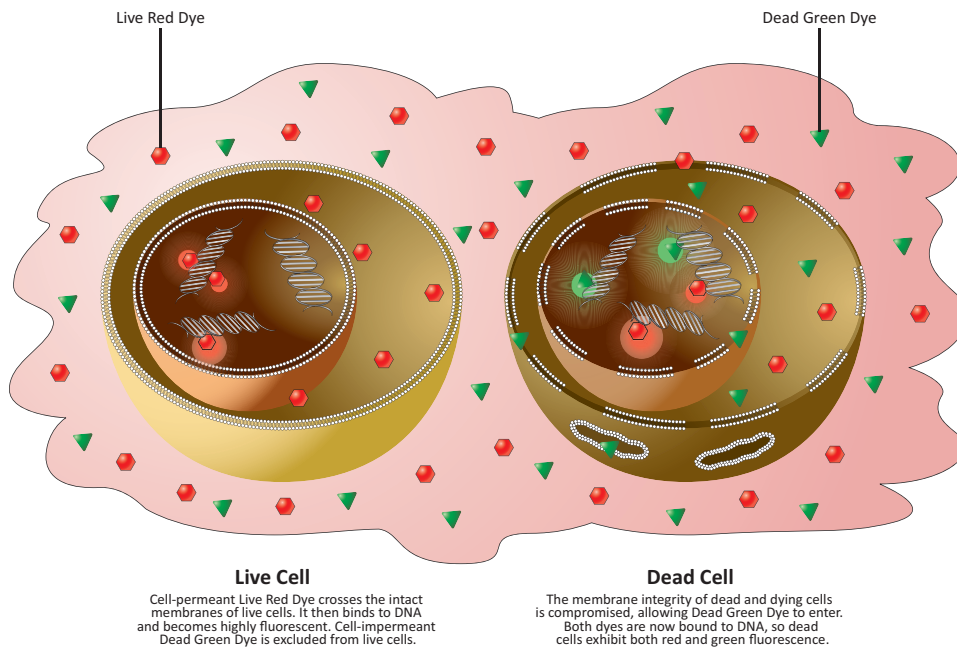
- Live Red Dye is cell permeant and marks both live and dead cells (Excitation: 622 nm/Emission: 645 nm).
- Dead Green Dye is cell impermeant and stains only cells with damaged outer membranes (Excitation: 503 nm/Emission: 526 nm).



Note: Use the appropriate wavelength setting for your instrument. See [Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer on page 9](#) or [Setup for Imaging Instruments on page 11](#).

The kit also contains a proprietary EarlyTox Masking Reagent to increase assay performance in a homogeneous format. Each kit enables a simple mix-and-read procedure whereby cells are incubated with reagents for 30 minutes and then transferred directly to an imaging instrument for evaluation. This eliminates time-consuming wash steps.

Assay Principle



EarlyTox Cell Integrity Kit Assay Principle



Kit Components

Components of the EarlyTox Cell Integrity Kits

Item	Explorer Kit (R8213)	Bulk Kit (R8214)
EarlyTox Live Red Dye	1 vial	1 vial
EarlyTox Dead Green Dye	1 vial	1 vial
EarlyTox Masking Reagent	1 bottle	3 bottles

- The entire EarlyTox Cell Integrity Explorer Kit (R8213) is sufficient for two (2) 96-well or 384-well plates.
- The entire EarlyTox Cell Integrity Bulk Kit (R8214) is sufficient for ten (10) 96-well or 384-well plates.

Materials Required But Not Provided

Reagents and Supplies

Item	Suggested Vendor
DPBS Buffer (1x Dulbecco's Phosphate Buffered Saline) pH 7.4	Corning Cellgro or equivalent

Storage and Handling

On receipt of the EarlyTox Cell Integrity Kit, store the contents at -20°C . Under these conditions the reagents are stable for six months in the original packaging.



WARNING! Reagents can contain chemicals that are harmful. Exercise care when handling reagents as described in the related safety data sheet (SDS or MSDS). The safety data sheet is available in the Knowledge Base on the Molecular Devices support web site: support.moleculardevices.com/



Standard Protocol

To run the standard protocol:

1. Plate the cells at the applicable plating densities (generally overnight) and treat the cells with compounds according to the specific experiment (generally for 24 to 48 hours).



Note: For imaging applications, use black-well, clear-bottom plates.

2. Prepare 1:2000 dilutions of the dyes (add 10 μ L of each reagent to 20 mL of DPBS) immediately before cell staining.
 3. Gently remove the cell media from the plates by pipetting or aspirating (do not wash the cells).
 4. Add the staining solution directly into the wells.
 - Add 100 μ L per well for 96-well plates.
 - Add 25 μ L per well for 384-well plates.
 5. Return the plates to the incubator and incubate for 15 to 30 minutes at 37°C and 5% CO₂.
 6. Proceed with image acquisition using the appropriate wavelengths for the Live Red Dye and Dead Green Dye. See [Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer on page 9](#) or [Setup for Imaging Instruments on page 11](#).
 - Live Red Dye can be detected using the 713 channel (Ex: 625 nm/Em: 713 nm).
 - Dead Green Dye can be detected by using the 541 channel (Ex: 456 nm/Em: 541 nm).
- Typical exposure times are 10 to 20 ms.



Note: Adjust the exposure times for the 713 and 541 channels to achieve the correct color balance between channels. Generally, 541 exposure time is 2x greater than 713 exposure time.

Homogeneous Assay Protocol



Note: In this protocol the culture medium is not removed from the wells. Use the optional Masking Reagent to reduce the background fluorescence typically created by the medium.

To run the Homogeneous Assay protocol:

1. Grow cells in 96-well or 384-well plates with the appropriate treatments. The final volume of media should be approximately 100 μ L for 96-well plates and 25 μ L for 384-well plates.
2. Prepare 2x (1:1000) solution of staining reagents in DPBS (for example, 10 μ L of each reagent in 10 mL of DPBS) immediately before cell staining.
3. Add a volume of 2x staining solution (equal to the volume already in the well) directly to the wells.
 - Add 100 μ L for 96-well plates.
 - Add 25 μ L for 384-well plates.
4. Return the plates to the incubator and incubate for 30 minutes at 37°C and 5% CO₂.
5. Optional:
 - Prepare a 5x solution of Masking Reagent (add 10 mL of DPBS to the bottle and vortex for ~1 minute).
 - Add the 5x solution of Masking Reagent directly into the media (25 μ L per 96-well plate, 6.3 μ L per 384-well plate).
 - Image the plate immediately.



Note: Do not add Masking Reagent simultaneously with staining solution.

6. Proceed with image acquisition using the appropriate wavelengths for the Live Red Dye and Dead Green Dye. See [Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer on page 9](#) or [Setup for Imaging Instruments on page 11](#).

Typical exposure time 10 to 20 ms. If Masking Reagent is used, it is necessary to increase imaging exposure times (100 to 200 ms).



Note: For optimal results it is important to use the recommended reagent volumes and concentrations. However, standard procedures vary across laboratories, and Molecular Devices recognizes that a variety of cell-handling conditions might be adopted at the discretion of the user. Assay development might be required for optimization.

Suggested Plating Volumes and Seeding Densities

Cell Type	96-well plate (100 μ L growth medium with compounds)	384-well plate (25 μ L growth medium with compounds)
Adherent cells	4000 to 20000 (cells per well)	1000 to 5000 (cells per well)
Non-adherent cells	20000 to 100000 (cells per well)	5000 to 20000 (cells per well)

Setup for SpectraMax i3x Reader With the Spectra Max MiniMax 300 Imaging Cytometer

Acquire images from the prepared plate using the SpectraMax i3x Multi-Mode Microplate Reader with the SpectraMax MiniMax 300 Imaging Cytometer and the SoftMax® Pro Data Acquisition and Analysis Software from Molecular Devices with the following suggested settings.

Wavelength:

- Excitation: 456 nm / Emission: 541 nm for Dead Green Dye
- Excitation: 625 nm / Emission: 713 nm for Live Red Dye

Image Acquisition Settings:

- Select control and treated wells for the **Max Image** and **Min Image** wells
- For Live Red Dye (713 nm), use a typical exposure time of 10 to 20 ms
- For Dead Green Dye (541 nm), use a typical exposure time of 20 ms

Image Analysis Settings:

- Use **Discrete Object Analysis**
- **Wavelength for Finding Objects:** 713 nm
- **Find Objects**
 - **Finding Method:** Draw on Images, or **Set Size and Intensity**



Note: If you are using Masking Reagent, select **Set Size and Intensity**.

- **Draw on Images** settings: **Nuclei** preset drawing analysis
Set Size and Intensity settings:
Min. Width: 5 µm
Max. Width: 30 µm
Intensity Above Background: 100–200 (adjust as needed for best results)
 - Click **Apply**
- **Classification**
 - **Settings:** Create New Classification
 - **Type A:** Rename to **Live**
Select the objects that are red but not green.
 - **Type B:** Rename to **Dead**
Select the objects that are both green and red.
 - Click **Apply**
- **Measurements**
 - Default (all parameters selected)

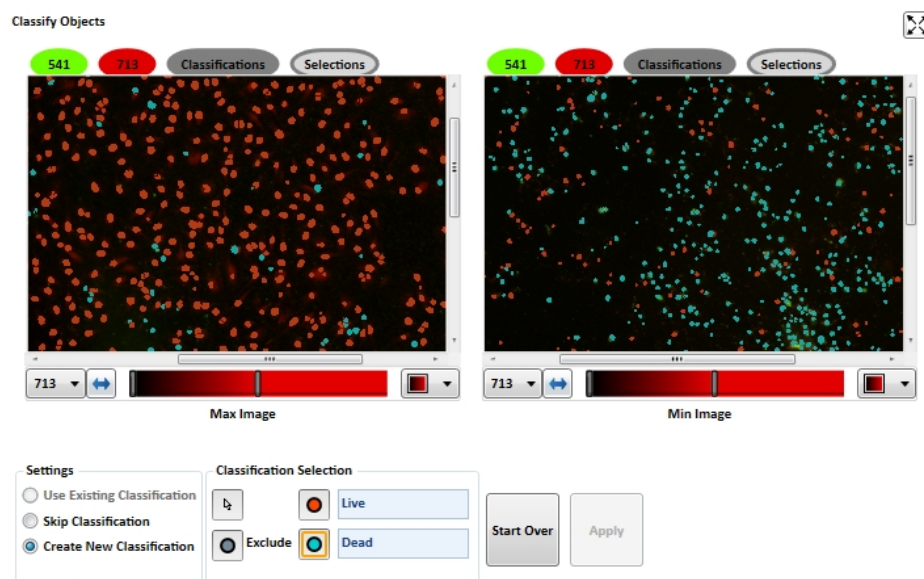
Click **OK** to save the acquisition and analysis settings.

Click **Read**  to acquire the images with the defined analysis settings.

In the **Plate** section, define a template for the tested compounds and concentrations.

Create appropriate tables and graphs using the desired output parameters.

Use SoftMax Pro Software or another program for additional analysis.



Example of Results of Classification for EarlyTox Cell Integrity Assay Data Analysis Using the SoftMax Pro Software

Setup for Imaging Instruments

Acquire images from the prepared plate using the ImageXpress Micro XLS Widefield High Content Analysis System, ImageXpress Micro 4 Automated Imaging System, and ImageXpress Nano Automated Imaging System. Use the following suggested configurations and settings in the MetaXpress High-Content Image Acquisition and Analysis Software from Molecular Devices.

- 20x, 10x, or 4x objectives
- Cy5 filter cube (Ex: 628 nm, Em: 692 nm) for Live Red Dye with a typical exposure time of 10 ms
- FITC filter cube (Ex: 485 nm, Em: 530 nm) for Dead Green Dye staining with a typical exposure time of 20 ms
- For 384-well plates with the 10x or 20x objectives, take one (1) or more images per well
- For 96-well plates with the 10x or 20x objectives, take two (2) to four (4) images per well
- For the 4x objective, take one (1) image per well

For image analysis use the MetaXpress Software with the Live-Dead image processing module. The images are first segmented into individual cells using the Live Red nuclear stain and designated Live or Dead based on the presence of the Dead Green stain. Typical settings for image analysis include the following:

Wavelength 1 Parameters (for Live Red Dye):

- **Wavelength 1 Source:** Cy5
- **Approximate min width:** 7 μm
- **Approximate max width:** 30 μm
- **Intensity above local background:** ~500 graylevels

Wavelength 2: Parameters (for Dead Green Dye):

- **Wavelength 1 Source:** FITC
- **Approximate min width:** 5 μm
- **Approximate max width:** 30 μm
- **Intensity above local background:** ~500 graylevels



Note: Nuclear sizes and brightness of staining can vary depending on specific cells and between experiments. The suggested settings should be adjusted for the specific experiment.

Wavelength 1 Parameters

Wavelength 1 Source:

Display result image:

Adaptive Background Correction™ system

Algorithm:

Stained cell type:

Stained area:

Approximate min width: $\mu\text{m} = 11$ pixels

Approximate max width: $\mu\text{m} = 46$ pixels

Intensity above local background: graylevels

Split touching objects

Wavelength 2 Parameters

Wavelength 2 Source:

Display result image:

Stained cell type:

Stained area:

Approximate min width: $\mu\text{m} = 8$ pixels

Approximate max width: $\mu\text{m} = 46$ pixels

Intensity above local background: graylevels

Split touching objects

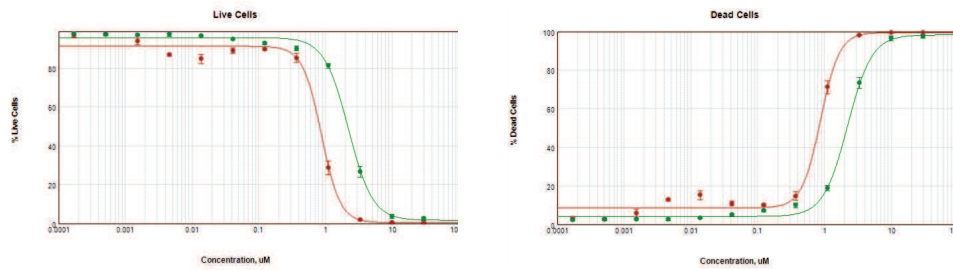
Suggested Settings for EarlyTox Cell Integrity Assay Data Analysis Using the MetaXpress Software and the Live-Dead Application Module

In the **Configure Summary Log** dialog, select the following parameters for toxicity evaluation:

- Total Cells (number of all cells, nuclear count)
- Live Cells
- Dead Cells
- % Live Cells
- % Dead Cells

Chapter 4: Data Analysis Examples

4



Examples of 4-parametric curve fits for concentration dependence of % of intact (left) or damaged (right) cells. Concentration responses are presented for staurosporine (red), IC_{50} $0.2 \mu\text{M}$; and mitomycin C (green), IC_{50} $3.4 \mu\text{M}$.



Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

Our Support website—support.moleculardevices.com/—describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to the Molecular Devices Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance, you can submit a request to Molecular Devices Technical Support.

Please have the instrument serial number (on the rear of the instrument), or reagent kit lot number, and any related sample data files available when you call.



Contact Us

Phone: [+1-800-635-5577](tel:+18006355577)
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.