

# ER-ID<sup>®</sup> Red assay kit (GFP-CERTIFIED<sup>®</sup>)

Widely cited Endoplasmic Reticulum staining  
with minimal toxicity to cells

Enzo Life Sciences ER-ID<sup>®</sup> Red assay kit (GFP-CERTIFIED<sup>®</sup>) contains a red-emitting, cell-permeable small molecule organic probe to stain Endoplasmic Reticulum (ER) in live cell, or fixed cells. The dye emits in the Texas Red region of the visible light spectrum and can be readily used in combination with other common UV and visible light excitable organic fluorescent dyes/fluorescent proteins in multi-color imaging and detection applications. The kit includes also the Hoechst 33342 dye for the nuclear staining.

Citations: 31

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-51026-K500	1Kit
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

- Specifically designed for use with GFP-expressing cell lines, as well as cells expressing blue, cyan or yellow fluorescent proteins (BFPs, CFPs, YFPs)
- Suitable for use with live or post-fixed cells
- Highly resistant to photo-bleaching, concentration quenching and photoconversion
- Suitable for use in conjunction with fluorescein- or coumarin-labeled antibodies

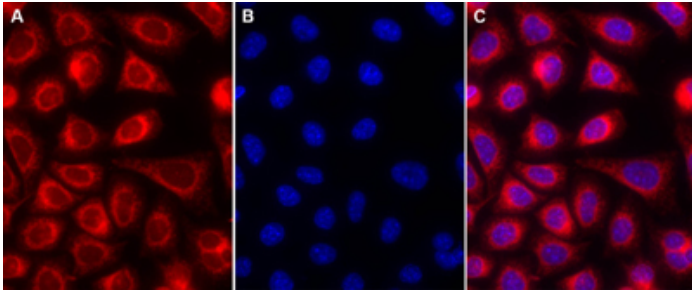


Figure 1: Live HeLa cells stained with ER-ID® Red dye (A), Hoechst dye (B) and resulting composite image (C).

# Handling & Storage

Use/Stability	With proper storage, the kit components are stable up to the date noted on the product label. Store kit at -20°C in a non-frost free freezer, or -80°C for longer term storage.
Handling	Protect from light. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Endoplasmic reticulum / Organelle
Application	Fluorescence microscopy
Application Notes	Enzo Life Sciences' ER-ID <sup>®</sup> Red assay kit (GFP-CERTIFIED <sup>®</sup> ) contains a novel endoplasmic reticulum-selective dye suitable for live cell, or detergent-permeabilized aldehyde-fixed cell staining.
Contents	<b>ER-ID<sup>®</sup> Red detection reagent:</b> 50µl <b>Hoechst 33342 nuclear stain:</b> 50µl <b>10X assay buffer:</b> 15ml
Quality Control	A sample from each lot of ER-ID <sup>®</sup> Red assay kit is used to stain HeLa cells, using the procedures described in the user manual. The selectivity of the ER-ID <sup>®</sup> Red dye is evident.
Quantity	500 assays

The ER-ID<sup>®</sup> Red assay kit (GFP-CERTIFIED<sup>®</sup>) 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: 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)