

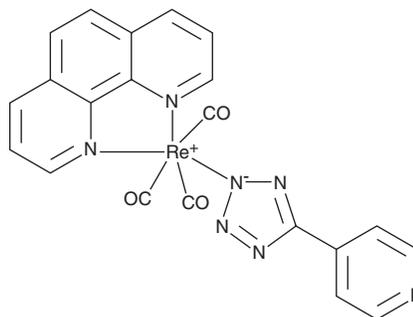
PRODUCT INFORMATION



ReZolve-ER™

Item No. 25908

CAS Registry No.: 1404104-40-0
Formal Name: (OC-6-33)-tricarbonyl
(1,10-phenanthroline-κN¹,κN¹⁰)
[4-(2H-tetrazol-5-yl-κN²)
pyridinato]-rhenium
MF: C₂₁H₁₂N₇O₃Re
FW: 596.6
Ex./Em. Max: 405/570 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

ReZolve-ER™ is a fluorogenic probe that can be used to label the endoplasmic reticulum in live and fixed cells. It easily crosses live cell membranes and accumulates in the endoplasmic reticulum, and its fluorescence can be used to assess endoplasmic reticulum structures using epifluorescent, confocal, and two-photon microscopy. ReZolve-ER™ can be washed from live cells and has low cytotoxicity making it suitable for intermittent monitoring. ReZolve-ER™ displays excitation/emission maxima of 405/570 nm, respectively, and can be used for live and fixed cell applications.

Assay Protocol

The amount of ReZolve-ER™ provided is sufficient to label 150-300 slides or 1,440-2,880 individual wells of cells when utilized in a 96-well plate format, depending on the protocol and application used.

1. Prepare ReZolve-ER™ solution

- Reconstitute ReZolve-ER™ with 300 µl of DMSO and mix thoroughly to prepare a 10 mM ReZolve-ER™ stock solution.
- Store ReZolve-ER™ stock solution at room temperature protected from light.

Note 1: ReZolve-ER™ should not be reconstituted in aqueous solutions such as PBS or cell culture media.

2. Prepare and stain the cells

a. Live adherent cells:

- Dilute ReZolve-ER™ stock solution in cell culture media to a final concentration of 50-100 µM to prepare a ReZolve-ER™ working solution.
- Remove cell culture medium from adherent cells and replace with ReZolve-ER™ working solution.
- Incubate cells at 37°C, 5% CO₂ for 15 minutes. Do not wash.
- Observe cells using fluorescence technique of choice.
- After imaging, aspirate ReZolve-ER™ working solution and wash cells with PBS. Cells may require several wash steps.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/31/2018

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



b. Live suspended cells:

- i. Dilute ReZolve-ER™ stock solution in cell culture media to a final concentration of 50-100 μM to prepare a ReZolve-ER™ working solution.
- ii. Adhere cells to a poly-L-lysine-coated coverslip and allow the cells to settle for 2-5 minutes.
- iii. Aspirate cell culture media and add ReZolve-ER™ working solution.
- iv. Incubate for 10 minutes.
- v. Wet mount coverslip.
- vi. Observe cells using fluorescence technique of choice.

c. Fixed cells:

- i. Fix cells in 4% paraformaldehyde for 20 minutes at room temperature.
- ii. Wash samples 3 x 10 minutes in PBS.
- iii. Dilute ReZolve-ER™ stock solution in PBS to a final concentration of 50-100 μM to prepare a ReZolve-ER™ working solution.
- iv. Incubate cells for 15 minutes at room temperature prior to imaging. Do not wash.
- v. Mount coverslips using ReZolve-ER™ working solution.
- vi. Observe samples using fluorescence technique of choice.

Note 2: Optimal staining may vary between cell lines. Staining conditions may be modified according to cell type.

Note 3: For epifluorescence applications, ReZolve-ER™ can be excited at approximately 365 nm (UV) or 405 nm. For confocal and two-photon applications, it can be excited at 400 nm and 800-830 nm, respectively.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM