

# MitoBright IM Red for Immunostaining *Technical Manual*

## General Information

Cells have intracellular organelles that function to produce energy, synthesize and degrade proteins. Typical organelles include the nucleus, mitochondria, endoplasmic reticulum, and lysosomes. Although research on the functions of these organelles has been actively conducted for some time, recent developments in imaging technology, such as super-resolution imaging, have revealed that not only the functions of each organelle but also the coordination among organelles is important. One such example is the mitochondria-associated endoplasmic reticulum membrane, which is the interface between mitochondria and the endoplasmic reticulum (ER). Visualization of mitochondria is necessary to observe the interaction between mitochondria and other organelles. Fluorescent protein labeling, immunostaining, and mitochondria-selective small molecule fluorochromes are commonly used for this purpose, although these methods have problems. The fluorescent protein labeling method requires gene transfer, and multiple staining may not be able to perform in the immunostain method depending on the animal species of the primary antibody. In addition, commercially available mitochondria-selective small molecule fluorescent dyes do not have good retention properties and are difficult to use in immunostaining.

To overcome the problems above, Dojindo has developed a new mitochondria-selective small-molecule fluorescent dye, MitoBright IM Red, which covalently binds to proteins and other biomolecules. MitoBright IM Red, therefore, is able to simultaneously perform the staining mitochondria and the immunostaining.

## Contents

	1 Tube	3 Tubes
MitoBright IM Red for Immunostaining	20 $\mu$ l x 1	20 $\mu$ l x 3

## Storage Condition

Store at -20°C

## Required Equipment and Materials

- Growth medium or HBSS (Hanks balanced salt solution)
- Micropipettes

## Preparation of Solutions

### Preparation of MitoBright IM Red working solution

Dilute the 1 mmol/l MitoBright IM Red solution with growth medium (serum-free) to prepare a 1  $\mu$ mol/l MitoBright IM Red working solution.

**Note: The MitoBright IM Red working solution cannot be stored and must be freshly prepared each day and used within the day.**

## General protocol

1. Seed cells on a dish. Culture the cells at 37°C overnight in a 5% CO<sub>2</sub> incubator.
2. Discard the supernatant and wash the cells once with growth medium.
3. Add the 1  $\mu$ mol/l MitoBright IM Red working solution to the dish and incubate at 37°C for 30 minutes in a 5% CO<sub>2</sub> incubator.
4. Discard the supernatant, and wash the cells once with growth medium.
5. Perform immunofluorescence staining.
6. Observe the cells under a fluorescence microscope.

## Usage Examples

### Multiple staining of MitoBright IM Red with Tom20 as a mitochondrial marker in HeLa cells.

1. HeLa cells (2.4 × 10<sup>4</sup> cells/well, 200  $\mu$ l) in MEM (10% FBS) were seeded in a  $\mu$ -slide 8 well plate and cultured at 37°C overnight in a 5% CO<sub>2</sub> incubator.
2. After removing the supernatant, the cells were washed once with 200  $\mu$ l of MEM.
3. MitoBright IM Red working solution (1  $\mu$ mol/l, 200  $\mu$ l) in MEM was added, and the cells were incubated at 37°C for 30 minutes in a 5% CO<sub>2</sub> incubator.
4. After removing the supernatant, the cells were washed twice with 200  $\mu$ l of MEM.
5. Paraformaldehyde (4% v/v) in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 15 minutes.
6. After removing the supernatant, the cells were washed three times with 200  $\mu$ l of PBS.
7. Triton X-100 (0.1% v/v) in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 10 minutes.
8. After removing the supernatant, 10% (v/v) Blocking One-P in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 1 hour.
9. After removing the supernatant, anti-Tom20 antibody (1  $\mu$ g/ml, 200  $\mu$ l) was added, and the cells were incubated at room temperature for 1 hour.
10. After removing the supernatant, the cells were washed three times with 200  $\mu$ l of PBS.
11. Alexa647 labeled anti-mouse IgG (2  $\mu$ g/ml, 200  $\mu$ l) was added, and the cells were incubated at room temperature for 1 hour.
12. After removing the supernatant, the cells were washed twice with 200  $\mu$ l of PBS.
13. The cells were observed under a confocal laser scanning microscope (LSM 800, Zeiss).

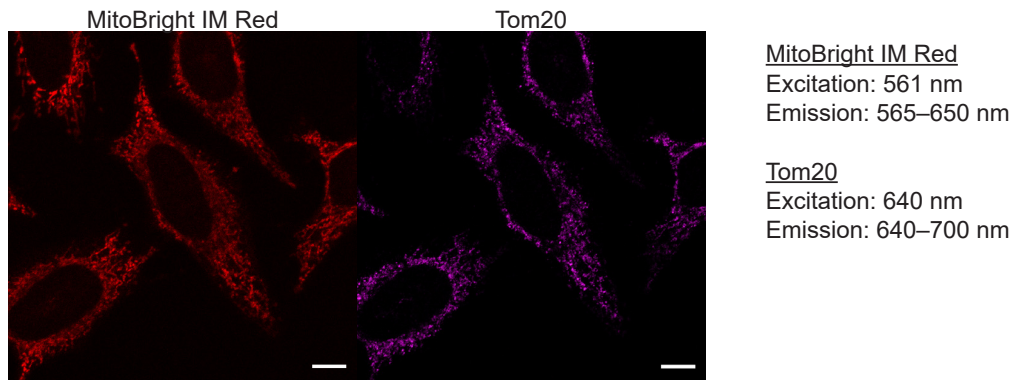


Figure 1 Fluorescence images using MitoBright IM Red and Tom20.

#### Multiple staining of MitoBright IM Red with KDEL as an ER marker in HeLa cells.

1. HeLa cells ( $2.4 \times 10^4$  cells/well, 200  $\mu$ l) in MEM (10% FBS) were seeded in a  $\mu$ -slide 8 well plate and cultured at 37°C overnight in a 5% CO<sub>2</sub> incubator.
2. After removing the supernatant, the cells were washed once with 200  $\mu$ l of MEM.
3. MitoBright IM Red working solution (1  $\mu$ mol/l, 200  $\mu$ l) in MEM was added, and the cells were incubated at 37°C for 30 minutes in a 5% CO<sub>2</sub> incubator.
4. After removing the supernatant, the cells were washed twice with 200  $\mu$ l of MEM.
5. Paraformaldehyde (4% v/v) in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 15 minutes.
6. After removing the supernatant, the cells were washed three times with 200  $\mu$ l of PBS.
7. Triton X-100 (0.1% v/v) in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 10 minutes.
8. After removing the supernatant, 10% (v/v) Blocking One-P in PBS (200  $\mu$ l) was added, and the cells were incubated at room temperature for 1 hour.
9. After removing the supernatant, anti-KDEL antibody (1  $\mu$ g/ml, 200  $\mu$ l) was added, and the cells were incubated at 4°C for overnight.
10. After removing the supernatant, the cells were washed three times with 200  $\mu$ l of PBS.
11. Alexa488 labeled anti-mouse IgG solution (2  $\mu$ g/ml, 200  $\mu$ l) was added, and the cells were incubated at room temperature for 1 hour.
12. After removing the supernatant, the cells were washed twice with 200  $\mu$ l of PBS.
13. The cells were observed under a confocal laser scanning microscope (LSM 800, Zeiss).

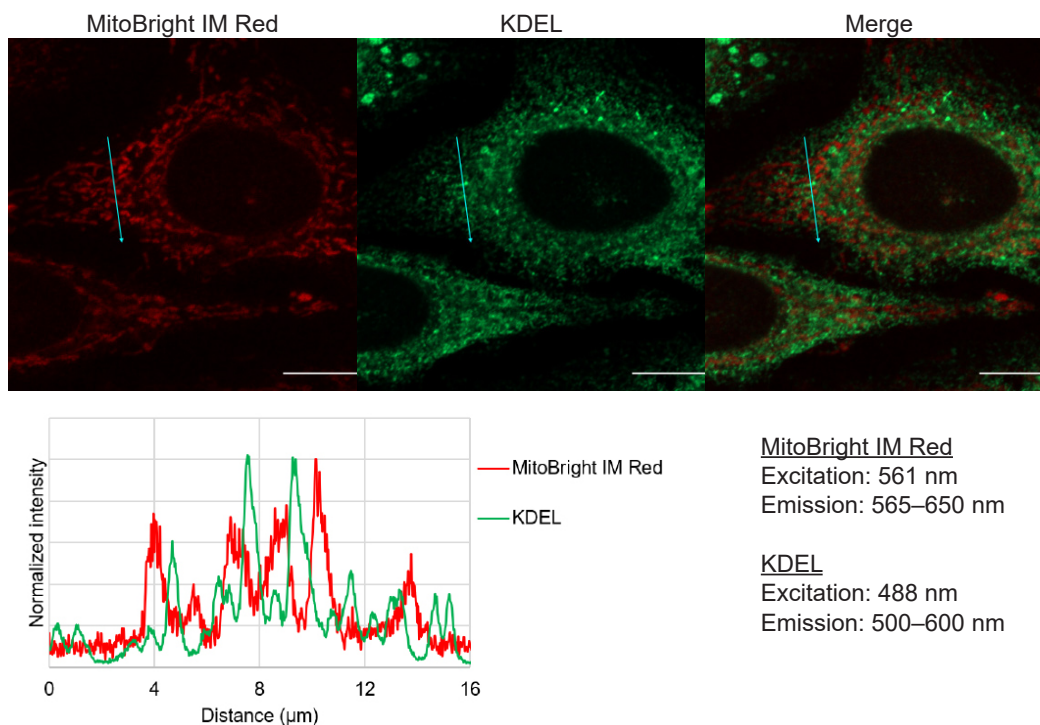


Figure 2 Multiple staining images and line profiles of MitoBright IM Red and KDEL.

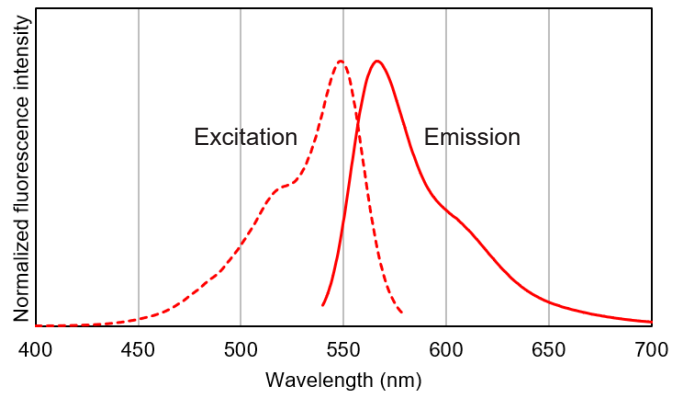


Figure 3 Excitation and emission spectra of MitoBright IM Red.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories  
2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto  
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525  
E-mail: [info@dojindo.co.jp](mailto:info@dojindo.co.jp) Web: [www.dojindo.co.jp](http://www.dojindo.co.jp)

Dojindo Molecular Technologies, Inc.  
Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>  
Dojindo EU GmbH  
Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>  
Dojindo China Co., Ltd  
Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>  
Tel: +86-10-8225-1765