

München, 1. August. 2019

Customer Notification

Dear Valued Customer,

As we continue to develop new products for offering better products and experience to customers, we have decided to discontinue the sales of MitoBright Dyes (MitoBright Green (Code: MT06), MitoBright Red (Code: MT07), MitoBright Deep Red (Code: MT08)) as of 1. January 2020.

However, we will continue to offer these products until 31. December 2019, with the service to support customer's experiment.

Discontinued Product		
Code	Name	Unit
MT06-10	MitoBright Green	50 µg x 3
MT07-10	MitoBright Red	50 µg x 3
MT08-10	MitoBright Deep Red	50 µg x 3

We are also glad to announce the new Mitochondria Staining Dyes, which are direct and even better functional replacements for the discontinued products have been released.

For detailed information of these new products, please visit the product page at www.dojindo.eu.com

Also, for further inquiries and technical questions are anytime welcome at info@dojindo.eu.com

Alternative Product		
Code	Name	Unit
MT10-10	MitoBright LT Green	20µL
MT10-12		400µL
MT10-14		400µL x 3
MT11-10	MitoBright LT Red	20µL
MT11-12		400µL
MT11-14		400µL x 3
MT12-10	MitoBright LT Deep Red	20µL
MT12-12		400µL
MT12-14		400µL x 3

We appreciate in advance your understanding.

Yours sincerely,

Dojindo EU GmbH