

# Analysis for Cellular Function

**Mitochondrial Analysis**

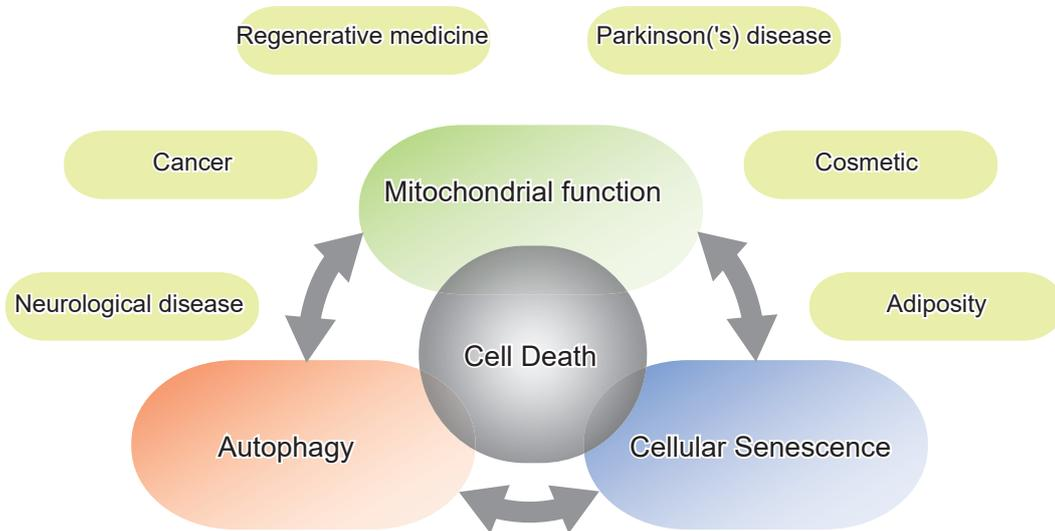
**Autophagy**

**Cellular Senescence**

**Cell Proliferation Cytotoxicity**

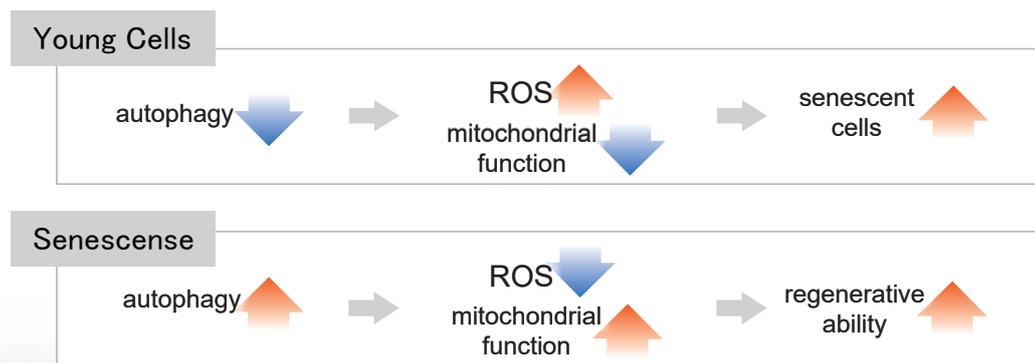
## Analysis for Cellular Function

Understanding for cellular functions are important to reveal the mechanism of diseases and to develop a drug-discovery. Especially autophagy, cellular senescence and mitochondrial function have been found to be closely related each other, and have been attracting attention in the medical research and cosmetics industry.



## Suppression of Cellular Senescence by Autophagy

As an example where autophagy, cellular senescence and mitochondrial function are closely related, there is an example using mouse stem cells. Muscle stem cells that remain resting stage of the cell cycle is dependent on autophagy. When the function of autophagy declines, the reactive oxygen species (ROS) increases due to mitochondrial dysfunction. As a result, it becomes cellular senescence. Even autophagy is inhibited in young cells, the cellular senescence is induced, conversely by inducing autophagy in senescent cells, the mitochondrial function is restored and recovery of regenerative ability has been confirmed.

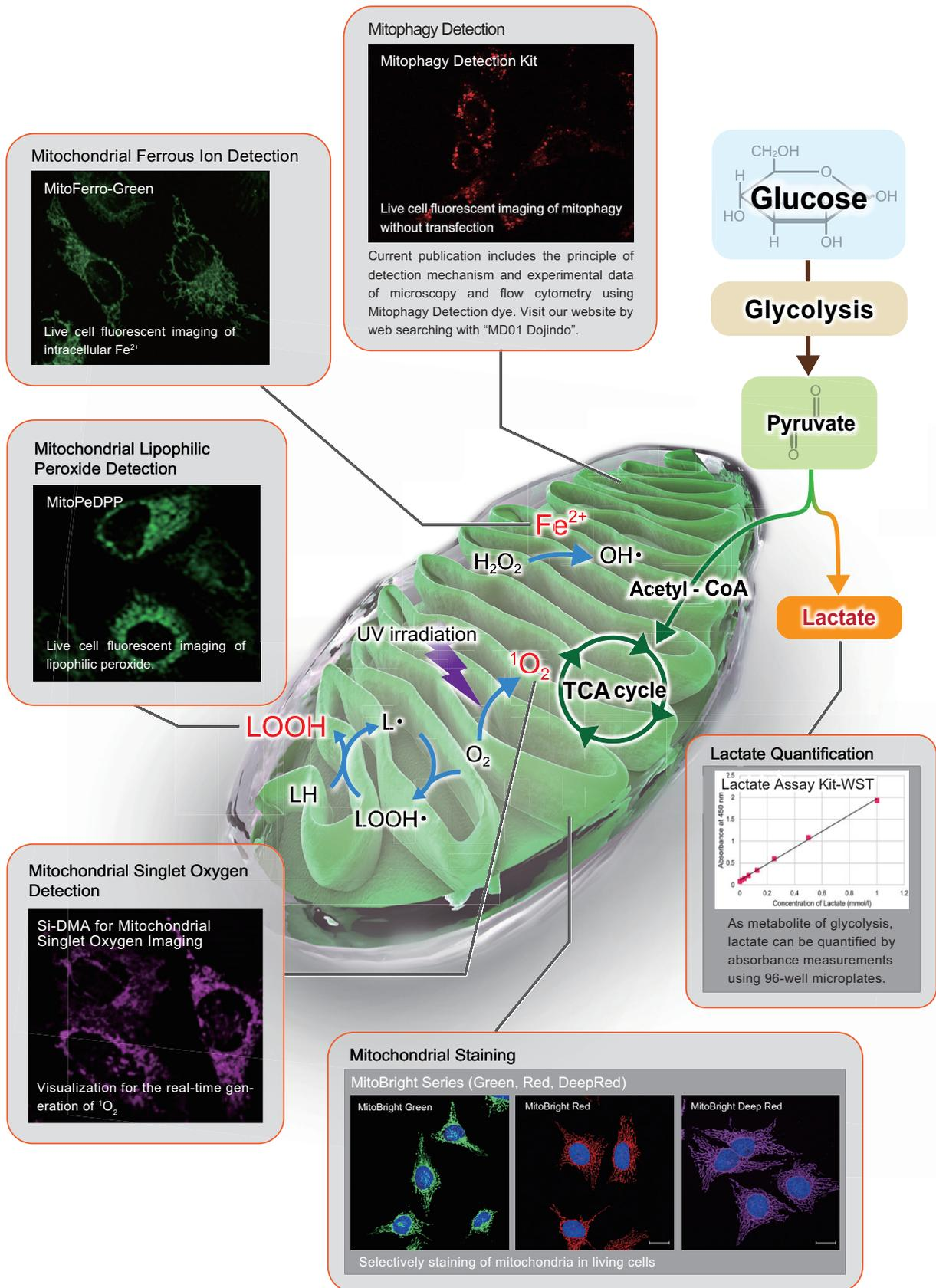


L. García-Prat, M. Martínez-Vicente, E. Perdiguerro, L. Ortet, J. Rodríguez-Ubreva, E. Rebollo, V. Ruiz-Bonilla, S. Gutarra, E. Ballestar, A. L. Serrano, M. Sandri & P. Muñoz-Cánoves, "Autophagy maintains stemness by preventing senescence." *Nature*, 2016, 529 (7584), 37

## Table of Contents

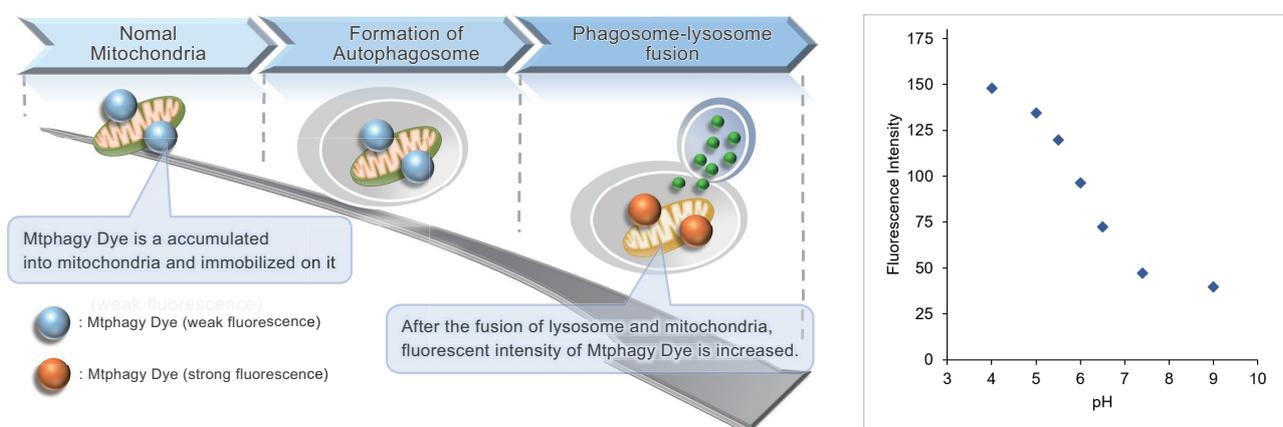
<b>Reagents for Mitochondrial Research</b>	...	3
<b>Mitophagy Detection Kit</b> Mitophagy Detection Kit	...	4
<b>Mitochondrial Iron Detection Reagent</b> Mito-FerroGreen	...	6
<b>Mitochondrial Lipophilic Peroxide Detection Reagent</b> MitoPeDPP	...	7
<b>Mitochondrial Singlet Oxygen Detection Reagent</b> Si-DMA for Mitochondrial Singlet Oxygen Imaging	...	8
<b>Mitochondrial Staining Reagent</b> MitoBright Series (Green / Red / DeepRed)	...	9
<b>Lactate Quantification Kit</b> Lactate Assay Kit	...	10
<b>Autophagy Detection Reagent</b> Autophagy Detection Kit	...	11
<b>Cellular Senescence Assay Kit</b> Cellular Senescence Detection Kit - SPiDER-βGal	...	15
<b>Cell Proliferation / Cytotoxicity</b> Cell Counting Kit-8	...	18
<b>Cytotoxicity Assay Kit</b> Cytotoxicity LDH Assay Kit-WST	...	19

# Reagents for Mitochondrial Research



# Mitophagy Detection Kit

This kit is composed of Mtpagy Dye, reagent for detection of mitophagy, and Lyso Dye. Mtpagy Dye accumulates in intact mitochondria, is immobilized on it with chemical bond and exhibits a weak fluorescence from the influence of surrounding condition. When Mitophagy is induced, the damaged mitochondria fuses to lysosome and then Mtpagy Dye emits a high fluorescence. To confirm the fusion of Mtpagy Dye-labeled mitochondria and lysosome, Lyso Dye included in this kit can be used.



The fluorescent intensity of Mtpagy Dye is increased at pH4-5.

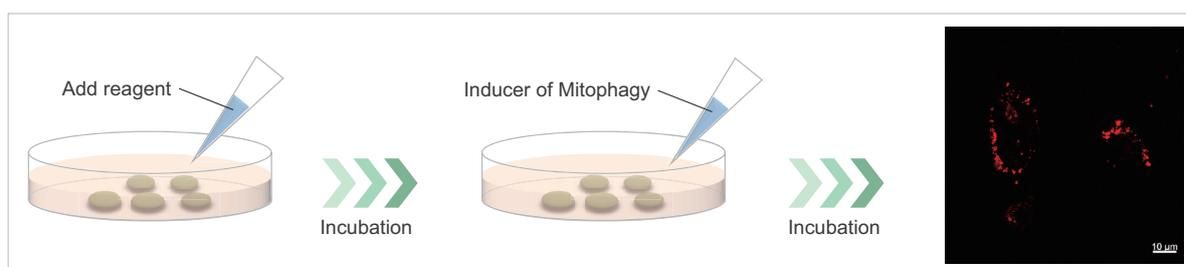
The publication of Mitophagy Detection Kit includes the detection principle and experimental data.

H. Iwashita, S. Torii, N. Nagahora, M. Ishiyama, K. Shioji, K. Sasamoto, S. Shimizu and K. Okuma, "Live Cell Imaging of Mitochondrial Autophagy with a Novel Fluorescent Small Molecule", *ACS Chem. Biol.*, **2017**, *12*, 2546.

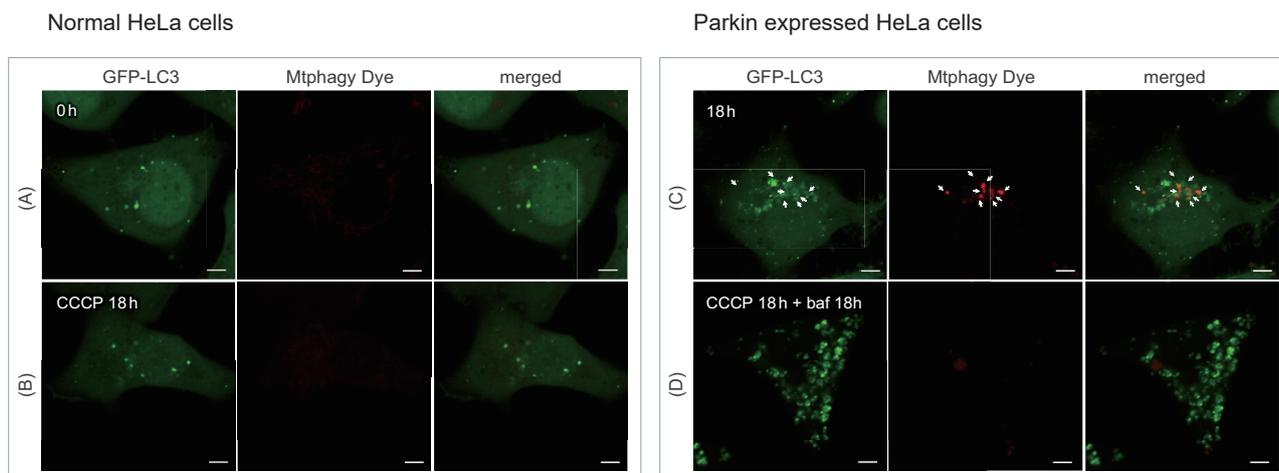
※ Other publications are listed online. Please visit at [www.dojindo.com](http://www.dojindo.com).

## Simple procedure

Transfection procedure is not need for mitophagy imaging.



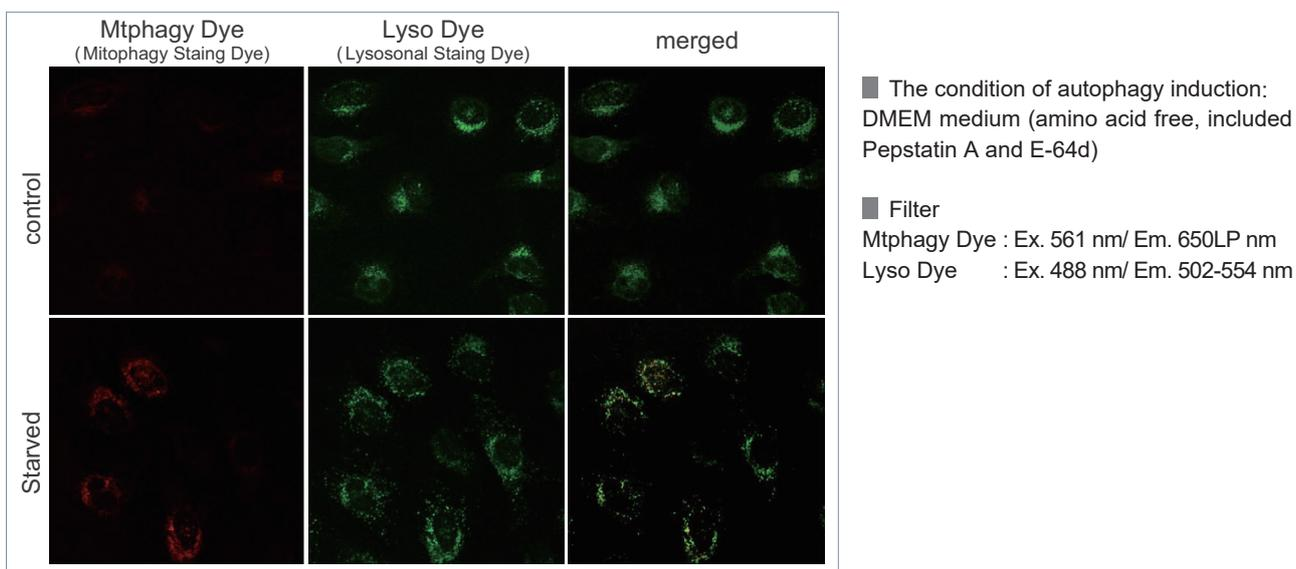
## Comparison with autophagy marker



CCCP (carbonyl cyanide m-chlorophenyl hydrazone) has been added to normal and Parkin expressed cells. The strong fluorescence was not observed in normal HeLa cells(A)(B). On the other hand, the strong fluorescence was shown in Parkin expressed cells in 18 hours after addition of CCCP(C). Some of the puncta are co-localized with the autophagy marker(GFP-LC3). In addition, suppressed fluorescence of Mtpahgy dye was observed when autophagy inhibitor, bafilomycin was added to Parkin expressed cells(D). Because lysosomal pH was increased by the addition of bafilomycin.

## Mitophagy detection in starved cells

Living HeLa cells are co-stained with Mtpahgy Dye and Lyso Dye under mitophagy induced condition.



Mtpahgy Dye was added to HeLa cells and the cells were incubated for 6 hours under starved condition. Before observation of the cells, lysosomal staining dye, Lyso Dye was added to HeLa cells. The fluorescent intensity of Mtpahgy Dye was increased in the starved HeLa cells but not in normal cells. In addition, Mtpahgy Dye was co-localized with Lyso Dye in the starved cells.

Discription	Unit	Code
Mitophagy Detection Kit	1 set	MD01

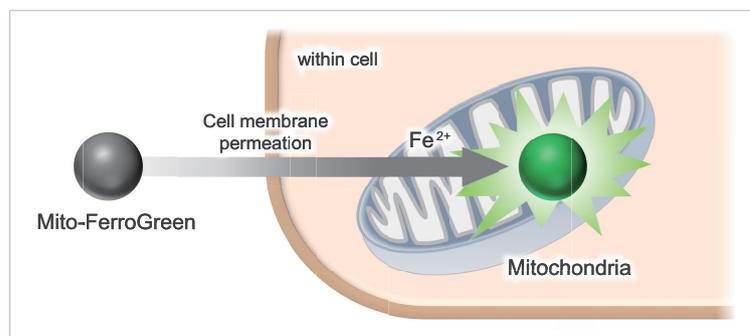
# Mitochondrial Iron Detection Reagent

## Mito-FerroGreen

### Cell permeability and Fe<sup>2+</sup> selectivity

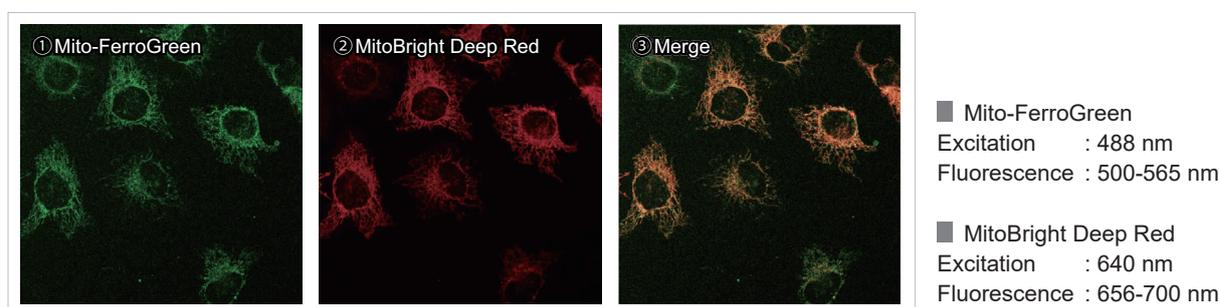
It is reported that iron is the most abundant transition metal element within an organism and shows various physiological activities. Recently, free iron in living cells is getting attention because its high reactivity is suggested to be related to cellular damage or death. Free iron exists in its stable redox states, ferrous ion (Fe<sup>2+</sup>) and ferric ion (Fe<sup>3+</sup>). In living cells, it is considered that understanding the behavior of Fe<sup>2+</sup> is more important than that of Fe<sup>3+</sup> because of the intracellular reductive environment, metal transporters and water solubility of Fe<sup>2+</sup>. Mito-FerroGreen is a novel fluorescent probe for the detection of ferrous ion (Fe<sup>2+</sup>) in mitochondria where Fe-S clusters and heme proteins are synthesized, and enables live cell fluorescent imaging of intracellular Fe<sup>2+</sup>.

This product was commercialized under the advisory of Dr. Hideko Nagasawa and Dr. Tasuku Hirayama (Gifu Pharmaceutical University).



### Double staining with mitochondrial staining probe

Co-localized Mito-FerroGreen and mitochondrial staining dye, MitoBright Deep Red (Dojindo, Code: MT08) were observed after addition of iron(II) in living HeLa cells.



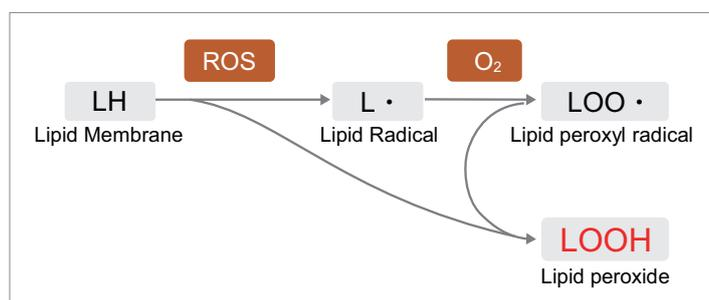
The final concentration of 5 μmol/l Mito-FerroGreen and 200 nmol/l MitoBright Deep Red were added to HeLa cells and the cells were incubated in a 5% CO<sub>2</sub> incubator for 30 minutes. Then, the final concentration of 100 μmol/l ammonium iron(II) sulfate was added to the cells and the cells were incubated in a 5% CO<sub>2</sub> incubator for 1 hour. After incubation, the cells were observed by a confocal fluorescent microscopy.

Discription	Unit	Code
Mito-FerroGreen	50 μg × 2	M489

# Mitochondrial Lipophilic Peroxide Detection Reagent

## MitoPeDPP

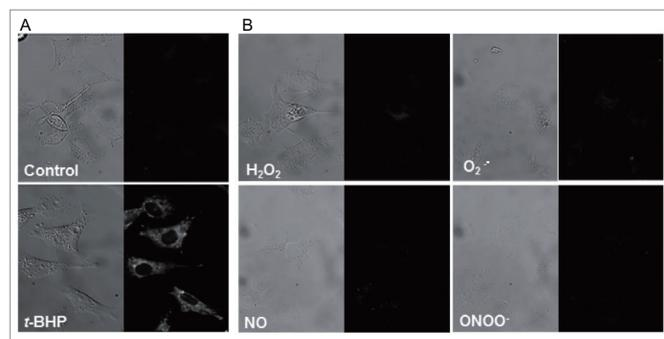
MitoPeDPP is a newly developed fluorescent dye which can penetrate cell membranes and is accumulated in the mitochondria. MitoPeDPP accumulated in mitochondrial inner membranes is oxidized by lipophilic peroxide and then emits strong fluorescence. Since the excitation and emission wavelengths of the oxidized MitoPeDPP (Ox-MitoPeDPP) are 452 nm and 470 nm, respectively, photodamage and autofluorescence of the samples can be minimized. Therefore, MitoPeDPP can be applicable for imaging of the lipophilic peroxide in living cells under a fluorescence microscope. This product was developed by Dr. Shioji, Department of Chemistry, Fukuoka University.



Lipid peroxidation is the metabolic process under the oxidative stress such as the presence of reactive oxygen species (ROS). Lipid peroxide is an important compound which formed by the chain reaction of lipid radicals and is a marker of oxidative damage in organelle membranes.

## Lipophilic peroxide Selectivity

Even though MitoPeDPP reacts with various peroxides (H<sub>2</sub>O<sub>2</sub>, *t*-BHP, ONOO<sup>-</sup>) in homogeneous systems (data is not shown), the MitoPeDPP is specifically-oxidized by *t*-BHP in mitochondria (A) but not with ROS and RNS (B).



A: MitoPeDPP stained cells with *t*-BHP treatment (*t*-BHP) and without (control).

B: MitoPeDPP stained cells with ROS or RNS exposure.

\*As ROS generators, PMA (O<sub>2</sub><sup>-</sup>), NOC7(NO), and SIN-1(ONOO<sup>-</sup>) were used in the experiments.

\*Abbreviations

*t*-BHP : *tert*-butylhydroperoxide

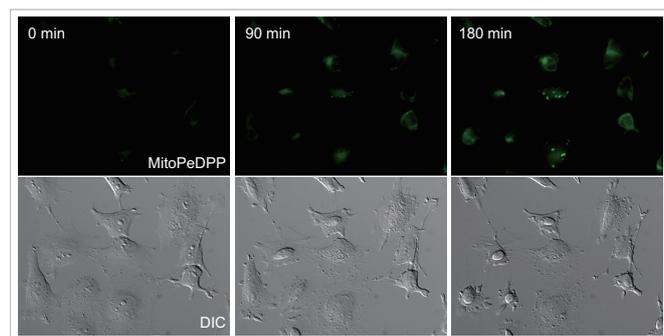
PMA : Phorbol Myristate Acetate

NOC7 : 1-Hydroxy-2-oxo-3-(N-methyl-3-aminopropyl)-3-methyl-1-triazine

SIN-1 : 3-(Morpholinyl)sydnonimine, hydrochloride

Emission : 470/40 nm, Excitation: 525/50 nm

## Detection of lipophilic peroxide generated with rotenone



The final concentration of 0.1 μmol/l MitoPeDPP was added to HeLa cells. The cells were then incubated in a 5% CO<sub>2</sub> incubator for 15 minutes. Then, the final concentration of 1 μmol/l rotenone was added to the cells and time-dependent fluorescence change was observed under a fluorescence microscope for 3 hours.

Emission : 470/40 nm

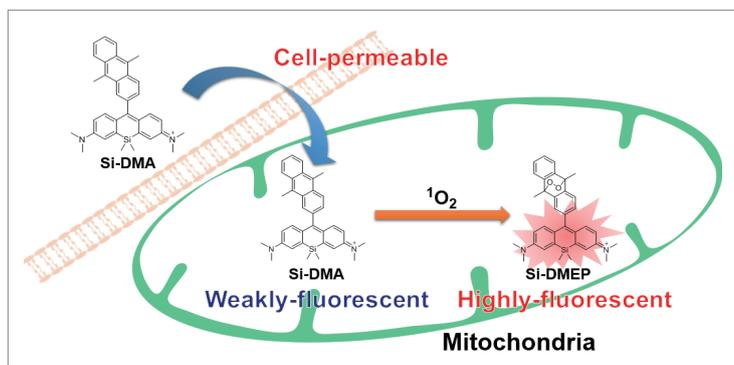
Excitation : 525/50 nm

Discription	Unit	Code
MitoPeDPP	1 set (5 μg×3)	M466

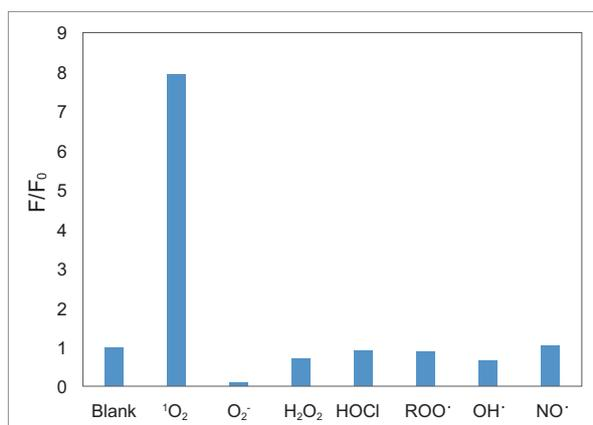
# Mitochondrial Singlet Oxygen Detection Reagent

## Si-DMA for Mitochondrial Singlet Oxygen Imaging

Majima *et al.* synthesized a new far-red fluorescence probe composed of silicon-containing rhodamine and anthracene moieties, namely Si-DMA, as a chromophore and a  $^1\text{O}_2$  reactive site, respectively. In the presence of  $^1\text{O}_2$ , fluorescence of Si-DMA increases 17 times due to endoperoxide formation at the anthracene moiety. Among seven different ROS, Si-DMA is able to selectively detect the  $^1\text{O}_2$ . In addition, Si-DMA is able to visualize the real-time generation of  $^1\text{O}_2$  from protoporphyrin IX in mitochondria with 5-aminolevulinic acid (5-ALA), a precursor of heme.



### Singlet oxygen selectivity

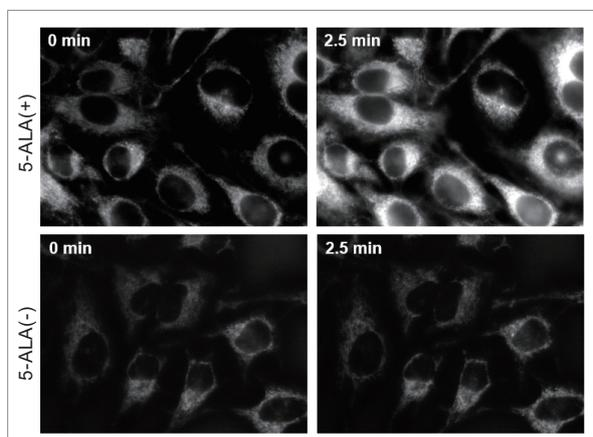


Si-DMA showed significant signal to singlet oxygen but not to other ROS and RNS.

#### Experimental condition

PBS(pH7.4) MeOH = 1 : 1  
 Si-DMA 1  $\mu\text{mol/l}$   
 ROS 10 mmol/l  
 RNS 10 mmol/l

### Singlet oxygen detection after the addition of 5-ALA



By the addition of 5-ALA(5-aminolevulinic acid) to HeLa cells, protoporphyrin IX was produced in mitochondria. Once the cells were irradiated under the excitation laser, singlet oxygen was generated. In the experiment, the generated singlet oxygen was detected by Si-DMA under a fluorescence microscope.

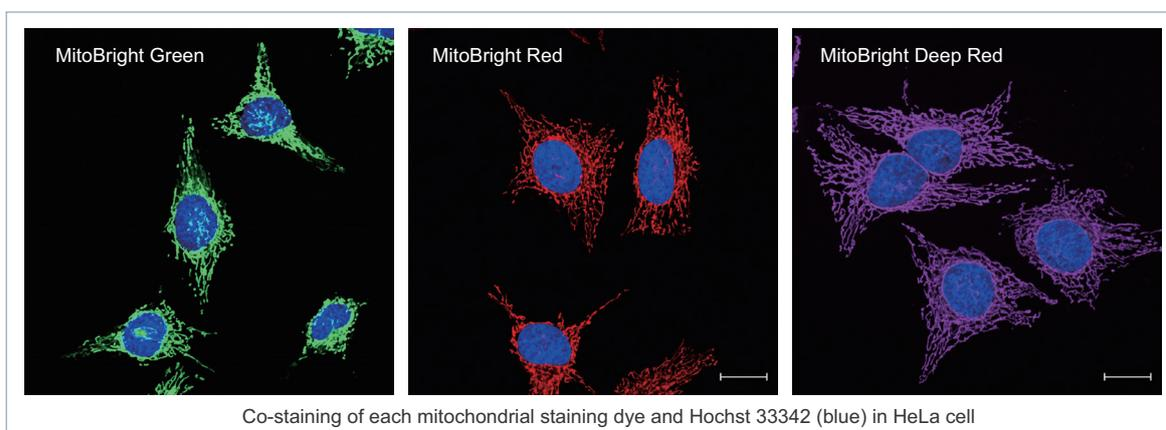
Emission : 575-625 nm  
 Excitation : 660-710 nm

Discription	Unit	Code
Si-DMA for Mitochondrial Singlet Oxygen Imaging	2 $\mu\text{g}$	MT05

# Mitochondrial Staining Reagent

## MitoBright Series Green / Red / DeepRed

There are various organelles playing an important role in cells. Mitochondria is not only a principal site of oxidative phosphorylation to produce ATP, but also an important organelle whose activity and dysfunction are relevant to cancer, cell senescent and neurodegenerative diseases such as Alzheimer's and Parkinson's. "MitoBright" is a fluorescent probe for selectively staining of mitochondria in living cells. It accumulates in healthy mitochondria due to the membrane potential dependency. In addition, MitoBright is well retained in mitochondria because of its covalent bond ability to proteins.



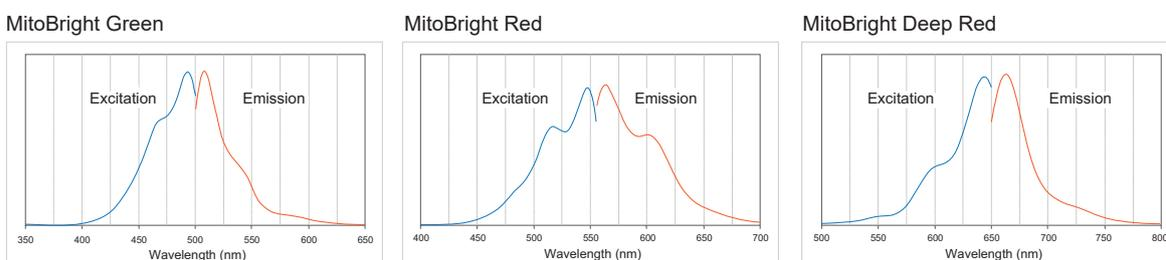
Co-staining of each mitochondrial staining dye and Hoechst 33342 (blue) in HeLa cell

### MitoBright series line-up

Description	$\lambda_{ex}$	$\lambda_{em}$	live imaging	staining before fixation*	staining after fixation
MitoBright Green	493 nm	508 nm	○	○	×
MitoBright Red	547 nm	563 nm	○	○	×
MitoBright Deep Red	643 nm	663 nm	○	○	×

※ PFA (paraformaldehyde) fixation

### Fluorescent Property



### Suitable unit size

The one tube (50  $\mu$ g) of each MitoBright series is corresponded more than 300 staining experiments when 35 mm dish is used.

Description	Unit	Code
MitoBright Green	50 $\mu$ g $\times$ 3	MT06
MitoBright Red	50 $\mu$ g $\times$ 3	MT07
MitoBright Deep Red	50 $\mu$ g $\times$ 3	MT08

# Lactate Quantification Kit

## Lactate Assay Kit-WST

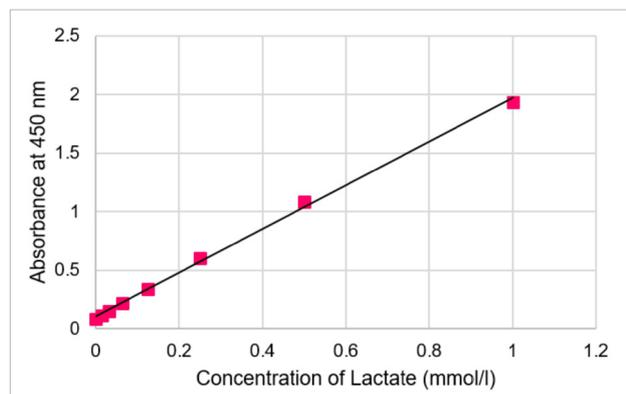
Lactate is a metabolite of glycolysis that is one of the main metabolic pathways in cells, and is known to be a biomarker for muscular fatigue and hyperlactacidemia. It also serves as a marker for monitoring the changes of intracellular metabolic pathways. In addition, recent metabolomic study suggests that lactate contributes as a major carbon source in the TCA cycle of tissues and cancer cells. Lactate Assay Kit-WST enables quantitation of lactate produced by glycolysis. This kit is formatted for 96-well microplate assays with a detection sensitivity limit of 0.02 mmol/l lactate.

### Simple procedure

After transferring the supernatant to a microplate, only the mixing with reagent and incubation procedures are needed.



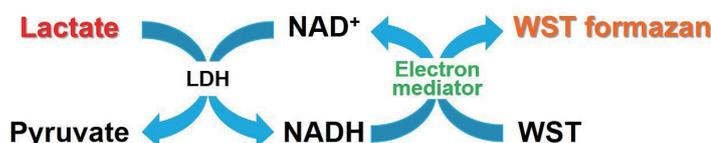
### Quantify through standard curve



The concentration of lactate in a sample is quantified with a standard curve which is prepared from the lactate standard in a kit. In case the concentration of lactate is over 1 mmol/l, the sample can be measured after a dilution procedure.

### Assay Principle

The kit has been optimized to quantitate lactate in cell culture supernatant by measuring the absorption derived from a colorimetric reaction of WST.



Discription	Unit	Code
Lactate Assay Kit-WST	50 tests	L256
	200 tests	

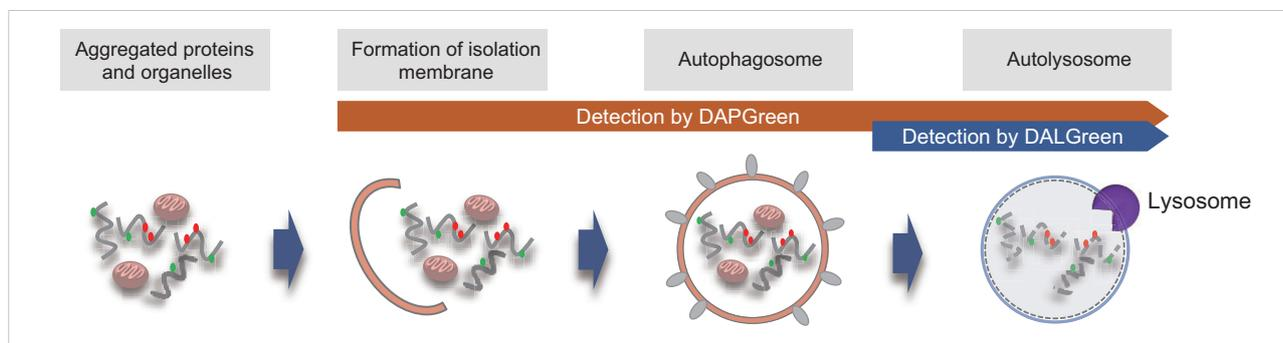
# Autophagy Detection Reagent

Detection of Autophagosome  
DAPGreen - Autophagy Detection

Detection of Autolysosome  
DALGreen - Autophagy Detection

## How they work

DAPGreen has been developed to monitor early-phase to late-phase autophagy, autophagosome and autolysosome. DALGreen is favorable for monitoring late-phase autophagy, autolysosome. With these probes that stain autophagosomes as they are being formed, the real-time change of autophagic phenomena of live cells may be traced.



**Autophagosome Detection Reagent DAPGreen**

Fluorescence enhanced DAPGreen

Autophagosome membrane      Autolysosome membrane

When an autophagosome membrane is formed, DAPGreen is incorporated inside of the membrane. The fluorescence of incorporated DAPGreen is enhanced under lipophilic condition. The analysis of DAPGreen also has a high correlation with that of LC3 which is a well-known autophagy marker. For details, please refer to the experimental data of DAPGreen.

**Autolysosome Detection Reagent DALGreen**

Fluorescence enhanced DAPGreen

Autophagosome membrane      Autolysosome membrane

In the same way as DAPGreen, DALGreen is incorporated inside of the autophagosome membrane when the membrane is formed. The fluorescence of DALGreen is enhanced under acidic condition after the autophagosome is fused with the lysosome.

## References

The following paper includes the detail principle of DAPGreen and DALGreen. The co-staining results with a lysosomal staining dye and the staining data in ULK1/2\* double-knockout cells are also described in this article.

\*ULK1/2: Uncoordinated 51-like kinases 1 and 2 which plays for homologous function of yeast Atg1.

H. Iwashita, H. T. Sakurai, N. Nagahora, M. Ishiyama, K. Shioji, K. Sasamoto, K. Okuma, S. Shimizu, and Y. Ueno, "Small fluorescent molecules for monitoring autophagic flux", *FEBS Lett.*, **2018**, 592(4), 559.

## General information

Below table guides you to choose the most suitable application for you.

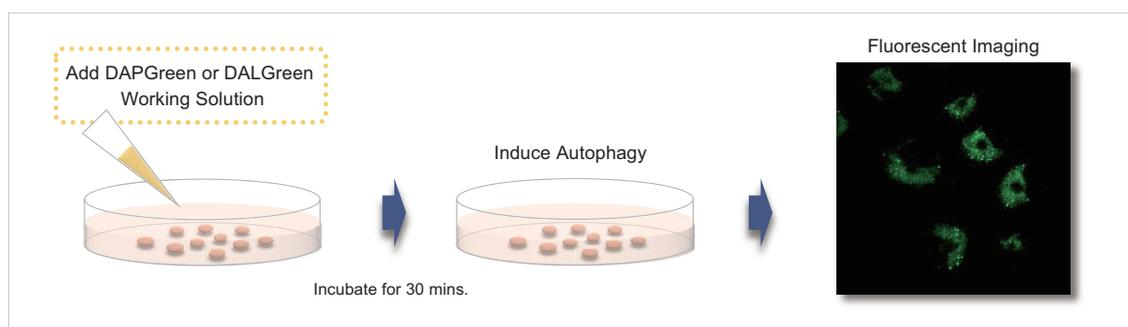
	Applicable instruments			Fluorescent property	Volume / the number of usable assays	Existing methods
	Fluorescent Microscope	Flow cytometer	Microplate reader			
DAPGreen	○	○	○	Ex. 425-475 Em. 500-560 *For confocal microscope, the sample can be excited at 488 nm	5 nmol x 1 / 35 mm dish x 25 (when used with max. concentration)	LC3-GFP MDC Cyto-ID etc.
DALGreen	○	○	×	Ex. 350-450 Em. 500-560 *For confocal microscope, the sample can be excited at 488 nm	20 nmol x 1 / 35 mm dish x 10 (when used with max. concentration)	LC3-GFP-RFP etc.

\*Fluorescent property is available on Dojindo HP.

\*Double staining with DAPGreen and DALGreen is not possible, unfortunately. If you like to double stain autophagosome and autolysosome in different color, please contact us.

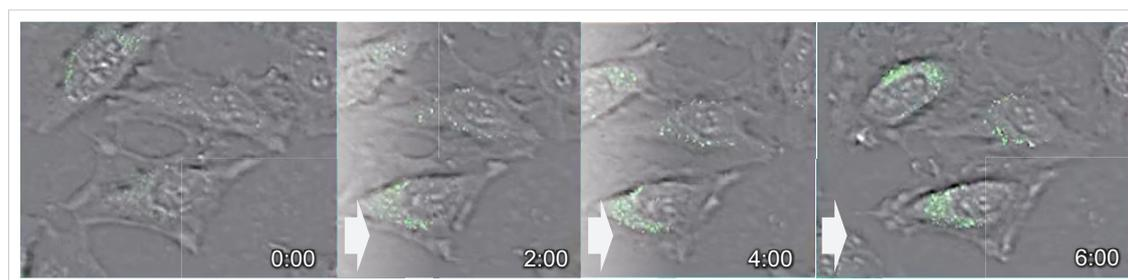
## Without genetic engineering

DAPGreen and DALGreen do not require any transfection. You only need to add the reagent to cultured cells.



## Time-lapse imaging of autophagy

Six hours of time-lapse imaging was taken using DALGreen and starved HeLa cells.



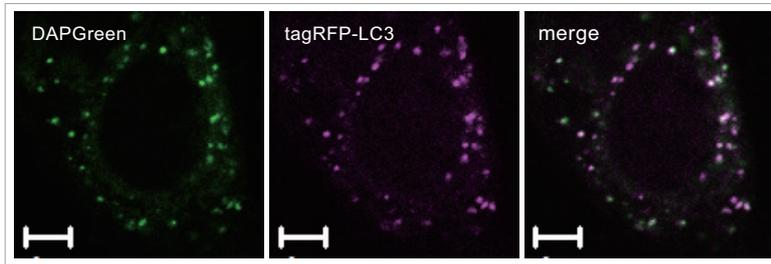
### Imaging condition

Medium : Amino acids-free culture medium  
 Machine : Confocal Quantitative Image Cytometer ( YOKOGAWA Electric Corporation : CQ1 )  
 Fluorescence Filters : Ex. 405 / Em. 525/50,  
 Magnification : 20X

The imaging movie has been uploaded on Dojindo HP.

## Good correlation with LC3

The HeLa cells were double stained by DAPGreen and tagRFP-LC3 to determine their colocalization.



### Result

Almost all of DALGreen signals were colocalized with LC3 reasonably.

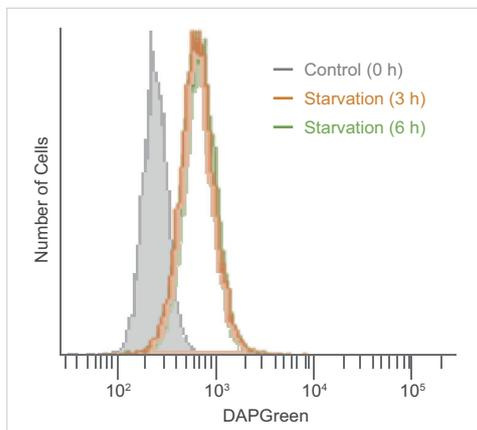
### Imaging condition

DAPGreen : Ex. 488 nm / Em. 500-563 nm  
Scale bar : 10  $\mu$ m

### Autophagy induced condition

After adding DAPGreen, RFP-LC3 expressed HeLa cells were treated with rapamycin. The images were obtained after 4 hours of incubation.

## Quantitative analysis using flow cytometer



### Result

Starved HeLa cells showed strong fluorescent signal after 3 hours of incubation.

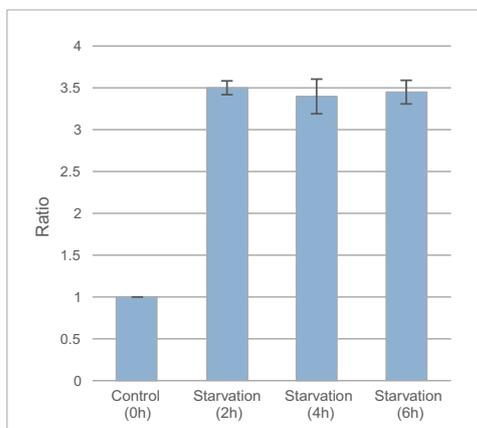
### Detection

Wavelengths: Ex. 488 nm / Em. 500-560 nm

### The condition of autophagy induction

After staining with DAPGreen, HeLa cells were incubated for 0, 3, 6 hours with amino acid-free medium.

## Quantitative analysis using microplate reader



### Result

After 2 hours of incubation under starved condition, enhanced fluorescence was observed. (3.5 times stronger than "Control")

### Detection

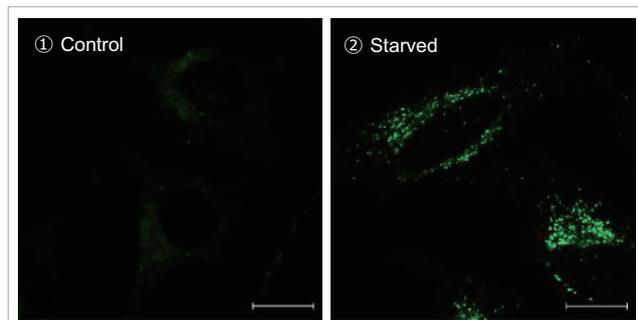
Wavelengths: Ex. 450 nm / Em. 535 nm

### The condition of autophagy induction

After staining with DAPGreen, HeLa cells were incubated for 0, 2, 4, 6 hours with amino acid-free medium.

## Autolysosome analysis by microscopy

Once DALGreen is added, HeLa cells were incubated either with a growth medium(1) or an amino acid-free medium(2).



### Result

The strong fluorescence of DALGreen was observed as puncta in starved HeLa cells.

### Imaging condition

Detection : Ex.488 nm / Em.500-563 nm

Scale bar : 20  $\mu$ m

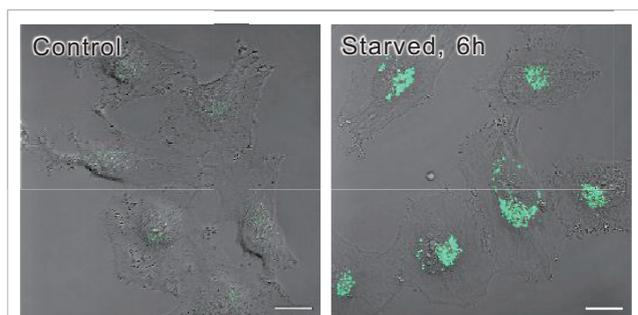
### The condition of autophagy induction

① Control: 6 hours incubation in growth medium

② Starved: 6 hours incubation in amino acid-free growth medium

## Comparison with MDC

Autophagy in starved HeLa cells was observed with DALGreen and MDC(Monodansylcadaverine).



Dye concentration : 1  $\mu$ mol/l  
 Incubation time : 30 min  
 Detection : Confocal fluorescence microscope  
 Ex. 488 nm/ Em. 500-563 nm  
 Scale bar : 40  $\mu$ m  
 The reagent was added **BEFORE** starvation for 6 hours

### Results

The fluorescent signal of DALGreen was increased in the starved HeLa cells, whereas MDC showed low S/N ratio.

### Advantage

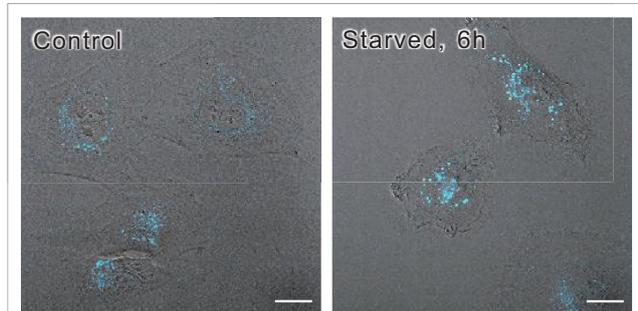
MDC requires excitation by ultraviolet which can damage living cells. On the other hand, DALGreen can be excited at a longer wavelength.

### The condition of autophagy induction

DALGreen : Addition of reagent  $\Rightarrow$  Induction of starvation

MDC : Induction of starvation  $\Rightarrow$  Addition of reagent

DALGreen can be used for monitoring autophagy because the reagent is added prior to induction of autophagy unlike MDC.



Dye concentration : 1  $\mu$ mol/l  
 Incubation time : 30 min  
 Detection : Confocal fluorescence microscope  
 Ex. 405 nm/ Em. 450-546 nm  
 Scale bar : 40  $\mu$ m  
 The reagent was added **AFTER** starvation for 6 hours

Discription	Unit	Code
 DALGreen - Autophagy Detection	20 nmol	D675
 DAPGreen - Autophagy Detection	5 nmol	D676

# Cellular Senescence Assay Kit

## Cellular Senescence Detection Kit - SPiDER-βGal

### Comparison with conventional method, X-Gal

Since X-Gal requires a counting by colorimetric visualization under a microscope, it is not suitable as a quantitative detection. Dojindo's β-galactosidase substrate, SPiDER-βGal allows fluorescent co-staining and quantitative analysis through flow cytometer, image cytometer and so on.

#### X-Gal

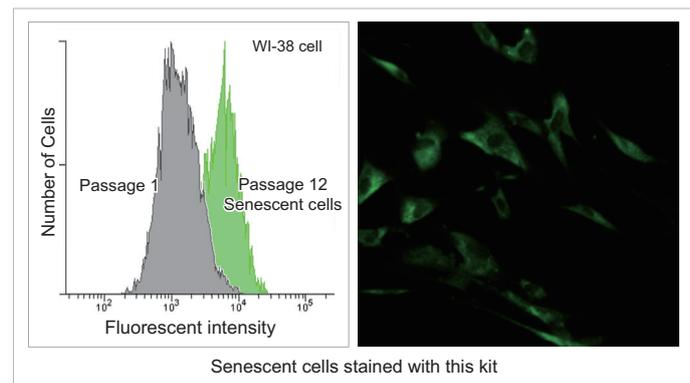
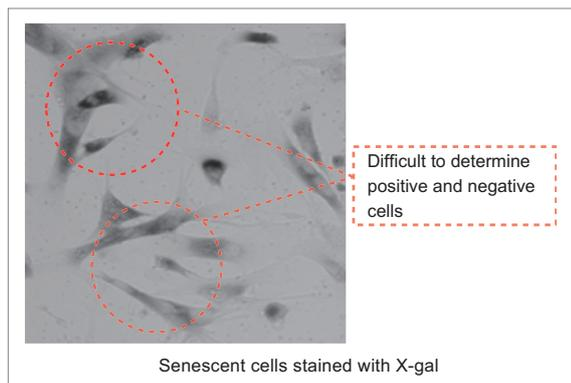
Visual counting required

- Disadvantages : - Personal error
- Time consuming procedure

#### SPiDER-βGal

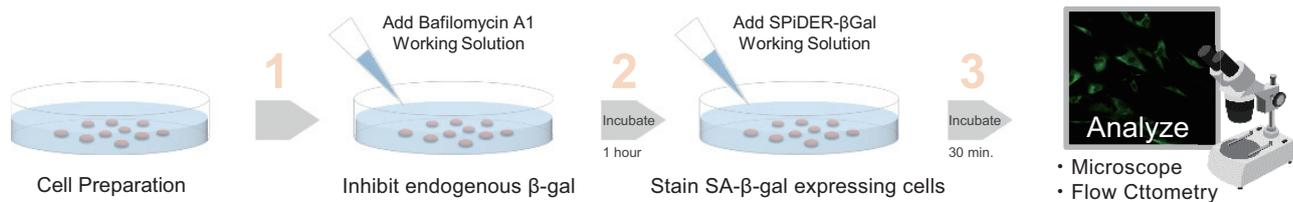
Reliable analysis by flowcytometry

- Advantages : - Quantitative method
- Simple procedure



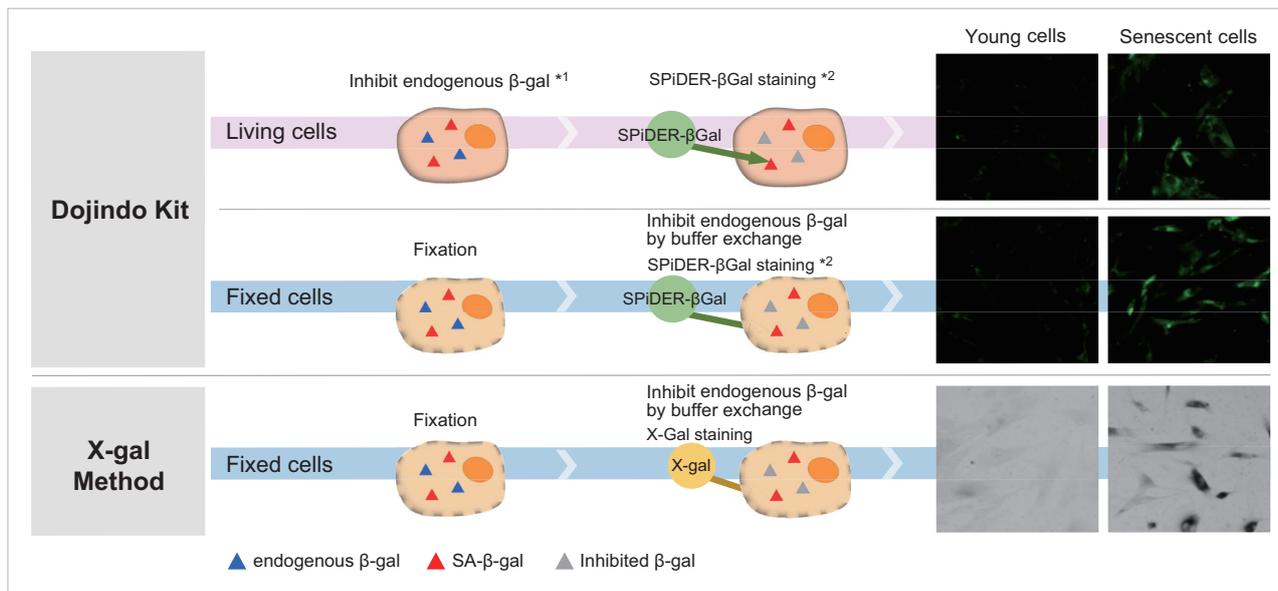
### Simple Procedure

The kit is applicable for both living and fixed cells. The SA-β-gal staining can be done in only 30 minutes. For living cells, just need to add 2 types of reagent included in the kit to cultured cells. If fixation is needed, it can be performed before staining.



**This kit is applicable for both living and fixed cells.**

The staining principles are shown in below.



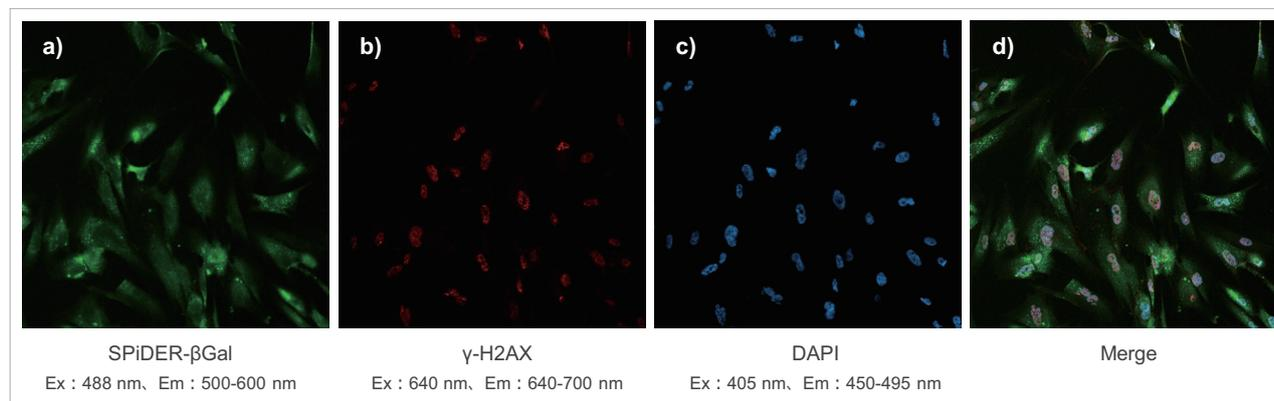
**\*1 Why endogenous β-galactosidase suppressed?**

Normally endogenous β-galactosidase exist in lysosome where acid pH enviloment. To detect acculate SA-β-gal selectively, Bafilomycin A1(included in the kit) works to supress such endogenous β-galactosidase by pH neutralization.

**\*2 How SPiDER-βGal beneficial compared with other small molecules?**

SPiDER-βGal has better cell permeability and retention ability than other commercially available dyes such as FDG and C12FDG. Therefore, it enables high sensitive fluorescence imaging.

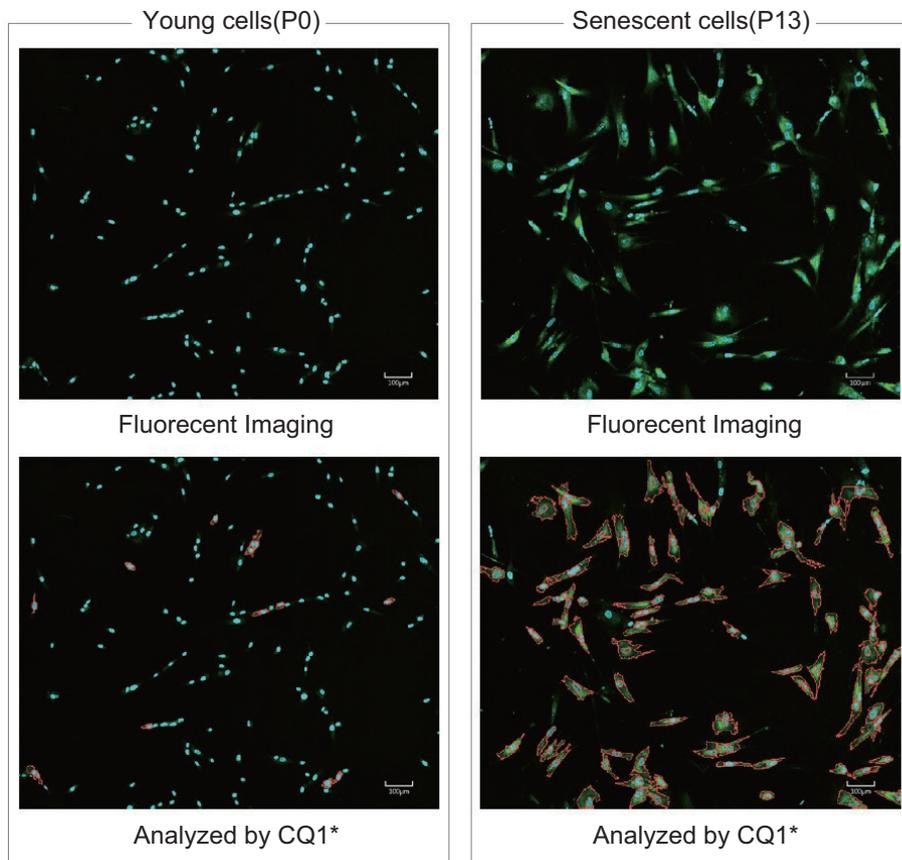
**Co-staining with other cellular senescence markers**



In this experiment, WI-38 cells were used after ten times of subculture (Passage 10) as the model of senescent cells. The cells were stained with a) this kit for SA-β-gal detection, b) anti γ-H2AX antibody for DNA damage detection, and DAPI for nuclear staining. In the results, both senescent markers, SA-β-gal and γ-H2AX staining were observed. The experimental procedure is informed at the product HP.

## Quantification with confocal quantitative image cytometer

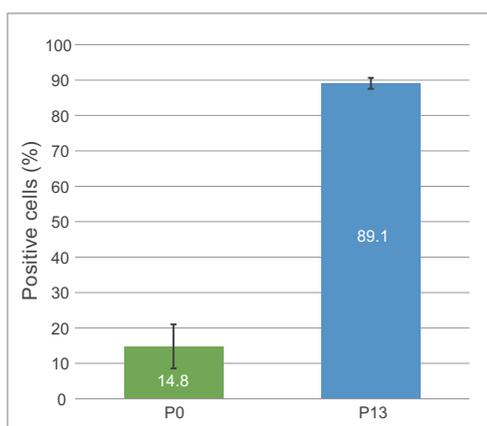
In the conventional method of X-gal, SA-β-gal-positive cells are counted under microscope and calculate the percent of the senescent cells by compared with total cells. The SA-β-gal-positive cells were stained with this kit and analyzed using confocal quantitative image cytometer CQ1(Yokogawa Electric Corporation).



### Imaging condition

Culture plate : 96 well plate  
 Objective lens : 10 times  
 Emission Detector  
 405 nm (Hoechst 33342) : cyan  
 488 nm (SPiDER-βGal) : green

\* SA-β-gal positive cells are marked with red circles.



### Quantification for SA-β-gal-positive cells

The total cells were stained with DNA staining dye, Hoechst 33342. The SA-β-gal-positive cells and total were counted and calculate the ratio of senescent cells.

The difference of SA-β-gal-positive cells ratio were shown in WI-38 cells depending on the number of passage. The data was quickly analysed with the confocal quantitative image cytometer compared with the manually counting procedure with X-gal staining method.

Discription	Unit	Code
Cellular Senescence Detection Kit - SPiDER-βGal	-	SG03 or SG04

# Cell Proliferation / Cytotoxicity Assay Kit

## Cell Counting Kit - 8

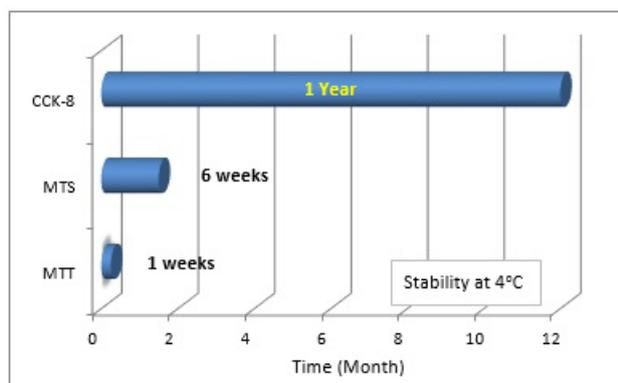
Cell Counting Kit-8 (CCK-8) allows sensitive colorimetric assays for the determination of cell viability in cell proliferation and cytotoxicity assays. Dojindo's highly water-soluble tetrazolium salt, WST-8, is reduced by dehydrogenase activities in cells to give a yellow-color formazan dye, which is soluble in the tissue culture media. The amount of the formazan dye, generated by the activities of dehydrogenases in cells, is directly proportional to the number of living cells. The detection sensitivity of CCK-8 is higher than the other tetrazolium salts such as MTT, XTT, MTS or WST-1.

### Simple procedure

Cell Counting Kit-8 requires 3 simple steps.



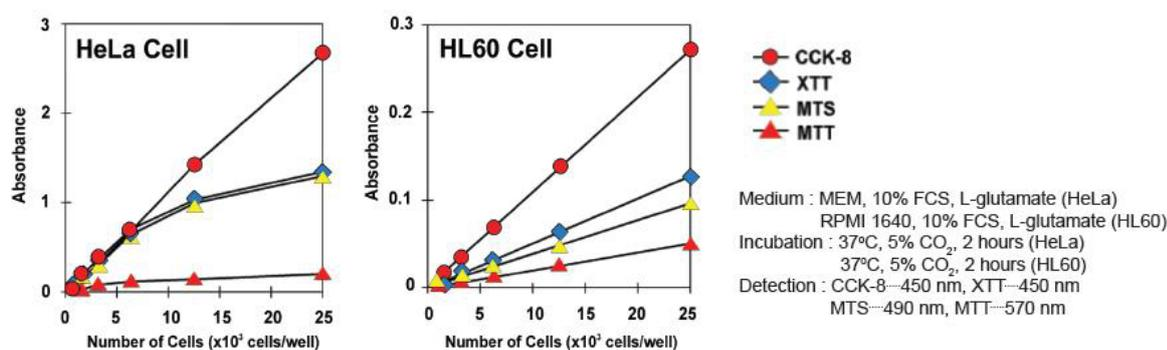
### Longer stability



Cell Counting Kit-8 is a ready-to-use solution. It is also stable at 4°C for 1 year. Your assay can be done anytime without thaw and freeze.

### Higher sensitivity

CCK-8 (WST-8) is the highest sensitive kit for cell based assay.



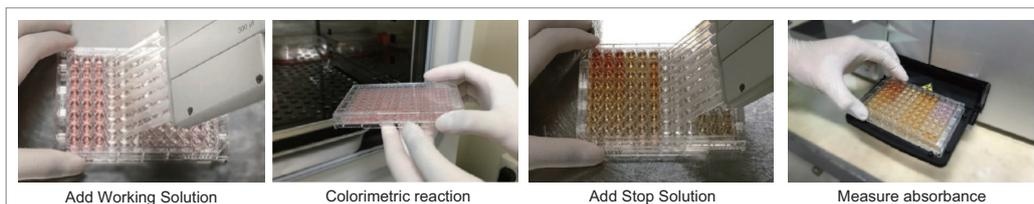
Discription	Unit	Code
Cell Counting Kit-8	100 - 10,000 tests	CK04

# Cytotoxicity Assay Kit

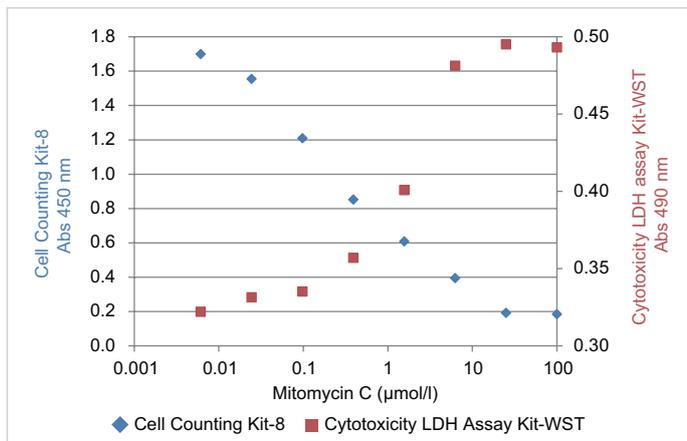
## Cytotoxicity LDH Assay Kit-WST

Lactate dehydrogenase(LDH) is an enzyme that presents in almost cell types and it catalyzes the oxidation of lactate to pyruvate in the presence of co-enzyme NAD+. Once Cells are impaired by stress, injuries, chemicals, or intercellular signals, LDH is rapidly released from the cell membrane. Thus, the measurement of the amount of released LDH from cells is one of the major methods to assess the cell death. Cytotoxicity LDH Assay Kit-WST is designed for use with 96-well micro plates formats for high-throughput screening.

### Simple procedure



### Combination use



To obtain the accurate result in cytotoxicity assay, the samples are measured by different principles. CCK-8 and LDH assay measure NADH as living cells and released LDH as dead cells, respectively. According to these results, the living cells were reduced and dead cells were increased with the high concentration of toxicant.

Discription	Unit	Code
Cytotoxicity LDH Assay Kit-WST	100 tests	CK12
	500 tests	
	2,500 tests	

For the detail information of each product, put the product code with "dojindo" and visit our web site.



If you have any questions or need any assistances, please feel free to contact us.

Japan (Headquarters)  
Dojindo Laboratories  
Tel: 096-286-1515  
E-mail: info@dojindo.co.jp  
URL: www.dojindo.co.jp

USA/Canada  
Dojindo Molecular Technologies, Inc.  
Tel: +1-301-987-2667  
E-mail: info@dojindo.com  
URL: www.dojindo.com

Europe  
Dojindo EU GmbH  
Tel: +49-89-3540-4805  
E-mail: info@dojindo.eu.com  
URL: www.dojindo.eu.com

China  
Dojindo China Co., Ltd  
Tel: +86-21-6427-2302  
E-mail: info@dojindo.cn  
URL: www.dojindo.cn