

Reagent for Cellular Function Analysis

Autophagy

- Autophagic Flux Assay Kit
- DALGreen-Autophagy Detection
- DAPGreen-Autophagy Detection
- DAPRed-Autophagy Detection

Senescence

- Cellular Senescence Detection Kit
-SPiDER-βGal
- Cellular Senescence Plate Assay Kit
-SPiDER-βGal

Neurodegenerative Diseases

Cancer

Senescence

Mitochondria

- Mitophagy Detection Kit
- JC-1 MitoMP Detection Kit
- MitoBright LT Series
- MT-1 MitoMP Detection Kit
- MitoBright ROS Deep Red
- Extracellular OCR Plate Assay Kit

Cellular Metabolism

- Glycolysis/OXPHOS Assay Kit
- ATP Assay Kit-Luminescence
- Lactate Assay Kit-WST

Ferroptosis

- FerroOrange
- Liperfluo
- Mito-FerroGreen
- MitoPeDPP
- Cystine Uptake Assay Kit
- MDA Assay Kit
- Lipid Peroxidation Probe
-BDP 581/591 C11-

Senescence

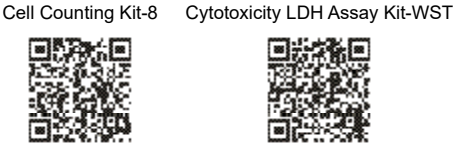
| Code | Product Name | Page |
|------------------------|--|------|
| SG03 | Cellular Senescence Detection Kit - SPiDER-βGal | 8 |
| SG05 | Cellular Senescence Plate Assay Kit - SPiDER-βGal | 9 |
| MT09 | JC-1 MitoMP Detection Kit | 27 |
| MT10, MT11, MT12 | MitoBright LT Series | 31 |
| MT13 | MT-1 MitoMP Detection Kit | 27 |
| MT16 | MitoBright ROS Deep Red - Mitochondrial Superoxide Detection | 29 |
| MD01 | Mitophagy Detection Kit | 28 |
| MT02 | Mtphagy Dye | 28 |
| E297 | Extracellular OCR Plate Assay Kit | 26 |
| A562 | Autophagic Flux Assay Kit | 11 |
| D675 | DALGreen - Autophagy Detection | 10 |
| D676 | DAPGreen - Autophagy Detection | 10 |
| D677 | DAPRed - Autophagy Detection | 10 |
| F374 | FerroOrange | 16 |
| L248 | Liperfluo | 14 |
| M489 | Mito-FerroGreen | 30 |
| UP05 | Cystine Uptake Assay Kit | 23 |
| L267 | Lipid Peroxidation Probe -BDP 581/591 C11- | 15 |
| G270 | Glycolysis/OXPHOS Assay Kit | 21 |
| A550 | ATP Assay Kit-Luminescence | 19 |
| L256 | Lactate Assay Kit-WST | 19 |
| R252 | ROS Assay Kit -Highly Sensitive DCFH-DA- | 12 |
| R253 | ROS Assay Kit -Photo-oxidation Resistant DCFH-DA- | 13 |
| G257 | GSSG/GSH Quantification Kit | 17 |
| LD01, LD02, LD03, LD04 | Lipi-Blue, Lipi-Green, Lipi-Red, Lipi-Deep Red | 40 |
| LD05 | Lipid Droplet Assay Kit - Blue | 41 |
| LD06 | Lipid Droplet Assay Kit - Deep Red | 41 |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

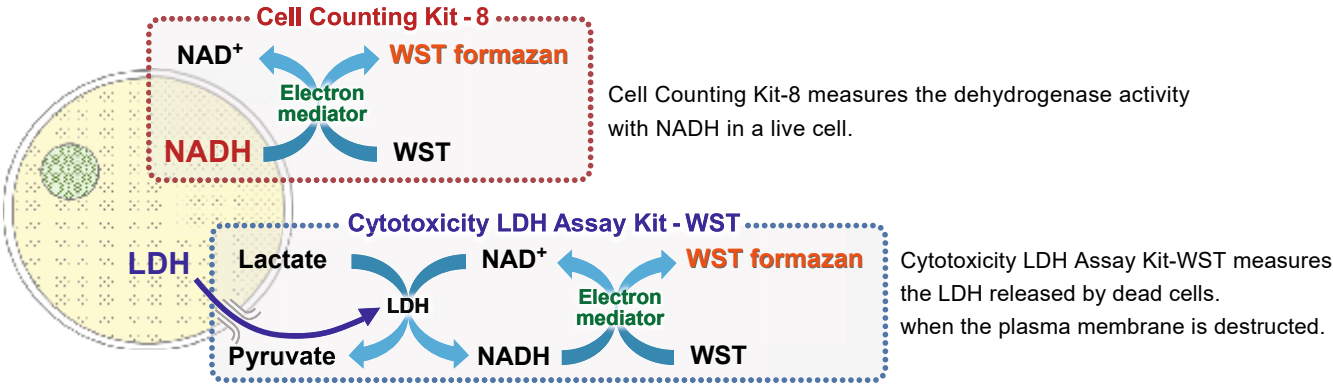
Cell Proliferation / Cytotoxicity Assay

Cell Counting Kit-8

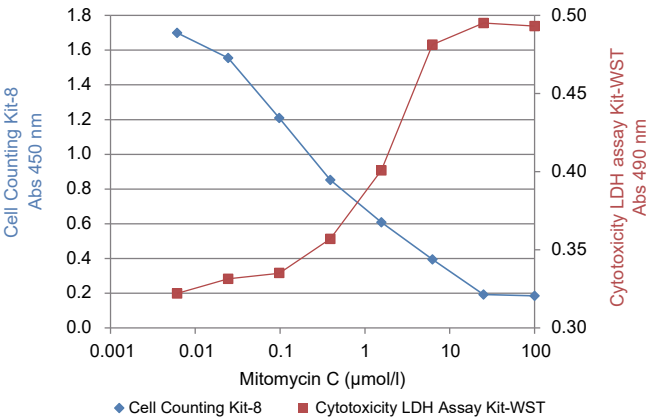
Cytotoxicity LDH Assay Kit-WST



Detection Principle



Simultaneous Usage of CCK-8 and Cytotoxicity LDH Assay Kit-WST



Drug: Mitomycin C

Cell Line: HeLa

Media: MEM, 10% FBS

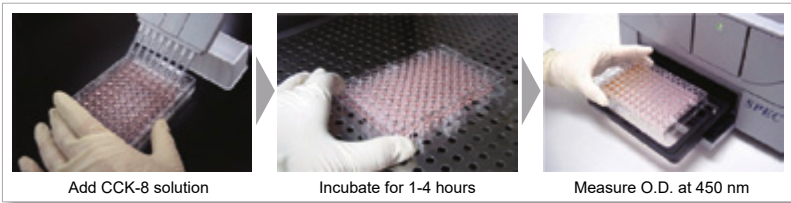
Incubation: 37°C, 5% CO₂ for 48 hours

Measuring Condition: Cell Counting Kit-8 (450 nm)

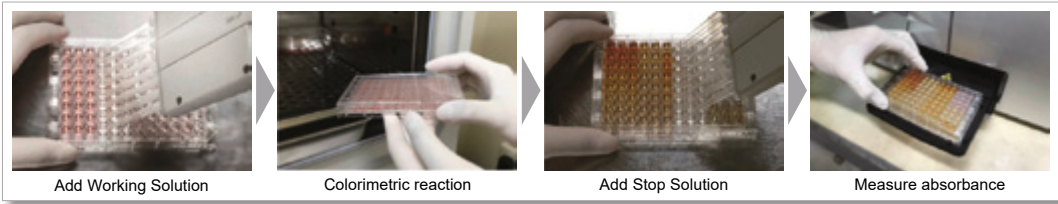
Cytotoxicity LDH Assay Kit-WST (490 nm)

Simple Procedure

• Cell Counting Kit-8

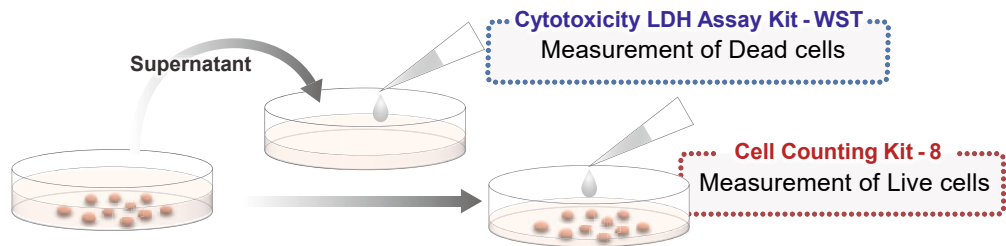


• Cytotoxicity LDH Assay Kit-WST



Same Samples can be used

Since same samples can be used for Cell Counting Kit-8 and Cytotoxicity LDH Assay Kit-WST, the method is convenient and time efficient.



| Description | Unit | Code |
|--------------------------------|-------------|---------|
| Cell Counting Kit-8 | 500 tests | CK04-05 |
| | 1000 tests | CK04-11 |
| | 3000 tests | CK04-13 |
| | 10000 tests | CK04-20 |
| Cytotoxicity LDH Assay Kit-WST | 100 tests | CK12-01 |
| | 500 tests | CK12-05 |
| | 2000 tests | CK12-20 |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Senescence Detection

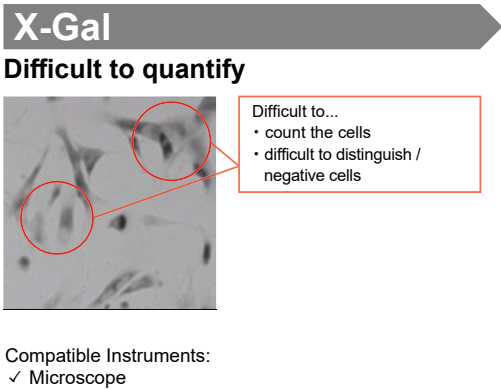
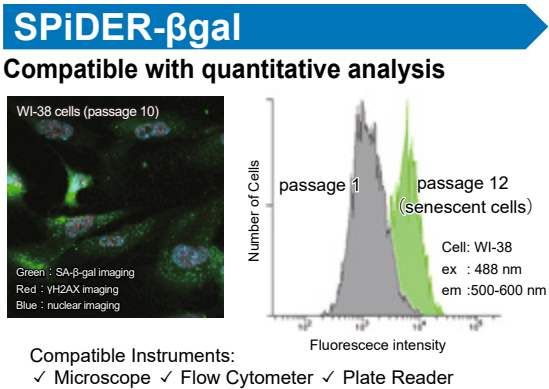
Cellular Senescence Detection Kit - SPiDER-βGal

Cellular Senescence Detection Kit - SPiDER Blue

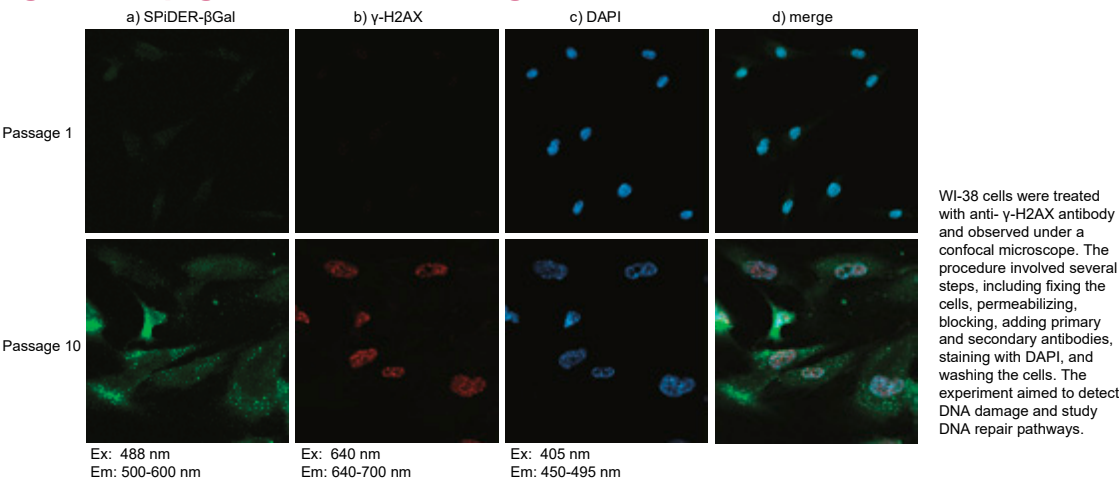
NEW



Cellular Senescence Detection Kit – SPiDER-βGal allows to detect SA-β-gal with high sensitivity and ease of use. SPiDER-βGal is a new reagent to detect β-galactosidase which possesses a high cell-permeability and a high retentivity inside cells. SA-β-gal are detected specifically not only in living cells but also fixed cells by using a reagent (Bafilomycin A1) to inhibit endogenous β-galactosidase activity. Therefore, SPiDER-βGal can be applied to quantitative analysis by flow cytometry.



Co-staining of SA- β-gal and DNA Damage marker in WI-38 cells



| Description | Unit | Code |
|---|-----------|---------|
| Cellular Senescence Detection Kit - SPiDER-βGal | 10 assays | SG03-10 |
| Cellular Senescence Detection Kit - SPiDER Blue | 1 Plate | SG07-10 |

Cellular Senescence Plate Assay Kit - SPiDER-βGal



This product is a simple detection kit by plate assay for senescence-associated β-galactosidase (SA-β-gal) activity which is used as a marker for senescent cells. By simply adding SPiDER-βGal, a reagent for detection of β-galactosidase, to 96 well plates, this kit allows you to quantify SA-β-gal activity and makes it possible to evaluate multiple samples. When normalization is done by the results obtained by counting cells, quantifying nucleic acids (a relevant product), or quantifying proteins, the measured values obtained using this kit become available for evaluating SA-β-gal activity according to cell number.

Correlation with Imaging Data

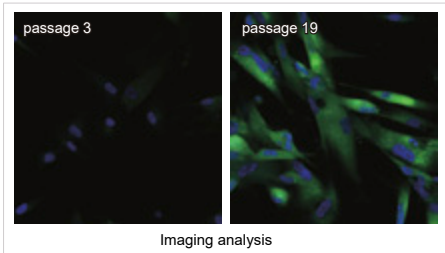
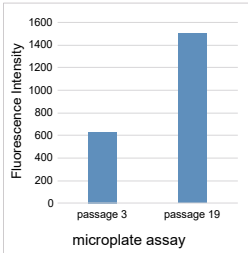


Plate Assay
Ex. 535nm / Em. 580nm

Imaging data
Green: Ex. 488nm / Em. 500-600nm (SA-β-Gal staining with Cellular Senescence Detection Kit – SPiDER-βGal(Code SG04))
Blue: Ex. 405nm / Em. 450-495nm (Nuclear staining with -Cellstain- DAPI solution(Code D523))

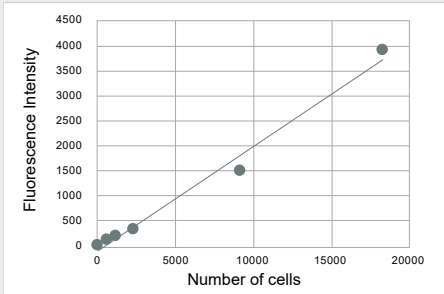
As a result, it was confirmed that in both kits, SA-β-gal staining increased in the high-passage WI-38 cells. Bear in mind that although initial cell seeding densities are the same, cell densities at the time of plate assay differ due to low proliferation rate of senescent cells at higher passage levels. Therefore, in this experiment, we used SA-β-Gal activity values normalized by the results obtained using the Cell Count Normalization Kit in which cell number is determined by a nuclear marker.

Cell Count Normalization Kit

Cell Count Normalization Kit includes nucleic acid staining dye, Hoechst 33342 which binds to nuclear DNA to emit blue fluorescence. By measuring this blue fluorescence, correction of the measured value can easily be carried out in simple steps whereas the visual cell counting method requires complicated procedure. Moreover, unlike the correction by protein or ATP amount, the kit requires no lysis procedure. In addition, Quenching Buffer included in the kit enables a direct measuring of fluorescence signal without any background.



Highly correlated to cell number



| Description | Unit | Code |
|---|------------|---------|
| Cellular Senescence Plate Assay Kit - SPiDER-βGal | 20 tests | SG05-01 |
| | 100 tests | SG05-05 |
| Cell Count Normalization Kit | 200 tests | C544-02 |
| | 1000 tests | C544-10 |

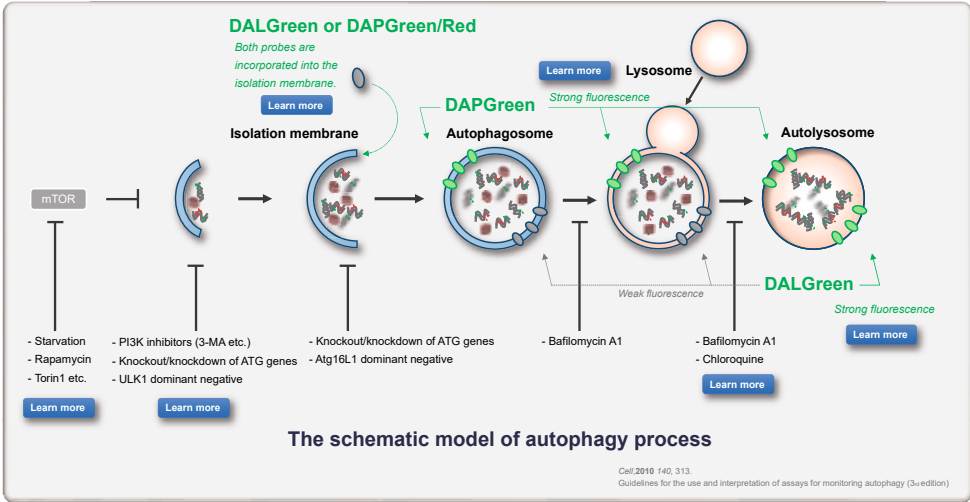
Autophagy

DAPGreen / Red - Autophagy Detection

DALGreen - Autophagy Detection



DAPGreen and DAPRed detect autophagosomes, while DALGreen detects autolysosomes. These dyes are permeable to cells and enables live cell imaging with fluorescence microscopy, and DAPGreen and DALGreen allow for quantitative assay by flow cytometry. Autophagy is an intracellular degradation system involving autophagosome formation, detected by DAPGreen and DAPRed, and lysosome fusion, detected by DALGreen, which fluoresces intensity increases in acidic conditions.



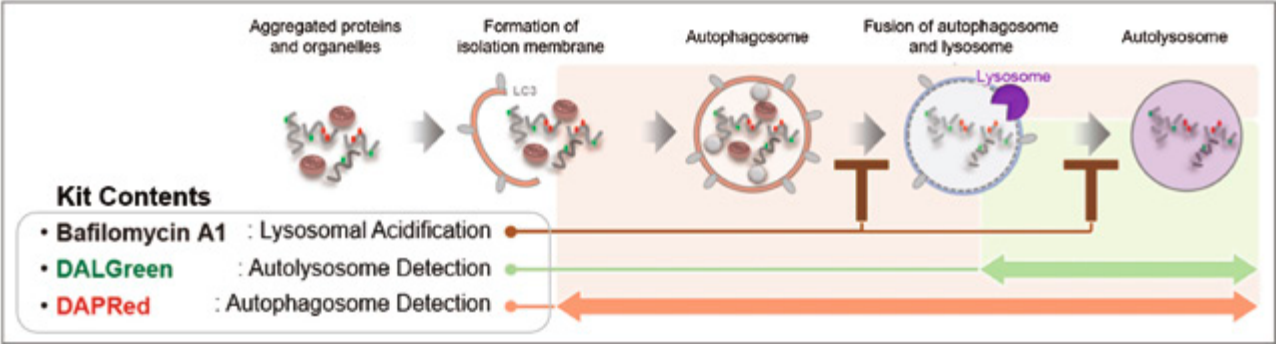
Feature of Each Dye

| | Applicable instruments | | | Fluorescent properties | Volume / the number of usable assays | Existing methods |
|----------|------------------------|----------------|-------------------|---|---|--------------------------------|
| | Fluorescent Microscope | Flow cytometer | Microplate reader | | | |
| DAPGreen | ✓ | ✓ | ✓ | Ex = 425-475 nm Em = 500-560 nm <small>* For confocal microscope, the sample can be excited at 488 nm</small> | 5 nmol x 1 / 35 mm dish: 25 (when used in 1.0 μmol/l) | LC3-GFP MDC Cyto-ID etc. |
| DAPRed | ✓ | | | Ex = 500-560 nm Em = 690-750 nm | 5 nmol x 1 / 35 mm dish: 25 (when used in 1.0 μmol/l) | |
| DALGreen | ✓ | ✓ | | Ex = 350-450 nm Em = 500-560 nm <small>* For confocal microscope, the sample can be excited at 488 nm</small> | 20 nmol x 1 / 35 mm dish: 10 (when used in 1.0 μmol/l) | LC3-GFP-RFP etc. |

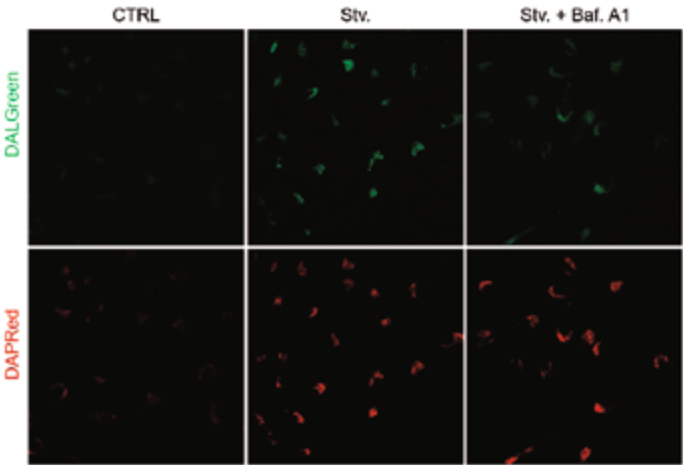
*Double staining imaging by DAPGreen and DALGreen is not possible

Autophagy

Autophagic Flux Assay Kit

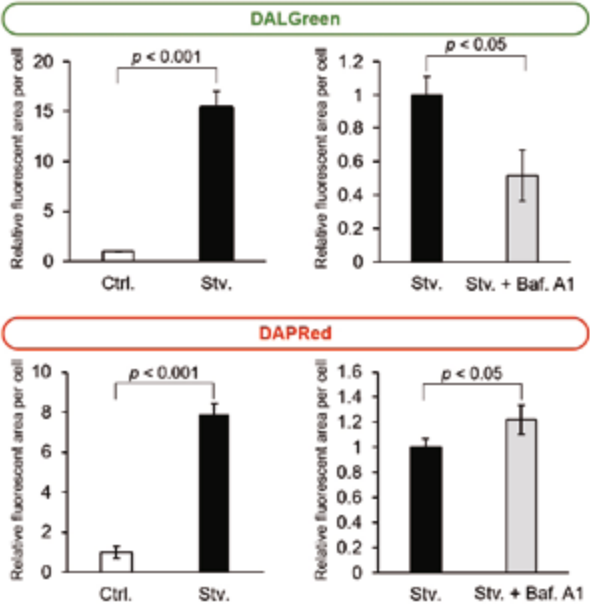


Experimental Example: Autophagy Flux Analysis



By culturing HeLa cells in HBSS with starvation, autophagy was induced and DAPRed and DALGreen fluorescence increased. Addition of Baf. A1 decreased DALGreen fluorescence, indicating that autolysosomes were reduced and Autophagy Flux was inhibited.

Quantification method: Fluorescence values (area) were obtained in Image J and normalized by the number of cells per field of view*. Number of samples: n=3
*Please obtain images with the same number of cells per field of view as possible.



| Description | Unit | Code |
|--------------------------------|---------|---------|
| Autophagic Flux Assay Kit | 1 set* | A562-10 |
| DALGreen - Autophagy Detection | 20 nmol | D675-10 |
| DAPGreen - Autophagy Detection | 5 nmol | D676-10 |
| DAPRed - Autophagy Detection | 5 nmol | D677-10 |

*Equivalent to 5 dishes (35 mm dish)

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Oxidative Stress

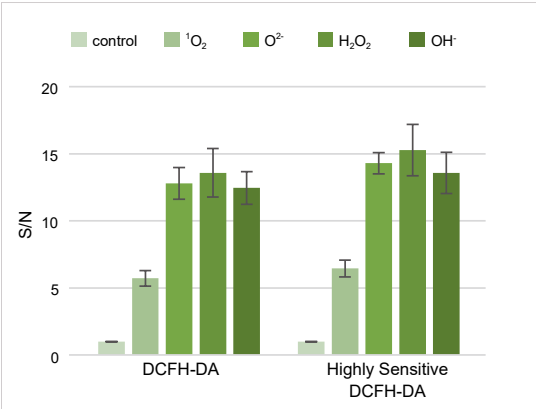
ROS Assay Kit -Highly Sensitive DCFH-DA-



ROS Assay Kit -Highly Sensitive DCFH-DA- overcomes these limitations. The dye allows ROS detection with higher sensitivity than DCFH-DA. Moreover, the Loading Buffer included in this kit maintains cellular health during assays.

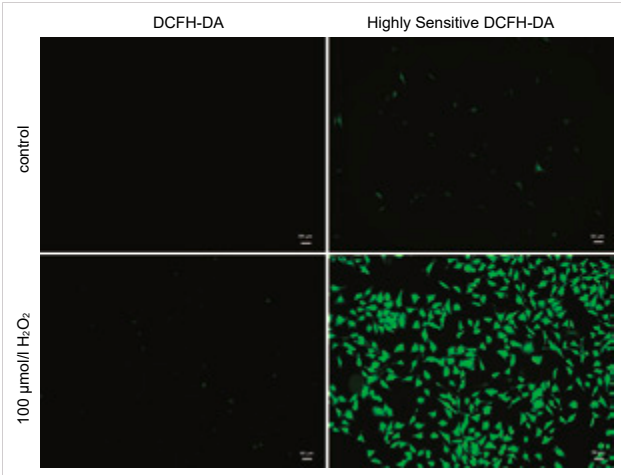
The reactivity of the Highly Sensitive DCFH-DA for ROS is similar to the reactivity of 2'-7' dichlorofluorescein diacetate (DCFH-DA). The Highly Sensitive DCFH-DA also has similar fluorescence characteristics (λ_{ex} : 505 nm, λ_{em} : 525 nm) to DCFH-DA. Therefore, ROS is detectable at the same excitation/fluorescence wavelength.

The selectivity for ROS

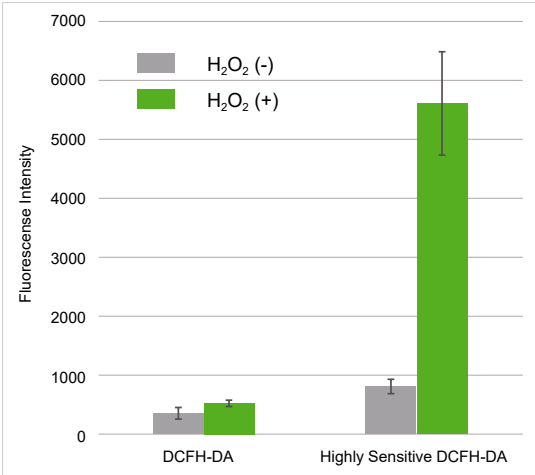


High Sensitive Detection Compared with DCFH-DA

Detection using fluorescent microscope



Detection using microplate reader



Hydrogen peroxide (H_2O_2)-treated HeLa cells (1×10^4 cells/ml) were stained with DCFH-DA or the ROS Assay Kit-Highly Sensitive DCFH-DA, and the fluorescence intensity of intracellular ROS was compared between two detection kits. As a result, the ROS Assay Kit-Highly Sensitive DCFH-DA in high-sensitivity detection of intracellular ROS was better than DCFH-DA.

| Description | Unit | Code |
|--|-----------|---------|
| ROS Assay Kit -Highly Sensitive DCFH-DA- | 100 tests | R252-10 |

Oxidative Stress

ROS Assay Kit -Photo-oxidation Resistant DCFH-DA-



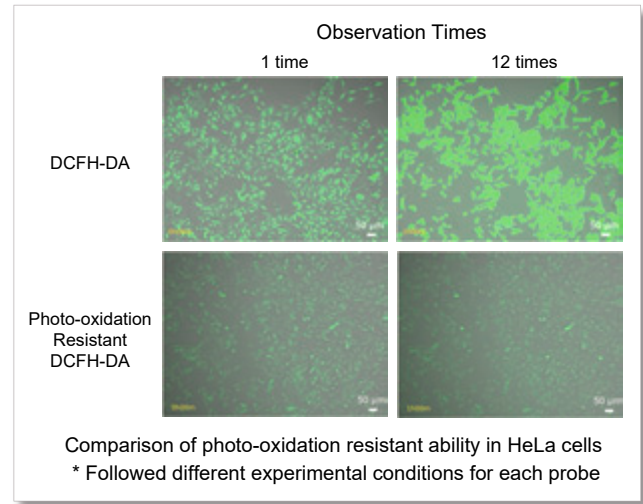
The dye that is employed in this kit allows ROS detection with higher sensitivity than DCFH-DA; It does not leak from cells because the fluorescent dye can immobilize protein via a chemical bond, and it is resistant to photo-oxidation compared with DCFH-DA. Moreover, the Loading Buffer in the kit maintains cellular health during assays.

Time-lapse imaging movie Available

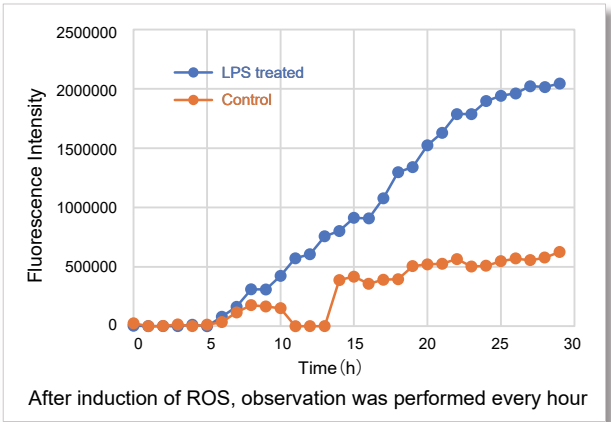
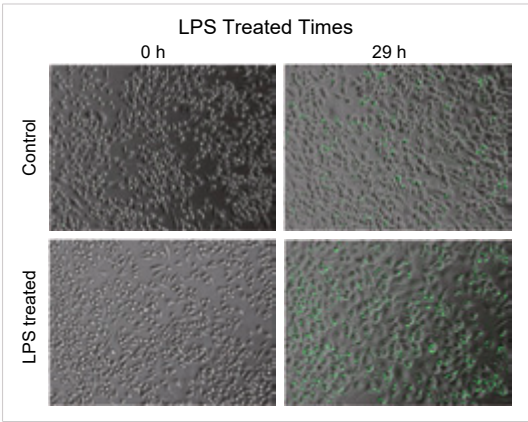
Other Probe

Photo-oxidation Resistant DCFH-DA-

Resistant to Photo-oxidation



Simultaneous Detection of ROS in LPS-treated macrophages



In Lipopolysaccharide (LPS) treated RAW 264.7 cells, after being stained with regular DCFH-DA, Highly Sensitive DCFH-DA, or Photo-oxidation Resistant DCFH-DA, the intracellular ROS level was compared. The results showed that the Dojindo Laboratories' probes could detect intracellular ROS with higher sensitivity.

| Description | Unit | Code |
|---|-----------|---------|
| ROS Assay Kit -Photo-oxidation Resistant DCFH-DA- | 100 tests | R253-10 |

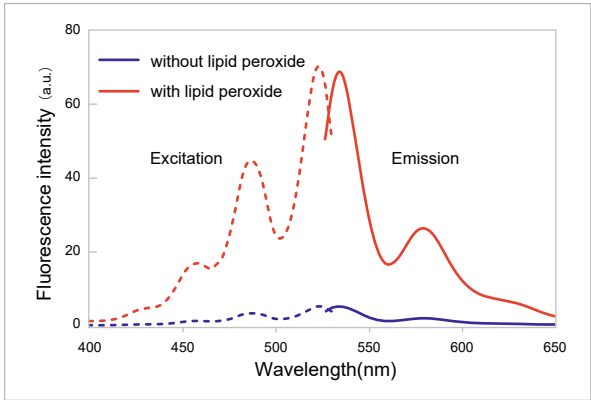
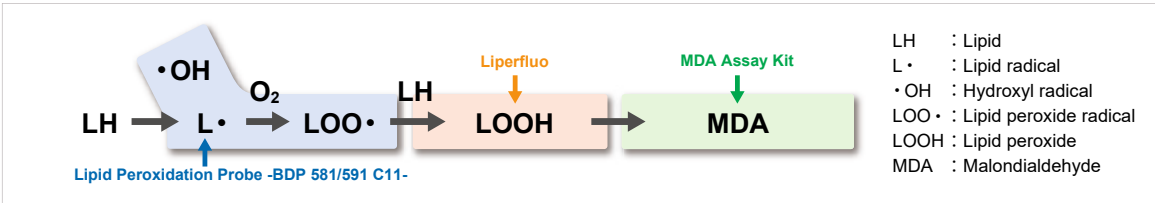
| |
|--|
| Proliferation |
| Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Lipid Peroxide Detection

Liperfluo

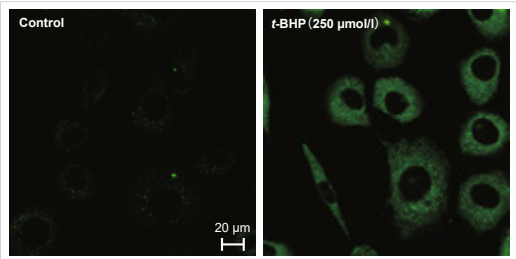


Liperfluo is a Dojindo-developed fluorescence probe to specifically detect lipid peroxides with minimal photo-damage or auto-fluorescence. It emits intense fluorescence in organic solvents and is nearly non-fluorescent in aqueous media. Liperfluo's tetraethyleneglycol group increases its solubility and makes it suitable for imaging lipid peroxides in cell membranes. It's used to monitor lipid peroxidation in ferroptosis research through fluorescence microscopy and flow cytometry.



Excitation and emission without lipid peroxide spectra of Liperfluo with or without lipid peroxide in ethanol.

Lipid Peroxide Detection in Living Cells



Liperfluo added to cells, t-BHP induced lipid peroxidation and cells were observed under confocal microscope to study ferroptosis.

Cell line: L929
 Microscope: Zeiss LSM510META
 Filter type: FITC (GFP, Alexa488) wide filter
 HFT UV/488
 NFT490
 BP505-550

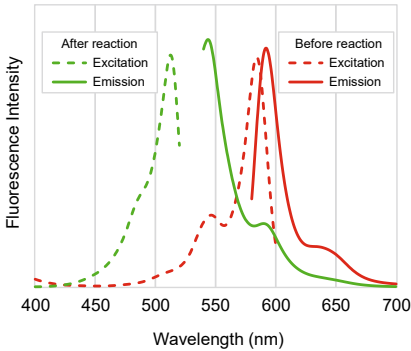
| | Description | Unit | Code |
|-----------|-------------|-------------------|---------|
| Liperfluo | | 1 set (50 μg × 5) | L248-10 |

Lipid Peroxidation Detection

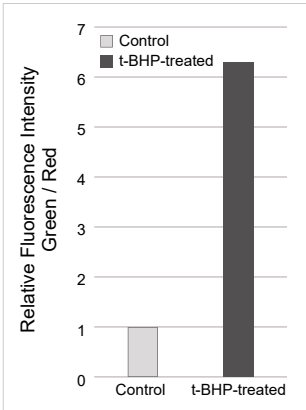
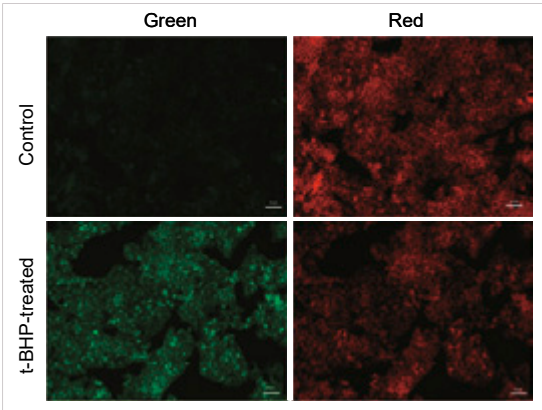
Lipid Peroxidation Probe -BDP 581/591 C11-



Lipid Peroxidation Probe -BDP 581/591 C11- is a fluorescent probe for detecting lipid peroxidation. This fluorescent probe does not react with lipid peroxides but reacts with lipid radicals generated when lipids are peroxidized, resulting in the detection of lipid peroxidation. The unreacted probe emits red fluorescence, but after reacting with radicals around lipids, it changes its fluorescence from red to green. Thus, lipid peroxidation can be detected with high sensitivity because it is detected by the ratio of red to green fluorescence intensity.



Lipid Peroxidation Assay



HepG2 cells stained with this probe were stimulated with HBSS solution containing 200 $\mu\text{mol/l}$ *t*-BHP for 2 hours, and the fluorescence intensity was compared with control cells. As a result, a decrease in red fluorescence and an increase in green fluorescence were observed with high sensitivity in *t*-BHP-treated cells compared to untreated cells. The cells were detected using a plate reader, and the values obtained were calculated as the intensity ratio of green/red fluorescence, which allowed quantified lipid peroxidation. Furthermore, an increase in the histogram of green fluorescence was observed when the cells were detected using a flow cytometer. Which improves that this dye is three different instruments.

<Experimental Conditions>
Fluorescent Microscope
Green: GFP filter (Ex = 450-490 nm, Em = 500-550 nm)
Red: TexasRed filter (Ex = 540-580 nm, Em = 600-660 nm)
Scale bar: 50 μm

Fluorescent Plate Reader
Green: Ex = 490 nm, Em = 520-540 nm
Red: Ex = 570 nm, Em = 600-620 nm

| Description | Unit | Code |
|--|-----------|---------|
| Lipid Peroxidation Probe -BDP 581/591 C11- | 200 tests | L267-10 |

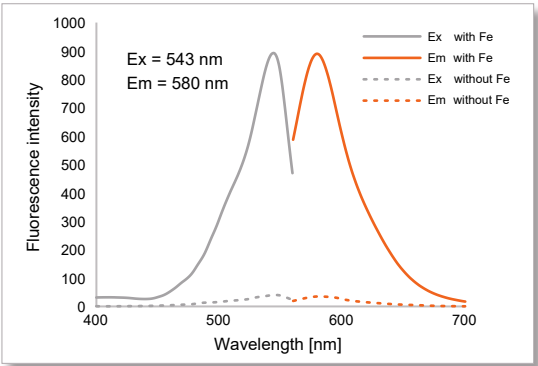
| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Intracellular Iron Ion Measurement

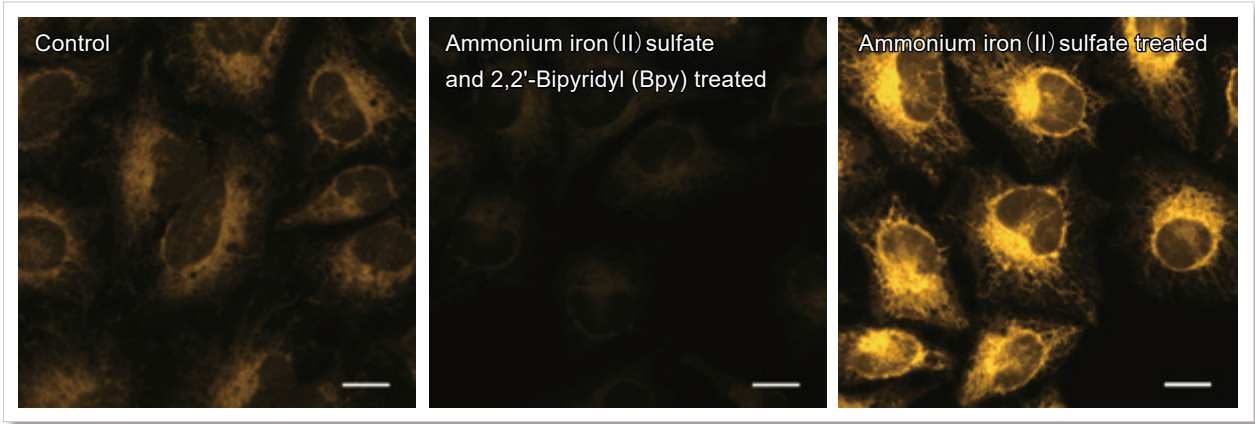
FerroOrange



Liperfluo is a Dojindo-developed fluorescence probe to specifically detect lipid peroxides with minimal photo-damage or auto-fluorescence. It emits intense fluorescence in organic solvents and is nearly non-fluorescent in aqueous media. Liperfluo's tetraethyleneglycol group increases its solubility and makes it suitable for imaging lipid peroxides in cell membranes. It's used to monitor lipid peroxidation in ferroptosis research through fluorescence microscopy and flow cytometry.



Experimental Example



HeLa cells treated with chelator of iron 2,2'-bipyridyl (Bpy) (100 μmol/l) or Ammonium iron (II) sulfate (100 μmol/l) were prepared. The change of intracellular Fe²⁺ in HeLa cells was detected by the FerroOrange.
Ex = 561 nm, Em = 570-620 nm, Scale bars 20 μm

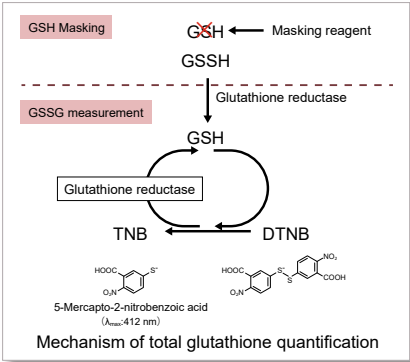
| Description | Unit | Code |
|-------------|--------|---------|
| FerroOrange | 1 tube | F374-10 |
| | 3 tube | F374-12 |

Quantification of Reduced (GSH) and Oxidized (GSSG) Glutathione

GSSG/GSH Quantification Kit



The GSSG/GSH Quantification kit contains Masking Reagent of GSH. GSH will be deactivated in the sample by simply adding the Masking Reagent. Then, using the enzymatic recycling system, only the GSSG will be detected by measuring the absorbance ($\lambda_{\text{max}} = 412 \text{ nm}$) of DTNB (5,5-dithio-bis- (2-nitrobenzoic acid)). The quantity of GSH can also be determined, by subtracting GSSG from the total amount of glutathione. With this kit, GSH/ GSSG concentrations from $0.5 \mu\text{mol/l}$ to $50 \mu\text{mol/l}$ and GSSG concentrations from $0.5 \mu\text{mol/l}$ to $25 \mu\text{mol/l}$ can be quantified.



Assay Procedure

1) GSSG/GSH Standard Solution and add Sample A or Sample B to each well.

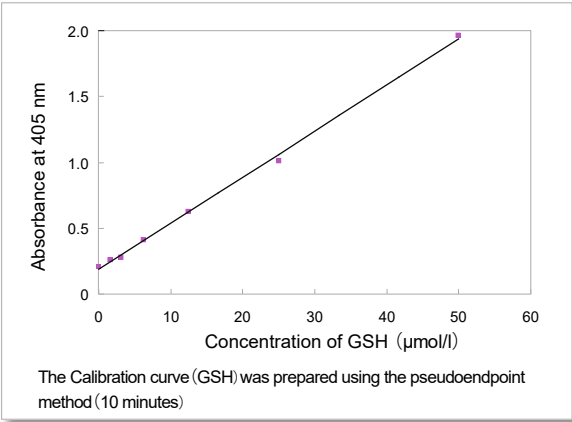
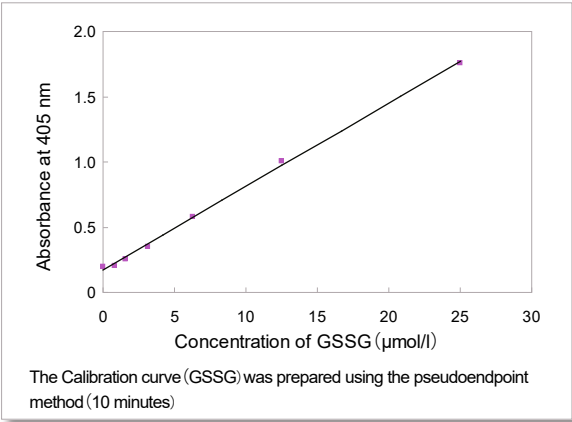
2) Add Buffer solution to each well

3) Incubate at 37°C for 1 h.

4) -5) Add substrate working solution and Enzyme/ Coenzyme working solution to each well.

6)-7) After incubating at 37°C for 10 minutes, measure the absorbance of each well with a microplate.

Calibration Curve



| Description | Unit | Code |
|-----------------------------|-----------|---------|
| GSSG/GSH Quantification Kit | 200 tests | G257-10 |

| |
|------------------------------|
| Proliferation |
| Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles |
| Exosome, Lipid Droplet, etc. |

Proliferation
Cytotoxicity

Senescence

Autophagy

Oxidative
Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Other Organelles
Exosome, Lipid Droplet, etc.

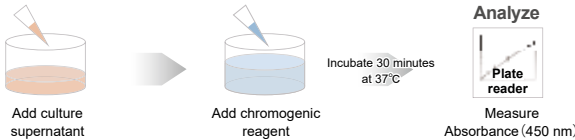
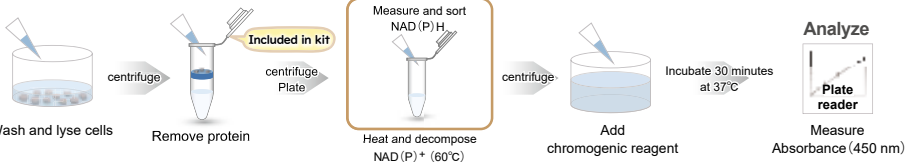
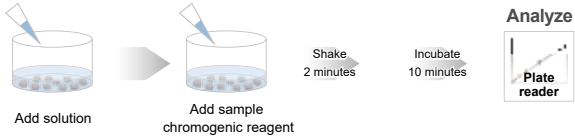
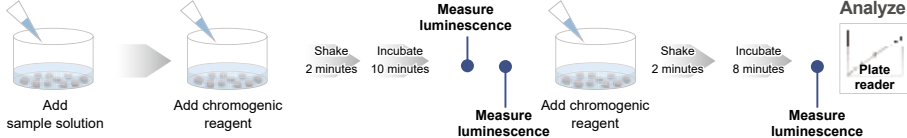
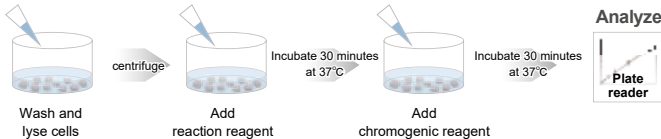
Measurements of Intracellular Metabolism



| Description | Unit | Code |
|---|-----------|---------|
| Starter Kit | | |
| Glycolysis/OXPHOS Assay Kit | 50 tests | G270-10 |
| Glycolysis/JC-1 MitoMP Assay Kit | 50 tests | G272-10 |
| Quantification for Intracellular Metabolism | | |
| ATP Assay Kit-Luminescence | 50 tests | A550-10 |
| | 200 tests | A550-12 |
| ADP/ATP Ratio Assay Kit-Luminescence | 100 tests | A552-10 |
| Glucose Assay Kit-WST | 50 tests | G264-05 |
| | 200 tests | G264-20 |
| Glutamine Assay Kit-WST | 100 tests | G268-10 |
| Glutamate Assay Kit-WST | 100 tests | G269-10 |
| α-Ketoglutarate Assay Kit-Fluorometric | 100 tests | K261-10 |
| | 50 tests | L256-10 |
| Lactate Assay Kit-WST | 200 tests | L256-20 |
| | 100 tests | N509-10 |
| NADP/NADPH Assay Kit-WST | 100 tests | N510-10 |
| Uptake Assay Kit | | |
| Glucose Uptake Assay Kit-Blue | 1 set | UP01-10 |
| Glucose Uptake Assay Kit-Green | 1 set | UP02-10 |
| Glucose Uptake Assay Kit-Red | 1 set | UP03-10 |
| Amino Acid Uptake Assay | 20 tests | UP04-10 |
| | 100 tests | UP04-12 |
| Cystine Uptake Assay Kit | 20 tests | UP05-10 |
| | 100 tests | UP05-12 |
| Fatty Acid Uptake Assay Kit | 100 tests | UP07-10 |

Simple Procedure for First Time User

For a first-time user, the kit includes the reagents and components necessary for measuring samples. You'll soon realize how easy it is to use.

| Determination index | Detection | Operation |
|---|--------------|---|
| <div>Glucose</div> <div>Lactate</div> <div>Glutamine</div> <div>Glutamate</div> <div>NAD/NADH</div> <div>NADP/NADPH</div> | Colorimetric | <p>Simply transfer the culture supernatant to a plate and mix it with the chromogenic reagent</p> <div></div> |
| | | <div></div> |
| <div>ATP</div> <div>ADP/ATP</div> | Luminescent | <p>Kit includes ATP standard - very easy to use</p> <div></div> |
| | | <div></div> |
| <div>α-ketoglutaric acid</div> | Fluorescent | <p>Less variable results than existing assays</p> <div></div> |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

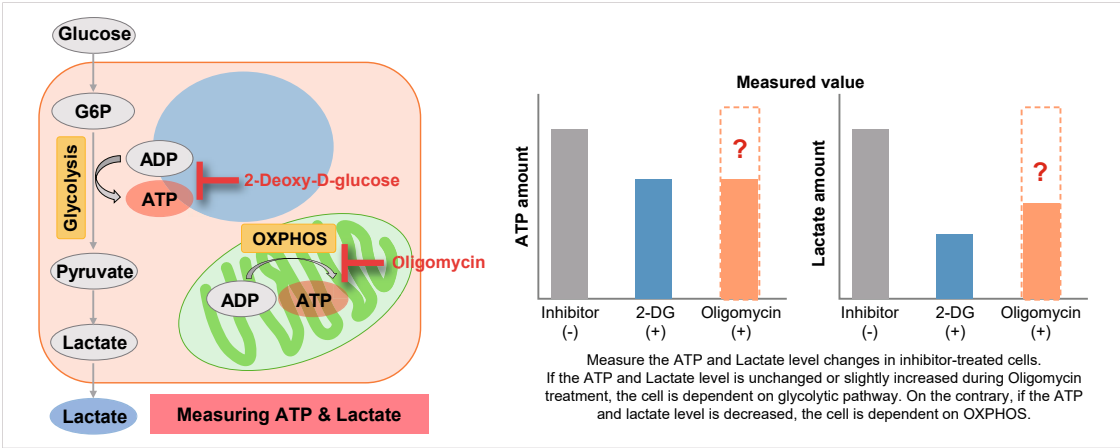
Intracellular Metabolism

Glycolysis/OXPHOS Assay Kit

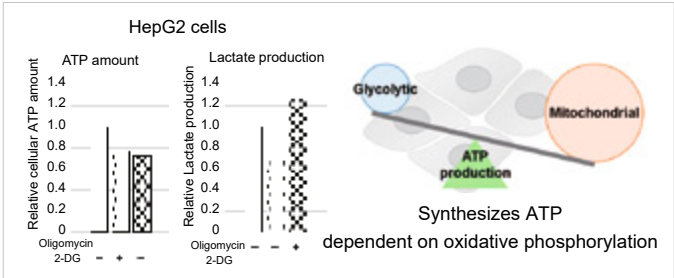
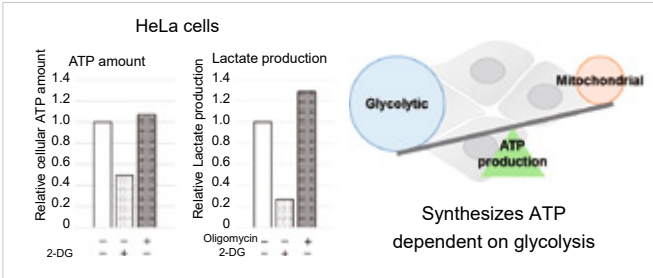


- Easy test via plate reader, no need for expensive equipment
- All reagent acquired is included, ready to use kit
- Easy-to-understand detailed protocol

Combining methods (1) and (2) can be used to measure the metabolic pathway dependency of cells. Cells are treated with oligomycin or 2-DG to inhibit OXPHOS or ATP synthesis in the glycolytic pathway, and the amounts of ATP and lactate production are measured, respectively. Changes in the amount of ATP can be used to determine the efficiency of energy production, and changes in the amount of lactate produced can be used to determine changes in glycolytic capacity and evaluate whether cells are dependent on glycolysis or OXPHOS.



Experimental Example: Comparison of metabolic pathway dependence in different cell line



| Description | Unit | Code |
|-----------------------------|----------|---------|
| Glycolysis/OXPHOS Assay Kit | 50 tests | G270-10 |

Proliferation
Cytotoxicity

Senescence

Autophagy

Oxidative
Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Other Organelles
Exosome, Lipid Droplet, etc.

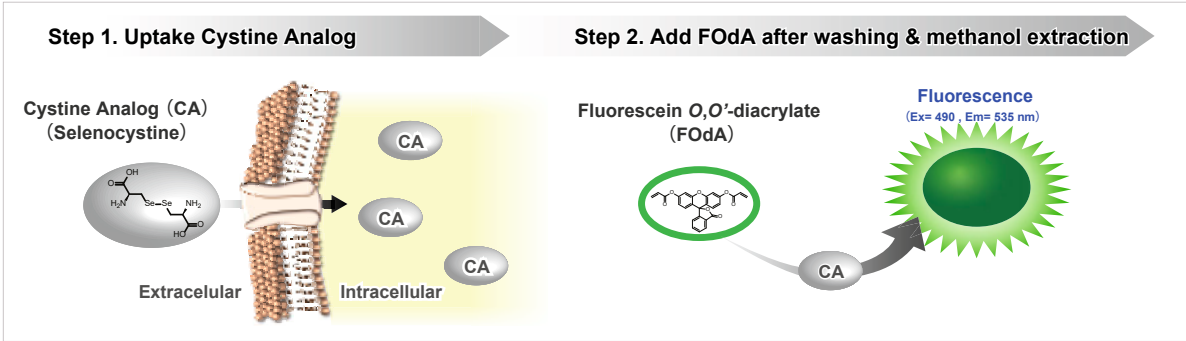
Intracellular Metabolism

Cystine Uptake Assay Kit



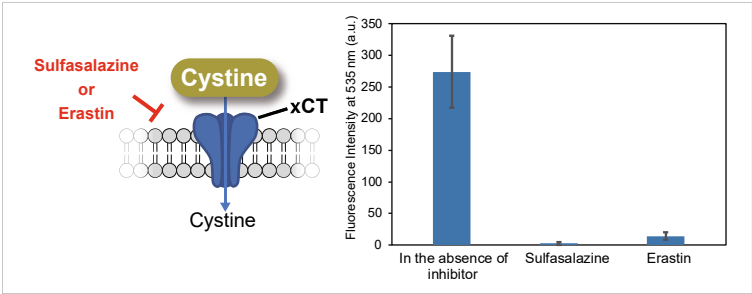
- Easier way to cystine uptake assay
- Applied for plate assay

The Cystine Analog (CA) in this kit can be taken up into cells via xCT, and the incorporated CA can be specifically detected using the Fluorescent Probe and Reducing Agent. Thus, the xCT activity can be measured easily.[Patent applied]



Evaluation of xCT inhibitor Sulfasalazine or Erastin

Using this kit, we measured the inhibitory effect of sulfasalazine and erastin on cystine uptake by HeLa cells. The fluorescence intensity of the sulfasalazine and elastin groups decreased significantly, indicating that both reagents inhibit cystine uptake.



Experiment Condiitons
Cell Line: HeLa cells
Pretreatment: DMEM (cystine-free, serum-free), 37°C, 5 min
Uptake conditions: 0.5 mmol/l sulfasalazine or 2 µmol/l erastin / Cystine Analog / DMEM (cystine-free, serum-free), 37°C, 30 min
Instrument: Fluorescent Plate Reader
Filter: Ex=485 nm, Em=535 nm

| Description | Unit | Code |
|--------------------------|-----------|---------|
| Cystine Uptake Assay Kit | 20 tests | UP05-10 |
| | 100 tests | UP05-12 |

| |
|--|
| Proliferation |
| Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |



Mitochondrial Research

Proliferation
Cytotoxicity
Senescence
Autophagy
Oxidative
Stress
Metabolism
Mitochondria
Lysosome
Endocytosis
Other Organelles
Exosome, Lipid Droplet, etc.

Mitochondrial Superoxide Detection

MitoBright ROS Deep Red - Mitochondrial Superoxide Detection

Allow to detecting mitochondrial superoxide with a long wavelength (Deep Red)

Singlet Oxygen Detection

Si-DMA for Mitochondrial Singlet Oxygen Imaging

Real-time visualization of $^1\text{O}_2$ generation

Lipophilic Peroxide Detection

MitoPeDPP

Live-cell fluorescent imaging of lipophilic peroxide

Ferrous Ion Detection

Mito-FerroGreen

Live-cell fluorescent imaging of intracellular Fe^{2+}

Mitophagy Detection

Mitophagy Detection Kit

Live-cell fluorescent imaging of mitophagy without transfection

Measurement of Glucose

Glucose Assay Kit-WST

Measurement of intracellular glucose concentrations via fluorescence

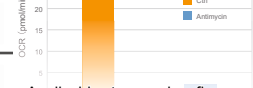
Measurement of Lactate

Lactate Assay Kit-WST

Measuring lactate to infer glycolytic activity

Oxygen consumption rate (OCR) Detection

Extracellular OCR Plate Assay Kit



Applicable to regular fluorescent plate reader with temperature-controlled incubation

Mitochondria Fluorescent Probe for Immunostaining

MitoBright IM Red for Immunostaining

Capable of co-stained with immunostaining. Higher retention in mitochondria after fixation & membrane permeabilization

Membrane Potential Detection

MT-1 MitoMP Detection Kit

Monitoring and observation even after fixation, with more sensitive detection than JC-1

Total ROS Detection

ROS Assay Kit - Highly Sensitive DCFH-DA

Detection with higher sensitivity than the original DCFH-DA

Mitochondrial Staining

MitoBright LT Series (Green / Red / DeepRed)

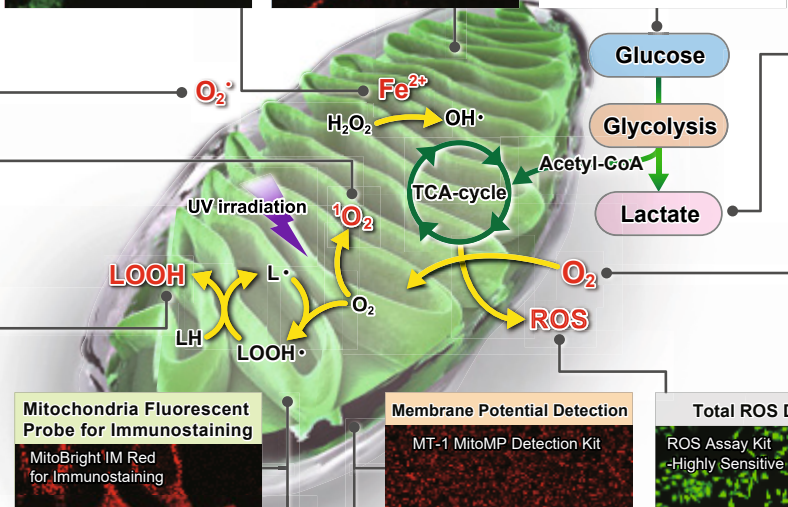
Green Red Deep Red

Selective staining of mitochondria in living cells

Membrane Potential Detection

JC-1 MitoMP Detection Kit

Analysis of mitochondrial membrane potential through fluorescence color ratios via microscopy, FCM, or microplate reader



| Description | Unit | Code |
|--|------------------------------------|---------|
| Metabolism | | |
| Extracellular OCR Plate Assay Kit | 100 tests | E297-10 |
| Glucose Assay Kit-WST | 50 tests | G264-05 |
| | 200 tests | G264-20 |
| Lactate Assay Kit-WST | 50 tests | L256-10 |
| | 200 tests | L256-20 |
| Mitochondrial Membrane Potential | | |
| MT-1 MitoMP Detection Kit | 1 set | MT13-10 |
| JC-1 MitoMP Detection Kit | 1 set | MT09-10 |
| Mitophagy | | |
| Mitophagy Detection Kit | 1 set | MD01-10 |
| Mtphagy Dye | 5 $\mu\text{g} \times 3$ | MT02-10 |
| Mitochondrial Staining | | |
| MitoBright LT Green | 400 μl | MT10-12 |
| MitoBright LT Red | 400 μl | MT11-12 |
| MitoBright LT Deep Red | 400 μl | MT12-12 |
| MitoBright IM Red for Immunostaining | 20 $\mu\text{l} \times 1$ | MT15-10 |
| | 20 $\mu\text{l} \times 3$ | MT15-12 |
| Oxidative Stress | | |
| MitoBright ROS Deep Red - Mitochondrial Superoxide Detection | 100 nmol $\times 1$ | MT16-10 |
| | 100 nmol $\times 3$ | MT16-12 |
| Mito-FerroGreen | 1 set (50 $\mu\text{g} \times 2$) | M489-10 |
| Si-DMA for Mitochondrial Singlet Oxygen Imaging | 2 μg | MT05-10 |
| MitoPeDPP | 5 $\mu\text{g} \times 3$ | M466-10 |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Proliferation
Cytotoxicity
Senescence
Autophagy
Oxidative
Stress
Metabolism
Mitochondria
Lysosome
Endocytosis
Other Organelles
Exosome, Lipid Droplet, etc.

Mitochondrial Research

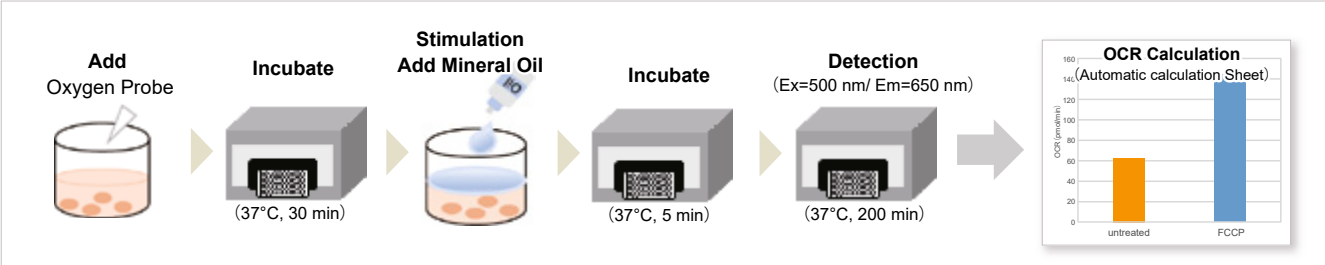
Extracellular OCR Plate Assay Kit



- Applicable to regular fluorescent plate reader with temperature-controlled incubation
- No need for an expensive instrument, special medium, and plates
- All-in-One Kit with OCR calculation Sheets



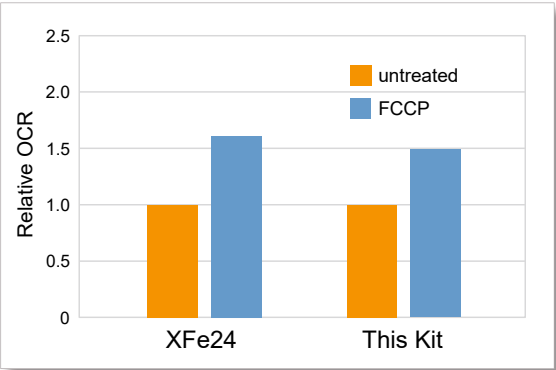
Procedure



Comparison with Flux Analyzer

Flux Analyzer (XFe24) and this kit were measured on the same day under the same conditions (cell type, cell number, and FCCP concentration). As a result, correlated data of oxygen consumption rate changes were obtained for XFe24 and this kit.

Cells: HepG2
Cell Number: 5×10^4 cells/well
Stimulation: FCCP (Carbonyl cyanide 4-(trifluoromethoxy) phenylhydrazone)
FCCP Concentration: 2 $\mu\text{mol/l}$



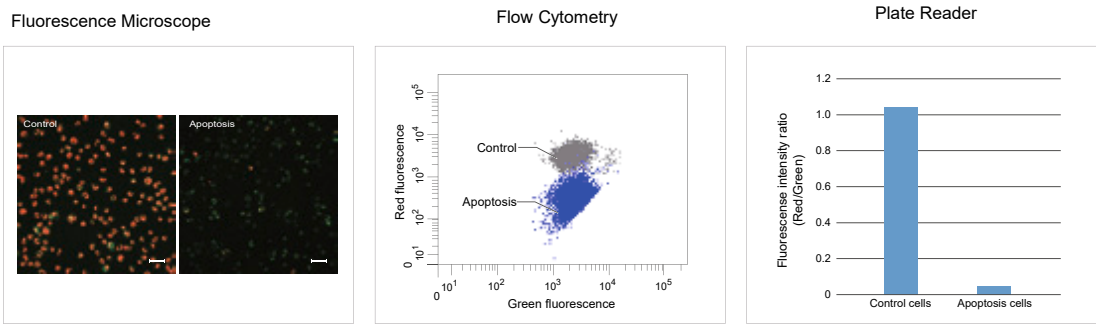
| Description | Unit | Code |
|-----------------------------------|-----------|---------|
| Extracellular OCR Plate Assay Kit | 100 tests | E297-10 |

Mitochondrial Membrane Potential Detection

JC-1 MitoMP Detection Kit



JC-1 forms aggregate (in healthy mitochondria) with red fluorescence. As membrane potential decreases, JC-1 becomes monomers, which shows in green fluorescence. The change in ratio of red to green fluorescence is used as an indicator of mitochondrial condition.



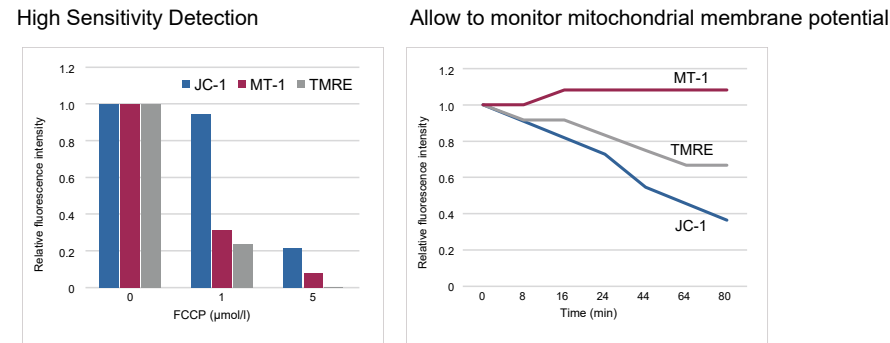
| Description | Unit | Code |
|---------------------------|-------|---------|
| JC-1 MitoMP Detection Kit | 1 set | MT09-10 |

Mitochondrial Membrane Potential Detection

MT-1 MitoMP Detection Kit



JC-1 dye, TMRE, and TMRM are widely used to monitor MMP, however, these dyes have some limitations, such as low photostability and poor retention after aldehyde fixation. These limitations result in poor reproducibility of experiments. Dojindo's MT-1 MitoMP Detection Kit overcomes these limitations. In addition, the Imaging Buffer included in this kit minimizes background fluorescence and maintains cell vitality while the assay is being performed.



| Description | Unit | Code |
|---------------------------|-------|---------|
| MT-1 MitoMP Detection Kit | 1 set | MT13-10 |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Proliferation
Cytotoxicity

Senescence

Autophagy

Oxidative
Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

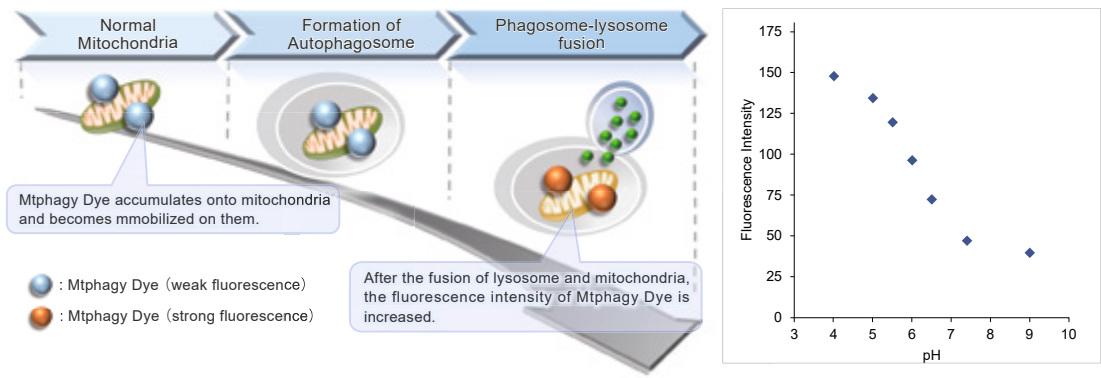
Other Organelles
Exosome, Lipid Droplet, etc.

Mitochondrial Research

Mitophagy Detection Kit

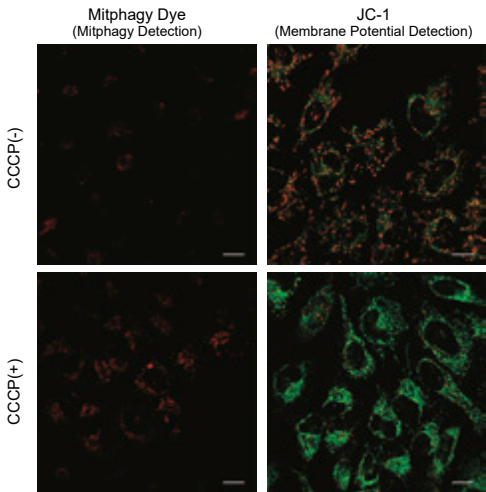


This kit is composed of Mtphagy Dye, reagent for detection of mitophagy, and Lyso Dye. Mtphagy Dye accumulates in intact mitochondria, is immobilized on it with chemical bond and exhibits a weak fluorescence from the influence of surrounding condition. When Mitophagy is induced, the damaged mitochondria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



The fluorescent intensity of Mtphagy Dye is increased at pH 4-5.

Mitophagy Induction and Mitochondrial Membrane Potential Changes



Mitochondrial condition in the carbonyl cyanide m-chlorophenyl hydrazine (CCCP) treated Parkin-expressing HeLa cells was compared with untreated cells using Mitophagy Detection Kit (MD01, MT02) and JC-1 MitoMP Detection Kit (MT09).

Result:

As a result, mitophagy was hardly detected in the CCCP-untreated cells, and mitochondrial membrane potential was maintained normally. On the other hand, in CCCP-treated cells, we observed a decrease in mitochondrial membrane potential (decrease in red fluorescence of JC-1) and induction of mitophagy (increase in fluorescence of Mtphagy Dye).

| Description | Unit | Code |
|-------------------------|----------|---------|
| Mitophagy Detection Kit | 1 set | MD01-10 |
| Mtphagy Dye | 5 μg × 3 | MT02-10 |

Mitochondrial Superoxide Detection

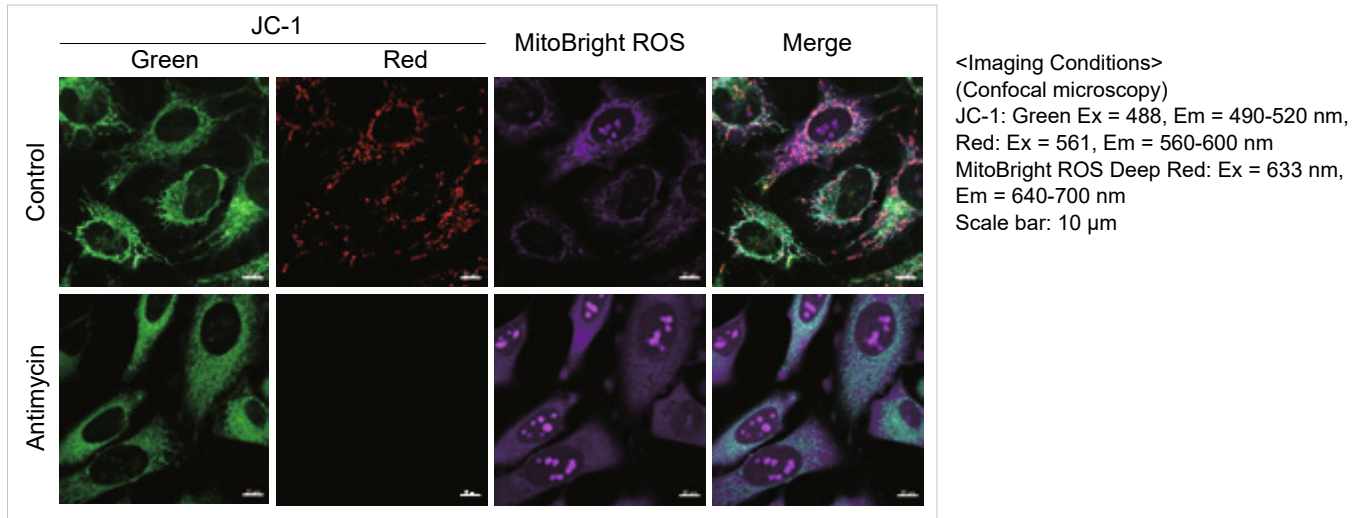
MitoBright ROS Deep Red - Mitochondrial Superoxide Detection



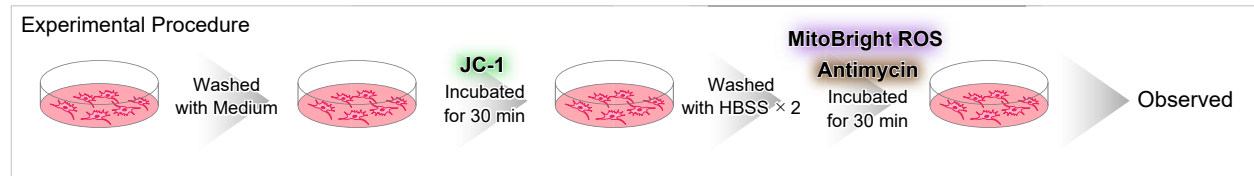
This dye emits deep red fluorescence; its fluorescence does not overlap with emission wavelengths that other red fluorescent markers use. Furthermore, the MitoBright ROS Deep Red is better able to selectively detect superoxide, compared to Company T's product Red.

Experimental Example

Simultaneously Evaluation of Mitochondrial Superoxide and Membrane Potential



After HeLa cells were washed with HBSS, co-stained with MitoBright ROS Deep Red and mitochondrial membrane potential staining dye (JC-1: code MT09), and the generated mitochondrial ROS and membrane potential were observed simultaneously. As a result, the decrease in mitochondrial membrane potential and the generation of mitochondrial ROS are simultaneously observed.



| Description | Unit | Code |
|--|--------------|---------|
| MitoBright ROS Deep Red - Mitochondrial Superoxide Detection | 100 nmol × 1 | MT16-10 |
| | 100 nmol × 3 | MT16-12 |

| |
|--|
| Proliferation Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

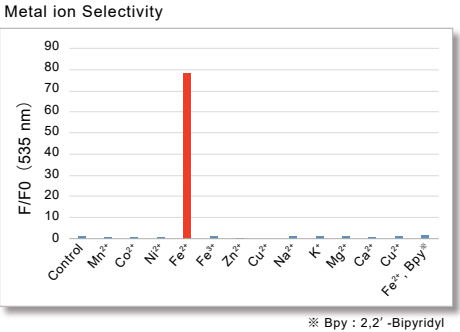
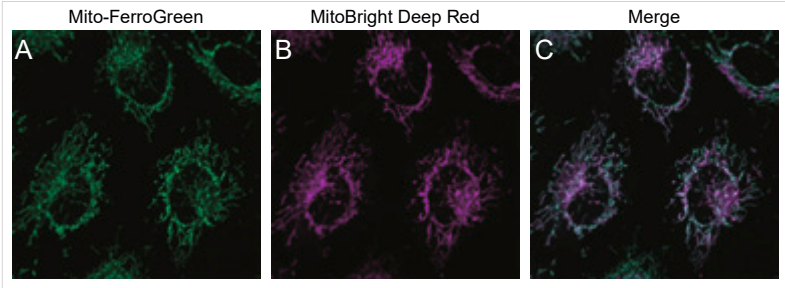
Mitochondrial Superoxide Detection

Mito-FerroGreen

Mito-FerroGreen is a novel fluorescent probe for the detection of ferrous ion (Fe²⁺) in mitochondria where Fe-S clusters and heme proteins are synthesized, and enables live cell fluorescent imaging of intracellular Fe²⁺. Mito-FerroGreen has no no chelating ability. Mito-FerroGreen and Fe²⁺ react irreversibly, which is different from the detection principle of calcium-iron probes such as Fluo-3.

Double staining with mitochondrial staining probe

HeLa cells incubated with Mito-FerroGreen and MitoBright Deep Red, treated with ammonium iron(II) sulfate, were observed by fluorescence microscopy.



Double staining with mitochondrial staining probe

Mito-FerroGreen (5 μmol/l) Ex/Em = 488 nm/ 500-550 nm

MitoBright Deep Red (200 nmol/l) Ex/Em = 640 nm/ 656-700 nm

A Mito-FerroGreen

B MitoBright Deep Red

C Merge

Iron Detection Dyes

| | Mito-FerroGreen (M489) | FerroOrange (F374) |
|----------------------|--|---|
| Localization | Mitochondria | Intracellular |
| Fluorescent Property | lex 505 nm, lem 535 nm | lex 543 nm, lem 580 nm |
| Instrument (filter) | Fluorescence microscope (FITC, GFP) | Fluorescence microscope, plate reader (Cy3) |
| Sample | Live Cell | Live cell |
| The number of assays | 1 set (50 μg x 2) 10 assays at 35 mm dish (final concentration 5 μmol/l) | 1 tube (24 μg) 17 assays at 35 mm dish (final concentration 1 μmol/l) |

| Description | Unit | Code |
|-----------------|-------------------|---------|
| Mito-FerroGreen | 1 set (50 μg × 2) | M489-10 |
| FerroOrange | 1 tube | F374-10 |
| | 3 tube | F374-12 |



Mitochondrial Staining

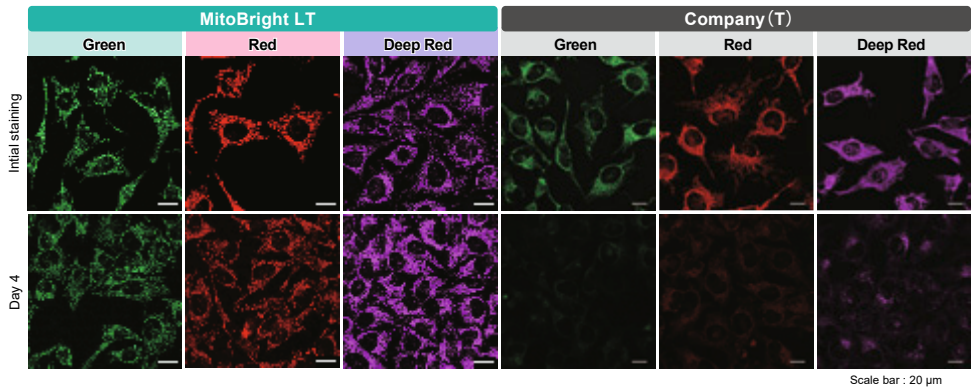
MitoBright LT Series



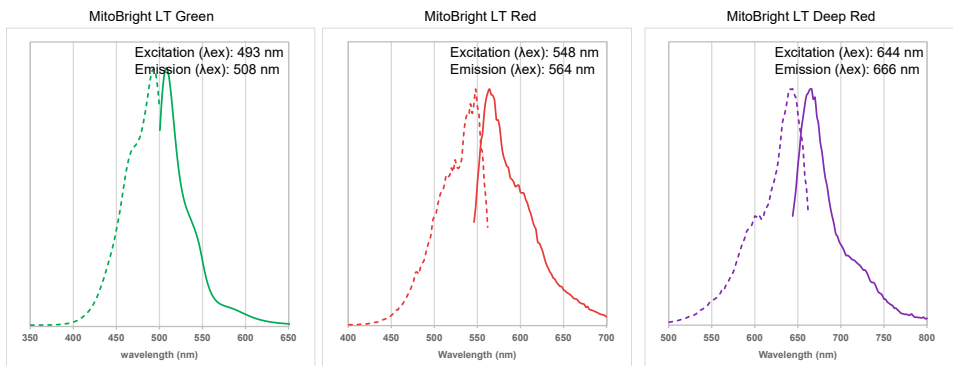
MitoBright LT dyes are designed to exhibit mitochondria retention for long-term visualization. In addition, the MitoBright LT dyes show stronger fluorescence signals compared with other commercially available dyes that contain the chloromethyl moiety. The MitoBright LT dyes offer three different color options (Green, Red and Deep Red), and are provided as a ready-to-use DMSO solution. A working solution can easily be prepared in a single dilution step with growth medium or HBSS.

Stained in serum-contained media

HeLa cells were stained with MitoBright LTs or an existing reagent and observed after 4 days. MitoBright LT remained unchanged and observable even after 7 days, while the existing reagent's intensity decreased.



Fluorescence Properties



| Description | Unit | Code |
|------------------------|--------|---------|
| MitoBright LT Green | 400 μl | MT10-12 |
| MitoBright LT Red | 400 μl | MT11-12 |
| MitoBright LT Deep Red | 400 μl | MT12-12 |

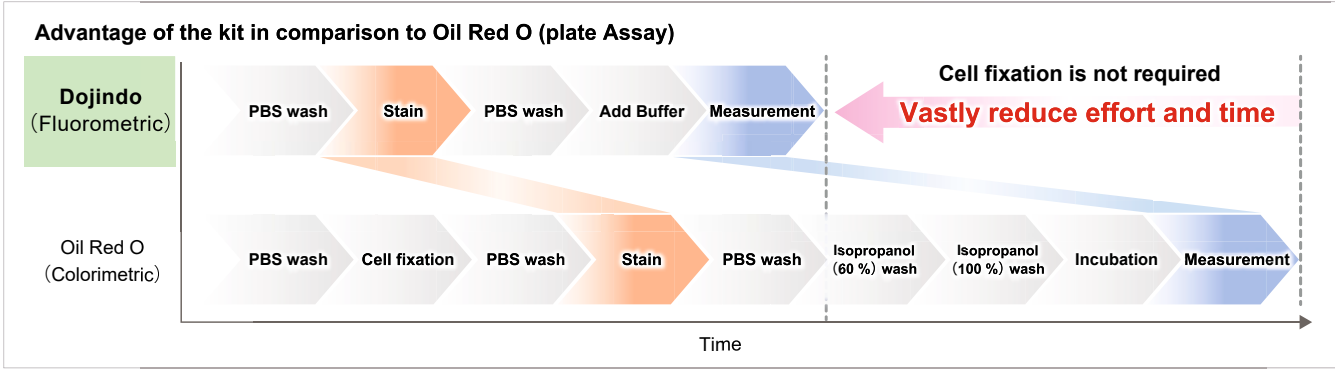
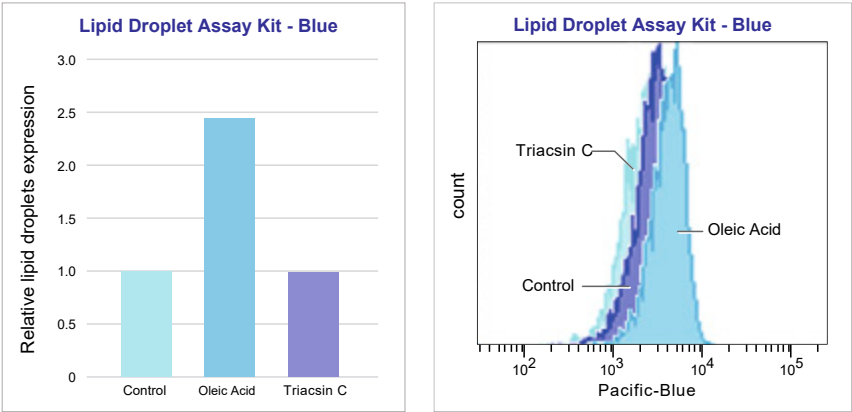
| |
|--|
| Proliferation |
| Cytotoxicity |
| Senescence |
| Autophagy |
| Oxidative Stress |
| Metabolism |
| Mitochondria |
| Lysosome |
| Endocytosis |
| Other Organelles Exosome, Lipid Droplet, etc. |

Lipid Droplet Staining

Lipid Droplet Assay Kit - Blue / Deep Red



The Lipid Droplet Assay Kit simplifies the quantification of fat droplets with provided protocols and buffers. It works for live cells, and its fluorescent dye is suitable for both live and fixed cells. Compared to colorimetric reagents, it reduces measuring time and increases experiment repeatability by avoiding dye deposition in the plate.



| Description | Unit | Code |
|----------------------------------|-------|---------|
| Lipid Droplet Assay Kit-Blue | 1 set | LD05-10 |
| Lipid Droplet Assay Kit-Deep Red | 1 set | LD06-10 |

Proliferation
Cytotoxicity
Senescence
Autophagy
Oxidative Stress
Metabolism
Mitochondria
Lysosome
Endocytosis
Other Organelles
Exosome, Lipid Droplet, etc.

Follow us on **LinkedIn**

