# **Reagent for Cellular Function Analysis**



7th Edition

# **Reagent for Cellular Function Analysis**

#### Autophagy

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

#### Senescence

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

## 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-

	Cancer	
Code	Product Name	Page
MT09	JC-1 MitoMP Detection Kit	31
MT10, MT11, MT12	MitoBright LT Series	35
MT13	MT-1 MitoMP Detection Kit	31
MT16	MtoBright ROS Deep Red - Mitochondrial Superoxide Detection	33
MD01	Mitophagy Detection Kit	32
MT02	Mtphagy Dye	32
E297	Extracellular OCR Plate Assay Kit	30
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	34
UP05	Cystine Uptake Assay Kit	27
L267	Lipid Peroxidation Probe -BDP 581/591 C11-	15
G270	Glycolysis/OXPHOS Assay Kit	22
A550	ATP Assay Kit-Luminescence	24
L256	Lactate Assay Kit-WST	23
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
G264	Glucose Assay Kit-WST	23
G268	Glutamate Assay Kit-WST	21

	Neurodegenerative Diseases	
Code	Product Name	Page
MT09	JC-1 MitoMP Detection Kit	31
MT10, MT11, MT12	MitoBright LT Series	35
MT13	MT-1 MitoMP Detection Kit	31
MT16	MtoBright ROS Deep Red - Mitochondrial Superoxide Detection	33
MD01	Mitophagy Detection Kit	32
MT02	Mtphagy Dye	32
E297	Extracellular OCR Plate Assay Kit	30
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	34
UP05	Cystine Uptake Assay Kit	27
L267	Lipid Peroxidation Probe -BDP 581/591 C11-	15
G270	Glycolysis/OXPHOS Assay Kit	22
A550	ATP Assay Kit-Luminescence	24
L256	Lactate Assay Kit-WST	23
L261	LysoPrime Green - High Specificity and pH Resistance	36
L264	LysoPrime Deep Red - High Specificity and pH Resistance	36
L265	pHLys Red - Lysosomal Acidic pH Detection	36
L266	Lysosomal Acidic pH Detection Kit	37
L268	Lysosomal Acidic pH Detection Kit-Green/Deep Red	37
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

	Senescence	
Code	Product Name	Page
SG03	Cellular Senescence Detection Kit - SPiDER-βGal	8
SG07	Cellular Senescence Detection Kit - SPiDER Blue	8
SG05	Cellular Senescence Plate Assay Kit - SPiDER-βGal	9
MT09	JC-1 MitoMP Detection Kit	31
MT10, MT11, MT12	MitoBright LT Series	35
MT13	MT-1 MitoMP Detection Kit	31
MT16	MtoBright ROS Deep Red - Mitochondrial Superoxide Detection	33
MD01	Mitophagy Detection Kit	32
MT02	Mtphagy Dye	32
E297	Extracellular OCR Plate Assay Kit	30
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	34
UP05	Cystine Uptake Assay Kit	27
L267	Lipid Peroxidation Probe -BDP 581/591 C11-	15
G270	Glycolysis/OXPHOS Assay Kit	22
A550	ATP Assay Kit-Luminescence	24
L256	Lactate Assay Kit-WST	23
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

#### Cell Proliferation / Cytotoxicity Assay Cell Counting Kit-8 Cytotoxicity LDH Assay Kit-WST Annexin V Apoptosis Plate Assay Kit

#### **Detection Principle**



Cell Counting Kit-8 measures the dehydrogenase activity with NADH in a live cell.



CK04

CK12

AD12

Annexin V Apoptosis Plate Assay Kit measures apoptosis using FITC-labeled Annexin V for phosphatidylserine binding.

#### **Procedure**





Annexin V Apoptosis Plate Assay Kit



The Annexin V Apoptosis Plate Assay Kit contains a reagent that quenches the fluorescence of Annexin V not bound to phosphatidylserine, allowing rapid detection of multiple samples using a plate reader without washing procedures.

#### **Experimental Example:** Changes in various indicators due to Staurosporine

HepG2 cells were treated with staurosporine to induce apoptosis, and phosphatidylserine, extracellular LDH and cell proliferation were detected. Phosphatidylserine was measured as an apoptosis marker using the Annexin V Apoptosis Plate Assay Kit, extracellular LDH was measured as an indicator of dead cells using the Cytotoxicity LDH Assay Kit-WST, and cell proliferation was measured using the Cell Counting Kit-8. The results showed that staurosporine treatment increased phosphatidylserine and extracellular LDH, and decreased cell proliferation.

#### Same Samples can be used

In addition, cell samples are separated into cells and culture media and measured with different indicators (using a combination of our kits), allowing more detailed analysis of cell death.



Description	Unit	Code
	500 tests	CK04-05
Coll Counting Kit 9	1000 tests	CK04-11
Cell Counting Kit-6	3000 tests	CK04-13
	10000 tests	CK04-20
	100 tests	CK12-01
Cytotoxicity LDH Assay Kit-WST	500 tests	CK12-05
	2000 tests	CK12-20
Annexin V Apoptosis Plate Assay Kit	100 tests	AD12-10

Senescence

Endocytosis

#### Senescence Detection Cellular Senescence Detection Kit - SPiDER-ßGal Cellular Senescence Detection Kit - SPiDER Blue

Cellular Senescence Detection Kit – SPiDER- $\beta$ Gal allows to detect SA- $\beta$ -gal with high sensitivity and ease of use. SPiDER- $\beta$ Gal is a new reagent to detect  $\beta$ -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  $\beta$ -galactosidase activity. Therefore, SPiDER- $\beta$ Gal can be applied to guantitative analysis by flow cytometry.

X-Gal



Cellular Senescence Detection Kit - SPiDER Blue

SPiDER-ßgal

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

8



Code

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



Proliferation Cytotoxicity

Senescence

Autophagy

This kit allows you to quantify SA- $\beta$ -gal activity and evaluate multiple samples in a 96-well plate by simply adding SPiDER- $\beta$ Gal, a reagent that can detect  $\beta$ -galactosidase.

#### 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-&Gal(Code SG04)) Blue: Ex. 405nm / Em. 450-495nm (Nuclear staining with -Cellstain- DAPI solution(Code D523))

#### Experimental Procedure



## **Cell Count Normalization Kit**

Combined Cellular Senescence Plate Assay protocol available online

#### When normalized to the results obtained by quantifying nucleic acids using the Cell Count Normalization Kit, the measured values of SA-Bgal activity become available for evaluation of SA-β-gal activity according to cell number.



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

Metabolism

Mitochondria

#### Autophagy DAPGreen / Red - Autophagy Detection DALGreen - Autophagy Detection

D675	D676	D677

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

	App Fluorescent	licable instrum	ents Microplate	Fluorescent properties	Volume / the number of usable assays	Existing methods	
DAPGreen	Microscope			Ex = 425-475 nm Em = 500-560 nm * For confocal microscope,the sample can be excited at 488 nm	5 nmol x 1 / 35 mm dish: 25 (when used in 1.0 μmol/l)	LC3-GFP	
DAPRed	0	×	×	Ex = 500-560 nm Em = 690-750 nm	5 nmol x 1 / 35 mm dish: 25 (when used in 1.0 µmol/l)	MDC Cyto-ID etc.	
DALGreen	0	0	×	Ex = 350-450 nm Em = 500-560 nm * For confocal microscope,the sample can be excited at 488 nm	20 nmol x 1 / 35 mm dish: 10 (when used in 1.0 µmol/l)	LC3-GFP-RFP etc.	

10

\*Double staining imaging by DAPGreen and DALGreen is not possible

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

## Autophagy Autophagic Flux Kit



Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Other Organelles Exosome, Lipid Droplet, etc.



#### 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

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### **Oxidative Stress ROS Assay Kit -Highly Sensitive DCFH-DA-**



#### The selectivity for ROS

DCFH-DA is widely used for ROS detection, but it has some limitations like weak fluorescence signals and high background. Dojindo's ROS Assay Kit -Highly Sensitive DCFH-DA- allows ROS detection with higher sensitivity than DCFH-DA with the similar ROS selectivity.



#### **Experimental Example:** High Sensitive Detection Compared with DCFH-DA



Ш Х			
Other Organelle psome, Lipid Drop	Hydrogen peroxide (H <sub>2</sub> O <sub>2</sub> )-treated HeLa cells ( $1 \times 10^4$ cells/ml) were stained with DC Highly Sensitive DCFH-DA, and the fluorescence intensity of intracellular ROS detection kits. As a result, the ROS Assay Kit-Highly Sensitive DCFH-DA in high-sense ROS was better than DCFH-DA.	FH-DA or the RC was compared I sitivity detection c	DS Assay Kit- between two f intracellular
let, e	Description	Unit	Code
etc.	ROS Assay Kit -Highly Sensitive DCFH-DA-	100 tests	R252-10

Endocytosis

## **Oxidative Stress ROS Assay Kit -Photo-oxidation Resistant DCFH-DA-**



Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

The dye used in this kit allows ROS detection with higher sensitivity than DCFH-DA and long-term observation of live cells due to its resistance to photooxidation.

#### Resistant to Photo-oxidation



Comparison of photo-oxidation resistant ability in HeLa cells



#### Experimental Example: Simultaneous Detection of ROS in LPS-treated macrophages



After induction of ROS, observation	was performed every ho	our	is et, etc.
In Lipopolysaccharide (LPS) treated RAW 264.7 cells, after being stained with Sensitive DCFH-DA, or Photo-oxidation Resistant DCFH-DA, the intracellular RC	regular DCFH-DA S level was comp	A, Highly bared.	ther Organelle ne, Lipid Dropl
Description	Unit	Code	
ROS Assay Kit -Photo-oxidation Resistant DCFH-DA-	100 tests	R253-10	Ŭ Ě

# Lipid Peroxide Detection



Liperfluo is a Dojindo-developed fluorescence probe to specifically detect lipid peroxides with minimal photodamage 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:** 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 $\mu$ g $ imes$ 5)	L248-10

14

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### 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.



#### **Experimental Example:** Lipid Peroxidation Assay



HepG2 cells stained with this probe were stimulated with HBSS solution containing 200 µmol/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.

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



#### Intracellular Iron Ion Measurement FerroOrange

FerroOrange is a specific probe for ferrous ions ( $Fe^{2+}$ ) and reacts irreversibly, which differs from the detection principle of chelating ability. Fluorescence imaging or quantification of intracellular  $Fe^{2+}$  is often used to evaluate iron-related non-apoptotic cell death, ferroptosis.



#### **Experimental Example**



HeLa cells treated with chelator of iron 2,2'-bipyridyl (Bpy) (100  $\mu$ mol/l) or Ammonium iron (II) sulfate (100  $\mu$ mol/l) were prepared. The change of intracellular Fe<sup>2+</sup> in HeLa cells was detected by the FerroOrange. Ex = 561 nm, Em = 570-620 nm, Scale bars 20  $\mu$ m

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

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

#### 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 ( $\lambda$ max = 412 nm) of DTNB (5,5-dithio-bis- (2-nitrobenzoic acid). The guantity of GSH can also be determined, by substracting GSSG from the total amount of glutathione. With this kit, GSH/GSSG concentrations from 0.5 µmol/l to 50 µmol/l and GSSG concentrations from 0.5 µmol/l to 25 µmol/l can be quantified.

#### GSH Masking GXH ← Masking reagent GSSH Glutathione reductase GSSG measurement GSH Glutathione reductase TNB DTNB 5-Mercapto-2-nitrobenzoic acid (λmax:412 nm) Mechanism of total glutathione quantification

#### **Experimental Procedure:** Experimental time is only 1-2 hours



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℃ for 1 h.



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



6)-7) After incubating at 37℃ for 10 minutes, measure the absorbane of each well with a microplate

#### Calibration Curve



GSSG/GSH Quantification Kit



## Related Products

## **Ferroptosis research**

"Ferroptosis" was coined by Stockwell et al. at Columbia University in 2012 and described as a form of irondependent cell death. \* It was reported to be a form of programmed cell death by the Nomenclature Committee on Cell Death (NCCD) in 2018.

Ferroptosis is a form of programmed cell death caused by iron ion-dependent accumulation of lipid peroxides. Ferroptosis has been shown to follow a different cell death pathway from apoptosis and thus is attracting attention as a new target for cancer therapy. It has also been found to be associated with various diseases, such as neurodegenerative diseases, cerebral apoplexy, and hepatitis (NASH).

#### Ferroptosis Pathway



#### Experimental Example: Evaluating Intracellular Uptake and Redox Balance

We investigated the transition of cellular metabolisms in A549 cells treated with Erastin, a known ferroptosis inducer. Our results revealed the following.



Endocytosis

## Measurements of Intracellular Metabolism



Description	Unit	Code
tarter Kit		
Glycolysis/OXPHOS Assay Kit	50 tests	G270-10
Glycolysis/JC-1 MitoMP Assay Kit	50 tests	G272-10
uantification for Intracellular Metabolism		
	50 tests	A550-10
ATP Assay Kit-Luminescence	200 tests	A550-12
ADP/ATP Ratio Assay Kit-Luminescence	100 tests	A552-10
	50 tests	G264-05
Glucose Assay Kit-WST	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-WS I	200 tests	L256-20
NAD/NADH Assay Kit-WST	100 tests	N509-10
NADP/NADPH Assay Kit-WST	100 tests	N510-10
ptake 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
Amine Asid Unteke Assess	20 tests	UP04-10
Amino Acid Optake Assay	100 tests	UP04-12
	20 tests	UP05-10
Cystine Optake Assay Kit	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.

Proliferation Cytotoxicity



#### Intracellular Metabolism **Glycolysis/OXPHOS Assay Kit**

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Exosome, Lipid Droplet, etc



- 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



Other Organelle	$\begin{array}{c ccccccccccccccccccccccccccccccccccc$	Synthesizes ATP dependent on glycolysis	0 0 0 0 0 0 0 0 0 0 0 0 0 0	9 0.4 9 0.2 0 0 0 0 0 0 0 0 0 0 0 0 0 0 0	Synthesizes ATP dependent on oxidative phosphorylat		
S,		Description			Unit	Code	
	Glycolysis/OXPHOS Assay Kit				50 tests	G270-10	

## Intracellular Metabolism Glucose Assay Kit-WST Lactate Assay Kit-WST





Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

**Metabolism** 

Mitochondria

Lysosome

Endocytosis

1256

The Glucose Assay Kit-WST and Lactate Assay Kit-WST are colourimetric kits for the quantification of glucose and lactate, respectively, both with a lower detection limit of 0.02 mmol/L. These two indicators are crucial for understanding glycolytic metabolism and are among the most commonly measured parameters in metabolic studies. In addition, the use of Dojindo's proprietary WST dye in the assay systems allows for easy and highly accurate detection.



#### **Experimental Example:** Evaluation of culture supernatant using two indicators

Phloretin, the glucose transporter inhibitor, was added to Jurkat cells and the intracellular metabolism change was evaluated using Glucose Assay Kit-WST and Lactate Assay Kit-WST.







#### <Description>

Glucose consumtion has decreased due to inhibition of glucose uptake by Phloretin, resulting in increase of glucose and decrease in lactate in culture supernatant.

glucose and decrease in lactate in culture supernatant.			elles oplet, etc.	
Description	Unit	Code	rgan bid D	
Glucose Assay Kit-WST	50 tests 200 tests	G264-05 G264-20	Other O me, Lip	
Lactate Assay Kit-WST	50 tests 200 tests	L256-10 L256-20	Exosol	

#### Intracellular Metabolism **ATP Assay Kit-Luminescence**

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism



ATP Assay Kit-Luminescence is a luciferase luminescence assay kit for quantification of intracellular ATP. ATP is an important energy source of living cells that is synthesized in both glycolysis and mitochondrial oxidative phosphorylation. Mitochondrial dysfunction reduces ATP levels in cells and the decreased ATP levels are known to be associated with cancer, aging, and neurodegenerative diseases. Therefore, ATP level is used as an indicator for mitochondrial activity. ATP level is a lso focused in Cancer research since the recent studies have revealed that although cancer cells were known to rely on glycolysis for ATP synthesis, a shift from glycolysis to oxidative phosphorylation occurs when glycolysis is suppressed.

#### **Experimental Example:** Change in intracellular metabolism of rotenone-treated cells

Rotenone, which is known to inhibit the mitochondrial electron transport chain, was added to Jurkat cells, followed by measurement of intracellular ATP using the ATP Assay Kit-Luminescence. As a result, ATP production in the mitochondrial respiratory chain (the electron transport chain) was inhibited, and ATP concentrations were lower than those in the control cells.



Other Organ			
elles	Description	Unit	Code
0,	ATP Assay Kit-Luminescence	50 tests 200 tests	A550-10 A550-12

#### Intracellular Metabolism Glucose Uptake Assay Kit



Proliferation Cytotoxicity

•	Highly	sensitive	and simple	measurement	of gluco	se uptake	capacity

- Applicable for microscopy & FCM
- Reduces dye leakage after staining

#### **Principle**



#### Experimental Example

Glucose uptake enhancement by insulin



Detailed results and other data are available online.

#### Comparison with Existing Method

The comparison of the Glucose Uptake Probe Series and the existing method(2-NBDG) is as below.

product name	Fluorescence microscope	Plate reader detection	FCM detection	Retention ability	Fluorescence characteristics
Glucose Uptake Assay Kit-Blue	0	×	0	1 hour *	λ <sub>ex</sub> :386 nm λ <sub>em</sub> :474 nm
Glucose Uptake Assay Kit-Green	0	0	0	1 hour *	λ <sub>ex</sub> :507 nm λ <sub>em</sub> :518 nm
Glucose Uptake Assay Kit- <mark>Red</mark>	0	0	0	1 hour *	λ <sub>ex</sub> :560 nm λ <sub>em</sub> :572 nm
2-NBDG	0	×	0	30 minutes or less *	λ <sub>ex</sub> :465 nm λ <sub>em</sub> :540 nm

\*Result of A549 cells, the retention time for other cell lines may be different.

Description	Unit*	Code
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

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### Intracellular Metabolism Amino Acid Uptake Assay Kit

**Principle** 



After amino acid analogs (BPA) are taken up into cells via amino acid transporters, the fluorescent probe permeates the cell membrane and binds to the amino acid analogs, emitting fluorescence

#### **Experimental Example:** Evaluation of BCH (Inhibitor of amino acids transporter)



Experiment Condiitons Cell Line: HeLa cells Medium: MEM (5.5 mmol/l Glucose) Incubation: 1 mmol/l BCH/HBSS (Hanks' Balanced Salt Solution), 37°C, 30 min Instrument: Fluorescent Microscopy (Ex=340-380 nm, Em: 435-485 nm) Instrument: Plate Reader (Ex=360 nm, Em: 460 nm)

:	Description	Unit	Code
	Amino Acid Uptake Assay Kit	20 tests 100 tests	UP04-10 UP04-12



Proliferation Cytotoxicity

#### Intracellular Metabolism Cystine Uptake Assay Kit

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

### **Principle**



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]

#### Experimental Example: 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.



Description	Unit
Cystine Uptake Assay Kit	20 tests 100 tests



UP05-10 UP05-12

## Mitochondrial Research





## Proliferation Cytotoxicity

Description	Unit	Code	ion
Metabolism			liferat
Extracellular OCR Plate Assay Kit	100 tests	E297-10	Cyt
	50 tests	G264-05	Sence
Giucose Assay Kit-WST	200 tests	G264-20	Senesc
	50 tests	L256-10	
Laciale Assay Kil-WST	200 tests	L256-20	tophag
Mitochondrial Membrane Potential			Aut
MT-1 MitoMP Detection Kit	1 set	MT13-10	tive
JC-1 MitoMP Detection Kit	1 set	MT09-10	Oxida
Mitophagy			E
Mitophagy Detection Kit	1 set	MD01-10	abolis
Mtphagy Dye	$5\mu g imes 3$	MT02-10	Met
Mitochondrial Staining			ndria
MitoBright LT Green	400 µl	MT10-12	itocho
MitoBright LT Red	400 µl	MT11-12	້ <b>E</b>
MitoBright LT Deep Red	400 µl	MT12-12	osome
MitePright IM Ded for Immunosteining	20 µl×1	MT15-10	Lys
	20 µl×3	MT15-12	tosis
Oxidative Stress			ndocy
MiteDright DOC Deep Ded. Mitechandriel Currensuide Detection	100 nmol × 1	MT16-10	ш ig
MitoBright ROS Deep Red - Mitochondrial Superoxide Detection	100 nmol × 3	MT16-12	lles plet, e
Mito-FerroGreen	1 set (50 $\mu$ g $ imes$ 2)	M489-10	Drgane oid Drc
Si-DMA for Mitochondrial Singlet Oxygen Imaging	2 µg	MT05-10	Other C me, Lij
MitoPeDPP	$5~\mu{ m g} imes3$	M466-10	Exoso

#### 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

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Exosome, Lipid Droplet, etc



#### 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<sup>4</sup> cells/well Stimulation: FCCP (Carbonyl cyanide 4-(trifluoromethoxy) phenylhydrazone) FCCP Concentration: 2 µmol/l



Other Organelles Description E297-10 100 tests Extracellular OCR Plate Assay Kit E297-12 300 tests

30

#### Mitochondrial Membrane Potential Detection JC-1 MitoMP Detection Kit MT-1 MitoMP Detection Kit



Proliferation Cytotoxicity

Senescence

JC-1 indicates mitochondrial condition by changing from red to green fluorescence as mitochondrial membrane
potential (MMP) decreases. While JC-1 and similar dyes such as TMRE and TMRM are popular for MMP detection,
they suffer from low photostability and poor retention. Dojindo's MT-1 MitoMP Detection Kit overcomes these
limitations and improves experimental reproducibility.

#### Comparison of Reagents

	Features	Sensitivity	Fixation	Monitoring	Fluorescence change (upon loss of mitochondrial membrane potential)	Detection (ex/em)
<b>JC-1</b> (JC-1 MitoMP Detection Kit)	Recomended for starting-up	\$			Color change from red to green	Green: 450-490 nm / 500-550 nm Red: 530-560 nm / 570-640 nm
MT-1 (MT-1 MitoMP Detection Kit)	Recommended for more detailed analysis	✓ (High)	1	1	Decrease in fluorescence intensity	530-560 nm / 570-640 nm
TMRE	Widely used	✓ (High)			Decrease in fluorescence intensity	530-560 nm / 570-640 nm

## Experimental Example

JC-1

Detection of MMP in Apoptotic Cells







Allow to monitor MMP

**MT-1** 



Description	Unit*	Code
JC-1 MitoMP Detection Kit	1 set	MT09-10
MT-1 MitoMP Detection Kit	1 set	MT13-10

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### 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 incresased at pH 4-5.

9 10

#### Experimental Example

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).

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	$5~\mu{ m g} imes3$	MT02-10

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

Autophagy

#### Mitochondrial Superoxide Detection MitoBright ROS Deep Red - Mitochondrial Superoxide Detection

fluorescent markers use. Furthermore, the MitoBright ROS Deep Red is better able to selectively detect

superoxide, compared to Company T's product Red.

This dye emits deep red fluorescence; its fluorescence does not overlap with emission wavelengths that other red



Proliferation Cytotoxicity



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.

Experimental Procedure Washed with Medium Washed for 30 min Washed MitoBright RC Mathematical MitoBright RC Antimycin Incubated for 30 min MitoBright RC	Observed	I
Description	Unit*	Code
MitoBright ROS Deep Red - Mitochondrial Superoxide Detection	100 nmol $ imes$ 1 100 nmol $ imes$ 3	MT16-10 MT16-12

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### Mitochondrial Superoxide Detection Mito-FerroGreen

Mito-FerroGreen is a novel fluorescent probe for the detection of ferrous ion (Fe<sup>2+</sup>) in mitochondria. Mito-FerroGreen has no chelating ability. Mito-FerroGreen and Fe<sup>2+</sup> react irreversibly, which is different from the detection principle of calcium-iron probes such as Fluo-3.

#### **Experimental Example** 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

Double staining with mitocondrial staining probe

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

## Selection Guide of Iron Detection Dyes

	Mito-FerroGreen (M489)	FerroOrange (F374)
Localization	Mitochondria	Intracellular
Fluorescent Property	λex 505 nm, λem 535 nm	λex 543 nm, λem 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*

	Description	Unit*	Code
Mito-FerroGreen		1 set (50 $\mu$ g $ imes$ 2)	M489-10
FerroOrange		1 tube 3 tube	F374-10 F374-12

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.





※ Bpy : 2,2' -Bipyridyl

Autophagy

#### Mitochondrial Staining MitoBright LT Series



Proliferation Cytotoxicity

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.

#### **Experimental Example:** 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 4 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

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### Lysosomal Analysis LysoPrime Green / Deep Red - High Specificity and pH Resistance pHLys Red - Lysosomal Acidic pH Detection





#### Resistance to pH changes



pHLys Red is highly specific to lysosomes and shows pH-dependent changes in fluorescence, and pH-resistant LysoPrime Green/Deep Red is retained in lysosomes even after adding Bafilomycin A1, a lysosomal acidity inhibitor. The lysosomal pH and mass of the same sample can be measured using these two dyes for a detailed analysis of lysosomal function.

#### pH-dependent Probe pHLys Red



#### pH dependence of pHLys Red



The fluorescence intensity of pHLys Red at each pH was confirmed in vitro, and it was confirmed that the fluorescence intensity changed sensitively within the range of lysosomal pH (pH 4.0-5.5).

Proliferation Cytotoxicity

Stress

Mitochondria

# Lysosomal Analysis Lysosomal Acidic pH Detection Kit

L266



The kit includes lysosome staining dyes, pHLys Red/Green (pH dependent), and LysoPrime Green/Deep Red (pHindependent). The pHLys and LysoPrime dyes accumulate in the intact lysosomes. The fluorescence intensity of pHLys dyes are enhanced as the acidity increases, and weak fluorescence is observed when lysosomes are neutralized due to the lysosomal dysfunction. On the other hand, LysoPrime dyes gives stable emissions even lysosomes are neutralized. Lysosomal pH and lysosomal mass can be measured by combining these pHLys and LysoPrime dyes.



#### Imaging Analysis: Green/Red (#L266-10)



<Experimental Conditions> LysoPrime Green: Ex = 488 nm, Em = 490 – 550 nm pHLys Red: Ex = 561 nm, Em = 560 – 620 nm

## FCM Analysis: Green/Deep Red (#L268-10)



#### <Experimental Conditions> pHLys Green: FITC Filter (Ex = 488 nm, Em = 515 – 545 nm) LysoPrime Deep Red: APC Filter (Ex = 640 nm, Em = 650 – 670 nm)

Description	Unit*	Code
Lysosomal Acidic pH Detection Kit – Green/Red *1	1 set	L266-10
Lysosomal Acidic pH Detection Kit – Green/Deep Red *2	1 set	L268-10
LysoPrime Green – High Specificity and pH Resistance	$10~\mu$ l $ imes$ 1 10 $\mu$ l $ imes$ 3	L261-10 L261-12
LysoPrime Deep Red - High Specificity and pH Resistance	1 tube 3 tube	L264-10 L264-12
pHLys Red - Lysosomal Acidic pH Detection	1 tube 3 tube	L265-10 L265-12

<sup>\*1</sup> Green/Red: combination of LysoPrime Green and pHLys Red, <sup>\*2</sup> Green/Deep Red: combination of pHLys Green and LysoPrime Deep Red \*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### Endocytosis ECGreen-Endocytosis Detection



ECGreen-Endocytosis Detection is a pH dependent fluorescence dye that localizes to vesicle membrane. The visualization of endocytosis using the ECGreen is a more direct method than fluorescent analogs and allows visualization endocytosis from the stage of early endosomes.

#### The detection mechanism of endocytosis



#### Clear visualization of intracellular vesicular trafficking

It has been known that Wortmannin inhibits the recycling of endosomes or transition to lysosomes and causes enlargement of endosomes. To evaluate these changes caused by Wortmannin, early endosomes were co-stained by ECGreen and Rab5-RFP (marker protein of early endosomes), and lysosomes were co-stained by ECGreen and lysosome staining reagent. In adding Wortmannin, ECGreen was colocalized with enlarged endosomes (Rab5-RFP). On the other hand, ECGreen wasn't colocalized with lysosomes.



Description Unit\* Code ECGreen-Endocytosis Detection 40 ul F296-10

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

Autophagy

### Endocytosis AcidSensor Labeling Kit – Endocytic Internalization Assay



This kit is an all-in-one kit that allows visualization of the endocytosis uptake of a target substance. The NH<sub>2</sub>-Reactive AcidSensor (fluorescent probe) included in the kit has an intramolecular active ester group that forms a stable covalent bond when mixed with an amino group-containing target substance (protein). The AcidSensor label can be excited at 633 nm, allowing for multiple staining with green or red fluorescence (Figure 1). The AcidSensor label shows little fluorescence in neutral conditions and fluoresces when acidified in the cells where it is taken up by endocytosis (Figure 2).



This kit includes a filtration tube necessary to remove the unreacted dye, and allows you to perform everything from labeling to purification operations.\* In addition, even first-time users can easily label AcidSensor by conducting experiments according to the instruction manual. \* Protein/Antibody is not included.



39

Proliferation

Senescence

Autophagy

Other Organelles Exosome, Lipid Droplet, etc.

#### Lipid Droplet Staining Lipi-Blue / Green / Red / Deep Red



Lipi probes are small molecules that emit strong fluorescence in a hydrophobic environment such as LDs, which can be observed without any washing steps after staining with Lipi probes.



A medium that contained oleic acid (200  $\mu$ mol/l) was added and incubated overnight. Then, the supernatant was removed and the cells were washed with PBS. Each Lipi product series (1  $\mu$ mol/l) was added and the cells were incubated for 15 minutes.

Lipi-Blue: Ex. 405 nm / Em. 450 – 500 nm, Lipi-Green: Ex. 488 nm / Em. 500 – 550 nm, Lipi-Red: Ex. 561 nm / Em. 565 – 650 nm, Lipi-Deep Red: Ex.640 nm / Em.650-700 nm

#### **Comparison of Reagents**

	Dojindo		Dojindo Other Products		6	
	Lipi-Blue	Lipi-Green	Lipi-Red	Oil Red O	Nile Red	Reagent B
Live Cells	1	1	1		1	1
Fixed Cells	1	1	1	1	1	1
Selectivity towards Lipid Droplet (Level of Background)	1	1	1			
General Filter Accommondation*1	1	1	✓*2	n.d.	*3	1
Retention in Live Cells	1	1		n.d.		

\*1 Please refer to our website for the co-staining filter.

 $^{'2}$  When co-staining with a green fluorescent dye, a green fluorescent emission filter less than 550 nm is recommended.  $^{'3}$  Leaks in GFP filter (500  $\sim$  540 nm)

Description	Unit*	Code
Lipi-Blue	10 nmol $ imes$ 1	LD01-10
Lipi-Green	10 nmol $ imes$ 1	LD02-10
Lipi-Red	100 nmol $ imes$ 1	LD03-10
Lipi-Deep Red	10 nmol $ imes$ 1	LD04-10

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

Endocytosis

## Lipid Droplet Staining Lipid Droplet Assay Kit - Blue / Deep Red

Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

Metabolism

Mitochondria

Lysosome

Endocytosis

Other Organelles Exosome, Lipid Droplet, etc.

The Lipid Droplet Assay Kit simplifies the quantification of fat droplets with provided protocols and buffers. It works 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.

#### Experimental Example of plate assay and flow cytometry



Changes in lipid droplets were examined after the addition of oleic acid or Triacsin C (acyl-CoA synthetase inhibitor) to the A549 cell culture medium.

#### Advantage of the kit in comparison to Oil Red O (Plate Assay)



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

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

105

#### Cell Membrane Staining PlasMem Bright Green / Red



PlasMem Bright dyes are designed to stain plasma membrane for over a day. Furthermore, the PlasMem Bright dyes are more water-soluble compared with other commercially available dyes and can be diluted with culture medium. The PlasMem Bright dyes offer two different color options (green and red) and are provided as ready-to-use DMSO solutions.

#### **Experimental Example:** High retentivity on plasma membrane



#### Fluorescence Properties



Description	Unit*	Code
PlasMem Bright Green	100 µl	P504-10
PlasMem Bright Red	100 µl	P505-10

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

Endocytosis

#### Nucleolus Staining Nucleolus Bright Green / Red

#### Detection Principle

Nucleolus Bright reacts to RNAs present besides nucleolus, but it shows strong fluorescence in nucleolus, which is the site of rRNA production. We recommend to co-stain with DAPI in order to image nucleolus clearly.



Nuclear

Nucleolus

rRNA

	Maximum Excitation Wavelength	Maximum Emission Wavelength	Fluorescence of MeOH fixed cells	Fluorescence of PFA fixed cells
Nucleolus Bright Green	513 nm	538 nm	0	0
Nucleolus Bright Red	537 nm	605 nm	0	0

#### Nucleolus Localization



Description	Unit*	Code
Nucleolus Bright Green	60 nmol	N511-10
Nucleolus Bright Red	60 nmol	N512-10

Proliferation Cytotoxicity

Nucleolus Bright Dye

Autophagy

\*Approximate usage depends on the experiment. Please refer to the Q&A section on our product webpage.

#### Exosome Staining Exosome Labeling Kits



The ExoSparkler series can be used to stain purified exosomal membrane or protein and allows imaging of labeled exosomes taken up by cells.



#### Labelling Procedure



ExoSparkler series contains filtration tubes available for the removal of dyes unreacted after fluorescence labeling, as well as an optimized protocol for labeling exosomes. Our ExoSparkler series makes it possible to prepare fluorescence labeling of exosomes using the simple procedure.

Description	Unit	Code
ExoSparkler Exosome Membrane Labeling Kit-Green	5 samples	EX01-10
ExoSparkler Exosome Membrane Labeling Kit-Red	5 samples	EX02-10
ExoSparkler Exosome Membrane Labeling Kit-Deep Red	5 samples	EX03-10
Exosparkler Exosome Protein Labeling Dye-Green	5 samples	EX04-10
Exosparkler Exosome Protein Labeling Dye-Red	5 samples	EX05-10
Exosparkler Exosome Protein Labeling Dye-Deep Red	5 samples	EX06-10

Proliferation Cytotoxicity



Proliferation Cytotoxicity

Senescence

Autophagy

Oxidative Stress

#### Easy to Use No Special Technique Required



Exo*lsolator* Exosome Isolation Kit can collect exosomes from cell supernatants with a recovery rate equivalent to the ultracentrifugation(UC) method. Science Exo*lsolator* Exosome Isolation Kit requires only the filtration procedure, unlike the UC, exosomes are obtained quickly without any complicated operations.

n.s.

Ultracentrifugation Exolsolator

Fig. 2a The number of particles

#### Recovery Rate Equivalent to Ultracentrifugation



Fig. 1 Particle size distribution



We isolated the exosomes from the supernatant of HEK293S using both the ultracentrifugation method and the Exo*lsolator* method. The results showed that the Exo*lsolator* recovered exosomes with the same particle size distribution and the number of particles as the ultracentrifugation method, and the amount of exosome marker expressions per protein were higher, indicating that Exo*lsolator* recovered exosomes with higher purity than the ultracentrifugation method.

Fig. 2b Expression level of exosome markers

Description	Unit	Code
Exolsolator Exosome Isolation Kit	3 tests	EX10-10
Exolsolator Isolation Filter	10 pieces	EX11-10

# Product line up for in vitro Diagnostics

For inquiries, please contact Dojindo Europe GmbH.

# Follow us on Linked in





European Headquarters

#### **Dojindo Europe GmbH**

Leopoldstr. 254, 80807, Munich, Germany Email: info@dojindo.eu.com Web: www.dojindo.com/EUROPE/