

2020 No.172 ISSN 0385-1516

マイグラソーム

Migrasomes and migracytosis, an emerging field in cell biology 清華大学 Li Yu (俞立) 北京大学 Yang Chen (陈扬) 清華大学 Dongju Wang (王东锔)

一注目の研究一

二重機能性蛍光色素(H-V)を用いた

フェロトーシスの解明

株式会社同仁化学研究所 北村 怜奈

新製品 エクソソーム蛍光染色キット 🖽 細胞周期測定試薬 💁 脂肪滴染色試薬・キット 📭 ミトコンドリア染色試薬 📭



表紙撮影:熊本県天草 「戦国時代に、この地の武将が山頂で 舞いを楽しんだ」という言い伝えから、 「高舞登山」と命名された。 photo:永島俊介氏

CONTENTS

Review Migrasomes and migracytosis, an emerging field in cell biology ● [1] 清華大学 Li Yu (俞立) 北京大学 Yang Chen (陈扬) 清華大学 Dongju Wang (王东锔)

Topics on Chemistry

二重機能性蛍光色素(H-V)を用いたフェロトーシスの解明 • [6] 株式会社同仁化学研究所 北村 怜奈

Commercial

細胞内鉄検出試薬 ●────[6] _____[**7**] 細胞内過酸化脂質検出試薬 ●──── グルタチオン測定キット ● [7] ミトコンドリア研究用試薬 ● [8] 細胞内代謝測定キット •-----[10] 実験例 ミトコンドリア電子伝達系の阻害実験 ●-----[11] 抗酸化能測定キット -DPPH ラジカル消去活性 ●-----[12] 抗酸化能測定キット -SOD 様活性 ● [12] 新製品 エクソソーム蛍光染色キット ● [4] 細胞周期測定試薬 •-----[4] 脂肪滴染色試薬・キット •------_____[5] ミトコンドリア染色試薬 ● [9] - 関連アプリケーションの紹介-フェロトーシス研究 ● [7] FAQ DPPH Antioxidant Assay Kit •----[13]

お知らせ

食品機能性評価・微生物検出試薬冊子 第二版 ●───[12] 細胞機能解析冊子 第3版 ●───[13] 販売中止製品のお知らせ ●────[13]

※希望納入価格には消費税等は含まれておりません。



Migrasomes and migracytosis, an emerging field in cell biology



Li Yu (俞立)

The State Key Laboratory of Membrane Biology, Tsinghua University-Peking University Joint Center for Life Sciences, School of Life Sciences, Tsinghua University.



Yang Chen (陈扬)

Center for Precision Medicine Multi-Omics Research, Peking University Health Science Center, Peking University.



Dongju Wang (王东锔)

The State Key Laboratory of Membrane Biology, Tsinghua University-Peking University Joint Center for Life Sciences, School of Life Sciences, Tsinghua University.

Abstract:

Migrasomes are newly discovered cellular organelles generated through migracytosis. They function in active intracellular communications which are essential for multiple core biological processes. In this review, we summarize the current research progress on migrasomes, including the molecular mechanisms involved in migrasome biogenesis and the physiological significance of migrasomes. We also propose the potential functions of migrasomes and the future directions of this emerging field.

Main text:

Migrasomes are newly discovered cellular organelles, first described in 2015 by Liang Ma et al^{1, 2} (Figure 1). Migrasomes are vesicles with diameters on the micron scale which contain numerous small vesicles inside. They are generated on the tips or intersections of retraction fibers (RFs), the long membrane projections left by cells when they migrate away. During biogenesis, cytosolic components are actively transported to migrasomes. Once the generation process is completed, migrasomes are either broken to release their contents into the extracellular space or taken up by incoming cells. The process of migrasome release by migrating cells is called migracytosis. Thus, migrasomes and migracytosis are proposed as mechanisms for spatial and temporal cell-cell communications.

1. Molecular mechanisms of migrasome biogenesis

1.1. Integrins provide the adhesion force for migrasome formation

Generation of migrasomes is dependent on cell migration. The number of migrasomes increases when migration is enhanced and decreases when migration is inhibited¹. In cultured cells, migrasomes adhere to the extracellular matrix (ECM) where they are formed, which indicates that adherent molecules play a role in the process of migrasome formation. Mass spectrometry analysis revealed that integrin $\alpha 5\beta 1$ is enriched on migrasomes. Integrin $\alpha 5\beta 1$ appears on retraction fibers before the actual growth of migrasome vesicles. Moreover, integrin $\alpha 5\beta 1$ is located at the bottom of migrasomes³. These spatial and temporal distribution features suggest that integrin $\alpha 5\beta 1$ specifies the site of migrasome formation. There are 18α and 8β integrins in mammals. Different integrins bind to different ECM proteins. Interestingly, we found that the pairing of integrins with their specific ECM partners is a determinant for migrasome formation³. This



Figure 1. TSPAN4 is enriched on migrasomes. L929 cells stably expressing TSPAN4-mcherry were plated on a glass dish coated with fibronectin and observed by confocal microscopy. Scale bar: 50 µm.

regulatable specificity provides great potential for migrasomes to function *in vivo* as signal transmitters.

1.2. Tetraspanins regulate migrasome formation

Tetraspanin 4 was found to be enriched on migrasomes and has been used as a marker for migrasomes in previous studies^{1,4}. There are 33 tetraspanins in the mouse genome. Overexpression of 14 of these tetraspanins can enhance migrasome formation in Normal Rat Kidney (NRK) cells, while knockout of the highest expressed migrasome-forming tetraspanins blocks this process⁵. In addition, tetraspanins and cholesterol form microdomains on migrasomal membranes. Reducing the cellular cholesterol level significantly impairs migrasome formation. Thus, tetraspanin and cholesterol are required for migrasome formation. To more accurately study the process of migrasome generation, we designed an in vitro system. Using purified tetraspanin 4-GFP (TSPAN4-GFP), cholesterol and other lipids, we generated giant unilamellar vesicles (GUVs). To mimic cell adhesion and migration, we first attached GUVs with bioengineered biotin to the bottom of a flow chamber coated with streptavidin. Then, by applying directional liquid flow in the chamber, we provided a mechanical force to move GUVs away from adherent spot. In this way, we successfully reconstituted the formation of migrasomes, and we found that both tetraspanin 4 and cholesterol colocalized on the reconstituted migrasomes. Without tetraspanin 4 or cholesterol, migrasomes could not be reconstituted; instead, small clusters of tetraspanin 4 or cholesterol were observed to be randomly distributed⁵. To study the dynamics of migrasome formation in vitro, we applied a direct pulling force using a glass needle attached to the GUV instead of directional liquid flow. By pulling the GUV away with a needle, we generated retraction fiber-like structures. In this way, we observed that the initially evenly-distributed small tetraspanin 4 clusters self-organized into macrodomains, which eventually swelled to form migrasomes. Thus, the mechanical force drove the formation of tetraspanin 4 macrodomains, which in turn shaped the migrasomes. We concluded that tetraspanin 4 and cholesterol are sufficient and necessary for migrasome formation⁵. Moreover, the *in vitro* reconstitution system provided us with an easily monitored platform to study migrasome formation with single components and made it possible to measure and calculate biophysical parameters. This made it possible to model migrasome formation. Our theoretical model explained that the shaping of migrasomes by tetraspanin 4 and cholesterol-enriched macrodomains may be explained by two biophysical properties: a high degree of bending stiffness of the migrasomal membrane and the line tension that acts at the boundary between the migrasome and the retraction fiber. The experimental parameters fit well with the calculated parameters from the modeling⁵. Thus, migrasome formation is mediated by assembly of micron-scale tetraspanin macrodomains. This work established tetraspanins as the key regulators of migrasome formation.

2. Physiological significance of migrasomes

2.1. Migrasomes are essential for zebrafish organ morphogenesis

The zebrafish embrvo is the first in vivo model system documented to have migrasomes⁶. We found that during zebrafish gastrulation, migrasomes are formed in the extracellular pockets of space between mesendodermal cells and in the pockets between the blastodermal margin and the yolk syncytial layer. Tetraspanin 4a, tetraspanin 7, and integrin β 1b were found to regulate migrasome biogenesis in zebrafish. Knockout of tetraspanins 4a and 7 resulted in organ laterality defects including left-right reversal and bilateral duplication. Injection of exogenous migrasomes partially rescues organ morphogenesis in tspan4a and tspan7 mutants⁶. Quantitative mass-spectrometry analysis showed that signaling molecules, including chemokines, morphogens, growth factors and cytokines, are enriched in migrasomes. Among these, Cxcl12 plays an important role in organ morphogenesis. We found that migrasomes were enriched around the embryonic shield and provided a source of Cxcl12a. The regional chemoattractants provided by migrasomes attract and hold dorsal forerunner cells (DFCs) at the front of the embryonic shield. The correct migration of DFCs ensures the formation of Kupffer's vesicle (KV), which is essential in establishing the left-right body axis⁶. The study of migrasomes in zebrafish embryonic development provides a new mechanism for establishing signaling cues, in which signaling molecules are packed into membrane-bound compartments for release. The spatial and temporal control of signaling release adds another layer of regulation to the coordination of embryonic development².

2.2. Identification of migrasomes in human serum through specific protein markers

We employed tandem mass tag (TMT) labeling followed by quantitative mass spectrometry to identify proteins enriched on migrasomes. By comparing the mass spectrometry data from migrasomes and exosomes, we identified and verified a set of specific protein markers that are enriched on migrasomes, but not on exosomes. These markers allowed us to analyze human serum samples biochemically⁷. We fractionated the serum and found that a certain fraction is positive for all the migrasome markers identified. Further electron microscopy analysis confirmed the presence of migrasomes in the fraction. Thus, migrasomes are present in human serum and can be detected by western blotting using a set of protein markers⁷. The identification of migrasome-specific protein markers will potentially be important in disease-related studies.

3. Future perspectives on migrasome research

Migrasomes are emerging as a new field in cell biology. So far, we have dissected the role of integrins, TSPAN4 and cholesterol

in migrasome biogenesis. Moreover, we demonstrated the essential role of migrasomes in organ morphogenesis in zebrafish embryonic development, which is the first example of the physiological significance of migrasomes. Furthermore, migrasomes were detected in human serum, which suggests their potential as vehicles for signal transduction in the circulatory system. We have already developed protein markers and a specific dye, WGA, for migrasome detection by imaging and biochemical approaches⁸. However, we are still at an early stage of migrasome research. The molecular mechanisms involved in migrasome biogenesis, for example the detailed regulatory mechanisms of migrasome biogenesis, the mechanisms for sorting the enriched contents of migrasomes, and the mechanisms of migrasome uptake by recipient cells, await dissection. Moreover, we are still far from fully appreciating the biological significance of migrasomes. More extensive research to establish a deeper understanding the physiological and pathophysiological relevance of migrasomes is urgently needed (Figure 2). One direction is to establish disease models in migrasome-deficient mice to study the different phenotypes caused by defective migrasome biogenesis. On the other hand, multi-omics approaches, including proteomics, transcriptomics, lipidomics, metabolomics, and post-translationomics analysis of human biofluid samples - for example, serum samples from large disease cohorts accompanied by data mining - is another way to unravel the link between migrasomes and disease. Last but not least, more accurate migrasome detection methods are required for both mechanistic research and translational research in the migrasome field. Development of a highly specific, easily manipulated dye with robust signals will be very much appreciated.

Migrasomes, which are generated through migracytosis, can be considered as signal vehicles for several core biological processes, including the regulation of immune responses, embryonic cell migration, and establishment of the tumor microenvironment.



Figure 2. Potential roles of migrasomes in normal physiology and tumor pathogenesis.

[References]

- L. Ma, Y. Li, J. Peng, D. Wu, X. Zhao, Y. Cui, L. Chen, X. Yan, Y. Du and L. Yu, *Cell Res.*, **2015**, 25(1), 24-38.
- 2. B. da Rocha-Azevedo and S. L. Schmid, Cell Res., 2015, 25(1), 1-2.
- D. Wu, Y. Xu, T. Ding, Y. Zu, C. Yang and L. Yu, *Cell Res.*, 2017, 27(11), 1397-1400.
- 4. Y. Chen, Y. Li, L. Ma and L. Yu, Methods Mol. Biol., 2018, 1749, 43-49.
- Y. Huang, B. Zucker, S. Zhang, S. Elias, Y. Zhu, H. Chen, T. Ding, Y. Li, Y. Sun, J. Lou, M. M. Kozlov and L. Yu, *Nat. Cell Biol.*, **2019**, *21*(8), 991-1002.
- D. Jiang, Z. Jiang, D. Lu, X. Wang, H. Liang, J. Zhang, Y. Meng, Y. Li, D. Wu, Y. Huang, Y. Chen, H. Deng, Q. Wu, J. Xiong, A. Meng and L. Yu, *Nat. Cell Biol.*, **2019**, *21*(8), 966-977.
- X. Zhao, Y. Lei, J. Zheng, J. Peng, Y. Li, L. Yu and Y. Chen, *Cell Discov.*, 2019, 5, 27.
- 8. L. Chen, L. Ma and L. Yu, Cell Discov., 2019, 5, 13.

[Contact]

Corresponding author:

Li Yu (俞立), School of Life Sciences, Tsinghua University, Beijing 100084, China. Tel: +86-(0)10-62792880 Email: liyulab@mail.tsinghua.edu.cn Graduation: Peking University: Ph.D Current position: Professor in the School of Life Sciences, Tsinghua University

Current interests: Migrasomes, a new organelle discovered in the Yu lab

First author:

Yang Chen (陈扬), Center for Precision Medicine Multi-Omics Research, Peking University Health Science Center, Beijing 100191, China Tel: +86-(0)10-82805521 Email: chenyang1816185048@bjmu.edu.cn Graduation: Peking University: Ph.D Current position: Tenure Track Investigator in the Health Science Center, Peking University Current interests: Multi-omics of extracellular vesicles

Second author:

Dongju Wang (王东锔), School of Life Sciences, Tsinghua University, Beijing 100084, China. Tel: +86-(0)10-62794552 Email: wangdj16@mails.tsinghua.edu.cn Graduation: Tsinghua University: B.S. Current position: Ph.D candidate in the School of Life Sciences, Tsinghua University Current interests: Migrasomes, a new organelle discovered in the Yu lab

新製品

エクソソーム蛍光染色キット

ExoSparkler Exosome

Membrane Labeling Kit-Green, Red, Deep Red Protein Labeling Kit-Green, Red, Deep Red

<特長>

- ・色素自身が凝集しない
- ・未反応色素を除くフィルトレーションチューブ同梱

ExoSparkler シリーズは、精製されたエクソソームの膜または タンパク質を染色し、細胞に取り込まれるエク 既存製品との比較

ソソームをイメージングすることができます。 また、一般的なエクソソーム染色試薬の課題 である色素自体の凝集を起こさず、染色前後で のエクソソームのサイズ分布や表面電荷状態を 変化させないため、より正確にエクソソームの 動態を可視化できます。







<時間依存的なエクソソームの局在変化>

超遠心法で精製したエクソソームを Membrane Labeling Kit-Deep Red(メーカーコード:EX03)で染色し、リソソーム染色 試薬で染色した HeLa 細胞に添加、1、4 時間後の蛍光画像を観 察しました。



<検出条件> EX03(紫): Ex 640 nm / Em 640-760 nm リソソーム染色試薬(緑): Ex 488 nm / Em 490-540 nm

その結果、時間経過に伴い EX03 の蛍光輝点(紫) はリソソームの局在(緑)と重なり(白)、時間依存的にエクソソームの局在が変化している様子が確認されました。

品名	容量*	希望納入価格(¥)	メーカーコード
ExoSparkler Exosome Membrane			
Labeling Kit -Green	5 samples	25,000	EX01
-Red	5 samples	25,000	EX02
-Deep Red	5 samples	25,000	EX03
ExoSparkler Exosome Protein			
Labeling Kit-Green	5 samples	20,000	EX04
-Red	5 samples	20,000	EX05
-Deep Red	5 samples	20,000	EX06

[※]精製済エクソソーム(超遠心法)として、

タンパク質: 1-10 μg/sample、粒子数: 10-100 × 10⁸ 個 /sample)

新製品

細胞周期測定試薬

Cell Cycle Assay Solution Blue Cell Cycle Assay Solution Deep Red

<特長>

- ・細胞の固定化、RNase 処理が不要で試薬を添加するだけ
- ・405 nm、633 nm のレーザーで励起が可能

一般的に使用される Propidium lodide (PI)を用いた手法と比較して、本製品は高い細胞膜透過性と DNA 選択性を有する色素を使用しているため、細胞懸濁液に試薬を添加するだけで測定が可能です。



Cell Cycle Assay Solution Deep Red	50 tests	13,000	C548
Cell Cycle Assay Solution Blue	50 tests	13,000	C549

細胞老化誘導に伴う細胞周期と ミトコンドリア膜電位の変化

細胞周期の G2/M 期に作用し細胞増殖を停止させ、細胞老化 を誘導することが知られる Doxorubicin (DOX) を A549 細胞へ添 加後、本製品で細胞周期の変化と、細胞老化 [Cellular Senescence Detection Kit - SPiDER- β Gal (メーカーコード: SG03)] およびミトコンドリア膜電位の低下 [JC-1 MitoMP Detection Kit (メーカーコード: MT09)] を確認しました。



新製品

脂肪滴染色試薬・キット

Lipi シリーズは、脂肪親和性の高い低分子蛍光試薬であり、疎水性環境下で蛍光が増強します。また、試薬を添加するだけで生細胞および固定化細胞中の脂肪滴を明瞭に観察することができます。

Lipi シリーズを用いて脂肪滴の量的変動を数値化できるキットもご用意しております。

蛍光顕微鏡用 Lipi-Blue, Green, Red, Deep Red

<測定例>

オレイン酸を添加した HeLa 細胞を生細胞の状態で、Lipi シリーズの各色素にて染色しました。



フローサイトメーター、プレートリーダー用 Lipid Droplet Assay Kit - Blue, Deep Red

<測定例>

細胞にオレイン酸または脂肪滴の形成阻害剤である Triacsin C を添加し、Lipid Droplet Assay Kit を用いて脂肪滴の量的変動を数値 化しました。その結果、コントロールならびに Triacsin C を加えた細胞と比較して、オレイン酸を添加した細胞で脂肪滴の量が増加 していることを確認しました。



Deep Red : Ex 623 - 633 nm、Em 649 - 669 nm

品名	容量	希望納入価格(¥)	メーカーコード
Lipi-Blue	10 nmol * 1	18,000	LD01
Lipi-Green	10 nmol * 1	18,000	LD02
Lipi-Red	100 nmol * 1	18,000	LD03
Lipi-Deep Red	10 nmol * 1	18,000	LD04
Lipid Droplet Assay Kit-Blue	1 set * 2	25,000	LD05
Lipid Droplet Assay Kit-Deep Red	1 set * 2	25,000	LD06

* 1 Lipi シリーズ色素: 35 mm dish 10 ~ 50 枚分

* 2 Assay Kit: 96 well plate 1 枚分、フローサイトメトリー 40 アッセイ分



細胞:HeLa

Blue \therefore Ex 405 nm $\$ Em 425 - 475 nm Deep Red \therefore Ex 640 nm $\$ Em 650 - 670 nm

Lipi シリーズを使用した論文は 小社 HP よりご覧いただけます。



Topics on Chemistry

二重機能性蛍光色素(H-V)を用いたフェロトーシスの解明

フェロトーシスは、アポトーシスともネクローシスとも異なる 経路で起こる新たな細胞死である。2012年にStockwellらは、 鉄依存的な細胞死としてフェロトーシスを提唱し、その経路とし て活性酸素種(Reactive Oxygen Species: ROS)と二価鉄の反 応により脂質の過酸化が増加することで細胞死へと誘導されるこ とを報告している(図1)¹⁾。しかしながら、現在フェロトーシ スの実行因子やその機構についてはまだ不明な点が多く、その解 明のため非常に活発な研究が行われている。

本稿では、Liらが発見した ROS の一つであるヒドロキシラジ カル及び細胞質の粘度変化に伴うフェロトーシスの研究について 紹介する²。

Li らは、細胞内のヒドロキシラジカル並びに細胞質内の粘度変 化を可視化するため、アニソールにビニル基を介しインドリウム 誘導体を導入した低分子蛍光プローブ(H-V)を開発した。H-V は、細胞質の粘度上昇依存的に緑色蛍光(Am = 520 nm)が増 加し、一方でヒドロキシラジカルにより酸化を受けた際には赤色 蛍光(λ_{em} = 652 nm)を発する特徴も持つ(図 2)。グリセロー ルを含むメタノールを試料として用いた In vitro 実験において、 H-V は検体の粘度増加に伴い約 50 倍の蛍光増大を示し、フェン トン反応により生成させたヒドロキシラジカルを約450倍の高 感度で検出した。また Li らは、HT-1080 細胞にエラスチンを用 いてフェロトーシスを誘導すると、細胞質の粘度上昇に伴う 520 nm 蛍光増加(約7倍)と、ヒドロキシラジカル産生に伴う 652 nm 蛍光増加(約26倍)が起きることを確認している。この結 果は、フェロトーシスにおいてヒドロキシラジカルの生成と共に 細胞質の粘度上昇が生じることを示している。興味深いことに、 細胞質内粘度の増加に伴い脂肪滴の形成が加速されることも脂肪 滴染色試薬を用いた実験で示唆されている。これは、鶴崎らが報 告した NASH(非アルコール性脂肪肝炎)とフェロトーシスの関 係性にも通ずる点である³⁾。

今回は、Liらの研究成果をもとにフェロトーシスを紹介した が、前述したようにまだまだ未知な点が多い研究分野である。よ り詳細な細胞内鉄(総量/二価鉄/三価鉄)の動態や量変化を検 出できる技術が開発されることで更なる研究の発展が期待される。

[参考文献]

- 1) S. J. Dixon *et al.*, *Cell*, **2012**, *149*, 1060.
- 2) H. Li et al., J. Am. Chem. Soc., 2019, 141, 18301.
- 3) S. Tsurusaki et al., Cell Death and Disease, 2019, 10, 449.



図1 フェロトーシスの分子機構

株式会社同仁化学研究所 北村 怜奈



図2 ヒドロキシラジカルと粘度に対する H-V の反応メカニズム

細胞内鉄検出試薬

FerroOrange

く特長>

- ・生細胞中の Fe²⁺を高感度に検出
- ・96 穴マイクロプレートアッセイに対応

本製品は、細胞内の二価鉄 (Fe²⁺) を選択的に検出するための 蛍光プローブです。培養細胞に添加するだけで細胞膜を透過し、 細胞内の Fe²⁺と反応して強い蛍光 (λ_{ex} :543 nm、 λ_{em} :580 nm)を発します。

<検出例>

細胞内に存在する Fe²⁺ を確認するため、鉄キレート試薬 Bpy (2,2'-bipyridine) または鉄(硫酸アンモニウム鉄(Ⅱ))を添加し た HeLa 細胞を用い、細胞内の Fe²⁺ を検出しました。

イメージング解析



マイクロプレートアッセイによる数値化



3 tubes

32,000

※ 1 tube あたり 24μg の FerroOrange が含まれます

- 関連アプリケーションの紹介-

フェロトーシス研究

アミノ酸トランスポーター(xCT)阻害

Erastin 刺激により、シスチン / グルタミン酸交換輸送体(xCT)を阻害すると、細胞内へのシスチンの取り込みおよび、細胞外へのグルタミン酸の放出が阻害されます。シスチンの細胞内取り込みが阻害されることにより、グルタチオンの生成が低下し、さらに、鉄に依存した脂質過酸化によるフェロトーシスが引き起こされます。

<実験例>

xCT が阻害された A549 細胞において、Glutamate 放出量および細胞内 GSH が減少し過酸化脂質が増大する結果が得られました。



細胞内過酸化脂質検出試薬

Liperfluo

<特長>

- ・細胞内の過酸化脂質の検出ができる
- ・過酸化脂質特異性が高い

Liperfluo は生細胞の過酸化脂質のイメージングやフローサイトメトリーによる過酸化脂質量の分析に使用することができます。

<測定例>

生細胞を用いたライブセルイメージング



装置: Zeiss LSM510META フィルター: FITC (GFP, Alexa488) wide filter HFT UV/488、NFT490、 BP505-550

細胞:L929

フェロトーシス研究にて Liperfluo を用いた 論文を続々公開! 詳細は小社 HP をご覧ください。

	L248	同仁	検索		
品名		容量	希望	望納入価格(¥)	メーカーコード
Liperfluo		50 μg × 9	5	21,200	L248

グルタチオン測定キット

GSSG/GSH Quantification Kit

<特長>

- ・酸化型(GSSG)、還元型(GSH) グルタチオンの分別定量が 可能
- ・比色によるマイクロプレートアッセイで多検体の測定が可能

キットには、GSH のマスキング剤が含まれており、酵素リサ イクリング法を用いた DTNB [5,5'-dithiobis (2-nitrobenzoic acid)] 発色 (λ max = 412 nm)を測定することで、GSSG の みを定量でき、別途測定した総グルタチオン量から GSSG 量を 差し引くことで GSH 量を求めることができます。



③酵素溶液および基質を

加える



② Buffer solution を加え37℃
 1時間インキュベート

	-	-			1
1	1		in the second	1	
2	VE.		12	1.4	
h.,	1			月	۰.
			-	4	1

 ④ 37℃で 10 分間インキュベー ト後 412 nm の吸光度を測定

品名	容量	希望納入価格(¥)	メーカーコード
GSSG/GSH Quantification Kit	200 tests	54,800	G257

ミトコンドリア研究用試薬



品名	容量	希望納入価格(¥)	メーカーコード
Mitophagy Detection Kit	1 set	36,000	MD01
Mito-FerroGreen	50 μ g $ imes$ 2	25,000	M489
MitoPeDPP	$5\mu{ m g} imes3$	19,200	M466
Si-DMA for Mitochondrial Singlet Oxygen Imaging	2µg	20,600	MT05
Lastata Assaul Kit WST	50 tests	29,000	1.056
	200 tests	68,000	L200
Chucoso Accov Kit WST	50 tests	18,000	C264
Glucose Assay Rit-W31	200 tests	38,000	G204
JC-1 MitoMP Detection Kit	1 set	23,000	MT09

新製品

DOIINDO MOLECULAR TECHNOLOGIES, INC.

ミトコンドリア染色試薬

MitoBright LT Green, Red, Deep Red

<特長>

- ・ミトコンドリアを長期間観察可能
- ・血清入り培地中で染色可能

・コラーゲンコートプレート上で培養した細胞のミトコンドリアを低バックグラウンドで検出可能

ミトコンドリアを長期間観察

HeLa 細胞を各 MitoBright LT または既存試薬で染色し、4 日間培養後ミトコンドリアを観察しました。その結果、既存試薬の蛍光 強度は大きく低下したのに対し、MitoBright LT は蛍光強度が維持され、ミトコンドリアを明瞭に観察できました。



タイムラプスイメージング

HeLa 細胞を CCCP 処理し、マイトファジー検出 試薬 (MtphagyDye メーカーコード: MT02) 及びミトコンドリア染色 試薬(MitoBright LT Green)にて共染色しタイムラプス撮影(6 時間)しました。詳細は HP にてご覧いただけます。

マイトファジーのリアルタイム観察

< 検出条件> ■ 緑:MitoBright LT Green



装置: LSM-700 Laser scanning confocal microscope (LSCM) (Carl Zeiss, Oberkochen, Germany) 励起波長: MitoBright LT Green 488 nm MtphagyDye 555 nm 対物レンズ:63倍 拡大:1.5倍 撮影インターバル:15秒



データ提供 タイムラプスビジョン

品名	容量	希望納入価格(¥)	メーカーコード
	20 μ I $ imes$ 1	6,000	
MitoBright LT Green	400 μ I $ imes$ 1	12,000	MT10
	400 μ I $ imes$ 3	30,000	
	20 μ I $ imes$ 1	6,000	
MitoBright LT Red	400 μ I $ imes$ 1	12,000	MT11
	400 μ I $ imes$ 3	30,000	
	20 μ I $ imes$ 1	6,000	
MitoBright LT Deep Red	400 μ I $ imes$ 1	12,000	MT12
	400 μ I $ imes$ 3	30,000	

よくある質問

- Q. MitoBright LT で染色した後、脱分極をすると染色の度合いに 影響はありますか?
- A. 脱分極条件や細胞種などにより影響は確認されます。また、 MitoBright LT の各色素で影響の度合は異なります。参考とし て小社では HeLa 細胞を各 MitoBright LT で染色し、下記条件 にて脱分極処理を行って蛍光染色の変化を観察しました。



<染色条件> HeLa 細胞を MitoBright LT(100 µmol/1、30 分インキュベーショ . ン)にて染色し HBSS にて洗浄 後 FCCP 処理(100 µmol/I、60 分 インキュベーション) し HBSS にて2回洗浄後観察した。

<検出条件> MitoBright LT Green : Ex 488 nm/Em 500-560 nm MitoBright LT Red : Ex 561 nm/Em 560-620 nm MitoBright LT Deep Red : Ex 640 nm/Em 650-700 nm

さらに詳細な情報は小社 HP で! 図各試薬の蛍光特性 ☑タイムラプス動画 ☑フローサイトメーターの検出例

MitoBright LT 同仁

検索

細胞内代謝測定キット

細胞内の代謝システムである解糖系や TCA 回路、ペントース - リン酸経路の解析は、細胞状態を理解する上で重要であり、グルコース、乳酸、NAD(P)+/NAD(P)H、グルタミン、グルタミン酸などの代謝産物を指標に評価されています。





1 キットあたり測定可能なサンプル数 (n = 3 で標準サンプル 8 点で検量線を作成した場合) グルタミン、NAD(P) / NAD (P) H: 12 サンプル グルタミン酸: 24 サンプル

品名	容量	希望納入価格(¥)	メーカーコード
Glutamine Assay Kit-WST	100 tests	55,000	G268
Glutamate Assay Kit-WST	100 tests	50,000	G269
Glucosa Assay Kit WST	50 tests	18,000	6264
Glucose Assay Kit-WST	200 tests	38,000	6204
Lactato Assav Kit-W/ST	50 tests	29,000	1 256
	200 tests	68,000	L250
NAD/NADH Assay Kit-WST	100 tests	54,000	N509
NADP/NADPH Assay Kit-WST	100 tests	54,000	N510

Antimycin

実験例

ミトコンドリア電子伝達系の阻害実験

Antimycin 刺激によりミトコンドリア電子伝達系を阻害すると、解糖系を強制的に維持しようとしピルビン酸から乳酸への代謝が亢 進し、NAD/NADH 比が低下します。また電子伝達系の阻害では、ミトコンドリア膜電位が低下し、活性酸素種(ROS)の増加に伴う GSH の枯渇を引き起こすことが知られています。

下記の実験では、Jurkat 細胞に Antimycin 50 μmol/l を添加し 24 時間インキュベート後の培養上清または細胞を用いて小社キット により測定を行いました。



抗酸化能測定キット -DPPH ラジカル消去活性

DPPH Antioxidant Assay Kit

<特長>

- ・必要な試薬が揃ったキット
- ・試薬調製の手間を大幅に削減
- ・マイクロプレートアッセイによる多検体分析が可能

高知大学の島村らは DPPH(2,2-Diphenyl-1-picrylhydrazyl)を 用いた施設間差の少ない抗酸化活性測定法を報告しています¹⁾。 本製品は島村らの方法をマイクロプレートアッセイに適用し、多 検体分析を可能にしました。また ready-to-use 仕様であるため 試薬調製の煩雑さを低減できるとともに、ヒト間および施設間に よるデータのバラツキを抑えることが可能になりました。

1) T. Shimamura et al., Anal. Sci., 2014, 30, 717-721.

操作の手間を大幅に削減



試薬を添加するだけの簡便な操作





 ①標準物質(Troiox)又はサンプル
 ②バッファー及び発色試薬(DPPH) を添加
 を添加



③ 25℃で 30 分間インキュベート

④マイクロプレートリーダーで測定 (517 nm)

本キットを使用した論文が公開されました。 詳細は小社 HP をご覧ください。

DPPH	同仁	検索

よくある問い合わせは P13 をご覧ください。

品名	容量	希望納入価格(¥)	メーカーコード
DPPH Antioxidant Assay Kit	100 tests	6,400	D679
	500 tests	19,000	D070

*本品1セットあたり、100 tests:1-3サンプル、500 tests: 8-15サンプル測定できます。(n = 3、8段階希釈の場合)

DOJINDO MOLECULAR TECHNOLOGIES, INC.

抗酸化能測定キット-SOD 様活性

SOD Assay Kit-WST

<特長>

- ・試薬添加だけの簡便な操作
- ・マイクロプレートアッセイによる多検体処理が可能

本キットでは、活性酸素種の一つであるスーパーオキシドの消 去能を持つ SOD(superoxide dismutase)様活性を測定するこ とができます。

<操作>

96 穴マイクロプレートを用い、サンプル調製から測定まで約 1 時間で完了します。



サンフル調製



高知大学の島村らは碁 石茶(ゴイシチャ)の製 造工程毎に SOD 様活性 をキットで測定していま す。 参照元:島村,日本食品科学 工学会誌,55(12),640.

本キットを使用した研究報告を小冊子に掲載しております。

	はじめてプロトコル	同仁	検	索		
	品名	容量		希望納	入価格(¥)	メーカーコード
	anav Kit MCT	100	tests		8,800	6211
SOD Assay NI-WST		500 tests		2	2,900	5311



FAQ

DPPH Antioxidant Assay Kit

- Q. <u>このキットでどのようなことができますか?</u>
- A. 農作物や食品、物質のスクリーニング測定を行うことで、機 能性のある農産物や食品、物質などを見つけ出すことができ ます。
- Q. 本キットを使用する利点を教えてください。
- A. 本キットは、必要な試薬が揃っており、課題であったサンプ ル溶液の pH や有機溶媒の濃度によるデータのばらつきの影 響を最小限に抑えるためのプロトコルと解析法を採用してい ます。また、時間と手間を大幅に削減することができます。
- Q. DPPH ラジカル消去活性だけで抗酸化活性を評価していいで しょうか?
- A. 複数のサンプルの抗酸化活性を比較するには、同一の方法で 測定する必要があり、他の方法で測定した値と比較すること はできません。また、他の方法で測定すると同じサンプルで も抗酸化活性の程度が異なる場合があります。幾つかの測定 方法で抗酸化活性を比較することをお勧めします。
- **Q**. どんなサンプルでもこのキットで測定できますか?
- A. 本キットはエタノールが反応溶媒の 50%程度を占めるため、 タンパク質を豊富に含む試料(例えば、牛乳など)はタンパ ク質の凝集を起こします。サンプルを 50%エタノール水溶液 に溶解し沈殿が生じる場合は、沈殿が生じない濃度まで希釈 してください。希釈できない場合は、反応溶媒が水系である SOD Assay Kit-WST での SOD 様活性での評価をおすすめし ます。

詳細は、右の選択ガイドを参照下さい。

- Q. DPPH Antioxidant Assay Kitの測定実績を教えてください。
- A. 農作物や食品抽出物、食品添加物などの測定例があります。 前処理方法等は HP に掲載しております。







ホームページアドレス URL:https://www.dojindo.co.jp/ E-mail:info@dojindo.co.jp



0120-021557 0120-489548 (受付時間:平日9:00~17:00 土・日・祝日は除く)



ドージンニュース No.172 令和2年3月6日発行 株式会社同仁化学研究所 DOJINDO LABORATORIES 熊本県上益城郡益城町田原2025-5 〒861-2202 発行責任者 松野寛明 編集責任者 金子詩織 年4回発行 許可なくコピーを禁ず