



オルガネラゾーン - 総 説-小胞体-ミトコンドリア接触領域の形成因子と 細胞機能 ^{熊本大学 山本 真寿・山口 知也} - 注目の研究--低分子蛍光色素によるミトコンドリア-リソソーム間相互作用の観察

株式会社同仁化学研究所 田尻 智巳

新製品 細胞膜染色試薬 🛤

エンドサイトーシス検出試薬 P3 エクソソーム染色キット 210 細菌二重染色キット 212 バイオフィルム形成阻害測定キット・ バイオフィルム薬剤効果測定キット 213



表紙撮影:熊本県宇土市住吉町 海苔の種付け 毎年10月末に海苔の種付けが行わ れ、漁業者によって異なる赤や青、 緑色などのノリ網が浮かび上がり、 海面を鮮やかに彩っている。 photo:永島俊介氏

CONTENTS

Review

小胞体-ミトコンドリア接触領域の形成因子と細胞機能 • [1] 熊本大学 山本 真寿 · 山口 知也

Topics on Chemistry

低分子蛍光色素による ミトコンドリア-リソソーム間相互作用の観察 • [6] 株式会社同仁化学研究所 田尻 智巳

Commercial

ミトコンドリア染色試薬 ●	[5]
微生物増殖アッセイキット ●	[13]

新製品

細胞膜染色試薬 ● [8]
エンドサイトーシス検出試薬 ● [9]
エクソソーム染色キット ● [10]
細菌二重染色キット ● [12]
バイオフィルム形成阻害測定キット・バイオフィルム薬剤効果測定キット ● [13]

お役立ち 用途で選ぶ細胞膜関連試薬 ●────[7] フェロトーシスの検出 ●───[11]

フォーラム・イン・ドージン延期のご連絡 2020年11月に開催を予定しておりました第31回フォーラム・イン・ドージンは、 新型コロナウイルス感染症(COVID-19)の感染拡大の防止という観点から、開催の 延期を決定いたしましたことをお知らせいたします。 今後の日程につきましては、決まり次第ホームページ等でご案内いたします。 何卒ご理解をいただきますよう、よろしくお願いいたします。

※希望納入価格には消費税等は含まれておりません。



小胞体-ミトコンドリア接触領域の形成因子と細胞機能 The composition and cellular function of ER-mitochondria contact sites



山本 真寿

熊本大学 大学院生命科学研究部 がん生物学講座 助教



山口知也

熊本大学 大学院生命科学研究部 がん生物学講座 准教授

Abstract

Eukaryotic cells contain various organelles which are compartmentalized by the intracellular membrane to maintain their identity and specialize biological process. These organelles are not only individual entities but also communicated between other organelles to perform their appropriate function. Such inter-organelle communication is achieved by canonical vesicular trafficking and physical associations between multiple organellar membranes. Here, we briefly review recent findings on the composition of ER-mitochondria contact sites and their cellular function.

1. はじめに

真核細胞内のオルガネラは、生体膜や相分離によって細胞質か ら区切られ個々の機能に特化した組成・構造を維持している細胞 内コンパートメントである。このことは、オルガネラ間の物質輸 送には大きな物理化学的障壁が存在することを意味しており、効 率的な物質輸送には自由拡散に依存しない物質輸送経路が必要で あると考えられる。このような細胞内輸送経路として「小胞輸送 経路」が知られているが、ミトコンドリアなど明確な小胞輸送経 路を持たないオルガネラの存在から、それ以外の経路の存在が考 えられてきた。この小胞輸送に依存しない輸送経路のひとつとし て考えられているのが「オルガネラ接触領域」を介した物質輸送 が近年注目されている。二つのオルガネラが物理的に近接した接 触領域では、イオンなど親水性分子が効率的に輸送されるだけで なく、リン脂質の様な疎水性分子も接触領域の繋留タンパク質を 介して輸送されていると考えられている。さらに近年、オルガネ ラ接触領域は物質輸送だけでなく、オルガネラの形成やオルガネ ラの形態・動態を制御していることも明らかになってきた。本稿 ではこれらオルガネラ接触領域の中でも特に研究の進んでいる 「小胞体-ミトコンドリア接触領域」に注目し、現在までに明らか にされた形成機構や生理機能について概説したい。

2. オルガネラ接触領域

オルガネラ接触領域の定義については議論が進んでいるものの ¹⁾、厳密な定義は定まっていない。現時点における多くの研究者 の認識は「繋留タンパク質を介して2つのオルガネラ膜が10~ 30 nm 程度に近接した領域」といえるだろう。オルガネラ接触 領域は「接触」や「Contact」という言葉が用いられているが、 オルガネラ同士の接触は繋留タンパク質を介した相互作用に限定



図1 細胞内の小胞体接触領域とミトコンドリア接触領域

されており、オルガネラ膜の融合は通常見られない。このことは 2つのオルガネラが混ざり合うのではなく、個を維持したまま物 理的に近接し、効率的に物質の輸送を行う「ゾーン」を形成する ことに接触領域の生理的な意義があるものと考えられる。

近年、細胞膜を含め様々なオルガネラ膜間に接触領域が存在す ることが相次いで報告されている。その多くは小胞体と他のオル ガネラとの接触であり、細胞内の膜を有するオルガネラとの接触 のほか²⁾⁻⁵⁾、ストレス顆粒やP体(processing body)など相分 離によって形成される膜を持たないオルガネラとの接触も報告さ れている⁶⁾(図1左)。このことは小胞体が膜を構成するリン脂 質分子を合成する場であることから、様々なオルガネラへ生体膜 の材料となる脂質分子を効率的に輸送することを反映しているの かもしれない。また、ミトコンドリアも小胞体^{71,8)}やペルオキ シソーム⁹⁾、エンドソーム¹⁰⁾、脂肪滴¹¹⁾など様々なオルガネラ との接触が報告されており、これらは脂肪酸やトランスフェリン を介した鉄イオンのミトコンドリアへの取り込みなどに関与して いると考えられている(図1右)。

このように様々なオルガネラ接触領域が報告されるに伴い、そのオルガネラ接触領域を形成する因子群の同定も進められてきた。後述する小胞体-ミトコンドリア接触領域に関しては複数の 繁留分子や調節因子の関与が見出されており、構成分子の異なる 接触領域間での機能的な違いなどにも関心が向けられつつある。 その一方で、他のオルガネラ間接触領域の形成因子の同定はあま り進んでおらず、詳細がほとんど不明なままの接触領域もある。 これらの新たに発見された接触領域については、単離精製法の開 発と形成因子の同定、細胞・個体レベルでの生理的機能の解析が 待たれる。

3. 小胞体-ミトコンドリア接触領域 (ER-Mitochondria Contact Sites, EMCSs)

小胞体とミトコンドリアの接触領域は歴史的に最も早く見出された接触領域であり、1950年代にラット肝細胞の電子顕微鏡解析において小胞体膜とミトコンドリア外膜の近接が報告された¹²(図 2)。

しかし当初は、この小胞体とミトコンドリア膜同士の近接に何 らかの機能があるとはあまり考えられておらず、接触領域を単離 精製する技術的な困難もあり長らく形態学的な解析に留まってい た。転機となったのは 1990 年の Vance による報告であり、ラッ トの肝臓から精製した粗ミトコンドリア画分に小胞体を主成分と するミクロソーム画分が含まれていること、この画分を Percoll 密度勾配遠心によって単離精製できることを示した"。この画分 は現在 MAM (Mitochondria-Associated Membrane) 画分と呼ば れるミトコンドリアと近接した小胞体膜画分であり、この発見に より EMCSs の生化学的な解析が可能となった。この MAM 画分 にはリン脂質¹³⁾やトリアシルグリセロール¹⁴⁾、ステロール¹⁵⁾ など様々な脂質合成酵素群が多く含まれていることから、MAM はミトコンドリアと接触した滑面小胞体画分であり、小胞体から ミトコンドリアへの脂質輸送を担っていることが予想された。他 方、粗面小胞体も EMCSs を形成することが電子顕微鏡で観察さ れており16),17)、リボソームを含むためか膜間の近接距離は50 nm 以上と滑面小胞体の場合よりも大きいということはわかって いるが、機能については不明な点が多く残されている。



HeLa細胞を共焦点顕微鏡及び透過型電子顕微鏡にて撮影

図2 小胞体-ミトコンドリア接触領域 (EMCSs)

4. EMCSs の形成機構と細胞機能

細胞内に存在する EMCSs は細胞種により多少の違いがある が、小胞体とミトコンドリアのそれぞれ 10%程度が接触領域を 形成している場合が多い。ミトコンドリアを基準に見てみると、 HeLa 細胞では表面積の 10 ~ 20%が^{8),18)}、マウス肝臓では 4 ~ 10%程度が小胞体と近接していると報告されている¹⁹⁾。逆に小 胞体は全体の 15%程度がミトコンドリアと近接していることが HeLa 細胞において報告されている¹⁸⁾。このように細胞内では、 想像以上に多くの小胞体とミトコンドリアが接触領域を形成して おり、様々な細胞機能にとって重要な場となっていることが予想 される。

EMCSsの形成機構は依然として不明な点が多いが、2つの膜に存在するタンパク質間の相互作用によって形成されていると考えられている。これまでにイオンチャンネルやミトコンドリア融合に関与するタンパク質など、様々な分子が小胞体とミトコンドリア間の相互作用を担う形成因子として同定されている。哺乳動物ではこれら形成因子の1つを破壊・抑制してもEMCSsの個数や面積は部分的な減少に留まることが多いため、EMCSsは一様な構造ではなく、複数の繋留複合体によって形成された多様性のある構造であることが予想される。以下では細胞内においてEMCSsの関与が報告されている機能毎に形成分子を概説したい。

4.1. 小胞体-ミトコンドリア間のCa²⁺イオン輸送

細胞内における Ca²⁺イオンは非常に重要なセカンドメッセン ジャーであり、Ca²⁺イオン濃度の変動は細胞のエネルギー代謝 や細胞死、筋肉の収縮や神経伝達物質の放出など様々な生命現象 に重要である。小胞体とミトコンドリアはともに細胞内における Ca²⁺イオン貯蔵庫であり、刺激に応じて小胞体から放出された Ca²⁺イオンをミトコンドリアが取り込むことで局所的な ATP 合 成を促すとともに、過剰な Ca²⁺イオンの流入は細胞死を誘導す ることが知られている²⁰⁾。この小胞体からミトコンドリアへの Ca²⁺イオンの効率的な輸送が EMCSs の生理的意義の一つであ ると考えられている⁸⁾。

そのことを裏付けるように、小胞体から Ca²⁺イオンを放出す るチャネルであるイノシトール三リン酸受容体(IP3R)自体が、 ミトコンドリア外膜のチャネルである VDAC と細胞質の GRP75 を介して 2 つのオルガネラを繋留する複合体として機能するこ とが見出されている²¹⁾(図 3)。さらに IP3R のシャペロンとして 安定化に寄与する Sigma-1 受容体も MAM に濃縮されており ²²⁾、間接的に EMCSs の形成に寄与している。Mitofusin2 (MFN2) はミトコンドリアの融合に関与する分子であるが、EMCSsの形 成因子としても広く知られている²³⁾。MFN2 はその一部が小胞 体に局在し、ミトコンドリアに局在する MFN1 及び MFN2 と結 合し2つのオルガネラを繋留するとされている²⁴⁾。Mfn^{-/-}ノッ クアウトの MEF では EMCSs の間隔が拡がり、ミトコンドリア への Ca²⁺イオン流入の減少がみられたことから Ca²⁺イオンの 輸送を担う EMCSs の形成因子と考えられる^{24), 25)} (MFN2 に関 しては EMCSs の形成に負に作用するという相反する報告もあり ^{26), 27)}、今後の詳細な検証が待たれる)。このほかBAP31-FIS1²⁸⁾、VAPB-PTPIP51²⁹⁾、PDZD8¹⁸⁾など様々な分子が EMCSs の形成に関与し、ミトコンドリアへの Ca²⁺イオン流入を促進す ることが示されている。一方、EMCSsの形成に阻害的に働く因 子として FATE1 が知られている。FATE1 は EMD や MIC60 と結 合し小胞体とミトコンドリアの距離を拡げることにより EMCSs の形成を阻害することで、ミトコンドリアへの Ca²⁺イオン流入 を減弱させることが報告されている³⁰⁾。

このように小胞体とミトコンドリアの Ca²⁺イオン輸送能については、多くの形成因子の関与が報告されており、水溶性の小分子である Ca²⁺イオンの放出や吸収を制御する仕組みとして膜同士の物理的な近接は重要な役割を担っているのかもしれない。



図3 Ca²⁺輸送に関与する EMCSs 形成因子

4.2. 小胞体-ミトコンドリア間の脂質輸送

リン脂質合成は主に小胞体で行われるが、いくつかのリン脂質 合成酵素はミトコンドリアにも存在しており、例えばホスファチ ジルエタノールアミン(PE)の合成経路の一つであるホスファ チジルセリン脱炭酸酵素(PSD)はミトコンドリア内膜に局在し ている³¹⁾。そのため小胞体で合成したホスファチジルセリン (PS)をミトコンドリアへと輸送し、ミトコンドリアで合成した PEを小胞体へ送り返していることになるが、前述の通り小胞体 とミトコンドリア間には小胞輸送経路がなく、EMCSsを介して 脂質輸送が行われていることが予想されていた⁷⁾。

その分子メカニズムの一端が明らかとなったのは、小胞体-ミ トコンドリアを人工的に繋留するリンカー分子を利用した出芽酵 母の遺伝学的スクリーニングにより、ERMES (ER-Mitochondria Encounter Structure) と呼ばれるタンパク質複合体が EMCSs の 繋留因子として同定されたことによる³²⁾。ERMES 複合体の欠損 株では、PS → PE → PC という変換反応が有意に減少すること が報告されており³²⁾、in vitro においても ERMES 複合体は、 ATP などのエネルギーを必要とせずに 2 つのオルガネラ膜間の 脂質輸送を担うことが明らかにされている³³⁾。これら ERMES 複合体を構成する4つのタンパク質のうち3つ(Mdm12, Mmm1, Mdm34)はSMP(Synaptotagmin-like Mitochondrial lipid-binding Profein)ドメインを有しており、疎水性度が高いポケットにリ ン脂質の炭化水素鎖が結合することが結晶構造解析から推定され ている^{34),35)}。このことからリン脂質はERMES 複合体によって 構成された疎水性環境を介して2つのオルガネラ間を行き来し ているものと考えられている(図4)。

ERMES 複合体は長らく哺乳動物など高等真核生物でのホモロ グ分子が見つかっていなかったが、近年、ERMES 複合体の1つ である Mmm1の哺乳類ホモログ分子として PDZD8 が見出され た¹⁸⁾。PDZD8 は EMCSs の形成と Ca²⁺イオン輸送能への関与 は明らかにされているものの¹⁸⁾、脂質輸送能については明らか ではなく、他の ERMES コンポーネントのホモログ分子の同定な どを含めて今後の研究が待たれる。

また、酵母から哺乳類まで遺伝的に保存された EMC(ERmembrane protein complex) 複合体も EMCSs における脂質輸 送への関与が報告されている^{36),37)}。しかし、EMC 複合体には脂 質輸送に関与するドメインを持つ構成タンパク質がないことか ら、その詳細なメカニズムについては不明な点が多く残されてい る(図 4)。



図4 リン脂質輸送に関する EMCSs 形成因子

4.3. ミトコンドリアダイナミクス

ミトコンドリアは融合と分裂を繰り返す動的なオルガネラであ るが、チューブ状の小胞体がミトコンドリアに絡みつくようにし て収縮させることで、ミトコンドリアを分裂させていることが報 告された³⁸⁾。このミトコンドリア分裂開始点に一過的に形成さ れる EMCSs では、まず小胞体に局在する Formin タンパク質の 一種 INF2 が集積し³⁹⁾、ミトコンドリアに存在する SPIRE1C と 結合することでアクチンの重合が促進されることが分かってきた ⁴⁰⁾。このアクチン繊維が収縮力となり、ミトコンドリアがくびれ た部分にダイナミン様タンパク質の DRP1 や DYN2 が重合する ことで、最終的にミトコンドリアが分裂するというモデルが考え られている^{38),41)}。

4.4. その他の機能

小胞体とミトコンドリアは ER ストレスや ATP 合成などの反応時に ROS を産生することが知られている。FRB-FKBP 二量体化法を用いて ROS センサーを EMCSs に標的化したところ、EMCSs 周辺は H₂O₂ が局所的に高濃度になっていることが見出

された⁴²⁾。これは EMCSs において Ca²⁺が小胞体からミトコン ドリアへ流入することが引き金となり、ミトコンドリアに蓄積された H₂O₂ が細胞質へ放出されることによって形成される⁴²⁾。またこのようにして EMCSs 周辺に形成された酸化的環境は、局所的に p38 α を活性化し、MIRO/TRAK 複合体を介したミトコンドリアの移動を阻害することが報告されている⁴³⁾。これによりミトコンドリアを EMCSs に留めることで、EMCSs を安定化させるという仮説が提唱されている。

さらに EMCSs はオートファジーとの関連についても報告され ている。前述の通り MAM は生体膜の主成分であるリン脂質合成 の場であり、特にオートファゴソーム膜や隔離膜のソースとして の役割が注目されている。飢餓刺激時にオートファゴソームの形 成に重要な ATG14 や ATG5 が、SNARE タンパク質 STX17 によっ てEMCSsに集積されることや⁴⁴⁾、EMCSs形成因子である MFN2 や PACS2⁴⁵⁾ の発現抑制によって ATG14 の集積が阻害さ れることから、EMCSs がオートファゴソーム形成の場であるこ とが示唆された⁴⁴⁾。また、出芽酵母においても ERMES 複合体 の欠損によって隔離膜の形成が阻害され、マイトファジーが不全 になることが報告されている⁴⁶⁾。その一方で、Ca²⁺輸送に関与 する繋留因子である VAPB と PTPIP51 の結合形成は、恒常的オー トファジーに対して阻害的に作用することが報告されている ⁴⁷⁾。小胞体からミトコンドリアへの Ca²⁺の輸送を阻害すると、 この効果は見られないことから、ミトコンドリアの Ca²⁺レベル が関与していると思われるが、飢餓誘導によるオートファジーと の違いなど詳細は明らかにされていない。

5. 今後の課題

近年 EMCSs の形成や制御に関わる因子の同定や、様々な細胞 機能との関連についての研究が大きく進んだ。しかしその一方 で、解決すべき問題も多く残されている。EMCSs をはじめとし たオルガネラ接触領域の研究に言えるのは、オルガネラ全体に与 える影響を最小限にしつつ接触領域のみを操作することの困難さ である。現在同定されている形成因子の多くは小胞体やミトコン ドリアに広く分布しており、これらの遺伝子破壊や発現抑制に よって得られた表現型が、接触領域の形成不全だけに起因すると 断言することは難しい。この問題に対するアプローチとして、人 エリンカーの発現によるレスキュー実験が行われることもあるが その解釈には議論の余地がある。今後、EMCSs 形成能や EMCSs への局在能だけを欠損した変異体を探索する必要がある と思われる。

また、EMCSsの生成や崩壊の過程などのダイナミクスも興味深 い点であるが不明な点が多い。数十nmのEMCSsを正確に解析 するには超解像ライブイメージングが必要と考えられるが、現在 は一般的な研究室での実施は困難である。別のアプローチとして EMCSsのマーカー分子の探索や構築が考えられる。その一例とし て、最近報告されたBRETベースのEMCSsプローブはEMCSs の定量や動態解析に有効なツールとなる可能性がある⁴⁸⁾。このプ ローブは、単分子リンカーや二量体化リンカーのように2つの膜 を人工的に繋留し続けることはなく、より生理的なEMCSsの動 態を解析できると考えられる。ただ現時点では光学顕微鏡での観 察は感度的に難しく、同報告にある顕微鏡観察用のFRETベース のプローブは S/N 比が低く今後の改良が期待される⁴⁸⁾。

最後に、疾患など個体レベルの生命現象における EMCSs の意義の解明も今後の大きな課題である。これまでにアルツハイマー病の患者やモデルマウス由来の線維芽細胞においては野生型より

も2倍以上のEMCSsが形成されることが報告されているほか ^{49),50)}、Mfn2欠損によってEMCSsを減少させた細胞ではγ-セ クレターゼの活性が半分程度にまで減少する⁴⁹⁾。また、PTEN ⁵¹⁾やAKT⁵²⁾、p53⁵³⁾などのがん関連タンパク質がEMCSsに局 在していることも知られており、特に小胞体からミトコンドリア へのCa²⁺の輸送を制御することでがん細胞の増殖や細胞死を制 御していることも報告されている。このほかにもパーキンソン病 など様々な疾患との関連が報告されているが、その病態において 果たしてEMCSsがどこまで重要な役割を担っているかはよく分 かっていない。これまでに蓄積されたEMCSs形成の分子メカニ ズムや細胞機能の知見を、モデル生物や患者由来サンプルを用い た解析に取り入れることによりEMCSsの生理的な意義について 解き明かしていく必要があるだろう。

多様な分野の研究者同士の「接触」がオルガネラ接触領域の謎 を解き明かし、新たな発見がなされることを願いたい。

[参考文献]

- 1) L. Scorrano et al., Nat. Commun., 2019, 10, 1-11.
- 2) M. Wu et al., J. Cell Biol., 2006, 174, 803-813.
- 3) D. Peretti et al., Mol. Biol. Cell, 2008, 19, 3871-3884.
- 4) J. L. Costello et al., J. Cell Biol., 2017, 216, 331-342.
- 5) R. Hua et al., J. Cell Biol., 2017, 216, 367-377.
- 6) J. E. Lee et al., Science, 2020, 367, eaay7108.
- 7) J. E. Vance, J. Biol. Chem., 1990, 265, 7248-7256.
- 8) R. Rizzuto et al., Science, **1998**, 280, 1763-1766.
- 9) N. Shai et al., Nat. Commun., 2018, 9, 1761.
- 10) A. Das et al., J. Cell Biol., **2016**, 214, 831-845.
- 11) I. Y. Benador et al., Cell Metab., 2018, 27, 869-885.
- 12) W. Bernhard and C. J. Rouiller, Biophys. Biochem. Cytol., 1956, 2, 73-78.
- 13) S. J. Stone et al., J. Biol. Chem., 2000, 275, 34534-34540.
- 14) S. J. Stone et al., J. Biol. Chem., 2009, 284, 5352-5361.
- 15) A. E. Rusiñol et al., J. Biol. Chem., **1994**, 269, 27494-27502.
- 16) D. F. Montisano et al., Anat. Rec., **1982**, 203, 441-450.
- 17) P. T. Wang et al., J. Cell Sci., 2015, 128, 2759-2765.
- 18) Y. Hirabayashi et al., Science, 2017, 358, 623-630.
- 19) A. Sood et al., Proc. Natl. Acad. Sci. USA, 2014, 111, 16017-16022.
- 20) C. Cárdenas et al., Cell, 2010, 142, 270-283.
- 21) G. Szabadkai et al., J. Cell Biol., 2006, 175, 901-911.
- 22) T. Hayashi et al., Cell, 2007, 131, 596-610.
- 23) A. Santel and M. T. Fuller, J. Cell Sci., 2001, 114, 867-874.
- 24) O. M. de Brito and L. Scorrano, Nature, 2008, 456, 605-610.
- 25) D. Naon et al., Proc. Natl. Acad. Sci. USA, 2016, 113, 11249-11254.
- 26) R. Filadi et al., Proc. Natl. Acad. Sci. USA, 2015, 112, E2174-E2181.
- 27) P. Cosson et al., PLoS One, **2012**, 7, e46293.
- 28) R. Iwasawa et al., EMBO J., 2011, 30, 556-568.
- 29) K. J. De Vos et al., Hum. Mol. Genet., 2012, 21, 1299-1311.
- 30) M. Doghman-Bouguerra et al., EMBO Rep., 2016, 17, 1264-1280.
- 31) J. Zborowski et al., FEBS Lett., **1983**, *1*57, 179-182.
- 32) B. Kornmann et al., Science, 2009, 325, 477-481.
- 33) R. Kojima et al., Sci. Rep., **2016**, 6, 30777.
- 34) H. Jeong et al., EMBO Rep., 2016, 17, 1857-1871.
- 35) H. Jeong et al., Proc. Natl. Acad. Sci. USA, 2017, 114, E9502-E9511.
- 36) S. Lahiri et al., PLoS Biol., 2014, 12, e1001969.
- 37) A. Janer et al., EMBO Mol. Med., 2016, 8, 1019-1038.
- 38) J. R. Friedman et al., Science, **2011**, 334, 358-362.
- 39) F. Korobova et al., Science, **2013**, 339, 464-467.
- 40) U. Manor *et al.*, *elife*, **2015**, *4*, e08828.
- 41) J. E. Lee et al., Nature, **2016**, 540, 139-143.
- 42) D. M. Booth *et al.*, *Mol. Cell.*, **2016**, *63*, 240-248.
 43) V. Debattisti *et al.*, *Cell Rep.*, **2017**, *21*, 1667-1680.
- 44) M. Hamasaki *et al.*, *Nature*, **2013**, *495*, 389-393.

- 45) T. Simmen et al., EMBO J., 2005, 24, 717-729.
- 46) S. Böckler et al., Dev. Cell, **2014**, 28, 450-458.
- 47) P. Gomez-Suaga et al., Curr. Biol., 2017, 27, 371-385.
- 48) V. Hertlein et al., Life Sci. Alliance, 2019, 3, e201900600.
- 49) E. Area-Gomez et al., EMBO J., 2012, 31, 4106-4123.
- 50) E. Area-Gomez et al., Am. J. Pathol., 2009, 175, 1810-1816.
- 51) A. Bononi et al., Cell Death Differ, 2013, 20, 1631-1643.
- 52) T. Szado et al., Proc. Natl. Acad. Sci. USA, 2008, 105, 2427-2432.
- 53) C. Giorgi et al., Proc. Natl. Acad. Sci. USA, 2015, 112, 1779-1784.

[著者プロフィール]

 氏名:山本 真寿(YAMAMOTO Masatoshi)
 所属:熊本大学 大学院生命科学研究部 がん生物学講座 〒 860-8556 熊本市中央区本荘 1-1-1 Tel:096-373-5107
 Email:myamamoto@kumamoto-u.ac.jp
 出身学校:京都大学 大学院生命科学研究科
 学位:博士(生命科学) 氏名:山口 知也 (YAMAGUCHI Tomoya) 所属:熊本大学 大学院生命科学研究部 がん生物学講座 〒 860-8556 熊本市中央区本荘 1-1-1 Tel:096-373-5107 FAX:096-373-5108 Email:tyamaguchi@kumamoto-u.ac.jp 出身学校:名古屋大学 大学院医学系研究科 学位:博士 (医学) 専門分野:がん生物学、細胞生物学

ミトコンドリア染色試薬

専門分野:細胞生物学 オルガネラ間相互作用

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。



Topics on Chemistry

低分子蛍光色素によるミトコンドリア-リソソーム間相互作用の観察

真核生物の細胞内には、膜構造を持つ細胞小器官(オルガネラ) が存在している。オルガネラは、それぞれの役割を担いながら他 のオルガネラと相互作用して、細胞内の様々な機能を維持してい る。その一例として、正常な細胞では、マイトファジーと呼ばれ る不良ミトコンドリア消去機構が維持されており、不良ミトコン ドリアが二重の膜構造により周囲から隔離され、リソソームとの 融合により消化される。ミトコンドリアの恒常性の破綻は、アル ツハイマー病やパーキンソン病などの神経変性疾患、がん、老化 の原因となる。このように、ミトコンドリアの品質と恒常性の維 持が疾患や老化と密接に関わっていることから、ミトコンドリア の動的プロセス(分裂、融合、オルガネラ間相互作用)を研究す るためのプローブは生物学的研究において重要である。

本稿では、Sánchez らが開発した、ミトコンドリア-リソソー ム間相互作用を観察できる低分子蛍光色素 "MitoBlue" について 紹介したい¹⁾。

MitoBlue は、蛍光ビスアミジン骨格を有するカチオン性の色 素である(図1)。この色素は既存のミトコンドリア染色色素 Rhodamine123 とは異なり、膜電位に依存することなくミトコン ドリアを染色し、細胞固定後もミトコンドリアに保持されるとい う特徴がある。また、ミトコンドリア膜タンパク質マーカーの Tom20 抗体や、ミトコンドリアマトリックスマーカーの PDH 抗 体との共染色を行うと、MitoBlue のシグナルは PDH と共局在す ることから、MitoBlue はミトコンドリアマトリックスを染色す る色素となっている。興味深いことに、MitoBlue で細胞を長時 間染色し続けると、小胞状に染色される様子が観察され、 Rhodamine123とは異なる染色パターンとなる。これは、 MitoBlue とリソソーム染色試薬である LysoTracker を共染色さ せた結果、MitoBlueとLysoTrackerの局在が一致してくること から、MitoBlue がミトコンドリアからリソソームに移動するこ とが示唆された。そこで、MitoBlue を既存のミトコンドリア染 色色素 MitoTracker Deep Red (MTDR) とリソソームマーカー の Lamp1 抗体で共染色し、時間変化を確認したところ、染色直 後のMitoBlueはMitoTracker Deep Redと局在が一致していたが、 30 分後には Lamp1 のシグナルと重なり始めることから、 MitoBlue は時間依存的にミトコンドリアからリソソームに移動 することが観察された(図2)。

MitoBlue がミトコンドリアからリソソームへ移動するメカニ ズムを解明するため、mRFP-GFP-LC3 を発現 A541 細胞を用い て解析を行った。MitoBlue で染色した細胞をオートファジー誘 導剤であるラパマイシンで処理すると、LC3 のシグナルと MitoBlue のシグナルが一部共局在することが分かった。このこ とから、MitoBlue はオートファジーを介してミトコンドリアか らリソソームに移動することが示唆された。また、MitoBlue で 染色した細胞を、ミトコンドリア電子伝達系を阻害し、ミトコン ドリア小胞(MDVs)の形成を誘導する Antimycin A で処理すると、 MitoBlue のシグナルはリソソームの局在と一致した。このこと から、MitoBlue はミトコンドリア小胞によってもリソソームに 移動することが確認された。

上記のように、MitoBlue は時間依存的にミトコンドリア小胞 やオートファジー、マイトファジーを介してリソソームに移動す る(図3)。これにより、ミトコンドリア-リソソーム間相互作用 だけでなく、ミトコンドリアの動的プロセスを観察することに成 功している。 株式会社同仁化学研究所 田尻 智巳



図3 MitoBlueのミトコンドリアからリソソームへの移動メカニズム¹⁾

イメージングで使用されるプローブの多くは、目的のオルガネ ラのみをターゲットとして染色するため、動的プロセスを直接分 析することは難しい。しかし、MitoBlue はミトコンドリアの状 態に依存することなくミトコンドリアを染色し、その後のミトコ ンドリアの動的プロセスを追跡することが可能である。このよう なミトコンドリアとリソソーム間の相互作用をモニタリングでき る化合物により、これまで実態の明らかとなっていないオルガネ ラ間相互作用を今後解明することが期待される。

[参考文献]

1) M. I. Sánchez et al., Sci. Rep., 2020, 10, 3528.

用途で選ぶ細胞膜関連試薬

細胞膜は細胞の内側と外側を区分する境界としての役割だけでなく、物質の選択的な透過や伝達系などの重要な役割も担っていま す。また、細胞膜はエンドサイトーシスや神経細胞の軸索など動的な変化も大きいことから、細胞膜の動態を捉えることは様々な研 究分野で有用です。

<細胞膜関連試薬のラインアップ>

同仁化学研究所が販売している細胞膜関連試薬と、期待される用途を紹介致します。



細胞膜染色試薬

PlasMem Bright Green PlasMem Bright Red

<特長> ・生細胞で使用できる、染色後に固定できる ・低毒性で色素の滞留性が高い ・培地に色素を加えるだけ

PlasMem Bright シリーズは低毒性で色素の細胞膜滞留性が高く、生細胞や固定化細胞を用いた様々な実験でご利用いただくことが可能です。既存の細胞膜染色試薬に対して特に多かった不満を PlasMem Bright シリーズが解決します。

製品名	細胞毒性*	滞留性	染色後の洗浄	血清入り培地の使用	染色後の固定化
PlasMem Bright (Green, Red)	低	24 時間	不要	可	可(PFA)
S 社 製品 P	_	0.5 時間以下	必要	不可	可
T 社 製品 D	_	0.5 時間以下	必要	可	不可
T 社 製品 C	高	1.5 時間	必要	可	可(PFA)

※各染色試薬で染色後、神経細胞の形態変化(凝集)を比較した場合

<細胞膜の鮮明な可視化:神経細胞の形態と軸索内のミトコンドリア局在観察>





緑:ミトコンドリア (MitoBright LT Green)

<細胞膜への長期滞留性>

各細胞膜染色試薬で染色した細胞を 24 時間培養したのち、それぞれの蛍光画像を比較しました。結果、PlasMem Bright シリーズ は既存試薬と比較して長時間、膜に滞留することが確認されました。



品名	容量	希望納入価格(¥)	メーカーコード
PlasMem Bright Green	100 μ I $ imes$ 1	25,000	P504
PlasMem Bright Red	100 μ I $ imes$ 1	25,000	P505

※〈使用回数の目安〉100 μ I あたり、35 mm dish 10 枚、 μ -Slide 8 well 10 枚

製品の詳細や他社製品と比較した実験例などは、小社 HP をご覧ください。

膜染色同仁 検索

観察条件 PlasMem Bright Green: Ex 488 nm / Em 500-600 nm

観察条件 PlasMem Bright Red: Ex 561 nm / Em 560-700 nm

エンドサイトーシス検出試薬

ECGreen-Endocytosis Detection

<特長>

- ・より正確にエンドサイトーシスを可視化できる
- ・生細胞でエンドサイトーシスを追跡できる
- ・pH 変化に対する応答性が高い

ECGreen-Endocytosis Detection は小胞膜に局在し、pH に依存して蛍光を発する試薬です。そのため、蛍光アナログよりも直接的 にエンドサイトーシスを可視化することができます。また、本品は従来用いられてきた蛍光標識デキストランと比較して、pH 変化に 対する応答性に優れています。そのため、初期エンドソームも高感度に検出することが可能です。



染色原理の模式図

各試薬の pH に対する蛍光強度変化

<実験例:細胞内小胞輸送系を鮮明に可視化>

Wortmannin はエンドソームのリサイクリングやリソソームへの移行を阻害し、エンドソームの肥大化を惹起することが知られています。Wortmannin によるこれらの変化を ECGreen と初期エンドソームマーカータンパク質 Rab5-RFP(蛍光蛋白質)との共染色、および ECGreen とリソソーム染色試薬との共染色で確認しました。その結果、Wortmannin 添加時には ECGreen は肥大化したエンドソーム(Rab5-RFP)と共局在(左図:Merge)し、リソソームとは共局在しない(右図:Merge)ことが確認できました。上記の結果から、ECGreen は細胞内小胞輸送系の変化を可視化することができます。



※<使用回数の目安> 40 µI あたり、35 mm dish 20 枚、µ-Slide 8 well 20 枚

製品の詳細や他社製品と比較した実験例などは、小社 HP をご覧ください。

エンドソーム 同仁 検索

エクソソーム染色キット

エクソソーム膜 蛍光染色キット

ExoSparkler Exosome Membrane Labeling Kit-Green, Red, Deep Red エクソソームタンパク質 蛍光染色キット

ExoSparkler Exosome Protein Labeling Kit-Green, Red, Deep Red

- <特長>
 - ・細胞外で凝集しない
 - ・キットだけで蛍光標識から精製まで可能
 - ・エクソソームの性質にほとんど影響しない

製品の詳細や他社製品と比較 小社 HP をご覧ください。	交した実験例などは、
エクソソーム 同仁	検索

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選択的に染色し、細胞に取り込まれるエクソソームを イメージングすることができます。

(1) ExoSparkler Exosome Membrane Labeling Kit-Deep Red またはS 社製品 P(緑または赤)で染色したエクソソームを HeLa 細胞へ添加し、細胞内へ取り込まれるエクソソームを蛍光顕微鏡で確認した結果、製品 P(緑または赤)で染色したエクソソームにおいては、色素の凝集が疑われる細胞外の蛍光輝点が確認されました。



■観察条件 ExoSparkler Exosome Membrane Labeling Kit-Deep Red(紫): Ex 640 nm / Em 640-760 nm S 社 製品 P(緑): Ex 488 nm / Em 490-540 nm

S 社 製品 P(赤): Ex 561 nm / Em 570-640 nm

(2) ExoSparkler シリーズは、エクソソームの標識に最適化したプロトコルに加え、蛍光標識後の未反応色素を除去できるフィルト レーションチューブを同梱しているため、簡単な操作で蛍光標識エクソソームを調製できます。

ExoSparkler シリーズにおける染色工程



精製手法(未反応色素の除去)と回収率の比較

	回収率※
フィルトレーションチューブ(本キット)	50% 程度
ゲルろ過法	10% 程度
※ 小社での実施例:	

精製前後のエクソソーム粒子数を NTA(ナノ粒子トラッキング解析)で比較

(3) ExoSparkler Exosome Membrane Labeling Kit (Green, Red, Deep Red) で染色する前後のエクソソームについて、NTA (ナノ 粒子トラッキング解析) とゼータ電位を測定した結果、小社製品は、エクソソームの性質にほとんど影響を及ぼさないことが確認されました。比較データは製品 HP に掲載しています。

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
ExoSparkler Exosome Membrane Labeling Kit-Red	5 samples	25,000	EX02
ExoSparkler Exosome Membrane Labeling Kit-Deep Red	5 samples	25,000	EX03
ExoSparkler Exosome Protein Labeling Kit-Green	5 samples	20,000	EX04
ExoSparkler Exosome Protein Labeling Kit-Red	5 samples	20,000	EX05
ExoSparkler Exosome Protein Labeling Kit-Deep Red	5 samples	20,000	EX06

フェロトーシスの検出

ネクローシスやアポトーシス、オートファジーは細胞死関連プロセスとして知られています。2012年にフェロトーシスは新しい細胞死の1つとして提案されました。フェロトーシスは、鉄イオン依存性脂質過酸化物の蓄積によって引き起こされる非アポトーシス細胞死として研究されています¹⁾。フェロトーシスに関わる指標と関連製品を紹介します。

1) 豊國 伸哉, DOJIN News, 2017, 162, 1-3



フェロトーシス研究に関連するさまざまな指標と小社製品群

品名	Liperfluo	MitoPeDPP	Mito-FerroGreen	FerroOrange
指標	脂質過酸化	脂質過酸化	鉄イオン(Fe ²⁺)	鉄イオン(Fe ²⁺)
検出	蛍光	蛍光	蛍光	蛍光
(Ex/Em)	524 nm / 535 nm	452 nm / 470 nm	505 nm / 580 nm	543 nm / 580 nm
局在	細胞内	ミトコンドリア	ミトコンドリア	細胞内
装置	蛍光顕微鏡 フローサイトメーター	蛍光顕微鏡	蛍光顕微鏡 プレートリーダー	蛍光顕微鏡
サンプル	生細胞	生細胞	生細胞	生細胞
メーカーコード	L248	M466	M489	F374

	GSSG/GSH	Glutamine Assay Kit-	Glutamate Assay Kit-
	Quantification Kit	WST	WST
指標	GSSG/GSH	グルタミン	グルタミン酸
		比色	比色
19,11	412 nm	450 nm	450 nm
装置	装置 プレートリーダー プレートリーダー		プレートリーダー
	細胞		
++ヽ,プu.	組織	細胞	細胞
	血漿	細胞培養液	細胞培養液
	赤血球		
メーカーコード	G257	G268	G269

掲載した製品の詳細や文献などの情報は、 小社 HP をご覧ください。

メーカーコード 同仁 検索

細菌二重染色キット

-Bacstain- Bacterial Viability Detection Kit-DAPI/PI -Bacstain- Bacterial Viability Detection Kit-CTC/DAPI -Bacstain- Bacterial Viability Detection Kit-CFDA/PI

<特長>

- ・蛍光二重染色に最適化したプロトコルで試薬を添加するだけ
- ・複数の指標で薬剤効果や菌の状態を評価できる
- ・培地法と比較して評価にかかる時間を大幅に短縮できる

このような方におススメします!

☑ これから蛍光染色法による実験を始めたい!
 ☑ 培養法で判定している検査を迅速化したい!
 ☑ 損傷菌や VNC 菌を検出したい!
 ☑ 菌に対する薬剤効果を正確かつ簡便に評価したい!

試薬を添加するだけの簡単操作

評価にかかる時間を大幅に短縮



※微生物の増殖能力によって時間は異なります。

蛍光染色法による細菌の生死判定などでよく用いられる膜損傷の有無を指標とした染め分けに加え、-Bacstain-Bacterial Viability Detection Kit シリーズでは複数の指標を選んで染め分けすることができるため、細菌の状態を様々な視点から評価できます。

検出対象	全菌	膜損傷	呼吸活性	酵素活性
-Bacstain- Bacterial Viability Deteciton Kit-DAPI/PI	\checkmark	\checkmark		
-Bacstain- Bacterial Viability Deteciton Kit-CTC/DAPI	\checkmark		\checkmark	
-Bacstain- Bacterial Viability Deteciton Kit-CFDA/PI		\checkmark		\checkmark
T 社 製品 L	\checkmark	\checkmark		

現在、菌に対する抗菌剤等の効果を評価する場合、培養法によるコロニー形成単位(CFU)の確認や蛍光染色試薬を用いた多 重染色法による評価が広く用いられています。近年では特に感染症における薬剤耐性菌の出現や損傷菌が引き起こす食中毒等の 問題を受けて、抗菌性評価の信頼性向上を目的にこれまで一般的に行われてきた単一指標ではなく、複数の指標で薬剤等の抗菌 性を評価する必要性が高まっています。

また、日本国内では 2020 年の 6 月から食品を扱う全事業者に対して HACCP(危害分析重要管理点)に沿った衛生管理が制度 化されることから、特に飲食料品の製造分野では菌検査の簡素化ならびに迅速化の需要が高まっています。

呼吸活性 CTC formazan CFDA	orescein	全菌 DAPI PI 膜損傷	
品名	容量	希望納入価格(¥)	メーカーコード
-Bacstain- Bacterial Viability Deteciton Kit-DAPI/PI	1 set	20,000	BS08
-Bacstain- Bacterial Viability Deteciton Kit-CTC/DAPI	1 set	36,000	BS09
-Bacstain- Bacterial Viability Deteciton Kit-CFDA/PI	1 set	36,000	BS10

※<使用回数の目安> 1 set あたり、約 100 検体分

バイオフィルム形成阻害測定キット・バイオフィルム薬剤効果測定キット

Biofilm Formation Assay Kit Biofilm Viability Assay Kit <特長>

- 測定の手間を大幅に低減
- ・バラツキを抑えることが可能



B601

既存法はマイクロプレートの底にバイオフィルムを形成するため、菌の培養に伴う培地交換や、染色工程前後の洗浄作業に多くの 手間を要し、バラツキの問題がありました。本キットは蓋に固定されたピン上にバイオフィルムを形成させるため、培地交換や染色 工程が蓋を移すだけで完了し、操作が非常に簡便です。



微生物増殖アッセイキット

Microbial Viability Assay Kit-WST

<特長>

- ・評価にかかる時間を大幅に短縮
- ・生菌の活性評価を濁度から色の変化へ

本キットに含まれる小社の開発品である還元発色剤 WST-8 は、 電子メディエーターを介して微生物内で生成する NAD(P)H に よって還元され発色します。還元された WST-8 ホルマザン量は 微生物の代謝活性に比例しており、吸光度を測定することで、微 生物の生存率や活性度合いを簡便に測定することができます。

<操作>

96 well プレートに菌懸濁液を準備し、試薬を添加する。



<従来法との比較>*菌種によって最適条件は異なります。

M439

同仁

	従来法		本方法
	コロニー形成法	濁度法	WST 法
所要 時間	24 時間以上	18-24 時間	約8時間* (培養6時間) (発色2時間)
特徴	・標準法 ・操作が煩雑	 ・マイクロプレート アッセイが可能 ・代謝活性は見れない 	 ・短時間でのマイク ロブレートアッセ イが可能 ・代謝活性を指標と する

品名	容量	希望納入価格(¥)	メーカーコード
Microbial Viability Assay Kit-WST	100 tests	6,000	M420
	500 tests	21,300	101439

※本ページの上記3製品は、福岡県工業技術センター生物食品研究所との共同開発品です。

検索



膜がなが~く染まる

細胞膜染色試薬

PlasMem Bright Green PlasMem Bright Red

SH-SY5Y より分化誘導した神経細胞 (核とミトコンドリアを多重染色)

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





ドージンニュース No.174 令和2年9月16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 岩永竜弥 年4回発行 許可なくコピーを禁ず