

Dharmacon™ Trans-Lentiviral packaging kitの使用法

製品説明

Dharmacon[™] Trans-Lentiviral[™] packaging systemは、複製能の ないレンチウイルス粒子を効率的に作製するための試薬です。作製 したレンチウイルス粒子は、分裂中または非分裂中の哺乳動物細胞 に、shRNAまたはORF(オープンリーディングフレーム)を発現す るコンストラクトを導入するために使用されます(文献1)。 Trans-Lentiviral Packaging Systemは、パッケージングコンポー ネントが5つのプラスミドに分離されているため、市販のレンチウ イルスベクターシステムの中でも、優れた安全性プロファイルを提 供します。Trans-Lentiviral Packaging Systemの詳細な説明は文 献2にあります。10回のパッケージング反応用のキット (Dharmacon HEK293Tウイルス産生用細胞付き、または無し)、 および50反応用と100反応用のより大きなバルクサイズが利用可能 です。

キットの詳細な説明、コンポーネントの保管および出荷条件につい ては、Appendix A (7ページ)を参照してください。

Dharmacon Trans-Lentiviral packaging systemのコンポーネント

1. Trans-Lentiviral packaging mix

パッケージングミックスには、HEK293Tウイルス産生用細胞 へのコトランスフェクション後のトランスファーベクターの ウイルスパッケージングを容易にするために、5つのパッケー ジングプラスミドの最適化された混合物が含まれています (文献3、4)。これらのプラスミドは、レンチウイルス粒子 を作製するために必要なトランスのヘルパー機能と構造タン パク質および酵素タンパク質を提供します。パッケージング プラスミドのコンポーネントの詳細については、Trans-Lentiviral packaging systemのバイオセーフティ機能という タイトルのセクションを参照してください。 2. Calcium Phosphate Transfection Reagent

塩化カルシウムは、DNAと複合体を形成し、リン酸緩衝液を 添加して共沈すると、形質転換された接着細胞へのDNAの取 り込みを促進します。このキットには、塩化カルシウムと 2xHEPES緩衝生理食塩水(2x HBSS)が含まれており、最適 なpHと安定性について厳密にテストされています。トランス ファーおよびパッケージングベクターの高効率リン酸カルシ ウムトランスフェクションは、HEK293T細胞で特異的に得ら れます。

3. Dharmacon HEK293T Packaging Cell Line (細胞付き10反 応キットとして提供、または個別購入も可能)

HEK293T細胞は、レンチウイルス粒子のパッケージングに理 想的であり、Trans-Lentiviral packaging mixとDharmacon™ のshRNA/ORFトランスファープラスミドをコトランスフェク ションすると、高力価のレンチウイルス粒子を作製すること ができます。これらの細胞はSV40ラージT抗原(Simian Vacuolating Virus 40 TAg)を安定して発現し、SV40複製起点 を含むベクターからタンパク質を高レベルで発現させること ができます。

4. Control Vector(10反応用キットにのみ含まれる)

Dharmacon[™] GIPZ[™] Non-silencing Control は、shRNA packaging kitで提供されます。このNon-silencing Controlは、 TurboGFP[™] (Evrogen、モスクワ、ロシア)レポーター遺伝 子と、既知の哺乳類遺伝子と相同性のないshRNA配列 (Cold Spring Harbor miR-30 expression scaffold内にある)を発現 します。ORF packaging kitでは、Precision LentiORF RFP Controlが提供されます。Precision LentiORF RFP Controlが 提供されます。Precision LentiORF RFP Controlk、 ヒトORF配列の代わりにTurboRFP (Evrogen、モスクワ、ロ シア)レポーター遺伝子を発現します。

Cat #	10 Reaction Kit	10 Reaction Kit (with Dharmcon HEK293T cells)	50 Reaction Kit	100 Reaction Kit
Dharmacon Trans-Lentiviral shRNA Packaging Kit with Calcium Phosphate Transfection Reagent	TLP5912	TLP5917	TLP5913	TLP5914
Dharmacon Trans-Lentiviral ORF Packaging Kit with Calcium Phosphate Transfection Reagent	TLP5916	TLP5918	TLP5919	TLP5920

Dharmacon HEK293T 細胞は単体で購入可能です(Cat#: HCL4517)。

horizondiscovery.com

Dharmacon Trans-Lentiviral packaging systemのバイオセーフティ機能

Dharmacon Trans-Lentiviral packaging systemは、他の市販のレ ンチウイルスベクターシステムと比較して優れた安全性プロファイ ルを提供します。パッケージングコンポーネントは5つのプラスミ ドに分離されています(Figure1)。さらに、gag-proおよびtatrevの発現は、条件付きテトラサイクリン応答性プロモーターエレ メント(TRE)の制御下にあり、これらのウイルス成分の発現は、 テトラサイクリントランス活性化因子を発現する細胞に厳密に制限 されます。トランスレンチウイルスパッケージングシステムの詳細 な説明は文献2にあります。



Figure1 Dharmacon Trans-Lentiviral packaging systemに含まれるプ ラスミドベクター

Dharmacon Trans-Lentiviral packaging systemに含まれる5つのプラスミドベクターは以下の通り:

1. pTLA1-PAK

gag-proパッケージングコンストラクトには、HIV-1分子クローン SG3 (座標258~8384)のDNAフラグメントが含まれています。こ れは、パッケージングシグナル(Ψ)の39塩基の欠失と、env遺伝 子の1,357塩基の欠失(座標5,827から7,184)により改変されてい ます。RTおよびIN遺伝子のpolオープンリーディングフレーム (ORF)は、それぞれの最初のコドンを終止コドン(TAA)に置き 換えることによって削除されました。Sall制限酵素切断部位での平 滑末端ライゲーションによってフレームシフト変異がVprをコード する配列に導入されました。これらの変更により、RT、IN、およ びアクセサリータンパク質であるVpr、Vif、Vpu、およびNefの発 現が削除されます。Gag-Proの発現はTREプロモーターによって制 御されています。

2. pTLA1-ENZ

RTおよびINタンパク質は、アクセサリータンパク質Vprとのインフレーム融合によってウイルス粒子にパッケージ化されます。 Vpr-RT-IN融合ポリタンパク質の発現は、Tatタンパク質によってトランス活性化されるHIV-2 LTRの制御下にあります。

3. pTLA1-ENV

VSV-G発現ベクターには、ヒトサイトメガロウイルス(CMV)の前 初期プロモーターの制御化でVSV-gを発現するコンストラクトが含 まれています。さらに、SV40ポリアデニル化シグナルがVSV-gの下 流にクローン化されています。

4. pTLA1-TOFF

Clontech[™] Tet-Off[™]発現ベクターには、hCMV前初期プロモー ターの制御下にあるテトラサイクリン制御トランス活性化因子 (tTA) が含まれています。 Tet-Offは、Gag-Pro発現(pTLA1-PAK)およびTAT / REV発現(pTLA1-TAT / REV)の転写活性化に 必要です。

5. pTLA1-TAT / REV

HIV-1アクセサリータンパク質TatおよびRevは、内部リボソーム進入部位(IRES)を含むバイシストロン性ベクターから発現します。 その発現はTREプロモーターの制御下にあります。Tatの発現は、 GIPZ shRNAやPrecision LentiORF、pTLA1-ENZウイルスベクター 等に見られる5'LTRのトランス活性化に必要です。Revは、gag-pro およびvpr-RT-IN(pol)遺伝子、および完全長のウイルスゲノムを コードするmRNA転写物の核外輸送および翻訳に必要です。

一般的な封じ込めに関する考慮事項

Dharmacon Trans-Lentiviral packaging systemを使用してレンチ ウイルス粒子を作製する場合、研究者は、適切な実験室の安全ガイ ドラインの下で安全に利用して、さまざまな哺乳類細胞で遺伝子発 現を抑制または促進できるウイルス粒子を作製していることになり ます。レンチウイルスベクターの使用に関連する主なリスクは、 (1) 複製能力のあるレンチウイルス粒子を産生する可能性、およ び(2) 主要な調節遺伝子の挿入活性化/不活性化による発癌の可能 性です。例えば、癌遺伝子の活性化または腫瘍抑制因子の不活性化 です。

これらのリスクを軽減するには、レンチウイルスベクター粒子を取り扱う際にBSL-2または強化されたBSL-2封じ込めのいずれかが必要です。レンチウイルスベクターの封じ込めに関するガイダンスについては、組換えDNA諮問委員会(RAC)のガイドラインを参照してください。

米国では、米国保健福祉省、米国疾病予防管理センター、および米 国国立衛生研究所のBiosafety in Microbiological and Biomedical Laboratoriesをダウンロードしてください。 https://www.cdc.gov/biosafety/publications/bmbl5/

こちらも参照:組換えDNA分子を含む研究のためのNIHガイドライン (NIHガイドライン) もご参考ください。 https://osp.od.nih.gov/biotechnology/nih-guidelines/

References

- 1. Naldini, L. et al., In Vivo Gene Delivery and Stable Transduction of Nondividing Cells by a Lentiviral Vector. Science 272, 263–267 (1996).
- 2. Wu, X. et al., Development of a Novel Trans-Lentiviral Vector That Affords Predictable Safety. Molecular Therapy 2, 47–55 (2000).
- 3. Kappes J.C., Wu X, Safety considerations in vector development. Somat Cell Mol Genet 26, 147-58 (2001).
- 4. Kappes J.C. et al, Production of trans-lentiviral vector with predictable safety. Methods Mol Med 76, 449-65 (2003).

horizondiscovery.com

プロトコール

HEK293T細胞でのレンチウイルス粒子の作製

以下は、Trans-Lentiviral Packaging Kitを使用してレンチウイルス 粒子を作製するためのプロトコールです。レンチウイルス粒子の作 製は、リン酸カルシウム試薬を使用して、Dharmacon Trans-Lentiviral packaging mixを、shRNAまたはORFトランスファーベク ターとともにHEK293Tパッケージング細胞にコトランスファク ションすることから始まります。コトランスフェクションすると、 複製能力のないビリオンが培地に放出されるため、それらを収集し 使用します。以下に示す例は、6ウェルまたは100 mmの細胞培養 プレートにおいて、平均力価(約10⁶ TU / mL)のレンチウイルス 粒子の作製に最適化されており、超遠心分離による濃縮により高力 価(約10⁸ TU / mL)レンチウイルス粒子となります。

トランスファーベクターとパッケージングプラスミドDNAのコト ランスフェクションを成功させるためのいくつかの一般的なガイド ラインは次のとおりです。

- 定期的に継代され、活発に分裂している指数関数的に増殖する 細胞を使用することが重要です。トランスフェクションの前日 に細胞をプレーティングすると、最高のトランスフェクション 効率が得られます。細胞が継代から回復するのに十分な時間を 与える必要があります(通常 12時間以上)。トランスフェク ションの前に、細胞培養ディッシュで細胞を約70%のコンフル エンシーまで増殖させ、1:2の比率で少なくとも2日間連続継 代することにより、Dharmacon HEK293T細胞が迅速な複製状 態にあることを確認します。この方法で細胞を準備することは、 再現性良く最大のレンチウイルス粒子収量を確保するのに役立 ちます。
- 低純度のプラスミドDNAは、夾雑物の抑制効果によりトランス フェクションが不十分になる場合があります。プラスミドDNA を注意深く調製および精製することにより、細胞毒性の影響を 回避します。260 nm / 280nmの吸光度比を元にDNAの品質を評 価します。一般に、1.8~2.0の比率が許容できると見なされま す。

細胞のプレーティング

レンチウイルス粒子にパッケージする各トランスファーベクターに ついて、6ウェルプレートの少なくとも1つのウェルまたは100 mm 細胞培養プレートに細胞を準備します。

- HEK293T細胞をトリプシン処理後、細胞を数えます。
- 通常の増殖培地(DMEM高グルコース、ピルビン酸ナトリウム、 10%FBS、1%Pen / Strep)で細胞を希釈し、選択した細胞培養 プレートに最適な細胞密度にします(トランスフェクションの 時点での最適な細胞コンフルエンシーは85-95%でなければな りません)。
 - a. 6ウェルプレートの各ウェルについて、2 mL中で1.2×10⁶ 個の細胞を準備します。
 - b. 100 mmの細胞培養プレートごとに、14 mL中で5.5×10⁶ 個の細胞を準備します。
- ──37℃、二酸化炭素5%にて細胞をオーバーナイト培養します。

トランスフェクション

CaCl₂と2x HBSSを37℃のウオーターバス中で短時間にて解凍しま す。解凍後、両方の試薬を4℃で数週間保存しても、機能が失われ ることはありません。トランスフェクションを進める前に、試薬を 常温に戻す必要があります。

 6ウェルプレートの各ウェルまたは100 mm細胞培養プレート について、5 mL (Fisher Scientific Cat#14-959-1A) または50 mL (Fisher Scientific Cat#14-432-22) のポリスチレンチュー ブ中で、表に示した量のトランスファーベクターDNAおよび Trans-Lentiviral Packaging Mixを調製します。滅菌水を使用し て、DNA混合物を指定の総液量にします。

horizondiscovery.com

	Lentiviral transfer vector DNA (shRNA or ORF)	Trans-Lentiviral packaging mix	Total volume (with sterile water)
One well of a 6-well plate	6 µg	4.3 µL	135 µL
100 mm plate	42 µg	30 µL	945 μL
	RしたDNA溶液に、	表に示した量のCa	aCl₂を加えます。
	Ca	CI2	

One well of a 6-well plate	15 μL
100 mm plate	105 µL

 こぼすことなく試薬を完全に混合するのに十分な速度で チューブをボルテックスします。ボルテックスしながら、表 に示した量の2x HBSSを滴下(drop-wise)します。

	2x HBSS
One well of a 6-well plate	150 µL
100 mm plate	1050 µL

- 室温で3分間インキュベートします。このインキュベーション 中に、軽いチョークのような白色沈殿物が現れるはずです (沈殿物は必ずしも明白ではない場合があります)。
- トランスフェクションミックスの総量 (300 μLまたは2.1 mL)を細胞に滴下 (drop-wise) します。注:ピペッティング ロスのために実際の量がわずかに少なくなる場合があります が、これはトランスフェクション効率に悪影響を与えること はありません。
- 6. 37℃、二酸化炭素5%にて細胞を10~16時間培養します(この時間を延長しないでください)。
- 7. 16時間以上の培養後、レンチウイルストランスファーベク ターがEvrogen TurboGFPなどの蛍光レポーターを有する場合 は、それらの蛍光レポータータンパク質を細胞が発現してい るか顕微鏡を用いて調べ、トランスフェクション効率の指標 とします。注:培地の色はオレンジまたはオレンジ/黄色の場 合があります。これはウイルスの産生には影響しません。
- 8. Reduced Serum Mediumを調製します。
 - a. High Glucose DMEM (Fisher Scientific Cat #SH30243.LS)b. 5% Fetal Bovine Serum (Fisher Scientific Cat
 - #SH30070.03)
 - c. 2 mM L-glutamine (Fisher Scientific Cat #SH30034.01)
 - d. 1x Penicillin/Streptomycin (Fisher Scientific Cat #SV30010)
- 9. リン酸カルシウム含有培地を細胞から除去し、表に示した量のReduced Serum Mediumと交換します。

	Reduced Serum Medium
One well of a 6-well plate	2 µL
100 mm plate	14 μL

10. 37℃、二酸化炭素5%にて細胞をさらに48時間培養します。

ウイルス粒子の回収と濃縮

- 培地交換の48時間後に、培地を15 mLの滅菌済みのキャップ 付きコニカルチューブに移して、ウイルス粒子を含む上清を 回収します。
 注意:この段階では、感染性のウイルス粒子(複製能を持た ない)を扱っていることを忘れないでください。BSL-2生物を 扱うための推奨ガイドラインに従ってください。
- 2. 1,600×g、4℃で10分間遠心分離して、付着していない細胞や 細胞の破片をペレット化します。

- 6ウェルまたは100 mmプレートのいずれかで作製された平均 的な力価の未濃縮ウイルス粒子の場合、上清をゆっくりと除 去し、新しいチューブに移します。ピペッティングにより混 合し、必要に応じてレンチウイルス粒子を分注します。力価 測定のために1つの小さなアリコート(~20µL)を準備します。 レンチウイルス粒子は常に-80℃で保管してください。
- 4. 濃縮された高力価ウイルス粒子を作製するには、低速での遠 心分離により残りの細胞破片を除去した後、上清を滅菌済み 0.22~0.45 µmタンパク質低結合フィルター(Millipore Millex-HV 0.45 µm PVDFフィルター等)を用いてろ過するこ とをお勧めします。スイングバケット超遠心ローターでの超 遠心分離により濃縮します。ろ過した上清を滅菌超遠心 チューブに移します。血清を含まないDMEMを使用して、 ローターバケットの必要な最小容量まで容量を調整します。 SW28ローターの場合、23,000 rpmで1.5~2時間、4°Cで遠心 分離します。
- 5. 遠心分離後、上清を注意深く吸引し捨てます。
- チューブの底にあるペレットに、必要量のDMEM(血清なし)をピペットで加えます。
- 目に見えるペレットは、トランスフェクトされた細胞の培地 からの血清タンパク質で主に構成されています。ウイルス粒 子は、このタンパク質ペレットから取り外す必要があります。 DMEMをペレットに加えた後、4℃で5~10分間インキュベー トします。次に、泡の形成を避けつつ、約30回ゆっくりと上 下にピペッティングします。
- 再懸濁したペレットを滅菌マイクロフュージチューブに移し、 フルスピードで3~4分間遠心分離します。この遠心分離によ り、チューブの底に付着した血清タンパク質がペレット化さ れます。遠心分離後、上清を新しいマイクロフュージチュー ブに移し、必要に応じて分注します。力価測定のために1つの 小さなアリコート(~20 μL)を準備します。レンチウイルス 粒子は常に-80℃で保管してください。

レンチウイス粒子ストックの力価測定

ウイルス産生用細胞株(HEK293Tなど)だけでなく、実際の実験 で使用する細胞株でもレンチウイルス粒子ストックの力価を決定す ることが重要です。このプロトコールは、EvrogenTurboGFPや TurboRFPなどのレポーター遺伝子を発現するレンチウイルスベク ターの力価を決定する際に使用できます。Dharmacon[™] TRIPZ[™] shRNA製品のように、レポーターがTet誘導性プロモーターシステ ムの制御下にある場合は、細胞培養液にドキシサイクリンを必ず含 めてください。

- 形質導入の前日、24ウェル組織培養プレートにHEK293T細胞 をウェルあたり5×10⁴の密度でDMEM(10%FBS、1%Pen / Strep)中にて播種します。翌日、ウェルは40~50%のコンフ ルエントになります。細胞プレーティング密度は、他の細胞 株については経験的に決定する必要があります。
- 無血清培地を使用して、丸底96ウェルプレートでウイルスス トックを希釈します。テストするウイルス粒子ストックごと に1行を使用して、Figure 2に示すようにプレートを利用しま す。ウイルスストックの希釈は、以下の手順(ステップ4から 開始)で行います。目標は、一連の5倍希釈液を作製して、 390,625倍の最終希釈に到達することです。
- 3. 各ウェルに80 µLの無血清培地を加えます。
- 20 μLの解凍したウイルス粒子ストックを1列目の対応する各 ウェルに加えます(5倍希釈)。ウェルの内容物を10~15回ピ ペットで上下させます。ピペットチップを廃棄します。
- 新しいピペットを使用して、1列目の各ウェルから2列目の対応するウェルに20 µLを移します。ウェルの内容物を上下に10~15回ピペッティングにすることで混合後、ピペットチップを廃棄します。
- 新しいピペットチップを使用して、2列目の各ウェルから3列 目の対応するウェルに20 μLを移します。ウェルの内容物を10 ~15回ピペッティングすることで混合後、ピペットチップを 廃棄します。

horizondiscovery.com

- 3列目から8列目まで20 μLの溶液の移動を繰り返し、10~15回 上下にピペッティングし、各希釈間でピペットチップを交換 します。高品質のマルチチャネルピペッターを使用すること を強くお勧めします。ウイルス粒子ストックの希釈液を室温 で5分間プレインキュベートします。
- 8. Figure 3に示すように、テストするウイルス粒子ストックごと に1行を使用して、24ウェルプレートにラベルを付けます。
- 9. 24ウェルプレートの細胞から培地を取り除きます。
- 10. 225 µLの無血清培地を各ウェルに加えます。
- 細胞を含む24ウェルデスティネーションプレート(Figure 3) の各ウェルに、96ウェルプレート(Figure2)から25μLの希 釈ウイルス粒子を加えて細胞を形質導入します。たとえば、 96ウェルプレートのウェルA2から24ウェルプレートのウェル A1に25 μLを移します(Table 1)。
- 12. 形質導入した培養物を37℃で4時間培養します。
- 形質導入ミックスを培養液から取り除き、1 mLのDMEM (10%FBS、1%Pen / Strep)を静かに加えます。TRIPZなど のTet誘導性レンチウイルストランスファーベクターの場合、 培地に1 μg/ mLのドキシサイクリンを加えます。
- 14. 細胞を72時間培養します。
- 15. TurboGFPまたはTurboRFPを発現している細胞のコロニーを 数えるために、適切な数の蛍光細胞を含むデスティネーショ ンプレートのウェルを選択します。細胞は48時間の培養期間 にわたって分裂するため、各マルチセルコロニーを1つの形質 導入細胞として数えます。Figure 4はこの原理を示しています。
- 1 mLあたりのトランスダクションユニット(TU / mL)は、次の式を使用して決定できます。カウントされたEvrogen TurboGFPまたはTurboRFP陽性コロニーの数×希釈係数× 40 =#TU / mL。
- ・ 例: 55個のEvrogen TurboGFP陽性コロニーをウェルA3でカウント
- 55(Evrogen TurboGFP陽性コロニー)×625(希釈係数)×40 = 1.38×10⁶ TU / mL

力価の単位は1 mLあたりのトランスダクションユニット(TU)で 示されるため、40という数字は、力価測定に使用した25 μL(「使 用される希釈ウイルスの量」、Table 1)を1ミリリットルに変換す るために使用されます。



Figure 2. ウイルスストックの5倍段階希釈



Figure 3. ウイルスの力価を測定するために使用されるDharmacon HEK293T細胞を播種した24ウェル組織培養プレート

Table 1.ウイルス粒子の希釈系列の作成

Well				
Originating (96-well plate)	Destination (24-well plate)	Volume of diluted viral particles used	Dilution factor	
A1		25 µL	5*	
A2	A1	25 µL	25	
A3	A2	25 µL	125	
A4	A3	25 µL	625	
A5	A4	25 µL	3125	
A6	A5	25 µL	15625	
A7	A6	25 µL	78125	
A8		25 µL	390625*	

*非常に低い力価または非常に高い力価を期待する場合は、それぞれ ウェル1またはウェル8のいずれかを含めることをお勧めします。



Figure 4.コロニーカウントによるウイルス力価の決定 HEK293T細胞に、段階希釈したGIPZレンチウイルス粒子を形質導入 しました。形質導入の72時間後に細胞を蛍光顕微鏡で観察しました。 各円は個々のコロニーを表します。

よくある質問

リン酸カルシウム試薬を使用してトランスファーベクターと Dharmacon Trans-Lentiviral packaging mixをコトランスフェクト すると、トランスフェクション後にすべてのHEK293T細胞が死滅 しました。何が起こったのでしょうか?

細胞死は、トランスフェクションの36~48時間後に一般的に観察 されることがあります。ウイルスの回収時近くで細胞が死ぬのは、 実際にはレンチウイルス粒子が活発に生産されていることの肯定的 な兆候です。Trans-Lentiviral packaging mixコンポネントの高発現 が原因である可能性があります。逆に、トランスフェクション後 24時間以内に高度の細胞死が観察されるべきではありません。こ れは多くの場合、トランスフェクション中の取り扱いエラーが原因 です。トランスフェクション試薬の添加および培地の交換中に細胞 が除去される可能性があります。試薬は必ずゆっくりと穏やかに細 胞に加えてください。トランスフェクション前のコンフルエンシー が85~95%であることを確認し、細胞培養フードからインキュ ベーターに輸送する際はプレートを慎重に取り扱ってください。

トランスフェクション後にHEK293T細胞の形態が変化するのはな ぜですか?

VSV-g糖タンパク質の発現により、HEK293T細胞が融合し、多核融 合細胞が出現します。この形態学的変化は正常であり、レンチウイ ルス粒子の産生には影響しません。

リン酸カルシウム-DNA複合体を形成するために各試薬を追加する 順序に違いはありますか?

示された順序で試薬を追加することにより、トランスフェクショ ンプロトコールを最適化しました。試薬の添加順序を変更してもト ランスフェクションは成功する可能性がありますが、プロトコール に示されている順序でトランスフェクション反応を組み立てること をお勧めします。

HBSSを追加しながらサンプルをボルテックスすることは本当に必要ですか?

いいえ、DNA反応ミックスをボルテックスすることは必須ではあ りませんが、可能な限り最高のトランスフェクション効率を達成す るのに役立つはずです。

リン酸カルシウム-DNA複合体は室温でどのくらいの時間インキュ ベートできますか?

リン酸カルシウムによるDNAの沈殿は、時間と温度に依存します [M. Jordan, A. Schallhorn, Transfecting mammalian cells: optimization of critical parameters affecting calcium-phosphate precipitate formation. Nucleic Acids Res. 24, 596-601 (1996)]。私 たちの経験では、最適なトランスフェクション効率は、HBSSの添 加後約3分で得られます。

リン酸カルシウムを使用したHEK293T細胞へのDNAのトランス フェクションにはどの程度の効率が期待できますか?

通常、90~95%のトランスフェクション効率が得られます。

回収した上清において、どの程度のウイルス力価を期待することが できますか?

達成可能なウイルス力価は、トランスファーベクター間で異なりま す。GIPZ shRNAコンストラクトをパッケージ化する場合、通常、 106 TU / mL以上の力価が得られます。レンチウイルス粒子を濃縮 するためにここで説明したプロトコールに従うと、通常、108 TU / mLの範囲の力価を得ることができます。多くの場合、トランス フェクトされたプレートの数を増やしたり、再懸濁量を減らしたり することで、109 TU / mL以上の力価を達成できます。

ウイルス粒子の濃縮に備えてウイルス上清をろ過するために、どの ミリポアフィルターを使用しますか?

ろ過する容量に応じて、さまざまなフィルターを使用します。

- ≤ 50 mL: Millipore Cat #SCGP00525, SteriflipTM-GP Filter Unit, 0.22 µM PES membraneOr, Millipore Cat #SE1M003M00, Steriflip-HV Filter Unit, 0.45 µM PVDF membrane
- ≤ 150 mL: Millipore Cat #SCGPU01RE, Stericup-GP Filter Unit, 0.22 µM PES membrane
- ≤ 500 mL: Millipore Cat #SCGPU05RE, Stericup-GP Filter Unit, 0.22 µM PES membrane
- ≤ 1000 mL: Millipore Cat #SCGPU10RE, Stericup-GP Filter Unit, 0.22 µM PES membrane

ウイルスを濃縮するためのプロトコールに記載されているのと同じ 種類のローターがありません。私のローターが機能することをどの ように知ることができますか?

SW28ローターシステムの使用は必要ありません。他のローターサ イズも使用てきます。別のローターを使用する場合は、ローターの G力と実行時間を相互参照します。さらに、次のジャーナル記事は、 レンチウイルスを濃縮するためのSW32Tiローター(19,500 rpmで 2時間)の使用について言及しています:[F. Di Nunzio, B. Piovani, Transduction of human hematopoietic stem cells by lentiviral vectors pseudotyped with the RD114-TR chimeric envelope glycoprotein. Hum. Gene Ther., 18, 811-820 (2007)]。以下のURLの ウェブサイトでrpmからrcfへの変換ができます: http://insilico.ehu.es/mini_tools/rcf_rpm.php

horizondiscovery.com

Use of Product and Label Licenses

The Products, use and applications, are covered by pending and issued patents. The following Label Licenses govern all use of the Products. It is each Purchaser's responsibility to determine if patent or other intellectual property rights held by third parties may restrict the use of Products for a particular application.

The Broad Institute Inc. and Massachusetts Institute of Technology

Limited License to Use. The purchase of this product from Dharmacon or its Affiliates conveys to the final purchaser ("Limited Licensee") a non-transferable right to use the CRISPR Products solely for research conducted by Limited Licensee in accordance with all of the following requirements: (i) The Limited Licensee shall not sell or otherwise transfer products (including without limitation any material that contains a CRISPR Product in whole or part) to any other person or entity, or use CRISPR Products to perform services for the benefit of any other person or entity, provided however, that a Limited Licensee that is an academic, nonprofit, or government entity or person may perform core facility services for its academic and teaching purposes; (ii) the Limited Licensee shall use only the amount of the CRISPR Products purchased from Dharmacon, and components of such CRISPR Products, and shall use any materials made using the CRISPR Products by such Limited Licensee, only for its internal research as a research tool for research purposes, and not for any commercial purposes, (iii) the Limited Licensee shall use CRISPR Products in compliance with all applicable laws and regulations, including without limitation applicable human health and animal welfare laws and regulations, (iv) Broad, MIT and Harvard shall provide no warranties of any kind to the Limited Licensee (statutory or implied concerning certain patent rights owned by Broad, MIT and Harvard or the CRISPR Products), including without limitation, as to product quality, condition, description, merchantability, fitness for a particular purpose, noninfringement of intellectual property rights or the absence of latent or other defects, and all such warranties are hereby expressly disclaimed, (v) Broad, MIT and Harvard shall also expressly disclaim any warranty regarding results obtained through the use of the CRISPR Products, including without limitation any claim of inaccurate, invalid or incomplete results, (vi) Broad, MIT and Harvard, and their directors, trustees, officers, employees, agents, faculty, affiliated investigators, and students, shall have no liability to the Limited Licensee, including, without limitation, for any loss of use or profits, business interruption or any consequential, incidental, special or other indirect damages of any kind, regardless of how caused and regardless of whether an action in contract, tort, strict product liability or otherwise, (vii) the Limited Licensee shall indemnify, defend and hold harmless MIT, Harvard and Broad against any liability, damage, loss, or expense (including without limitation reasonable attorneys' fees and expenses) incurred by or imposed upon MIT, Harvard and Broad in connection with any claims, suits, investigations, actions, demands or judgments arising out of or related to the exercise of any rights granted to the Limited Licensee under the Limited License or any breach of the Limited License by such Limited Licensee; provided, however, that the foregoing shall not apply to a Limited Licensee that is an academic, non-profit, or government entity or person, and such Limited Licensee instead agrees that it, and not MIT, Harvard and Broad, shall be responsible for any liability, damage, loss or expense arising out of or related to the exercise of any rights granted to the Limited Licensee under the Limited License or any breach of the Limited License by the Limited Licensee, and (viii) the CRISPR Product and its use may be the subject of one or more issued patents and/or pending patent applications owned by Broad, MIT and Harvard and the purchase of the CRISPR Product does not convey a license under any claims in the foregoing patents or patent applications directed to the CRISPR Product or use, production or commercialization thereof.

Salk Institute Restricted Use Label License THE SALK INSTITUTE PURCHASER NOTIFICATION

1. The Nature of the LICENSEE License: Licensee has a license to sell the Product containing WPRE, under the terms described below. Any use of WPRE outside of Licensee's Product or the Product's intended use, requires a license as detailed below. Before using the Product containing WPRE, please read the following license agreement. If you do not agree to be bound by its terms, contact Licensee within 10 days for authorization to return the unused Product containing WPRE and to receive a full credit.

- Patents: The WPRE technology is covered by patents issued to The Salk Institute for Biological Studies.
- Individual License: The WPRE technology is covered by patents issued to The Salk Institute for Biological Studies.
- 4. Agreement: Licensee grants you a non-exclusive license to use the enclosed Product containing WPRE in its entirety for its intended use. The Product containing WPRE is being transferred to you in furtherance of, and reliance on, such license. Any use of WPRE outside of Licensee's Product or the Product's intended use, requires a license from the Salk Institute for Biological Studies.
- Termination of License: This license agreement is effective until terminated. You may terminate it at any time by destroying all Products containing WPRE in your control. It will also terminate automatically if you fail to

horizondiscovery.com

comply with the terms and conditions of the license agreement. You shall, upon termination of the license agreement, destroy all Products containing WPRE in your control, and so notify Licensee in writing. This License shall be governed in its interpretation and enforcement by the laws of the State of California.

6. Contact for WPRE Licensing: The Salk Institute for Biological Studies, 10010 North Torrey

Pines Road, La Jolla, CA 92037 Attn.: Office of Technology Management Phone: (858) 453-4100 extension 1703 Fax: (858) 546-8093

Sigma-Aldrich Restricted Use Label License

These Products are For Research Use Only - Not for any Clinical, Therapeutic or Diagnostic use in Humans. These Products may be subject to the following patents (issued or pending):

Oxford technology

PCT Application (filing date)	Priority Application(s) (filing date)	National Application	Issued Patents
	US 07/586,603 (21 Sept 1990)	US 08/361,839	5,817,491
PCT/US91/05699 (9 Aug 1991)	US 07/658,632 (19 Feb 1991)	EP 91915104.3 JP 3-514518 AU 84302/91 CA 2,104,396	0572401 3547129 663470 2,104,396
	US 07/1 70,515 (21 March 1988) US 07/1 70,515 (21 March 1988)	US Continuation 08/156,789 US Continuation 08/462,492 US Continuation 10/205,179	5,591,624 5,716,832
PCT/US91/06852 (20 Aug 1991) WO 92/05266	US 07/586,603 (21 Sept 1990)	AU 88424/91 AU Divisional 47984/96	665176 690427
PCT/GB97/02857 (17 Oct 1997) WO 98/1 7815	GB 9621680.9 (17 Oct 1996) GB 962445739 (25 Nov 1996)	EP 97909436.4 EP Divisional 00202432.1 US 09/224.014 US Divisional 09/91 5,169 US CIP 10/661,761 US Continuation 11/646,041 JP 10-519086 AU 47122/97 CN97198767.X NZ 334860	0904392 6,312,682 6,669,936 7,198,784 725143 2L97198767.X 334860
PCT/GB97/02858 (17 Oct 1997) WO 98/17816	GB 9621680.9 (17 Oct 1996)	GB 9903117.1 EP 97909437.2 US 09/284,011 JP 10-519087 AU 47123/97 CN 97198883.8 NZ 334522	2331522 6,235,522 737801 ZL971 98883.8 334522
PCT/GB97/02696 28 Oct 1997) WO 98/1 8934	GB 9622500.8 (29 Oct 1996)	US CIP 10/324,616 US Continuation 11/155,043 US Continuation 11/726,679 JP 10-520197	6,924,123 7,056,699

Benitec technology licensed patents

Name Inventor	Title	Patent No.	Application No.
Graham/Rice	Control of Gene Express	ZA 2000/4507* AU2001100608* SG75542* US 6,573,099* GB2353282* AU 743316* NZ 506648* US 10/346853* PCT/AU/99/00195* BR P19908967-0* CA 2323726* CB 99804255-2* CZ PE2000-3346* EP 99910039.9* US10/646,070 US 10/759,841* EP 04015041.9* AU2005211538	HK 01105904.3* HU P0101225* IN 2000/00 169/DEL* JP P2000-537990 KR 7010419/2000* MX008631* PL P.343064* SK PV 1372-2000* AU 35647/02* SG200205122-5* US 09/646807* PP2492/98* AUPP2499/98* US 10/821,710* US 10/821,726* AU 2005209648
Graham/Rice/M/R	Genetic Silencing	WO 01/70949 GB 237722 AU PQ6363	SG 91678 ZA2002/7428 AU2001240
Grahm, et al.	Double-Stranded Nucleic Acid	AU 2003906281 AU 2003906281 US 10/861,191	AU 2004902279 PCT/AU04/000759

*Indicates a "Co-Owned Patents" are jointly owned by Benitec and CSIRO.

Appendix A

Dharmacon Trans-Lentiviral ORF Packaging Kit with Calcium Phosphate Transfection Reagent

Kit contents	TLP5916 (10 rxn)	TLP5918 (10 rxn + cells)	TLP5919 (50 rxn)	TLP5920 (100 rxn)	Shipping condition	Storage
Dharmacon Trans- Lentiviral Packaging Mix	285 µg	285 µg	285 µg x 5	285 µg x 10	Wet ice	-20 °C
Dharmacon HEK293T Packaging Cell Line	not included	1.0 mL	not included	not included	Dry ice	Liquid N ₂
CaCl ₂ Reagent	1.2 mL	1.2 mL	6 mL	6 mL x 2	Wet ice	-20 °C
2x HBSS Reagent	12 mL	12 mL	60 mL	60 mL x 2	Wet ice	-20 °C
Dharmacon Precision LentiORF RFP Control DNA	45 μg (0.45 μg/μL)	45 μg (0.45 μg/μL)	not included	not included	Wet ice	-20 °C

Dharmacon Trans-Lentiviral shRNA Packaging Kit with Calcium Phosphate Transfection Reagent

Kit contents	TLP5912 (10 rxn)	TLP5917 (10 rxn + cells)	TLP5913 (50 rxn)	TLP5914 (100 rxn)	Shipping condition	Storage
Dharmacon Trans- Lentiviral Packaging Mix	285 µg	285 µg	285 µg x 5	285 µg x 10	Wet ice	-20 °C
Dharmacon HEK293T Packaging Cell Line	not included	1.0 mL	not included	not included	Dry ice	Liquid N ₂
CaCl ₂ Reagent	1.2 mL	1.2 mL	6 mL	6 mL x 2	Wet ice	-20 °C
2x HBSS Reagent	12 mL	12 mL	60 mL	60 mL x 2	Wet ice	-20 °C
Dharmacon GIPZ Non- Silencing Control DNA	45 μg (0.45 μg/μL)	45 μg (0.45 μg/μL)	not included	not included	Wet ice	-20 °C

For more information



©2023 Horizon Discovery Group Ltd and its affiliated companies. All rights reserved. Revvity is a trademark of Revvity, Inc., registered in the United States and other countries. Horizon Discovery is a trademark of Horizon Discovery Ltd. Dharmacon is a trademark of Dharmacon, Inc. All Revvity, Inc., Horizon Discovery Ltd and Dharmacon, Inc. trademarks are used with permission. Other product names and brand names may be trademarks or registered trademarks of their respective owners.



本文書は英語版プロトコルを元にした日本語版です。内容は英語版プロトコルの最新バージョンとは異なることがあります。

B048-2106-v3