



Instruction Manual

PureLink™ 384 Plasmid Purification System

Catalog no. 12263-026

Version B

20 April 2005

25-0451

Table of Contents

Introduction 1

Preparing 96- or 384-Well Cell Cultures 3

Protocol A, Direct Load Method 6

Protocol B, Harvested Cells Method 9

Troubleshooting Guide 12

Centrifuge Information 14

Related Products 15

Technical Service 16

Introduction

Introduction

PureLink™ 384 Plasmid Purification Plates are part of a unique system for high-throughput plasmid DNA isolation. This solid-phase lysis system replaces the numerous reagents and mixing steps associated with traditional alkaline lysis. Since the lysis occurs within a solid-phase matrix, there is no mixing or transfer of lysate to create cross-contamination. Harvesting and resuspending of cells are avoided by growing cultures in rich media.

System Overview

To use PureLink™ 384 Plasmid Purification Plates as part of a high-throughput plasmid purification system, first you prepare cell cultures from rich media in 96- or 384-well growth blocks (see page 3). Since proper growth of bacterial cultures is a key element in this system, we strongly recommend that you grow and analyze test cultures before attempting purification.

Then, depending on the desired yield of plasmid DNA, proceed to Protocol A, Direct Load Method (2–3 µg DNA), or Protocol B, Harvested Cells Method (3–5 µg DNA). Protocol A is the simpler, less time-consuming method and should provide ample yield for most users.

In Protocol A (page 6), you stack the 384-well filter plate on top of a receiver plate containing isopropanol and load pre-lysis buffer containing lysozyme onto the filters. Then you load the cells directly into the wells of the filter plate, followed by lysis buffer. Centrifuge the stacked plates to recover the plasmid DNA and precipitate it in one step. Genomic DNA and other cell components will remain within the filter matrix.

In Protocol B (page 9), you stack the 384-well filter plate on top of a receiver plate containing isopropanol and load pre-lysis buffer containing lysozyme onto the filters. You then resuspend harvested cells to the desired concentration and load them into the wells of the filter plate, followed by lysis buffer. Centrifuge the stacked plates to recover the plasmid DNA and precipitate it in one step. Genomic DNA and other cell components will remain within the filter matrix.

System Features

This system has the following features:

- Protocols are compatible with a variety of automated devices.
- Approximately 66 plates (>25,000 samples) can be processed in one shift using a robot with a 96-well head, a 96/384 well dispenser, and one centrifuge with a 6-place rotor.
- Yields of 2–3 µg/well for Protocol A (Direct Load Method) and 3–5 µg/well for Protocol B (Harvested Cells Method) are typical for high copy number plasmids.
- Filter plates can be processed one at a time or in stacks of up to three filter plates per centrifuge plate carrier.

Plasmid DNA isolated using this method is suitable for use in automated fluorescent DNA sequencing, PCR, or restriction digestion.

Continued on next page

Introduction, Continued

Contents

PureLink™ 384 Plasmid Purification Plates are sold in packages of 25 filter plates (Cat. no. 12263-026).

The plates are shipped at room temperature and should be stored at room temperature.

Reagents Needed

The following reagents are needed before starting the procedures.

Component	Composition	Amount	Cat. No.
Cell suspension buffer	50 mM Tris-HCl (pH 8.0), 10 mM EDTA	500 ml	12056-016
Pre-lysis buffer	Proprietary	125 ml	12263-034
Lysis buffer	Proprietary	1 liter	12057-022
RNase A solution	20 mg RNase A/ml in 50 mM Tris-HCl (pH 8.0), 10 mM EDTA	10 ml	12091-021
		25 ml	12091-039
TE buffer	10 mM Tris-HCl (pH 8.0), 0.1 mM EDTA	100 ml	12090-015
<u>Media</u>			
Terrific Broth	See page 3 for recipe	500 g	22711-022
MR2001	See page 3 for source		—
Antibiotic	See page 3 for concentrations		—
Lysozyme			—
Isopropanol			—
70% ethanol			—

Materials Needed

The following equipment and materials are needed before starting the procedures.

- PureLink™ Air-Porous Tape (Catalog no. 12262-010)
- PureLink™ Foil Tape (Catalog no. 12261-012)
- PureLink™ 96-well Growth Blocks (Catalog no. 12256-020) **or** 384-well growth blocks (required volume: 200 µl/well)
- 384-well polypropylene receiver plates (required volume: 200 µl/well)
Contact technical service (see page 16) for assistance with obtaining appropriate receiver plates
- Shaking incubator at 37°C
- Refrigerated centrifuge with swinging bucket rotor capable of reaching 3000 × g and with a plate height clearance of 3.6 cm (see page 14 for additional information)

Preparing 96- or 384-Well Cell Cultures

Procedural Overview

In this procedure, you prepare cell cultures from rich media in 96- or 384-well growth blocks.



Note

For direct loading of cultures, it may be easier to attain optimal concentrations of cells using 96-well growth blocks.

We strongly recommend that you grow and analyze some test cultures before attempting purification.

Recommended Cell Type

While this system is compatible with various strains of *E. coli*, DH10B cells are recommended for optimal performance and plasmid yield. This strain is compatible with direct cloning of genomic DNA and is available as both chemically competent and electrocompetent cells.

Media Recipes and Sources

The following table shows the recipe or source for recommended rich media:

Media	Recipe for 1 liter or Source
Terrific Broth	Available from Invitrogen (500 g, Cat. no. 22711-022) Recipe: 11.8 g peptone 23.6 g yeast extract 9.4 g dipotassium hydrogen phosphate (K_2HPO_4) 2.2 g potassium dihydrogen phosphate (KH_2PO_4) 4 ml glycerol Bring to 1 liter with deionized water and autoclave for 15 minutes at 121° C.
MR2001	MacConnell Research, San Diego, CA

Recommended Antibiotic Concentrations

The following table shows the recommended concentrations for different types of antibiotics:

Antibiotic	Concentration in culture medium
Ampicillin (Cat. no. 11593-019)	100 µg/ml
Chloramphenicol	30 µg/ml
Gentamicin (Cat. no. 15750-060)	10 µg/ml
Kanamycin (Cat. no. 15160-054)	50 µg/ml
Streptomycin (Cat. no. 11860-038)	100 µg/ml
Tetracycline	25 µg/ml

Continued on next page

Preparing 96- or 384-Well Cell Cultures, Continued

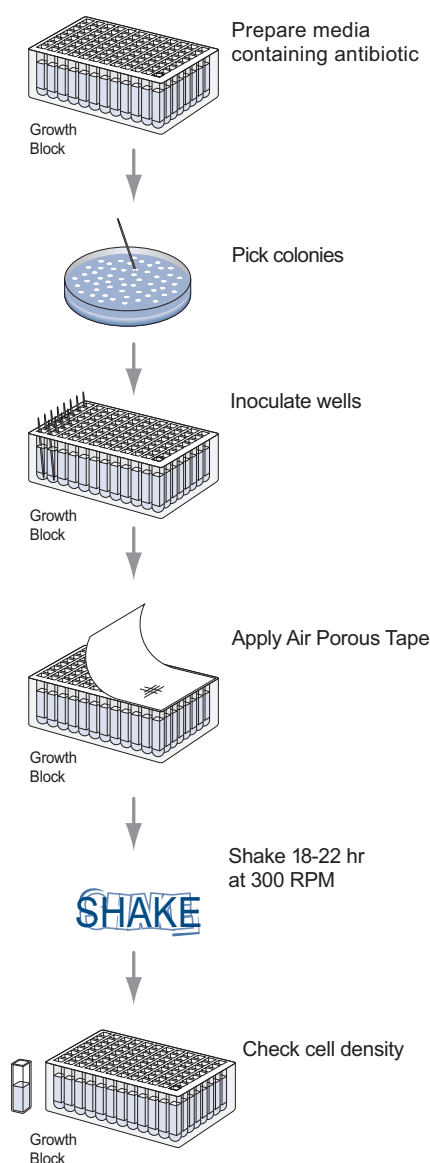
Critical Parameters

For the following procedure, the parameters listed below are critical:

- Antibiotic: Verify concentration and freshness (see previous page).
- Age of colonies: <1 week (the fresher the better).
- Recommended media: Terrific Broth or MR2001 (check for clarity).
- Incubation temperature: 37°C.
- Shaking speed: 300–320 RPM.
- Incubation time: 18–22 hours (20–24 hours for glycerol stock inoculation).

Preparing Cultures

Follow the steps below to prepare your 96- or 384-well cell cultures:



1. Add an appropriate concentration of antibiotic to rich media (Terrific Broth or MR2001 is recommended; see previous page for recipe/source).
2. Add 0.5 ml of antibiotic-containing media to each well of a 96-well growth block
OR
Add 0.15 ml of antibiotic-containing media to each well of a 384-well growth block (well capacity: 200 µl/well).
3. Inoculate all wells of the growth block with *fresh* colonies (<1 week old).
Note: Glycerol stocks should be used with great care. Old glycerol stocks can produce low plasmid copy numbers, even though the cells grow adequately. Cell growth will be delayed out of glycerol stocks (see increased culture times in step 5).
4. Cover the growth block with air porous tape. Seal cover securely.
5. Incubate in a 37°C shaker at 300–320 RPM for 18–22 hours (20–24 hours for cells from glycerol stocks).
6. Check growth of a few random wells in the block by measuring OD₆₀₀ of a 1:33 dilution in TE. (Note that different dilutions will result in different calculated undiluted OD₆₀₀.) Be certain that cells are well suspended before removing aliquot. Use 30 µl media plus 970 µl TE as a blank. See recommended cell densities for the different protocols on the next page.

Continued on next page

Preparing 96- or 384-Well Cell Cultures, Continued

Recommended Cell Densities

For optimal performance, the OD₆₀₀ of a 1:33 dilution should be **0.18–0.33** (undiluted OD₆₀₀ 6–10) for Protocol A, Direct Load, or **>0.12** (undiluted OD₆₀₀ >4) for Protocol B, Harvested Cells.

Storing Cultures

Cultures can be refrigerated or frozen, depending on how soon you want to use them:

- Cultures can be stored up to 3 days at 4°C before use. Resuspend cultures thoroughly just before using.
 - Cultures can be stored up to one month frozen at -20°C. Air porous tape should be removed and replaced with foil tape before storing.
-

Thawing Cultures

To thaw frozen cultures:

1. Store block at 4°C for at least 18 hours.
 2. Resuspend cells thoroughly before use.
-

Protocol A, Direct Load Method

Protocol Overview

In this protocol, you stack the 384-well filter plate on top of a receiver plate containing isopropanol and load pre-lysis buffer containing lysozyme onto the filters. Then you load the cells directly into the wells of the filter plate, followed by lysis buffer. Centrifuge the stacked plates to recover the plasmid DNA and precipitate it in one step. Genomic DNA and other cell components will remain within the filter matrix.

Required Cell Density

For optimal performance, this protocol requires an OD₆₀₀ of **0.18–0.33** for a 1:33 dilution (undiluted OD₆₀₀ 6–10). If the density of the 1:33 dilution is lower, optimize growth conditions as described in the previous section or use Protocol B, Harvested Cells (see page 9).

Preparing the Pre-lysis Buffer with Lysozyme

Immediately before using this protocol, follow the steps below to prepare the pre-lysis buffer with lysozyme:

1. Calculate the necessary volume to prepare using the following formula:
$$(\text{total vol pre-lysis buffer}) = [(\# \text{ plates}) \times (2 \text{ ml/plate})] + (\text{waste vol})$$

The waste volume is the volume required to fill the lines or troughs of the robot or pipetting device to be used. The waste volume should be a minimum of 2 ml.
 2. Calculate the necessary amount of lysozyme to weigh using the following formula:
$$(\text{total mg lysozyme}) = (\text{total vol pre-lysis buffer}) \times (5 \text{ mg/ml})$$
 3. Add the lysozyme to the pre-lysis buffer and invert 30 or more times to thoroughly dissolve.
Note: This solution should be prepared fresh daily. Alternatively, a large batch can be prepared and stored aliquoted for single use at –20°C. Aliquots should be frozen and thawed only once.
-

Preparing the Lysis Buffer with RNase A

Immediately before using this protocol, prepare the lysis buffer/RNase A mixture according to the directions below:

1. Calculate the necessary volume of lysis buffer and RNase A using the following formulas:
$$[(\# \text{ plates}) \times (12 \text{ ml/plate})] + (\text{waste volume}) = (\text{total vol lysis buffer})$$
$$(\text{total vol lysis buffer}) \times (0.05) = (\text{total vol RNase A})$$
 2. Mix the lysis buffer and RNase A by inverting 20 times to mix thoroughly. Use this solution within 24 hours.
-

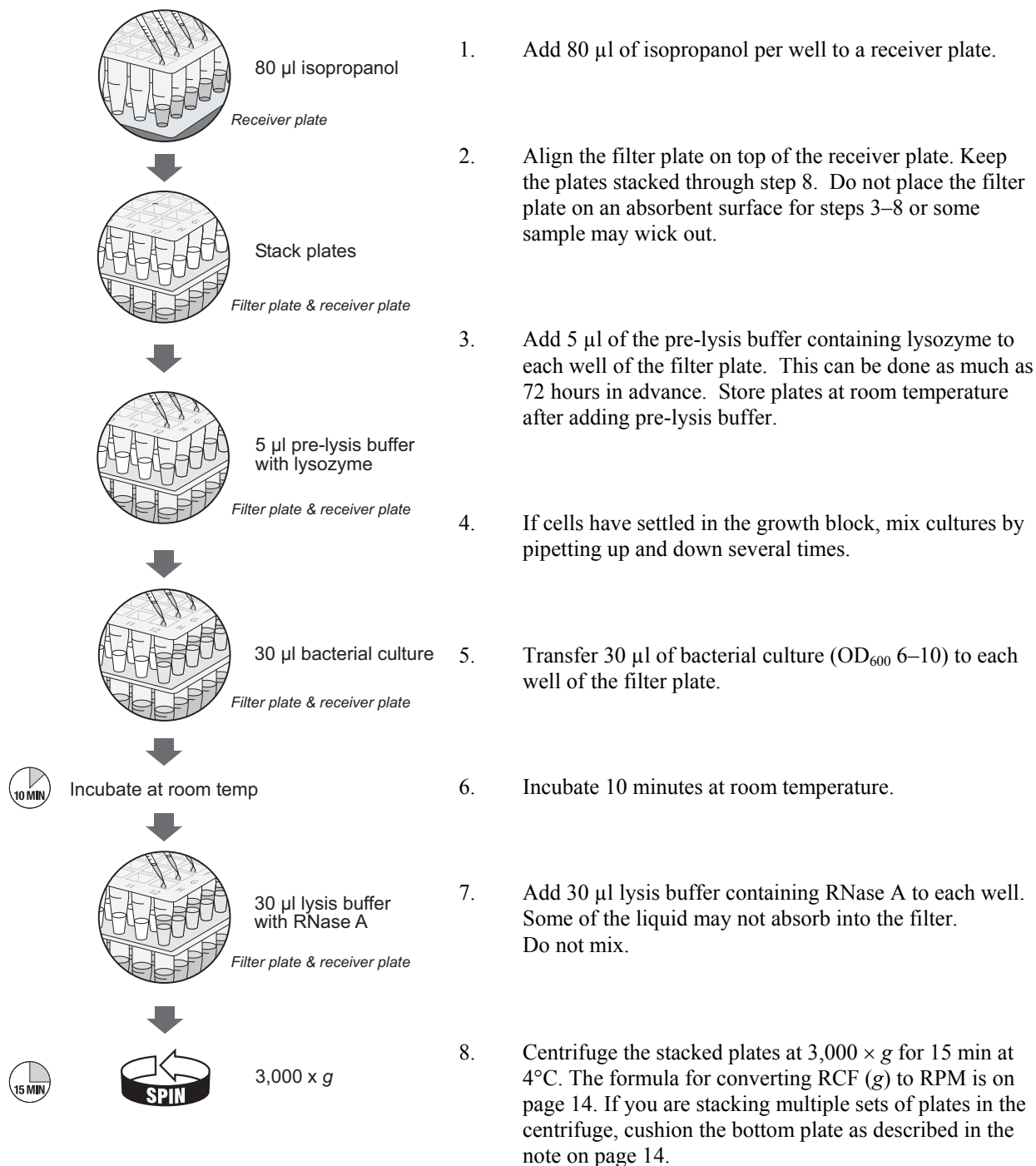
Expected Amount

The expected amount of recovered DNA for Protocol A is **2–3 µg DNA per well**.

Continued on next page

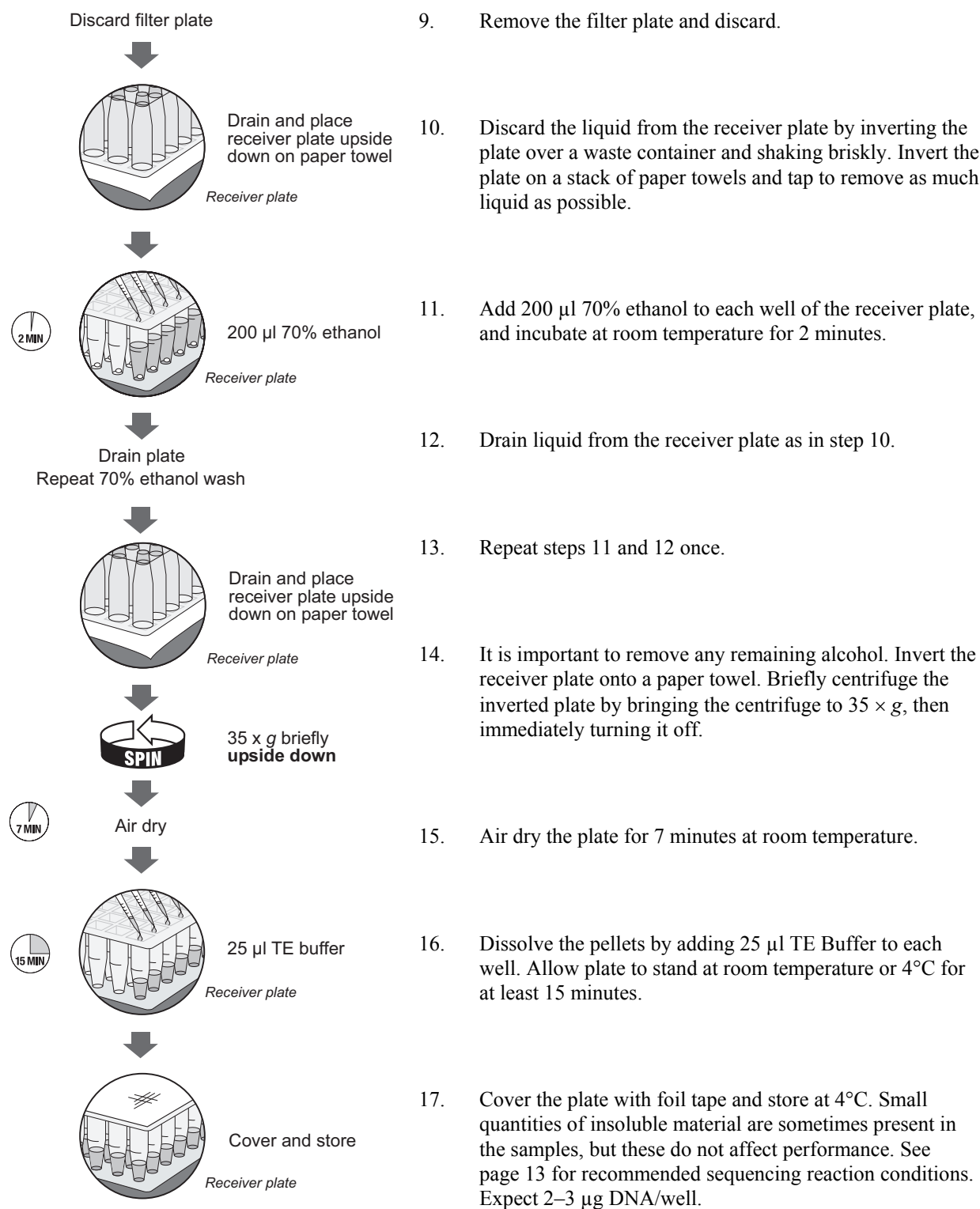
Protocol A, Direct Load Method, Continued

Protocol Procedure



Continued on next page

Protocol A, Direct Load Method, Continued



Protocol B, Harvested Cells Method

Protocol B Overview

In this protocol, you stack the 384-well filter plate on top of a receiver plate containing isopropanol and load pre-lysis buffer containing lysozyme onto the filters. You then resuspend harvested cells to the desired concentration and load them into the wells of the filter plate, followed by lysis buffer. Centrifuge the stacked plates to recover the plasmid DNA and precipitate it in one step. Genomic DNA and other cell components will remain within the filter matrix.

Required Cell Density

For optimal performance, this protocol requires an OD₆₀₀ of **>0.12** for a 1:33 dilution (undiluted OD₆₀₀ >4).

Preparing the Pre-lysis Buffer with Lysozyme

Immediately before using this protocol, follow the steps below to prepare the pre-lysis buffer with lysozyme:

1. Calculate the necessary volume to prepare using the following formula:

$$(\text{total vol pre-lysis buffer}) = [(\# \text{ plates}) \times (2 \text{ ml/plate})] + (\text{waste vol})$$

The waste volume is the volume required to fill the lines or troughs of the robot or pipetting device to be used. The waste volume should be a minimum of 2 ml.

2. Calculate the necessary amount of lysozyme to weigh using the formula below:

$$(\text{total mg lysozyme}) = (\text{total vol pre-lysis buffer}) \times (5 \text{ mg/ml})$$

3. Add the lysozyme to the pre-lysis buffer and invert 30 or more times to thoroughly dissolve.

Note: This solution should be prepared fresh daily. Alternatively, a large batch can be prepared and stored aliquoted for single use at -20°C . Aliquots should be frozen and thawed only once.

Preparing the Lysis Buffer with RNase A

Immediately before using this protocol, prepare the lysis buffer/RNase A mixture according to the directions below:

1. Calculate the necessary volume of lysis buffer and RNase A using the following formulas:

$$[(\# \text{ plates}) \times (10 \text{ ml/plate})] + (\text{waste volume}) = (\text{total vol lysis buffer})$$

$$(\text{total vol lysis buffer}) \times (0.05) = (\text{total vol RNase A})$$

2. Mix the lysis buffer and RNase A by inverting 20 times to mix thoroughly. Use this solution within 24 hours.
-

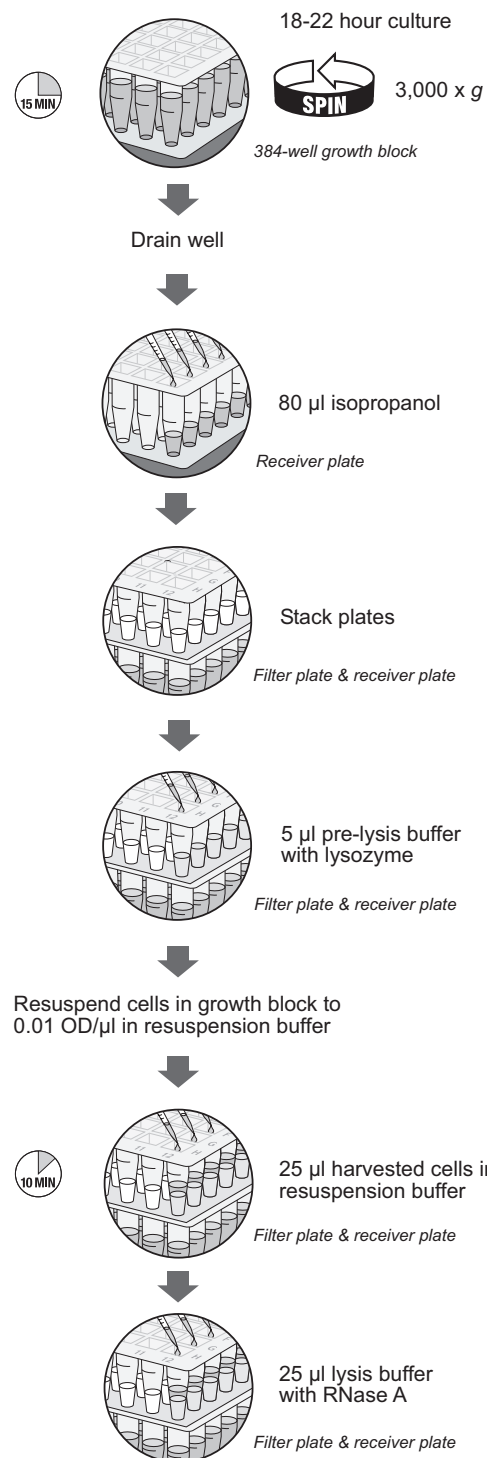
Expected Amount

The expected amount of recovered DNA for Protocol B is **3–5 μg DNA per well**.

Continued on next page

Protocol B, Harvested Cells Method, Continued

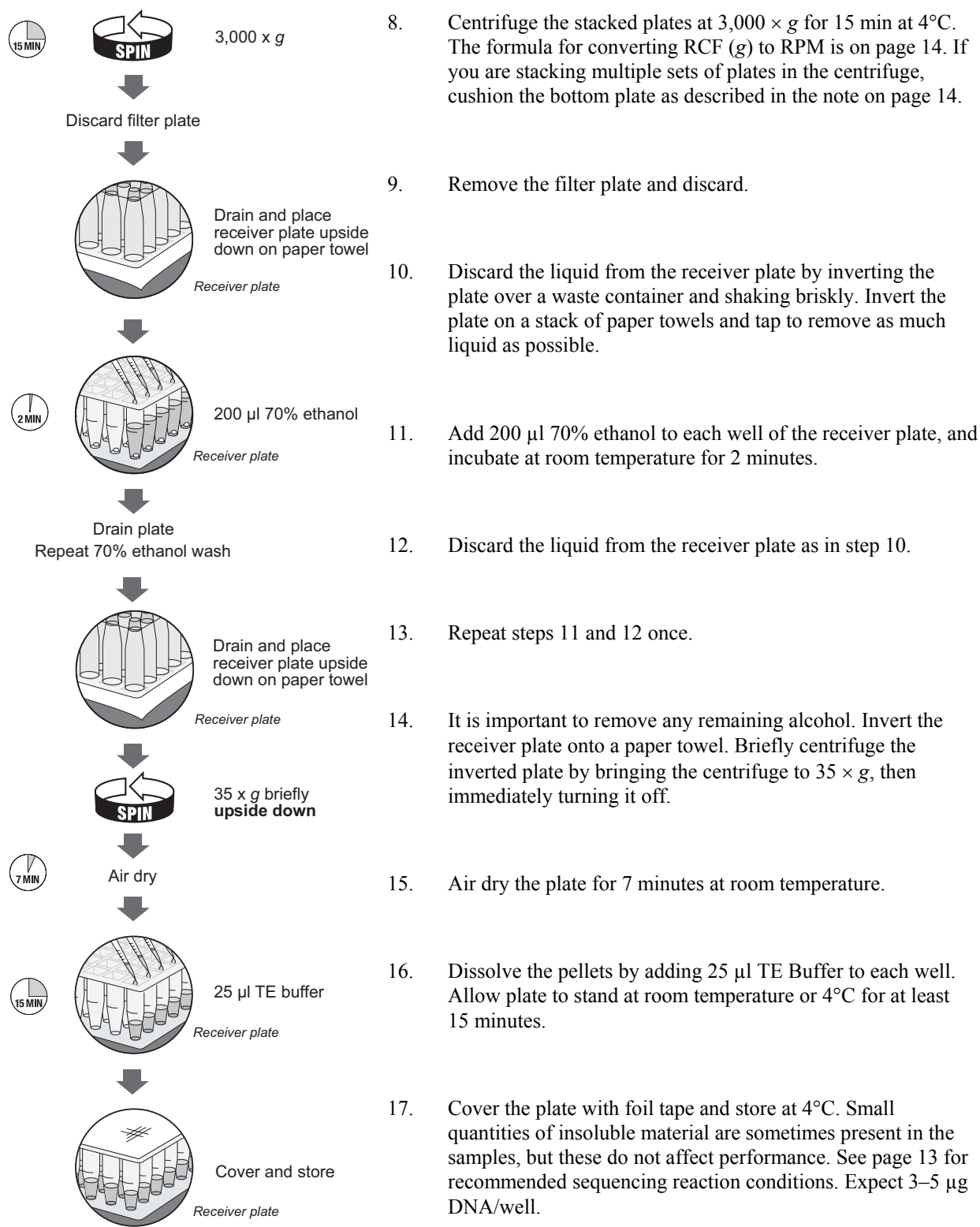
Protocol Procedure



1. Harvest cells from overnight growth by centrifuging growth block at 3,000 $\times$ g for 15 minutes at 4–25°C. Drain well.
2. While cells are pelleting, add 80 μ l of isopropanol to each well of a receiver plate.
3. Align filter plate on top of receiver plate. Keep the plates stacked through step 8. Do not place the filter plate on absorbent surface for steps 3–8 or sample may wick out.
4. Pipet 5 μ l of the pre-lysis buffer with lysozyme onto each filter in the plate. This can be done as much as 72 hours in advance. Store plates at room temperature after adding pre-lysis buffer.
5. Thoroughly resuspend cells in each well of the growth block in cell suspension buffer so that the final cell density will be 0.01 OD/ μ l. Complete resuspension is critical, as clumps of cells will lead to reduced yield. Pipet cells up and down repeatedly or vortex on a plate vortexer to resuspend thoroughly. (**Example:** If an OD₆₀₀ reading of a 1:33 dilution of bacterial culture is 0.21 (7 OD/ml) and there is ~100 μ l of culture in each well (150 μ l media in a 384-well growth block will evaporate to ~100–110 μ l overnight), then there are $7 \times 0.100 = 0.7$ OD of cells in each well. To obtain a cell suspension containing 0.01 OD/ μ l, resuspend cells in 70 μ l of cell suspension buffer.)
6. Transfer 25 μ l of cells into each well of the filter plate (0.25 OD/well). Incubate 10 minutes at room temperature.
7. Add 25 μ l lysis buffer containing RNase A to each well of the filter plate. Some of the liquid may not absorb into the filter. Do not mix.

Continued on next page

Protocol B, Harvested Cells Method, Continued



Troubleshooting Guide

Problem	Possible Cause	Suggested Solutions
Low yield of plasmid DNA	Culture not dense enough	Use fresh colonies, NOT glycerol stocks. Use fresh Terrific Broth or MR2001 media. Use fresh antibiotic at the concentration recommended on page 3. Ensure that the incubator is at 37°C. Ensure that the incubator is shaking at 300–320 RPM. Increase incubation time to as much as 22 hours.
	High copy number plasmid behaves like low copy number plasmid	Use fresh antibiotic at the concentration recommended on page 3 Use DH10B cells to ensure high plasmid DNA production. Ensure that the incubator is at 37°C. Low incubation temperature can reduce plasmid copy number. Ensure that all residual detergent has been rinsed from vessels.
	Insufficient centrifugation	Centrifuge at $3000 \times g$ to recover and precipitate DNA. If your centrifuge cannot reach $3000 \times g$, longer centrifugation times can be used, as specified on page 14. Be sure to calculate g -force using the radius of the rotor with the microplate carrier in place and fully extended.
	Loss of pellet	Centrifuge at 4°C for best adherence of pellet to wall.
Excessive insoluble material	Too many cells used	For Protocol A, Direct Load Method, culture should be grown to an OD_{600} of 6–10. If cultures are growing to a higher density, reduce the incubation time or shaking speed. For Protocol B, Harvested Cells Method, check the density of a few random samples to confirm that the correct number of cells is being used (OD_{600} of 0.25 per well).
	Temperature too high during centrifugation	Ensure that the centrifuge is operating at 4°C during the run, or place the centrifuge in a cold room.
	Excessive centrifugation	Centrifuge at $3,000 \times g$ to recover and precipitate DNA. Do not use higher g -force or longer times.
Weak fluorescence signal in sequencing	DNA pellet insufficiently washed	Two 200 μ l washes with 70% ethanol are recommended in the protocol. Be sure to thoroughly remove the supernatant before each wash to minimize the salt carryover.
	Suboptimal sequencing reaction conditions	See the recommended sequencing reaction conditions on page 13.
Weak PCR Amplification	DNA pellet insufficiently washed	Two 200 μ l washes with 70% ethanol are recommended in the protocol. Be sure to thoroughly remove the supernatant before each wash to minimize the salt carryover.

Continued on next page

Troubleshooting Guide, Continued

Problem	Possible Cause	Suggested Solutions
Chromosomal DNA contamination	Too many cells used	For Protocol A, Direct Load Method, culture should be grown to OD ₆₀₀ of 6-10. If cultures are growing to a higher density, reduce the incubation time or shaking speed. For Protocol B, Harvested Cells Method, check the density of a few random samples to confirm that the correct number of cells are being used (OD ₆₀₀ of 0.25 per well).
RNA contamination	Excessive centrifugation	Centrifuge at 3,000 × g to recover and precipitate DNA. Do not use a higher g-force or longer times.
	Insufficient RNase A	Add RNase A to the lysis buffer immediately before use. Do not store reagent mix for more than one week.

Recommended Sequencing Reaction Conditions

The conditions below are provided as an example; other conditions may be suitable.

1/4 × sequencing reaction:

2.5 µl plasmid DNA from PureLink™ 384 Plasmid Purification System

2.0 µl BigDye Terminator Ready Reaction Mix (ABI)

0.8 pmole primer

1 µl sequencing buffer (see below)

Water sufficient to bring volume to 10.5 µl

Sequencing Buffer:

400 mM Tris-HCl, pH 9.0

10 mM MgCl₂

Note: Omission of the sequencing buffer can result in low fluorescent signal strength.

Centrifuge Information

Compatible Centrifuges and Rotors

The protocols described in this manual are compatible with a wide variety of centrifuges and rotors. Centrifuge and rotor combinations must be able to accommodate a 3.6-cm microtiter plate stack and provide refrigeration.

The protocols also specify a centrifugation speed of $3000 \times g$; however, if your centrifuge is not capable of reaching $3000 \times g$, longer centrifugation times can be used. See Centrifugation at Lower g -Force below for details.



Note

If two or three sets of filter and receiver plates are stacked in one plate carrier in the rotor, a cushion should be placed under the bottom receiver plate to prevent breakage of the outer plate walls. The cushion should be a firm silicon or rubber mat, fit inside the bottom of the receiver plate, and be about 0.5 cm thick.

Conversion Formula

The formula for converting RCF (g) to RPM is:

$$\text{RCF (g)} = 1.12 \times \text{radius in mm} \times (\text{RPM}/1000)^2$$

Where:

RPM is the centrifuge speed in revolutions per minute

RCF is the relative centrifugal force (g)

Radius is the distance (in mm) from the center of the centrifuge spindle to the bottom of the plate in the rotor when the microplate carrier is fully extended as during centrifugation.

Centrifugation at Lower g -Force

If your centrifuge is not capable of reaching $3000 \times g$, longer centrifugation times can be used. See the following table for guidelines:

g-Force	Centrifugation time
1500	30 min
2000	25 min
2500	20 min
3000	15 min

Related Products

Related Products

The following table lists products that support PureLink™ 384 Plasmid Purification Plates:

Product	Size	Cat. No.
PureLink™ 96 Plasmid Purification System	4 plates	12263-018
PureLink™ 96 Plasmid Components:		
PureLink™ 96 Filter Plates	Pkg. of 50	12192-035
PureLink™ 96 Receiver Plates	Pkg. of 50	12193-025
Media:		
Terrific Broth	500 g	22711-022
Glycerol	500 ml	15514-011
Transformation and Testing:		
ElectroMAX™ DH10B™ Cells	5 × 0.1 ml	18290-015
ElectroMAX™ DH10B™ T1 Phage Resistant Competent Cells	5 × 0.1 ml	12033-015
MAX Efficiency® DH5α™ Competent Cells	5 × 0.2 ml	18258-012
MAX Efficiency® DH5α™ T1 Phage Resistant Competent Cells	5 × 0.2 ml	12034-013
CloneChecker System	100 reactions	11666-013
Sequencing:		
Custom Primers (see www.invitrogen.com for information)		

Technical Service

World Wide Web



Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

<http://www.invitrogen.com>

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

Contact Us

For more information or technical assistance, please call, write, fax, or email. Additional international offices are listed on our Web page (www.invitrogen.com).

Corporate Headquarters:

Invitrogen Corporation
1600 Faraday Avenue
Carlsbad, CA 92008 USA
Tel: 1 760 603 7200
Tel (Toll Free): 1 800 955 6288
Fax: 1 760 602 6500
E-mail:
tech_service@invitrogen.com

Japanese Headquarters:

Invitrogen Japan K.K.
Nihonbashi Hama-Cho Park Bldg. 4F
2-35-4, Hama-Cho, Nihonbashi
Tel: 81 3 3663 7972
Fax: 81 3 3663 8242
E-mail: jpinfo@invitrogen.com

European Headquarters:

Invitrogen Ltd
Inchinnan Business Park
3 Fountain Drive
Paisley PA4 9RF, UK
Tel: +44 (0) 141 814 6100
Tech Fax: +44 (0) 141 814 6117
E-mail: eurotech@invitrogen.com

MSDSs

MSDSs are available from our web site: www.invitrogen.com.

Continued on next page



Corporate Headquarters:

*Invitrogen Corporation
1600 Faraday Avenue
Carlsbad, California 92008
Tel: 1 760 603 7200
Tel (Toll Free): 1 800 955 6288
Fax: 1 760 603 7229
Email: tech_service@invitrogen.com*

European Headquarters:

*Invitrogen Ltd
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Tel (Free Phone Orders): 0800 269 210
Tel (General Enquiries): 0800 5345 5345
Fax: +44 (0) 141 814 6287
Email: eurotech@invitrogen.com*

International Offices:

*Argentina 5411 4556 0844
Australia 1 800 331 627
Austria 0800 20 1087
Belgium 0800 14894
Brazil 0800 11 0575
Canada 800 263 6236
China 10 6849 2578
Denmark 80 30 17 40*

*France 0800 23 20 79
Germany 0800 083 0902
Hong Kong 2407 8450
India 11 577 3282
Italy 02 98 22 201
Japan 03 3663 7974
The Netherlands 0800 099 3310
New Zealand 0800 600 200
Norway 00800 5456 5456*

*Spain & Portugal 900 181 461
Sweden 020 26 34 52
Switzerland 0800 848 800
Taiwan 2 2651 6156
UK 0800 838 380
For other countries see our Web site*

www.invitrogen.com