



Instruction Manual

Relay™ 96 Protein Screen

For high-throughput isolation of recombinant proteins from bacterial cells

Catalog no. 12346-011

Version E

042602

25-0448

Table of Contents

General Information.....	1
Introduction	1
Contents	1
Shipping and Storage.....	2
Advantages of Relay™ 96 Protein Screen.....	2
Materials Supplied by the User	2
Product Qualification	2
Bacterial Cells	3
Introduction	3
Preparing Lysis Buffer with Lysozyme	3
Isolating Proteins from Bacterial Cells	4
Isolating Proteins with Low Expression Levels from Bacterial Cells	5
Insect Cells.....	6
Introduction	6
Isolating Proteins from Insect Cells.....	6
Phosphate Buffer	7
Isolating Proteins with Low Expression Levels from Insect Cells	7
Troubleshooting	8
License Information.....	9
Limited Use License Number 94: Protein Screen 96 System	9
Technical Service	10
World Wide Web	10
Contact Us	10
MSDS Requests.....	11
Emergency Information.....	11
Limited Warranty	12

General Information

Introduction

The Relay™ 96 Protein Screen provides a simple and rapid high-throughput method for isolating soluble and insoluble fractions of the recombinant proteins from various types of cell cultures. The protocol eliminates the need to harvest cells and does not require any special equipment.

The Relay™ 96 Protein Screen consists of a 96-well Filter Plate containing a solid phase lysis matrix. To isolate your recombinant protein from a cell culture, add the cell culture directly to the Filter Plate, add lysis buffer, and centrifuge the plate.

After centrifugation, the soluble protein is collected into the Receiver Plate. The cell debris, chromosomal DNA, and insoluble proteins remain trapped in the Filter Plate. The insoluble proteins can be isolated from the same sample using the Denaturing Buffer.

Contents

The components included in Relay™ 96 Protein Screen are listed below. Sufficient reagents are provided in the kit to perform 384 isolations.

Item	Amount	Catalog no.
Relay™ 96 Protein Screen System	Lysis Buffer – 25 ml Denaturing Buffer – 80 ml 96-Well Filter Plates – 4 96-Well Receiver Plates – 8	12346-011
Relay™ 96 Protein Screen Filter Plates	25	12346-045
Relay™ 96 Protein Screen Lysis Buffer	150 ml	12346-029
Relay™ 96 Protein Screen Denaturing Buffer	500 ml	12346-037
Relay™ 96 Protein Screen Receiver Plates	50	12346-052

Continued on next page

General Information, Continued

Shipping and Storage

All components of the Relay™ 96 Protein Screen are shipped at +4°C. Upon receipt, store the denaturing buffer at +4°C and all the other components at room temperature.

Advantages of Relay™ 96 Protein Screen

Using the Relay™ 96 Protein Screen to isolate your recombinant proteins from cell cultures has the following advantages:

- Rapid, inexpensive method designed for high-throughput application
 - Gentle detergent lysis method ensures minimal protein denaturation
 - Designed to isolate soluble and insoluble proteins from the same sample in 30 minutes
 - Minimal contamination of the protein sample with chromosomal DNA
 - Compatible with difficult mucoid strains such as BL21
 - Protein sample suitable for downstream applications such as enzyme assays, ELISA, and affinity chromatography
-

Materials Supplied by the User

- Cell culture containing your recombinant protein
 - Centrifuge capable of centrifuging 96-well plates
 - For bacterial cells, Lysozyme
 - Gels for analyzing protein samples. We recommend using NuPAGE® Novex 10% BisTris Gels (Cat. no. NP0301).
 - For proteins with low expression levels from insect cells, 50 mM phosphate and 100 mM NaCl to use as a buffer solution (see page 7).
-

Product Qualification

Protocols, plates, and buffers were tested using BL21 (DE3) *E. coli*. Presence of isolated proteins was confirmed on NuPAGE® Novex 10% BisTris Gels against Mark12™ Standards.

Bacterial Cells

Introduction

This section provides protocols for isolating soluble and insoluble proteins from bacterial cells.

Preparing Lysis Buffer with Lysozyme

To lyse bacterial cells, you must add Lysozyme to the Lysis Buffer supplied in the kit.

Prepare 6 ml of Lysis Buffer containing Lysozyme in a final concentration of 25 $\mu\text{g/ml}$ as described below for use with one 96-well plate. If you are using more than one 96-well plate, scale up accordingly.

To prepare the appropriate concentration of Lysozyme in Lysis Buffer:

1. Dissolve 25 mg of Lysozyme in 1 ml sterile water
 2. Add 6 μl of this solution to 6 ml Lysis Buffer.
 3. Mix well to achieve a final concentration of 25 $\mu\text{g/ml}$.
 4. Proceed to the following protocol.
-

Continued on next page

Bacterial Cells, Continued

Isolating Proteins from Bacterial Cells

Use the following protocol to recover soluble and insoluble proteins from bacterial cells. Before proceeding, prepare the Lysis Buffer with Lysozyme as described on the previous page.

Do not place the Filter Plate on any absorbent surface during the protocol to prevent the loss of sample

1. If you are using bacterial cells from a growth block, mix cultures by pipetting up and down several times.
2. Place the Filter Plate on top of a Receiver Plate. Transfer 150 μl of induced cell culture to each well of the Filter Plate.
3. Add 50 μl of Lysis Buffer containing Lysozyme to each well of the Filter Plate. All of the Lysis Buffer may not absorb into the filter. Do not mix by pipetting.
4. Incubate the stacked plates at room temperature for 10–15 minutes.
5. Centrifuge the stacked plates at $1000\text{--}2500 \times g$ for 5 minutes at room temperature. You can centrifuge at $+4^\circ\text{C}$ if your protein is sensitive to high temperature.
6. Soluble proteins will be in the Receiver Plate. Remove the Filter Plate and set aside. (Do not place it on an absorbent surface if you want to recover the insoluble proteins.) Analyze 15 μl of the soluble proteins by SDS-PAGE.

After you have isolated the soluble proteins, insoluble proteins will remain in the Filter Plate. Use the following procedure to recover these proteins.

7. Place the Filter Plate containing your insoluble protein on top of a new Receiver Plate.
8. Add 200 μl of Denaturing Buffer to each well of the Filter Plate. All of the Denaturing Buffer may not absorb into the filter.
9. Incubate the stacked plates at room temperature for 10 minutes.
10. Centrifuge the stacked plates at $1000\text{--}2500 \times g$ for 5 minutes at room temperature. You may centrifuge the plates at $+4^\circ\text{C}$ if your protein is sensitive to high temperature.
11. Insoluble proteins will be in the Receiver Plate. Discard the Filter Plate. Analyze 15 μl of the insoluble proteins by SDS-PAGE.

Continued on next page

Bacterial Cells, Continued

Isolating Proteins with Low Expression Levels from Bacterial Cells

The following protocol can be used to isolate proteins with low expression levels from bacterial cells.

Before proceeding, prepare the Lysis Buffer with Lysozyme as described on page 3.

Note: Do not place the Filter Plate on any absorbent surface during the protocol to prevent the loss of sample.

1. Centrifuge the cell culture at $3000 \times g$ or higher for 10 minutes to pellet the cells.
2. Remove some of the media to achieve the desired concentration. You can concentrate the cells up to $OD_{600} = 15$ (higher concentrations may clog the filter).
3. Resuspend the cells in the media by vortexing briefly or pipetting up and down.
4. Place the Filter Plate on top of a Receiver Plate. Add $150 \mu\text{l}$ of cell culture to each well of the Filter Plate.
5. Add $50 \mu\text{l}$ of Lysis Buffer (with Lysozyme) to each well of the Filter Plate and incubate for 10–15 minutes at room temperature. All of the Lysis Buffer may not be absorbed into the filter. Do not mix by pipetting.
6. Centrifuge the stacked plates at $2000 \times g$ for 5 minutes.
7. Soluble proteins will be in the Receiver Plate. Remove the Filter Plate and set aside. (Do not place it on an absorbent surface if you want to recover the insoluble proteins.) Analyze $15 \mu\text{l}$ of the soluble proteins by SDS-PAGE.

After you have isolated the soluble proteins, insoluble proteins will remain in the Filter Plate. Use the following steps to recover these proteins:

8. Place the Filter Plate containing your insoluble protein on top of a new Receiver Plate.
 9. Add $200 \mu\text{l}$ of Denaturing Buffer to each well of the Filter Plate and incubate for 10 minutes.
 10. Centrifuge the stacked plates at $2000 \times g$ for 5 minutes.
 11. Insoluble proteins will be in the Receiver Plate. Discard the Filter Plate. Analyze $15 \mu\text{l}$ of the insoluble proteins by SDS-PAGE.
-

Insect Cells

Introduction

This section provides protocols for isolating soluble and insoluble proteins from insect cells.

Isolating Proteins from Insect Cells

Use the following protocol to recover soluble proteins from insect cells. (Do not add Lysozyme to the Lysis Buffer for insect cells.)

Note: Do not place the Filter Plate on any absorbent surface during the protocol to prevent the loss of sample.

1. Place the Filter Plate on top of a Receiver Plate. Transfer 150 μl of cell culture to each well of the Filter Plate.
2. Add 50 μl of Lysis Buffer to each well of the Filter Plate. All of the Lysis Buffer may not absorb into the filter. Do not mix by pipetting.
3. Incubate the stacked plates at room temperature for 10–15 minutes.
4. Centrifuge the stacked plates at 1000–2500 $\times g$ for 5 minutes at room temperature. You may centrifuge the plates at +4° C if your protein is sensitive to high temperature.
5. Soluble proteins will be in the Receiver Plate. Remove the Filter Plate and set aside. (Do not place it on an absorbent surface if you want to recover the insoluble proteins.) Analyze 15 μl of the soluble proteins by SDS-PAGE.

After you have isolated the soluble proteins, insoluble proteins will remain in the Filter Plate. Use the following steps to recover these proteins:

6. Place the Filter Plate containing your insoluble protein on top of a new Receiver Plate.
7. Add 200 μl of Denaturing Buffer to each well of the Filter Plate. All of the Denaturing Buffer may not absorb into the filter.
8. Incubate the stacked plates at room temperature for 10 minutes.
9. Centrifuge the stacked plates at 1000–2500 $\times g$ for 5 minutes at room temperature. You may centrifuge the plates at +4° C if your protein is sensitive to high temperature.
10. Insoluble proteins will be in the Receiver Plate. Discard the Filter Plate. Analyze 15 μl of the insoluble proteins by SDS-PAGE.

Continued on next page

Insect Cells, Continued

Phosphate Buffer

For Step 3 of the following protocol, prepare a simple Phosphate Buffer of 50 mM phosphate and 100 mM NaCl at pH 8.0. Prepare 150 μ l of buffer per 1 ml of cell culture.

Isolating Proteins with Low Expression Levels from Insect Cells

The following protocol can be used to isolate soluble and insoluble proteins with low expression levels from insect cells.

The amounts in the protocol are per well. Scale up as needed.

Note: Do not place the Filter Plate on any absorbent surface during the protocol to prevent the loss of sample.

1. Centrifuge 1 ml of cell culture at $3000 \times g$ or higher for 10 minutes to pellet the cells.
2. Remove all of the media by aspiration or tapping the container upside down.
3. Add 150 μ l of the Phosphate Buffer described above to the cell pellet.
4. Resuspend the cells in the buffer by vortexing briefly or pipetting up and down.
5. Place the Filter Plate on top of a Receiver Plate. Add 150 μ l of cell solution to each well of the Filter Plate.
6. Add 50 μ l of Lysis Buffer to each well of the Filter Plate and incubate for 10–15 minutes at room temperature. All of the Lysis Buffer may not be absorbed into the filter. Do not mix by pipetting.
7. Centrifuge the stacked plates at $2000 \times g$ for 5 minutes.
8. Remove the Filter Plate. (Do not place it on an absorbent surface if you want to recover the insoluble proteins.) Analyze 15 μ l of the soluble protein in the Receiver Plate by SDS-PAGE.

After you have isolated your soluble proteins, insoluble proteins will remain in the Filter Plate. Use the following steps to recover these proteins:

9. Place the Filter Plate containing your insoluble protein on top of a new Receiver Plate.
 10. Add 200 μ l of Denaturing Buffer to each well of the Filter Plate and incubate for 10 minutes.
 11. Centrifuge the stacked plates at $2000 \times g$ for 5 minutes.
 12. Discard the filter plate. Analyze 15 μ l of the insoluble protein in the receiver plate by SDS-PAGE.
-

Troubleshooting

Problem	Cause	Solution
Clogged filter	Cell density is too high	Reduce the cell density. Avoid loading bacterial cultures with an $OD_{600} > 15$.
Low protein yield	Bacterial cells—poor induction of expression	If you are using a 96-well growth block, induction may be difficult and slow; we recommend growing bacterial cell cultures in 24- or 48-well growth blocks for optimal expression.
	Low protein expression	For proteins with low expression levels, use the Isolating Proteins with Low Expression Levels from Bacterial Cells protocol on page 5 or the Isolating Proteins with Low Expression Levels from Insect Cells protocol on page 7.
	Low protein expression or protein may be toxic to cells	Optimize growth conditions or try a different expression system to increase protein expression.

License Information

Limited Use License Number 94: Protein Screen 96 System

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for commercial purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for the commercial purposes of the buyer, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for commercial purposes. Commercial purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned by Invitrogen based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the products with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 1600 Faraday Avenue, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500.

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
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK

Tel: +44 (0) 141 814 6100

Tel (Toll Free): 0800 5345 5345

Fax: +44 (0) 141 814 6287

E-mail:

eurotech@invitrogen.com

Continued on next page

Technical Service, Continued

MSDS Requests

To request an MSDS, visit our Web site (www.invitrogen.com).

1. On the home page, go to the left-hand column under 'Technical Resources' and select 'MSDS Requests'.
 2. Follow instructions on the page and fill out all the required fields.
 3. To request additional MSDSs, click the 'Add Another' button.
 4. All requests will be faxed unless another method is selected.
 5. When you are finished entering information, click the 'Submit' button. Your MSDS will be sent within 24 hours.
-

Emergency Information

In the event of an emergency, customers of Invitrogen can call the 3E Company, 24 hours a day, 7 days a week for disposal or spill information. The 3E Company connects the customer with poison control or with the University of California at San Diego Medical Center doctors. 3E Company, Voice: 1-760-602-8700

Technical Service, Continued

Limited Warranty

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives.

Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

©2002 Invitrogen Corporation. All rights reserved.

Notes:

Notes:



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 website*

www.invitrogen.com

