

USER GUIDE

novex
by *life* technologies

E-PAGE™ 48 Protein Electrophoresis System

For electrophoresis of 48 protein samples

Catalog Number EP048-08

Publication Part Number 25-0735

Publication Number MAN0000374

Revision A.0

For Research Use Only. Not for use in diagnostic procedures.

life
technologies

For Research Use Only. Not for use in diagnostic procedures.

Information in this document is subject to change without notice.

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

IMPORTANT LICENSING INFORMATION

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

TRADEMARKS

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Contents

E-PAGE™ 48 Experienced Users Procedure.....	iv
Product Contents.....	1
Introduction.....	2
E-PAGE™ Protein Electrophoresis System.....	2
Methods.....	5
Guidelines for Sample Preparation.....	5
Electrophoresis of E-PAGE™ 48 Gels.....	10
Guidelines for Loading E-PAGE™ Gels by Automated Liquid Handling.....	15
Maintenance of the E-PAGE™ Device.....	16
Opening the E-PAGE™ Cassette.....	17
Visualizing and Staining of E-PAGE™ Gels.....	18
Visualizing Lumio™ Fusion Proteins.....	18
Coomassie Staining.....	19
Blotting of E-PAGE™ Gels.....	21
Semi-Dry Blotting of E-PAGE™ Gels.....	21
Expected Results.....	24
Using E-Editor™ 2.02 Software.....	26
Troubleshooting.....	27
Appendix.....	29
Product Specifications.....	29
Technical Support.....	32

Product Contents

Types of kits

This manual is supplied with the following product:

Product	Quantity	Catalog no.
E-PAGE™ 48 8% Gels	8-pack	EP048-08

Kit contents

Each pack of E-PAGE™ Gels contain the following contents:

E-PAGE™ Gel 8-Pack Contents	Quantity
E-PAGE™ Gels	8 gels
E-PAGE™ Loading Buffer 1 (4X)	4.5 mL
E-PAGE™ Blotting Pad	1

Shipping and storage

The E-PAGE™ 48 and E-PAGE™ Blotting Pad are shipped and stored at room temperature.

Do not allow the temperature to drop below 4°C or rise above 40°C when storing the gels.

The E-PAGE™ 48 8% Gels are guaranteed stable for 6 months from the date of production when stored properly. The expiration date is printed on the package.

The E-PAGE™ Loading Buffer is shipped at room temperature. Store the E-PAGE™ Loading Buffer 1 (4X) at room temperature or at 4°C.

Introduction

E-PAGE™ Protein Electrophoresis System

About the E-PAGE™ Protein Electrophoresis System

The E-PAGE™ Protein Electrophoresis System is designed for fast, medium-to-high-throughput protein electrophoresis in a horizontal format.

The E-PAGE™ 48 System consists of the following components:

- E-PAGE™ 48 8% Pre-cast Gels
 - E-Base™ Electrophoresis Device
 - E-PAGE™ Loading Buffer 1 (4X)
 - E-PAGE™ Blotting Pad
 - E-Editor™ 2.02 Software
-

Applications

The E-PAGE™ 48 Protein Electrophoresis System is ideal for protein analysis and screening including:

- Coomassie R-250 staining (page 19)
 - In-gel staining with Lumio™ Green Reagent (page 18)
 - Silver staining and SYPRO® Ruby staining techniques (protocols available in the E-PAGE™ Technical Guide)
 - Western blotting
 - Functional assays
-

E-PAGE™ 48 Gels

- **IMPORTANT!** E-PAGE™ 48 Gel are **not compatible** with the E-Gel® 96 base (Cat nos. G7100-01/G7200-01, previously available from Invitrogen). The older E-Gel® 96 bases do not have the 'E-Base™' inscription on the platform.
- E-PAGE™ 48 Gels are self-contained, pre-cast gels that include a buffered gel matrix and electrodes packaged inside a disposable, UV-transparent cassette.
- Each E-PAGE™ 48 Gel contains 48 sample lanes and 4 marker lanes. This configuration provides a 3.2-cm run length. The wells of E-PAGE™ 48 Gels are compatible for loading with a multichannel pipetter in alternating lanes or with an automated liquid handling system (see page 15 for automation specifications).
- After electrophoresis, the E-PAGE™ 48 cassette is easily opened with a Gel Knife to remove the gel for staining or blotting applications.
- Each E-PAGE™ 48 cassette is labeled with an individual barcode to facilitate identification of the gel using commercial barcode readers (page 10).

See page 29 for E-PAGE™ 48 8% Gel specifications.

E-PAGE™ Protein Electrophoresis System, Continued

Diagram of an E-PAGE™ 48 Cassette

A diagram of the E-PAGE™ 48 Gel cassette is shown below.



Separation range

E-PAGE™ 48 8% Gels have a unique separation profile, which gives protein resolution similar to that of a 4-12% Tris-Glycine gel. Specifically, the following table describes the separation range of the E-PAGE™ 48 8% Gels (see page 29):

10% Resolution Range	30% Lower Resolution Range	30% Upper Resolution Range
50-80 kDa	10-50 kDa	80-200 kDa

About the E-Base™ System

E-PAGE™ 48 Gels are used with a specially designed electrophoresis device that combines a base and a power supply. Two types of devices are available from Life Technologies:

- The **Mother E-Base™** Device (Cat. no. EB-M03) has an electrical plug that can be connected directly to an electrical outlet and is used for electrophoresis of one E-PAGE™ 48 Gel.
Note: The Mother E-Base™ Device has been tested for electrophoresis with up to three Daughter E-Bases™ connected at one time.
- The **Daughter E-Base™** Device (Cat. no. EB-D03) connects to the Mother E-Base™, and together they can be used for the independent electrophoresis of 2 or more E-PAGE™ 48 Gels.
Note: The Daughter E-Base™ Device does not have an electrical plug and cannot be used without a Mother E-Base™ Device.

E-PAGE™ Protein Electrophoresis System, Continued

E-PAGE™ Loading Buffer

E-PAGE™ Gels are supplied with E-PAGE™ Loading Buffer 1 (4X), which is optimized for E-PAGE™ Gels when performing routine SDS-PAGE and staining or blotting applications. **Do not use any other SDS-PAGE sample buffer.** See page 7 for more information about loading buffers.

E-PAGE™ Blotting Pad

E-PAGE™ Blotting Pads are supplied with the E-PAGE™ 48 Gels. It is necessary to use the pad during semi-dry blotting to ensure a good transfer. The pad is reusable as long as the pad retains porosity and liquid retaining capacity.

E-Editor™ 2.02 Software

The E-Editor™ 2.02 software allows you to quickly reconfigure digital images of E-PAGE™ Gel results for analysis and documentation.

Capture an image of the gel and use the E-Editor™ 2.02 software to:

- Align and arrange the lanes in the image.
- Save the reconfigured image for further analysis.
- Copy and paste selected lanes or the entire image into other applications for printing, saving, emailing, and/or publishing.

The E-Editor™ 2.02 software can be downloaded FREE at www.lifetechnologies/epage. Follow the instructions to download the software and user manual.

Methods

Guidelines for Sample Preparation

Introduction

For optimal results using the E-PAGE™ System, follow the guidelines for preparing your protein samples as described in this section.

Prepare your protein samples as described in this section for electrophoresis on E-PAGE™ 48 Gels.

We recommend that you read this section before preparing your samples. Use the sample preparation method and loading buffer appropriate for your detection method:

Application	Method	Loading Buffer
Routine staining and western blotting	1 (page 8)	Loading Buffer 1
Lumio Green Detection	2 (page 9)	Lumio™ Gel Sample Buffer

Materials needed

The following items are needed for sample preparation. See page 31 for ordering information.

- Protein sample
 - NuPAGE® Sample Reducing Agent (10X)
 - E-PAGE™ Loading Buffer 1 (4X) included in the kit
 - Deionized water
 - Heating block set at 70°C
 - Molecular weight markers (page 7)
 - (Optional) Lumio™ Green Detection Kit for detection of Lumio™ fusion proteins (page 9)
-



Note

E-PAGE™ Gels contain SDS and are designed for performing electrophoresis under denaturing conditions.

To obtain the best results, perform SDS-PAGE under reducing conditions.

If you need to perform SDS-PAGE under non-reducing conditions, **do not** add NuPAGE® Sample Reducing Agent (10X) during sample preparation.

Guidelines for Sample Preparation, Continued

Total sample volume

Avoid loading less than 5 μL of sample in wells, and maintain a uniform loading volume. If you do not have enough samples to load all the wells of the gel, load an equal volume of deionized water into any empty wells.

E-PAGE™ Gel Type	Recommended Loading Volume
E-PAGE™ 48 8% Gel	20 μL

We recommend loading 5–10 μL of deionized water into all wells of the E-PAGE™ Gel prior to loading samples or molecular weight markers.

Amount of protein

Load up to 20 μg of protein per well of the E-PAGE™ Gel. The amount of protein required depends on the staining or western detection method used for visualization. If you are unsure of how much protein to use, test a range of concentrations to determine the optimal concentration for your sample.

Note: To ensure a proper LDS (lithium dodecyl sulfate from Loading Buffer 1) to protein ratio, limit sample protein or lipid (from the sample) amount to 2 $\mu\text{g}/\mu\text{L}$ of the final sample volume. Excess protein will cause poor resolution.

High salt or detergent samples

Samples containing high salt or detergents result in loss of resolution on E-PAGE™ Gels. Dilute the samples such that the final concentration of the salt or detergent in the sample is as described below. NT= not tested.

Detergent or Salt	Final Concentration for E-PAGE™ 48Gels
Triton® X-100	<0.3%
Tween® 20	<0.3%
CHAPS	<0.3%
NP-40	<0.3%
RIPA	<0.25X
SDS	<2% (already in loading buffer)
Tris	<300 mM
NaCl	<300 mM
Ammonium sulfate	<100 mM
Sodium acetate	<200 mM
EDTA	<20 mM
MES	Not recommended
DTT, Glycine, Urea, Imidazole	No effect seen up to 500 mM

Guidelines for Sample Preparation, Continued

Loading buffer

Based on your application, use the appropriate loading buffer as described below:

- **SDS-PAGE and routine staining (Method 1, page 8)**
For SDS-PAGE and routine staining or blotting, use the E-PAGE™ Loading Buffer 1 (4X) included in the kit for preparing samples. The E-PAGE™ Loading Buffer 1 (4X) is optimized for E-PAGE™ Gels. **Do not** use any other types of SDS-PAGE sample buffer.
 - **SDS-PAGE and detection of Lumio™ Fusion Proteins (Method 2, page 9)**
For in-gel detection of Lumio™ fusion proteins with the Lumio™ Green Detection Kit, use the Lumio™ Gel Sample Buffer (4X) included with the Lumio™ Green Detection Kit. This buffer is specifically formulated to provide optimal results with the Lumio™ Green Detection Reagent. **Do not** use the E-PAGE™ Loading Buffer 1 (4X) or any other type of SDS-PAGE buffer to prepare samples for Lumio™ Green detection.
-

Molecular weight standards

The following protein molecular weight standards and loading volumes are recommended for E-PAGE™ Gels. Because molecular weight standards have different migration patterns in different gel systems, see page 24 for apparent molecular weight calibration on E-PAGE™ Gels. See page 31 for ordering information.

Gel Type	Standard	Amount	Application
E-PAGE™ 48 8% Gel	SeeBlue® Plus2 Pre-stained Standard	5 µL	Electrophoresis
	MagicMark™ XP Western Protein Standard	10 µL	Electrophoresis, followed by staining
	MagicMark™ XP Western Protein Standard	5 µL	Western blotting
	BenchMark™ Fluorescent Protein Standard	5 µL	Fluorescent Detection

Guidelines for Sample Preparation, Continued

Method 1: Routine staining and blotting

Use this protocol if you are performing SDS-PAGE followed by routine staining or blotting.

If the E-PAGE™ Loading Buffer 1 (4X) is stored at 4°C, bring the buffer to room temperature and mix briefly prior to use.

1. Prepare your samples in a total volume of 10 µL in the E-PAGE™ Loading Buffer 1 (4X) as described below. If you need to prepare samples in a volume of 5–15 µL, adjust the volumes accordingly.

Reagent	Reduced	Non-reduced
Protein Sample	x µL	x µL
E-PAGE™ Loading Buffer 1 (4X)	2.5 µL	2.5 µL
NuPAGE® Sample Reducing Agent (10X)	1 µL	—
Deionized Water	to 10 µL	to 10 µL

2. Incubate the samples at 70°C for 10 minutes.
 3. Proceed to **Loading E-PAGE™ Gels**, page 12.
-

Guidelines for Sample Preparation, Continued

Method 2: Lumio™ detection

A brief protocol to prepare samples for specific detection of Lumio™ fusion proteins using the Lumio™ Green Detection Kit is described below.

For details on the Lumio™ Green Detection Kit, refer to the manual available at www.lifetechnologies.com or contact Technical Support (page 32).

1. Refer to the Lumio™ Detection manual for details on each type of protein. Prepare protein samples as follows:

Protein Sample	Sample Volume	Lumio™ Gel Sample Buffer (4X) Volume
Bacterial samples	7.5 µL	2.5 µL
Mammalian lysate	7.5 µL	2.5 µL
Partially purified sample	7.5 µL	2.5 µL
Purified sample	7.5 µL	2.5 µL
<i>In vitro</i> expressed	10 µL	Not needed*

*There is no need to add Lumio™ Gel Sample Buffer (4X), as the sample is already prepared in this buffer.

2. Thaw the Lumio™ Green Detection Reagent and mix well.
3. Add 0.1 µL Lumio™ Green Detection Reagent to the protein samples from Step 1 in a fume hood. Mix well. Return the Lumio™ Green Detection Reagent to –20°C immediately after use.
4. Incubate the samples at 70°C for 10 minutes.
5. Allow samples to cool for 1–2 minutes and centrifuge briefly at maximum speed in a microcentrifuge.
6. Thaw the Lumio™ In-Gel Detection Enhancer and mix well. Add 1 µL of Lumio™ In-Gel Detection Enhancer to the samples.
7. Mix well and incubate the samples at room temperature for 5 minutes. Return the Lumio™ In-Gel Detection Enhancer to –20°C immediately after use.
8. Proceed to **Loading E-PAGE™ Gels**, page 10.



Note

When performing electrophoresis of Lumio™ fusion proteins, extending the run time of the gel for an additional 2 minutes can prevent the formation of a fluorescent dye front.

Electrophoresis of E-PAGE™ 48 Gels

Introduction

After preparing your samples, you are ready to load E-PAGE™ Gels. This section describes the procedure for loading protein samples and molecular weight standards.

The Mother E-Base™ and Daughter E-Base™ Electrophoresis Devices are designed to fit most robotic platforms allowing you to load and run E-PAGE™ Gels directly on the automated liquid handling system.

If you are using an automated liquid handling device, it is important to align the robotic tip loading assembly to the proper setting prior to loading samples on the E-PAGE™ 48 Gel. This ensures proper loading of samples into the wells. See page 15 for automation guidelines.

If you need to load multiple gels on a robotic platform while other gels are running on the E-Base™ Electrophoresis Devices, use an E-Holder™ Platform (page 15).

General guidelines

- Perform manual loading with a pipette or multichannel pipetter (load samples into alternate wells of the gel followed by a second round of loading into the remaining wells).
 - Use short, rigid tips for loading E-PAGE™ Gels.
 - **Always load 5–10 µL deionized water first into all wells prior to loading sample or molecular weight standard.**
 - E-PAGE™ Gels can only be used once. **Do not re-use.**
 - To obtain the best results, run the E-PAGE™ Gel immediately after removal from the pouch and loading.
 - Store and run E-PAGE™ Gels at room temperature.
 - For optimal results, do not run reduced and non-reduced samples on the same gel. If you do choose to run these samples on the same gel, avoid running reduced and non-reduced samples in adjacent lanes as the reducing agent may have a carry-over effect on the non-reduced samples.
 - Avoid running samples containing different salt or protein concentrations in adjacent lanes.
 - Avoid touching the gel during electrophoresis.
-

Using the barcode

Each E-PAGE™ 48 Gel cassette is labeled with an individual barcode. The barcode facilitates the identification of each gel cassette during electrophoresis of multiple gels. Each E-PAGE™ cassette contains an EAN 39 type of barcode, which is recognized by the majority of commercially available barcode readers. Refer to the manufacturer's instructions to set up the barcode reader.

Note: When capturing an image of an E-PAGE™ Gel, note that the barcode label is easily overexposed. To ensure that the barcode label is distinct and readable in the image, experiment with different shutter settings for your particular camera.

Electrophoresis of E-PAGE™ 48 Gels, Continued

Select a program on the E-Base™ Device

Use **program EP** for running E-PAGE™ 48 Gels. Select the program **prior** to inserting a gel into the base.

Brief instructions for using the E-Base™ Device are included in this manual. For details, refer to the E-Base™ manual.

1. Plug Mother E-Base™ Device into an electrical outlet using the electrical plug on the base. Connect a Daughter E-Base™ Device to a Mother E-Base™ Device or another Daughter E-Base™ Device connected to a Mother E-Base™ Device.
2. The display shows EP or last program used (EG or EP).
3. Press and release the **pwr/prg** button to select **program EP**.
4. Press and release the **time** button to view the time setting. The default run time for program EP is 14 minutes.
5. Press and hold the **time** button to increase the time to **25 minutes**. If the time button is not released, the time setting increases until it reaches 00. To begin cycling through the numbers again, starting from 00, press the time button again.

Do not run E-PAGE™ 48 Gels for more than 30 minutes.



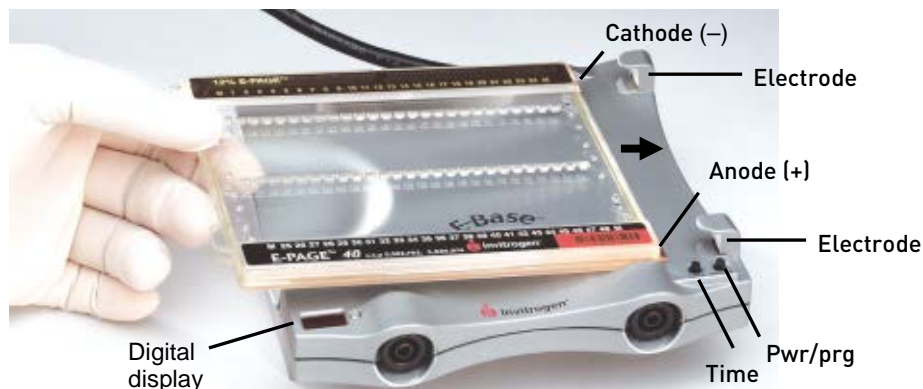
Note

- When performing electrophoresis of Lumio™ fusion proteins, extend the run time of the gel for an additional 2 minutes to prevent the formation of a fluorescent dye front.
 - If your sample contains high salt or detergent concentrations, you may need to manually increase the run time.
 - To increase the run time when a cassette is inserted, press and release the time button to increase the time setting by 1-minute intervals or press and hold the time button to increase the time continuously.
 - To increase the run time while a run is in progress, see next page. To manually interrupt or stop a run, see page 14.
-

Electrophoresis of E-PAGE™ 48 Gels, Continued

Insert the gel cassette

1. Open the package and remove the gel.
2. Remove the plastic combs from the gel.
3. Slide the gel into the Mother E-Base™ or Daughter E-Base™ Device. The two copper electrodes on the right side of the gel cassette must be in contact with the two electrode connections on the base, as shown below.



When the gel is properly inserted into the base, a fan in the base begins running, and a red light illuminates at the lower-left corner of the base. The digital display shows the appropriate time for a selected program or the last time setting (Ready Mode).

Note: If a cassette is inserted into the base before selecting program EP, remove the cassette, select program EP, and reinsert the cassette.

Load the gel

There should be a uniform volume of liquid in each well. The recommended total volume for each well is 20 μ L.

1. Load 5–10 μ L deionized water to each well of the E-PAGE™ Gel prior to loading your samples or protein molecular weight standard.
 2. Load 10–5 μ L of samples in loading buffer into the wells using a multichannel pipetter or an automated liquid handling system.
 3. Load the appropriate protein molecular weight standards in the marker wells of the gel. See page 7 for recommended molecular weight standards.
-

Electrophoresis of E-PAGE™ 48 Gels, Continued

Perform electrophoresis with the E-Base™ Device

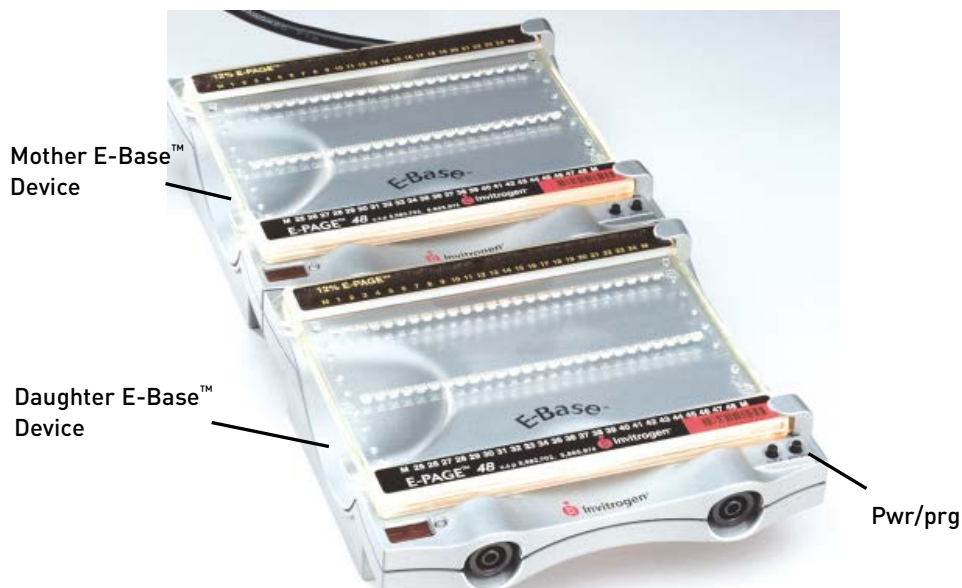
Note: It is not necessary to have a gel in the Mother E-Base™ Device if you are using a Daughter E-Base™ Device. However, the Mother E-Base™ Device must be plugged into an electrical outlet.

1. Press and release the **pwr/prg** button to begin electrophoresis.

The **red light** changes to **green** and the digital display shows the count down time while the run is in progress.



On a Daughter E-Base™ Device, press and release the **pwr/prg** button on the Daughter E-Base™ Device to begin electrophoresis.



To increase the run time while the run is in progress, press the time button to select the desired time and then release the time button.

Do not run an E-PAGE™ 48 Gel for more than 30 minutes.

To interrupt or stop a run in progress, see page 14.

Electrophoresis of E-PAGE™ 48 Gels, Continued

Perform electrophoresis with the E-Base™ Device, continued

2. The Mother E-Base™ or Daughter E-Base™ Device signals the end of the run with a flashing red light and an audible alarm that beeps rapidly for 2 minutes, followed by a single beep every minute.
The digital display shows the original time setting (it does not indicate time changes that were made during electrophoresis). The digital display also shows the elapsed time since the end of the run (up to 19 minutes).
3. **Press and release** the pwr/prg button to stop the beeping. The light turns to a steady red and the digital display shows the last time setting.
4. Remove the gel cassette from the Mother E-Base™ or Daughter E-Base™ Device.
5. For visualizing Lumio™ fusion proteins, see page 18. For opening the E-PAGE™ cassette, see page 17.

Note: The bands in the gel will diffuse within 40 minutes.

Interrupting electrophoresis

You can interrupt an electrophoresis run at any time by **pressing and releasing** the pwr/prg button to stop the current. The stopped current is indicated by a steady red light and the digital display will flash to indicate that the run was interrupted.

You can remove the gel from the E-Base™ Device to check the progress of the run. Then:

- To **continue** the run from the point at which it was stopped, reinsert the gel and press and release the pwr/prg button. The light changes to steady green and the digital display shows the count down time.
- To **cancel** the rest of the interrupted run, press and hold the pwr/prg button for a few seconds. The digital display will reset and the base will return to Ready Mode. If desired, you can then program a new run time as described on page 11 and rerun the gel.

In case of an **external power failure** (loss of electricity or the electrical cord is accidentally removed from the outlet), the run will continue when the power resumes. The Mother E-Base™ or Daughter E-Base™ Device signals the end of the run as described on the previous page, except the light will be an alternating red/green to indicate that an external power failure had occurred during the run.

Guidelines for Loading E-PAGE™ Gels by Automated Liquid Handling

Automated loading of E-PAGE™ 48 Gels

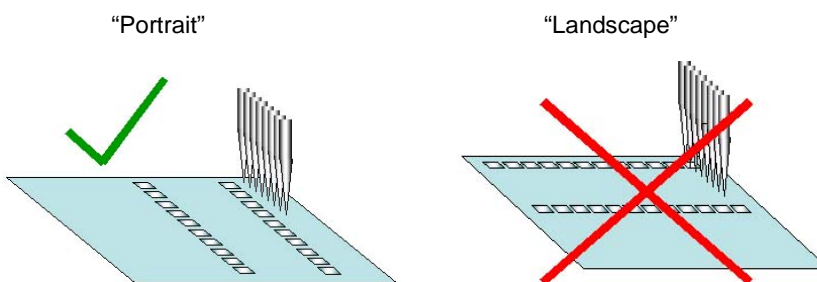
E-PAGE™ 48 Gels are compatible with any automated liquid handling system that has an 8-span loading head with either fixed (9-mm) or variable distance between the loading heads. These loading patterns are described below.

To download programming scripts for your automated liquid handling system, go to www.lifetechnologies.com/epage.



Important

For automated loading of E-PAGE™ 48 Gels, position the plate in the “Portrait” orientation rather than the “Landscape” position, as shown below:

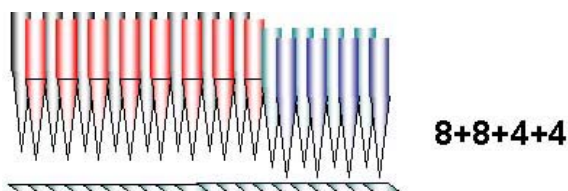


This orientation is available for some automation systems with the use of a 90° adapter. If your system does not have an adapter that allows the “Portrait” configuration, please contact Technical Support (page 32) to obtain a 90° adapter.

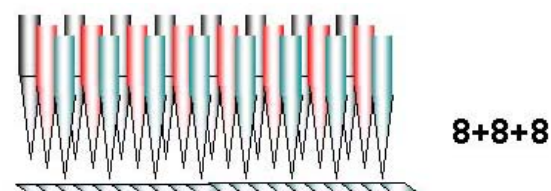
Loading patterns for E-PAGE™ 48 Gels

Loading E-PAGE™ 48 Gels requires one of the following loading patterns, depending on your machine:

Fixed Tip (4 movements per row: 8+8+4+4)



Variable Tip (3 movements per row: 8+8+8)



Continued on next page

Maintenance of the E-PAGE™ Device

Maintaining the E-Base™ Device

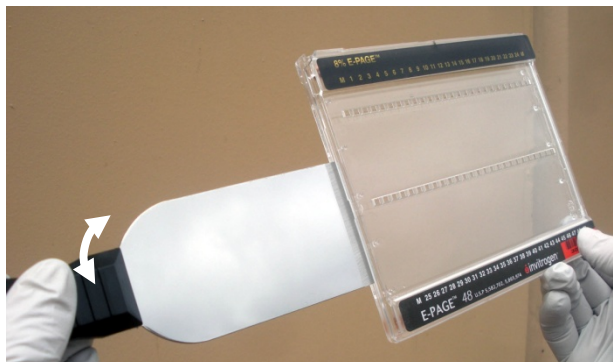
Keep the surfaces of the Mother E-Base™ Device and Daughter E-Base™ Device free of contaminants. To clean, disconnect bases from power source and wipe with a dry cloth. Do not attempt to open or service the bases. To honor the warranty, bases should only be opened and serviced by Life Technologies. Disconnect the Mother E-Base™ Device from the outlet when not in use for a prolonged period of time.

Opening the E-PAGE™ Cassette

Opening the E-PAGE™ 48 cassette with the Gel Knife

To remove the E-PAGE™ 48 Gel from the cassette for blotting or staining, use the Gel Knife to open the cassette.

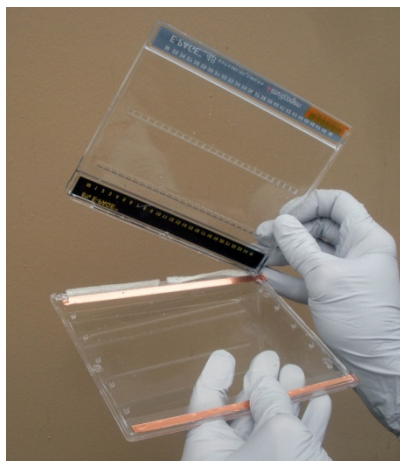
1. Separate each of the bonded sides of the E-PAGE™ 48 cassette by inserting the Gel Knife into the gap between the two plastic plates that make up the cassette. The side of the cassette with the wells should face up.



2. Lever the handle of the knife gently to separate the plates. Repeat on each side of the cassette until the plates are completely separated.

Caution: Use caution while inserting the Gel Knife between the two plates to avoid excessive pressure on the gel.

3. Gently pull apart the cassette halves with your hands until the cassette halves are completely separated and the gel is exposed.



4. Carefully remove the gel from the cassette.
5. Use the Gel Knife to trim the top and bottom electrode areas of the gel.
6. Proceed to blotting (page 21) or staining (pages 18).

Note: Small pieces of gel material may remain in the wells of an E-PAGE™ Gel after removal of the gel from the cassette. To obtain the best staining or blotting results, remove any small pieces of gel material in the wells of the gel by gently rubbing a gloved hand over the well side of the gel.

Visualizing and Staining of E-PAGE™ Gels

Visualizing Lumio™ Fusion Proteins

Introduction

The steps involved in detecting Lumio™ fusion proteins run on an E-PAGE™ Gel are described below.

To visualize Lumio™ fusion protein bands after electrophoresis, you will need a UV transilluminator or a laser-based scanner (see below). For further details on imaging Lumio™ fusion proteins, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 32)

Visualizing Lumio™ fusion proteins

After electrophoresis is complete, immediately visualize and image the gel as described below. **There is no need to remove the E-PAGE™ Gel from the cassette to visualize Lumio™ fusion proteins.**

1. Place the gel cassette on a UV transilluminator equipped with a camera and select the ethidium bromide or SYBR® Green filter on the camera.

You may also use a laser-based scanner with a laser line that falls within the excitation maxima of the stain (500 nm), and a 535 nm long pass filter or a band pass filter centered near the emission maxima of 535 nm.

Note: Adjust the settings on the camera prior to turning on the UV transilluminator. Avoid exposing the gel to UV light for long periods of time.

2. Image the gel with a suitable camera with the appropriate filters using a 4–10 second exposure. You may need to adjust the brightness and contrast to reduce any faint non-specific bands.

You should see fluorescent bands of Lumio™ fusion proteins and the gel should have minimal background, as shown on page 25.

Coomassie Staining

Introduction

Instructions for staining E-PAGE™ Gels using Coomassie R-250 are described in this section.

To obtain maximum sensitivity, the total staining time for Coomassie R-250 staining is 1.5 h plus overnight destaining. The use of the Coomassie R-250 microwave protocol reduces the amount of time needed for staining and destaining, however for minimal background, overnight destaining is recommended.

Materials needed

You will need the following items for staining one E-PAGE™ Gel:

- Clean staining containers or incubation trays (if using the Coomassie R-250 microwave protocol, make sure the container is **microwave safe**. Do NOT use the Incubation Tray cat no. LC2102)
- Rotary shaker

For Coomassie R-250 Staining

- Coomassie R-250 Stain (see note below)
 - Destaining solution (8% acetic acid in deionized water, see note below)
 - Methanol (regular protocol only)
 - Ethanol (microwave protocol only)
 - 2 pieces of nylon membrane (microwave protocol only)
-



The volume of fixing, staining and destaining solutions will depend on the volume of your staining container. To obtain good results, the volume of solution must be sufficient to cover the gel completely and to allow the gel to move freely during all of the steps.



When using the Coomassie R-250 microwave staining protocol, warm the Coomassie R-250 stain and destaining solution to 50°C without boiling. **It is important NOT to boil the solutions.**

Since microwave ovens differ significantly, we recommend testing various times (10 second intervals) and power settings of your microwave oven to achieve a temperature of 50°C in the volume of solution required for your particular staining container. Perform these steps without the gel.

Once you have optimized the time and settings for your microwave, use these settings for staining.

Coomassie Staining, Continued



Note

The staining solutions for the Coomassie R-250 staining protocol and the microwave staining protocol are different. Be sure to use the correct stain for the correct protocol.

Coomassie R-250 staining protocol

For all staining and destaining steps described below, be sure to use sufficient reagents to completely cover the gel using a suitable container such that the gel moves freely during the staining and destaining steps.

1. After electrophoresis, remove the gel from the cassette (page 17) and place the gel in a clean incubation tray.
 2. Prepare Coomassie stain (0.03% Coomassie R-250 in 30% methanol and 10% acetic acid). See note on previous page.
 3. Stain the gel in the prepared stain for 1.5 hours at room temperature with gentle shaking.
 4. Destain the gel in destaining solution (see previous page) at room temperature with gentle shaking with intermittent changes of solution until the bands are visible or overnight for maximum sensitivity and clear background.
-

Coomassie R-250 staining microwave protocol

For all staining and destaining steps described below, be sure to use sufficient reagents to completely cover the gel in a **microwave safe** container such that the gel moves freely during the staining and destaining steps.

1. After electrophoresis, remove the gel from the cassette (page 17) and place the gel in a clean **microwave safe** container.
2. Prepare Coomassie stain (0.015% Coomassie R-250 in 30% ethanol and 10% acetic acid). See note on previous page.
3. Add enough stain to completely cover the gel in the **microwave safe** container.
4. Warm the gel and solution to about 50°C in a microwave oven (see page 19). **Note:** Do NOT boil the solution.
5. Incubate the gel in the warmed staining solution for 30 minutes on an orbital shaker at room temperature.
6. Discard the stain, rinse the gel briefly with water and discard the water.
7. Add enough destaining solution (see previous page) to cover the gel during incubation
8. Place two pieces of positively charged nylon membrane on top of the destaining solution to speed up the destaining process.
9. Warm the destaining solution, gel and nylon membrane to 50°C in a microwave oven (see previous page). **Note:** Do NOT boil the solution.
10. Incubate the gel in the warm destaining solution on an orbital shaker at room temperature until the desired background is achieved.

Note: To obtain a clear background, perform destaining overnight. Results obtained with Coomassie stain are shown on page 24.

Blotting of E-PAGE™ Gels

Semi-Dry Blotting of E-PAGE™ Gels

Introduction

A semi-dry blotting procedure for blotting E-PAGE™ 48 Gels is described in this section. You will need a semi-dry apparatus that can accommodate the dimensions of an E-PAGE™ Gel.

Additional blotting protocols

Guidelines for semi-wet and other blotting methods can be found in the E-PAGE™ Technical Guide available at www.invitrogen.com or contact Technical Support (page 32).

Materials needed

For ordering information, see page 31.

- Semi-dry blotter
- Methanol
- NuPAGE® Transfer Buffer (20X)
- NuPAGE® Antioxidant
- Blotting membranes: Invitrolon™/Filter Paper Sandwich or Nitrocellulose/Filter Paper Sandwiches
- E-PAGE™ Blotting Pad (supplied with E-PAGE™ Gels or available separately)
- 4 pieces of 2.5 mm Blotting Filter Paper
- Blotting Roller
- Incubation Tray

IMPORTANT! To ensure a good transfer, it is necessary to use the E-PAGE™ Blotting Pad and 2.5 mm Blotting Filter Paper.

Preparing 2X transfer buffer

We recommend using 2X NuPAGE® Transfer Buffer with 15% methanol and NuPAGE® Antioxidant for optimal transfer of most proteins from E-PAGE™ Gels. When the complete transfer of higher molecular weight proteins (>150 kDa) is desired, reduce methanol to 10%.

For one gel, prepare 500 mL of 2X NuPAGE® Transfer Buffer:

Buffer Component	2X NuPAGE® Transfer Buffer with 15% Methanol	2X NuPAGE® Transfer Buffer with 10% Methanol
NuPAGE® Transfer Buffer (20X)	50 mL	50 mL
NuPAGE® Antioxidant	0.5 mL	0.5 mL
Methanol	75 mL	50 mL
Deionized Water	to 500 mL	to 500 mL

Semi-Dry Blotting of E-PAGE™ Gels, Continued

Equilibrating the gel

Equilibration of the gel in transfer buffer results in the removal of salts that may increase conductivity and heat during transfer. Perform equilibration for the recommended time, as longer equilibration results in protein diffusion.

1. After electrophoresis, remove the gel from the cassette as described on page 17.
 2. Using the Butterfly Opener or a gel knife, trim off the top and bottom electrode areas of the gel.
 3. Equilibrate the E-PAGE™ Gel in 200 mL 2X NuPAGE® Transfer Buffer (see previous page) for 30 minutes with shaking.
-

Preparing blotting membrane

Nitrocellulose

1. Use pre-cut Nitrocellulose/Filter Paper Sandwich or cut nitrocellulose membrane to the appropriate size (8.6 cm x 13.5 cm).
2. Soak the membrane in a 2X NuPAGE® Transfer Buffer (see previous page) for several minutes in the Incubation Tray.

PVDF

1. Use pre-cut Invitrolon™/Filter Paper Sandwich or cut PVDF membrane to the appropriate size (8.6 cm x 13.5 cm).
 2. Pre-wet the membrane for 30 seconds in methanol, ethanol, or isopropanol. Briefly rinse the membrane in deionized water.
 3. Soak the membrane in a 2X NuPAGE® Transfer Buffer (see previous page) for several minutes in the Incubation Tray.
-

Semi-Dry blotting protocol

1. In a clean container or Incubation Tray, soak 4 pieces of 2.5 mm Blotting Filter Paper (8.6 cm x 13.5 cm) in 2X NuPAGE® Transfer Buffer (see page 21). Remove any air bubbles trapped between filter paper sheets using the Blotting Roller while the paper is still submerged in buffer.
 2. In a clean container or Incubation Tray, soak the E-PAGE™ Blotting Pad in 2X NuPAGE® Transfer Buffer (see page 21). Press the pad to ensure the elimination of any visible air bubbles. Inspect both sides of the pad for air bubbles before use.
Note: The Blotting Pad has no specific orientation (either side may be facing toward the gel).
 3. Place 2 pieces of pre-soaked 2.5 mm Blotting Filter Paper from Step 1 on the anode plate of a semi-dry blotting apparatus. Ensure that all filter paper sheets are aligned properly and remove any air bubbles with the Blotting Roller.
 4. Place the pre-soaked blotting membrane on top of the filter paper stack and remove any air bubbles with the Blotting Roller.
 5. Remove the gel from the transfer buffer. Gently rub a gloved finger over the well side of the gel to remove small gel pieces from the gel surface. Re-submerge the gel in transfer buffer to remove any gel pieces from the gel, as they can cause air bubbles and field distortion during transfer.
-

Semi-Dry Blotting of E-PAGE™ Gels, Continued

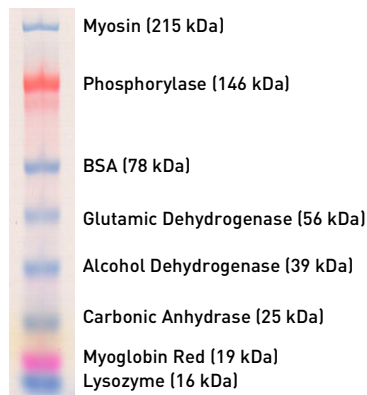
Semi-Dry blotting protocol, continued

6. Place the flat side of the gel on top of the blotting membrane such that the well side of the gel is facing up and remove any air bubbles with the Blotting Roller. Fill the wells of the gel with 2X NuPAGE® Transfer Buffer (see page 21).
 7. Place the pre-soaked E-PAGE™ Blotting Pad on the gel and gently roll out air bubbles using the Blotting Roller.
 8. Place 2 pieces of 2.5 mm Blotting Filter Paper from Step 1 on top of the Blotting Pad. Ensure that all filter paper sheets are aligned properly and flush with the gel/membrane sandwich. Remove any air bubbles with the Blotting Roller.
 9. Place the cathode plate on the stack without disturbing the blot sandwich. Follow the manufacturer's instructions to further assemble the semi-dry blotting apparatus.
 10. Transfer at 25 V for 1 h (~19 V/cm). You may need to optimize the transfer conditions for your specific proteins or semi-dry blot apparatus.
-

Expected Results

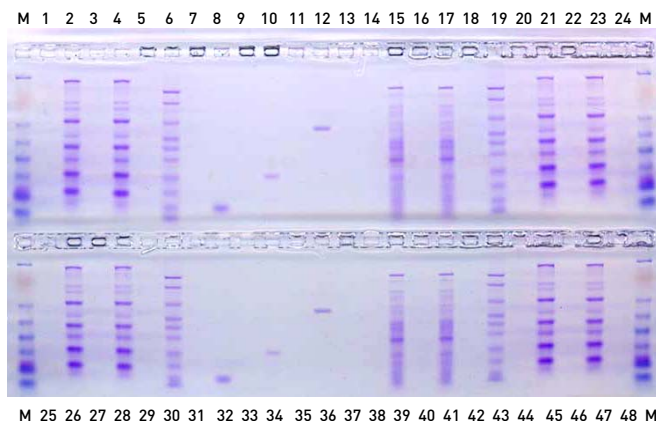
Molecular weight calibration

Apparent molecular weight values for SeeBlue® Plus2 Pre-stained Standard on an E-PAGE™ 48 8% Gel are shown below:



Results with Coomassie R-250 staining

Samples were run on an E-PAGE™ 48 8% Gel and stained with Coomassie R-250 using the microwave staining protocol as described in this manual (page 20). Results after staining are shown below.



The gel contains the following samples (rows not indicated are blank):

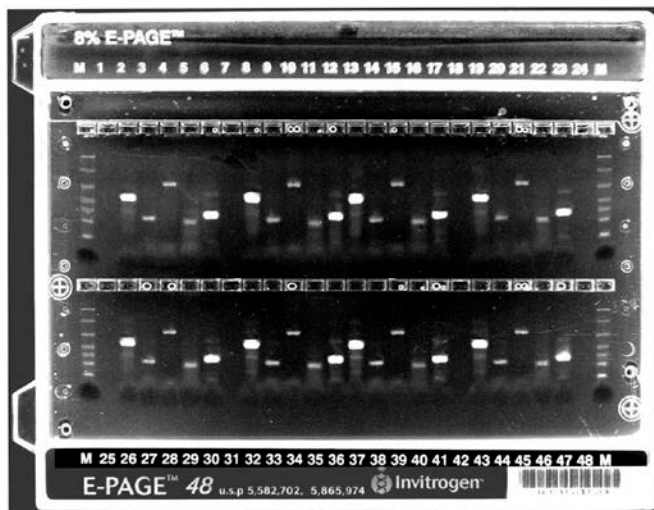
Lane	Sample
M	SeeBlue® Plus2 Pre-stained Standard (5 µL)
2, 4, 21, 23, 26, 28, 45, 47	MagicMark™ XP Western Protein Standard (10 µL)
6, 19, 30, 45	Benchmark™ His-tagged Protein Standard (10 µL, diluted 1:5)
8, 32	Lysozyme (200 ng)
10, 34	Carbonic Anhydrase (200 ng)
12, 36	BSA (100 ng)
15, 17, 39, 41	<i>E. coli</i> lysate (5 µL)

Expected Results, Continued

Lumio™ Green detection

Results obtained with an E-PAGE™ 48 8% Gel using the Lumio™ Green Detection Kit are shown below.

Various concentrations of a Lumio™ fusion protein were labeled with Lumio™ Green Detection Kit and run on an E-PAGE™ 48 8% Gel as described in this manual.



The gel contains the following samples (lanes not indicated are blank):

Lane	Sample
M	BenchMark™ Fluorescent Marker (5 µL)
2, 8, 13, 19, 26, 32, 37, 43	Human kinase Lumio™ fusion protein (10 µL)
3, 9, 14, 20, 27, 33, 38, 44	<i>E. coli</i> CAT Lumio™ fusion protein (10 µL)
4, 10, 15, 21, 28, 34, 39, 45	<i>E. coli</i> GUS Lumio™ fusion protein (10 µL)
5, 11, 16, 22, 29, 35, 40, 46	<i>E. coli</i> calmodulin Lumio™ fusion protein (10 µL)
6, 12, 17, 23, 30, 36, 41, 47	<i>E. coli</i> kinase D Lumio™ fusion protein (10 µL)

Using E-Editor™ 2.02 Software

Introduction

The E-Editor™ 2.02 software for Windows® allows you to reconfigure digital images of E-PAGE™ Gels for analysis and documentation. E-Editor™ 2.02 software reconfigures the wells of the E-PAGE™ Gels into a side-by-side format for easy comparison and analysis.

You can reconfigure gels that were scanned in the original gel cassette, or gels that were removed from the cassette. You can also group the images of multiple gels loaded from a 384-well microtiter plate into a single image with a layout corresponding to that of the original plate.

Capture an image of the gel as described below and then use the E-Editor™ 2.02 software to:

- Align and arrange the lanes in the image
 - Save the reconfigured image for further analysis
 - Copy and paste selected lanes or the entire reconfigured image into other applications for printing, saving, emailing, and/or publishing
-

Imaging the gel

Use an appropriate gel documentation system to capture a digital image of the gel. When imaging, the gel should be properly aligned (i.e., not at an angle) and gel features should be clear and distinct. Proceed to **Downloading Software**.

Downloading software

E-Editor™ 2.02 software can be downloaded FREE from our website. Go to www.lifetechnologies.com/epage and follow the instructions to download the software and user manual.

Troubleshooting

Troubleshooting

The table below provides some solutions to possible problems you might encounter during the electrophoresis of E-PAGE™ Gels.

Observation	Cause	Solution
No current	Daughter E-Base™ Device used without a Mother E-Base™ Device	Do not use the Daughter E-Base™ Device without a Mother E-Base™ Device. The Daughter E-Base™ Device does not have an electrical plug to connect to an electrical outlet.
No electric contact (no red light when cassette is inserted) or run does not start (no green light)	Copper contacts in the base are damaged due to improper use	Make sure that the copper contacts in the base are intact.
	Expired or defective gel cassette used	Use properly stored gels before the specified expiration date.
	E-PAGE™ cassette is not correctly inserted into base	Remove cassette and reinsert. When the cassette is correctly inserted and power is on, a fan in the base begins to run and a steady red light illuminates on the base (page 13).
Sample leaking from the wells	Sample is overloaded or wells are damaged	Be sure to load the recommended volume of sample per well (page 12). Remove the comb carefully without damaging the wells.
Poor resolution or smearing of bands	Sample is overloaded	Do not load more than 20 µg of protein sample per well.
	Very low volumes of sample were loaded	Do not load less than 5 µL of sample. Always load 10–20 µL deionized water in all wells prior to sample loading. For proper band separation, we recommend keeping sample volumes uniform and loading deionized water into empty wells.
	Incorrect loading buffer used	Make sure that protein sample is in one of Use the recommended loading buffers as described on page 7.
	Electrophoresis was not started immediately after sample loading	For best results, the gel should be run within 15 minutes of sample loading.

Continued on next page

Troubleshooting, Continued

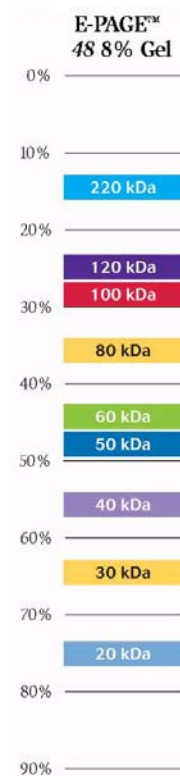
Observation	Cause	Solution
Poor resolution or smearing of bands	High salt or detergent concentration in samples	Be sure the final concentration of salt or detergent in the sample is as described on page 6. You may need to manually increase the run time for high salt or detergent samples to obtain optimal results.
	Expired gel used	Use properly stored gels before the specified expiration date.
Over-run the gel or need more time to run gel	Accidentally selected an incorrect program	Select program EP for E-PAGE™ Gels. If you accidentally selected an incorrect program and are at the beginning of the run, stop the run and select the desired program. If you are well into the run, check the gel to see where the loading dye is running. Estimate the amount of time remaining and then manually stop the run.
Protein bands distorted on membrane after semi-dry blotting	Non-uniform electric field created around wells	Be sure that the E-PAGE™ Blotting Pad is used correctly.
	Incorrect gel orientation	Be sure that the well side of the gel is not facing the membrane.
Weak transfer of high molecular weight samples during semi-dry blotting	Not enough SDS in sample	Reduce methanol concentration in transfer buffer from 15% to 10% if transferring E-PAGE™ 48.
Weak transfer of low molecular weight samples	Use of large pore membranes allow small proteins to “blow through”	Use 0.2 µm nitrocellulose membrane for optimal capture of small proteins.
Uneven transfer of proteins and edge lanes during semi-dry blotting of E-PAGE™ 48 Gel	No methanol in transfer buffer	Use 10–15% methanol in the transfer buffer.
Weak transfer of proteins during semi-wet blotting	No methanol in transfer buffer	Use 10% methanol in the transfer buffer.

Appendix

Product Specifications

E-PAGE™ gel separation range

The migration and resolution range of proteins run on E-PAGE™ 48 8% Gel.



E-PAGE™ 48 Gel specifications

Each E-PAGE™ 48 gel contains 48 sample wells and 4 marker wells (M).

Cassette Size: 13.5 cm (l) × 10.8 cm (w) × 0.67 cm (thick)

Gel Thickness: 3.7 mm

Gel Volume: 50 mL

Gel Formulation: Proprietary, operating at a neutral pH

Well Depth: 3 mm

Well Volume: 15 µL

Well Opening: 3.6 mm (l) × 2.2 mm (w)

Running Distance:
(one well to the next) 3.2 cm

Space between Wells: 4.5 mm

Note: E-PAGE™ 48 8% Gels have a unique separation profile, which gives protein resolution similar to that of a 4–12% Tris-Glycine gel.

Accessory Products

E-PAGE™ Gels

The following E-PAGE™ Gel kits are available from Life Technologies.

Product	Quantity	Catalog no.
E-PAGE™ 48 8% Gels	8-pack	EP048-08
E-PAGE™ 96 6% Gels	8-pack	EP096-06

Electrophoresis Bases

The following electrophoresis bases are available from Life Technologies for running E-PAGE™ gels.

- The Mother E-Base™ Device (Cat. no. EB-M03) is used for electrophoresis of one E-PAGE™ Gel.
 - The Daughter E-Base™ Device (Cat. no. EB-D03) attaches to the Mother E-Base™ Device and together are used for the independent electrophoresis of two or more E-PAGE™ Gels.
-

E-Holder™

The E-Holder™ Platform is used to hold an E-PAGE™ Gel in place while loading. Ordering information can be found on the following page.

E-Editor™ 2.02 Software

The E-Editor™ 2.02 software is available FREE with the purchase of any E-PAGE™ Gels or related equipment. The software may be downloaded from www.lifetechnologies.com/epage.

iBlot® Gel Transfer Device

The iBlot® Gel Transfer Device is available from Life Technologies for transfer of proteins from E-PAGE™ gels to nitrocellulose or PVDF membranes.

Product	Quantity	Catalog no.
iBlot® Gel Transfer Device	1 unit	IB1001, IB1001UK, IB1001EU
iBlot® Gel Transfer Stack, Nitrocellulose, Regular	10-pack	IB3010-01
iBlot® Gel Transfer Stack, PVDF, Regular	10-pack	IB4010-01

Continued on next page

Accessory Products, Continued

Additional products

The following products for use with E-PAGE™ gels are available separately from Life Technologies:

Product	Quantity	Catalog no.
SeeBlue® Plus2 Pre-Stained Standard	500 µL	LC5925
E-PAGE™ SeeBlue® Pre-stained Protein Standard	500 µL	LC5700
MagicMark™ XP Western Standard	250 µL	LC5602
E-PAGE™ MagicMark™ Unstained Protein Standard	250 µL	LC5701
BenchMark™ Fluorescent Protein Standard	125 µL	LC5928
Lumio™ Green Detection Kit	1 kit	LC6090
SYPRO® Ruby Protein Gel Stain	1 L	S-12000
SimplyBlue™ SafeStain	1 L	LC6060
SilverQuest™ Silver Staining Kit	1 kit	LC6070
SilverXpress® Silver Staining Kit	1 kit	LC6100
InVision™ His-tag In-gel Stain	500 mL	LC6030
NuPAGE® Transfer Buffer (20X)	125 mL	NP0006
NuPAGE® Antioxidant	15 mL	NP0005
NuPAGE® Sample Reducing Agent (10X)	10 mL	NP0009
Nitrocellulose/Filter Paper Sandwich 0.45 µm	16/pk	LC2006
Nitrocellulose/Filter Paper Sandwich 0.2 µm	16/pk	LC2009
Invitrolon™ PVDF/Filter Paper Sandwich 0.45 µm	16/pk	LC2007
Blotting Filter Paper (2.5 mm)	50/pk	LC2008
E-PAGE™ Blotting Pad	4/pk	LC2101
Blotting Roller	1	LC2100
Incubation Tray	8/pk	LC2102
Gel Knife	1	EI9010
Large Gel Drying Kit	1 kit	NI2207
Gel-Dry™ Drying Solution (1X)	500 mL	LC4025
WesternBreeze® Chromogenic Kit		
Anti-Mouse	1 kit	WB7103
Anti-Rabbit	1 kit	WB7105
Anti-Goat	1 kit	WB7107
WesternBreeze® Chemiluminescent Kit		
Anti-Mouse	1 kit	WB7104
Anti-Rabbit	1 kit	WB7106
Anti-Goat	1 kit	WB7108
Pro-Q® Diamond Phosphoprotein Gel Stain	1 L	P-33300
Pro-Q® Diamond Phosphoprotein Gel Destain Solution	1 L	P-33310
Pro-Q® Sapphire 532 Oligohistidine Gel Stain	500 mL	P-33354

Technical Support

Obtaining support For the latest services and support information for all locations, go to www.lifetechnologies.com

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
 - Search through frequently asked questions (FAQs)
 - Submit a question directly to Technical Support (techsupport@lifetech.com)
 - Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
 - Obtain information about customer training
 - Download software updates and patches
-

Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/support.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Limited warranty

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies General Terms and Conditions of Sale found on the Life Technologies web site at <http://www.lifetechnologies.com/termsandconditions>. If you have any questions, please contact Life Technologies.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

29 October 2014

