

E-PAGE™ Technical Guide

General information and protocols for using E-PAGE™ Gels

Catalog Number EP096-06

Revision B.0

Publication Number MAN0000374

Publication Part Number 25-0644



Manufacturer: Life Technologies Corporation | 29851 Willow Creek | Eugene, Oregon 97402 USA

Information in this document is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

© 2021 Thermo Fisher Scientific Inc. All rights reserved.

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

www.lifetechnologies.com

06 October 2021

ThermoFisher
S C I E N T I F I C

Contents

Product Contents	iv
Introduction	1
E-PAGE™ Protein Electrophoresis System	1
E-PAGE™ System Components	3
Methods	8
Electrophoresis of E-PAGE™ Gels.....	8
Guidelines for Sample Preparation.....	8
Electrophoresis of E-PAGE™ Gels	13
Running E-PAGE™ Gels.....	17
Guidelines for Loading E-PAGE™ Gels by Automated Liquid Handling	18
Using E-Holder™ Platform	20
Opening the E-PAGE™ Cassette	24
Blotting E-PAGE™ Gels.....	25
Blotting E-PAGE™ Gels with the iBlot® Device	25
Semi-Dry Blotting of E-PAGE™ Gels.....	31
Semi-Wet Blotting of E-PAGE™ Gels	35
Visualizing and Staining of E-PAGE™ Gels	38
Visualizing Lumio™ Fusion Proteins	39
SYPRO® Ruby Staining Protocol	40
Coomassie Staining	41
SimplyBlue™ SafeStain	43
Silver Staining Methods.....	45
InVision™ His-tag Protein Gel Stain.....	48
Drying of E-PAGE™ Gels	49
Gel Drying	49
Analysis and Troubleshooting	51
Molecular Weight Calibration	51
Expected Results.....	53
Using E-Editor™ 2.02 Software.....	60
Troubleshooting.....	61
Appendix.....	63
Additional Staining Protocols.....	63
Product Specifications.....	67
Accessory Products	70
Technical Support.....	72

Product Contents

Purpose of the Guide

The E-PAGE™ Technical Guide contains information about E-PAGE™ Gels and is intended to supplement the manuals supplied with E-PAGE™ Gels. Details for sample preparation, electrophoresis conditions, blotting, and staining are included in this guide.

Types of Kits

E-PAGE™ gels are designed for use with the E-Base™ Electrophoresis Device, and are available in 48-well or 96-well formats. See pages 3–4 for further specifications. See page 71 for ordering information.

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

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 96 Gels, E-PAGE™ Blotting Pad and Mother E-Base™ 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™ 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 self-contained E-PAGE™ 48 or 96 Pre-cast Gels consist of a buffered gel matrix and electrodes packaged inside a disposable, UV-transparent cassette. The gels can be loaded by multichannel pipetter or automated loading system.

Electrophoresis of the samples is performed with the E-Base™ Electrophoresis Device. The E-Base™ Electrophoresis Device consists of a base for electrophoresis, and an integral power supply. After electrophoresis is completed, the cassette can be opened with the Invitrogen™ Gel Knife (Cat. No. EI9010, sold separately) to perform downstream applications such as western blotting or staining.

The E-Editor™ 2.02 software allows digital images of the E-PAGE™ Gels to be reconfigured into a side-by-side format for easy comparison and analysis.

Applications

The E-PAGE™ Protein Electrophoresis System is ideal for screening protein samples using the following applications:

- In-gel staining with Lumio™ Green Reagent
 - Western blotting
 - Staining (Coomassie, silver, or fluorescent stains)
 - Functional assays
-

Advantages

Using E-PAGE™ Gels for protein analysis offers the following advantages:

- Produces fast, safe, and consistent electrophoresis
 - Eliminates the need to prepare gels and buffers
 - Compatible with most commercially available automated liquid handling systems for high-throughput electrophoresis
 - Available in two well formats to suit your applications
 - Easy to open gel cassettes facilitate downstream staining and blotting applications
 - Ability to track samples during electrophoresis of multiple gels using an individual barcode on each gel cassette
-

E-PAGE™ Protein Electrophoresis System, Continued

Choosing a Gel for your Application

To obtain the best results for your application it is important to choose the correct E-PAGE™ Gel. The table below describes the major features of each format. For specifications and separation range of each gel type, see page 67–68.

E-PAGE™ Gel	No. Rows	No. Wells	Well Volume	Run Length	Run Time
48 Gel	2*	48 sample 4 marker	15 µL 15 µL	3.2 cm	25 min.
96 Gel	8*	96 sample 8 marker	20 µL 20 µL	1.6 cm	14 min.

*Wells compatible for loading with a multichannel pipetter or with an automated liquid handling system.

Western Blotting of E-PAGE™ Gels

E-PAGE™ Gels are suitable for western blotting. The iBlot® Gel Transfer Device is recommended method for blotting E-PAGE™ Gels (page 25). Semi-dry blotting (page 25), and semi-wet blotting may also be performed (page 35). Tank blotting is not recommended, but can be performed if even pressure is maintained between the gel and the membrane, and cooling is provided.

Staining E-PAGE™ Gels

E-PAGE™ Gels are compatible with many standard Coomassie, silver, or fluorescent staining protocols. E-PAGE™ Gels are thicker than most SDS-PAGE mini-gels, so additional time may be required for staining and destaining steps. We recommend the following stains for E-PAGE™ Gels. Detailed staining protocols are described on pages 38–48.

Total Protein Stains:

- SYPRO® Ruby Protein Gel Stain (page 38)
- Coomassie Stain (page 41)
- SimplyBlue™ SafeStain (page 43)
- SilverQuest™ Silver Staining Kit (page 46)
- SilverXpress® Silver Staining Kit (page 47)

Specific Protein Stains:

- Lumio™ Green Detection Reagent for detecting Lumio™ fusion proteins (page 38)
- InVision™ His-tag In-Gel Stain for detecting 6X His-tagged proteins (blotting recommended first, page 48)

E-PAGE™ System Components

Introduction

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

- E-PAGE™ 48 or 96 Pre-cast Gels
- E-Base™ Electrophoresis Device
- E-Holder™ Platform
- E-PAGE™ Loading Buffer 1 (4X)
- E-PAGE™ Blotting Pad
- E-Editor™ 2.02 Software

Details on each system component are described in this section.

E-PAGE™ 48 Gels

- 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 17 for automation specifications).
- After electrophoresis, the E-PAGE™ 48 cassette is easily opened with the Invitrogen™ Gel Knife (Cat. No. EI9010, sold separately) 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 13).

See page 67 for E-PAGE™ 48 8% Gel specifications.

Diagram of an E-PAGE™ 48 Cassette

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



E-PAGE™ System Components, Continued

E-PAGE™ 96 Gels

- E-PAGE™ 96 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™ 96 Gel contains 96 sample lanes and 8 marker lanes in a patented staggered well format that provides a 1.6-cm run length. The wells of the E-PAGE™ 96 gels are compatible for loading with a multichannel pipetter or with an automated liquid handling system (See page 17 for automation specifications).
- After electrophoresis, the E-PAGE™ cassette is easily opened with the Invitrogen™ Gel Knife (Cat. No. EI9010, sold separately) to remove the gel from the cassette for staining or blotting applications.
- Each E-PAGE™ 96 cassette is labeled with an individual barcode to facilitate identification of the gel using commercial barcode readers (page 13).

See page 67 for E-PAGE™ 96 6% Gel specifications

Diagram of an E-PAGE™ 96 Cassette

A diagram of the E-PAGE™ 96 Gel cassette is shown below.



E-PAGE™ System Components, Continued

E-Base™ Electrophoresis Device

E-PAGE™ 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™** (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™ Gel. The Mother E-Base™ has been tested for electrophoresis with up to three Daughter E-Bases™ connected at one time.
 - **The Daughter E-Base™** (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™ Gels. The Daughter E-Base™ does not have an electrical plug and cannot be used without a Mother E-Base™.
-

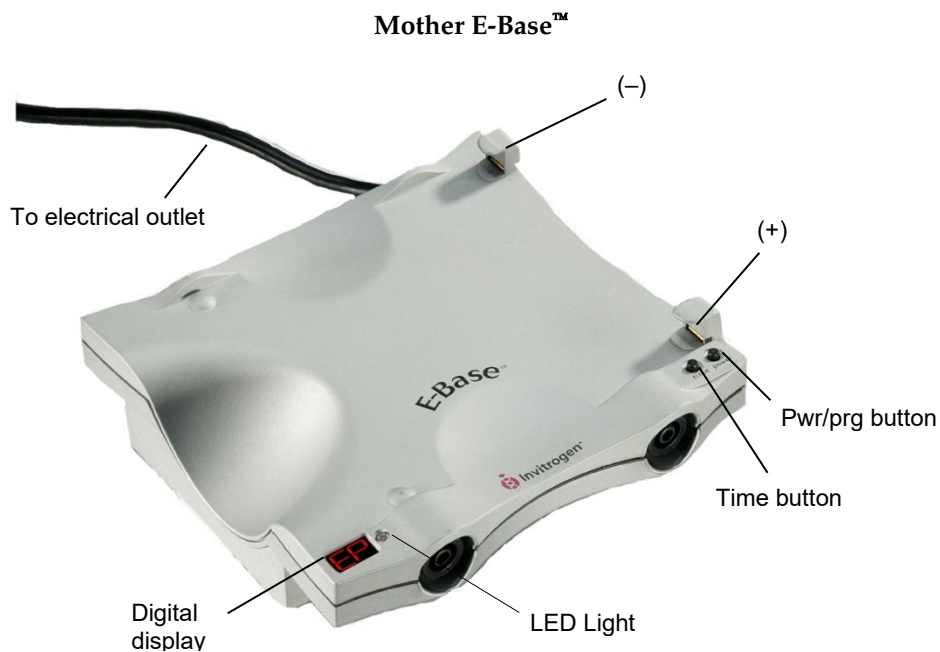


E-PAGE™ Gels are **not compatible** with the E-Gel® 96 bases (Cat nos. G7100-01/G7200-01) previously available from Life Technologies. The older E-Gel® 96 bases do not have the 'E-Base™' inscription on the platform.

Mother E-Base™

Each Mother E-Base™ has a pwr/prg (power/program) button (right side) and a time button (left side) on the lower-right side of the base. The lower-left side of each Mother E-Base™ contains a light LED and a digital display. The gel cassette is inserted into the two electrode connections. The Mother E-Base™ is connected to an electrical outlet with the electrical plug.

The E-Base™ EP program is set to 14 minutes default time. See page 15 to adjust time as recommended.



E-PAGE™ System Components, Continued

Daughter E-Base™ The Daughter E-Base™ is similar to the Mother E-Base™ except the Daughter E-Base™ does not have an electrical cord and cannot be connected to an electrical outlet.

The Daughter E-Base™ is connected to a Mother E-Base™ or to another Daughter E-Base™ (already connected to a Mother E-Base™). Once connected to a Mother E-Base™, each Daughter E-Base™ is designed to function independently of the Mother E-Base™ or other Daughter E-Bases™

Mother E-Base™/Daughter E-Base™



E-PAGE™ System Components, Continued

E-PAGE™ Loading Buffer

E-PAGE™ Gels are supplied with E-PAGE™ Loading Buffer 1 (4X), which is optimized for E-PAGE™ Gels, and is recommended for routine SDS-PAGE and staining or blotting applications. **It is not recommended to use any other SDS-PAGE sample buffer.** See page 10 for more information about loading buffers.

E-Holder™ Platform

The E-Holder™ Platform is designed to hold E-PAGE™ Gels during loading. Use the E-Holder™ when you need to load multiple gels while other gels are running on the E-Base™.

Note: The E-Holder™ is not a power supply unit, cannot be connected to an electrical outlet, and cannot be used to run gels.

See page 17 for more information about the E-Holder™ Platform.

E-PAGE™ Blotting Pad

The E-PAGE™ Blotting Pad is supplied with the E-PAGE™ Gels and E-PAGE™ Starter Kits. 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

Electrophoresis of E-PAGE™ Gels

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.



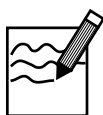
We recommend that you read this section before preparing your samples. Use the sample preparation method and loading buffer appropriate for your detection method:

Application	Method	Loading Buffer
Routine staining and western blotting	1 (page 11)	Loading Buffer 1
Lumio Green Detection	2 (page 12)	Lumio™ Gel Sample Buffer

Materials Needed

The following items are needed for sample preparation. See page 71 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 10)
 - (Optional) Lumio™ Green Detection Kit for detection of Lumio™ fusion proteins (page 12)
-



Note

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

To obtain the best results, we recommend performing 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

For best results, 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 Volume
E-PAGE™ 96 6% Gel	15 μL
E-PAGE™ 48 8% Gel	10 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™ 48	Final Concentration for E-PAGE™ 96
Triton® X-100	<0.3%	<0.5%
Tween® 20	<0.3%	<0.5%
CHAPS	<0.3%	NT
NP-40	<0.3%	NT
RIPA	<0.25X	NT
SDS	<2% (already in loading buffer)	<4%
Tris	<300 mM	<200 mM
NaCl	<300 mM	<250 mM
Ammonium sulfate	<100 mM	NT
Sodium acetate	<200 mM	NT
EDTA	<20 mM	NT
MES	Not recommended	NT
DTT, Glycine, Urea, Imidazole	No effect seen up to 500 mM	NT

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 11)**
 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 12)**
 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 51 for apparent molecular weight calibration on E-PAGE™ Gels. See page 71 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
E-PAGE™ 96 6% Gel	E-PAGE™ SeeBlue® Pre-stained Protein Standard	10 µL	Electrophoresis
	E-PAGE™ MagicMark™ Unstained Protein Standard	10 µL	Electrophoresis, followed by staining
	E-PAGE™ MagicMark™ Unstained 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 16.
-

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.

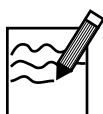
For details on the Lumio™ Green Detection Kit, refer to the manual available at www.lifetechnologies.com or contact Technical Support (page 72).

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 13.



Note

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

Electrophoresis of E-PAGE™ 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 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 17).

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 or 96 Gels. This ensures proper loading of samples into the wells. See page 17 for automation guidelines.



- Dispose of gels as hazardous waste.
 - Avoid touching the gel while the gel is running.
-

Using the Barcode

Each E-PAGE™ 48 and 96 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™ Gels, Continued



Important

- 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, we do not recommend running 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.
 - E-PAGE™ 48 and 96 Gels are not compatible with the E-Gel® 96 Mother and Daughter Bases (Cat nos. G7100-01/G7200-01) available previously from Life Technologies. The older E-Gel® 96 bases do not have the “E-Base™” inscription on the platform.
-

Method of Loading Samples

We recommend the following methods of sample loading based on the gel type. Use short, rigid tips for loading E-PAGE™ Gels.

Gel Type	Method of Loading
E-PAGE™ 48	Manual pipetter, multichannel pipetter (load samples into alternate wells of the gel followed by a second round of loading into the remaining wells), or automated liquid handling devices (8- or 12-tip). See page 17 for automated liquid handling system guidelines.
E-PAGE™ 96	Manual pipetter, multichannel pipetter, or automated liquid handling devices (8-, 12-, or 96-tip). See page 17 for automated liquid handling system guidelines.

Electrophoresis of E-PAGE™ Gels, Continued

Selecting a Program on the E-Base™ Device

The recommended program for E-PAGE™ Gels is EP. **Select program EP 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 the 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 (power/program) button to select the **program EP**. The default run time for program EP is 14 minutes.
4. To manually change the default run time, press and release the time button located on the lower-right corner of the base to view the time setting.
5. Press and hold the time button to increase the time to the desired run time. 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.

Gel Type	Time in Minutes
E-PAGE™ 48 8%	25 minutes
E-PAGE™ 96 6%	14 minutes (default run time for program EP)

Do not run an E-PAGE™ 96 Gel for more than 25 minutes, or an E-PAGE™ 48 Gel for more than 30 minutes.



Note

- When performing electrophoresis of Lumio™ fusion proteins, we recommend extending 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 21.
-

Electrophoresis of E-PAGE™ Gels, Continued

Loading Gels

Each E-PAGE™ Gel is individually wrapped and ready for use.

Refer to pages 17–19 for automated liquid handling system guidelines for loading E-PAGE™ Gels. To load your samples on the gels using the E-Holder™ Platform, refer to page 17 for detailed instructions.

To load E-PAGE™ Gels on the E-Base™ Device:

1. Open the package and remove the gel.
2. Remove the plastic comb from the gel.
3. Slide the gel into the two electrode connections on the Mother E-Base™ Device 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.

4. Load deionized water to each well of the E-PAGE™ Gel prior to loading your samples or protein molecular weight standard. There should be a uniform volume of liquid in each well. Load samples into the gels using a multichannel pipetter or an automated liquid handling system as described below:

E-PAGE™ Gel Type	Deionized Water	Sample in Loading Buffer
E-PAGE™ 48 Gel	5–10 µL	10–5 µL
E-PAGE™ 96 Gel	5–15 µL	15–5 µL

5. Load the appropriate protein molecular weight standards in the marker wells of the gel. See page 10 for recommended molecular weight standards.
6. Proceed immediately to **Running E-PAGE™ Gels**, page 17.

Running E-PAGE™ Gels

Introduction

After loading your protein samples on the E-PAGE™ Gels, proceed immediately to electrophoresis using the E-Base™ Device. Refer to page 22 for the E-Base™ Quick Reference Guide.

Using E-Base™

Instructions for running E-PAGE™ Gels on the Mother E-Base™ Device or Daughter E-Base™ Device are provided below.

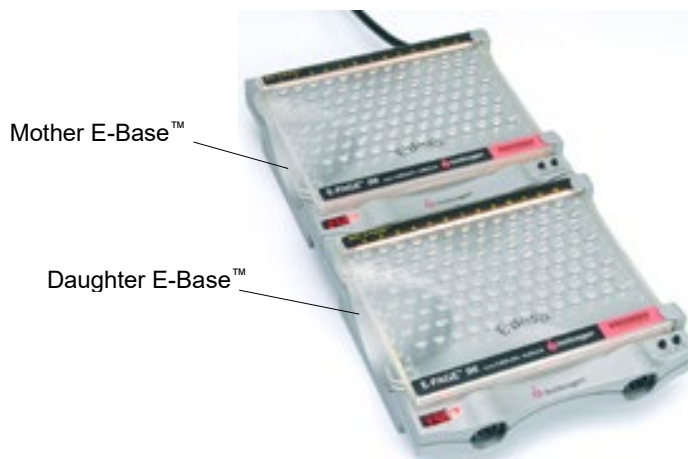
Note: It is not necessary to have a gel in the Mother E-Base™ Device if you are using a Daughter E-Base™ Device. However, the Mother E-Base™ Device must be plugged into an electrical outlet.

1. To begin electrophoresis, press and release the pwr/prg button located on the lower-right corner of the Mother E-Base™ Device.

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



2. If you are using a Daughter E-Base™ Device, press and release the pwr/prg button located on the lower-right corner of the Daughter E-Base™ Device.



To add to 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™ 96 Gel for more than 25 minutes or an E-PAGE™ 48 Gel for more than 30 minutes.

To interrupt or stop a run in progress, see page 21.

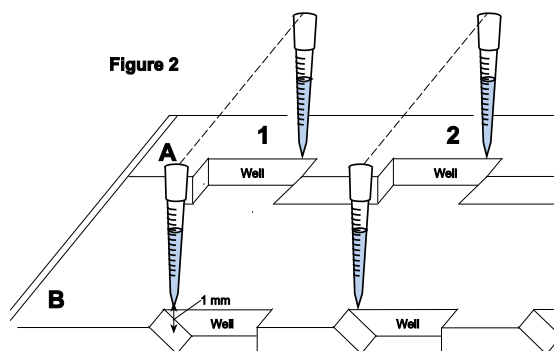
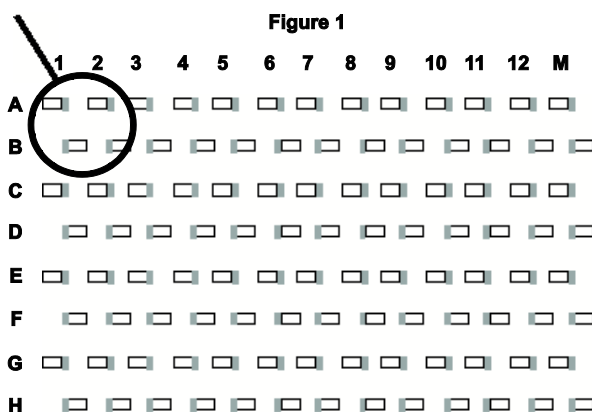
Guidelines for Loading E-PAGE™ Gels by Automated Liquid Handling

Introduction

E-PAGE™ 48 and 96 Gels are designed for automated loading using standard SBS (Society for Biomolecular Screening)-compatible liquid handling systems with an 8, 12, or 96-tip robotic loading apparatus. In general, follow the manufacturer instructions for your automated loading system. Specific guidelines for loading each gel format are described in the section below. To download programming scripts for your automated liquid handling system, go to www.lifetechnologies.com/epage.

Automated Loading of E-PAGE™ 96 Gels

The wells of the E-PAGE™ 96 Gel are staggered to provide maximum run length (see Figure 1, below). For proper loading of samples, it is important to program your automated liquid handling system to set the A1 tip of the 8, 12 or 96-tip robotic head over the E-PAGE™ 96 Gel cassette as described below. Set the position of the first tip approximately 1 mm above the slope of the A1 well (see Figure 2, below). This ensures that the remaining tips are aligned above the slopes of the remaining wells. Refer to the manufacturer's manual of your automated liquid handling system to program this setting. After programming the setting, load your samples. During loading, the samples drop onto the slopes of the wells and are drawn into the wells by capillary force.



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

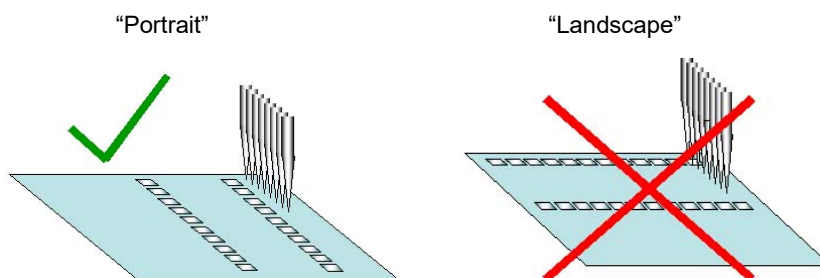
Automated Loading of E-PAGE™ 48 Gels

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:

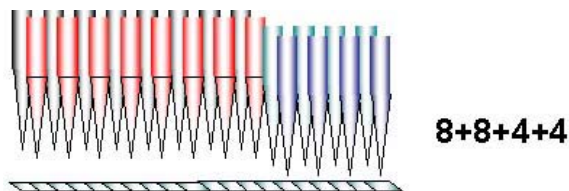


This orientation is available for some automation systems with the use of a 90° adapter. If your system does not have an adapter that allows the “Portrait” configuration, please contact Technical Support (page 72) to obtain a 90° adapter.

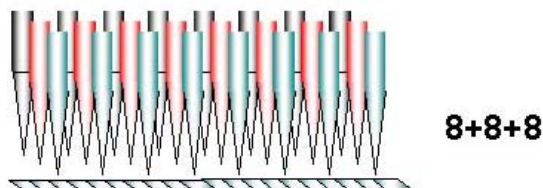
Loading Patterns for E-PAGE™ 48

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)



Using E-Holder™ Platform

Introduction

The E-Holder™ Platform is designed to hold E-PAGE™ Gels during loading. You can use the E-Holder™ Platform when you need to load multiple gels while other gels are running on the E-Base™

Note: The E-Holder™ Platform is not a power supply unit, cannot be connected to an electrical outlet, and cannot be used to run E-PAGE™ 48 or 96 Gels.

To obtain the best results, run E-PAGE™ Gels on the Mother E-Base™ Device or Daughter E-Base™ Device within 15 minutes after loading on E-Holder™ Platform.

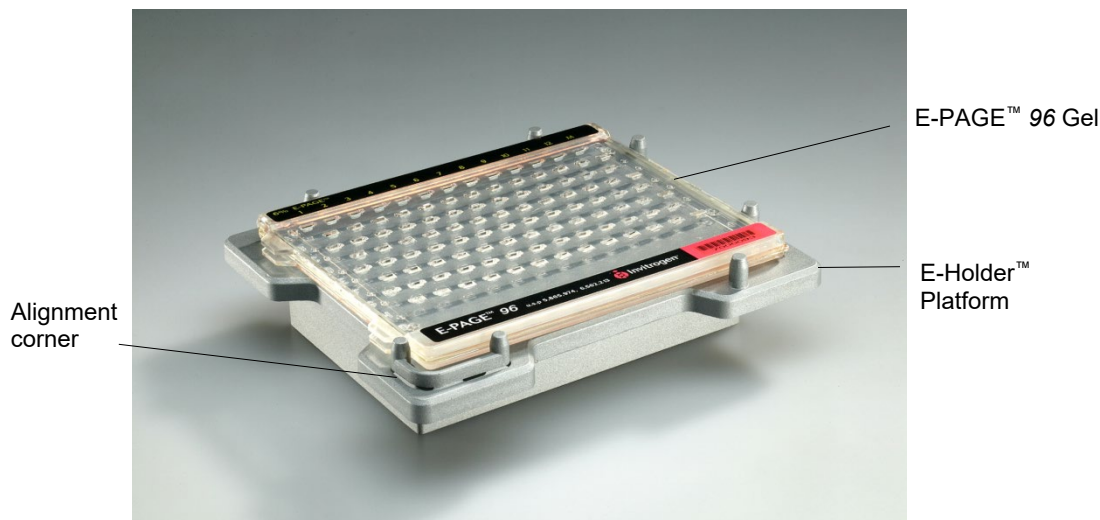


Important

The E-PAGE™ Gels will not fit on the E-Gel™ 96 Holder previously available from Life Technologies.

Using E-Holder™

1. Open the package and remove the E-PAGE™ Gel.
2. Remove comb from the E-PAGE™ cassette.
3. Place the E-PAGE™ cassette in the E-Holder™ Platform. Align the bottom-left end of the cassette in the lower-left alignment corner of the E-Holder™ Platform as shown in the figure below. Proceed to loading the gel or to Step 4 below for automated liquid handling system loading.



4. **Optional:** Set up your automated liquid handling system to load samples into the E-PAGE™ 48 or 96 Gel placed on an E-Holder™ Platform. Program your automated liquid handling system to load the samples approximately 5 minutes before the previous electrophoresis run is complete. This ensures that the loaded gel from the E-Holder™ Platform is placed onto an E-Base™ Device within the recommended time of 15 minutes.
-

Running E-PAGE™ Gels, Continued

Using E-Base™, Continued

3. The Mother E-Base™ Device or Daughter E-Base™ Device signals the end of the run with a flashing red light and rapid beeping for 2 minutes followed by a single beep every minute.

At the end of the run, the digital display shows the original time setting. The digital display does not any time change that was made during electrophoresis. The digital display also shows the elapsed time (up to 19 minutes with a negative sign) since the end of the run.

4. **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.
5. Remove the gel cassette from the Mother E-Base™ Device or Daughter E-Base™ Device.
6. For visualizing Lumio™ fusion proteins, see page 38. For opening the E-PAGE™ cassette, see page 24.

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

Interrupting an Electrophoresis Run

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 15 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™ Device 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.



Note

We recommend that you disconnect the Mother E-Base™ Device from the electrical outlet when not in use for a prolonged period of time.

Maintaining E-Base™

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.

E-Base™ Quick Reference Guide

Introduction

A quick reference guide for operating the Mother E-Base™ Device and Daughter E-Base™ Device is provided below. Operating modes and electrophoresis runs are described.

Mode	Action	Sound	Light	Digital Display
Base plugged in	Mother E-Base™ connected to an electrical outlet	1 beep	No light if a cassette is not inserted, or red light if a cassette is inserted	Without gel cassette: last program used (EP or EG) With gel cassette in: last time setting
Ready (with no current flowing through gel)	Gel cassette inserted into a base	—	Steady red	Default time setting (12 minutes for EG, 14 minutes for EP, or last time setting)
Run	Press and release the pwr/prg button	—	Steady green	Count down time
End of run	Automatic	Continuous beeping for 2 minutes followed by a single beep every minute	Flashing red until the time button is pressed	Negative time display (00 to -19 minutes)
Run ends after an external power failure	Automatic	Continuous beeping for 2 minutes followed by a single beep every minute	Alternating red and green	Negative time display (00 to -19 minutes)
Pause (manually end the run)	Press and release the pwr/prg button during the run	—	With gel cassette in: steady red Without gel cassette: no light	Flashing time display

E-Base™ Quick Reference Guide, Continued

Mode	Action	Sound	Light	Digital Display
Return to Ready mode after an automatic stop	Press and release the pwr/prg button	—	Steady red	Last time setting
Restart after a manual stop	Press and release the pwr/prg button	—	Steady green	Count down time
Return to Ready mode after a manual stop	Press and hold the pwr/prg button	—	With gel cassette in: steady red Without gel cassette: no light	With gel cassette in: last time setting Without gel cassette: last program setting
Failure	Press and hold pwr/prg button for 2 seconds and remove gel from the base	Continuous loud beeping	—	Flashing “ER”
No cassette	—	—	—	EP, last program used (EP or EG)
Time setting	With gel cassette in: Press and release the time button	—	With gel cassette: steady red	Time increases by 1 minute increments
	With and without gel cassette: Press and hold the time button	—	With gel cassette in: steady red Without gel cassette: no light	Time increases continuously and automatically stops at 00
Program setting	Press and release the pwr/prg button when no cassette is inserted into the E-Base™ to select the desired program	1 beep	No light	Selected program EP or EG

Opening the E-PAGE™ Cassette

Opening E-PAGE™ 48 or E-PAGE™ 96 Cassettes with the Gel Knife

To remove the E-PAGE™ 48 or E-PAGE™ 96 Gel from the cassette for blotting or staining, use the Gel Knife to open the cassette.

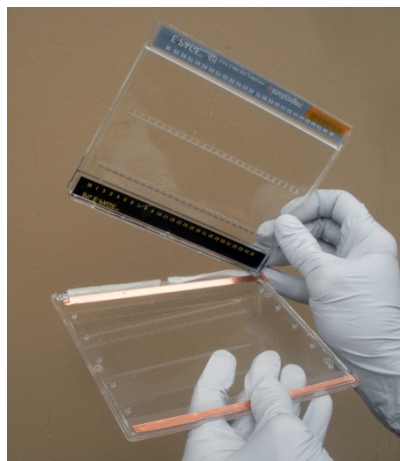
1. Separate each of the bonded sides of the E-PAGE™ cassette by inserting the Gel Knife into the gap between the two plastic plates that make up the cassette. The side of the cassette with the wells should face up.



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 (pages 25–37) or staining (pages 38–48).

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.

Blotting E-PAGE™ Gels

Blotting E-PAGE™ Gels with the iBlot® Device

Introduction

Instructions are provided in this section to assemble the iBlot® Gel Transfer Device with the De-Bubbling Roller for blotting E-PAGE™ Gels. The recommended program for blotting E-PAGE™ Gels is Program P3 at 20 volts for 7–8 minutes.

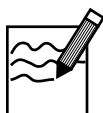
Materials Needed

You will need the following items:

- Pre-run E-PAGE™ gel containing your protein samples and standards
 - iBlot® Gel Transfer Stack, Regular (page 70)
-

Removing the Gel

Remove the gel from the cassette for transfer after completion of electrophoresis as described on page 25.



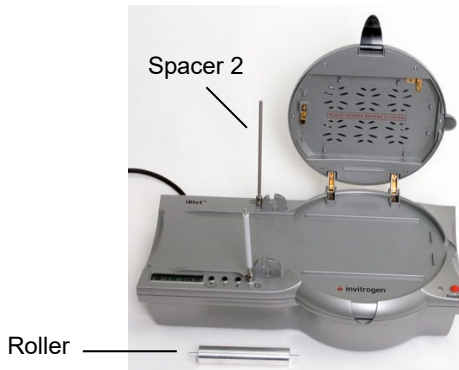
Note

- There is no need for any pretreatment of the gel after electrophoresis. Wash the E-PAGE™ gel briefly in deionized water to remove any small gel pieces attached to the gel.
 - The transfer membrane is supplied in a ready-to-use format in the stacks without any need for pretreatment. Do not treat the PVDF membrane with methanol as the PVDF membrane is preactivated prior to assembly with the transfer stack.
 - To obtain the best blotting results with the E-PAGE™ gels, we recommend that you use the De-bubbling Roller. For an alternative protocol using the Blotting Roller, refer to the manual for the iBlot® Dry Blotting System.
-

Using the iBlot[®] Device with the De-Bubbling Roller, Continued

Assembling the iBlot[®] Device

1. Open the lid of the device and pull up the Metal Spacers 1 and 2. If you have attached the De-bubbling Roller to the device, then remove the roller as shown in the figure below.



2. Remove the package labeled iBlot[®] Anode Stack, Bottom from the iBlot[®] Gel Transfer Stacks Box. Remove the laminated sealing of the iBlot[®] Anode Stack, Bottom and **keep the stack in the transparent plastic tray**.



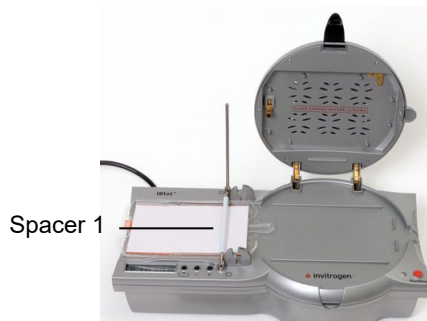
3. Place the iBlot[®] Anode Stack, Bottom stack **with the tray** to the left of the blotting surface area such that the tab on the tray is on the right side of the De-bubbling Roller, as shown below. Slide the bottom stack to the left until the stack is blocked by the Gel Barriers present on the left side of the device.
Note: Handle the iBlot[®] Anode Stack, Bottom using the plastic tray to avoid disturbing the gel and membrane layers in the stack. **Do not** touch the transfer membrane on the stack.



Using the iBlot® Device with the De-Bubbling Roller, Continued

Assembling the iBlot® Device, continued

4. Clean the Metal Spacer 1 with a damp cloth or tissue and place the spacer on the membrane as shown below.



5. Place the prerun gel containing your protein samples on Metal Spacer 1 such that the gel is aligned to the lower right corner of the bottom stack with the wells of the E-PAGE™ gel facing up.



6. Clean the Metal Spacer 2 with a damp cloth or tissue and place the spacer over the gel as shown below.

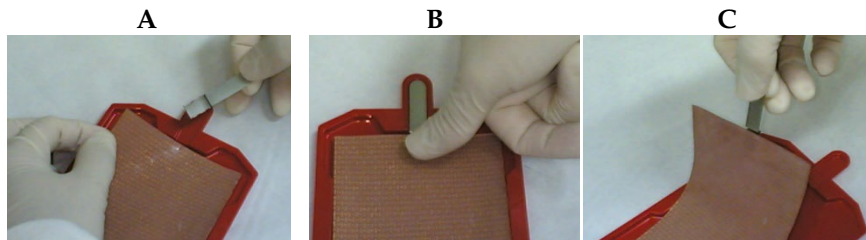


7. Remove the package labeled iBlot® Cathode Stack, Top from the iBlot® Gel Transfer Stacks Box. Remove the iBlot® Cathode Stack, Top from the package.
-

Using the iBlot® Device with the De-Bubbling Roller, Continued

Assembling the iBlot® Device, continued

8. Insert the steel iBlot® E-PAGE™ Tab in the plastic tray groove with the tab teeth facing up (figure A). Gently press the iBlot® Cathode Stack over the teeth to allow the teeth to penetrate into the copper electrode (figure B). Remove the iBlot® Cathode Stack, Top from the red plastic tray using the iBlot® E-PAGE™ Tab (figure C).



9. Place the iBlot® Cathode Stack, Top **without the tray** on top of Metal Spacer 2 with the copper electrode side facing up (and agarose side facing down). Ensure that all layers are aligned to the right to perform efficient de-bubbling.



10. Insert the De-bubbling Roller into the two grooves and lower the roller to its lowest location while holding the pull tab. The resulting assembly consists of the gel, and cathode and anode stacks placed between two Metal Spacers 1 and 2 with the De-bubbling Roller on top of the assembly as shown below.



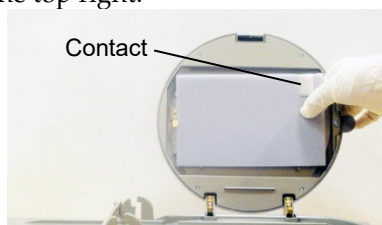
Using the iBlot® Device with the De-Bubbling Roller, Continued

Assembling the iBlot® Device, continued

11. Hold the iBlot® E-PAGE™ Tab and plastic tab on the iBlot® Anode Stack, Bottom together and pull the assembly (anode and cathode stacks, and gel) together through the De-bubbling Roller towards the blotting surface, in one smooth, uninterrupted movement until the assembly reaches the Gel Barriers on the blotting surface (figure A). At the end of de-bubbling, all layers are aligned to the right as shown below (figure B).



12. Place the iBlot® Disposable Sponge on the inner side of the lid (between the small protrusions on the lid that hold the sponge in its place) such that the metal contact is to the top right.



The sponge absorbs any excess liquid generated during blotting and exerts an even pressure on the stack surface.

Performing Blotting

After assembling the iBlot® Gel Transfer Device, perform blotting within 15 minutes of assembling the stacks with the gel as described below.

1. Close the iBlot® Lid and secure the latch. The red light is on indicating a closed circuit. Ensure the correct program is selected (Program P3 at 20 volts for 7–8 minutes).
2. Press the Start/Stop button to start the transfer. The red status light changes to green. The transfer continues using the programmed parameters.
3. At the end of the transfer, current automatically shuts off and the iBlot® Gel Transfer Device signals the end of transfer with repeated beeping sounds, and flashing red light and digital display.

Note: Previous versions of the iBlot® Gel Transfer Device (firmware versions prior to 2.7.9), signaled the end of transfer with repeated beeping sounds, and flashing **green** light (instead of red light) and digital display.

4. Press and release the Start/Stop button to stop the beeping. The light turns to a steady red light.
-

Using the iBlot® Device with the De-Bubbling Roller, Continued

Disassembling the iBlot® Gel Transfer Device

To obtain good transfer and detection results, disassemble the device and stacks within 30 minutes of ending the blotting procedure.

1. Open the lid of the iBlot® Device.
2. Remove the iBlot® E-PAGE™ Tab (used for blotting E-PAGE™ gels only). Rinse the tab with deionized water and store in a dry place for future use. **Do not discard the iBlot® E-PAGE™ Tab.**
3. Discard the iBlot® Disposable Sponge and iBlot® Cathode Stack, Top.
4. Carefully remove and discard the gel and filter paper (if used) as shown below. Remove the transfer membrane from the stack and proceed with the blocking procedure or stain the membrane (see next page for details).

Note: If you are using PVDF membranes, place the membrane immediately into water, as PVDF membranes dry quickly. If the PVDF membrane is dried, re-wet the membrane with methanol and rinse with deionized water a few times before use. Transfer the membrane to your blocking or staining solution only after you are sure that is completely wet, because reactivating after the membrane is exposed to the blocking solution may be problematic.



5. Discard the iBlot® Anode stack, Bottom.
6. At this point, the iBlot® Gel Transfer Device is ready for another run (no cooling period is required). If you are not using the device, turn off the power switch located on the back of the iBlot® Gel Transfer Device.



Important

Do not reuse the iBlot® Disposable Sponge, iBlot® Filter Paper, and iBlot® Cathode and Anode Stacks after blotting. Discard after each use.

Semi-Dry Blotting of E-PAGE™ Gels

Introduction

A semi-dry blotting procedure for blotting E-PAGE™ Gels is described in this section. You will need a semi-dry transfer apparatus that is capable of accommodating the dimensions of an E-PAGE™ Gel (8.6 cm x 13.5 cm) and a power supply.

Materials Needed

You will need the following items. See page 71 for ordering information.

- Semi-dry blotter
 - Methanol (for transfer of E-PAGE™ 48 Gels)
 - NuPAGE® Transfer Buffer (20X)
 - NuPAGE® Antioxidant
 - Blotting membranes: Invitrolon™ /Filter Paper Sandwiches or Nitrocellulose Membrane/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
-

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

E-PAGE™ 48 Transfer Buffer

We recommend using 2X NuPAGE® Transfer Buffer with **15% methanol** and NuPAGE® Antioxidant for the transfer of most proteins from E-PAGE™ 48 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 as follows:

Buffer Component	Transfer Buffer with 15% Methanol	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

E-PAGE™ 96 Transfer Buffer

We recommend using 2X NuPAGE® Transfer Buffer with NuPAGE® Antioxidant **without methanol** for the transfer of most proteins from E-PAGE™ 96 Gels.

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

Buffer Component	Transfer Buffer without Methanol
NuPAGE® Transfer Buffer (20X)	50 mL
NuPAGE® Antioxidant	0.5 mL
Deionized Water	to 500 mL

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

Equilibrating the Gel

Equilibration of the gel in 2X Transfer Buffer results in 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 24.
 2. Using 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. Perform this equilibration on a shaker.
-

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 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 in deionized water.
 3. Soak the membrane in 2X NuPAGE® Transfer Buffer (see previous page) for several minutes in the Incubation Tray.
-

Guidelines for Semi-Dry Blotting

- Use the correct 2X NuPAGE® Transfer Buffer as described on the previous page based on the E-PAGE™ Gel type.
 - Equilibrate the gel as described above.
 - To ensure a good transfer, it is necessary to use the E-PAGE™ Blotting Pad and 2.5-mm Blotting Filter Paper.
 - Remove any trapped bubbles by using the Blotting Roller at each step of creating the blot sandwich.
 - Since the E-PAGE™ Gel is thick and easily compressed, avoid applying any off-center pressure that may deform the blot sandwich, cause bubble formation, or break the contact between the gel and membrane resulting in poor transfer.
 - If the blot sandwich is disturbed during assembly, we recommend disassembling the sandwich and re-assembling it again.
-

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

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 32). 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 32). Inspect both sides of the pad for air bubbles. Press the pad to ensure the elimination of any visible air bubbles.
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 (see page 32) 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.
 6. Place the flat side of the gel on top of the blotting membrane (well side up), and remove any air bubbles with the Blotting Roller. Fill the wells of the gel with 2X NuPAGE® Transfer Buffer (see page 32).
 7. Place the pre-soaked E-PAGE™ Blotting Pad on the gel and gently roll out air bubbles with the Blotting Roller.
 8. Place 2 pieces of pre-soaked 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 trapped between filter paper sheets using 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. See page 53 for results obtained with semi-dry blotting of E-PAGE™ Gels.
-

Semi-Wet Blotting of E-PAGE™ Gels

Introduction

A semi-wet blotting procedure using an XCell II™ Blot Module for transfer of proteins from E-PAGE™ Gels is described below.

Semi-Wet Transfer Apparatus

Two XCell II™ Blot Module units are required to transfer one E-PAGE™ Gel. To obtain the best transfer, we recommend transferring only one half of an E-PAGE™ Gel in one XCell II™ Blot Module. Inefficient transfer may occur if you transfer both halves of the E-PAGE™ Gel in one blot module at the same time, as the E-PAGE™ Gel is thicker than regular mini-gels.



Important

If you are using the XCell II™ Blot Module for transferring one E-PAGE™ Gel, we recommend that you have two XCell II™ Blot Modules on hand before proceeding for transfer.

Other semi-wet transfer apparatus that can accommodate an E-PAGE™ Gel (8.6 cm x 13.5 cm) may be used. However you may get sub-optimal results if your blotting apparatus is prone to transfer distortion when accommodating thick gels. It may be necessary to optimize your semi-wet blotting procedure based on the protocol below, or it may be necessary to use a semi-dry blotting apparatus to obtain optimal results.

Note: Using the Criterion Blotter module from BioRad (BioRad Cat. no. 170-4070) is not recommended.

Materials Needed

You will need the following items. Ordering information is on page 71.

- XCell II™ Blot Module (2 units required for transfer of 1 E-PAGE™ Gel)
- XCell *SureLock*™ Mini-Cell
- Power supply (1–2 amps)
- Methanol
- NuPAGE® Transfer Buffer (20X)
- NuPAGE® Antioxidant
- Invitrolon™ or Nitrocellulose Membrane/Filter Paper Sandwiches
- Blotting Roller
- Incubation Tray

Note: Use the blotting pads that fit the XCell II™ Blot Module. **Do not** use the E-PAGE™ Blotting Pads for semi-wet transfer.

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

Preparing 1X Transfer Buffer

We recommend using 1X NuPAGE® Transfer Buffer with NuPAGE® Antioxidant and 10% methanol for semi-wet transfer of proteins from the E-PAGE™ Gel.

Note: If a large amount of heat is produced during the transfer, use a cooling unit or perform transfer with cold buffer.

For one blot module, prepare 1000 mL of 1X NuPAGE® Transfer Buffer as follows:

Buffer Component	Transfer Buffer with 10% Methanol
NuPAGE® Transfer Buffer (20X)	50 mL
NuPAGE® Antioxidant	1 mL
Methanol	100 mL
Deionized Water	849 mL

Preparing Blotting Pads, Membrane, and Gel

1. Use ~ 700 mL 1X NuPAGE® Transfer Buffer (see previous page) to soak the blotting pads of the blot module until saturated. Remove air bubbles by squeezing the blotting pads while they are submerged in buffer.
Note: Use the blotting pads that fit the XCell II™ Blot Module. **Do not** use the E-PAGE™ Blotting Pads for semi-wet transfer.
 2. Prepare pre-cut membrane/filter paper sandwiches as described below.
 - **PVDF membrane:** Pre-wet the PVDF membrane for 30 seconds in methanol, ethanol, or isopropanol. Briefly rinse in deionized water and place the membrane in a shallow dish containing 50–100 mL 1X NuPAGE® Transfer Buffer for several minutes.
 - **Nitrocellulose membrane:** Place the membrane directly into a tray containing 1X NuPAGE® Transfer Buffer for several minutes.
 - **Filter paper:** Soak briefly in 1X NuPAGE® Transfer Buffer before using.
 3. Transfer the E-PAGE™ Gel immediately following electrophoresis. You will need to cut the gel in half so it can fit into the XCell II™ Blot Module. **Do not** soak the E-PAGE™ Gel in transfer buffer.
-

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

Semi-Wet Blotting Procedure

1. Remove any small pieces of gel from the well areas by gently rubbing a gloved finger over the well side of the gel. The gel pieces on the gel may cause air bubbles and field distortion during transfer.
 2. Lay the gel on a flat surface, well side up, in a suitable container and pour sufficient 1X NuPAGE® Transfer Buffer (see previous page) over the gel to fill all wells.
 3. Place a piece of pre-soaked filter paper on top of the gel and remove any trapped air bubbles by gently rolling the Blotting Roller over the filter paper. Turn the gel over so the filter paper and gel are facing downwards over a gloved hand or clean flat surface
 4. Place a pre-soaked membrane on the gel and remove any trapped air bubbles with the Blotting Roller.
 5. Place another pre-soaked filter paper on top of the membrane and remove any trapped air bubbles with the Blotting Roller.
 6. Place two pre-soaked blotting pads into the cathode core of the blot module.
Note: Use the blotting pads that fit the XCell II™ Blot Module. **Do not** use the E-PAGE™ Blotting Pads for semi-wet transfer.
 7. Place the gel/membrane assembly on the blotting pads, such that the gel is closest to the cathode core.
 8. Add another pre-soaked blotting pad on top of the membrane assembly. Remove air bubbles with the Blotting Roller.
 9. Add enough pre-soaked blotting pads to rise to 0.5 cm over the rim of the cathode core. Place the anode (+) core on top of the pads.
 10. Hold the blot module together firmly and slide it into the guide rails on the lower buffer chamber.
 11. Insert the gel tension wedge into the lower buffer chamber and lock the wedge into position.
 12. Fill the blot module with 1X NuPAGE® Transfer Buffer (see page 36) until the gel/membrane assembly is just covered.
 13. Fill the outer buffer chamber with 650 mL deionized water.
 14. Place the lid on the unit and connect electrical leads to the power supply.
 15. Perform transfer for nitrocellulose or PVDF membranes using 35 V constant for 1 hour. The expected starting current is about 170 mA.
-

Visualizing and Staining of E-PAGE™ Gels

Choosing a Staining Protocol

A summary of the time needed and sensitivity obtained with E-PAGE™ gels using various staining protocols is described below. Based on the starting material and your requirements for sensitivity and background, choose an appropriate staining protocol.

For background: ++++ (very low background), +++ (low background), ++ (moderate background).

Protocol	Time	Sensitivity	Background
Lumio™ Green Detection Reagent	Pre-label proteins during sample preparation, run the gel, and visualize (total time: ~1 hour).	1–5 pmole	++++
SYPRO® Ruby Protein Gel Stain	Fix: 0.5 h Stain: 1.5 h Destain: 2 h-overnight	1–2.5 ng	++++
Coomassie R-250 Stain	Stain: 1.5 h Destain: overnight	20 ng	+++
Coomassie R-250 Stain Microwave Protocol	Stain: 0.5 h Destain: overnight	20 ng	+++
SimplyBlue™ SafeStain	Fix: 0.5 h Stain: 0.5 h–2 h Destain: 3 h-overnight	20 ng	+++
SilverQuest™ Silver Stain	Total: 4.5 h	10 ng	++++
SilverXpress® Silver Stain	Total: 6.5 h	1 ng	++
InVision™ His-tag In-Gel Stain	Blot: 7–8 min Rinse: 2 min Stain: 20 min	40 ng His-tagged protein	+++

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 72)

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 54.

SYPRO® Ruby Staining Protocol

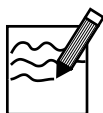
Introduction

Instructions for staining E-PAGE™ Gels using SYPRO® Ruby Protein Gel Stain is described in this section. To visualize protein bands after electrophoresis using SYPRO® Ruby, you will need a UV transilluminator or a laser-based scanner (see below). For further details on imaging SYPRO® Ruby proteins, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 72). **Total staining time is 4 hours, with optional destaining overnight.**

Materials Needed

You will need the following items for staining one E-PAGE™ Gel. See page 71 for ordering information.

- SYPRO® Ruby Protein Gel Stain (see note below)
 - Fixing solution (20% acetic acid, see note below)
 - Destaining solution (10% methanol, 7% acetic acid, see note below)
 - Incubation Trays
 - Rotary shaker
 - UV transilluminator equipped with a standard camera or an appropriate laser scanner (see note below)
-



Note

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.

Staining Protocol

1. After electrophoresis, remove the gel from the cassette (page 24) and place the gel in a clean Incubation Tray.
2. Fix the gel in 20% acetic acid for 30 minutes on an orbital shaker.
3. Stain the gel in undiluted SYPRO® Ruby Protein Gel Stain for 1.5 hours on an orbital shaker.
4. Transfer the gel to a clean Incubation Tray and destain in 10% methanol, 7% acetic acid for 2 hours. If complete removal of background is desired, perform the destaining step overnight.
5. Place the gel on a UV transilluminator equipped with a standard camera and select the ethidium bromide 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 (610 nm).
6. Image the gel with a suitable camera with the appropriate filters using a 1–4 second exposure. You may need to adjust the brightness and contrast to reduce any faint non-specific bands.

You should see fluorescent protein bands and the gel should have minimal background as shown on page 55.

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.

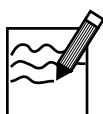
Materials Needed

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)
-



Note

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.



Important

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 24) 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 24) 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 41). **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 56.

SimplyBlue™ SafeStain

Introduction

Instructions for staining E-PAGE™ Gels using SimplyBlue™ SafeStain are described in this section. For further details, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 72). **The total staining time for SimplyBlue™ SafeStain is 0.5–3 hour plus overnight destaining.** Results obtained with SimplyBlue™ SafeStain are shown on page 57.

Materials Needed

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

- Clean Incubation Tray
 - Rotary shaker
 - SimplyBlue™ SafeStain (see note below)
 - Fixing solution: (see note below)
 - For E-PAGE™ 48: 10% acetic acid/30% methanol in deionized water
 - For E-PAGE™ 96: 20% acetic acid in deionized water
 - Wash solution: (see note below)
 - For E-PAGE™ 48: 2% NaCl in deionized water
 - For E-PAGE™ 96: Deionized water
-



Note

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.

E-PAGE™ 48 and E-PAGE™ 96 Gels require different staining protocols. Be sure to use the correct protocol for your gel.

E-PAGE™ 48 Gels

E-PAGE™ 48 Gel Staining Protocol:

1. After electrophoresis, remove the gel from the cassette (page 24) and place the gel in a clean incubation tray.
 2. Fix the gel in fixing solution (10% acetic acid/30% methanol in deionized water) for 1 hour with gentle shaking at room temperature. Decant the fixing solution.
 3. Stain the gel with sufficient SimplyBlue™ SafeStain to cover the gel. Incubate at room temperature for 2 hours with gentle shaking. Decant the stain.
 4. Wash the gel briefly with deionized water.
 5. Wash the gel with 2% NaCl solution with gentle shaking at room temperature with intermittent changes of wash solution until the bands are visible or overnight for maximum sensitivity and clear background.
-

SimplyBlue™ SafeStain, Continued

E-PAGE™ 96 Gels For routine staining of E-PAGE™ 96 Gels, use Protocol A. To obtain the clearest background for imaging, use Protocol B, which includes a 12–24 hour washing step to improve the background.

Protocol A

1. After electrophoresis, remove the gel from the cassette (page 24) and place the gel in a clean incubation tray.
2. Rinse the gel 3 times for 5 minutes each with deionized water to remove SDS and buffer salts.
3. Stain the gel with sufficient SimplyBlue™ SafeStain to cover the gel. Incubate at room temperature for 1.5 hours with gentle shaking. Decant the stain.
4. Wash the gel with deionized water for 3 hours with intermittent water changes.

Protocol B

1. After electrophoresis, remove the gel from the cassette (page 24) and place the gel in a clean incubation tray.
 2. Fix the gel in 20% acetic acid for 30 minutes at room temperature. Decant the solution.
 3. Stain the gel with sufficient SimplyBlue™ SafeStain to cover the gel. Incubate at room temperature for 30 minutes with gentle shaking. Decant the stain.
 4. Rinse the gel briefly with deionized water. Wash the gel in deionized water for 12–24 hours at room temperature, changing the water at least once.
-

Silver Staining Methods

Introduction

Instructions for staining E-PAGE™ Gels using SilverQuest™ Silver Staining Kit or SilverXpress® Silver Staining Kit are described in this section.

Total staining time with SilverQuest™ Silver Staining Kit is 4.5 hours and SilverXpress® Silver Staining Kit is 6.5 hours.

Materials Needed

You will need the following items for staining one E-PAGE™ Gel (see page 71 for ordering information):

- Ultrapure water (>18 megohm/cm resistance recommended, see below)
- Clean staining containers (if using the SilverQuest™ Silver Staining Kit, make sure the container is **microwave safe**. Do NOT use the Incubation Tray cat no. LC2102 for this purpose)
- Rotary shaker
- Clean glass bottles for reagent preparation

For SilverQuest™ Silver Staining:

- SilverQuest™ Silver Staining Kit
- Microwave oven (700–1000 watts)
- 30% ethanol (made with ultrapure water)
- 100% ethanol
- Fixative (40% ethanol, 10% acetic acid, made with ultrapure water)

For SilverXpress® Silver Staining:

- SilverXpress® Silver Staining Kit
 - Methanol
 - Acetic acid
-



- Always use ultrapure water of >18 megohm/cm resistance for preparing all solutions and rinsing gels and containers. Poor water quality may increase the background or impair band development.
 - To achieve best results, be sure to keep the volume of all solutions and incubation time of all steps exactly as given in the protocol. Changes in the protocol can result in high background or poor band development.
 - When using a microwave oven for staining E-PAGE™ Gels with the SilverQuest™ Silver Staining Kit, always use the microwave oven at full power setting. Always use a large volume of buffer (2-fold) and a microwave safe container. **Avoid microwaving the E-PAGE™ Gel for more than 45 seconds because gel decomposition can occur.**
-

Silver Staining Methods, Continued

SilverQuest™ Silver Staining Protocol

A brief protocol for staining one E-PAGE™ Gel with the SilverQuest™ Silver Staining Kit is described below. For details, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 72).

After electrophoresis, remove the gel from the cassette (page 24). For all staining and washing steps described below, be sure to use sufficient volume of reagent to completely cover the gel in a **microwave safe** container such that the gel moves freely during all staining and washing steps.

Results obtained with SilverQuest™ Silver Staining Kit are shown on page 58.

Step	Reagent	Fast Procedure
Fix	Ethanol	80 mL
	Acetic Acid	20 mL
	<u>Ultrapure water</u>	<u>100 mL</u>
	Final Volume	200 mL
Wash	Ethanol	60 mL
	<u>Ultrapure water</u>	<u>140 mL</u>
	Final Volume	200 mL
Sensitize	Ethanol	30 mL
	Sensitizer	10 mL
	<u>Ultrapure water</u>	<u>60 mL</u>
	Final Volume	100 mL
First Wash	Ultrapure water	150 mL
Second Wash	Ultrapure water	150 mL
Stain	Stainer	1 mL
	<u>Ultrapure water</u>	<u>99 mL</u>
	Final Volume	100 mL
Wash	Ultrapure water	100 mL
Develop	Developer	10 mL
	Developer Enhancer	1 drop
	<u>Ultrapure water</u>	<u>90 mL</u>
	Final Volume	100 mL
Stop	Stopper	10 mL
Wash	Ultrapure water	100 mL

Silver Staining Methods, Continued

SilverXpress® Silver Staining Protocol

A brief protocol for staining one E-PAGE™ Gel with the SilverXpress® Silver Staining Kit is described below. For details, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 72).

After electrophoresis, remove the gel from the cassette (page 24). For all staining and washing steps described below, be sure to use sufficient volume of reagent to completely cover the gel in a suitable container such that the gel moves freely during all staining and washing steps.

Results obtained with SilverXpress® Silver Staining Kit are shown on page 59.

Note: The final volumes of solutions containing both methanol and water reflect a volume shrinkage that occurs when these two reagents are mixed. Do not adjust volumes of components or final volume.

Step	Reagent	Procedure
Fix	Ultrapure water 270 mL Methanol 300 mL <u>Acetic Acid</u> 60 mL Final Volume 600 mL (see note above)	Fix the gel in 3 changes (~200 mL each) of fixing solution for 90 minutes. Change the fixing solution every 30 minutes.
Sensitize	Ultrapure water 105 mL Methanol 100 mL <u>Sensitizer</u> 5 mL Final Volume 200 mL (see note above)	Incubate the gel in 2 changes of Sensitizing Solution for 1 h, each. Decant the solution.
Wash	Ultrapure water 200 mL	Wash the gel twice with ultrapure water for 15 minutes each.
Stain	Stainer A 5 mL Stainer B 5 mL <u>Ultrapure water</u> 90 mL Final Volume 100 mL	Incubate the gel in Staining Solution for 30 minutes. Decant the solution.
Wash	Ultrapure water 200 mL	Wash the gel twice with ultrapure water for 10 minutes each.
Develop	Developer 5 mL <u>Ultrapure water</u> 95 mL Final Volume 100 mL	Incubate the gel in Developing Solution for 1–5 minutes.
Stop	Stopping Solution 5 mL Add the Stopping Solution directly to the Developing Solution when the desired staining intensity is reached.	Incubate the gel in Stopping solution for 20 minutes. Decant the solution.
Wash	Ultrapure water 200 mL	Wash the gel three times with ultrapure water for 20 minutes each.

InVision™ His-tag Protein Gel Stain

Introduction

For detection of His-tagged proteins, use the following protocol. Due to the thickness of E-PAGE™ Gels, we recommend transferring proteins onto nitrocellulose membrane (see pages 25-37) and then detecting them using InVision™ His-tag In-Gel Stain. For details, refer to the product manual available at www.lifetechnologies.com or contact Technical Support (page 72)

Note: This procedure is **not recommended** for staining PVDF membranes.

Materials Needed

You will need the following items for staining one E-PAGE™ Gel. See page 71 for ordering information.

- InVision™ His-tag In-gel Stain
 - Ultrapure water (>18 megohm/cm resistance recommended)
 - Incubation Tray
 - Rotary shaker
 - UV transilluminator equipped with a standard camera, or laser-based scanner.
-

Staining Procedure

1. After electrophoresis, remove the gel from the cassette (page 24) and blot proteins onto nitrocellulose membrane (page 25).
2. Rinse the nitrocellulose membrane (containing the transferred proteins) with deionized water for 2 minutes.
3. Stain the nitrocellulose membrane with 20 mL of ready-to-use InVision™ His-tag In-gel Stain for 20 minutes at room temperature.
4. Rinse the membrane briefly with ultrapure water.
5. Place the membrane on a UV transilluminator equipped with a camera. Visualize and image the membrane by exposing the membrane to UV light from the bottom or from the top (you may place the UV transilluminator on its side to illuminate the blot or use epi-illumination) for 4–8 seconds.

You may also use a laser-based scanner with the appropriate filters to visualize and image the membrane.

Drying of E-PAGE™ Gels

Gel Drying

Introduction

We recommend drying E-PAGE™ Gels using passive air-drying methods such as the Large Gel Drying Kit. Due to the thickness of E-PAGE™ Gels, vacuum drying is not recommended and may cause the gel to crack.

Materials Needed

You will need the following items for drying one E-PAGE™ Gel. See page 71 for ordering information.

- Large Gel Drying Kit
 - Gel-Dry™ Drying Solution (or prepare your own gel drying solution containing 30% methanol and 5% glycerol)
 - Rotary shaker
 - Staining Tray
-

Gel Drying Protocol

A brief gel drying protocol using the Large Gel Drying Kit is provided below. For more details on this system, refer to the Large Gel Drying Kit manual available at www.lifetechnologies.com or contact Technical Support (page 72).

1. At the end of all staining and destaining steps, wash the E-PAGE™ Gel three times for 2 minutes each time in deionized water (100 mL) on a rotary shaker.
 2. Decant the water and add fresh Gel-Dry™ Drying Solution (100 mL per E-PAGE™ gel).
 3. Equilibrate the gel in the Gel-Dry™ Drying Solution by shaking the gel for 15–20 minutes in the staining container. **Note:** Do not equilibrate gels stained with Coomassie R-250 in the Gel-Dry™ Drying Solution for more than 5 minutes to avoid losing band intensity.
 4. Cut any rough edges off the gel (including the wells and the gel foot) using a gel knife.
 5. Remove 2 pieces (per gel) of cellophane from the package.
 6. Immerse one sheet at a time in the Gel-Dry™ Drying Solution. Allow 10 seconds for complete wetting before adding additional sheets. Do not soak the cellophane for more than 2 minutes.
-

Gel Drying, Continued

Gel Drying Protocol, continued

7. Place the solid square of the Large Gel Drying Frame on a paper towel with the corner pins facing up.
 8. Lay a piece of pre-wetted cellophane from Step 5 over the frame carefully without trapping any air bubbles.
 9. Lay the E-PAGE™ Gel on the center of the cellophane sheet making sure no bubbles are trapped between the gel and the cellophane. Add some Gel-Dry™ Drying Solution to the surface of the cellophane, if necessary.
 10. Carefully lay the second sheet of cellophane over the gel so that no bubbles are trapped between the cellophane and the gel. Add some Gel-Dry™ Drying Solution if necessary. Gently smooth out any wrinkles in the assembly with a gloved hand or pipette.
 11. Place the plastic frame, beveled side up, on top of the cellophane. Push the plastic clamps onto three edges of the frame. Tilt the frame up on the remaining unclamped edge to drain excess solution, and then install the final clamp.
 12. Lay the gel dryer assembly flat on a bench top where it will be undisturbed. Be careful to avoid drafts as they can cause an uneven rate of drying which leads to cracking. Drying will take between 4–5 days.
 13. When the cellophane covering the gel is dry to the touch, remove the gel/cellophane sandwich from the drying frame. Trim off the excess cellophane.
 14. Press the dried gel(s) between the pages of a notebook under light pressure for approximately 2 days. Gels will then remain flat for scanning, photography, display, and overhead projection.
-

Analysis and Troubleshooting

Molecular Weight Calibration

Molecular Weight Calibration

The molecular weight of a protein can be determined based upon its relative mobility by constructing a standard curve with protein standards of known molecular weights.

The protein mobility in SDS gels is dependent on:

- Length of the protein in its fully denatured state
- Secondary structure of the protein
- Buffer system used

An identical molecular weight standard may have slightly different mobility resulting in different apparent molecular weight when run in different gel buffer systems.

If you are using molecular weight standards from Life Technologies, see the apparent molecular weights of these standards on E-PAGE™ Gels listed under **Assigned Apparent Molecular Weights** to determine the apparent molecular weight of your protein.

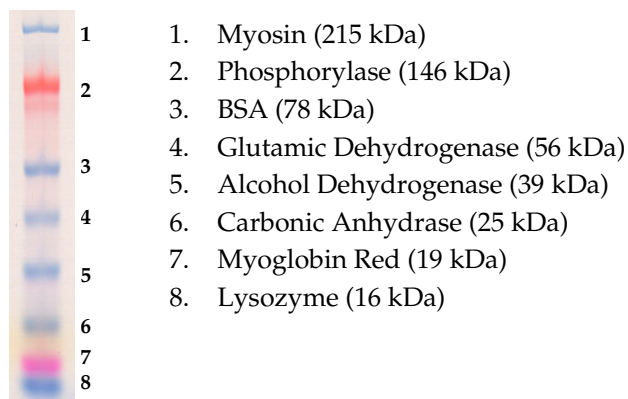
You may need to generate calibration curves in your lab with any other manufacturer's standards.

Assigned Apparent Molecular Weights

The apparent molecular weight values shown below are derived from the construction of a calibration curve in the E-PAGE™ 48 and 96 buffer systems. Use the values listed in this section for the most accurate calibration of your protein on an E-PAGE™ Gel.

SeeBlue® Plus2 Pre-stained Standard

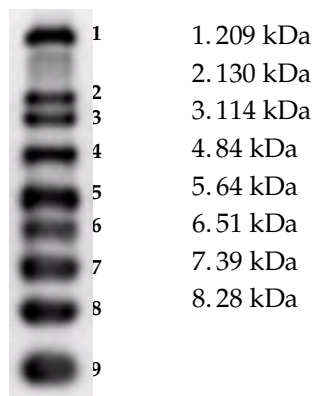
Molecular Weight on E-PAGE™ 48 8% Gel



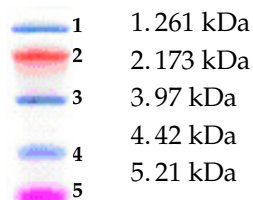
Molecular Weight Calibration, Continued

**Assigned Apparent
Molecular Weights,
Continued**

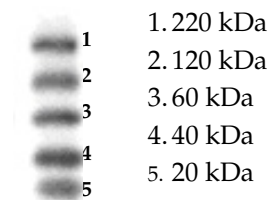
**MagicMark™XP Western Protein Standard
Molecular Weight on E-PAGE™ 48 8% Gel**



**E-PAGE™ SeeBlue® Pre-stained
Protein Standard Molecular Weight
on E-PAGE™ 96 6% Gel**



**E-PAGE™ MagicMark™ Western
Protein Standard Molecular Weight
on E-PAGE™ 96 6% Gel**



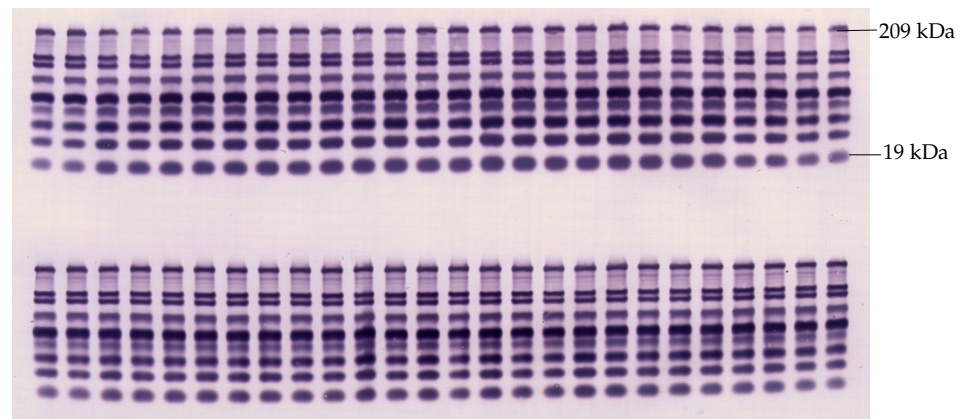
Expected Results

Western Blotting

Results obtained using an E-PAGE™ 48 8% Gel that was blotted using the semi-dry protocol described in this manual are shown below.

MagicMark™XP Western Protein Standard (5 µL) was loaded in all sample and marker wells. The gel was electrophoresed for 23 minutes using standard conditions. Proteins were transferred to a 0.45-µm nitrocellulose membrane using the semi-dry blotting protocol described in the manual.

Detection was performed with WesternBreeze® Anti-Mouse Immunodetection Kit using a 1:1000 dilution of Anti-Xpress™ primary antibody from Life Technologies. Both top and bottom rows are identical.

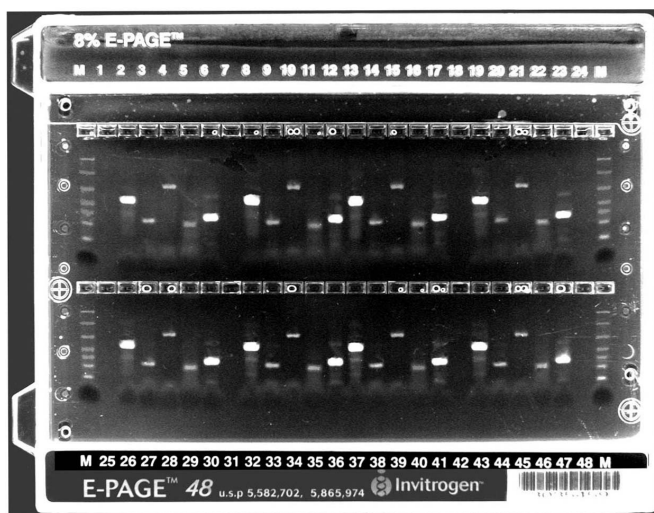


Expected Results, Continued

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):

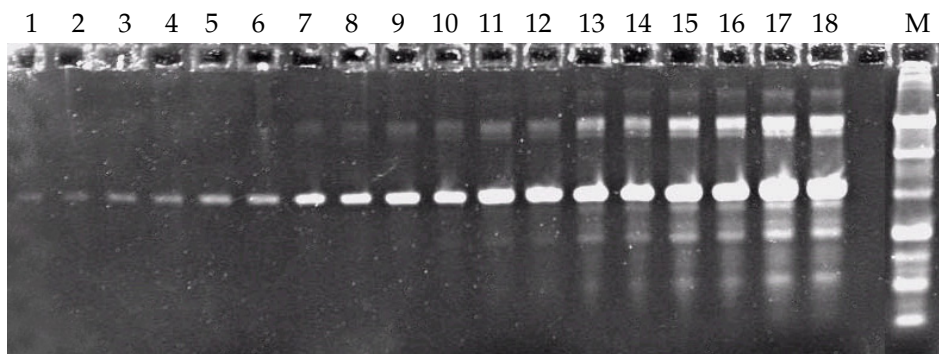
Lane	Sample
M	BenchMark™ Fluorescent Marker (5 µL)
2, 8, 13, 19, 26, 32, 37, 43	Human kinase Lumio™ fusion protein (10 µL)
3, 9, 14, 20, 27, 33, 38, 44	<i>E. coli</i> CAT Lumio™ fusion protein (10 µL)
4, 10, 15, 21, 28, 34, 39, 45	<i>E. coli</i> GUS Lumio™ fusion protein (10 µL)
5, 11, 16, 22, 29, 35,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

Expected Results, Continued

SYPRO® Ruby Protein Gel Stain

Results obtained using an E-PAGE™ 48 8% Gel stained with SYPRO® Ruby are shown below.

BSA (2.5 - 1000 ng) was run on an E-PAGE™ 48 8% Gel and stained with SYPRO® Ruby Protein Gel Stain as described in this manual.



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

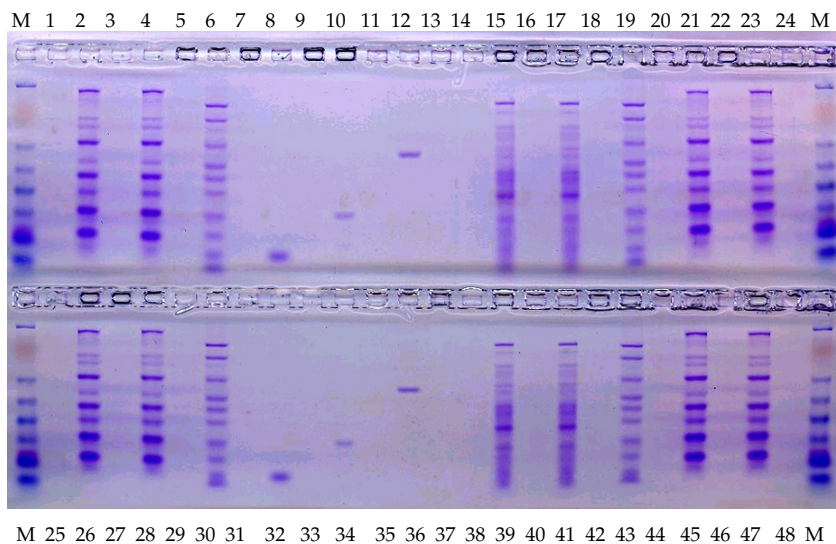
Lane	Amount of BSA
1, 2	2.5 ng
3, 4	5.0 ng
5, 6	10 ng
7, 8	25 ng
9, 10	50 ng
11, 12	100 ng
13, 14	250 ng
15, 16	500 ng
17, 18	1000 ng
M	BenchMark™ Fluorescent Marker (5 µL)

Expected Results, Continued

Coomassie® R-250 Stain

Results obtained using an E-PAGE™ 48 8% Gel stained with Coomassie R-250 Stain are shown below.

Samples were run on an E-PAGE™ 48 8% Gel and stained with Coomassie R-250 Stain using the microwave protocol as described in this manual.



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

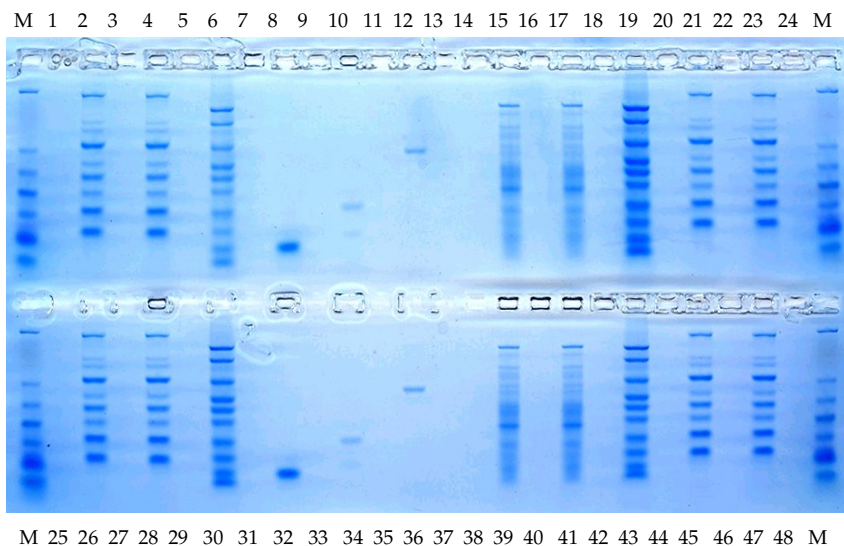
Lane	Sample
M	SeeBlue® Plus2 Pre-stained Standard (5 µL)
2, 4, 21, 23, 26, 28, 45, 47	MagicMark™ XP Western Protein Standard (10 µL)
6, 19, 30, 45	BenchMark™ His-tagged Protein Standard (10 µL, diluted 1:5)
8, 32	Lysozyme (200 ng)
10, 34	Carbonic anhydrase (200 ng)
12, 36	BSA (100 ng)
15, 17, 39, 41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

Expected Results, Continued

SimplyBlue™ SafeStain

Results obtained using an E-PAGE™ 48 8% Gel stained with SimplyBlue™ SafeStain are shown below.

Samples were run on an E-PAGE™ 48 8% Gel and stained with SimplyBlue™ SafeStain as described in this manual.



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

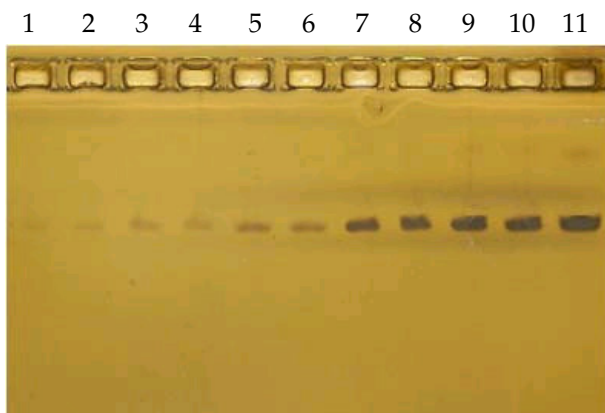
Lane	Sample
M	SeeBlue® Plus2 Pre-stained Standard (5 µL)
2, 4, 21, 23, 26, 28, 45, 47	MagicMark™ XP Western Protein Standard (10 µL)
6, 19, 30, 45	BenchMark™ His-tagged Protein Standard (10 µL, diluted 1:5)
8, 32	Lysozyme (200 ng)
10, 34	Carbonic anhydrase (200 ng)
12, 36	BSA (100 ng)
15, 17, 39, 41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

Expected Results, Continued

SilverQuest™ Silver Staining Kit

Results obtained using an E-PAGE™ 48 8% Gel stained with the SilverQuest™ Silver Staining Kit with are shown below.

BSA (0.5–1000 ng) was run on an E-PAGE™ 48 8% Gel and stained with the SilverQuest™ Silver Staining Kit as described in this manual.



The gel contains the following samples:

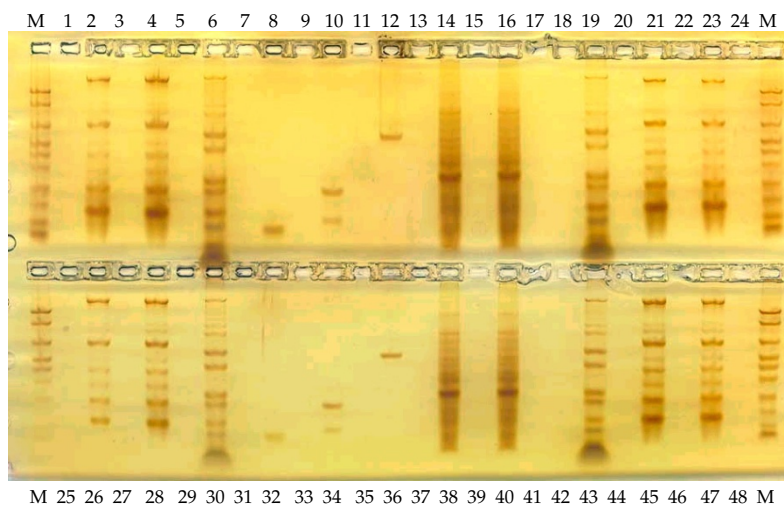
Lane	Amount of BSA
1	5 ng
2	5 ng
3	10 ng
4	10 ng
5	20 ng
6	20 ng
7	50 ng
8	50 ng
9	100 ng
10	100 ng
11	200 ng

Expected Results, Continued

SilverXpress™ Silver Staining Kit

Results obtained using an E-PAGE™ 48 8% Gel stained with the SilverXpress™ Silver Staining Kit with are shown below.

Samples were run on an E-PAGE™ 48 8% Gel and stained with the SilverXpress™ Silver Staining Kit as described in this manual.



The gel contains the following samples: (lanes 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, 38, 40	<i>E. coli</i> lysate (5 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. For issues with protein transfer, refer to the manual for the iBlot® Gel Transfer Device.

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 17).
Sample leaking from the wells	Sample is overloaded or wells are damaged	Be sure to load the recommended volume of sample per well (page 16). 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 10.
	Electrophoresis was not started immediately after sample loading	For best results, the gel should be run within 15 minutes of sample loading.

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 9. You may need to manually increase the run time for high salt or detergent samples to obtain optimal results.
	A1 tip not aligned	Be sure to align the A1 tip properly prior to automated loading of E-PAGE™ 96 Gel (page 18).
	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. Make sure transfer buffer contains no methanol if transferring E-PAGE™ 96.
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

Additional Staining Protocols

Pro-Q® Diamond Phosphoprotein Gel Stain

Instructions for specific staining of phosphoproteins on E-PAGE™ Gels using Pro-Q® Diamond Phosphoprotein Gel Stain is described in this section.

The total staining time is ~5 hours.



Note

Due to the thickness of E-PAGE™ Gels, the Pro-Q® Diamond Phosphoprotein Gel Stain **may not be optimally sensitive** for your particular staining application.

Materials Needed

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

- Pro-Q® Diamond Phosphoprotein Gel Stain
- Fixing solution (50% methanol, 10% acetic acid, made with deionized water)
- Pro-Q® Diamond Phosphoprotein Gel Destaining solution (or use 50 mM sodium acetate pH 4.0 containing 20% acetonitrile as destaining solution)
- Deionized water
- **Microwave safe** staining tray
- Rotary shaker
- Microwave oven (700–1000 watts)
- UV transilluminator equipped with a standard camera or laser scanners

See page 71 for ordering information.

Phosphoprotein Staining Protocol

For all steps described below, be sure to use sufficient volume of reagent to completely cover the gel using a suitable container such that the gel moves freely.

1. After electrophoresis, remove gel from the cassette (page 24) and place the gel in a clean **microwave safe** staining container.
 2. Fix the gel in fixing solution (50% methanol, 10% acetic acid in deionized water) for 45 minutes at room temperature with gentle shaking. Decant the fixing solution and add fresh fixing solution to the container and repeat the fixing step again for 45 minutes at room temperature. Decant the fixing solution.
 3. Add deionized water to the gel and microwave at high power for 45 seconds. Remove the gel from the microwave and wash the gel for 10 minutes at room temperature with gentle shaking. Decant the water and repeat the wash step again with deionized water. Decant the water.
 4. Stain the gel with Pro-Q® Diamond Phosphoprotein Gel Stain. Incubate the gel in the **DARK** for 1–2 hours at room temperature with gentle shaking. Decant the stain.
-

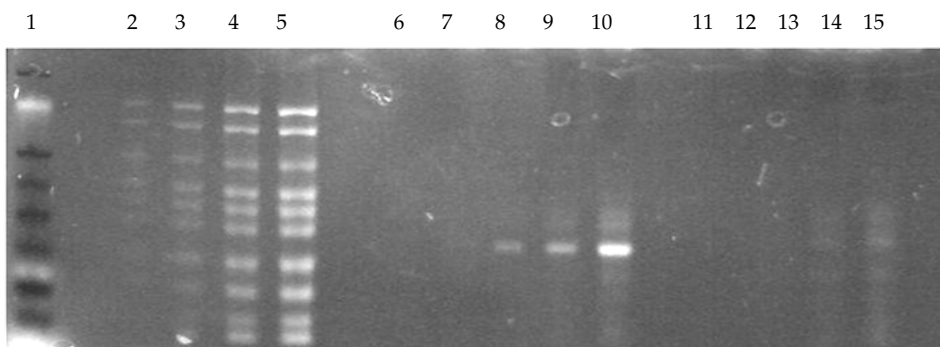
Additional Staining Protocols, Continued

Phosphoprotein Staining Protocol, Continued

5. Destain the gel with Pro-Q® Diamond Phosphoprotein Gel destaining solution twice for 30 minutes each time at room temperature with gentle shaking. You may wash the gel overnight to improve sensitivity and obtain clear background.
6. Place the gel on a UV transilluminator equipped with a standard camera and select an appropriate filter on the camera (stain emission maxima is 555–580 nm).
You may also use a laser-based scanner with a laser line that falls within the excitation maxima of the stain (580 nm).
7. Image the gel with a suitable camera with the appropriate filters using a 3–4 second exposure.

Results

Results obtained using an E-PAGE™ 48 8% Gel stained with Pro-Q® Diamond Phosphoprotein Gel Stain are shown below.



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

Lane	Sample
1	SeeBlue® Plus2 Pre-stained Standard (5 µL)
2, 3, 4, 5	1, 2.5, 5, and 10 µL, respectively, of BenchMark™ Fluorescent Protein Standard
6, 7, 8, 9, 10	<i>E. coli</i> lysate containing His-tagged urate oxidase protein, 35 kDa (0.05, 0.1, 0.25, 0.5, and 1 µg)
11, 12, 13, 14, 15	Negative control (<i>E. coli</i> lysate with non-His-tagged urate oxidase 0.05, 0.1, 0.25, 0.5, and 1 µg)

Additional Staining Protocols, Continued

Pro-Q® Sapphire 532 Oligohistidine Gel Stain

Instructions for specific staining His-tagged proteins on E-PAGE™ Gels using Pro-Q® Sapphire 532 Oligohistidine Gel Stain is described in this section.

The total staining time is ~5 hours.



Note

Due to the thickness of E-PAGE™ Gels, the Pro-Q® Sapphire 532 Oligohistidine Gel Stain **may not be optimally sensitive** for your particular staining application. **To visualize His-tagged proteins on E-PAGE™ Gels, we recommend the InVision™ His-tag In-Gel Stain (page 48).**

Materials Needed

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

- Pro-Q® Sapphire 532 Oligohistidine Gel Stain
 - Fixing solution (50% ethanol, 10% acetic acid, made with deionized water)
 - Deionized water
 - **Microwave safe** staining tray
 - Rotary shaker
 - Microwave oven (700–1000 watts)
 - UV transilluminator equipped with a standard camera or laser scanners
-

Oligohistidine Staining Procedure

For all steps described below, be sure to use sufficient volume of reagent to completely cover the gel using a suitable container such that the gel moves freely.

1. After electrophoresis, remove the gel from the cassette (page 24) and place the gel in a clean **microwave safe** staining container.
 2. Fix the gel in fixing solution (50% ethanol, 10% acetic acid in deionized water) for 45 minutes at room temperature with gentle shaking. Decant the fixing solution and add fresh fixing solution to the container and repeat the fixing step again for 45 minutes at room temperature. Decant the fixing solution.
 3. Add deionized water to the gel and microwave at high power for 45 seconds. Remove the gel from the microwave and wash the gel for 10 minutes at room temperature with gentle shaking. Decant the water and repeat the wash step for another 10 minutes again with deionized water. Decant the water.
 4. Stain the gel with Pro-Q® Sapphire 532 Oligohistidine Gel Stain. Incubate the gel in the **DARK** for 30–60 minutes at room temperature with gentle shaking. Decant the stain.
 5. Wash the gel with deionized water twice for 1 hour each time at room temperature with gentle shaking. You may wash the gel overnight to improve sensitivity and to obtain clear background.
-

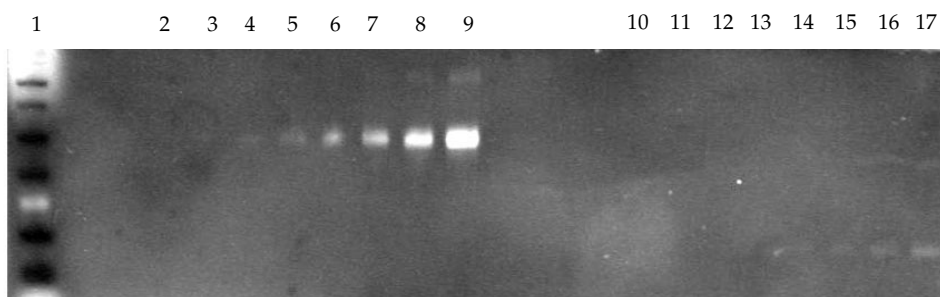
Additional Staining Protocols, Continued

Oligohistidine Staining Procedure, Continued

7. Place the gel on a UV transilluminator equipped with a standard camera and select an appropriate filter on the camera (stain emission maxima is 572 nm). You may also use a laser-based scanner with a laser line that falls within the excitation maxima of the stain (532 nm).
8. Image the gel with a suitable camera with the appropriate filters using a 1–4 second exposure.

Results

Results obtained using an E-PAGE™ 48 8% Gel stained with Pro-Q® Sapphire 532 Oligohistidine Gel Stain are shown below.



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

Lane	Sample
1	SeeBlue®Plus2 Pre-stained Standard (10 µL)
2, 3, 4, 5, 6, 7, 8, 9	Purified ovalbumin phosphoprotein, 35 kDa (5, 10, 25, 50, 100, 250, 500, 1000 ng)
10, 11, 12, 13, 14, 15, 16, 17	Negative control, lysozyme (5, 10, 25, 50, 100, 250, 500, 1000 ng)

Product Specifications

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: 3.2 cm

(one well to the next)

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.

E-PAGE™ 96 Gel Specifications

Each E-PAGE™ 96 gel contains 96 sample wells and 8 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: 20 µL

Well Opening: 3.8 mm × 1.8 mm

Running Distance: 1.6 cm

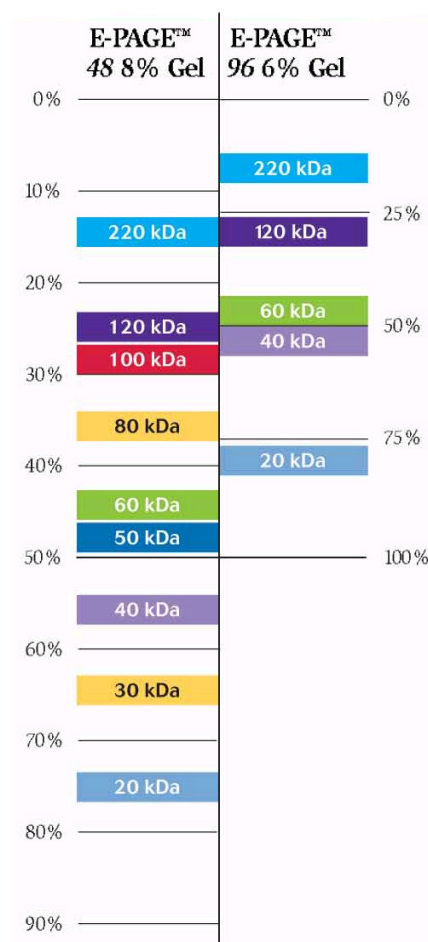
(one well to the next)

Space between Wells: 9 mm

Product Specifications, Continued

E-PAGE™ Gel Separation Range

The migration and resolution range of proteins run on E-PAGE™ 48 8% Gels and E-PAGE™ 96 6% Gels are shown.



E-Base™ Specifications

The specifications for the Mother E-Base™ and Daughter E-Base™ are listed below.

Dimensions:	14.6 cm (l) x 15 cm (w) x 5.3 cm (h)
Weight:	Mother E-Base™ - 370 g Daughter E-Base™ - 271 g
Safety:	Double Insulation, UL listed, and CE certified
Temperature:	Ambient 15°C to 40°C
Built-in Features:	Digital time display (00–99 minutes), alarm, light LED

The SBS (Society for Biomolecular Screening) standard 96-well plate format of the E-Base™ fits on most robotic platforms allowing the loading and electrophoresis of gels on the E-Base™ directly on the automated liquid handling system. See page 67 for an explanation of symbols and warnings used on the E-Base™.

Explanation of Symbols and Warnings

E-Base™



E189045

The Mother E-Base™ and Daughter E-Base™ comply with the Underwriters Laboratories Inc. regulation and the European Community Safety requirements. Operation of the E-PAGE™ bases is subject to the following conditions:

- Indoor use.
- Altitude below 2,000 meters.
- Temperature range: 5–40° C.
- Maximum relative humidity: 80%.
- Installation categories (over voltage categories) II; Pollution degree 2
- Mains supply voltage fluctuations not to exceed 10% of the nominal voltage (100–240V, 50/60Hz, 1500 mA).
- The Mother E-Base™ has been tested with up to 3 Daughter E-Bases™ connected at one time.
- Main plug is a disconnect device and must be easily accessible.
- Do not attempt to open E-Base™ devices. To honor the warranty, E-Base™ can only be opened and serviced by Life Technologies.
- The protection provided by the equipment may be impaired if the equipment is used in a manner not specified by Life Technologies.

Life Technologies Israel Ltd., is the manufacturer and owner of the UL file. For more information, contact:

Life Technologies Israel Ltd.
12 Hamada St.
P.O. Box 4035
Rehovot, Israel 74103



Caution

The **Caution** symbol denotes a risk of safety hazard. Refer to accompanying documentation.



Double Insulation

Class II product

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™ (Cat. no. EB-M03) is used for electrophoresis of one E-PAGE™ Gel
- The Daughter E-Base™ (Cat. no. EB-D03) attaches to the Mother E-Base™ 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

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

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.

Life Technologies and/or its affiliate(s) disclaim all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. 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.
