



Contents

Catalog Numbers EBM03, EBD03

Product	Cat. No.	Amount
Mother E-Base™ Device	EBM03	1 each
Daughter E-Base™ Device	EBD03	1 each



Product description

- The E-Base™ Device is an easy-to-use, programmable, automated small-footprint device combining a base and power supply to simplify electrophoresis of precast E-Gel™ 48, E-Gel™ 96, E-PAGE™ 48 and E-PAGE™ 96 gels.
- High-throughput and automation-compatible.
- Provides fast, safe, consistent, high-resolution electrophoresis.
- Eliminates the need to prepare agarose gels, buffers, and to stain gels.



Required materials

DNA analysis

- E-Gel™ 48 or 96 agarose gels (See [Gel selection guide](#))
- E-Gel™ DNA Ladder (See [Ladder selection guide](#))
- UltraPure™ DNase/RNase-Free Distilled Water (Cat. Nos. 10977015, 10977035)
- E-Gel™ Imager System with Blue Light Base (Cat. No. 4466612)
- (Optional) 1X E-Gel™ Sample Loading Buffer (Cat. No. 10482055)
- (Optional) Safe Imager™ 2.0 Blue-Light Transilluminator (Cat. No. G6600)

Protein Analysis

- E-PAGE™ 48 or 96 gels (See [Gel selection guide](#))
- E-PAGE™ SeeBlue™ Pre-stained Standard (Cat. No. LC5700)
- E-PAGE™ Loading Buffer 1 (Cat. No. EPBUF01)

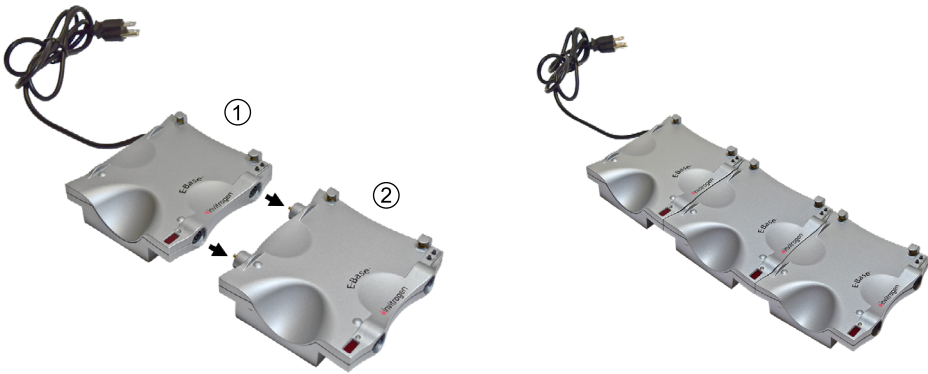


Online resources

- Visit our [product pages](#) for protocols, safety, and additional product information.
- Go online to view related [E-Gel™ products](#).
- For support, visit [thermofisher.com/support](#).

Connecting a Daughter E-Base™ Device

- For higher throughput, up to three E-Base™ Daughter units can be connected to an E-Base™ Mother Device.
- **IMPORTANT!** Ensure the Mother E-Base™ Device is unplugged before connecting any Daughter E-Base™ Devices.



① Mother E-Base™ Device with
② Daughter E-Base™ Device

Mother E-Base™ Device with
multiple Daughter E-Base™ Devices

i Troubleshooting

For detailed troubleshooting instructions see the E-Base™ Electrophoresis System User Guide at [thermofisher.com](#) or contact Technical Support.

i Limited product warranty and licensing information

Disclaimer: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288




©2019 Thermo Fisher Scientific Inc. All rights reserved. Triton is a Trademark of The Dow Chemical Company. Tween is a registered trademark of Croda International PLC. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

E-Base™ DNA electrophoresis protocol


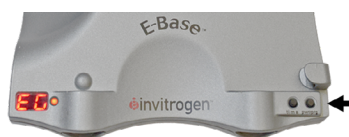
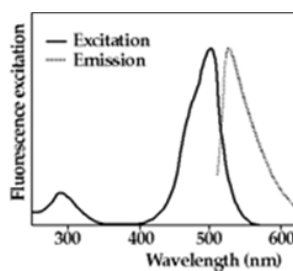


Important guidelines

- Use 10–200 ng DNA per band for samples with one unique band or up to 500 ng per lane for samples with multiple bands.
- Dilute samples with high salt concentrations (>50 mM NaCl, >100 mM KCl, >10 mM acetate ions, >10 mM EDTA) 2- to 5-fold in deionized water, TE, or 1X E-Gel™ Sample Loading Buffer, in a final volume of 15 µL (48-well gels) or 20 µL (96-well gels).
- Load E-Gel™ agarose gels within 30 minutes after opening the pouch; run gels within 1–3 minutes after loading samples.

Step		Action
1–5 min	1 	Prepare samples <p>Prepare DNA samples in deionized water OR 1X E-Gel™ Sample Loading Buffer.</p> <ul style="list-style-type: none"> ▪ For optimal separation use 20–100 ng of DNA per band for samples with one unique band or up to 500 ng per lane for samples with multiple bands. ▪ The total sample volume for 48-well gels is 15 µL. ▪ The total sample volume for 96-well gels is 20 µL.
5–10 min	2 	Prepare gel cassette <ol style="list-style-type: none"> Plug the Mother E-Base™ Device into an electrical outlet. Remove the gel from the package and gently remove the comb(s) from the E-Gel™ cassette. Insert the cassette into the E-Base™ Device, starting from the right edge. When properly inserted, the device indicates its initialized status with a steady red light. <p>Note: The protocol type on the display shows EG for E-Gel™ DNA cassettes, and EP for E PAGE™ cassettes.</p>
	3 	Load samples <p>Load samples with a multichannel pipettor.</p> <p>Load a volume of 15 µL in each well for 48-well gels. Load a volume of 20 µL in each well for 96-well gels.</p> <ol style="list-style-type: none"> Load prepared samples into sample wells. Keep all sample volumes uniform. Load prepared E-Gel™ DNA ladder into marker wells. Load 1X E-Gel™ Sample Loading Buffer or deionized water in all empty wells. The buffer for empty wells should have a similar salt concentration to adjacent sample wells.

E-Base™ DNA electrophoresis protocol




Step		Action						
12–30 min	<div>4</div> 	<div>Run the gel</div> <div><div><div>Gel type</div><div>Recommended run time</div><div>Maximum run time</div></div><table><tr><td>E-Gel™ 48 agarose gel</td><td>20 min</td><td>25 min</td></tr><tr><td>E-Gel™ 96 agarose gel</td><td>12 min</td><td>17 min</td></tr></table></div> <div>a. Select the EG program (default run time 12 minutes) for running E-Gel™ cassettes by pressing and releasing the “pwr/prg”.</div> <div>b. Select the recommended run time for a specific gel type by pressing and releasing the time button, then press and hold the time button to increase the time. Release time button when the desired run time for the gel is reached.</div> <div>c. Start the run by pressing and releasing the “pwr/prg” button. The red indicator light will change to green.</div>	E-Gel™ 48 agarose gel	20 min	25 min	E-Gel™ 96 agarose gel	12 min	17 min
	E-Gel™ 48 agarose gel	20 min	25 min					
E-Gel™ 96 agarose gel	12 min	17 min						
<div>5</div> 	<div>End the run</div> <div>a. A flashing red indicator light and rapid beeping indicates the end of the run. Press and release “pwr/prg” to stop the device.</div> <div>b. For better detection sensitivity, allow the gel to cool down for 10 minutes after the end of the run.</div>							
1–2 min	<div>6</div> 	<div>Analyze the gel</div> <div>a. Visualize the with a DNA imager using blue-light transillumination (e.g., with the E-Gel™ Imager System with Blue Light Base).</div> <div>▪ SYBR Safe™ DNA gel stain has an excitation maxima at 280 and 502 nm, and an emission maximum at 530 nm when bound to nucleic acid.</div> <div>▪ Use the E-Editor™ 2.0 software available at thermofisher.com/egel to analyze 96-well format digital images.</div>						

E-Base™ protein electrophoresis protocol



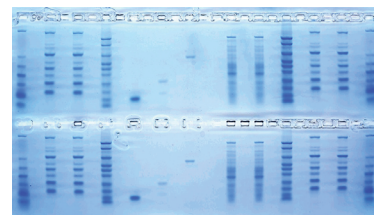


Important guidelines

- E-PAGE™ Gels contain SDS and are designed for electrophoresis under denaturing conditions.
- Dilute samples with high salt or detergent concentrations to prevent loss of resolution. ⓘ See **Table of recommended final concentrations**.
- For optimal separation use up to 20 µg of protein per well. Limit the protein (or lipid) amount in the sample to 2 µg/µL in the final sample volume for a proper LDS to protein ratio.

Step		Action															
1–5 min	<div>1</div> <div></div> <div>Prepare samples</div>	<div>a. Prepare protein samples in a total volume of 10 µL according to the following table. Scale volumes according the to the required sample volume.</div> <div><div><div>▪ The total sample volume for 48-well gels is 10 µL.</div><div>▪ The total sample volume for 96-well gels is 15 µL.</div></div><table><tr><th>Reagent</th><th>Reduced</th><th>Non-reduced</th></tr><tr><td>Protein sample</td><td>x µL</td><td>x µL</td></tr><tr><td>4X E-PAGE™ Loading Buffer 1</td><td>2.5 µL</td><td>2.5 µL</td></tr><tr><td>10X NuPAGE™ Sample Reducing Agent</td><td>1 µL</td><td>—</td></tr><tr><td>Deionized water</td><td>to 10 µL</td><td>to 10 µL</td></tr></table></div> <div>b. Incubate the samples at 70°C for 10 minutes.</div>	Reagent	Reduced	Non-reduced	Protein sample	x µL	x µL	4X E-PAGE™ Loading Buffer 1	2.5 µL	2.5 µL	10X NuPAGE™ Sample Reducing Agent	1 µL	—	Deionized water	to 10 µL	to 10 µL
	Reagent	Reduced	Non-reduced														
Protein sample	x µL	x µL															
4X E-PAGE™ Loading Buffer 1	2.5 µL	2.5 µL															
10X NuPAGE™ Sample Reducing Agent	1 µL	—															
Deionized water	to 10 µL	to 10 µL															
5–10 min	<div>2</div> <div></div> <div>Prepare gel cassette</div>	<div>a. Remove the gel from the package and gently remove the combs from the E-PAGE™ cassette.</div> <div>b. Insert the cassette into the E-Base™ Device, starting from the right edge. When properly inserted, the device indicates its initialized status with a steady red light.</div> <div>Note:The protocol type on the display shows EG for E-Gel™ DNA cassettes, and EP for E-PAGE™ cassettes.</div>															
	<div>3</div> <div></div> <div>Load samples</div>	<div>Load samples with a multichannel pipettor.</div> <div>Load a volume of 10 µL in each well for 48-well gels.</div> <div>Load a volume of 15 µL in each well for 96-well gels.</div> <div>a. Load 5–10 µL of deionized water into all wells prior to adding samples or standards.</div> <div>b. Load prepared samples into sample wells. Keep all sample volumes uniform.</div> <div>c. Load prepared E-PAGE™ standard into marker wells.</div> <div>d. Load deionized water in all empty wells.</div>															

E-Base™ protein electrophoresis protocol

Step		Action									
14–30 min	<div>4</div> <div></div> <div>Run the gel</div>	<div>a. Select the EP program (default run time 14 minutes) for running E-PAGE™ cassettes by pressing and releasing the “pwr/prg”.</div> <div>b. Select the recommended run time for a specific gel type by pressing and releasing the time button, then press and hold the time button to increase the time. Release time button when the desired run time for the gel is reached.</div> <table><thead><tr><th>Gel type</th><th>Recommended run time</th><th>Maximum run time</th></tr></thead><tbody><tr><td>E-PAGE™ 48 gel</td><td>25 min</td><td>30 min</td></tr><tr><td>E-PAGE™ 96 gel</td><td>14 min</td><td>25 min</td></tr></tbody></table> <div>c. Start the run by pressing and releasing the “pwr/prg” button. The red indicator light will change to green.</div>	Gel type	Recommended run time	Maximum run time	E-PAGE™ 48 gel	25 min	30 min	E-PAGE™ 96 gel	14 min	25 min
	Gel type	Recommended run time	Maximum run time								
E-PAGE™ 48 gel	25 min	30 min									
E-PAGE™ 96 gel	14 min	25 min									
	<div>5</div> <div></div> <div>End the run</div>	<div>a. A flashing red indicator light and rapid beeping indicates the end of the run. Press and release “pwr/prg” to stop the device.</div> <div>b. For better detection sensitivity, allow the gel to cool down for 10 minutes after the end of the run.</div>									
1–2 hr	<div>6</div> <div></div> <div>Stain the gel</div>	<div>a. Open the gel cassette.</div> <div>b. Visualize the protein by staining the gel using any of the following techniques (See the E-PAGE™ Technical Guide for details on staining and imaging).</div> <div><div>▪ SYPRO™ Ruby Protein Gel Stain protocol</div><div>▪ Coomassie R-250 protocol</div><div>▪ SimplyBlue™ SafeStain protocol</div><div>▪ SilverQuest™ Silver Stain protocol</div><div>▪ SilverXpress™ Silver Stain protocol</div></div>									

Gel selection guide

Application	Product	Gel %	Sample wells	In-gel stain ^[1]	Amount	Cat. No.
DNA sample analysis	E-Gel™ 48 Agarose Gels, 1%	1%	48 + 4 ladder lanes	SYBR Safe™	8 gels	G820801
					4 x 8 gels	G820841
	E-Gel™ 48 Agarose Gels, 2%	2%	48 + 4 ladder lanes	SYBR Safe™	8 gels	G820802
					4 x 8 gels	G820842
	E-Gel™ 96 Agarose Gels, 1%	1%	96 + 8 ladder lanes	SYBR Safe™	8 gels	G720801
					4 x 8 gels	G720841
Protein sample analysis	E-Gel™ 96 Agarose Gels, 2%	2%	96 + 8 ladder lanes	SYBR Safe™	8 gels	G720802
					4 x 8 gels	G720842
Protein sample analysis	E-PAGE™ 8% Protein Gels, 48-well	8%	48 + 4 ladder lanes	—	8 gels	EP4808
	E-PAGE™ 6% Protein Gels, 96-well	6%	96 + 8 ladder lanes	—	8 gels	EP9606

[1] For other stain options visit thermofisher.com/egel.

Ladder selection guide

Product	Recommended DNA ladder		
	E-Gel™ 96 High Range DNA Ladder (Cat. No. 12352019)	E-Gel™ 50 bp DNA Ladder (Cat. No. 10488099)	E-Gel™ Low Range Quantitative DNA Ladder (Cat. No. 12373031)
E-Gel™ 48 Agarose Gels, 1%	✓	—	—
E-Gel™ 48 Agarose Gels, 2%	—	✓	—
E-Gel™ 96 Agarose Gels, 1%	✓	—	—
E-Gel™ 96 Agarose Gels, 2%	—	—	✓

For more ladder options visit thermofisher.com/egelladders.