

NativePAGE™ Bis-Tris Gels

	Package Contents	<table border="0"> <thead> <tr> <th>Product</th> <th>Quantity</th> </tr> </thead> <tbody> <tr> <td>3–12% Bis-Tris Gels</td> <td>Box of 10 gels</td> </tr> <tr> <td>4–16% Bis-Tris Gels</td> <td>Box of 10 gels</td> </tr> </tbody> </table>	Product	Quantity	3–12% Bis-Tris Gels	Box of 10 gels	4–16% Bis-Tris Gels	Box of 10 gels
Product	Quantity							
3–12% Bis-Tris Gels	Box of 10 gels							
4–16% Bis-Tris Gels	Box of 10 gels							
	Storage Conditions	<ul style="list-style-type: none"> Store at 2–8°C for a 1-year shelf life. Do not freeze. 						
	Required Materials	<ul style="list-style-type: none"> Protein sample and standard NativePAGE™ Running Buffer Kit NativePAGE™ Sample Prep Kit Novex® Power Supply Adapters (Cat. no. ZA10001) if not using a Life Technologies™ power supply XCell SureLock® Mini-Cell gel running tank 						
	Timing	<p>Run Time: 90–115 minutes for the 3–12% gel 105–120 minutes for the 4–16% gel</p> <p>Voltage: 150 V constant</p>						
	Selection Guide	<p>Specialized Protein Gels Go online to view related products.</p>						
	Product Description	<p>NativePAGE® Bis-Tris Gels provide resolution for a wide range of proteins (15–10,000 kDa) under non-denaturing conditions.</p> <p>NativePAGE® Bis-Tris Gels are available in the following variations:</p> <ul style="list-style-type: none"> Polyacrylamide percentages: 3–12% and 4–16% Well formats: 10 and 15 wells Thickness: 1.0 mm 						
	Important Guidelines	<ul style="list-style-type: none"> This system is designed for use in the XCell SureLock® Mini-Cell gel running tank. Load samples into sample wells filled with 1X NativePAGE™ Dark Blue Cathode Buffer prior to filling the cathode chamber to better visualize the sample wells. For western blotting or two-dimensional (2D) electrophoresis applications, replace the Dark Blue Cathode Buffer with Light Blue Cathode Buffer during electrophoresis to improve protein transfer. Do not use SDS-PAGE samples for native gel electrophoresis. 						
	Online Resources	<p>Visit our product page for additional information and protocols. For support, visit www.lifetechnologies.com/support.</p>						



Protocol Outline

- Prepare samples, buffers, and gels.
- Assemble the gel apparatus.
- Load buffer, samples, and standards.
- Perform electrophoresis.

Electrophoresis Protocol

- See page 2 to view a procedure for preparing and running your electrophoresis experiment.

Choosing the Right Gel Type for Your Application

- Review the table in the pop-up to determine the best gel type for your experiment.

Choosing the Right Gel Percentage and Buffer

- Refer to the migration charts in the pop-up to find the gel best suited for your application. As a general rule, your proteins of interest should migrate through ~70% of the length of the gel for the best resolution. When protein molecular weights are wide ranging or unknown, gradient gels are usually the best choice.

Choosing a Well Format and Gel Thickness

- We offer NativePAGE® polyacrylamide gels in a choice of two well formats (10 or 15 wells) and one thickness (1.0 mm). When loading large samples, a gel with fewer wells is more appropriate.

Choosing a Protein Standard for your Application

Choose a Life Technologies™ standard based on your experiment:

- Unstained:** NativeMark™ Unstained Protein Standard
- Western:** NativeMark™ Unstained Protein Standard

For all other specialty standards, please view further information [here](#).

Limited Product Warranty and Disclaimer Details

NativePAGE™ Bis-Tris Mini Gel Electrophoresis Protocol

Follow the procedure below to prepare for and perform native gel electrophoresis using NativePAGE™ Bis-Tris Mini Gels.

Timeline	Steps	Procedure Details															
1	Prepare samples	<table border="1"> <thead> <tr> <th>Components</th> <th>Sample with Detergent</th> <th>Detergent-free Sample</th> </tr> </thead> <tbody> <tr> <td>Sample</td> <td>x μL</td> <td>x μL</td> </tr> <tr> <td>NativePAGE® Sample Buffer (4X)</td> <td>2.5 μL</td> <td>2.5 μL</td> </tr> <tr> <td>NativePAGE® 5% G-250 Sample Additive</td> <td>0.25–1 μL*</td> <td>optional</td> </tr> <tr> <td>Deionized Water</td> <td>to 10 μL</td> <td>to 10 μL**</td> </tr> </tbody> </table> <p>* Ensure that the final G-250 concentration is $\frac{1}{4}$th the detergent concentration. ** For additional sample preparation methods, follow instructions in the NativePAGE™ Sample Prep Kit. Do not heat samples for native gel electrophoresis. Prepare 1X Sample Buffer for dilutions of samples, if needed.</p>	Components	Sample with Detergent	Detergent-free Sample	Sample	x μ L	x μ L	NativePAGE® Sample Buffer (4X)	2.5 μ L	2.5 μ L	NativePAGE® 5% G-250 Sample Additive	0.25–1 μ L*	optional	Deionized Water	to 10 μ L	to 10 μ L**
Components	Sample with Detergent	Detergent-free Sample															
Sample	x μ L	x μ L															
NativePAGE® Sample Buffer (4X)	2.5 μ L	2.5 μ L															
NativePAGE® 5% G-250 Sample Additive	0.25–1 μ L*	optional															
Deionized Water	to 10 μ L	to 10 μ L**															
2	Prepare buffers	<p>1X NativePAGE™ Anode Buffer: Add 50 mL of 20X NativePAGE™ Running Buffer to 950 mL of deionized water.</p> <p>1X NativePAGE™ Dark Blue Cathode Buffer: Add 50 mL 20X NativePAGE™ Running Buffer and 50 mL 20X NativePAGE™ Cathode Additive to 900 mL deionized water.</p> <p>1X NativePAGE™ Light Blue Cathode Buffer: Add 50 mL 20X NativePAGE™ Running Buffer and 5 mL 20X NativePAGE™ Cathode Additive to 945 mL deionized water.</p>															
3	Prepare gels	<ol style="list-style-type: none"> Remove the comb, and rinse the gel wells three times with 1X NativePAGE™ Dark Blue Cathode Buffer. Remove the white tape near the bottom of the gel cassettes. Place the gels in the XCell SureLock® Mini-Cell gel running tank. Fill the gel wells with 1X NativePAGE™ Dark Blue Cathode Buffer. 															
4	Load samples and standards	<p>Load samples into sample wells filled with 1X NativePAGE™ Dark Blue Cathode Buffer prior to filling the cathode chamber to better visualize the sample wells.</p> <p>Load an appropriate volume and protein mass of samples on the gel. Then, load your standards.</p>															
5	Load buffers	<p>Fill the Upper Cathode Buffer Chamber with 200 mL 1X NativePAGE™ Dark Blue Cathode Buffer, and fill the Lower Anode Buffer Chamber with 550 mL NativePAGE™ Anode Buffer.</p>															
6	Run	<p>Note: If you are not using a Life Technologies™ power supply, install the Novex® Power Supply Adapters (Catalog number ZA10001).</p> <p>For western blotting or 2D electrophoresis applications, replace the Dark Blue Cathode Buffer with Light Blue Cathode Buffer after the dye front has migrated about one third of the way through the gel.</p> <p>When using the 3–12% gel, run for 90–115 minutes at 150 V constant. When using the 4–16% gel, run for 105–120 minutes at 150 V constant.</p>															