



Contents and storage

Gel type	Amount	Storage
NuPAGE™ Bis-Tris Gels	Box of 2 or 10 gels	Store at 4–25°C for up to 1 year. Do not freeze.



Product description

NuPAGE™ Bis-Tris Gels are precast polyacrylamide gels designed for optimal separation and resolution of small- to medium-sized proteins (1.5–300 kDa) under denaturing gel electrophoresis conditions.

NuPAGE™ Bis-Tris Mini Gels are available with the following specifications:

- **Polyacrylamide percentage:** 8%, 10%, 12%, and 4–12%
- **Well format:** 1, 9, 10, 12, 15, 17, IPG, and 2D wells
- **Thickness:** 1.0 mm and 1.5 mm



Required materials

- Protein sample and protein ladder
- NuPAGE™ MES or MOPS SDS Running Buffer (20X)
- NuPAGE™ LDS Sample Buffer (4X)
- NuPAGE™ Sample Reducing Agent (10X) (for reduced samples)
- NuPAGE™ Antioxidant (for reduced samples)
- Novex™ Power Supply Adapters (Cat. No. ZA10001) if not using a Thermo Fisher Scientific™ power supply
- Mini Gel Tank (Cat. No. A25977) or XCell SureLock™ Mini-Cell (Cat. No. EI0001)



Online resources

- Visit thermofisher.com/proteingels for additional information and protocols.
- For support, visit thermofisher.com/support.

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

Choosing a well format

Thicker 1.5 mm gels with fewer wells are recommended for large samples (>30 µL). Thinner 1 mm gels are recommended for blotting because of better protein transfer.

Well type	Maximum loading volume ^[1]		Maximum protein load
	1 mm thickness	1.5 mm thickness	
1-well	700 µL	—	12 µg/band
IPG-well	7-cm IPG strip	—	—
2D-well	400 µL	600 µL	12 µg/band
9-well	28 µL	—	0.5 µg/band
10-well	25 µL	37 µL	0.5 µg/band
12-well	20 µL	—	0.5 µg/band
15-well	15 µL	25 µL	0.5 µg/band
17-well	15 µL	—	0.5 µg/band

[1] Not every format is available for every gel type.

Choosing a protein ladder for your application

Type	Marker	Cat. No.
Pre-Stained	PageRuler™ Prestained Protein Ladder	26616
	PageRuler™ Plus Prestained Protein Ladder	26619
Unstained	PageRuler™ Unstained Protein Ladder	26614
	PageRuler™ Unstained Broad Range Protein Ladder	26630
Western blot	iBright™ Prestained Protein Ladder	LC5615
	MagicMark™ XP Western Protein Standard	LC5602

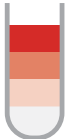

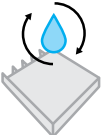

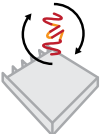

Go to thermofisher.com/proteinladders for more information on protein ladders.

Choosing buffers for your application

Buffer	Application	Cat. No.
NuPAGE™ MOPS SDS Running Buffer	Resolve mid-size proteins	NP0001
NuPAGE™ MES SDS Running Buffer	Resolve small molecular weight proteins	NP0002
NuPAGE™ Transfer Buffer	Wet transfer	NP0006

Limited product warranty and licensing information

Perform denaturing protein gel electrophoresis using NuPAGE™ Bis-Tris Mini Gels

Step		Action																		
<div data-bbox="69 363 107 395" style="border: 1px solid black; padding: 2px; display: inline-block;">1</div> 	<p>Prepare samples</p>	<p>Prepare 1X Sample Buffer for dilutions of samples if needed. Volumes are provided for a 10-μL sample size. Scale volumes proportionally for larger sample sizes.</p> <table border="1" data-bbox="672 245 1937 512"> <thead> <tr> <th data-bbox="672 245 1245 288">Components</th> <th data-bbox="1245 245 1592 288">Reduced sample</th> <th data-bbox="1592 245 1937 288">Non-reduced Sample</th> </tr> </thead> <tbody> <tr> <td data-bbox="672 288 1245 331">Sample</td> <td data-bbox="1245 288 1592 331">x μL</td> <td data-bbox="1592 288 1937 331">x μL</td> </tr> <tr> <td data-bbox="672 331 1245 375">NuPAGE™ LDS Sample Buffer (4X)</td> <td data-bbox="1245 331 1592 375">2.5 μL</td> <td data-bbox="1592 331 1937 375">2.5 μL</td> </tr> <tr> <td data-bbox="672 375 1245 418">NuPAGE™ Reducing Agent (10X)</td> <td data-bbox="1245 375 1592 418">1 μL</td> <td data-bbox="1592 375 1937 418">—</td> </tr> <tr> <td data-bbox="672 418 1245 461">Deionized Water</td> <td data-bbox="1245 418 1592 461">to 6.5 μL</td> <td data-bbox="1592 418 1937 461">to 7.5 μL</td> </tr> <tr> <td data-bbox="672 461 1245 512">Total Volume</td> <td data-bbox="1245 461 1592 512">10 μL ^[1]</td> <td data-bbox="1592 461 1937 512">10 μL ^[1]</td> </tr> </tbody> </table> <p>[1] See "Choosing a well format" for recommended loading volumes.</p> <p>Heat samples at 70°C for 10 minutes.</p>	Components	Reduced sample	Non-reduced Sample	Sample	x μ L	x μ L	NuPAGE™ LDS Sample Buffer (4X)	2.5 μ L	2.5 μ L	NuPAGE™ Reducing Agent (10X)	1 μ L	—	Deionized Water	to 6.5 μ L	to 7.5 μ L	Total Volume	10 μ L ^[1]	10 μ L ^[1]
Components	Reduced sample	Non-reduced Sample																		
Sample	x μ L	x μ L																		
NuPAGE™ LDS Sample Buffer (4X)	2.5 μ L	2.5 μ L																		
NuPAGE™ Reducing Agent (10X)	1 μ L	—																		
Deionized Water	to 6.5 μ L	to 7.5 μ L																		
Total Volume	10 μ L ^[1]	10 μ L ^[1]																		
<div data-bbox="69 671 107 703" style="border: 1px solid black; padding: 2px; display: inline-block;">2</div> 	<p>Prepare buffers</p>	<p>Add 50 mL of 20X NuPAGE™ MES or MOPS SDS Running Buffer to 950 mL of deionized water to prepare 1X SDS Running Buffer.</p> <p>For reduced samples, add 1 mL of NuPAGE™ Antioxidant to 400 mL 1X SDS Running Buffer.</p>																		
<div data-bbox="69 852 107 884" style="border: 1px solid black; padding: 2px; display: inline-block;">3</div> 	<p>Prepare gel</p>	<ol style="list-style-type: none"> Remove the comb, and rinse the gel wells three times using 1X Running Buffer. Remove the white tape near the bottom of the gel cassettes. Place the gels in the mini gel tank. 																		
<div data-bbox="69 1023 107 1054" style="border: 1px solid black; padding: 2px; display: inline-block;">4</div> 	<p>Load buffers</p>	<p>Fill the chambers with the appropriate 1X running buffer.</p> <p>Mini Tank: Add 400 mL of buffer to each chamber.</p> <p>XCell SureLock™ Mini-Cell: Add 600 mL of buffer to the lower chamber, and 200 mL to the upper chamber (for reduced samples, use running buffer with antioxidant in the upper chamber).</p>																		
<div data-bbox="69 1208 107 1240" style="border: 1px solid black; padding: 2px; display: inline-block;">5</div> 	<p>Load samples and ladders</p>	<ol style="list-style-type: none"> Load the appropriate volume of your samples in the appropriate wells. Load your protein ladder in the appropriate well. 																		
<div data-bbox="69 1374 107 1406" style="border: 1px solid black; padding: 2px; display: inline-block;">6</div> 	<p>Run the gel</p>	<p>Optimal run times vary depending on gel percentage and power supply used for electrophoresis.</p> <p>If using MES Running Buffer, run for 35 minutes at 200 V constant.</p> <p>If using MOPS Running Buffer, run for 50 minutes at 200 V constant.</p> <p>Note: If you are not using a Thermo Fisher Scientific™ power supply, install Novex™ Power Supply Adapters</p>																		

Buffer formulation

The following recipes are provided to allow preparation of buffers from scratch.

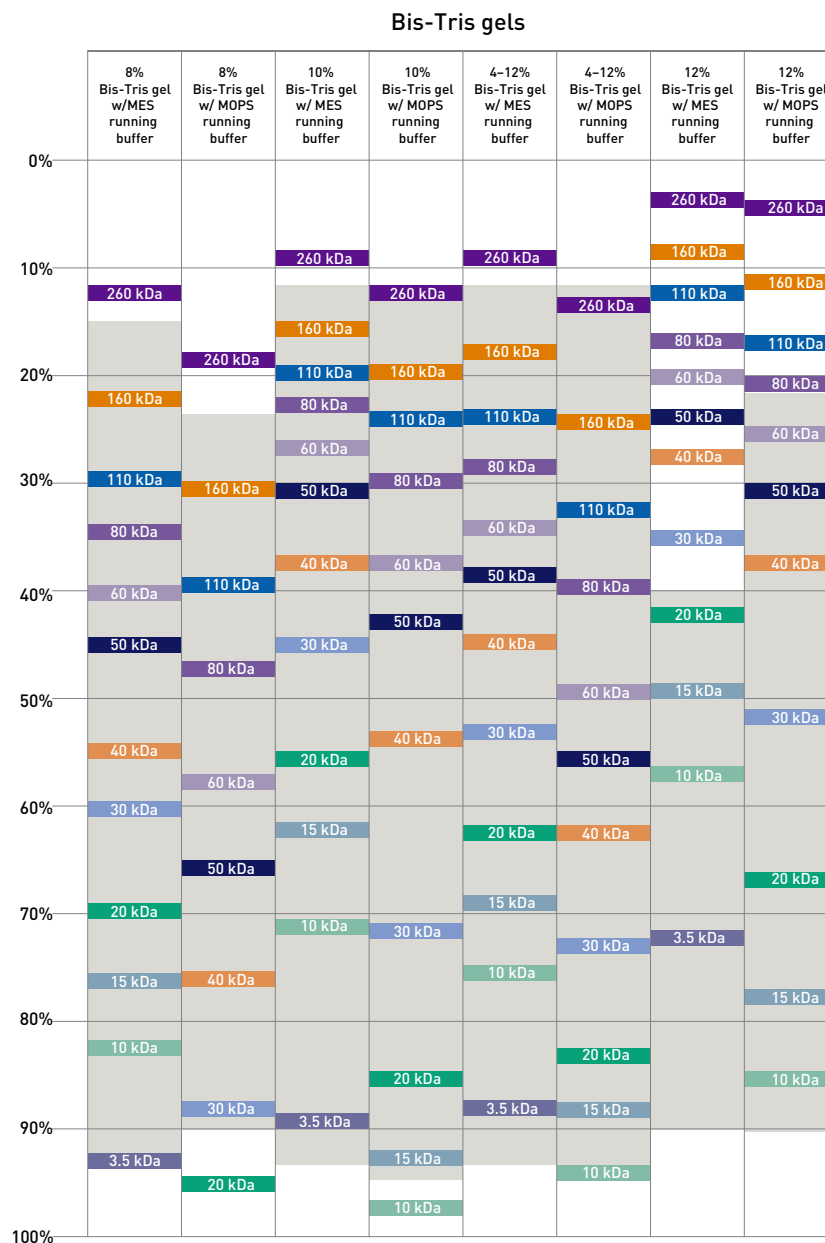
The pH listed for each buffer is for the 1X solution. **Do not use acid or base to adjust the pH.** Buffers are stable for 6 months when stored at 4°C.

Prepare 500 mL of 20X MES SDS Running Buffer	Prepare 500 mL of 20X MOPS SDS Running Buffer																				
50 mM MES, 50 mM Tris Base, 0.1% SDS, 1 mM EDTA, pH 7.3	50 mM MOPS, 50 mM Tris Base, 0.1% SDS, 1 mM EDTA, pH 7.7																				
1. Dissolve the following reagents in 400 mL ultrapure water.	1. Dissolve the following reagents in 400 mL ultrapure water.																				
<table border="1"> <thead> <tr> <th>Reagent</th> <th>Amount</th> </tr> </thead> <tbody> <tr> <td>MES</td> <td>97.6 g</td> </tr> <tr> <td>Tris Base</td> <td>60.6 g</td> </tr> <tr> <td>SDS</td> <td>10.0 g</td> </tr> <tr> <td>EDTA</td> <td>3.0 g</td> </tr> </tbody> </table>	Reagent	Amount	MES	97.6 g	Tris Base	60.6 g	SDS	10.0 g	EDTA	3.0 g	<table border="1"> <thead> <tr> <th>Reagent</th> <th>Amount</th> </tr> </thead> <tbody> <tr> <td>MOPS</td> <td>104.6 g</td> </tr> <tr> <td>Tris Base</td> <td>60.6 g</td> </tr> <tr> <td>SDS</td> <td>10.0 g</td> </tr> <tr> <td>EDTA</td> <td>3.0 g</td> </tr> </tbody> </table>	Reagent	Amount	MOPS	104.6 g	Tris Base	60.6 g	SDS	10.0 g	EDTA	3.0 g
Reagent	Amount																				
MES	97.6 g																				
Tris Base	60.6 g																				
SDS	10.0 g																				
EDTA	3.0 g																				
Reagent	Amount																				
MOPS	104.6 g																				
Tris Base	60.6 g																				
SDS	10.0 g																				
EDTA	3.0 g																				
2. Mix well and adjust the volume to 500 mL with ultrapure water.	2. Mix well and adjust the volume to 500 mL with ultrapure water.																				
3. Before electrophoresis, dilute buffer to 1X with water.	3. Before electrophoresis, dilute buffer to 1X with water.																				

Prepare 125 mL of 20X Bis-Tris Transfer Buffer								
25 mM Bicine, 25 mM Bis-Tris (free base), 1 mM EDTA, pH 7.2								
1. Dissolve the following reagents in 100 mL ultrapure water.								
<table border="1"> <thead> <tr> <th>Reagent</th> <th>Amount</th> </tr> </thead> <tbody> <tr> <td>Bicine</td> <td>10.2 g</td> </tr> <tr> <td>Bis-Tris (free base)</td> <td>13.1 g</td> </tr> <tr> <td>EDTA</td> <td>0.75 g</td> </tr> </tbody> </table>	Reagent	Amount	Bicine	10.2 g	Bis-Tris (free base)	13.1 g	EDTA	0.75 g
Reagent	Amount							
Bicine	10.2 g							
Bis-Tris (free base)	13.1 g							
EDTA	0.75 g							
2. Mix well and adjust the volume to 125 mL with ultrapure water.								
3. Before western transfer, dilute buffer to 1X with water.								

Migration patterns of protein standards on NuPAGE™ Bis-Tris gels

Refer to the migration chart to find the gel best suited for your application. Your proteins of interest should migrate through ~70% of the length of the gel for the best resolution.



Choosing the right gel type for your application

Thermo Fisher Scientific protein gels					
Gel type	Gel % available	Separation range	Shelf life	Average run time	Applications
Bolt™ Bis-Tris Plus	8%, 10%, 12%, 4–12%	6 to 400 kDa	up to 16 months	22–45 min	Best choice for separation of small- to medium-sized proteins. Neutral pH environment minimizes protein modifications. Wedge well design can accommodate large sample volumes. Ideal for Western blot transfer and analysis, and all techniques in which protein integrity is crucial.
NuPAGE™ Bis-Tris	8%, 10%, 12%, 4–12%	1.5 to 300 kDa	12 months	35–50 min	Separation of small- to medium-sized proteins. Neutral pH environment minimizes protein modifications.
NuPAGE™ Tris-Acetate	7%, 3–8%	36 to 500 kDa	6 months	60 min	Separation of larger proteins.
Novex™ Tricine	10%, 16%	2 to 200 kDa	1–2 months	90 min	Separation of small proteins and peptides.
NativePAGE™ Bis-Tris	3–12%, 4–16%	15 to 10,000 kDa	12 months	90–120 min	Separation of native proteins.
Novex™ Tris-Glycine	6%, 8%, 10%, 12%, 4–20%, 4–12%, 14%, 16%, 8–16%, 10–20%	8 to 250 kDa	up to 12 months	90 min	Separation of small- to medium-sized proteins using traditional Laemmli-style gels.
E-PAGE™	E-PAGE™ 96 6% E-PAGE™ 48 8%	10 to 220 kDa	6 months	14–25 min	High-throughput for recombinant production analysis and protein profiling.
IEF	pH 3–7, pH 3–10	pI 3.5 to pI 8.5	2 months	150 min	Separation of proteins based on isoelectric point.
Zymogram	10% (gelatin)	10 to 220 kDa	2 months	90 min	Separation of proteins based on size and activity.