

Precise™ Protein Gels

25200-25244

1472.9

Precise Protein Gel Selection Table*

Product Number	% Acrylamide	Running Buffer	Cassette Size	Wells	Well Volume	Separation Range (kDa)
25200	8	Tris-HEPES-SDS	100 × 85	10	50µL	205-45
25201	10	Tris-HEPES-SDS	× 4.5mm	10	50µL	205-24
25202	12	Tris-HEPES-SDS		10	50µL	205-14
25203	8-16	Tris-HEPES-SDS		10	50µL	205-14
25204	4-20	Tris-HEPES-SDS		10	50µL	205-6.5
25220	8	Tris-HEPES-SDS	100 × 85	12	30µL	205-45
25221	10	Tris-HEPES-SDS	× 4.5mm	12	30µL	205-24
25222	12	Tris-HEPES-SDS		12	30µL	205-14
25223	8-16	Tris-HEPES-SDS		12	30µL	205-14
25224	4-20	Tris-HEPES-SDS		12	30µL	205-6.5
25240	8	Tris-HEPES-SDS	100 × 85	15	25µL	205-45
25241	10	Tris-HEPES-SDS	× 4.5mm	15	25µL	205-24
25242	12	Tris-HEPES-SDS		15	25µL	205-14
25243	8-16	Tris-HEPES-SDS		15	25µL	205-14
25244	4-20	Tris-HEPES-SDS		15	25µL	205-6.5

*Choose a Precise Protein Gel equivalent to the gel that is used in the Laemmli system.

Important Note:

***Do not use Tris-glycine-SDS running buffer with Precise Protein Gels.
Use only Tris-HEPES-SDS running buffer as described in these instructions.***

Table of Contents

Introduction	2
Instructions for using Precise Protein Gels	3
A. Preparing the Gel Cassette and Gel Tank.....	3
B. SDS Sample Preparation.....	4
C. Sample Loading	4
D. Running Conditions	4
E. Removing a Gel from the Cassette.....	5
Staining and Drying Gels.....	5
Western Blotting Protocols for Precise Protein Gels	5
Buffer Recipes	6
Troubleshooting	7
Related Thermo Scientific Products	8

Introduction

SDS-PAGE (sodium dodecyl sulfate polyacrylamide gel electrophoresis) is used to determine the approximate molecular weights of proteins with reference to the mobility of standard proteins on the same gel. Gradient gels provide the appropriate gel pore size making molecular weight estimation more accurate by sharpening stained protein bands (Figure 1).

The most popular buffer system for SDS-PAGE is the Laemmli system [Nature 227, 680–686 (1970)]. This system, however, is limited by extended running times and gel instability from hydrolysis of polyacrylamide to acrylic acid in alkaline conditions. This chemical change alters gel conductivity, causing a constant change in the gel's migration pattern with time.

When the gel pH is neutral, hydrolysis does not occur. Thermo Scientific™ Precise™ Protein Gels are cast at pH 7, yielding a long shelf life and assured reproducibility of the migration pattern (Figure 2). The extra stability of Precise Protein Gels is combined with a new Tris-HEPES running buffer to give speed and excellent resolution with the same size ranges as the Laemmli system. Advantages of Precise Protein Gels are as follows:

- Sample wells reinforced with plastic eliminate damage when loading
- Sample well dividers do not deform or fall over
- Sample wells hold up to twice the volume of other gels
- Easy-to-open cassette
- Universally compatible cassette design
- 12-month shelf-life warranty from date of purchase
- 45-minute run time
- 90-minute transfer time

Please visit our web site or catalog for a complete list of electrophoresis reagents including MW markers, gel staining, Western blotting and products for sample preparation.

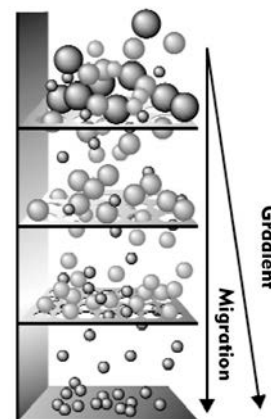


Figure 1. Gradient gel electrophoresis.

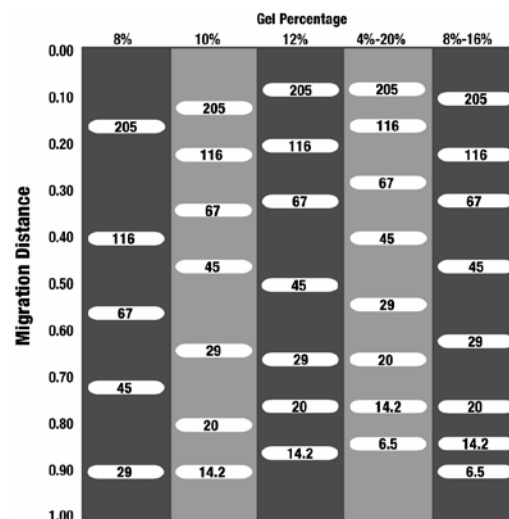


Figure 2. Protein migration table of the various gels.

Precise Protein Gels Specifications

Cassette Dimensions: 10cm × 8.5cm × 4.5mm

Gel Dimensions: 8cm × 5.8cm × 1mm

Storage Conditions: 4°C

Shelf Life: 12 months from date of purchase

Stacking Gel: 4%

Buffer System in Gel: Tris-HCl, pH 7

SDS: None

Required Running Buffer: Tris-HEPES-SDS

Recommended Sample Buffer: Tris-HCl-SDS

Compatible Gel Tanks

- Novex™ XCell I and II™
- Novex XCell II Surelock™
- Bio-Rad™ Mini-PROTEAN™ II and 3
- Hoefer™ Tall Mighty Small™ (SE 280)
- Hoefer Mighty Small II (SE 260/SE 250)
- IBI Universal Protein System
- Owl Road Runner, Penguin
- Owl Single Sided Vertical System

Instructions for using Precise Protein Gels

A. Preparing the Gel Cassette and Gel Tank

Important: Please see notes at the end of this section (Section A) concerning special instructions for using Bio-Rad Mini PROTEAN Cell and NOVEX Tanks.

1. Dissolve one packet of Thermo Scientific BupH Tris-HEPES-SDS Running Buffer (Product No. 28398) in 500mL of ultrapure water. This buffer volume of running buffer (500mL) is sufficient for one electrophoresis unit. (See Buffer Recipes on page 6 for instructions for preparing a 10X stock of the required Tris-HEPES-SDS Running Buffer.)

Note: Do not use Tris-glycine-SDS running buffer. This buffer formulation is not compatible with Precise Gels. Proteins will not migrate or resolve properly. Use only the required Tris-HEPES-SDS running buffer.

2. Remove Precise Protein Gel from the pouch and insert into the gel running apparatus. Refer to the apparatus manufacturer's instructions.
3. Add sufficient volume of Tris-HEPES-SDS running buffer into the inner tank of the gel running apparatus to cover the sample wells by 5-7mm.
4. Add the remaining volume of Tris-HEPES-SDS running buffer to the outer tank to ensure proper cooling. The buffer in the outer tank should be approximately level with the bottom of the sample wells.

Note: For best resolution, the buffer in the outer tank must reach the bottom of the sample wells to keep the gels cool.

5. Using a transfer (pasteur) pipette, rinse the sample wells thoroughly with Tris-HEPES-SDS running buffer to remove air bubbles and to displace any storage buffer. The gel may be pre-electrophoresed for 5-10 minutes.

Notes for using Bio-Rad Mini-PROTEAN Cell and NOVEX Tanks:

Using a Bio-Rad Mini-PROTEAN Cell

- To use a Bio-Rad Mini-PROTEAN Cell Apparatus, remove the gasket from the inner frame (Figure 3), turn it around so the flat side is facing outwards and re-insert into the inner frame.

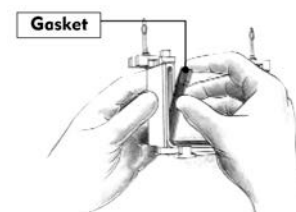


Figure 3. Removing the gasket from the cell.

Using a NOVEX Gel Running Apparatus

Note: A tight seal must be formed between the gel cassette and the gasket of the running frame to prevent leakage. In the NOVEX Tanks, adaptor plates must be used to form a tight seal. Below are instructions for adaptor plate placements when running only one gel and when running two gels.

Precise Protein Gels are 8.5cm high and are positioned lower in the Novex Tank than other brands of precast gels. However, the resolving portion of the gel is the same length as a Novex Gel.

Running One Precise Protein Gel in a NOVEX Tank

- Two adaptor plates are required when running just one gel.
- Place the gel onto the running frame.
- Place two adaptor plates onto the back of the running frame. (Ensure that the wells are facing away from the inner tank so that buffer does not leak.)
- Put the running frame into the tank.
- Using the wedge device, clamp the running frame and the gels into place (Figure 4). The running frame should be fixed tightly and the gels should not be able to move.

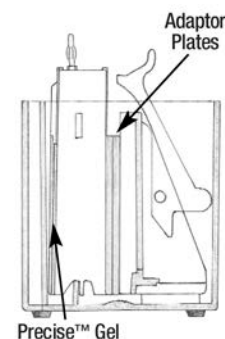


Figure 4. Adaptor plates placement when using one gel.

Running Two Precise Protein Gels in a NOVEX Tank

- One adaptor plate is required when running two gels.
- Place both gels onto the running frame.
- Place one adaptor plate behind the back gel. (Ensure that the cassettes are facing away from the inner tank so that the buffer does not leak.)
- Place the running frame into the tank.
- Using the wedge device, clamp the running frame and the gels into place (Figure 5). The running frame should be fixed tightly and the gels should not be able to move.

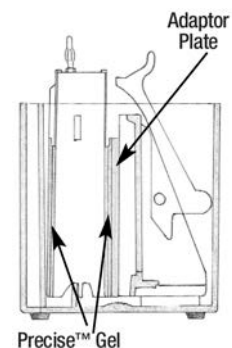


Figure 5. Adaptor plate placement when using two gels.

B. SDS Sample Preparation

Add 1 part Lane Marker Reducing or Non-Reducing 5X Sample Buffer (Product No. 39000 or 39001) to 4 parts sample. Alternatively, use the sample buffer recipe on page 6.

For solubilizing lyophilized samples, mix 100 μ L of Tris-HCl SDS Sample Buffer (1X) per mg of protein. Heat sample for 3-5 minutes at approximately 100°C. Clarify by centrifugation at 6000 rpm for 3 minutes and collect the supernatant.

Note: Tris-HEPES gels are compatible with Tris-Glycine SDS sample loading buffer (non-reducing and reducing). If the sample is thermally labile, add SDS to the sample as a 4% solution. Incubate sample for 1 hour at room temperature then add an equal volume of solution containing 0.05M HEPES, 1M sodium chloride, 10% glycerol and 0.05% Bromophenol blue. Dissolution may be helped by sonication.

C. Sample Loading

Note: These Precise Protein Gels are designed to open without any mechanical devices (e.g., keys, knives, etc.). However, this easy-to-open design requires that you verify the center pin at the top of the cassette is pinched tight to prevent leakage during sample loading and electrophoresis.

Apply 5-50 μ g (total protein) per sample well. Each sample well holds from 25-50 μ L. For a sample with a total protein concentration of 10mg/mL, apply 2-5 μ L per well. For best results, use pipette tips specifically designed for gel loading.

Caution: Inserting the pipette tip too far into the cassette may cause the cassette to separate.

Note: Optimal sample size must be established empirically. Overloading will cause smearing and distortion. Excessive loading of proteins with free carbohydrate may also result in band distortion or failure of the protein to penetrate into the gel (see Troubleshooting Section).

D. Running Conditions

Connect the gel rig leads to the power supply (Figure 6) and electrophorese according to Table 1.



Figure 6. Gel tank and power supply.

Table 1. Electrophoresis conditions for Precise Gels.

Voltage	Approximate Current		Run Time per Gel*
	Start	Finish	
100-120 V	100-130mA/gel	40-60mA/gel	~45 minutes

*Gel running time is dependent on the temperature in the laboratory. These run times are recommended at a laboratory temperature of 20°C.

E. Removing a Gel from the Cassette

1. Once the run is finished, remove the gel from the gel tank according to the manufacturer's instructions.
2. To open the cassette, insert a coin in one of the slots on the side and twist (Figure 7a).
3. Pull the top plate of the cassette away from the bottom plate (Figure 7b). The two halves will snap apart completely, exposing the gel.
4. Loosen the gel at the bottom with water and remove.



Figure 7a and 7b. Opening the cassette to expose the gel.

Staining and Drying Gels

All standard SDS staining procedures may be used with Precise Protein Gels.

Gels can be dried using standard drying techniques. When using commercially available gel drying reagents, follow the manufacturer's instructions.

Coomassie Staining

- To increase staining sensitivity and decrease staining time, wash gel three times for 5 minutes each in 200mL of water before staining. This wash will remove SDS from the gel.
- Commercially available stains, such as Thermo Scientific™ Imperial™ Protein Stain (Product No. 24615), Thermo Scientific™ GelCode™ Blue Stain (No. 24590), and homemade coomassie stains may be used. Best results are obtained with methanol concentrations of < 30%.

Silver Staining

- For best results, before staining, wash gels in ultrapure water for 10-15 minutes to remove SDS.
- Commercially available stains, such as Thermo Scientific™ Pierce™ Silver Stain Kit (Product No. 24612), Pierce Color Silver Stain Kit (Product No. 24597) and homemade silver stains may be used.

Western Blotting Protocols for Precise Protein Gels

Standard blotting procedures may be used with Precise Gels. Below are protocols for wet blotting and semi-dry blotting.

Wet Blotting Protocol

1. Cool the transfer buffer to 4°C.
2. Equilibrate the gels in transfer buffer for 5 minutes.
3. Soak filter papers (8 × 10cm) in transfer buffer.
4. Soak membrane(s) (8 × 10cm) in transfer buffer (PVDF membranes must be wetted in methanol first and then equilibrated in aqueous solution).
5. Soak the Scotch-Brite™ Pads in transfer buffer.
6. Assemble the transfer sandwich as follows:
 - Cathode (---)
 - Scotch-Brite Pad
 - 2x filter paper
 - Gel
 - Transfer Membrane
 - 2x filter paper
 - Scotch-Brite Pad
 - Anode (+++)

Note: Blotter should be firmly packed. If two membranes are to be blotted, repeat the above transfer sandwich. If only one gel is to be blotted, fill the space with more filter paper and another Scotch-Brite Pad.

7. Pour the transfer buffer through the sandwich and place it into the apparatus. Fill the apparatus with transfer buffer.
8. Transfer at 40V for 90 minutes (maintain buffer temperature at ~4°C).
9. Gently remove gel from sandwich and rinse with transfer buffer.
10. Use a cotton swab to remove any adhering gel from the membrane.

Semi-Dry Blotting Protocol

1. Cool the transfer buffer to 4°C.
2. Soak the filter paper, membrane and gel in Tris-Glycine Transfer Buffer (Product No. 28380) for 15 minutes.
3. Assemble the blotting sandwich in a semi-dry blotting apparatus as follows:
 - Anode (+++)
 - Filter paper
 - Membrane
 - Gel
 - Filter paper
 - Cathode (---)
4. Transfer the blot for 30 minutes at 20V.
5. Remove the gel from the sandwich and rinse with transfer buffer.
6. Use a cotton swab to remove any adhering gel from the membrane.

Staining Membranes

The Pierce Reversible Protein Stain Kits (Product No. 24580 for nitrocellulose; Product No. 24585 for PVDF) contain Reversible Protein Stain, Destain and Stain Eraser. Sensitivity is ~25ng of protein (~10 times the sensitivity of Ponceau S stain) and staining and destaining can be accomplished in 15 minutes.

Buffer Recipes

Use high-purity reagents and high-purity water when making buffers.

Tris-HEPES-SDS Running Buffer (10X)

Tris Base (MW = 121)	121g
HEPES (free acid MW = 238)	238g
SDS (MW = 288)	10g

Add ultrapure water to 1L

- Before use dilute 10-fold with water. The pH of the 1X buffer should be ~8.0; do not adjust pH.
- Final composition of the 1X buffer is 100mM Tris, 100mM HEPES, 0.1% (~3mM) SDS, pH 8.0.
- Do not use Tris-glycine-SDS running buffer with Precise Protein Gels.

Protein Transfer Buffer

Dissolve one BupH™ Tris-Glycine Transfer Buffer Pack (Product No. 28380) in 400mL ultrapure water

Add 100mL methanol (20%) and cool to 4°C

Note: The pH of the buffer should be ~8.0

Sample Buffer (2X)

10% (w/v) Sodium Dodecyl Sulfate (SDS) Electrophoresis Grade	4.0mL
Glycerol	2.0mL
0.1% (w/v) Bromophenol Blue	1.0mL
0.5M Tris•HCl, pH 6.8	2.5mL
2-β Mercaptoethanol or DTT*	2-5% v/v

Add ultrapure water to 10mL

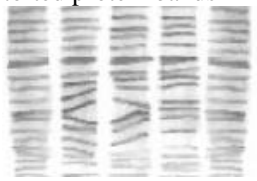
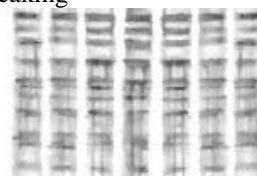


*Add if cleavage of disulfide bonds is desired.

Alternative Protein Transfer Buffer

Tris Base	3.00g
Bicine	4.08g
Ethanol or Methanol	100mL

Add ultrapure water to 1L

Troubleshooting

Problem	Cause	Solution
Distorted protein bands 	Air bubbles in the sample wells, or between gel and cassette, or at the bottom of the cassette	Use a transfer pipette to displace the air bubbles from the sample wells
	Sample contains appreciable carbohydrate	Remove the carbohydrate by enzymatic or chemical means
	Sample contains lipoproteins	Use a gel with a large pore size at top or try adding a non-ionic detergent
Streaking 	Poorly soluble or weakly charged particles (such as carbohydrates) in sample	Centrifuge samples
		Change pH of sample buffer
		Heat sample in the presence of SDS
Bands difficult to distinguish 	Incorrect gel selection, sample overloading and insufficient cooling buffer	Select a gel that separates in the desired molecular weight range
		Reduce sample size
		Increase buffer volume in the outer tank
		For proteins of similar molecular weight, a 2D separation may be required
Sample spreading across gel 	Excess salt in the sample	Reduce salt by dialysis or ultra-filtration
	Too much protein applied to the gel	Optimize the amount of protein applied to the gel
Protein denaturation and band inversion	Excessive heating	Start with chilled buffer (< 15°C)
Diffuse protein zones in the gel after staining	SDS present in the gel	Wash gel extensively (3 × 5 minutes) with ultrapure water and use 30% methanol to destain gel
	Protein bands are diffusing	Use 10% TCA to fix the proteins in the gel
Buffer front and proteins migrate only partly down the gel and are distorted	Incorrect running buffer used (e.g., Tris-glycine-SDS)	Use only Tris-HEPES-SDS Running Buffer (100mM Tris, 100mM HEPES, 1% SDS, pH 8.0)

Related Thermo Scientific Products

26616	PageRuler™ Prestained Protein Ladder, 10 to 170kDa
26619	PageRuler Plus Prestained Protein Ladder, 10 to 250kDa
26634	Spectra™ Multicolor Broad Range Protein Ladder, 10 to 260kDa
39000	Lane Marker Reducing Sample Buffer (5X), 5mL
77720	Bond-Breaker™ TCEP (Odorless reducing agent), 5mL
28398	BupH Tris-HEPES-SDS Running Buffer, 10 packs
28380	BupH Tris-Glycine Transfer Buffer, 40 packs
24615	Imperial™ Protein Stain, 1L
24594	GelCode Blue Safe Protein Stain, 1L
24590	GelCode Blue Stain Reagent, 500mL
24612	Pierce Silver Stain Kit
24597	Pierce Color Silver Stain Kit
24582	E-Zinc™ Reversible Stain Kit

Please see our web site or catalog for more information on our complete line of Western blotting products, such as the following:

- Nitrocellulose, PVDF and nylon transfer membranes
- Blocking buffers
- Labeled secondary antibodies
- X-ray film
- Chemiluminescent, chemifluorescent and colorimetric substrates
- Wash buffers and detergents
- Background Eliminator for film
- Western blot stripping buffer

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts (“Documentation”). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER’S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER’S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2013 Thermo Fisher Scientific Inc. All rights reserved. Bio-Rad, Mini-PROTEAN and Mini-Trans-Blot are trademarks of Bio-Rad Laboratories. Novex, Surelock & Xcell II are trademarks of Life Technologies Corp. Mighty Small is a trademark of Harvard Bioscience. Scotch-Brite is a trademark of 3M Company. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.