



critterion **XT**

Timeless Beauty.

With up to a 12-month shelf life, new Criterion™ XT gels give you beautiful results any time.



BIO-RAD

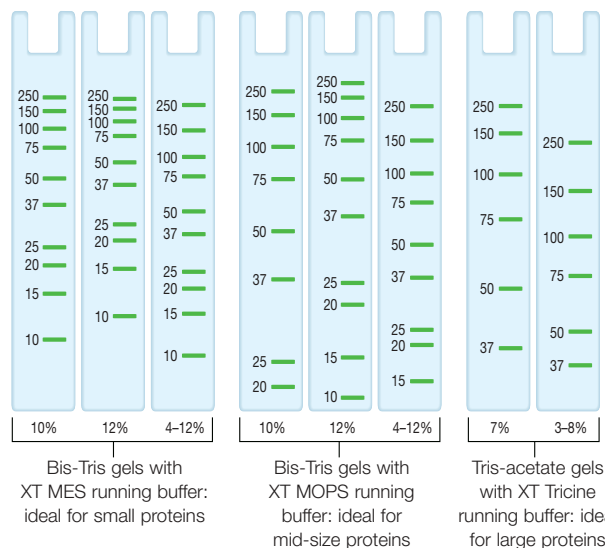
Criterion XT Precast Gels

With eXTended shelf life and room temperature storage!

Criterion XT gels are the highest-quality extended shelf-life gels available for protein electrophoresis with all the benefits of the Criterion system, including an integrated upper buffer chamber for ease-of-use, and gels that run up to 26 samples in an hour.

- Formulated at near-neutral pH to ensure:
 - Long shelf life (12 months for Bis-Tris, 8 months for Tris-acetate)
 - Better protein stability — Ideal for downstream applications such as protein sequencing and mass spectrometry
- Optimized sample and running buffers for sharp bands and minimal preparation time
- Choose from 3 buffer systems for flexibility in protein separations (see migration charts at right)

Visit criterion.bio-rad.com to schedule a demo.



Criterion XT Gels Separate More Samples: Demonstrated by Bio-Rad Precision Plus Protein™ standards separated on a Criterion XT 12% Bis-Tris gel (left) and the leading competitor's 12% Bis-Tris gel (right).

SDS-PAGE Gel to Criterion XT Gel Conversion Table

If you use ...	Then go with ...
10–20%, 16.5% peptide/Tricine gels	10%, 12%, 4–12% Bis-Tris gels with XT MES buffer
10% Tris-HCl gels	10% Bis-Tris gels with XT MOPS buffer
12% Tris-HCl gels	12% Bis-Tris gels with XT MOPS buffer
15%, 4–15%, 4–20% Tris-HCl gels	4–12% Bis-Tris gels with XT MOPS buffer
5% Tris-HCl gels	3–8% Tris-acetate gels with XT Tricine buffer
7.5% Tris-HCl gels	7% Tris-acetate gels with XT Tricine buffer

Ordering Information

Description	12+2 Well Comb* 45 µl Samples	18-Well Comb 30 µl Samples	26-Well Comb* 15 µl Samples	Prep+2 Well Comb 800 µl Samples	IPG+1 Well Comb 11 cm IPG Strip	IPG-Well Comb 11 cm IPG Strip
Criterion XT Bis-Tris Gels						
10% Resolving Gel	345-0111	345-0112	345-0113	345-0114	345-0115	345-0116
12% Resolving Gel	345-0117	345-0118	345-0119	345-0120	345-0121	345-0122
4–12% Resolving Gel	345-0123	345-0124	345-0125	345-0126	345-0127	345-0128
Criterion XT Tris-Acetate Gels						
3–8% Resolving Gel	345-0129	345-0130	345-0131	345-0132	345-0133	345-0134
7% Resolving Gel	345-0135	345-0136	345-0137	345-0138	345-0139	345-0140

* Multichannel pipet compatible.

Criterion Cell

165-6001 Criterion Cell, includes tank, lid with power cables, sample loading guides and instructions

Criterion XT Buffers and Reagents

161-0788	XT MOPS Running Buffer, 20x, 500 ml	161-0793	XT MOPS Buffer Kit, includes 500 ml 20x XT MOPS running buffer, 10 ml 4x sample buffer, 1 ml XT reducing agent
161-0789	XT MES Running Buffer, 20x, 500 ml		XT MES Buffer Kit, includes 500 ml 20x XT MES running buffer, 10 ml 4x sample buffer, 1 ml XT reducing agent
161-0790	XT Tricine Running Buffer, 20x, 500 ml	161-0796	XT Tricine Buffer Kit, includes 500 ml 20x XT Tricine running buffer, 10 ml 4x sample buffer, 1 ml XT reducing agent
161-0791	XT Sample Buffer, 4x, 10 ml		
161-0792	XT Reducing Agent, 1 ml	161-0797	

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191 Canada (905) 712-2771 Czech Republic + 420 2 41 43 05 32 China (86-21) 63052255 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 8126 672 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-62729877 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000