

New **Refresh** Kit Components (see pages 4-5) • **Free** PCR Kit Offer (see page 13)  
Up to **40%** Educator Savings • New **Free PCR Poster** Offer (see page 11)

# Biotechnology Explor<sup>™</sup>

CAPTIVATING SCIENCE EDUCATION



DNA > RNA > Protein > **Refresh!**

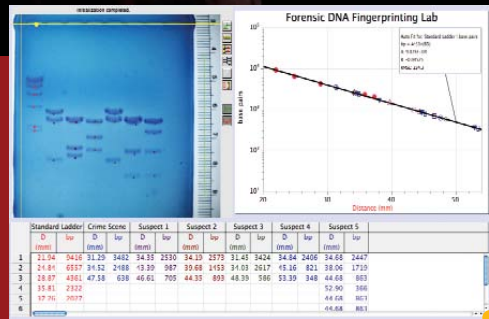
**BIO-RAD**

CURRICULA • EQUIPMENT • **KITS** • PROFESSIONAL DEVELOPMENT • TECHNICAL SUPPORT



# Collaboration

**Collaboratory.** At Bio-Rad our mission is to partner with you, our customers, so that we can deliver just what you want for your teaching laboratories. This year we are beginning the first in a series of initiatives to help you be green — by introducing biodegradable baggies for many of our products, reducing waste and providing purchasing flexibility at the same time. Our first step is the introduction of new packaging options for our classroom kits, helping you to reduce waste and maximize your space and budget for the key items you need. Read all about our new options on pages 4-5.



# laboratory



# laboratory

**Partnership.** The Biotechnology Explorer program was developed from a community of educators and researchers and we continue to learn from our colleagues. The core beliefs of the Biotechnology Explorer program are exemplified by Vernier Software & Technology. Bio-Rad and Vernier are teaming up to provide educators with the gold standard in hands-on science materials and technology. Vernier's expertise in data analysis complements Bio-Rad's lab activities for the life sciences. In this new partnership, Bio-Rad and Vernier are working together to broaden and enhance classroom experiences with joint experiments and curricula.

**Analysis of DNA Fingerprinting results using Vernier's Logger Pro® software.**

Logger Pro helps your students create a standard curve and calculate the number of base pairs (molecular weight) for each experimental band in only a few minutes. [www.vernier.com/biotech](http://www.vernier.com/biotech)



# Refresh.

**Refresh kit components** reduce packaging waste, reuse components, and refresh your kits while saving storage space by purchasing individual items. Bio-Rad now has many individual components for Biotechnology Explorer kits available for purchase.

Visit us at [explorer.bio-rad.com](http://explorer.bio-rad.com) to learn more. To request our kit components purchasing guide (bulletin 5872), please email us at [biotechnology\\_explorer@bio-rad.com](mailto:biotechnology_explorer@bio-rad.com).

Description	Catalog No.	Price
pGLO Bacterial Transformation Kit Refill Package	166-0555EDU	\$53.00
pGLO Plasmid, 20 µg, lyophilized	166-0405EDU	29.00
Arabinose, 600 mg, lyophilized	166-0406EDU	20.00
Ampicillin, 30 mg, lyophilized	166-0407EDU	15.00
<i>E. coli</i> Strain HB101 K-12, lyophilized	166-0408EDU	16.00
Transformation Solution, 15 ml	166-0409EDU	19.00
LB Broth, 10 ml	166-0421EDU	16.00
LB Nutrient Agar Powder, 20 g, makes forty 60 mm agar plates	166-0600EDU	10.00
LB Nutrient Agar Powder, 500 g, makes one thousand 60 mm agar plates	166-0472EDU	83.00
Petri Dishes, 60 mm, sterile, 500	166-0470EDU	119.00
Inoculation Loops, 10 µl, sterile, 80 — Limited time promotional price	166-0471EDU	9.99*
Jellyfish Foam Floating Racks, 8 racks, 12 microcentrifuge tube wells	166-0479EDU	29.00
Long-Wave UV Lamp, 1	166-0500EDU	32.00
Long-Wave UV Pen Light, 1	166-0530EDU	10.00
TAE, 50x, 1 L	161-0743EDU	56.00
TAE, 50x, 5 L cube	161-0773EDU	153.60
Sample Loading Dye, 5x, 10 ml	161-0767EDU	15.20
Certified™ Molecular Biology Agarose, 25 g	161-3100EDU	39.75
Certified Molecular Biology Agarose, 125 g	161-3101EDU	113.25
Certified Molecular Biology Agarose, 500 g	161-3102EDU	392.00
Fast Blast™ DNA Stain, 500x, 100 ml	166-0420EDU	22.00
Small Fast Blast DNA Electrophoresis Reagent Pack	166-0450EDU	48.00
Medium Fast Blast DNA Electrophoresis Reagent Pack	166-0455EDU	160.00
Large Fast Blast DNA Electrophoresis Reagent Pack	166-0460EDU	500.00
Gel Staining Trays, 4	166-0477EDU	14.00
Genes in a Bottle DNA Extraction Refill Package	166-2001EDU	37.00
Lysis Buffer, 150 ml	166-2002EDU	21.00
DNA Necklace Module, 18 amulet necklaces	166-2200EDU	33.00
MasterMix for PCR, 2x, 90 units, 1.2 ml	166-5009EDU	72.00
Disposable Plastic Transfer Pipets, sterile, 500	166-0474EDU	95.00
Disposable Plastic Transfer Pipets, nonsterile, 500	166-0480EDU	48.00
Colored 1.5 ml Microcentrifuge Tubes, 6 colors, 600	166-0473EDU	29.00
EZ Micro™ Test Tubes, 1.5 ml, natural, 500	223-9480EDU	21.60
EZ Micro Test Tubes, 2 ml, natural, 500	223-9430EDU	23.20
Conical Tubes, 1.5 ml, with installed O-ring screwcaps, sterile, 500	224-0110EDU	86.40
Conical Centrifuge Tubes, 15 ml, 50	166-0475EDU	24.00
Cell Culture Tubes, 17 x 100 mm, 14 ml, sterile, 25	166-0476EDU	14.00
0.2 ml Tubes With Domed Caps, natural, 1000	TWI-0201EDU	42.40
PCR Tube Capless Adapters, 500	223-9500EDU	17.60
96-Place PCR-Tube Rack and Cover, 5	TRC-0501EDU	28.80

\* Promotional price. Limited time offer expires January 31, 2010.

# pGLO™ Bacterial Transformation & GFP Extension Kits

Hands-on activities for AP LAB 6

## pGLO Bacterial Transformation and GFP Extension Kits

Free UV pen light included with each pGLO kit



Each kit contains sufficient materials for 8 student workstations. We recommend 2–4 students per workstation.

### pGLO Bacterial Transformation Kit

Catalog # .....EDU Price  
**166-0003EDU .....\$89.00**

### pGLO Transformation Lab Prep DVD

Catalog # .....EDU Price  
**166-0540EDU .....\$12.00**

pGLO bacterial transformation lab preparation and techniques instructional DVD contains everything you need to know to set up and run the lab like a pro.

### GFP Chromatography Kit

Catalog # .....EDU Price  
**166-0005EDU .....\$99.00**

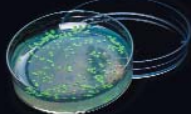
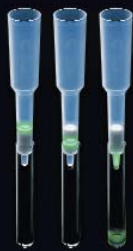
### pGLO Kit SDS-PAGE Extension

Catalog # .....EDU Price  
**166-0013EDU .....\$85.00**

Includes curriculum, 1 g DTT, 100 µl Precision Plus Protein™ Kaleidoscope™ protein standards, 100 ml Bio-Safe™ Coomassie stain, 1 L 10x TGS, 30 ml Laemmli sample buffer.

### Key Kit Features

- Aligns with AP Biology [AP LAB 6]
- Transform bacteria with jellyfish gene
- Capture the glow
- Analyze the proteome

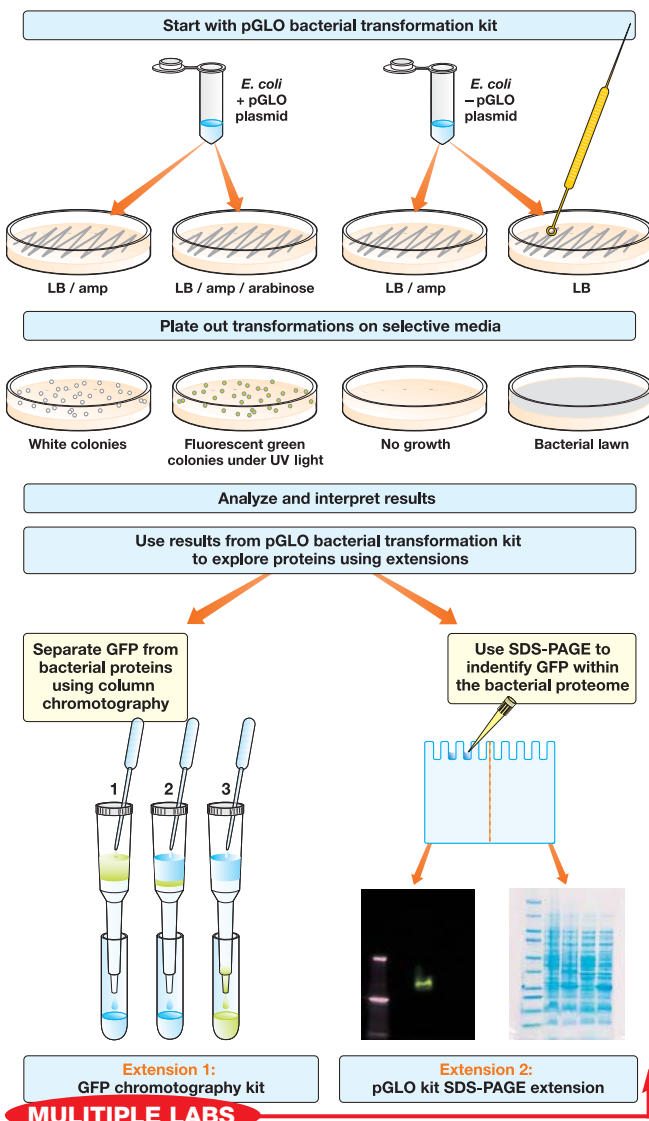


## New Refresh Kit Components

(see pages 4–5 and 18)



**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.



**MULTIPLE LABS**

Genetic engineering is the process of manipulating the genetic material of an organism — often to include the DNA from a foreign organism. In the **pGLO bacterial transformation kit**, students transform bacteria with a gene from a bioluminescent jellyfish to express a new protein, green fluorescent protein (GFP). This same procedure is used for creating “designer proteins” and has led to the explosion of new medicines, agricultural applications, and environmental solutions.

**What's next after the pGLO kit?** Transformation is only the first step in producing bioengineered proteins. Discover more about GFP with two great lab extensions that dive deeper into the proteins expressed in the pGLO bacterial transformation lab and allow you to find or purify the protein in the haystack. Use the **GFP chromatography kit** to purify GFP using a core technique in biomanufacturing: chromatography. Or, analyze the entire bacterial proteome using one of the most commonly cited molecular techniques with the **pGLO kit SDS-PAGE extension**.

Show your students the relevance of the science they learn in the classroom to science in their lives. When students genetically engineer bacteria with the genes from a bioluminescent jellyfish, they never forget the central mantra of molecular biology:

**DNA > RNA > Protein > Trait — Green Fluorescence!**

# Forensic DNA Fingerprinting Kit

Hands-on activities for AP LAB 6

## Forensic DNA Fingerprinting Kit



Each kit contains sufficient materials for 8 student workstations. We recommend 2-4 students per workstation.

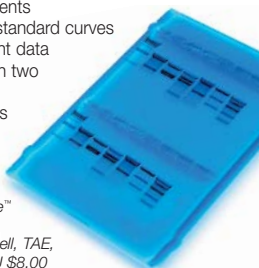
### Forensic DNA Fingerprinting Kit

Catalog # .....EDU Price  
**166-0007EDU .....\$130.00**

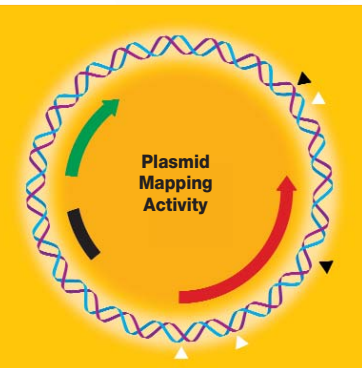
Convenient lyophilized reagents. Ships at room temperature. Immediately store temperature-sensitive reagents at -20°C.

#### Key Kit Features

- Aligns with AP Biology [AP LAB 6]
- Use electrophoresis to visualize DNA fragments
- Construct standard curves from student data
- Complete in two 45 minute lab sessions



ReadyAgarose™ precast gel  
 1.0%, 2 x 8 well, TAE,  
 161-3057EDU \$8.00  
 Find out more on page 18.



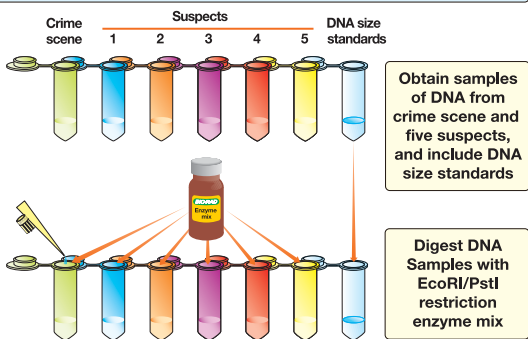
## New Jellyfish Foam Floating Racks

(see page 19)



**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.

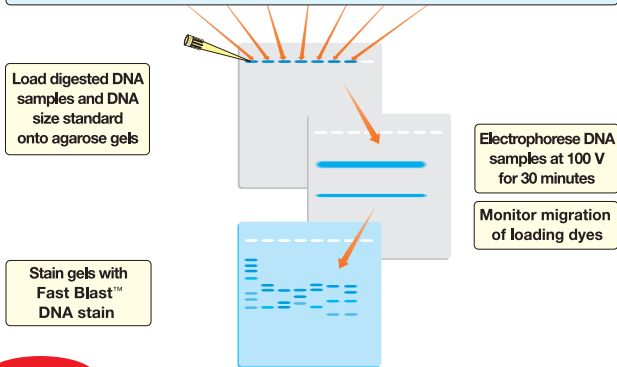
### Rehydrate DNA samples and restriction enzymes



### Incubate at 37°C for 45 minutes or overnight at room temperature

#### LAB 1

### Add loading dye to all digested DNA samples



#### LAB 2

### Match crime scene DNA with suspect's DNA. Who done it?

### Construct a standard curve using DNA size standards. Determine size of unknown fragments in DNA samples.

### Map the plasmids using restriction enzymes

#### LAB 3

**Using real DNA as evidence**, your students play the role of crime scene investigators. DNA evidence assists in criminal, missing person, mass disaster, and paternity cases. It can also be used to exonerate the innocent.

The six DNA samples in this kit are plasmids engineered to mimic the variations in DNA that exist between one human being and another. One DNA sample has been collected from a "crime scene" and five samples have been obtained from various "suspects". Each sample is digested using a mixture of two DNA restriction enzymes, generating a distinct pattern of DNA fragments. From their electrophoresis results, students construct standard curves, determine DNA fragment sizes, and place a suspect at the scene of the crime.

Cutting DNA with restriction enzymes is the first step in recombinant DNA and genetic engineering. Open the door to rich discussions about the scientific, ethical, and legal implications of DNA profiling and genetic engineering.

The ability to cut DNA with restriction enzymes makes modern molecular biology possible. Cut the DNA fingerprinting kit plasmids with the PstI and EcoRI restriction enzymes and try the new plasmid mapping activity now described in detail within this kit's curriculum manual which aligns with AP Biology Lab 6.

Get our free curriculum from the Web: [explorer.bio-rad.com](http://explorer.bio-rad.com).

# Genes in a Bottle™ Kit



Free Catalog!

CAN YOU SEE YOUR DNA?

## Genes in a Bottle Kit



Each kit contains sufficient materials for 36 students.

### Genes in a Bottle Kit

Catalog # .....EDU Price

**166-2300EDU\* ..... \$134.00**

Ships and stores at room temperature.

\* Includes one DNA extraction module and two DNA necklace modules

**166-2000EDU ..... 95.00**

DNA Extraction Module

**166-2200EDU ..... 33.00**

DNA Necklace Module (18 necklaces)

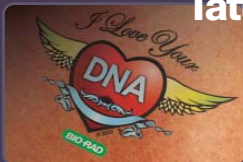
### Key Kit Features

- Use as introductory or capstone activity
- Perform real research techniques
- Extract, precipitate, and bottle your DNA
- Complete in one 45 minute lab session



DNA Necklace Extension Activity

## I Love Your DNA Tattoos



Pack of 200 temporary tattoos

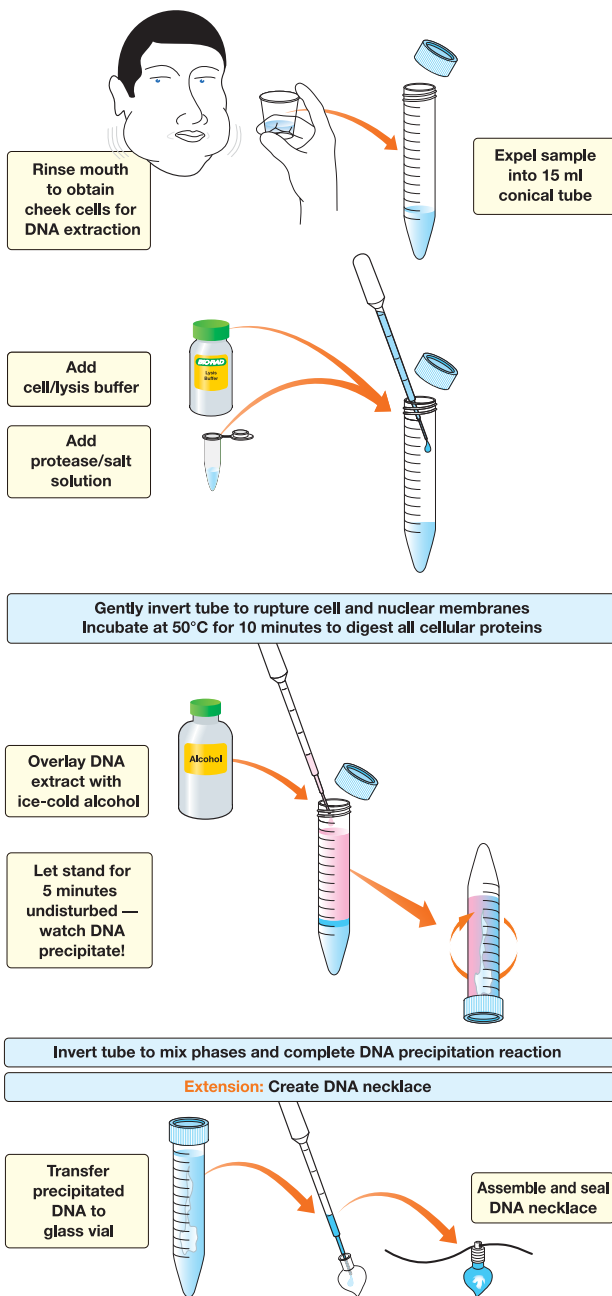
**166-2004EDU ..... 21.00**

Great for fundraising!

**email for ideas and a PDF fact sheet.**

[biotechnology\\_explorer@bio-rad.com](mailto:biotechnology_explorer@bio-rad.com)

**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.



## LAB 1

**Make your biology personal.** Introduce your students to molecular biology with their own DNA! Enable your students to see the normally invisible substance of life and begin to comprehend the meaning of their own genetic makeup. In this activity your students employ the same real-world laboratory procedure used to extract DNA from many different organisms for a variety of biotechnology research applications. Students extract genomic DNA from their own cheek cells, then precipitate and bottle it in a fabulously cool necklace.

From cell structure, to genetics, to the chemistry of life — this kit integrates multiple life science standards in a single lesson. Seeing DNA makes it real. Wearing it makes the lesson memorable!



# Free Life Sciences Workshops at a Conference Near You!

## Life Sciences Core Content Standards

**Scientific inquiry:** field and lab investigations, scientific problem solving, environmental considerations, impact of research on society, career options

**Genetics:** genes and heredity, chromosomes and alleles, mutations, variation; DNA structure, function, replication; central dogma

**Cell and molecular biology:** cell structures and processes; plant, microbe, and animal cell types; growth, development, and interaction

**Chemistry of life:** structure and function of cellular components and processes (proteins, carbohydrates, lipids, nucleic acids, etc.)

**Evolution:** change in heritable traits in populations; natural selection; allelic and species diversity, phylogeny, extinction, and fossil record

**Environmental and health science:** interactions in natural systems; biodiversity, natural resources, food webs, energy pyramid, human impact; disease and epidemiology

**Advanced concepts and techniques:** going beyond the textbook, applying the results to the real world

•	•					•		<b>Microbes and health kit</b> (166-5030EDU)	Introductory
•	•	•	•					<b>Genes in a Bottle™ kit</b> (166-2300EDU)	
•	•	•	•					<b>pGLO™ bacterial transformation kit</b> (166-0003EDU)	
•		•	•					<b>Green fluorescent protein (GFP) chromatography kit</b> (166-0005EDU)	
•	•	•	•			•		<b>Secrets of the Rainforest™ kit</b> (166-0006EDU)	
•			•					<b>Size exclusion chromatography kit</b> (166-0008EDU)	
•			•					<b>Got Protein?™ kit</b> (166-2900EDU)	Intermediate
•	•	•	•			•		<b>ELISA Immuno Explorer™ kit</b> (166-2400EDU)	
•	•		•	•				<b>Forensic DNA fingerprinting kit</b> (166-0007EDU)	
•	•	•	•					<b>Analysis of precut lambda DNA kit</b> (166-0001EDU)	
•	•	•	•					<b>Restriction digestion and analysis of lambda DNA kit</b> (166-0002EDU)	
•		•	•					<b>pGLO kit SDS-PAGE extension</b> (166-0013EDU)	
•	•		•	•				<b>Crime Scene Investigator PCR Basics™ kit</b> (166-2600EDU)	Advanced
•	•	•	•	•		•		<b>Comparative proteomics kit I: protein profiler module</b> (166-2700EDU)	
•	•	•	•	•				<b>Comparative proteomics kit II: western blot module</b> (166-2800EDU)	
•	•	•	•	•				<b>PV92 PCR informatics kit</b> (166-2100EDU)	
•	•	•	•	•				<b>GMO Investigator™ kit</b> (166-2500EDU)	
•	•		•	•		•		<b>Crime Scene Investigator PCR Basics real-time PCR starter kit</b> (166-2660EDU)	
•	•	•	•	•		•		<b>GMO Investigator real-time PCR starter kit</b> (166-2560EDU)	
•	•	•	•	•		•		<b>Cloning and Sequencing Explorer Series</b> (166-5000EDU)	

2009 Fall

## Teacher Conference and Workshop Schedule

### October

22-25	CSTA	Palm Springs	California
29-31	NSTA Midwestern Regional	Minneapolis	Minnesota
Oct 31-Nov 3	STANYS	Rochester	New York

### November

5-7	STAT-CAST	Galveston	Texas
11-14	NABT	Denver	Colorado
12-14	NSTA Eastern Regional	Fort Lauderdale	Florida
19-20	NCSTA	Greensboro	North Carolina

### December

3-5	NSTA Western Regional	Phoenix	Arizona
-----	-----------------------	---------	---------

A complete schedule can be found on the Web at [explorer.bio-rad.com](http://explorer.bio-rad.com).



## Educational Pricing Contract Application

As an educator you are entitled to discount pricing on all Bio-Rad products. To initiate setup of a new education account, simply fax or mail this page to Bio-Rad. If you are ready to place your first order, send this page and attach a purchase order with the numbers and descriptions of the items you wish to purchase.

## Biotechnology Explorer Program

Bio-Rad Laboratories, Inc.  
2000 Alfred Nobel Drive  
Hercules, CA 94547 USA

**Fax to: 1-800-879-2289  
or 1-510-741-5800**

## Re: Educational Discount

We are an educational facility interested in using Bio-Rad products to teach our students. We would like to receive a teaching lab discount and understand that the Bio-Rad education discount is only available to educators at the K-12 and undergraduate levels. This letter is our confirmation that the products we order through our educational account will be used only to educate students in a classroom or teaching lab environment and will not be used for basic scientific research. We are sending this letter in order to:

- Initiate setup of a new educational account with Bio-Rad Laboratories  
(If tax exempt, please include copy of the school's tax certificate.)
- Place a new order, with PO attached
- Confirm an order identification number (please fill in) \_\_\_\_\_

Name of contact: \_\_\_\_\_

Department: \_\_\_\_\_

School or institution: \_\_\_\_\_

We are a (please check one):

- High school     Community college     Undergraduate university     Other \_\_\_\_\_

**Bill to address:** \_\_\_\_\_

City/state/ZIP code: \_\_\_\_\_

Phone number/fax number: \_\_\_\_\_

**Ship to address:** \_\_\_\_\_

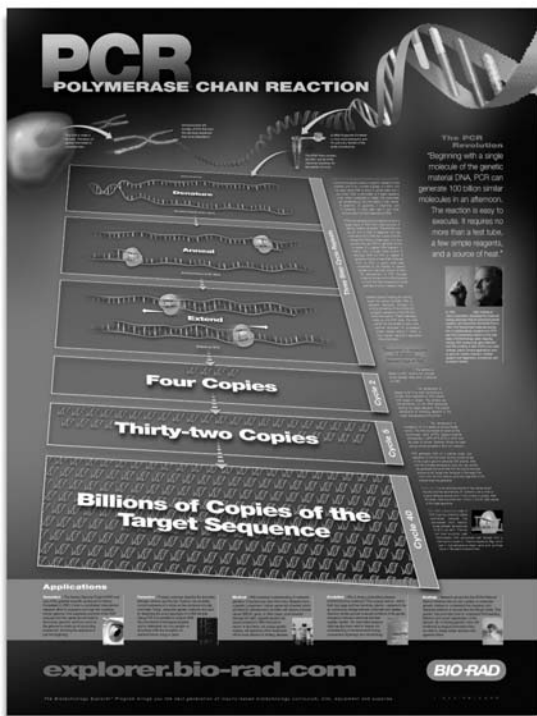
City/state/ZIP code: \_\_\_\_\_

Phone number/fax number: \_\_\_\_\_

Email address: \_\_\_\_\_

Signature: \_\_\_\_\_

# Get Your Free Biotechnology Explorer™ Stuff



## Free Giant 40" Full-Color PCR Poster

While supplies last, get your **free PCR poster** — see the detailed visual process of the polymerase chain reaction process: how it works, why it works, and what you need it for in the scientific research world.

- **Learn** about the PCR revolution
- **See** the process step by step
- **Learn** about the hot research areas using PCR now

To get your **free Biotechnology Explorer catalog**, call **1-800-424-6723**, fill out and mail the attached postage-paid card, or visit us on the Web at **explorer.bio-rad.com**.



Find out more about the **Biotechnology Explorer™ program professional development options & get some free stuff!**

20 to **40% off** list prices

Name \_\_\_\_\_ Title \_\_\_\_\_

Institution \_\_\_\_\_

Department \_\_\_\_\_ Bldg \_\_\_\_\_ Room no. \_\_\_\_\_

Address \_\_\_\_\_

City \_\_\_\_\_ State \_\_\_\_\_ ZIP code \_\_\_\_\_

( ) \_\_\_\_\_

Phone \_\_\_\_\_

( ) \_\_\_\_\_

Fax \_\_\_\_\_

Email (By providing my email address I agree to receive email about Bio-Rad products and events)

### We want to hear from you!

What are your professional development objectives? \_\_\_\_\_

\_\_\_\_\_

What challenges do you face in preparing students for biotech careers? \_\_\_\_\_

\_\_\_\_\_

Please have a Bio-Rad curriculum training specialist contact us about meeting our teaching goals and professional development needs in our own district today.

### Free stuff and more:

- Free PCR poster (bulletin 5886)
- Free Biotechnology Explorer catalog (bulletin 2112)
- Hey! Have my Bio-Rad account representative call me
- I would like to receive the Bio-Rad eFocus email newsletter

Fill out and return this card today, or for immediate information, call **1-800-4BIORAD (1-800-424-6723)**.

Bulletin 5873 09-0628 809

**Professional development — designed to fit your needs and surpass your expectations.** With over 30 years of classroom teaching experience, our curriculum training specialists know how to develop and implement curriculum within your classroom that will boost student achievement in science. Let them help you with all your training needs and make the most of your precious time and resources. **Can't travel?** They can come to your classroom, train the educators within your region or district, and help you find the resources you need to make teaching biotech exciting and engaging!

**Contact us today for:**

- Personalized training to meet your curriculum goals
- Laboratory activities aligned with your education standards
- Help setting up a biotech lab or curriculum

[professional\\_development@bio-rad.com](mailto:professional_development@bio-rad.com)



SYBR® is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR® Green I for use in real-time PCR, for research purposes only. Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

*Notice regarding Bio-Rad thermal cyclers and real-time systems:* Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160-163 and 167 only), and 6,703,236 (claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

This product is covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's United States Patent 6,814,934 B1 for use in research and for all other fields except veterinary diagnostics.

Coomassie is a trademark of BASF Aktiengesellschaft. GenBank is a trademark of United States Department of Health and Human Services. Geospiza is a trademark of Geospiza, Inc. Vernier and Logger Pro are trademarks of Vernier Software & Technology.



NO POSTAGE  
NECESSARY  
IF MAILED  
IN THE  
UNITED STATES

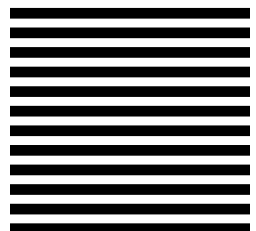
**BUSINESS REPLY MAIL**

FIRST CLASS PERMIT NO 42 RODEO, CALIF

POSTAGE WILL BE PAID BY ADDRESSEE



Bio-Rad Laboratories, Inc.  
2000 Alfred Nobel Drive  
Hercules, CA 94547-9980



# Get a Set of PCR-Tube Racks and One PCR Lab Kit Free!

## MyCycler™ Thermal Cycler Special Offer

### Free 96-Place PCR-Tube Rack and Cover Set

See page 18 for more information (TRC-0501EDU).

### Your Choice of One PCR Kit Free!

See below for details.

### MyCycler Thermal Cycler

With its 96-well sample capacity and rapid heating and cooling technology, this compact thermal cycler provides high sample throughput in less time. This classroom-scale instrument also features a large 12 cm screen, intuitive programming, and real-time graphical display of your PCR protocol, at an economical price.

### Specifications

Sample capacity	96 wells x 0.2 ml
Speed of ramping	Up to 2.5°C/sec
Temperature range	4–100°C
Heated lid	Adjustable to 110°C
Reaction volumes	15–100 µl
Display	12 cm high-resolution 1/4 VGA
Ports	1 USB
Memory	99 typical programs
Dimensions	25 x 44 x 21 cm (10 x 17 x 8") (W x D x H)

## MyCycler Thermal Cycler Special Offer

Purchase a MyCycler thermal cycler and get a free set of PCR-Tube racks and your choice of one PCR kit free — over \$188 value. Choose from the following kits: Crime Scene Investigator PCR Basics™ kit, PV92 PCR informatics kit, and GMO Investigator™ kit (Quote #09-Q25874).

Catalog #	Description	List Price	EDU price	Special Offer
170-9701EDU	MyCycler Thermal Cycler, 96 wells, 120/240V	4,750.00	3,575.00	<b>3,575.00</b>
TRC-0501EDU	96-Place PCR-Tube Rack and Cover, 5	36.00	28.80	<b>Free</b>



### Choose one of the following PCR kits:

166-2600EDU	Crime Scene Investigator PCR Basics kit	200.00	160.00	<b>One Free</b>
166-2100EDU	PV92 PCR informatics kit	248.75	199.00	
166-2500EDU	GMO Investigator kit	218.75	175.00	

Must reference quote #09-Q25874. Provide all catalog numbers including free PCR racks and PCR kit choice when placing order. Limited to two MyCycler thermal cyclers with free PCR racks and kit per customer. Limited-time offer expires January 31, 2010.

HAVE YOUR FAVORITE FOODS BEEN GENETICALLY MODIFIED?

## GMO Investigator Kit



Each kit contains sufficient materials for 8 student workstations. We recommend 2–4 students per workstation.

### GMO Investigator Kit'

Catalog # .....EDU Price  
**166-2500EDU ..... \$175.00**

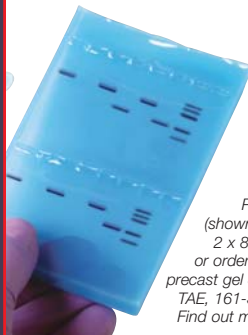
Required DNA electrophoresis reagents available separately; see p. 18. Ships at room temperature. Immediately store temperature-sensitive reagents as indicated.

### GMO Investigator Kit' Plus Small DNA Electrophoresis Reagent Pack

Catalog # .....EDU Price  
**166-2550EDU ..... \$219.00**

#### Key Kit Features

- Extract and amplify DNA from 8 food samples
- Perform genuine diagnostic procedures
- Use PCR and electrophoresis to find GM foods
- Complete in three 45 minute lab sessions



Pour your own gel (shown) 3.0% agarose, 2 x 8 well, TAE buffer, or order ReadyAgarose™ precast gel 3.0%, 1 x 8 well, TAE, 161-3017EDU \$6.40. Find out more on page 18.

## MJ Mini™ 48-Well Thermal Cycler (see page 19)

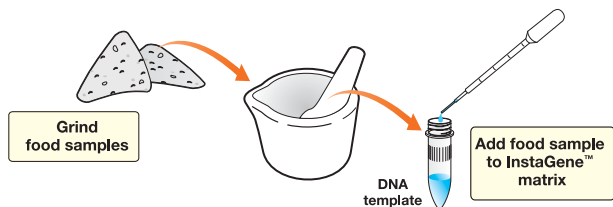


Hot real-time PCR kit extension  
 Free downloadable PDFs

**Web Check This Out!**

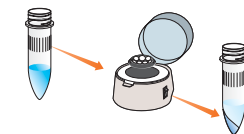
[explorer.bio-rad.com](http://explorer.bio-rad.com)

**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.



Incubate at 100°C for 5 minutes

**Centrifuge samples for 5 minutes to pellet matrix**

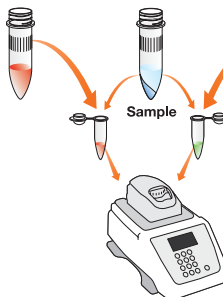


DNA template preparation

### LAB 1

**Red** master mix with GMO primers

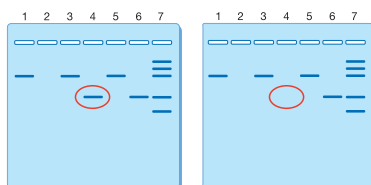
**Green** master mix with plant primers



**Place tubes in thermal cycler and amplify target DNA sequences**

Set up polymerase chain reaction and amplify DNA samples

### LAB 2



PCR sample lanes:  
 1 Non-GMO food with plant primers  
 2 Non-GMO food with GMO primers  
 3 Test food with plant primers  
 4 Test food with GMO primers  
 5 GMO-positive control DNA with plant primers  
 6 GMO-positive control DNA with GMO primers  
 7 PCR molecular weight ruler

Electrophorese PCR products and stain gels

### LAB 3

Dry gels and analyze results

**Genetically modified (GM) foods** do not require labeling as such in the U.S. Using state-of-the-art DNA extraction techniques, PCR, and electrophoresis, students determine whether their favorite foods have been genetically modified — or not!

Students test for the presence of two different genetically modified organism (GMO)-associated DNA sequences present in most (>85%) of the genetically modified crops that are approved for distribution worldwide. This kit incorporates multiple controls, allowing students to engage in a complete investigation and to assess the validity of their results.

Are GM crops a good thing? Some argue that there is a potential for “superweeds” to arise through cross-pollination or for “superbugs” to evolve. Proponents of GM crops argue that they increase crop yield and reduce the use of chemicals that are potentially toxic to the environment and human health. Regardless of where you stand in the GM debate, wouldn't it be interesting to know if the corn- or soy-based foods you eat are GM foods?

Download the free curriculum from the Web and see how at [explorer.bio-rad.com](http://explorer.bio-rad.com).

# Crime Scene Investigator PCR Basics™ Kit

WHICH HUMAN DNA SEQUENCES ARE USED IN FORENSIC INVESTIGATIONS?

## Crime Scene Investigator PCR Basics Kit



Each kit contains sufficient materials for 8 student workstations. We recommend 2-4 students per workstation.

### Crime Scene Investigator PCR Basics Kit'

Catalog # .....EDU Price  
**166-2600EDU ..... \$160.00**

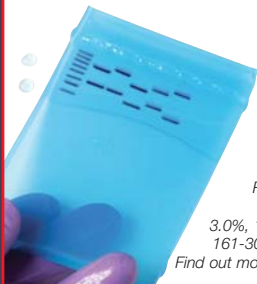
Required DNA electrophoresis reagents available separately. Ships at room temperature. Immediately store temperature-sensitive reagents at -20°C.

### Crime Scene Investigator PCR Basics Kit' Plus Small DNA Electrophoresis Reagent Pack

Catalog # .....EDU Price  
**166-2650EDU ..... \$205.00**

#### Key Kit Features

- Perform genuine DNA profiling
- Use PCR to amplify multiple DNA samples
- Use electrophoresis to visualize results
- Complete in two 45 minute lab sessions



ReadyAgarose™ precast gel  
 3.0%, 1 x 8 well, TAE,  
 161-3017EDU \$6.40.  
 Find out more on page 18.

## New Refresh Kit Components

(see pages 4-5 and 18)

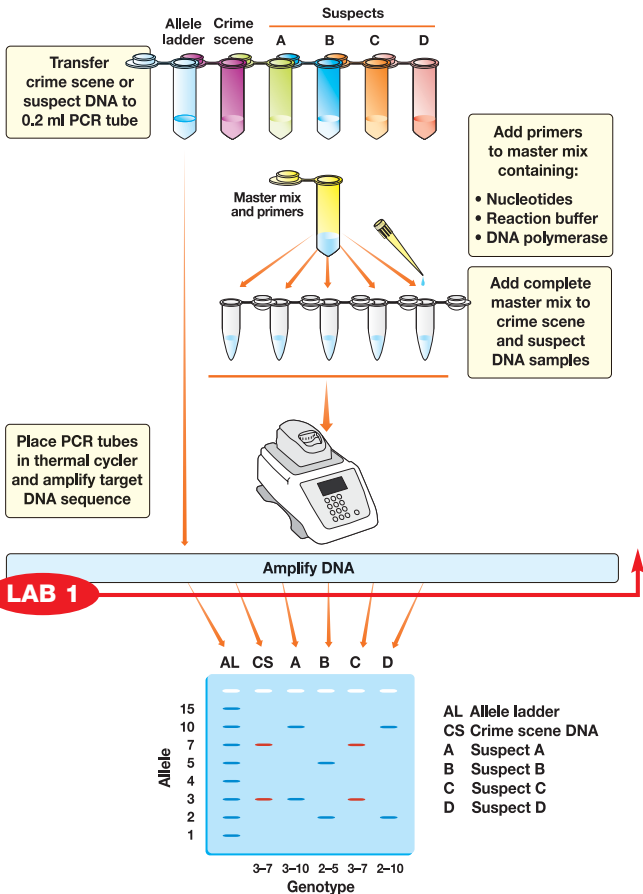


Hot real-time PCR kit extension  
 Free downloadable PDFs

**Web Check This Out!**

[explorer.bio-rad.com](http://explorer.bio-rad.com)

**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.



**LAB 2**

Electrophorese PCR samples in agarose gels at 100V for 30 minutes. Stain with Fast Blast™ DNA stain

Determine genotypes of samples from suspects and crime scene. Use the "power of discrimination" to verify the likelihood of a genotype match due to chance

**Extension:** Use web-based interactive animated tutorial

**Using real-world forensic lab techniques**, students learn which human DNA sequences are used for crime scene, missing person, mass disaster, and paternity investigations. Using the polymerase chain reaction (PCR) and DNA electrophoresis, students determine the genotypes of five DNA samples and experience how real-world crime labs solve mysteries.

DNA profiling determines the exact genotype of a DNA sample and distinguishes one human being from another by identifying a DNA "bar code" that is unique to every individual. There are over 3 billion bases in the human genetic blueprint, and more than 99.5% of them do not vary between human beings. Within the variant areas of the genome exist unique polymorphic sequences called short tandem repeats (STRs) of uncertain biological function. It is these DNA sequences that are used in forensic applications.

The DNA samples contained in this **introductory PCR kit** are plasmids that have been engineered to mimic the natural variations in human DNA that exist between one human being and another at a single STR locus. Students determine the genotypes of five samples and perform a simple statistical analysis on their own results to solve a forensic mystery.

# A Complete, Modular and Flexible Teaching Solution:



- Nested PCR
- Degenerate primers
- Exonuclease treatment

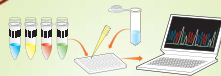


- Agarose gel analysis
- Visualize PCR products



- Genomic DNA extraction
- Cellular membrane disruption
- Isolation of genomic DNA

## It's Your Choice Complete Or Individual



- Directional sequencing primers
- Automated sequencing
- Web-based data analysis
- Contig assembly
- Intron/exon prediction



## Free Full-Color Printed Curriculum Manual Special Offer

**Purchase the (eight module) Cloning and Sequencing Explorer Series and get a free Full-Color Printed Cloning and Sequencing Explorer Series Curriculum Manual — a \$236 value. (Quote #09-Q25875)**

Catalog #	Description	List Price	EDU price
166-5000EDU	Cloning and Sequencing Explorer Series	\$1,868.75	\$1,495.00
166-5002EDU	Full-Color Printed Cloning and Sequencing Explorer Series Curriculum Manual	236.25	<b>Free</b>

Must reference quote #09-Q25875. Provide all catalog numbers when placing order. Limited-time offer expires January 31, 2010.

**Prepare your students for the real world of scientific research — publish novel research in GenBank!** Engage them with the opportunity to perform novel, relevant research that can actually contribute to the body of scientific knowledge. In this unique **modular lab series**, students are guided through an innovative research workflow identical to those performed in genomics labs worldwide. Over a multiple-week lab course, students will combine traditional and cutting-edge molecular biology techniques and bioinformatics. Your students will clone, sequence, and analyze a housekeeping gene from a plant of your choice allowing each class to produce unique and novel data.



# The Cloning and Sequencing Explorer Series

## Series Modules



- PCR product purification
- Size exclusion chromatography



- Blunting of PCR products
- Ligation
- Create competent bacteria
- Bacterial transformation



- Grow transformed cells
- Cell selection
- Ampicillin resistance



- Plasmid isolation
- Restriction enzyme digestion

### Complete Cloning & Sequencing Explorer Series Eight Modules

Catalog #                      List Price      . . . . . EDU Price  
**166-5000EDU**    \$1,868.75    . . . **\$1,495.00**

Each module contains materials sufficient for 12 student workstations.

Ships with both temperature-sensitive and room-temperature components. Immediately store temperature-sensitive items at 4°C or -20°C as indicated.

#### Series Contents:

Nucleic acid extraction module	1
GAPDH PCR module	1
Electrophoresis module	1
PCR Kleen™ spin purification module	1
Ligation and transformation module	1
Microbial culturing module	1
Aurum™ plasmid mini purification module	1
Sequencing and bioinformatics module	1
EZ Micro™ test tubes	1
Curriculum resource CD	1
(includes curriculum manual, overview webinar presentation, and master planning guide)	

### Individual Cloning & Sequencing Explorer Series Modules and Supplies

Catalog #	List Price	. . . . .	EDU Price
<b>Nucleic Acid Extraction Module</b>			
<b>166-5005EDU</b>	\$206.25	. . . . .	<b>\$165.00</b>
<b>GAPDH PCR Module</b>			
<b>166-5010EDU</b>	483.75	. . . . .	<b>387.00</b>
<b>Small Ethidium Bromide DNA Electrophoresis Reagent Pack</b>			
<b>166-0451EDU</b>	115.00	. . . . .	<b>58.00</b>
<b>Small Fast Blast™ DNA Electrophoresis Reagent Pack</b>			
<b>166-0450EDU</b>	92.00	. . . . .	<b>48.00</b>
<b>PCR Kleen™ Spin Purification Module</b>			
<b>732-6300EDU</b>	77.14	. . . . .	<b>54.00</b>
<b>Ligation and Transformation Module</b>			
<b>166-5015EDU</b>	377.50	. . . . .	<b>302.00</b>
<b>Microbial Culturing Module</b>			
<b>166-5020EDU</b>	118.75	. . . . .	<b>95.00</b>
<b>Aurum™ Plasmid Mini Purification Module</b>			
<b>732-6400EDU</b>	212.86	. . . . .	<b>149.00</b>
<b>Sequencing and Bioinformatics Module</b>			
<b>166-5025EDU</b>	312.50	. . . . .	<b>250.00</b>
<b>EZ Micro™ Test Tubes</b>			
<b>223-9503EDU</b>	39.00	. . . . .	<b>31.20</b>
<b>Curriculum Resource CD</b>			
<b>166-5001EDU</b>	36.25	. . . . .	<b>29.00</b>
<b>Full-Color Printed Cloning and Sequencing Explorer Series Curriculum Manual</b>			
<b>166-5002EDU</b>	236.25	. . . . .	<b>189.00</b>

**Educational discounts** apply only to items ordered with an EDU suffix. EDU price discounts are for qualified educational institutions and educators only.

## New Refresh Kit Components



- LB Nutrient Agar Powder**  
500 g, makes 1,000 plates (60 mm)  
**166-0472EDU \$83.00**



- Petri Dishes (500)**  
60 mm, sterile  
**166-0470EDU \$119.00**



- Inoculation Loops (80)**  
10 ul, sterile  
**166-0471EDU \$9.99\***  
*\* Limited-time offer expires January 31, 2010.*



- Gel Staining Trays (4)**  
**166-0477EDU \$14.00**



- Disposable Plastic Transfer Pipets (500) Sterile**  
**166-0474EDU \$95.00**



- Disposable Plastic Transfer Pipets (500) Nonsterile**  
**166-0480EDU \$48.00**



- 1.5 ml Microcentrifuge Tubes (600) 6 colors**  
**166-0473EDU \$29.00**



- Conical Centrifuge Tubes (50) 15 ml**  
**166-0475EDU \$24.00**



- Cell Culture Tubes (25)**  
17 x 100 mm, 14 ml, sterile  
**166-0476EDU \$14.00**



ReadyAgarose gels ship and store at room temperature.

- ReadyAgarose™ Precast Gels**  
Individually packaged precast agarose gels provide resolution of nucleic acids from 20 to 20,000 base pairs.

- UV-transparent trays
- ReadyAgarose gels lock into Mini-Sub® cell GT or ReadySub-Cell™ GT cells electrophoresis chambers.

Visit us on the Web for a complete gel listing at [discover.bio-rad.com](http://discover.bio-rad.com).



- 96-Place PCR-Tube Rack and Cover**  
Stackable storage units for 0.2 ml PCR tubes, 5  
**TRC-0501EDU \$28.00**



- PowerPac™ Mini Power Supply**  
• 100 V or 200 V constant  
• 200 mA maximum current  
• 40 W maximum power  
• Constant voltage output  
**165-5048EDU \$175.00**



EN-61010 safety compliant

- PowerPac™ Basic Power Supply**  
Constant voltage or current output, real-time monitoring, and pause/resume features  
• Runs 4 cells simultaneously  
• 400 mA maximum current  
• 10–300 V in 1 V steps  
• Timer controlled  
• Fully programmable  
**164-5050EDU \$341.25**



- Mini-PROTEAN® Tetra Cell**  
for 1-D vertical gel electrophoresis.  
• Runs one to four mini gels  
• Easy assembly  
• Leak-free electrophoresis  
• Run precast or handcast gels



- 165-8004EDU \$390.75**  
4-gel system



- 165-8005EDU 267.00**  
2-gel system



- Mini-PROTEAN Tetra Cell and Mini Trans-Blot Module**  
for Ready Gel® precast gels includes 165-8004EDU and 170-3935EDU  
**165-8030EDU \$628.50**

- Mini Trans-Blot Module**  
Includes 2 gel holder cassettes, 4 fiber pads, modular electrode assembly, blue cooling unit only (no tank or lid)  
**170-3935EDU \$285.75**



- DNA Electrophoresis Systems**

- Mini-Sub® Cell GT Cell**  
Includes 7 x 10 cm gel tray and two 8-well combs  
**166-4000EDU \$225.00**

- Mini ReadySub-Cell™ GT Cell**  
for ReadyAgarose™ precast gels.  
**170-4487EDU 175.00**

SYBR® is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR® Green I for use in real-time PCR, for research purposes only. Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Notice regarding Bio-Rad thermal cyclers and real-time systems: Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents 5,656,493, 5,333,673, 5,475,610 (claims 1, 44, 158, 160-163 and 167 only), and 6,703,236 (claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

This product is covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's United States Patent 6,814,934 B1 for use in research and for all other fields except veterinary diagnostics.

Coomassie is a trademark of BASF Aktiengesellschaft. GenBank is a trademark of United States Department of Health and Human Services. Geospiza is a trademark of Geospiza, Inc. Vernier and Logger Pro are trademarks of Vernier Software & Technology.



Free Catalog!

20-40% off

### Cool **New** Floaty



- Jellyfish Foam Floating Racks**  
Package of 8 racks,  
12 microcentrifuge tube wells  
**166-0479EDU** **29.00**



- Water Bath**  
Temperature-controlled,  
dependable, and affordable
    - Stainless-steel tank and lid
    - Temperature range up to 100°C
    - Over-temperature protection
    - Includes thermometer
- 166-0504EDU** **\$590.00**



- MJ Mini™ Thermal Cycler**  
**PTC-1148EDU** **\$3,196.00**



- Mini Incubation Oven**  
80-dish capacity, up to 60°C,  
includes thermometer  
**166-0501EDU** **\$367.00**



- MiniOpticon™ Two-Color Real-Time PCR Detection System**  
Precise thermal control,  
temperature gradient capability,  
performs applications including  
quantitative real-time PCR,  
relative gene expression  
analysis, and allelic discrimination



- Mini Centrifuge** (2,000 x g)  
Includes microtube and PCR  
strip tube rotors, and 0.4 ml and  
0.5 ml tube adaptors  
• Max. speed 6,000 rpm  
**166-0603EDU** **\$299.00**

- Compact, portable system
  - 48-well plate or 6 x 8-tube strips
  - 10–100 µl sample volumes
- CFB-3120EDU** **\$15,196.00**



- Model 16K** (16,000 x g)  
**Microcentrifuge**
    - 1.5 ml or 2.0 ml tubes
    - Quick-spin feature
    - Max. speed 14,000 rpm
- 166-0602EDU** **\$1,900.00**



- Professional Adjustable-Volume Digital Micropipets**

- Slender, contoured grip
  - Ergonomic tip ejector
  - Autoclavable
- 166-0505EDU** (0.5–10 µl)  
**166-0506EDU** (2–20 µl)  
**166-0507EDU** (20–200 µl)  
**166-0508EDU** (100–1,000 µl)  
Each pipet only **\$219.00**



- PCR Tube Adaptor**  
(for #166-0602EDU)  
Holds two PCR 8-tube strips  
or 16 individual 0.2 ml tubes  
**166-0620EDU** **\$105.00**



- BR-2000 Vortexer**  
General-purpose vortex mixer  
**166-0610EDU** **\$275.00**



- Classroom Adjustable-Volume Digital Micropipets**

- 166-0550EDU** (0.5–10 µl)  
**166-0551EDU** (2–20 µl)  
**166-0552EDU** (20–200 µl)  
**166-0553EDU** (100–1,000 µl)  
Each pipet only **\$110.00**



- Flathead Dimpled Adaptor**  
(for #166-0610EDU)  
**166-0622EDU** **\$45.00**



- Fixed-Volume Micropipets**

- 166-0511EDU** (5 µl)  
**166-0512EDU** (10 µl)  
**166-0513EDU** (20 µl)  
**166-0515EDU** (50 µl)  
Each pipet only **\$24.00**



- Long-Wave UV Pen Light**  
Small, disposable ultraviolet light  
**166-0530EDU** **\$10.00**



- Long-Wave UV Lamp**  
Portable, mini ultraviolet lamp  
**166-0500EDU** **\$32.00**

**Discount EDU prices.**  
Get free quotes.

Or, set up a **Bio-Rad account today!**

Use your Visa, MasterCard, or American Express card.



Prices shown reflect education discount and are limited to qualified educational institutions and educators only. \* Limited-time offers expire January 31, 2010.

Email: Send your request for a price quote to: [lsg.quotes.us@bio-rad.com](mailto:lsg.quotes.us@bio-rad.com).

Phone: 1-800-424-6723 x6757. Always include catalog number(s), your name, address, and phone and fax numbers. (Quote requests may also be left on voicemail.)

**explorer.bio-rad.com**

For more laboratory equipment or for immediate information, call 1-800-4BIORAD (1-800-424-6723).

PRESORTED STD  
U.S. POSTAGE  
**PAID**  
Lincoln, NE  
Permit No. 365

**BIO-RAD**

**Bio-Rad Laboratories, Inc.**  
2000 Alfred Nobel Drive  
Hercules, CA 94547

**Address service requested**

# Get a Set of PCR-Tube Racks and One PCR Lab Kit Free!

See page 13 for more information.



**Free 96-Place  
PCR-Tube Rack  
and Cover Set**



**MyCycler  
Thermal Cycler**

**MyCycler™  
Thermal Cycler**  
Special Offer

**Your Choice of  
One PCR Kit Free!**



## New Giant-Sized 40" Full-Color **PCR Poster**

See page 11 to order yours today — **it's free!**



For a free catalog or to order, call: 1-800-4BIORAD (1-800-424-6723). Use your Visa, MasterCard, or American Express card. Discounts and specials may not be combined with other promotional offers. Limited-time offers expire January 31, 2010.