
Biotechnology Explorer™

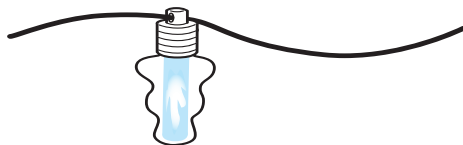
Genes in a Bottle™ Kit DNA Helix Necklace Module

Instruction Manual
Catalog #166-2250EDU

(Module contains 36 helix necklaces)

Duplication of any part of this document is permitted for classroom use only. Please visit explorer.bio-rad.com to access our selection of language translations for Biotechnology Explorer kit curricula.

For technical support, call your local Bio-Rad office, or in the U.S., call 1-800-424-6723



BIO-RAD

Genes in a Bottle: Capture Your Unique Essence!

Once your students have extracted genomic DNA from their cheek cells using the DNA extraction module (catalog #166-2000EDU), the DNA strands can be collected and transferred to a helix vial. The helix vial is then fashioned into a necklace that can be worn with pride, kept for posterity, or shared with a loved one!

Be the first on your block to wear DNA!

Read more: explorer.bio-rad.com.

Learning opportunities for all levels of instruction.

This activity is designed for any classroom environment and requires no specialized equipment or stains. For middle school students, it's a perfect introduction to the exciting world of DNA science. For secondary and college level instruction, lessons on DNA structure and function, cell structure, and enzyme function can be introduced or reinforced with this laboratory activity.

We welcome your comments and suggestions. Have fun!

Biotechnology Explorer Team
Bio-Rad Laboratories
6000 James Watson Drive
Hercules, CA 94547
explorer.bio-rad.com

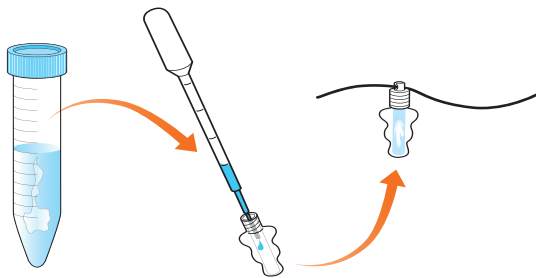
The DNA helix necklace module contains enough material to prepare 36 DNA helix necklaces.

Inventory Check List	Amount Provided
Helix vials	36
Screw caps	36
Cords	36

Genes in a Bottle: Capture Your Unique Essence!

Instructions

1. Using a disposable plastic transfer pipet, carefully transfer an appropriate portion of the DNA in alcohol into the helix vial. You do not have to fill the entire vial with alcohol. (Note that students can share plastic transfer pipets for transferring their DNA into the helix vial.)
2. Place the screw cap onto the top of the helix vial and tighten.
3. Slip the cord through the screw cap and tie the cord.



Optional: To permanently seal your DNA necklace, add super glue to the outside threads of the helix vial (do not allow super glue to contact the alcohol solution or hazing may occur), and then tighten the cap. Allow super glue to dry with cap tightened.

Congratulations, you've created your very own DNA helix necklace!

BIO-RAD

**Bio-Rad
Laboratories, Inc.**



*Life Science
Group*

Web site www.bio-rad.com **USA** 800 424 6723
Australia 61 2 9914 2800 **Austria** 01 877 89 01
Belgium 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532
Denmark 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0
Greece 30 210 9532 220
Hong Kong 852 2789 3300
Hungary 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091
Japan 03 6361 7000 **Korea** 82 2 3473 4460
Mexico 52 555 488 7670
The Netherlands 0318 540666
New Zealand 64 9 415 2280
Norway 23 38 41 30 **Poland** 48 22 331 99 99
Portugal 351 21 472 7700
Russia 7 495 721 14 04
Singapore 65 6415 3188
South Africa 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700
Switzerland 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000