

12/9/cl 12/8/98

GenePath Program Upgrade Kit

Instruction Manual

Catalog Number

310-0075



Table of Contents

	Page
Section 1 Introduction	1
1.1 Components	1
1.2 Additional Items Required	1
1.3 Precautions	1
Section 2 Instructions	1
2.1 Installing the RS-232 Cable	1
2.2 Download the GenePath Programs	1
Section 3 Programs	3
3.1 Programs Parameters	3
Section 4 Troubleshooting	5

Caution: Please read thoroughly and familiarize yourself with the entire contents of this instruction manual before using the product for the first time.

Section 1 Introduction

The GenePath instrument is based on the pulsed field method of CHEF (Contoured Clamped Homogeneous Electric Field). The system has been preprogrammed to analyze a variety of organisms (*E. coli*, *Pseudomonas*, *Enterobacter*, *Enterococcus*, *Staphylococcus*, and *Candida*). The system was also designed to easily accept new programs. The GenePath Program Upgrade Kit contains 15 new programs that are based on the separation of a particular size range.

1.1 Components

Item	Quantity
GenePath Program Upgrade disk	1
RS-232 Cable	1
Manual	1

1.2 Required Items

Item

PC Computer with the following minimum requirements:

- 486 co-processor
- DOS 6.x Operating System
- 8 Mb of RAM memory
- 3.25 inch disk drive
- 9-pin serial or COM port

1.3 Precautions

The program is to be used only with the GenePath System.

Section 2 Instructions

2.1 Installing the RS-232 Cable

1. Connect the female end of the 9-pin RS-232 cable to either the COM1 or COM2 port (serial port) on the PC computer. For most PC's this port is located on the back panel.
2. Connect the male end of the 9-pin RS-232 cable to the port on the back of the GenePath.

2.2 Downloading the GenePath Programs

1. Turn on the PC computer. Turn on the GenePath and use the INCREASE or DECREASE keys on the front panel to select the data load mode (display shows dLd).
2. Insert the GenePath Upgrade Program disk into drive A on the computer.
3. In DOS, select the A drive by typing in **A:**. Type in **gpsend** to start the GenePath program.

4. The main menu of the GenePath Download Program will be displayed. To view an explanation of a menu item, use the Up and Down keys to highlight the desired menu item and press the ENTER key. When you are finished reading the explanation, press ENTER to return to the main menu.
5. Select **Setup the RS-232 parameters** and press ENTER. Enter the number of the COM port (1 or 2) to which the RS-232 cable was connected as directed in Section 2.1. Press ENTER to return to the main menu.
6. Select **Check GenePath Status** and press ENTER. The message "The GenePath is Ready" should be displayed. Press ENTER to return to the main menu. If the message "The GenePath is not ready" is displayed, check that the GenePath is in the "dLd" program, that the RS-232 cable is firmly connected, and that the correct COM port was selected.
7. Select **Download to GenePath** and press ENTER. The program will download to the GenePath. During this process the bytes sent and percent completion are displayed. An irregular clicking sound will be made by the GenePath, indicating the download is in progress. After about 35 seconds, the clicking will stop and a message will be displayed indicating the download is complete. Press ENTER to return to the main screen.
Note: If the program needs to be stopped during the download process, press the escape key (ESC). Press ENTER to return to the main screen. To re-install the new programs, turn the GenePath system off and then back on. Repeat the downloading process from step 2.2.6.
8. To leave the GenePath Upgrade Program, select **Exit to DOS** and press ENTER.
9. Check that the new programs have been downloaded to the GenePath. All of the programs listed in Section 3 should be accessible.

Section 3 Programs

3.1 Program Parameters

Program 1: PSU

This program is to separate *Pseudomonas*.

Run time: 19.5 hours

Program 2: Eco

This program is to separate *E.coli*.

Run time: 20.0 hours

Program 3: Enb

This program is to separate *Enterobacter*.

Run time: 17.0 hours

Program 4: CAn

This program is to separate *Candida*.

Run time: 48.0 hours

Program 5: StA

This program is to separate *Staphylococcus*.

Run time: 20.0 hours

Program 6: Enc

This program is to separate *Enterococcus*.

Run time: 20.0 hours

Program 7: SCREEN

This program is to quickly check your sample plugs for preliminary analysis.

Run time: 4.0 hours

Program 8: 1-50

This program is to separate DNA in the size range of 1-50 Kb.

Run time: 7.0 hours

Program 9: 1-100

This program is to separate DNA in the size range of 1-100 Kb.

Run time: 12.0 hours

Program 10: 1-200

This program is to separate DNA in the size range of 1-200 Kb.

Run time: 12.7 hours

Program 11: 25-250

This program is to separate DNA in the size range of 25-250 Kb.

Run time: 18.5 hours

BIO-RAD**Bio-Rad
Laboratories****Diagnostics
Group**

4000 Alfred Nobel Drive
Hercules, California 94547
Telephone 510 724-7000
FAX 510 741-5824

www.bio-rad.com

AUSTRALIA, Bio-Rad Laboratories Pty. Ltd. Unit 1 Block Y, 391 Park Road, Regents Park NSW 2143 • Phone 61-2-9914-2800 • Telefax 61-2-9914-2888
AUSTRIA, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, A-1130 Vienna • Phone 43-1-877-8901 • Telefax 43-1-876-5629
BELGIUM, Bio-Rad S.A.-N.V. Begoniastraat 5, B-9810 Nazareth Eke • Phone 32-9-385-5511 • Telefax 32-9-385-8554
CANADA, Bio-Rad Laboratories, Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone 1-905-712-2771 • Telefax 1-905-712-2990
CHINA, Bio-Rad China Limited, 14 Zhi Chun Road, Haidian District, Beijing 100088 • Phone 86-10-6205-1850 • Telefax 86-10-6205-1876
DENMARK, Bio-Rad Laboratories, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 45-39-17-99-47 • Telefax 45-39-27-16-98
FINLAND, Bio-Rad Laboratories, Pihatorma 1 A, FIN-02240 Espoo • Phone 358-9-804-22-00 • Telefax 358-9-804-11-10
FRANCE, Bio-Rad S.A., 94-96 rue Victor Hugo B.P.220, 94203 Ivry sur Seine Cedex • Phone 33-1-43-90-4690 • Telefax 33-1-46-71-2467
GERMANY, Bio-Rad Laboratories GmbH, Heidemannstrasse 164, D-80939 Munich • Phone 49-89-318840 • Telefax 49-89-318-84100
HONG KONG, Bio-Rad Pacific Ltd., Unit 1111, 11F, New Kowloon Plaza 38, Tai Kok Tsui Rd., Tai Kok Tsui, Kowloon • Phone 852-2789-3300 • Telefax 852-2789-1257
INDIA, Bio-Rad Laboratories (India) Pte. Ltd., C-248 Defence Colony, New Delhi 110024 • Phone 91-11-460-3676 • Telefax 91-11-461-0765
ISRAEL, Bio-Rad Laboratories Ltd., 14 Homa Street, New Industrial Area, Rishon Le Zion 75150 • Phone 972-3-9514127 • Telefax 972-3-9514129
ITALY, Bio-Rad Laboratories S.r.l., Via Cellini 18/A, 20090 Segrate, Milan • Phone 39-2-216091 • Telefax 39-2-21609-399
JAPAN, Nippon Bio-Rad Laboratories K.K., 7-18 Higashi-Nippori 5-Chome, Arakawa-ku, Tokyo 116 • Phone 81-3-5811-6290 • Telefax 81-3-5811-6262
KOREA, Bio-Rad Korea Ltd., 6F Hang Woo Building, 1342-3 Seocho-Dong, Seocho-Ku, Seoul • Phone 82-2-3473-4460 • Telefax 82-2-3472-7003
THE NETHERLANDS, Bio-Rad Laboratories BV, Fokkerstraat 10, 3905 KV Veenendaal • Phone 31-318-540666 • Telefax 31-318-542216
NEW ZEALAND, Bio-Rad New Zealand, Unit B 156 Bush Road, Albany, Auckland • Phone 64-9-415-2280 • Telefax 64-9-415-2284
SINGAPORE, Bio-Rad Laboratories (Singapore) Pte. Ltd., 221 Henderson Road #05-19 Henderson Building, Singapore 159557 • Phone 65-272-9877 • Telefax 65-273-4835
SPAIN, Bio-Rad Laboratories S.A., Avda Dr. Severo Ochoa, Parque Casablanca Edificio B2, P3, Pol. Ind. Alcobendas, 28100 Alcobendas, Madrid • Phone 34-1-661-7085 • Telefax 34-1-661-9698
SWEDEN, Bio-Rad Laboratories A.B., Ekensbergsv., 128, Vintergatan 1, Box 1097, S-172 22, Sundbyberg • Phone 46-8-627-50-00 • Telefax 46-8-627-54-00
SWITZERLAND, Bio-Rad Laboratories AG, Kanalstrasse 9, CH-8152 Glattbrugg • Phone 41-1-809-55-55 • Telefax 41-1-809-55-00
UNITED KINGDOM, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Ave., Hemel Hempstead, Herts HP2 7TD • Phone 44-1442-232552 • Telefax 44-1442-259118