



Model GS-800[®] Calibrated Imaging Densitometer

User's Guide

For Catalog Numbers

170-7980 GS-800 Densitometer, PC

170-7981 GS-800 Densitometer, Mac

For Technical Service
Call Your Local Bio-Rad Office or
In the U.S. Call **1-800-4BIORAD**
(1-800-424-6723)

WWW.BIO-RAD.COM

Warranty Statement

Bio-Rad Model GS-800 Calibrated Imaging Densitometer

This warranty may vary outside of the United States. Please contact your local Bio-Rad office for the exact terms of your warranty.

Bio-Rad Laboratories warrants to the customer that the Model GS-800 Calibrated Imaging Densitometer (Catalog Number 170-7980 or 170-7981) will be free from defects in material and workmanship, and will meet all performance specifications for the period of one year from the date of shipment. This warranty covers all parts and labor.

In the event that the instrument must be returned for repair under warranty, the instrument must be packed for return in the original packaging.

Bio-Rad shall not be liable for any incidental, special or consequential loss, damage, or expense directly or indirectly arising from the use of the Model GS-800 Calibrated Imaging Densitometer. Bio-Rad makes no warranty whatsoever in regard to products or parts furnished by third parties, such being subject to the warranty of their respective manufacturers. Service under this warranty shall be requested by contacting the customer's nearest Bio-Rad office.

This warranty does not extend to any instruments or parts thereof that have been subject to misuse, neglect or accident, or that have been modified by anyone other than Bio-Rad or Bio-Rad's authorized agent, or that have been used in violation of Bio-Rad instructions.

For any inquiry or request for repair service, contact Bio-Rad Laboratories. Inform Bio-Rad of the model and serial number of your instrument.

Table of Contents

| | |
|---|----|
| Section 1 General Information | 6 |
| 1.1 Introduction..... | 6 |
| 1.2 Installation Qualification and Operational Qualification(IQ/OQ)..... | 6 |
| 1.3 Safety..... | 7 |
| 1.4 Regulatory Compliance | 7 |
| Section 2 Setting Up..... | 8 |
| 2.1 Checking the Contents | 8 |
| 2.2 Unpacking the Contents | 8 |
| 2.3 Unlocking the Densitometer and the Transparency Module..... | 9 |
| 2.4 Locking the Densitometer and the Transparency Module..... | 10 |
| 2.5 Taking a Closer Look..... | 10 |
| Section 3 Installation..... | 12 |
| 3.1 About SCSI Devices | 12 |
| 3.2 Changing the SCSI ID Number..... | 12 |
| 3.3 Checking the Terminators | 13 |
| 3.4 Installation for the Macintosh or PC Computer | 14 |
| 3.5 Testing the Densitometer..... | 15 |
| Section 4 Operation..... | 16 |
| 4.1 Overview of Operational Components | 16 |
| 4.2 Start-Up | 18 |
| 4.3 Reflective Scanning..... | 18 |
| 4.4 Transmittance Scanning..... | 18 |
| Section 5 Maintenance..... | 20 |
| 5.1 Cleaning the Image Window Glass..... | 20 |
| 5.2 Cleaning the Cabinet Exterior..... | 20 |
| Section 6 Troubleshooting | 21 |
| 6.1 Troubleshooting..... | 21 |
| 6.2 Technical Service..... | 23 |
| Section 7 Technical Specifications..... | 24 |

Section 1 General Information

1.1 Introduction

The Model GS-800 Calibrated Imaging Densitometer is a high performance, calibrated imaging densitometer that converts transparent and opaque electrophoretic samples into digital data. These data are analyzed with Quantity One® software operating in either Microsoft Windows® or Macintosh® computer systems. For a step-by-step guide to scanning and analyzing images, please refer to your Quantity One software user manual.

The main features of the Model GS-800 Calibrated Imaging Densitometer include:

- Adjustable resolution from 36.3 microns to 127.0 microns;
- Reflectance and transmittance
- Scanning area of 29 cm by 40 cm
- Sampling rate of up to 700 dots per inch (dpi).
- Stationary platen for higher reliability
- Variable wavelength lamp/filter architecture for improved color discrimination
- Analog to digital conversion at full 12-bit accuracy
- Calibrated transmission tablet to ensure accurate OD readings
- Reflective gray scale tablet to ensure reproducible reflectance values
- Supported by Quantity One software, which provides versatile image display, optimization, and quantitation features

1.2 Installation Qualification and Operational Qualification (IQ/OQ)

The calibration of the Model GS-800 Calibrated Imaging Densitometer can be validated by following the GS-800 Calibrated Densitometer Installation Qualification and Operational Qualification protocols. These protocols are available as an accessory to the GS-800 Calibrated Imaging Densitometer (Catalog number 170-7956). These straightforward and simple to follow IQ/OQ protocols include instructions for the installation of the hardware and software, and instructions for the verification of the reflectance and transmittance calibration functions.

Following proper installation, a linear dynamic range of 3.0 OD for the Transmissive scanning and 2.0 OD for Reflectance scanning can be confirmed with greater than 95% confidence. For additional details, contact your local Bio-Rad Representative.

1.3 Safety

The Model GS-800 Calibrated Imaging Densitometer system uses high voltage and high current and should be operated with care at all times. To avoid shock, set up the GS-800 Densitometer in a dry area. Immediately wipe up any spilled solutions.

- Turn off the Densitometer, and then disconnect the power plug when you want to clean the Densitometer case or glass plate, or when the Densitometer needs service or repair.
- Place the Densitometer on a level surface.
- To ensure proper ventilation, allow a minimum of 15 cm free space around each side of the Densitometer.
- Do not leave photographs, gels, or film on the glass plate for excessive periods of time. The heat from the light source in the Densitometer may cause them to deteriorate.
- Do not operate the Densitometer when the environmental temperature falls below 5° C or rises above 40° C.
- Do not operate the Densitometer when the environmental humidity falls below 25% or rises above 85%.

1.4 Regulatory Compliance*

Declaration of Compliance

The party responsible for product compliance

Corporate Name: UMAX Technologies, Inc.

Address: 3561 Gateway Blvd. Fremont, CA 94538, U.S.A.

Telephone No.: 510-651-4000

*See UMAX PowerLook 2100XL Operation Manual

Section 2 Setting Up

2.1 Checking the Contents

Your Model GS-800 Calibrated Imaging Densitometer system should arrive complete with one of each of the following:

- GS-800 Calibrated Imaging Densitometer
- GS-800 Calibrated Imaging Densitometer User Guide
- Quantity One™ Image Analysis Software
- UMAX PowerLook 2100XL Operation Manual
- 25/50-pin SCSI Cable
- Power Cord
- Photoperfect and Adobe Photoshop LE
- MagicScan CD and manual

These components may arrive in several boxes. If your system is missing any of these items, immediately contact your local Bio-Rad office for a prompt replacement. **Please retain all packaging material.** Additional charges will be assessed if packaging is not available for instrument warranty service shipping.

Caution: Use care when unpacking your equipment. The equipment is delicate and damage can occur if it is dropped or otherwise mistreated.

2.2 Unpacking the Contents

Unpacking the Model GS-800 Calibrated Imaging Densitometer shipping container:

Warning: The Model GS-800 Calibrated Imaging Densitometer weighs approximately 45.8 lbs. (20.8 kg). To avoid personal injury, two people may be needed to lift the Densitometer out of the shipping container. Lifting the unit just by the lid may cause damage to the instrument

1. Carefully lift the Model GS-800 Calibrated Imaging Densitometer out of the shipping carton and place it on a stable, dry, flat surface. The chassis of the Model GS-800 Calibrated Imaging Densitometer attempts to conform to any uneven surface. This may result in a poor-quality scan.
2. Remove the plastic wrapping and the packing materials from the Densitometer.
3. The Model GS-800 Calibrated Imaging Densitometer should be located close to the computer system to which it will be connected. Select a flat, dry surface that provides adequate ventilation on all sides to prevent overheating.

4. The Model GS-800 Calibrated Imaging Densitometer can operate with 100V AC to 264V AC. Check the power cord to insure that it is suitable for your area.

Caution: The power cord plug provided has a third grounding pin. This plug fits in a grounding-type outlet only. This is a safety feature. If you are unable to plug the power cord into a wall outlet, locate another outlet that can accommodate the plug or contact your electrician to change the outlet. Do not change the plug.

2.3 Unlocking the Densitometer and the Transparency Module

The optical assembly of the Densitometer's transparency module (located in the lid) is held in place during shipment by a carriage lock (Figure 2.1) on the underside of the transparency module. The carriage lock must be unlocked before the instrument will work properly. Additionally, if the carriage lock is not unlocked prior to use, the optical system may be damaged while scanning. Insert a coin into the Carriage Lock. Turn it until the mark points align with the unlocked mark.

Note: Leave the locking screw in place so that it can be re-locked if you have to move the Densitometer over long distances.

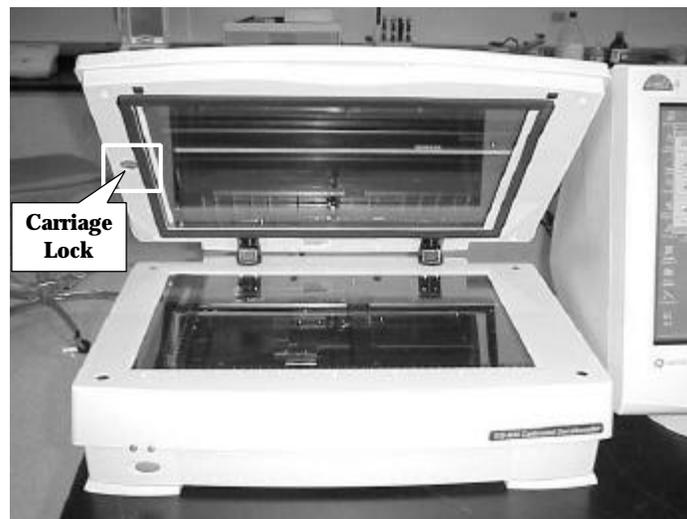


Fig. 2.1 Front view of Densitometer. Unlock the carriage lock by turning with a coin.

Warning: Always lower the transparency module slowly and release it only when it lies flat on the body of the Densitometer.

2.4 Locking the Densitometer and the Transparency Module

To transport or ship the Densitometer, be sure to re-lock the optical assembly to avoid damage to it. Before re-locking the assembly, make sure the optical assembly of the transparency adapter is in the Home Position by doing the following:

1. If the Densitometer is powered on, turn the Densitometer off.
2. Turn the Densitometer's power back on.
3. Wait for the Densitometer's Ready and Power indicators to light up.
4. Turn the Densitometer off again. The transparency assembly should be in the Home position.

To lock the optical assembly, insert a coin in the Carriage Lock. Turn it until the marked points align with the locked marks.

2.5 Taking a Closer Look

Now that you have the Model GS-800 Calibrated Imaging Densitometer out of the box, take a closer look to familiarize yourself with the parts. Figures 2.2 and 2.3 illustrate the locations of the different parts of your Model GS-800 Calibrated Imaging Densitometer.

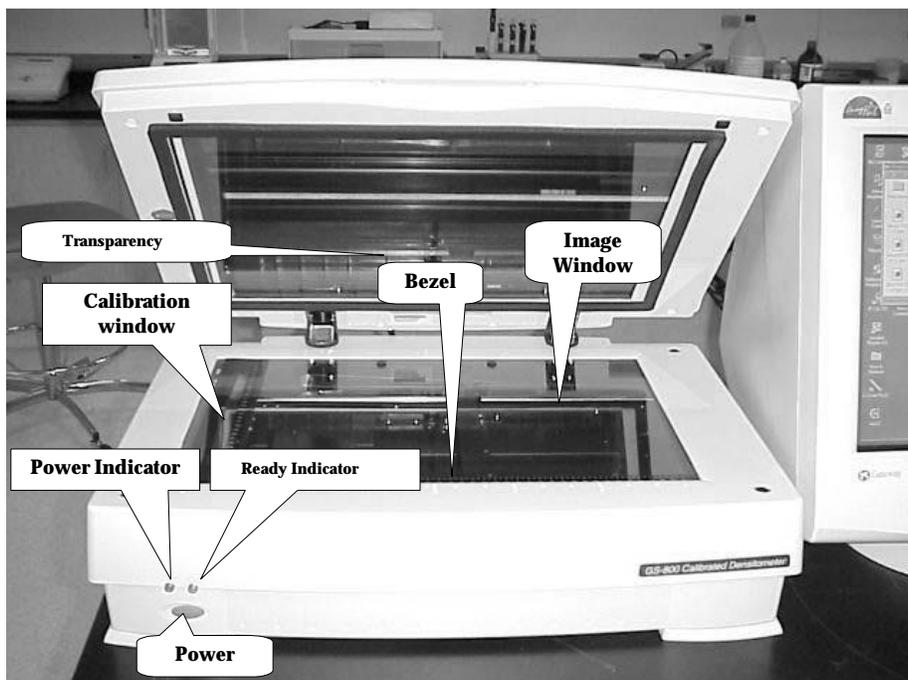


Fig. 2.2 The front of Model GS-800 Calibrated Imaging Densitometer.

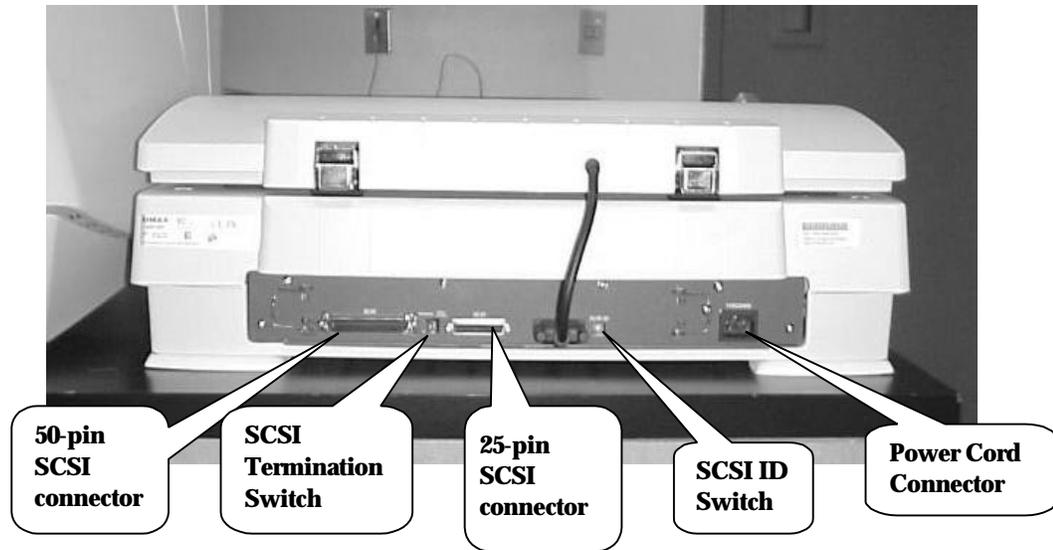


Fig. 2.3 The back of the Model GS-800 Calibrated Imaging Densitometer.

Section 3 Installation

3.1 About SCSI Devices

The Model GS-800 Calibrated Imaging Densitometer is a Small Computer System Interface (SCSI) device. It communicates with your computer by using the SCSI-2 standard. The SCSI communication standard allows you to connect more than one peripheral device to the same port of your computer in chain fashion.

A unique SCSI ID number is assigned to each device in the SCSI chain enabling your computer to identify the device with which it is communicating and the priority of each device.

Warning: If two SCSI devices have the same ID number, your system will not work properly and you may damage your SCSI devices.

A SCSI chain requires an electronic component called a 'terminator' which absorbs old signals traveling along the cables and keeps the path open for new signals. The chain should never have more than two terminators, one at each end. It is important to remember that using too many or too few terminators may damage your SCSI devices. Some SCSI devices have built-in terminators and must therefore be placed at the beginning or end of your SCSI chain.

Note: The Model GS-800 Calibrated Imaging Densitometer has a built-in terminator.

3.2 Changing the SCSI ID Number

Your Densitometer's SCSI ID setting is factory preset at #6. Check to see if this ID setting is used by another device connected to your computer's SCSI port.

If SCSI ID #6 is not used, you do not need to change your Densitometer's SCSI ID number. You can directly proceed to hardware connection and software installation. For installation instructions, proceed to Section 3.4 if you are connecting to a Macintosh computer or to Section 3.5 if you are connecting to a PC.

If you find however, that another connected device is already using SCSI ID #6, then you must reset the SCSI ID on your Densitometer.

To reset the SCSI ID, do the following:

1. Make sure the Densitometer power is off.
2. Use a screwdriver to rotate the notch on the SCSI ID switch until an arrowhead points to an unused number.

Note: Do not use SCSI ID setting 7 through 9 on your Densitometer. They are for factory use only.

3.3 Checking the Terminators

There should be two terminators in a SCSI chain. It is best to place the terminators at each end of the SCSI chain. The GS-800 Calibrated Densitometer has a built-in active terminator switch for you to turn on and off. Number 1 represents **active terminator** and number 2 represents **terminator power**. Turning both the **active terminator** and **terminator power** switch on by pressing them down will enable the built-in “active terminator” function.

The simplest configurations for using the Densitometer with a SCSI adapter card are as follows:

Case 1: Connect the Densitometer to an interface card that does not have another SCSI device attached to it. e. g. Densitometer is the **only** SCSI device connected to your computer.

*In this situation, the card will have a built-in terminator. The card forms one end of the SCSI chain, the Densitometer forms the other end. Turn **on** the **active terminator** and **terminator power** switch by pressing them down to enable the built-in “active terminator” function.*

Case 2: Connect the Densitometer to a SCSI card that has another SCSI device attached to it.

*In this case, turn **off** the **active terminator** on your Densitometer. If the SCSI device next to Densitometer is the last SCSI device of the SCSI chain, please refer to the instruction manual of this device to set its terminator on. If you experience unreliable SCSI operation and suspect terminator problems, contact Bio-Rad Technical Service.*

Case 3: Densitometer is the **last** SCSI device in the SCSI Chain

The setting of active terminator is the same with Case 1. Simply turning on the active terminator of your Densitometer.

3.4 Installation for the Macintosh or PC Computer

This section describes setting up the Model GS-800 Calibrated Imaging Densitometer with your Macintosh or PC computer. You must first choose and set a SCSI ID number (Section 3.2), then connect the Densitometer to your computer, and lastly test the connection.

SCSI Adapter Card

The Model GS-800 Calibrated Imaging Densitometer requires a SCSI adapter card to work with your computer. The Densitometer driver software supplied with the Densitometer supports most Adaptec adapter cards.

Please check the following documentation:

- The Read Me file on the Quantity One software disk for up-to-date information.
- The documentation supplied with your Adaptec SCSI adapter card for instructions on card installation.

SCSI Adapter Card Installation

Before you can use your Densitometer with your computer, you need to install the interface card into your computer. Contact your local Bio-Rad Sales Representative or Technical Service for the recommended Adaptec SCSI adapter. To install the card into one of the computer's expansion slots, observe the following procedures:

Note: The interface card is sensitive to static electricity. Handle the card by its mounting bracket, particularly when removing the card from its anti-static packaging.

1. Turn the computer power off and unplug the power cord.
2. Remove the housing cover of the computer. Follow the instructions provided in your computer's reference manual.
3. Remove the metal cover corresponding to your chosen slot. Keep the removed screw so that it can be used to fasten the interface card.
4. Gently insert the interface card into the slot until it is firmly seated in the slot.
5. Secure the card in place with the screw removed from the expansion slot cover in step 3 above.
6. Replace the housing cover following the instructions provided in the computer's reference manual.

Connect the Densitometer to the Computer

With settings on the Densitometer and card correctly set and the interface card properly installed in your computer, you can now connect the Densitometer and the computer, as follows:

1. Ensure that the Densitometer's SCSI ID is properly set. Determine which SCSI ID numbers are already assigned and which numbers are free. Refer to "Changing the SCSI ID Number" in Section 3.2 for instructions on setting the Densitometer's SCSI ID.
2. Turn on Active terminator and termination power (please refer to Section 3.3 for more details on setting termination), connect the SCSI cable to SCSI port and connect the other end of the SCSI cable to the Densitometer.
3. Connect the power cord to the Densitometer.
4. Turn on the Densitometer power.
5. Turn on the computer power.

3.5 Testing the Densitometer

Prepare your Densitometer for installation through the following steps:

1. Check and reset (if necessary) the Densitometer's SCSI ID
2. Run the automatic Densitometer self-test (see below)

Test the Densitometer

The Densitometer automatically performs a simple self-test each time it is turned on. The self-test checks the status of certain Densitometer devices.

Start the Densitometer self-test by following the steps below:

1. Connect the power cord to the Densitometer.
2. Connect the other end of the power cord to a wall outlet.
3. Turn on the power of the Densitometer. The on/off control is a push-button switch located on the left front of the unit.

At power-on, the front panel indicators flash once. The power indicator then glows and the ready indicator blinks. When the test is completed, the power and the ready indicators glow steadily.

The Densitometer performs a self-test for approximately 35 seconds after which the Densitometer indicator stabilizes. The Transparency Module indicator and the scanning lamp stay on. If a problem occurs during the self-test, refer to Section 6, Troubleshooting.

Section 4 Operation

4.1 Overview of Operational Components

This chapter provides a summary of the Model GS-800 Calibrated Imaging Densitometer components used during operation and explains the startup procedure.

Transparency Module

The Transparency Module (Figure 2.2) is secured to the Model GS-800 Calibrated Imaging Densitometer by two elongated hinges at the rear of the Densitometer. This Transparency Module should be lowered onto the Densitometer base for reflectance and transmittance scanning. The thick border gasket on the transparency module allows you to place samples up to 3.0 mm thick on the image window for scanning.

Warning: Always lower the transparency module slowly and release it only when it lies flat on the body of the Densitometer.

Image Window

The image window (Figure 4.2) is the area where the sample to be scanned is positioned. Light collected from the sample is focused through a lens onto a charge coupled device (CCD) array that captures the image.

Caution: The image window is glass. To avoid breakage, do not press firmly or place anything other than a sample for scanning on the image window.

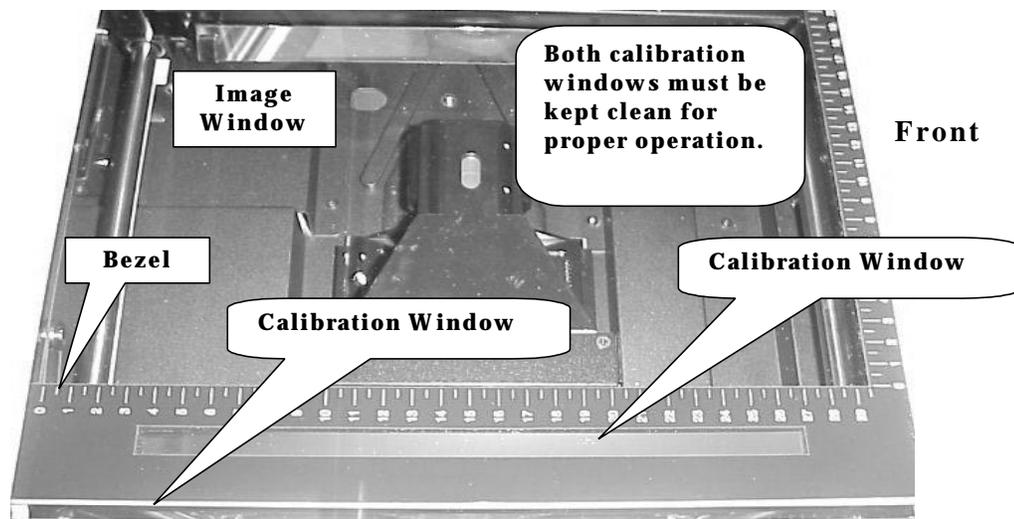


Figure 4.2 The image window and bezel on the Densitometer

Calibration Tablet Areas/Calibration Windows

The GS-800 Densitometer is equipped with reflectance and transmission step wedges. The transmission step wedge is located on the left hand side of the image window. The reflectance step wedge is internal. When you scan a transparent sample, place the sample on the Densitometer's image window Do not cover the calibration windows (Figure 4.2) or a calibration error will occur.

Validation and Documentation the calibration performance can be accomplished by the following GS-800 IQ/OQ Protocols.

See section 1.2

Bezel

The bezel (Figure 4.2) is the raised frame around the image window. It has marks in inches and centimeters to help you correctly position the sample on the image window for scanning.

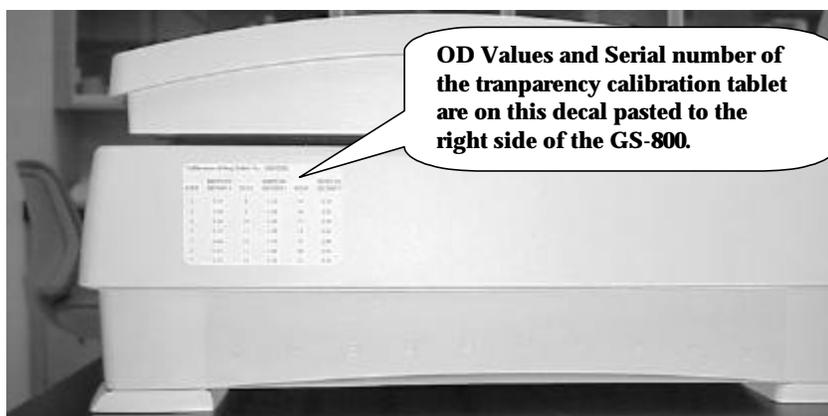
Indicator Lights

The indicator lights (Figure 2.2) are located on the front left side of the Model GS-800 Calibrated Imaging Densitometer cabinet. This panel consists of two lights that indicate the status of the unit.

| Light | Description |
|-----------------|--|
| Power Indicator | Is lit when the power switch is in the on position (green light). |
| Ready Indicator | Flashes during self-test(start-up) and during scans (green light). |

Calibration Strip Values

The OD values of the built in Calibration Strip can be found either on the decal posted on the side of the GS-800 (see below) or on the enclosed OD value sheet .



4.2 Start-Up

4.2 Start-Up

Turn on the Densitometer. The on/off control is a push-button switch located on the left front of the unit. The Power indicator lights up. The Densitometer performs a self-test that lasts approximately 35 seconds during which the Ready indicator flashes briefly. The scanning lamp turns on and both indicator lights will stay on once the Densitometer is powered up.

For optimum performance, allow the Densitometer to warm up for five minutes after power-up.

The Model GS-800 Calibrated Imaging Densitometer is now ready for use.

4.3 Reflective Scanning

To scan a reflective sample:

1. Raise the Transparency Module.
2. Place the sample to be scanned facedown on the image window.
3. Guide the top corner of the sample to the lower left (0,29) corner of the image window.
4. Lower the Transparency Module

Warning: Always lower the transparency module slowly and release it only when it lies flat on the body of the Densitometer.

5. For more information about scanning operations and options, refer to the Quantity One software manual.

Note: The Reflective density values for the Reflectance Calibration strip are already included as default values in the software so you do not need to enter them to calibrate the instrument. Keep the image window clean and free of oil and fingerprints. See Section 5, Maintenance, for cleaning procedures.

4.4 Transmittance Scanning

To scan a transparent sample:

1. Raise the Transparency Module.
2. Place the sample face down on the Densitometer's image window.
3. Lower the Transparency module.

Warning: Always lower the transparency module slowly and release it only when it lies flat on the body of the Densitometer.

4. For more information about scanning operations and options, refer to the Quantity One software manual.

Note: The OD values for the Transparency step tablet are included on a separate sheet with this Manual and are also on the Decal pasted on the right side of the instrument. If for any reason these values are lost contact Bio-Rad Technical Service. Keep the image window clean and free of oil and fingerprints. See Section 5, Maintenance, for cleaning procedures.

Section 5 Maintenance

The Model GS-800 Calibrated Imaging Densitometer requires easy periodic cleaning for continued optimum performance.

Warning: To prevent personal injury, always turn the power switch off and unplug the power cord from the wall outlet before performing any maintenance on the Model GS-800 Imaging Densitometer.

Caution: Do not use ammonia-based cleaning products to clean the Model GS-800 Imaging Densitometer. Do not disassemble the Model GS-800 Calibrated Imaging Densitometer or lubricate any parts.

5.1 Cleaning the Image Window Glass

To clean the glass:

1. Turn off the power switch.
2. Unplug the power cord from the wall outlet.
3. Use a slightly damp cloth with a small amount of mild detergent or alcohol to wipe the glass.
4. Dry the glass with a clean, lint-free cloth.

Caution: To avoid breakage, do not rub heavily on the glass. Do not spray liquids directly onto the image window. If spots remain on the glass, carefully use a razor blade to remove them.

5.2 Cleaning the Cabinet Exterior

To clean the cabinet exterior:

- 1 Turn off the power switch.
- 2 Unplug the power cord from the wall outlet.
- 3 Use a slightly damp cloth to wipe the exterior.
- 4 Dry the cabinet with a clean, lint-free cloth.

Caution: If the cabinet is stained and a damp cloth does not remove the stain, a mild detergent solution can be used.

Section 6 Troubleshooting

Though the Model GS-800 Calibrated Imaging Densitometer is very reliable, problems can occur during operation. This section lists some problems and the recommended corrective actions. Most problems are easily remedied.

6.1 Troubleshooting

Electronic Failure

Call Bio-Rad Technical Service.

Will not scan again, stays in home position, or grinds.

Make sure that the shipping lock down screw for the Transparency module has been properly unlocked (See Section 2.3). If this still happens then it is probably related to a mechanical failure somewhere inside the Densitometer. A part may have come loose and is obstructing the path of the lamp assembly. Contact Bio-Rad Technical Service.

The Power indicator fails to light up.

1. Verify the power connection to the Densitometer.
2. Check that the power switch is turned on.
3. If you have confirmed that there is power to the Densitometer, it is likely that the Densitometer fuse needs to be replaced. Contact Bio-Rad Technical Service.

The scanned image is marred by specks or horizontal lines.

Hardware failure. Contact Bio-Rad Technical Service.

The scanned image is showing random dark specks or vertical L bands.

Clean the image window glass. Refer to Section 5.1 on Maintenance.

The scanned image is marred by Moiré patterns

Moiré patterns are often seen when scanning a halftone. A halftone is a picture in which the gradations of light are simulated by dots of varying sizes or densities and produced by superimposing a screen over an image (e.g., pictures in newspapers and magazines). It is best to avoid scanning halftones because of the moiré, but if it is unavoidable, scan at a lower resolution.

The scanned image is too dark.

1. If you are scanning a transparent sample, make sure you are using the transparency selection.
2. If the problem persists, contact Bio-Rad Technical Service.

The computer does not boot.

Your computer cannot find its hard disk due to a conflict with the SCSI ID numbers of the devices you have attached. Turn off the power to the Densitometer and unplug the cord.

Disconnect all SCSI devices and connect them one by one, beginning with the Densitometer, to identify the device that causes the problem.

The Quantity One software returns a "Calibration error " or similar message when attempting to calibrate.

1. Check if the OD values for the Transparency target in the EDIT CALIBRATION window are entered properly.
2. Make sure the calibration windows are clean and free of dust, dirt, fluid, and other contaminants that can obstruct the path of light.
3. Check for correct offset value. For the GS-800 Calibrated Imaging Densitometer the correct offset value is 00 for transmissive scanning and 0.05 for reflectance scanning.
4. Make sure that the transparency scan image shows up correctly at the bottom of the Quantity One acquisition window and that during the calibration the software identifies each step wedge with a number inside a yellow rectangle in the sequence shown in the photo below:



The Quantity One software returns a "Densitometer not ready" or similar message.

1. Verify the power connection to the Densitometer.
2. Check that the power switch is turned on.
3. Check the Installation procedure (Section 3), to see if you followed the instructions. Pay special attention to the setting of the SCSI ID number.
4. Verify the integrity of the terminators and the cables. If there is a problem with a cable or with the SCSI board of the Densitometer, contact Bio-Rad Technical Service.
5. Disconnect all SCSI devices and connect them one by one, beginning with the Densitometer, to identify the device that causes the problem.
6. Verify that the Densitometer imaging device is selected in the FILE menu of the Quantity One software.

ASPI error message appears.

Call Bio-Rad Technical Service for assistance.

6.2 Technical Service

Bio-Rad provides complete technical support on all features of the GS-800 Calibrated Imaging Densitometer system, including all hardware, software, or spare parts issues. If you have any problems or any questions on the use of the features, contact your local Bio-Rad office, or in the U.S. call 1-800-424-6723.

Section 7 Technical Specifications

| | |
|-------------------------|---|
| Color Separation | 3 linear CCDs and 3 color (RGB) filter layers in conjunction with fluorescent lamps. |
| Scanning Resolution | Variable resolution from 200 to 700 dpi |
| Gray Scales | 12-bit A/D conversion |
| Effective Scanning Area | Reflective Mode 29 cm x 40 cm (8.5 in. x 14 in.) Transmittance Mode 29 cm x 40 cm (8 in. x 10 in.) |
| Interface | Built-in SCSI II, one 25-pin and one 50-pin connector |
| Transmission Speed | 1 Mbyte/second |
| Lamp | 8W daylight fluorescent light |
| Warm up time | 35 seconds, 5 minute for stable operation |
| Voltage | Line voltage: 100 to 240 Volts Frequency: 50 to 60 Hz Power: 30 Watts maximum |
| Dimensions | Length: 47.3cm (18.6 in.) Width: 62.6 cm (24.6 in.) Height: 19.3 cm (7.6 in.) |
| Weight | Approximately 20.8 kg (45.8 lb.) |
| Environment | Operating temperature 5°C to 40° C (40° F to 104° F) Relative Humidity 20% to 85% Surrounding space 15 cm on each side |

GS-800 and Quantity One are trademarks of Bio-Rad Laboratories. Macintosh is a trademark of Apple Computer, Inc. Windows is a trademark of Microsoft Corporation.



**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 **China** (86-21) 63052255 **Denmark** 45 44 52-1000 **Finland** 358 (0)9 804 2200 **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** 64-9-4152280
Norway 47-23-38-41-30 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-2729877 **South Africa** 00 27 11 4428508 **Spain** 590 5200
Sweden 46 (0)8-55 51 27 00 **Switzerland** 061 717-9555 **United Kingdom** 0800-181134
