

# The **Discovery Series**<sup>™</sup>

**Quantity One**<sup>®</sup> 1-D Analysis Software

**PDQuest**<sup>™</sup> 2-D Analysis Software

**BIO-RAD**

# Simply the Best Software for Gel Analysis

**Bio-Rad Laboratories is the world's largest supplier of imaging products for life science research. Our unmatched expertise in gel electrophoresis is the driving force behind our success with imaging instruments and image analysis software. Bio-Rad is proud to present a range of software packages for gel and blot analysis in The Discovery Series software.**

**The Discovery Series software has a long tradition of reliability in gel imaging and analysis. For more than 15 years, scientists have used Quantity One 1-D analysis software and PDQuest 2-D analysis software to evaluate and interpret the results of 1-D and 2-D electrophoretic separation of DNA and proteins in agarose and polyacrylamide gels.**

## Research tools evolve —

biomolecule separation techniques improve, the capacity and speed of computers increases, imaging technologies advance, and information technology tools become available. The Discovery Series software has been continually developed to meet the evolving requirements of the modern research lab. Bio-Rad has a proven track record of adding new, sophisticated analysis and automation features, and updating our imaging software for compatibility with the latest computer technology and operating systems.

As a leader in imaging, Bio-Rad will be updating and supporting The Discovery Series software for many years to come.

## Integrated Image Analysis Systems

From gel documentation to isotopic and chemiluminescent analysis, Bio-Rad brings you the most advanced technologies and the largest range of innovative imaging systems to match your needs. Patent-pending technologies guarantee clean and beautiful images from high-resolution and quantitative imaging systems.

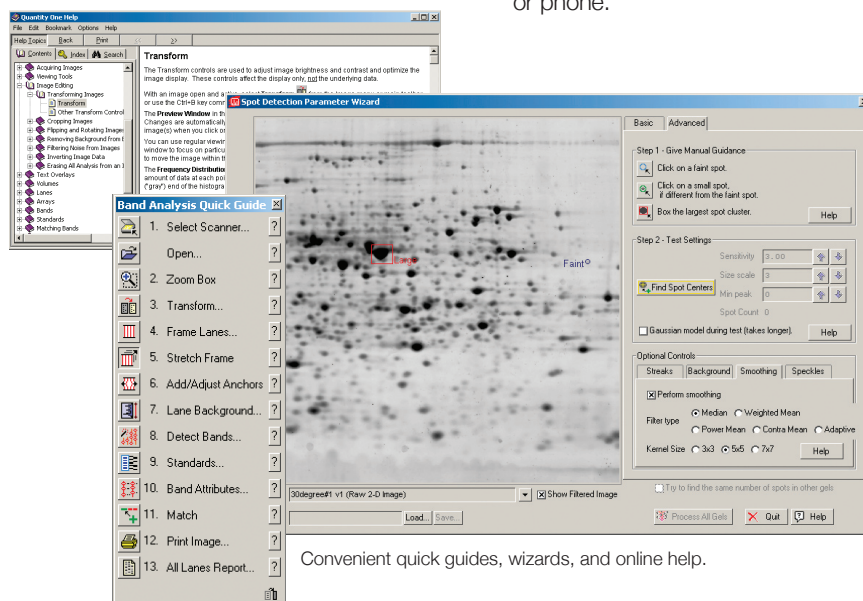
Integrated image analysis systems that combine image capture and analysis eliminate the expense and training time needed for other complex, application-specific software packages. The Discovery Series software interface is

application based, and it automatically configures your imaging system with the appropriate filters, lasers, etc., for your research sample. Our integrated systems take you rapidly from a sample to an optimized image and printout.

You can connect one or more of Bio-Rad's imaging systems to a single PC or Macintosh computer to control the scanning and analysis of all of your laboratory's samples. The Discovery Series software can also be easily networked for rapid data acquisition and flexible analysis at multiple remote stations. This is useful for multiuser core facilities, large institutions, and collaborating laboratories.

## Ease of Use

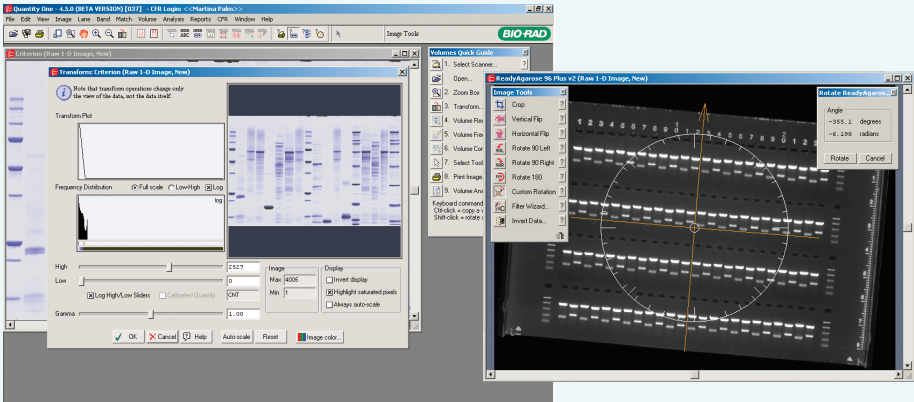
Simple built-in tools and quick guides let you start immediately. The Discovery Series offers an easy-to-use graphical interface with familiar pull-down menus, toolbars, and keyboard commands. File formats and menu commands are shared across Quantity One and PDQuest, allowing you to switch easily between 1-D and 2-D analysis. Onscreen, context-sensitive help is available at your fingertips. Still have a question? Contact Bio-Rad's knowledgeable team of technical support representatives via e-mail or phone.



Convenient quick guides, wizards, and online help.

# The Perfect Image

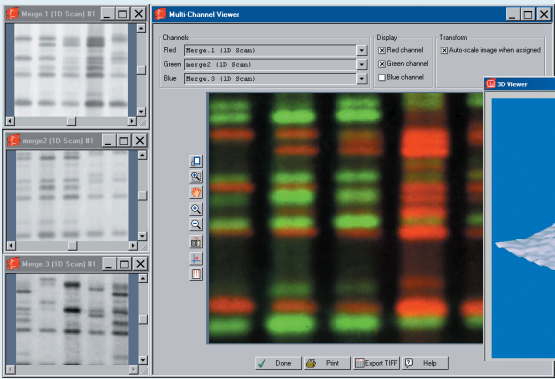
The Discovery Series includes tools to help you get perfect images. Publication-quality images are easily generated with a comprehensive range of image optimization tools. You can rapidly crop your images, filter noise, adjust brightness, improve contrast, and rotate skewed images incrementally. The color palette allows realistic color representation of your Coomassie-, silver-, or fluorescently stained samples.



Improving contrast and rotating images to capture the perfect image.

# Visualizing Data

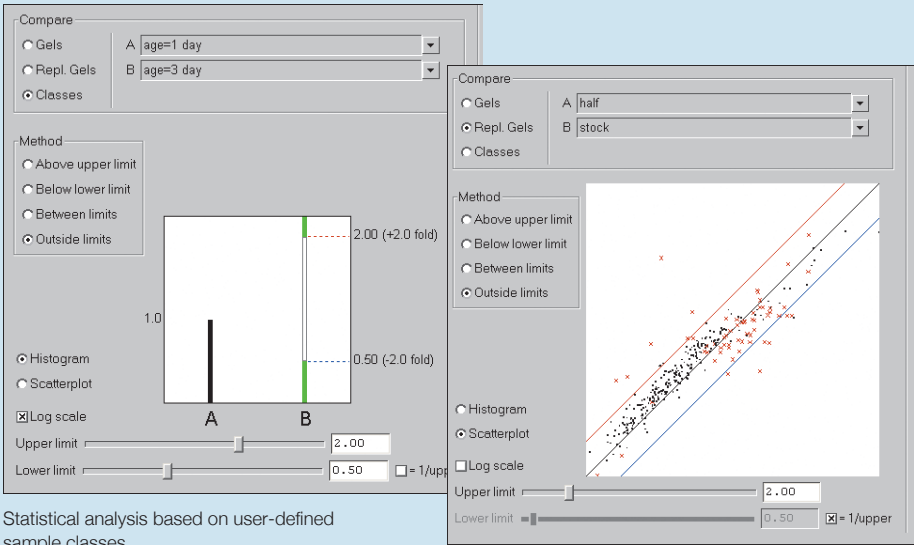
In addition to standard graphs and data reports, The Discovery Series offers sophisticated ways to visualize your data. To confirm that your data are interpreted correctly, use 3D Viewer for 3-D modeling of any user-defined area of the gel, and the Image Stacking tool for aligned viewing of multiple gel images. To conveniently view multicolor data, the Multichannel Viewer allows merging of three independent color channels into one image.



Merging and viewing up to three images in different color channels.

# Sophisticated Analysis

Whether you are quantitating a DNA gel, looking for the absence or presence of a band, determining purity of a protein sample, calculating molecular weight on bands in an unknown sample, or analyzing protein expression, The Discovery Series offers comprehensive and sophisticated analysis for 1-D and 2-D gels and blots. You will also find tools for colony counting and analysis of arrays and microplates.



Statistical analysis based on user-defined sample classes.

## Annotations

The Discovery Series features versatile annotation tools to add text comments, arrows, and lines to your images. For publication-quality images, annotations can easily be configured with the desired properties (font, type size, and color) and automatically aligned.

## Data Security

Regulated organizations where users require data security and integrity can add an optional security module that provides features for compliance with US Food and Drug Administration 21 CFR Part 11 regulations. The Discovery Series software confirms the accuracy of electronic records using a secure checksum. Secure time-stamped audit trails are implemented to record the date and time of operator actions that create, modify, or delete electronic records. To allow previewing and approval of data files prior to further processing, electronic signature capability is implemented.

## Sharing Data

Many features of The Discovery Series are available free in the basic mode of the software. You can share images with colleagues and collaborators and submit TIFF or JPEG images to journals for publication. The Discovery Series software can be loaded onto an unlimited number of computers, allowing everyone in your lab free access to the basic functionality.

## Software Updates

Bio-Rad continually updates the software with new functionalities. The latest software and user guides can be downloaded from the Web. Check for new releases at **www.bio-rad.com/softwaredownloads/**

### Ordering Information

Catalog #	Description
170-9650	The Discovery Series CD-ROM
<b>Quantity One 1-D Analysis Software</b>	
170-9600	The Discovery Series Quantity One 1-D Analysis Software
170-9601	The Discovery Series Quantity One 1-User Network License
170-9602	The Discovery Series Quantity One 2-User Network License
170-9603	The Discovery Series Quantity One 3-User Network License
170-9604	The Discovery Series Quantity One 4-User Network License
170-9605	The Discovery Series Quantity One 5-User Network License
170-9606	The Discovery Series Quantity One 10-User Network License
170-9607	The Discovery Series Quantity One 20-User Network License
170-9608	The Discovery Series Quantity One Add 1 User to Network License
170-9610	The Discovery Series Quantity One Version Upgrade
170-9612	The Discovery Series Quantity One User Guide
170-9615	The Discovery Series Quantity One CFR Module
<b>PDQuest 2-D Analysis Software</b>	
170-9630	The Discovery Series PDQuest 2-D Analysis Software
170-9631	The Discovery Series PDQuest 2-D 1-User Network License
170-9632	The Discovery Series PDQuest 2-D 2-User Network License
170-9633	The Discovery Series PDQuest 2-D 3-User Network License
170-9634	The Discovery Series PDQuest 2-D 4-User Network License
170-9635	The Discovery Series PDQuest 2-D 5-User Network License
170-9636	The Discovery Series PDQuest 2-D 10-User Network License
170-9637	The Discovery Series PDQuest 2-D 20-User Network License
170-9638	The Discovery Series PDQuest 2-D Add 1 User to Network License
170-9640	The Discovery Series PDQuest 2-D Version Upgrade
170-9642	The Discovery Series PDQuest 2-D User Guide
170-9645	The Discovery Series PDQuest 2-D CFR Module

Coomassie is a trademark of BASF Aktiengesellschaft.

Information in this tech note was current as of the date of the writing (2008) and not necessarily the date this version (Rev A, 2008) was published.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400  
**Canada** 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65  
**Germany** 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **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** 0508 805 500  
**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** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000