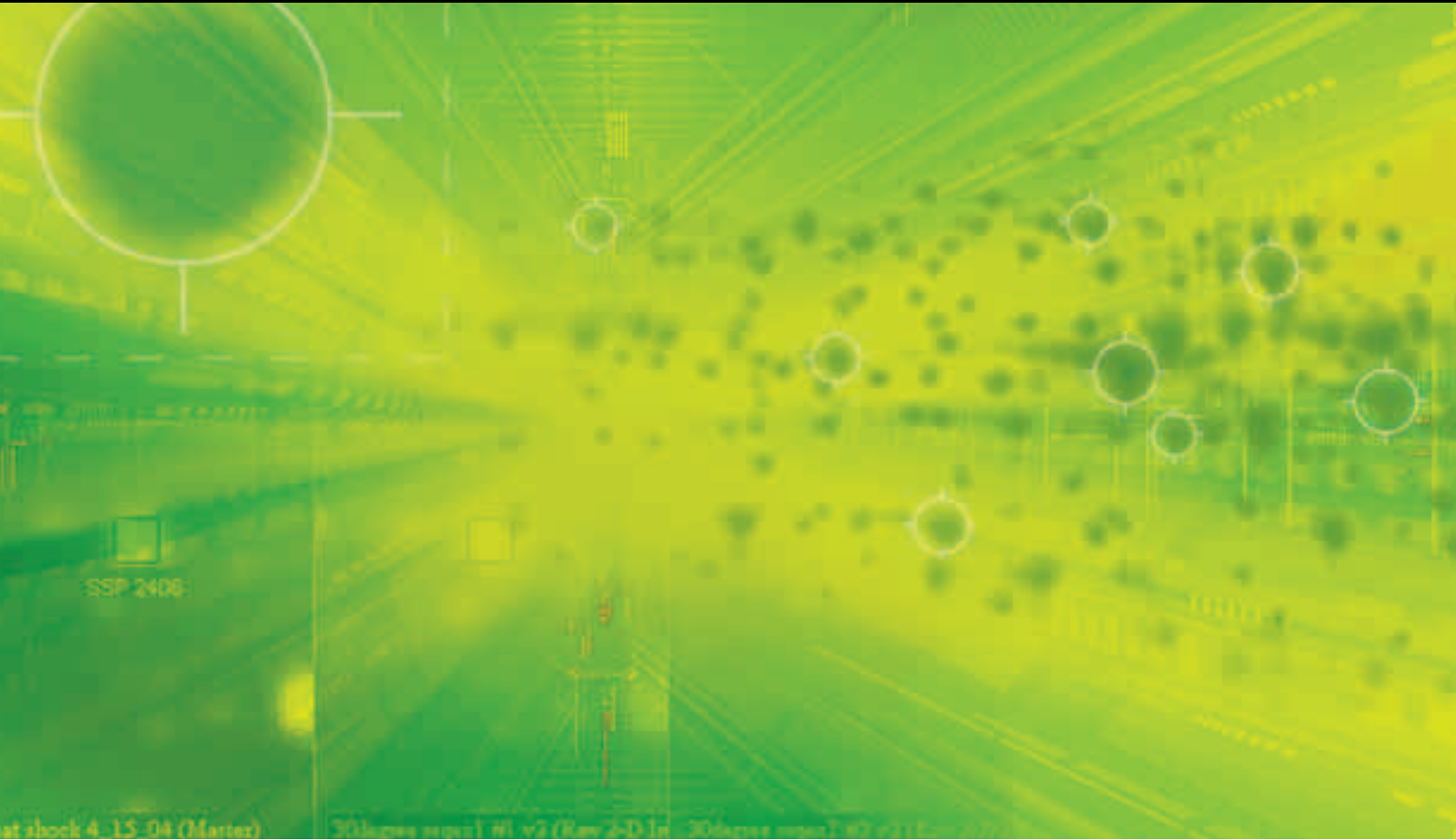


EXQuest™ Spot Cutter



Automated Gel Excision
Expertise for Every Lab



Bio-Rad Expression Proteomics

Bio-Rad's approach to expression proteomics focuses on four technologies: sample preparation, two-dimensional (2-D) electrophoresis, imaging and analysis, and protein identification. Each technology area is ideally suited to a 2-D gel electrophoresis approach for analysis of sample proteins. By providing choices in methodology, protocols, and products, Bio-Rad's expression proteomics tools help you achieve optimal results in any 2-D experiment.

Integrated Workflow Expertise

The EXQuest spot cutter is the latest addition to the expression proteomics suite of products designed by our team of experts. EXQuest provides the accuracy and precision required for reproducible mass spectrometric protein identification after gel-based separation. Continue to discover with confidence — delivered by Bio-Rad's 40+ years of experience developing research tools. EXQuest provides EXcision EXPertise for every laboratory.

Cut With Confidence

Precision, Accuracy, and Speed in Gel Excision

Gel Excision for Every Lab

Studying complex biological relationships requires reproducible methods and tools for purifying, separating, imaging, analyzing, and excising interesting proteins before they can be identified through mass spectroscopy.

The EXQuest spot cutter offers fast, hands-free gel excision with unparalleled accuracy and reliability. Bio-Rad Laboratories has reinvented gel excision with convenience and precision in mind, and has made it accessible to all types of laboratories. Whether you require occasional gel cutting or high-throughput gel cutting on a daily basis, EXQuest will improve your laboratory workflow and increase the accuracy of your results.



Expression Proteomics Workflow

● Sample preparation

● Electrophoresis

● Imaging and analysis

Science

Intelligent Workflow

- Gel excision is integrated with gel analysis software tools, allowing easy transition from gel analysis to gel cutting
- Gel image is matched to the original analysis image on the day of gel cutting, to increase precision by accounting for tearing, shrinking, swelling, or fading of protein spots
- Easily review spots to be cut, with real-time imaging of gel on cutter alongside original analysis image
- A variety of editing tools allow you to create multiple cuts on large spots (automatically or manually), move cuts, select different gels within a set, or pool cuts from multiple gels

Hands-Free Multi-Gel Imaging and Plate Processing

- Imaging and cutting of up to 4 Criterion™ gels at a time
- High-throughput capacity of up to 4 microplates

Sophisticated Liquid Handling

- Vacuum pickup and liquid-facilitated delivery for the most accurate gel recovery available, >99.5%
- Automatic calculation of dispensing volumes based on gel thickness and well volume
- Automated gel hydration to ensure gel integrity
- Multiple wash/rinse stations with “flow-through” tip rinsing for the most accurate spot identifications



Auto alignment of analysis set and cutter images.

Lane volume excision tool.

High-Speed Precision Robotics

- “New-generation” Cartesian robotics for reliable delivery of up to 600 spots per hour
- Resolution of 100 µm for unbeatable precision
- Smooth operation for any benchtop

Application and Format Flexibility

- Visible light- or UV-based applications, including SYPRO Ruby and Orange, silver, and Coomassie Blue
- Flexible spot placement options, including single or multiple spots per well in 384-well, 96-well, or 96-tube plates
- Cut gels from 0.5 to 2.0 mm thick of any gel format, including mini, Criterion, and large format (24 cm)
- Automated excision of both 1-D and 2-D gels
- Optional membrane cutting tool for cutting blots
- Excision from backed or free gels
- Import cut lists from any 2-D software package



Analysis set excision control.



Analysis set image and cutter image alignment setup.

Specify excision options.

● Protein identification

Spot cutting

● Image gel on cutter

● Align gel to original analysis image

● Review cut list

● Cut gels

● ID proteins



Versatile applications for UV- and fluorescently stained gels.

Ordering Information

Catalog # Description

EXQuest Spot Cutter

165-7200 EXQuest Spot Cutter
165-7201 EXQuest Spot Cutter With PC

Accessories

165-7202 Cutting Tip, 1.0 mm
165-7203 Cutting Tip, 1.5 mm
165-7204 Glass Bottle, 1 L
165-7205 Calibration Pucks, 10
165-7206 Membrane Cutting Head, with 1.0 mm tip
165-7207 Membrane Cutting Tip, 1.0 mm
165-7208 Gel Cutting Sheets, 15
165-7209 Gel Holding Clips, 2
165-7210 Calibration Target
165-7212 Micro Tubes, 1.5 ml, 20
165-7214 Bottle Holder
165-7215 Gel Tray
165-7216 Transilluminator Lamp
165-7217 Round-Bottomed Microplates, 96-well, 20
165-7218 Ferrule, 10–32, 1/16" OD, 10
165-7219 Bar Code Reader
165-7220 Microplate Holder

Related Products

Catalog # Description

PDQuest™ 2-D Analysis Software

170-9630 PDQuest 2-D Analysis Software
170-9640 PDQuest Version Upgrade
170-9645 PDQuest CFR Module
170-9642 PDQuest User Guide
170-9631 PDQuest 1-User Network License
170-9632 PDQuest 2-User Network License
170-9633 PDQuest 3-User Network License
170-9634 PDQuest 4-User Network License
170-9635 PDQuest 5-User Network License
170-9636 PDQuest 10-User Network License
170-9637 PDQuest 20-User Network License
170-9638 PDQuest Add 1 User to Network License

Quantity One® 1-D Analysis Software

170-9600 Quantity One 1-D Analysis Software
170-9610 Quantity One Version Upgrade
170-9615 Quantity One CFR Module
170-9612 Quantity One User Guide
170-9601 Quantity One 1-User Network License
170-9602 Quantity One 2-User Network License
170-9603 Quantity One 3-User Network License
170-9604 Quantity One 4-User Network License
170-9605 Quantity One 5-User Network License
170-9606 Quantity One 10-User Network License
170-9607 Quantity One 20-User Network License
170-9608 Quantity One Add 1 User to Network License

Coomassie is a trademark of BASF Aktiengesellschaft.
SYPRO is a trademark of Molecular Probes, Inc.

BIO-RAD

**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 2527 3454
Canada (905) 712-2771 China (86 21) 6426 0808 Czech Republic + 420 2 41 43 05 32 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)-2398112/3/4, 5018111, 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 55-52-00-05-20 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 7 095 721 1404 Singapore 65-64153188 South Africa 00 27 11 4428508
Spain 34 91 590 52 00 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan (886 2) 2578 7189/2578 7241 United Kingdom 020 8328 2000