



# The **Proteome Works™** Spot Cutter Fluorescent Enclosure

Highly accurate  
spot cutting from  
fluorescent and visibly  
stained protein gels  
and membranes

The  
**Proteome**WORKS  
System™

[proteomeworkssystem.com](http://proteomeworkssystem.com)

# The ProteomeWorks<sup>™</sup> Spot Cutter with Fluorescent Enclosure

The ProteomeWorks spot cutter with fluorescent enclosure provides automated spot cutting with expanded capabilities and versatility. Simply adjust a few settings and the spot cutter is ready to cut spots from gels or membranes stained with fluorescent SYPRO Ruby protein stains, as well as with traditional visible stains like Coomassie Blue and silver. Switching modes is nearly effortless and requires no recalibration, so lab workflow is unaffected by frequent imaging changes.

The ProteomeWorks spot cutter with fluorescent enclosure offers:

- Highly accurate spot cutting from gels or membranes stained with visible or fluorescent SYPRO Ruby protein stain
- Quick-change settings to switch between UV and visible light modes
- A small bench footprint (57 x 73 x 72 cm) that fits into any proteome lab
- Compatibility with existing ProteomeWorks spot cutters already in use worldwide

## Flexibility without Compromise

The ProteomeWorks spot cutter with fluorescent enclosure has 2 light sources. White lights under the cutting platform are used for visible stains like Coomassie Blue and silver, while the UV lights mounted above the platform are used to image fluorescent SYPRO Ruby protein stain. The enclosure creates a darkroom that produces quality images in either light mode.

## Contamination Control

The ProteomeWorks spot cutter provides excellent contamination control by drastically reducing keratin contamination compared with manual spot cutting (request bulletin 2503). The addition of the fluorescent enclosure adds another layer of protection from ambient



contamination and makes the system even more valuable to a busy proteome lab. The contamination control features built into the spot cutter include:

- Significantly reduced human interaction during spot cutting
- PDQuest<sup>™</sup> software sorts the cut list according to the increasing amount of protein in each spot
- Stainless-steel cutting tip
- 4 wash wells to wash the cutting tip
- Fluorescent enclosure that creates a physical barrier to ambient contamination

## Ordering Information

Catalog #	Description
165-7045	ProteomeWorks Fluorescent Enclosure, to add to previously purchased ProteomeWorks spot cutter (PDQuest software not included)
165-7042	ProteomeWorks Spot Cutter with Fluorescent Enclosure, PC (PDQuest software not included)
165-7043	ProteomeWorks Spot Cutter with Fluorescent Enclosure, Macintosh (PDQuest software not included)
165-7009	ProteomeWorks Spot Cutter, PC, without fluorescent enclosure (PDQuest software not included)
165-7039	ProteomeWorks Spot Cutter, Macintosh, without fluorescent enclosure (PDQuest software not included)
170-8603	PDQuest 2-D software, Windows
170-8611	PDQuest 2-D software, Macintosh

Coomassie is a trademark of Imperial Chemical Industries PLC. Macintosh is a trademark of Apple Computer. SYPRO is a trademark of Molecular Probes, Inc. Windows is a trademark of Microsoft Corp.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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** 86-10-8201-1366/68 **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 **Israel** 03 951 4124 **Italy** 34 91 590 5200 **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  
**Russia** 7 095 979 98 00 **Singapore** 65-2729877 **Spain** 34-91-590-5200 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 061-717-9555 **United Kingdom** 0800-181134