

Spot Cutter Template Instruction Guide

The Spot Cutter Template is a convenient tool for trimming large gels (>23 x 23 cm) to fit the ProteomeWorks spot cutter platform (figure 1). The ProteomeWorks spot cutter platform is 25 x 25 cm. Within this area, the functional cutting tip range is 22 cm starting from the right edge measured to the left, and 20.5 cm centered front to back within the platform. The inner viewing window of the Spot Cutter Template corresponds to the functional cutting area of the spot cutter's platform. The Spot Cutter Template allows the use of larger gels for optimal protein separation and resolution by easily and quickly trimming the gels to produce gels that fit perfectly on the platform.



Fig. 1. The Spot Cutter Template.

Note: Use caution when handling the Spot Cutter Template, the plastic cutting blade is sharp.

Trimming Procedure

1. Place the gel on a clean, flat surface.
Note: Keep in mind the orientation of the gel.
2. Position the Spot Cutter Template over the gel so the desired spot cutting area appears through the Spot Cutter Template's inner viewing window. See Figure 2 for illustration.

Note: Position the Spot Cutter Template over the gel without touching the gel surface or the gel may be improperly cut or torn.

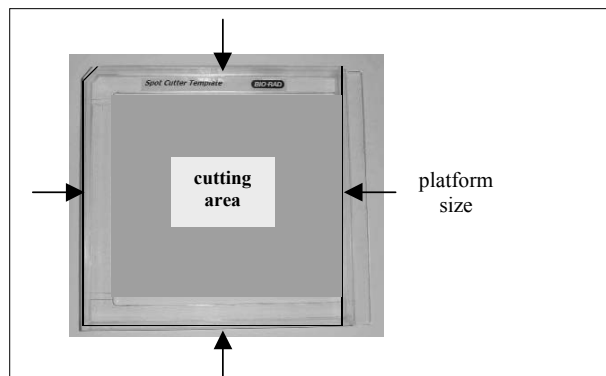


Fig. 2. The outside line corresponds to the ProteomeWorks Spot Cutter platform size. The inside line is the Spot Cutter Template's inner viewing window, which corresponds to the ProteomeWorks Spot Cutter's cutting area.

3. Gently place the Spot Cutter Template onto the gel surface.
4. Press down firmly on the top edges of the Spot Cutter Template to cut the gel.

Note: Make sure you apply pressure on all sides and corners of the Spot Cutter Template to make a clean cut through the gel's entire thickness.

Note: The cutting blade has a notch in the upper left corner to help distinguish the gel orientation.
5. Remove the excess gel edges that were trimmed from the gel.
6. Gently lift the Spot Cutter Template up from the gel.

Note: The gel may become attached to the cutting blade during trimming. Gently tap on the gel through the window to release it from the blade.
7. Rinse the Spot Cutter Template off with water.
8. Store the template in a secure place to protect the blade from damage.

Chemical Compatibility and Maintenance

The Spot Cutter Template is made of polyurethane. Several chemicals are not compatible with polyurethane: acetaldehyde, acetic acid, acetone, alcohol (including ethanol and methanol), benzene, chloroform, ethyl acetate, hydrazine, and methylene chloride. This list is representative, but not exhaustive. Some detergents contain strong chemicals, especially very basic detergents, which may cause damage to the product. Test the detergent on a small area of the product to determine if it is suitable to use.

Rinse the Spot Cutter Template with water after each use. Store the template in a secure place to protect the blade from damage.

Ordering Information

Catalog

Number Description

Spot Cutter Template

165-7050 Spot Cutter Template, 1 each

Other ProteomeWorks Spot Cutter Accessories

165-7011 1.5 mm Cutting Tip, 2 each
 165-7010 1.0 mm Cutting Tip, 2 each
 165-7018 Gel Cutting Sheets, 25 x 25 cm, 15 each



**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 **Czech Republic** (420) 2-4141 0532 **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 **Poland** (48) 22-8126 672 **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