



Instructions for Multi-Sample Tray II (170-7819)

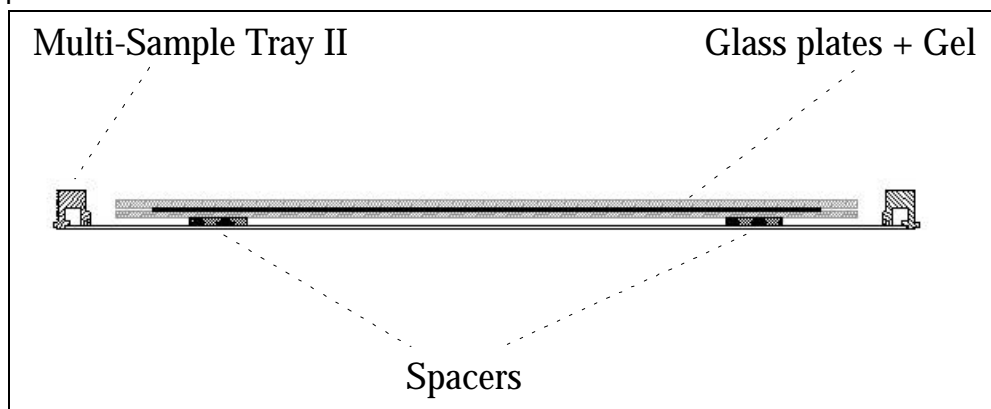
Usage:

The multi-Sample Tray II is designed specifically for imaging gel or samples that are bound to or are on top of a glass plate. It is designed for optimal sensitivity and resolution for differential display gels, VNTR, and mini-satellite gels. It also works well with fluorescent tissue sections, cell samples and fluorescent microarrays on glass slides.

For gels in between glass plates

1. Wipe glass plate - dry on both sides.
2. Place glass plate on top of an appropriate spacer(see table 1) inside the Multi-Sample Tray II. See figure 1.
3. Insert tray into the scanner and scan.

Figure 1



Hints:

For best sensitivity and resolution, remove the top plate. As much as a 30% increase in signal may be seen if the top glass plate is removed. Make sure to scan the gel immediately after removing the top plate since the gel may dry out. The gel can be kept moist by spray fine mist of water on top of the gel. Avoid pouring water directly on top of the gel since the sample may diffuse out - especially for thin gels.

For tissues, cells, DNA or protein on a glass slide

1. Place sample on appropriate spacer as described on table 1. Make sure the sample is facing upward.
2. Tape the glass slide firmly to the Multi-Sample tray II using clear scotch tape. See figure 2.

Table 1

Recommended spacers for samples bound on glass plate with different thickness.

Estimated glass plate thickness	Spacer use	Bio-Rad Part No.
4.5 mm (3/16 in.)	2.5mm	921-0117
3 mm (1/8 in.)	4.0mm	921-0118
1 mm	6.0mm	921-0119