



IMAGING AND ANALYSIS SOFTWARE

Gel/Blot Alignment Templates

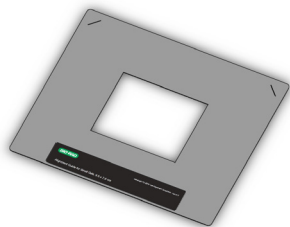
ChemiDoc™ Touch Imaging System

Quick Start Guide

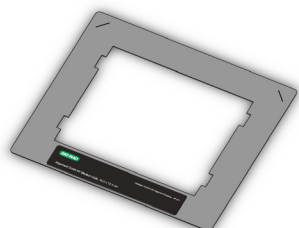
The Bio-Rad Gel/Blot Alignment Templates enable you to quickly and consistently place gels and blots on ChemiDoc™ Touch sample trays.

The Gel/Blot Alignment Templates (catalog #170-8376) contain the following:

- Alignment Template for Small Gels, 9.5 x 7.6 cm



- Alignment Template for Medium Gels, 15.5 x 12.4 cm



You can use the gel/blot alignment templates with all supported tray types. For more information about supported tray types, see the ChemiDoc Touch Imaging System with Image Lab Touch Software User Guide (#10042202).

Prepare the Gel/Blot Alignment Template

1. Select the gel/blot alignment template that matches the size of your sample.
2. Bend a corner at the pre-cut slit to make a tab.
Use the tab to remove the template from the tray after use.
3. Place the gel/blot alignment template on the sample tray.
4. Place the gel or gel tray on the open area of the template.



Alignment template for medium gels

Clean and Store Templates

It is important to keep the templates clean and to store them flat.

1. After using a gel/blot alignment template, rinse it with water and allow it to air dry.
2. Store the gel/blot alignment templates flat to prevent warping.



**Bio-Rad
Laboratories, Inc.**



Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **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** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000
