

Safe Band Removal with the Gel Excision Kit

Quick Guide

The Gel Excision Kit from Bio-Rad allows you to safely excise bands of interest from agarose or acrylamide gels for downstream applications such as DNA cloning or mass spectrometry. By using single-use plastic gel cutters instead of razor blades, you can avoid harming yourself or the transilluminator tray while minimizing cross-contamination between samples.

The procedure for excising bands depends on the type of sample tray you use. For specific instructions for safely excising bands on UV/Stain-Free, White, or Blue Trays for the GelDoc Go Imaging System, please refer to the Band Excision Using the GelDoc Go Imaging System Quick Guide (10000122134).

Excising Bands on a Tray

- 1 Pull out the transilluminator drawer and place a tray on the transilluminator. If using a UV/Stain-Free Tray, put on the required protective gear and ensure any ultraviolet (UV) shields are installed.
- 2 Place a gel sample in the center of the tray. Place the ruler next to the gel such that it is oriented along the direction of separation (Figure 1A).
- 3 Turn on the transilluminator to illuminate the gel with the appropriate type of light. Use the ruler to note the centimeter position of the gel band of interest.
- 4 Turn off the transilluminator and excise the bands.
 - a. If the bands are not readily visible by eye, use the ruler and noted location of the band to determine where to excise.
 - b. Carefully cut around the gel band of interest using a chopping (rather than sawing) motion with one of the disposable gel cutters (Figure 1B). For best results, remove as much gel as possible around the relevant band.
 - c. Pick up the gel band using the cutter and place in a centrifuge tube (Figure 1C).
- 5 When you have finished excising bands, acquire an image of the gel to confirm each band of interest has been excised.
- 6 Turn off the transilluminator and remove the gel and ruler from the tray. Slide the drawer in to close.



Caution: Sharp cutting tools can easily damage the surface of the tray. **Using the provided plastic disposable gel cutters with a chopping motion can reduce the chance of damage.**

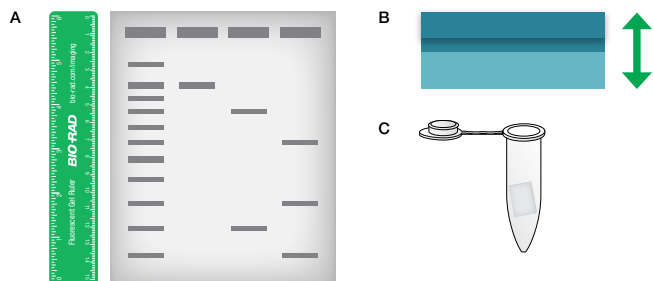


Fig. 1. Components of the Gel Excision Kit. The kit comes with a ruler (A) for determining band position, a disposable plastic gel cutter (B) for safe excision, and centrifuge tubes (C).

Complete Solutions for Your Needs with the GelDoc Go Imaging System

The GelDoc Go Imaging System gives you a benchtop imaging solution that is powerful enough to move your research forward, yet small enough to fit anywhere. Acquire high-resolution, publication-quality images of both nucleic acid and protein gels (Figure 2). Image up to four mini gels at once with the large imaging area provided by the system's advanced optical design. Easily set up and control the GelDoc Go Imaging System with Image Lab Touch Software.

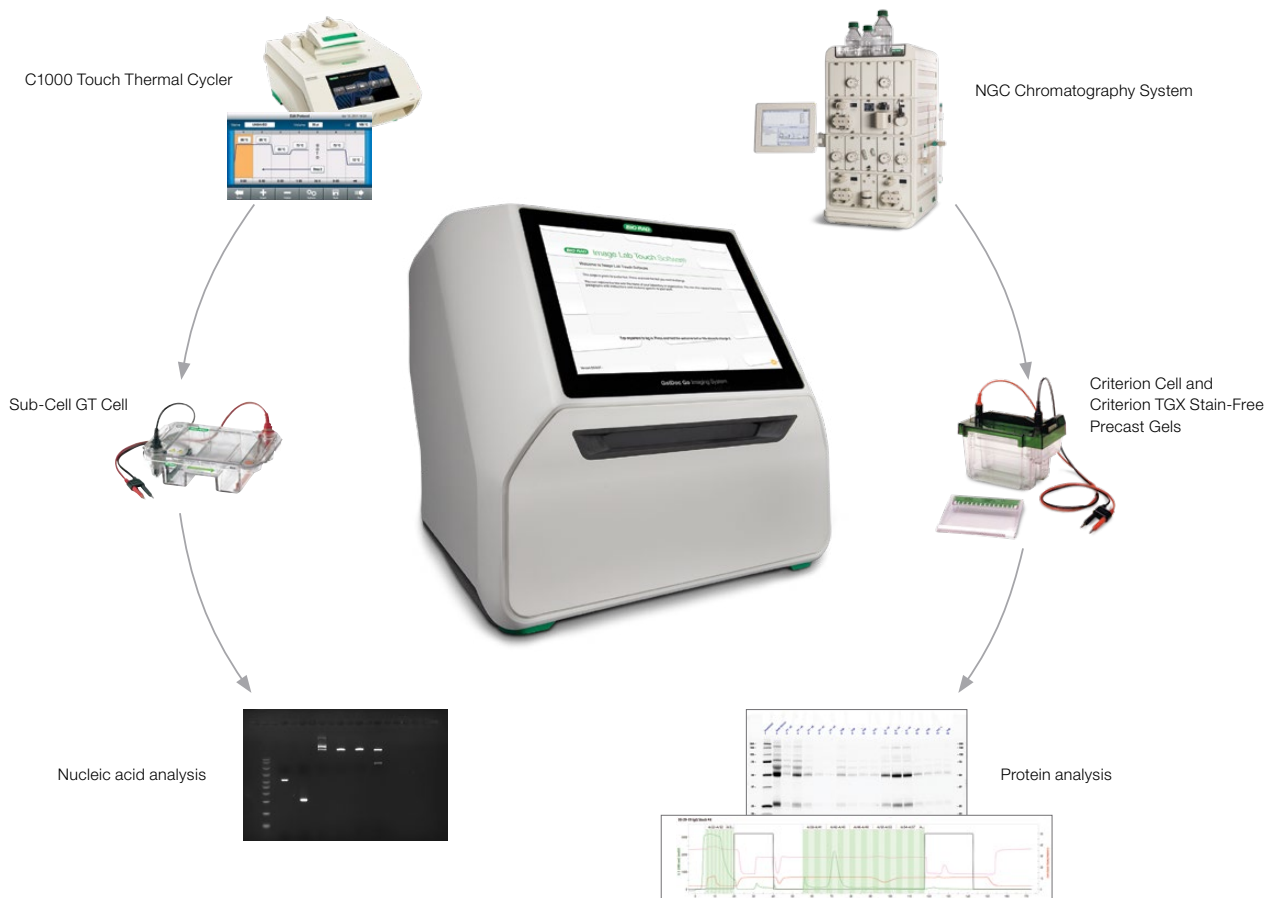


Fig. 2. Seamless integration into nucleic acid and protein workflows. The GelDoc Go Imaging System works with the Bio-Rad products you trust to provide complete solutions for nucleic acid and protein research. Use the C1000 Touch Thermal Cycler and the Sub-Cell GT Cell with the GelDoc Go Imaging System to generate and analyze PCR products. For protein workflows, pair the GelDoc Go Imaging System with the Criterion Cell and Criterion TGX Stain-Free Precast Gels to assess fractions from the NGC Chromatography System.

Ordering Information

Catalog #	Description
12009077	GelDoc Go Gel Imaging System with Image Lab Touch Software
2239480	EZ Micro Test Tubes, 1.5 ml

Visit bio-rad.com/SeeGelDocGo for more information.

BIO-RAD and SUB-CELL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

