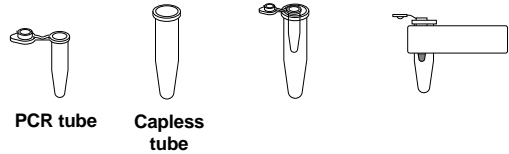


Quick Guide

Lesson 1: Setting up the PCR Reactions

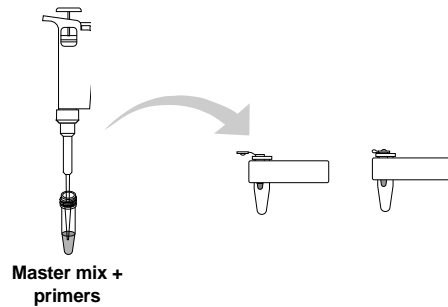
1. Label 5 PCR tubes CS, A, B, C, or D, and include your group name or initials as well. Place each PCR tube into a capless micro centrifuge tube in the foam float on ice.



2. Using the chart below as a guide, transfer 20 μ l of the appropriate template DNA into the correctly labeled tube. **Important: use a fresh aerosol barrier pipet tip for each DNA sample.**

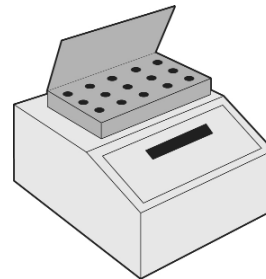
Label PCR tubes	Add DNA template	Add Master mix + primers
CS + your initials	20 μ l Crime Scene DNA	20 μ l MMP (blue)
A + your initials	20 μ l Suspect A DNA	20 μ l MMP (blue)
B + your initials	20 μ l Suspect B DNA	20 μ l MMP (blue)
C + your initials	20 μ l Suspect C DNA	20 μ l MMP (blue)
D + your initials	20 μ l Suspect D DNA	20 μ l MMP (blue)

3. Transfer 20 μ l of the blue MMP (master mix + primers) into each of the 5 PCR tubes containing template DNA. Pipet up and down to mix. Cap each tube after adding blue MMP. **Important: use a fresh aerosol barrier pipet tip each time. Immediately cap each tube after adding MMP.**



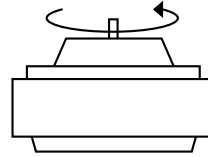
4. Place your capped PCR tubes in their adaptors on ice.

5. When instructed to do so, place your tubes in the thermal cycler. Your instructor will program the thermal cycler for PCR.

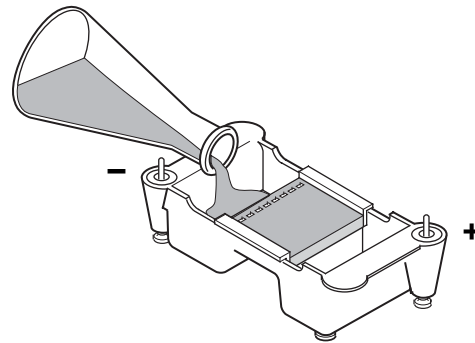


Lesson 2: Electrophoresis of PCR Products

1. Set up your gel electrophoresis equipment as instructed.
2. Obtain your 5 PCR tubes from the previous lesson. Place your PCR tubes in capless tubes and pulse-spin in a balanced microcentrifuge for a few seconds to collect all liquid to the bottom of the tube.
3. Transfer 10 μl of Orange G loading dye (from the tube labeled 'LD') into each of your PCR tubes. Pipet up and down to mix, and pulse-spin to collect liquid in the bottom of the tube.
4. Place an agarose gel in the electrophoresis apparatus. Check that the wells of the agarose gel are near the black (-) electrode and the base of the gel is near the red (+) electrode.
5. Fill the electrophoresis chamber with enough 1x TAE buffer to cover the gel. This will require ~275 ml of 1x TAE buffer.
6. Using a clean tip for each sample, load 20 μl of the samples into 6 wells of the gel in the following order:



Centrifuge

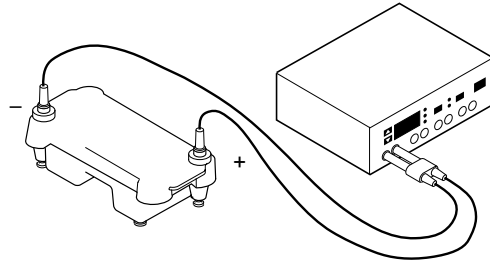


Lane	Sample	Load volume
1	Allele Ladder	20 μl
2	Crime Scene	20 μl
3	Suspect A	20 μl
4	Suspect B	20 μl
5	Suspect C	20 μl
6	Suspect D	20 μl

7. Secure the lid on the gel box. The lid will attach to the base in only one orientation: red to red and black to black. Connect the electrical leads to the power supply.

8. Turn on the power supply and electrophorese your samples at 100 V for 30 minutes.

9. Stain in Fast Blast DNA stain. Refer to the Student Manual for specific instructions.





**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 2527 3454
Canada (905) 712-2771 **China** (86 21) 6426 0808 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 318 84-0 **Greece** 30 210 777 4396 **Hong Kong** (852) 2789 3300 **Hungary** 36 1 455 8800
India (91-124)-2398112/3/4, 5018111, 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** 55-52-00-05-20 **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** 7 095 721 1404 **Singapore** 65-64153188 **South Africa** 00 27 11 4428508
Spain 34 91 590 52 00 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** (886 2) 2578 7189/2578 7241 **United Kingdom** 020 8328 2000