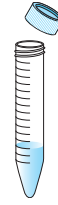


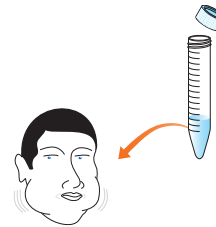
## Quick Guide for DNA Extraction and Precipitation

1. Obtain 15 ml tube containing 3 ml water from your instructor. Label the tube with your initials.



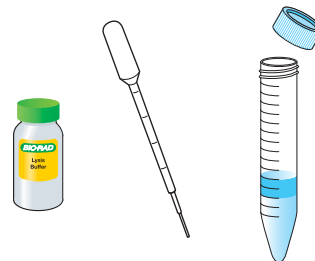
2. Gently chew the insides of your cheeks for 30 seconds. It is NOT helpful to draw blood!

3. Take the water from the 15 ml tube into your mouth, and swish the water around vigorously for 30 seconds.



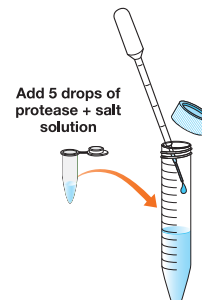
4. Carefully expel the liquid back into the 15 ml tube.

5. Obtain the tube of lysis buffer from your workstation, and add 2 ml of lysis buffer to your tube.

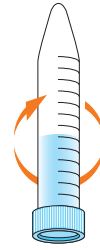


6. Place the cap on the tube, and gently invert the tube 5 times (don't shake your tube!). Observe your tube — do you notice any changes? If you do, write them down.

7. Obtain the tube of protease (**prot**) at your workstation. Add 5 drops of protease to your tube.



8. Place the cap on your tube, and gently invert it a few times.



9. Place your tube in a test tube rack or beaker in the water bath and incubate at 50°C for 10 minutes. Remove your tubes from the water bath.



50°C for 10 min

10. Obtain the tube of cold alcohol from your instructor or at the common workstation. Holding your tube at a 45° angle, fill your tube with cold alcohol, by adding approximately 10 mls to your tube. It will take repeated additions to add 10 ml of the cold alcohol using the disposable plastic transfer pipet.

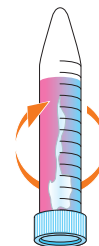


11. Place your cap on your tube, and let it sit undisturbed for 5 minutes. Write down anything you observe happening in the tube.



12. After 5 minutes, slowly invert the tube 5 times to help the DNA, which has begun to precipitate, to aggregate.

13. With a disposable plastic transfer pipet, carefully transfer the precipitated DNA along with approximately 750  $\mu$ l to 1 ml of the alcohol solution into a small glass vial provided in the DNA necklace kit (166-2200EDU). If you are not going to make a DNA necklace, save your DNA in a flip-top tube provided in this kit.





**Bio-Rad  
Laboratories, Inc.**

*Life Science  
Group*

**Web site** [www.bio-rad.com](http://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