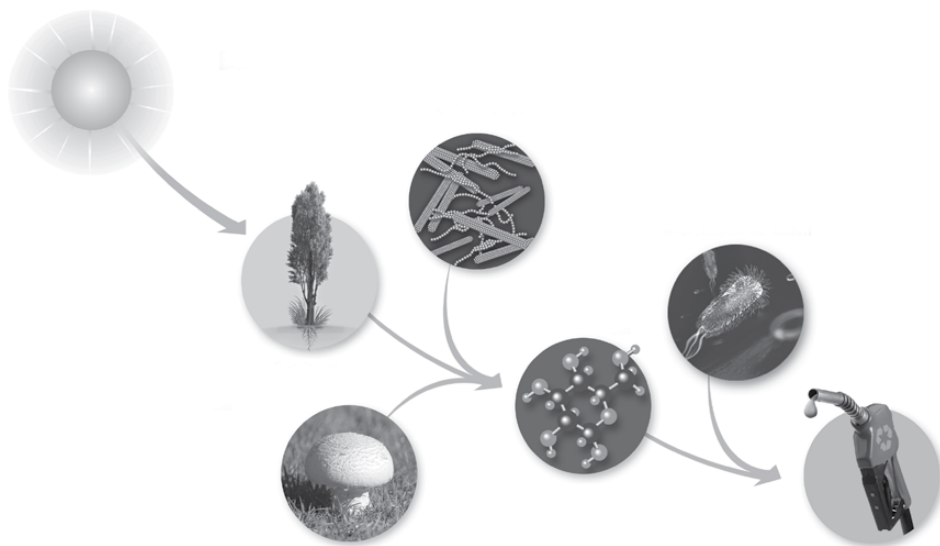

Biofuel Enzyme Reactions Kit for AP Biology: A ThINQ!™ Investigation

Quick Guide
explorer.bio-rad.com

Catalog #17001235EDU



Duplication of any part of this document is permitted for classroom use only.

Please visit explorer.bio-rad.com to access our selection of language translations for Bio-Rad Explorer kit curricula.

BIO-RAD

Quick Guide

Investigation 1: Determine How Mushroom Extracts Compare in Terms of Cellobiase Activity

1. Write down your mushroom type

2. Carefully remove the stem of your mushroom if you are using one with a woody stem like white button or shiitake. You will not need to remove the stem if you are using oyster or other soft-stemmed mushrooms.

3. Using a razor or knife, cut through the center of the cap of the mushroom. For mushrooms that do not have a distinguishable cap, cut through the meatiest section of the mushroom.



Mushroom skin
Mushroom gills

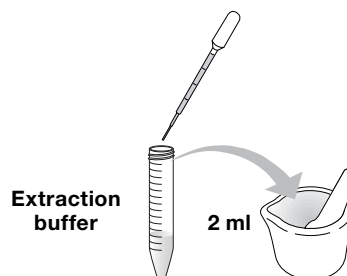
4. Using a cutting implement such as a razor cut out approximately 1 g (1,000 mg) of the internal flesh of the cap. Be careful to avoid including skin and gills in your sample.



Cut the meaty part of the cap, being careful to avoid skin and gills

5. Weigh out your sample and place it in a mortar: _____ mg

6. Add 2 ml of extraction buffer to the mortar for every gram of mushroom.



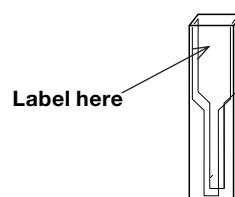
7. Using a pestle, grind your mushroom to produce a slurry — a semiliquid mixture.

- Strain the solid particles out of your slurry. If you have a centrifuge, scoop the slurry into a 1.5 ml microcentrifuge tube and then pellet the solid particles by spinning at top speed for 2 min. If you do not have a centrifuge, use a piece of cheesecloth to squeeze the extract into a 1.5 ml microcentrifuge tube. If you use the cheesecloth method, you will need to use 2 g mushroom and 4 ml of extraction buffer.

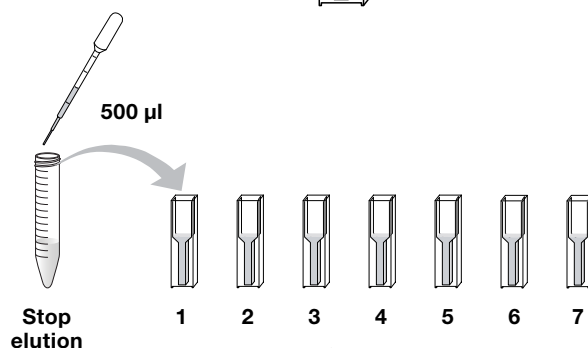


Note: You will need at least 250 μ l of extract to perform the enzymatic reaction portion of this protocol.

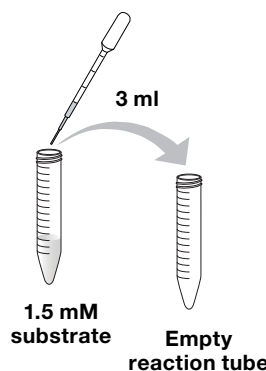
- Label your cuvettes 1–7. Write only on the upper part of the cuvette face so that your writing won't interfere with spectroscopy.



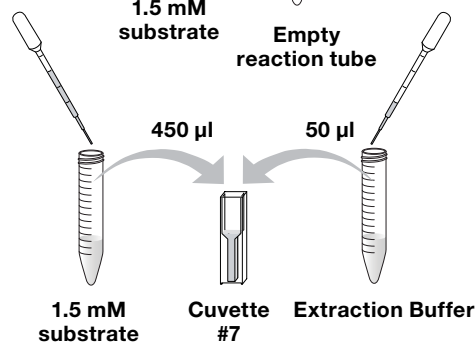
- Using a clean DPTP, pipet 500 μ l of stop solution into cuvettes 1–7. Rinse the DPTP thoroughly with water.



- Label a 15 ml conical tube with the type of mushroom you are using. Using a clean DPTP, pipet 3 ml of substrate into the tube.

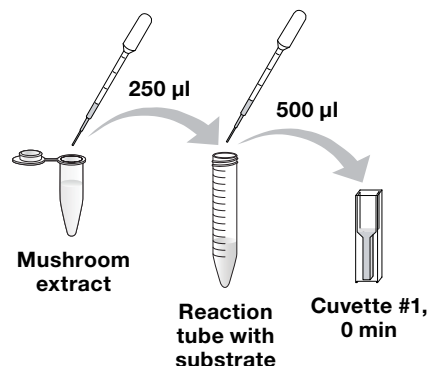


- Using the same DPTP from the previous step, add 450 μ l of substrate and 50 μ l of extraction buffer to cuvette #7 and set it aside for analysis later on. This cuvette will act as a control during your investigation because it does not contain mushroom extract.



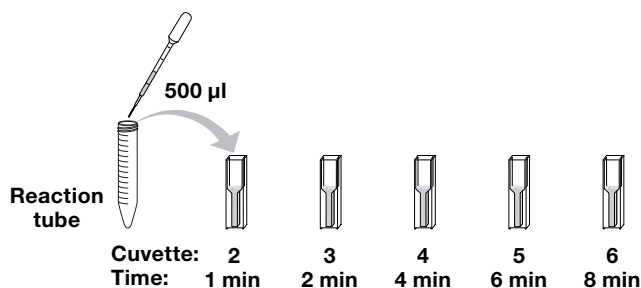
Please read and understand steps 13 and 14 fully before proceeding. These steps are time sensitive!

13. Using a clean DPTP, pipet 250 µl of your mushroom extract into the 15 ml conical tube containing 3 ml of substrate. This will now be referred to as the reaction tube. Pipet the liquid up and down to mix. Then pipet 500 µl from the 15 ml conical (reaction) tube to cuvette #1. **START YOUR TIMER.**



14. At the times indicated in the table remove 500 µl of mushroom extract/substrate mixture from the reaction tube and add it to the appropriately labeled cuvette that already contains stop solution.

Time	Cuvette
0 min	1
1 min	2
2 min	3
4 min	4
6 min	5
8 min	6



15. Rinse all DPTPs with copious amounts of water and save them for later investigations. After you have finished your analysis, rinse your reaction (conical) tubes and cuvettes with copious amounts of water and save them for later investigations.

Note: Do not discard the unused stock solutions or cuvettes containing standards. They will be used for the next investigation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

