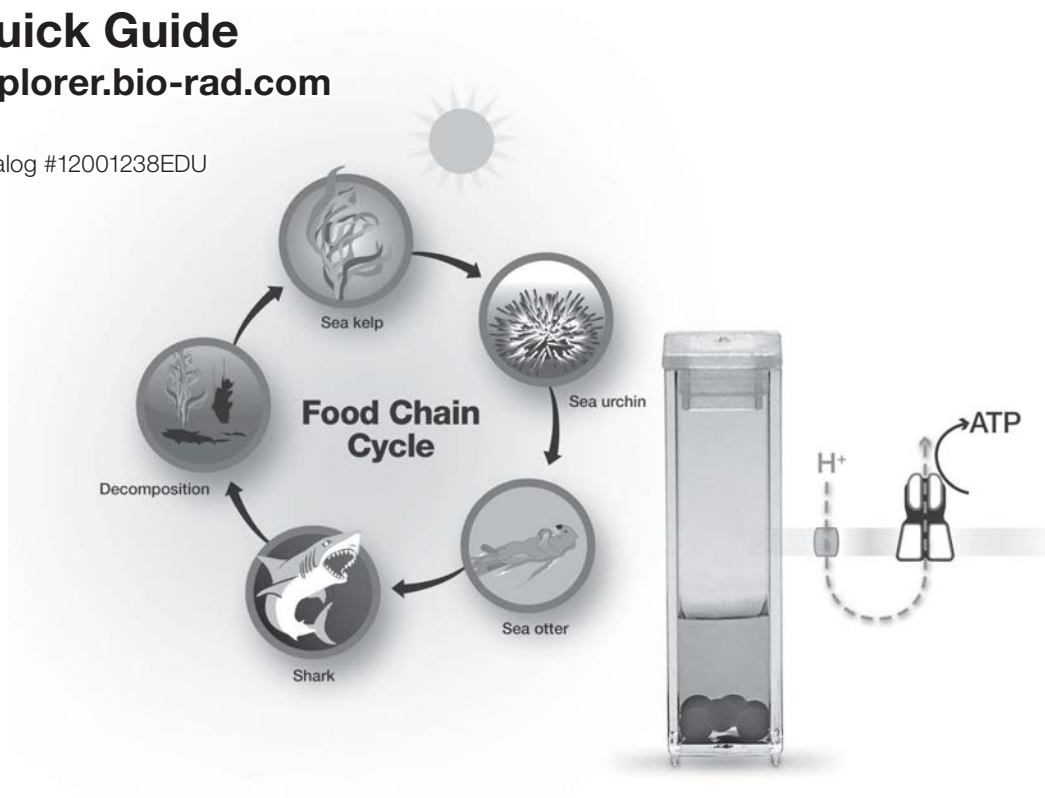


---

# Photosynthesis and Cellular Respiration Kit for AP Biology: A ThINQ!™ Investigation

Quick Guide  
[explorer.bio-rad.com](http://explorer.bio-rad.com)

Catalog #12001238EDU



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

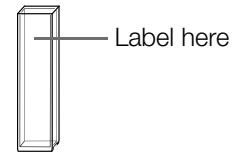
Please visit [explorer.bio-rad.com](http://explorer.bio-rad.com) to access our selection of language translations for Bio-Rad Explorer kit curricula.

**BIO-RAD**

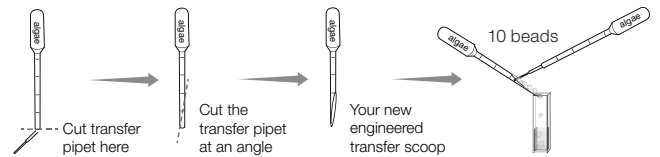
## Quick Guide

### Investigation #2: Photosynthesis and Cellular Respiration Core Lab

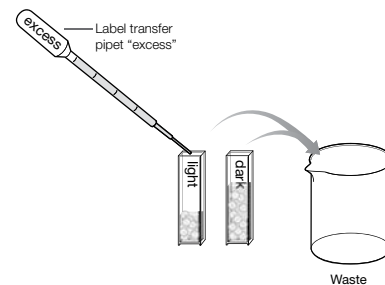
1. Label one empty cuvette **light**, and the other cuvette **dark**. Label each cuvette so that it does not obstruct light reaching the algae beads.



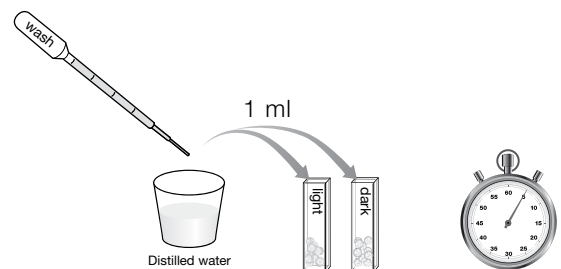
2. Label a transfer pipet **algae** and convert it into a scoop by cutting the transfer pipet at the 100  $\mu$ l mark diagonally. Use the **algae** transfer pipet to transfer 10 algae beads into each of the **light** and **dark** cuvettes.



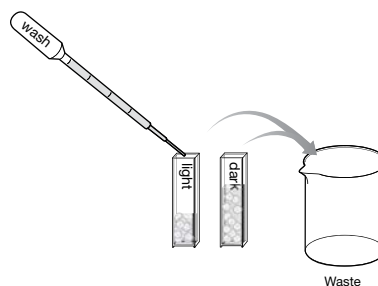
3. Label a new transfer pipet **excess** and use it to remove and discard the liquid that transferred along with the beads.



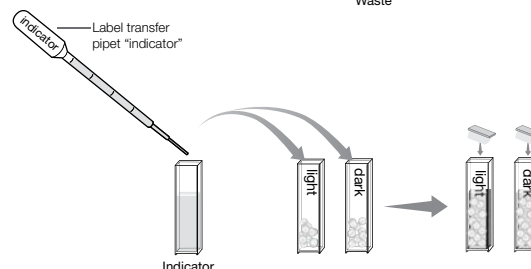
4. Label a new transfer pipet **wash** and use it to add 1 ml of distilled water to each of the cuvettes. Let the algae beads incubate in the water for 5 min to allow indicator within the bead to wash out.



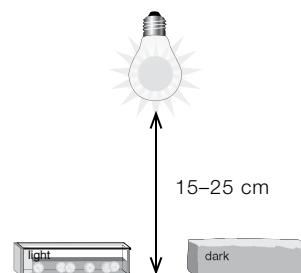
5. Use the wash transfer pipet to remove the water from the cuvette. Discard the water into the waste container.



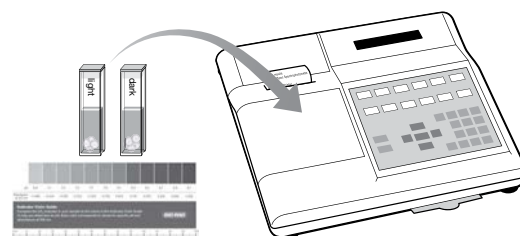
6. Label a new transfer pipet **indicator** and use it to transfer 1 ml of CO<sub>2</sub> indicator to each cuvette. Cap cuvettes tightly.



7. Wrap the cuvette labeled **dark** in aluminum foil being sure to cover both ends as well as all sides. Place both the cuvettes labeled **light** and **dark** on their sides 15–25 cm from the lamp. Ensure that the beads are distributed evenly throughout the cuvette and the clear side of the **light** cuvette faces the light.



8. Collect data starting at time = 0 min. Every 5 min, thoroughly mix the CO<sub>2</sub> indicator in the cuvettes and determine the color. This can be done by comparing the color of the CO<sub>2</sub> indicator in your cuvette to the provided Indicator Color Guide, or by reading the absorbance at 550 nm ( $A_{550}$ ) in a spectrophotometer (make sure your teacher has zeroed the machine). Be quick about taking this reading and immediately return the cuvettes to the experimental conditions.



9. If enough time remains after the last time point, switch the light and dark cuvettes. Place the cuvette labeled **light** in the dark and the cuvette labeled **dark** in the light. Continue to record pH or  $A_{550}$  every 5 min.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

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

