

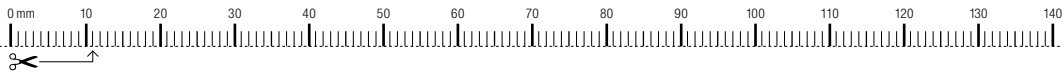


pH 6.9 7.1 7.3 7.5 7.7 7.9 8.1 8.3 8.5 8.7 8.9 9.1

Absorbance at 550 nm -0.485 -0.434 -0.402 -0.353 -0.223 -0.126 0.000 0.178 0.400 0.616 0.849 1.009

Indicator Color Guide

Compare the CO₂ indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.

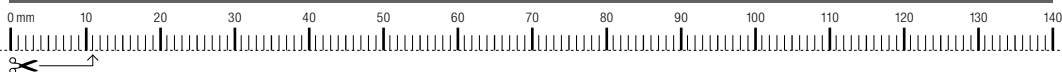


pH 6.9 7.1 7.3 7.5 7.7 7.9 8.1 8.3 8.5 8.7 8.9 9.1

Absorbance at 550 nm -0.485 -0.434 -0.402 -0.353 -0.223 -0.126 0.000 0.178 0.400 0.616 0.849 1.009

Indicator Color Guide

Compare the CO₂ indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.

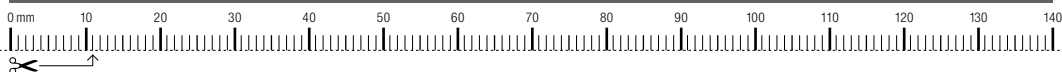


pH 6.9 7.1 7.3 7.5 7.7 7.9 8.1 8.3 8.5 8.7 8.9 9.1

Absorbance at 550 nm -0.485 -0.434 -0.402 -0.353 -0.223 -0.126 0.000 0.178 0.400 0.616 0.849 1.009

Indicator Color Guide

Compare the CO₂ indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.

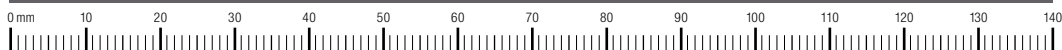


pH 6.9 7.1 7.3 7.5 7.7 7.9 8.1 8.3 8.5 8.7 8.9 9.1

Absorbance at 550 nm -0.485 -0.434 -0.402 -0.353 -0.223 -0.126 0.000 0.178 0.400 0.616 0.849 1.009

Indicator Color Guide

Compare the CO₂ indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.



Equilibrated CO₂ indicator color

pH 6.9 7.1 7.3 7.5 7.7 7.9 8.1 8.3 8.5 8.7 8.9 9.1

Absorbance at 550 nm -0.485 -0.434 -0.402 -0.353 -0.223 -0.126 0.000 0.178 0.400 0.616 0.849 1.009

Instructor's Indicator Color Guide

The color range of equilibrated CO₂ indicator is shown above. The color of your equilibrated CO₂ indicator may differ slightly depending on several variables, including the water you use and altitude. The absorbance values at 550 nm correspond to readings that were taken with a spectrophotometer that was zeroed with CO₂ indicator that equilibrated to pH 8.1.

Online Resources

A PDF of the Indicator Color Guide is available at bio-rad.com/algae.

You may use it to print additional copies. However, depending on your printer, the colors may vary significantly from the Indicator Color Guide provided in the kit.

Additional resources for the Photosynthesis and Cellular Respiration Kit are available at bio-rad.com/algae. These include:

- Teacher model process
- Student manual
- Science case study
- And many more teaching resources



Follow us [@BioRadEducation](https://twitter.com/BioRadEducation) to get the latest in science teaching news, stories, and events.