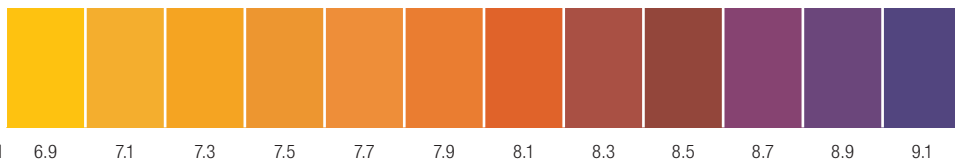
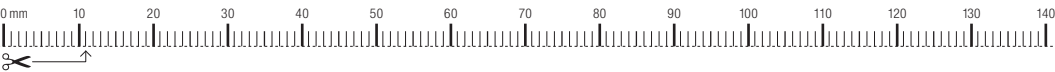


Absorbance at 550 nm	-0.485	-0.434	-0.402	-0.353	-0.223	-0.126	0.000	0.178	0.400	0.616	0.849	1.009
----------------------	--------	--------	--------	--------	--------	--------	-------	-------	-------	-------	-------	-------

**Indicator Color Guide**

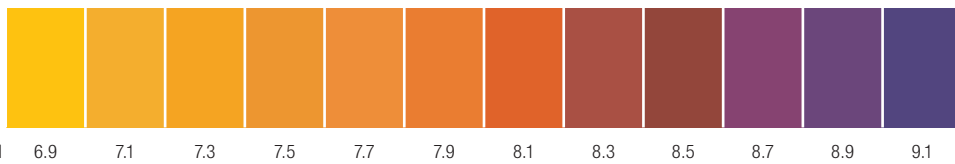
Compare the CO<sub>2</sub> indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.



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**

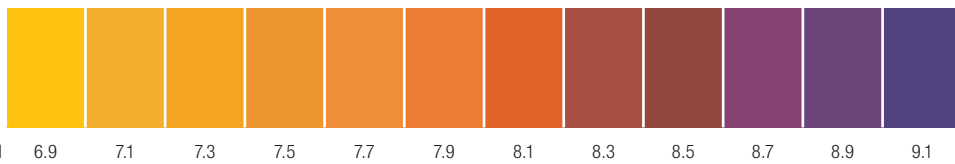
Compare the CO<sub>2</sub> indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.



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**

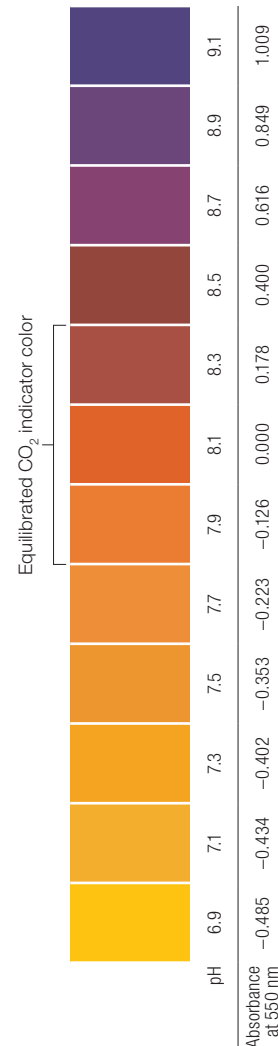
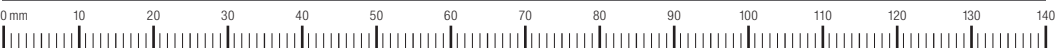
Compare the CO<sub>2</sub> indicator in your sample to the colors in this Indicator Color Guide to help you determine its pH. Each color corresponds to values for specific pH and absorbance at 550 nm.



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 CO<sub>2</sub> 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.



**Instructor's Indicator Color Guide**

The color range of equilibrated CO<sub>2</sub> indicator is shown above. The color of your equilibrated CO<sub>2</sub> 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 CO<sub>2</sub> indicator that equilibrated to pH 8.1.



**Online Resources**

A PDF of this guide is available at [bio-rad.com/ColorGuide](http://bio-rad.com/ColorGuide). 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.

Visit [explorer.bio-rad.com](http://explorer.bio-rad.com) for other great teaching resources.



Follow us @BioRadEducation to get the latest in science teaching news, stories, and events.