

<b>Date:</b>	2022-11-08
<b>Reference:</b>	ImageXpress Pico Installation Guide (5063666 M) ImageXpress Pico User Guide (5063667 L)
<b>Overview:</b>	This Update Information Sheet describes changes to the gas supply tubing and corresponding labels for the ImageXpress Pico Environmental Control System. These changes are not shown in the current Installation Guide (5063666 M) and User Guide (5063667 L). Both guides will be updated in the next revision. The packing list included with the instrument corresponds to the updated gas tubing.
<b>Description:</b>	The current guides show BLUE tubing used for all gas supply connections. The ImageXpress Pico Environmental Control System now includes color-coded tubing (BLUE, BLACK, and GRAY). In addition, the labels on the rear of the instrument are also color-coded. This improvement will help you clearly identify each gas supply port and the corresponding tubing.

## Documentation Changes

All page references refer to Installation Guide (5063666 M) and the User Guide (5063667 L).



Pages 20 (Installation) and 19 (User):  
Updated rear view of ImageXpress Pico with the Environmental Control System showing the updated labels for gas supply connections.



Pages 53 (Installation) and 41 (User):  
Updated labels for gas supply connections with gray, blue, and black markings corresponding to the colors of the gas supply tubing.



Pages 55 (Installation) and 43 (User):  
Updated labels next to Environmental Control Tray connector.



Pages 57 (Installation) and 45 (User):  
Compressed air (blue) and humidifier column  
connected. The CO<sub>2</sub> (gray) and N<sub>2</sub> (black) ports  
are closed with blind plugs.



Pages 57 (Installation) and 45 (User):  
Cover of Environmental Control Tray installed.



Pages 53 and 54 (User only):  
Compressed Air (blue) connected. The CO<sub>2</sub>  
(gray) and N<sub>2</sub> (black) ports closed with blind  
plugs.



New Photo:  
All gas supply ports connected: CO<sub>2</sub> (gray),  
Compressed air (blue) and N<sub>2</sub> (black).