



August 25, 2021

From: Eric Giesing, Director

Re: HIAC ROC Instrument Obsolescence

Dear HIAC ROC Customer,

This is an obsolescence notice for the HIAC ROC remote industrial online particle counter. Due to a combination of parts obsolescence and a strategic business pivot, the instrument will become obsolete and no longer available for purchase as of August 25, 2021. We will honor and ship instruments for purchase orders received up to the obsolescence date of August 25, 2021.

There is not currently, nor is there projected to be, a replacement or alternative product available from Beckman Coulter Life Sciences for the public market. Please note that after the date of obsolescence, additional software and hardware improvements or updates will not be made.

We will continue to fully service and support the HIAC ROC for a period of five (5) years after the obsolescence date up to August 25, 2026. During this time, calibrations, repairs, and reagents required for routine service will be available, however, breakdown repairs cannot be guaranteed if the required repair component is made obsolete by our supplier.

Instrument part numbers affected by the HIAC ROC obsolescence include:

PC9001-01	ROC, PC9001 LP FLOW WITH DISPLAY
PC9001-11	ROC, PC9001 HP FLOW WITHOUT DISPLAY
PC9001-21	ROC, PC9001 BYPASS WITHOUT DISPLAY
PC9001-31	ROC, PC9001 NO CNTRL WITHOUT DISPLAY
PC9001-41	ROC, PC9001 HP FLOW WITH DISPLAY
PC9001-51	ROC, PC9001 BYPASS WITH DISPLAY
PC9001-61	ROC, PC9001 NO CNTRL WITH DISPLAY
PC9001-71	ROC, PC9001 MP FLOW WITHOUT DISPLAY
PC9001-81	ROC, PC9001 MP FLOW WITH DISPLAY
ROC-01	ROC, HIAC LP FLOW WITH DISPLAY
ROC-11	ROC, HIAC HP FLOW WITHOUT DISPLAY
ROC-21	ROC, HIAC BYPASS WITHOUT DISPLAY
ROC-31	ROC, HIAC NO CNTRL WITHOUT DISPLAY
ROC-41	ROC, HIAC HP FLOW WITH DISPLAY
ROC-41-ISO	ROC, HIAC HP FLOW W DISP ISO-MTD
ROC-51	ROC, HIAC BYPASS WITH DISPLAY
ROC-51-ISO	ROC, HIAC BYPASS W DISP ISO-MTD

ROC-61	ROC, HIAC NO CNTRL WITH DISPLAY
ROC-71	ROC, HIAC MP FLOW WITHOUT DISPLAY
ROC-81	ROC, HIAC MP FLOW WITH DISPLAY
ROC-91	ROC, HIAC LP FLOW WITHOUT DISPLAY
C49712	HIAC ROC-51 with CD

We apologize for the abrupt nature of this decision and are happy to help answer questions you may have moving forward. Please contact your local sales representative to find out how this may impact a current or planned order. For technical questions or details regarding continued service and support, please contact Customer Technical Support (CTS) at (800) 369-0333.

Thank you,



Eric Giesing  
Director, Biotechnology Business Unit, Beckman Coulter Life Sciences