



August 25, 2021

From: Eric Giesing, Director
Re: HIAC PODS+ Instrument Obsolescence

Dear HIAC PODS+ Customer,

This is an obsolescence notice for HIAC PODS+ portable industrial liquid particle counters. 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 PODS+ instrument 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 PODS+ obsolescence include:

B95061	PODS+ SDM, 1.1
B95062	PODS+ SDM, 1.2, R
B95063	PODS+ SDM, 1.3
B95064	PODS+ SDM, 1.4, R
B95065	PODS+ SDM, 1.5
B95066	PODS+ STD, 2.1
B95067	PODS+ STD, 2.2, R
B95068	PODS+ STD, 2.3
B95069	PODS+ STD, 2.4, R
B95070	PODS+ STD, 2.5
B95071	PODS+ STD, 2.6, R
B95072	PODS+ STD, 2.7
B95073	PODS+ STD, 2.8, R
B95074	PODS+ STD, 2.9
B95075	PODS+ STD, 2.10, R
B95077	PODS+ STD, 2.12, R
B95078	PODS+ STD, 3.1

B95079	PODS+ STD, 3.2, R
B95080	PODS+ STD, 3.3
B95081	PODS+ STD, 3.4, R
B95082	PODS+ STD, 3.5
B95083	PODS+ STD, 3.6, R
B95084	PODS+ STD, 3.7
B95085	PODS+ STD, 3.8
B95086	PODS+ STD, 3.9
B95087	PODS+ STD, 3.10
C29888	PODS+ F-35

Accessory part numbers affected by the HIAC PODS+ obsolescence include:

B88730	Bottle Cap, GL45 QC Insert Simriz
B92123	Transit Case, PODS+ & Accessories
B93458	Dust Cover, HIAC PODS+
B95366	On-Line Adapter Kit, PODS+
B96273	Quick Start Guide, PODS+ English
B96581	Ship Kit, PODS+ STD
B96582	Ship Kit, PODS+ STD with Online
C29889	Ship Kit, PODS+ F-35
C29906	Transit Case, PODS+ F-35 and Accessories
C34596	Document, CE, PODS+

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