



August 25, 2021

From: Eric Giesing, Director
Re: LS 13 320 XR Instrument Obsolescence

Dear LS 13 320 XR Customer,

This is an obsolescence notice for the LS 13 320 XR particle size analyzer. Due to 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. 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 LS 13 320 XR 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 LS 13 320 XR obsolescence include:

B98100	LS 13 320 XR Optical Bench MW
B98101	LS 13 320 XR Optical Bench SW
C27180	LS 13 320 XR Optical Bench MW w/Workstation
C27181	LS 13 320 XR Optical Bench SW w/Workstation
C53018	LS 13320 XR (MW) and (ULM) Module
C53019	LS 13320 XR (MW) and (DPS) Module
C53020	LS 13320 XR (MW) with (ULM) and (DPS) Modules
C53021	LS 13320 XR (SW) and (ULM) Module
C53022	LS 13320 XR (SW) and (DPS) Module
C53023	LS 13320 XR (SW) with (ULM) and (DPS) Modules

Instrument accessory part numbers affected by the LS 13 320 XR obsolescence include:

B95435	Upgrade Kit, Sonicator
B98103	Dry Powder System
B98105	Universal Liquids Module
C06826	HiCap HEPA Vacuum (110 Volt, Model CFM: 99)
C20773	Filter, RFI, 250AC 60Hz, 10A, IEC320
C20930	Workstation Assy LS13320XR, Configured with Software

C24287	Bench Dust Cover
C32506	Hose, Vac, ESD., 1.26OD., 1.00ID., PVC., 10FT

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