



August 25, 2021

From: Eric Giesing, Director
Re: HIAC 8011+ Instrument Obsolescence

Dear HIAC 8011+ Customer,

This is an obsolescence notice for HIAC 8011+ industrial laboratory liquid particle counters. Due to a combination of parts obsolescence and a strategic business pivot, the 8011+ 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 8011+ 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 8011+ obsolescence include:

2089800-01	8011+ HIAC LIQUID PARTICLE COUNTER
8011+051	KTO: 8011+, HRLD-100, PSL
8011+053	KTO: 8011+, HRLD-100, MTD
8011+054	KTO: 8011+, HRLD-100, 11171
8011+061	KTO: 8011+, MC-05, PSL
8011+071	KTO: 8011+, HRLD-150, PSL
8011+073	KTO: 8011+, HRLD-150, MTD
8011+074	KTO: 8011+, HRLD-150, 11171
8011+081	KTO: 8011+, HRLD-150JA, PSL
8011+083	KTO: 8011+, HRLD-150JA, MTD
8011+084	KTO: 8011+, HRLD-150JA, 11171
8011+091	KTO: 8011+, HRLD-400, PSL
8011+0A1	KTO: 8011+, HRLD-400HC, PSL
8011+0B1	KTO: 8011+, HRLD-600JS, PSL
8011+0C1	KTO: 8011+, HRLD-100HC, PSL
8011+0C3	KTO: 8011+, HRLD-100HC, MTD
8011+0C4	KTO: 8011+, HRLD-100HC, 11171

8011-3833	8011, HRLD-150JA, 110V, MTD, Corrosive
C18729	Nalco 8011+, HRLD-150JA
C18800	KTO: Nalco 8011+ with HRLD-100

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