



November 16, 2022

To: Beckman Coulter Life Sciences Customers

From: Bhagya Wijayawardena, Vi-CELL Global Commercial Product Manager

Re: Vi-CELL XR Cell Viability Analyzer Product Life Cycle Update

Dear Valued Customer,

This letter is to inform you of our product life cycle plan for the Vi-CELL XR Cell Viability Analyzer.

First, an explanation of the terminology used in our product life cycle management process:

- **Obsolescence Date** - Date when new instruments can no longer be sold
- **Discontinuance** – Service of the instruments is no longer supported
- **Best Effort** – Full support cannot be guaranteed; Beckman Coulter will continue to provide support and maintenance to the extent that the necessary parts and software components are available.

The planned obsolescence date for Vi-CELL XR instruments is **December 31st, 2022**, with Service and Support anticipated to continue through **December 2027**. Our Service Organization will assess parts availability throughout the post-obsolescence period to determine if/when support will transition to Best Effort and/or Discontinuance.

Additional notification will be provided once an official discontinuation date has been established for ongoing support.

We hope this information will help you plan for the pending obsolescence of the Vi-CELL XR analyzer. For your convenience, your Beckman Coulter Life Sciences representative is here to help you with forecasting your needs for Vi-CELL XR analyzers and/or evaluating the Vi-CELL BLU cell viability analyzer as a replacement.

Thank you,
Bhagya Wijayawardena
Vi-CELL Global Commercial Product Manager

Product is not verified or validated for use in diagnostic procedures.

©2022 Beckman Coulter, Inc. All rights reserved.

Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

2022-GBL-EN-100556-v1