

In this issue: Identify the purity grade needed for your workflow

Issue | 001

What's the difference between sterile & Certified Free consumables, and which purity grade do I need?

Protecting your sample from degradation and contamination throughout the workflow is critical to consistently achieving a high-quality final product. Because every workflow and product is different, with differing requirements for various purity grades, it's important to select consumables supporting these needs. Sterile and Certified Free* consumables from Beckman Coulter are available to meet these varying needs.

Senior Applications Scientist Ross VerHeul, PhD

Sterile vs. Certified Free

Sterile – Absent of Living Organisms

Sterility is defined as an aseptic state where living organisms, such as bacteria and fungi, do not exist. Beckman sterile consumables are sterilized with a validated sterilization process using gamma irradiation or ethylene oxide. Sterilization procedures such as these and autoclaving eliminate viable organisms, though nonviable contaminants may remain.





Certified Free – Free of Nonviable Biologics

Certified Free consumables exhibit no detectable DNA, DNase, RNase, PCR inhibitors, or endotoxins. These nonviable contaminants are not removed via standard sterilization procedures, such as autoclaving, and the absence of such contaminants is critical for a variety of applications from nucleic acid sequencing to gene therapy.

Determining the Necessary Purity Grade

Each application and workflow is unique, with varying acceptable levels of viable and nonviable biologic materials. Accordingly, samples may require consumables verified as sterile, Certified Free, or both. Choosing the right tube, bottle, or liner provides your sample or experiment the best possible outcome.

Purity Grade Guidance by Application



Summary

Selecting the appropriate consumable purity grade is an important step in maintaining the integrity of your product, and Beckman Coulter offers a variety of sterile and Certified Free consumables designed to protect your samples.



*Based on sample results below detectable limit.

© 2018 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service names mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com CENT-4393FLY11.18