



The Particle Analyzers Advancing Life Science Research

Scientists share how automated cell counting is revolutionizing research

Technology that automates cell counting provides scientists with the ability to expand analyses to detect thousands of cells in one reading, saving precious time and preventing the risk of human error associated with laborious manual hemocytometer-based methods.

Beckman Coulter is a leader in technology for this purpose, and offers a range of cell counters and particle analyzers suitable for accurate analysis of any cell type --as well as the analysis of other particles-- for a broad range of applications including cell biology, biopharmaceuticals, the brewing industry, material analysis, and environmental monitoring.

Based on Coulter Principle technology, instruments in the Multisizer line of analyzers quantitatively count and size cells and particles between 0.2 μm to 1,600 μm at very high resolution in a single measurement. This enables fast analysis of cell concentration and viability to deliver consistent, accurate data.

Here, we share the opinions of scientists working in the life sciences and pharmaceutical industry, to learn how these analyzers are improving their research.



"Reproducible, fast and reliable. This has saved our lab a lot of time."

Mark Parkinson, of Covance Laboratories, uses the **Multisizer 4e Coulter Counter** in pharmaceutical analysis.





"This product saves lots of time and is less laborious. It is more efficient than manual cell counting. This is an example of evolution in the cell culture field."

Dr. Yuling Wu, of Medimmune, uses the **Vi-CELL® XR Cell Viability Analyzer** for translational research using cell cultures



"This is a very high-end, easy-to-use product that makes counting cells so much easier than a hemocytometer! One of the best pieces of equipment in our lab."

Dr. Elizabeth Tovar, of Wayne State University, uses the **Vi-CELL® XR Cell Viability Analyzer** in tissue culture research and for downstream assays.



Find out more about the **Vi-CELL® XR Cell Viability Analyzer** and the **Multisizer 4e Coulter Counter**



© 2018 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. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com

PART-336401.18