

FEATURES IN THIS BULLETIN

- Bacteria
Aggregation
- Disinfection
Efficacy
- Reporting the
Results

Evaluating Contact Lens Disinfection Efficacy Using the Beckman Coulter™ Multisizer™ 3 COULTER COUNTER®

Beckman Coulter Application Bulletin

Multisizer™ 3 COULTER COUNTER®

The percentage of aggregates in bacteria cultures may be determined by employing the Coulter Principle, also known as Electrical Sensing Zone (ESZ) technique. The ISO International Standard 13319 describes this technique. A suitable electrolyte solution is required to perform the analysis, in many cases the Phosphate Buffered Saline (PBS) solution used to suspend the cells may be used as the electrolyte. The sample is prepared and then analyzed using a Beckman Coulter Multisizer 3 to determine the size distribution of the cells. The results are reported as number percent of aggregates.

The use of the Multisizer 3 provides a fast, easy, accurate and automatic method to determine cell aggregation. This instrument also provides reliable results not dependent of the operator's judgment as opposed to results reported when using a hemocytometer.



EVALUATING CONTACT LENS DISINFECTION EFFICACY

SIGNIFICANCE

Contact lens care disinfecting products must be evaluated for their safety and effectiveness. Procedures used to evaluate the efficacy of lens care disinfection solutions routinely employ a specific panel of potentially pathogenic challenge organism suspensions. Microscopic examination of challenge organism suspensions has demonstrated the presence of cell clumps. These aggregates prevent uniform exposure to the test disinfectant and result in underestimation of the true number of challenge organisms. In order to reduce the number of clumps, and therefore produce a more uniform cell suspension, some methods suggest filtration of bacterial challenge organism suspensions prior to use as test inoculate. The method also allows the use of suspension mediums containing 0.05% Tween 80 as a dispersant. The Multisizer 3 Coulter Counter may be used to conduct analysis on cell suspensions pre- and post-filtration to evaluate filtration effectiveness by determining the size distribution and relative number of cell clumps in the bacterial cell suspension. The effectiveness of the Tween 80 may also be evaluated.

INSTRUMENT SET UP AND CALIBRATION

A 30 μm and a 100 μm aperture tubes should be used for the analysis. The linear dynamic range for any aperture is 2% to 60% of its diameter. A 30 μm aperture tube will be capable of analyzing the particle concentration and size distribution from 0.6 μm to 18 μm and a 100 μm from 2 μm to 60 μm . The two apertures combined yields an analysis range of 0.6 μm to 60 μm . Set up and calibrate the instrument according to the Multisizer 3 Operator's Manual. For determining number percent size distribution the control mode for the instrument may be Time Mode selecting 60 seconds as the run time.

PROCEDURE

1. *Electrolyte Preparation*

Isoton[®] II is a balanced electrolytic solution manufactured by Beckman Coulter. It is approximately 1% total salt concentration, (sodium and potassium chloride plus a phosphate buffer system). Isoton II may be used to suspend the cells for analysis on the Multisizer 3. If a different saline solution is used, there will be two possibilities:

- The custom made saline solution contains approximately 1% total salt concentration. In this case the electrolyte in the external container may be Isoton[®] II or the custom made saline solution.
- The custom made saline solution total salt concentration is different from 1%. In this case the electrolyte in the external container must be the custom made saline solution.

In both cases the cells may be analyzed using the custom made saline solution as electrolyte for the beaker containing the sample.

2. *Sample Analysis.*

Use a small round bottom beaker to analyze the cells using a 100 μm aperture. Place the beaker containing the sample into the analyzer, use the Multisizer stirrer to ensure a uniform distribution of the sample in the beaker. Flush the aperture tube before each analysis. After each run rinse the aperture and electrode before proceeding to the next sample. When all the samples have been analyzed, filter each using a 10 μm filter. Run the filtrates using a 30 μm aperture tube.

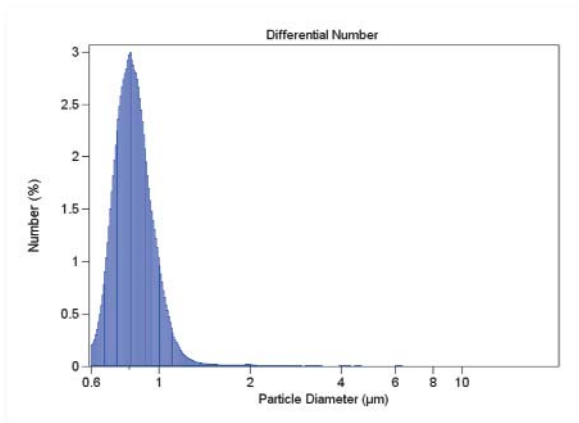
3. *Data Presentation.*

The results from the analyses using the 30 μm and 100 μm apertures for each sample may be presented as a single size distribution covering the combined range for both apertures. Refer to the Multisizer 3 Operator's Manual for information about the "Multi-tube Overlap" function.

EVALUATING CONTACT LENS DISINFECTION EFFICACY

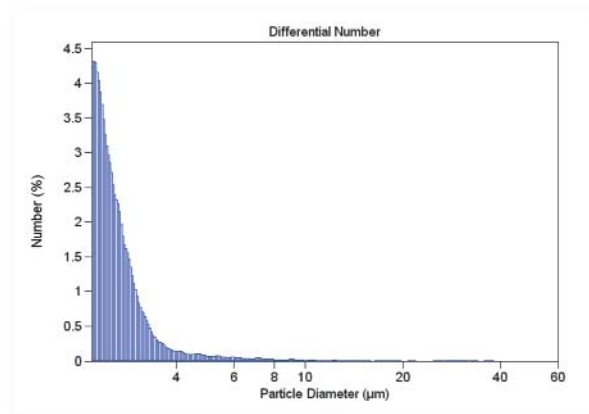
REPORTING THE RESULTS

Results may be represented as the number percentage of cells above selected size categories in the range of the size distribution by using the Interpolation feature in the Multisizer 3 Software.

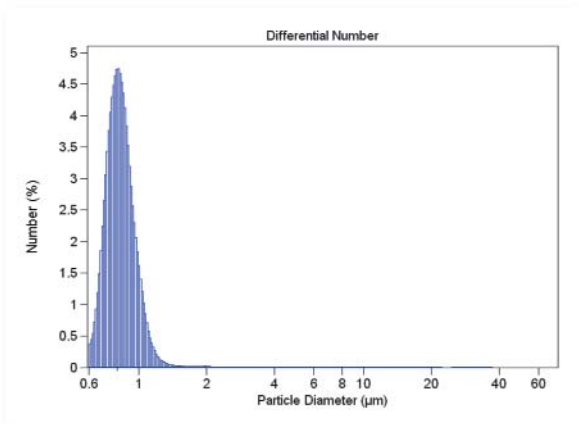


Results using a 30 µm aperture.

Defining which one is the larger size for the distribution of the normal cell population, it is possible to calculate the percentage of agglomerates. In the following example it was considered that above 3µm the cells are present in agglomerates. From this assumption the cell aggregation is 1.37% of the total population.



Results using a 100 µm aperture.



Combined results from the 30 µm and 100 µm aperture.

Particle Diameter µm	Number % >
0.6	100.00
1	23.76
2	3.23
3	1.37
5	1.18
10	0.84
15	0.63
25	0.49

By comparing the percentage of agglomerates present in a sample it is possible to control the filtration efficiency and therefore the potential effectiveness of the antimicrobial agent.

It is also possible to refine filtration process and the effectiveness of detergents in preventing cell clumps.

Visit us at our website: www.beckmancoulter.com

WORLD-CLASS GLOBAL SUPPORT: AUSTRALIA (61) 2 9844 6000 CANADA (1) 905 819 1234 CHINA (86) 106 515 6028 EASTERN EUROPE, MIDDLE EAST, AFRICA (41) 22 994 0707 FRANCE (33) 1 49 909083 GERMANY (49) 2151 333 625 HONG KONG (852) 2814 7431 ITALY (39) 2 25221 JAPAN (81) 3 5404 8359 LATIN AMERICA (1) (305) 380 3810 MEXICO (52) 55 5605 7770 NETHERLAND (31) 2972 30630 SINGAPORE (65) 63393633 SOUTH AFRICA (27) 11 805 2014 / 5 SPAIN (34) 91 728 7900 SWEDEN (46) 8 564 85900 SWITZERLAND 0800 050 810 TAIWAN (886) 2 2378 3456 U.K. (44) 1494 429162 U.S.A. (1) 800 523 3713

