

M I C R O B I A L C Y T O M E T R Y R E F E R E N C E S

General Microbiology

- Alvarez AM, Ibanez M, Rotger R. Beta-galactosidase activity in bacteria measured by flow cytometry. *Biotechniques*. 1993;15:974-976.
- Conrads G, Herrler A, Moonen I, Lampert F, Schnitzler N. Flow cytometry to monitor phagocytosis and oxidative burst of anaerobic periodontopathogenic bacteria by human polymorphonuclear leukocytes. *Journal of Periodontal Research*. 1999;34:136-144.
- Lebaron P, Catala P, Parthuisot N. Effectiveness of SYTOX green stain for bacterial viability assessment. *Applied and Environmental Microbiology*. 1998;64:2697-2700.
- Nebe-von-Caron G, Stephens PJ, Hewitt CJ, Powell JR, Badley RA. Analysis of bacterial function by multi-colour fluorescence flow cytometry and single cell sorting. *Journal of Microbiological Methods*. 2000;42:97-114.
- Porter J, Pickup R. Nucleic acid-based fluorescent probes in microbial ecology: application of flow cytometry. *Journal of Microbiological Methods*. 2000;42:75-79.
- Robinson JP. Microbiological Applications. *Current Protocols in Cytometry*. New York, NY: John Wiley and Sons; 2001:11.0.1-11.14.9.
- Roth BL, Poot M, Yue ST, Millard P. Bacterial viability and antibiotic susceptibility testing with SYTOX green nucleic acid stain. *Applied and Environmental Microbiology*. 1997;63: 2421-2431.
- Servais P, Courties C, Lebaron P, Troussellier M. Coupling bacterial activity measurements with cell sorting by flow cytometry. *Microbial Ecology*. 1999;38:180-189.
- Shapiro HN. Microbial analysis at the single-cell level: tasks and techniques. *Journal of Microbiological Methods*. 2000;42:3-16.
- Steen HB. Flow cytometric studies of microorganisms. *Flow Cytometry and Sorting*. 2nd Ed. New York, NY: Wiley-Liss, Inc; 1990: 605-622.
- Valdivia RH, Falkow S. Bacterial genetics by flow cytometry: rapid isolation of *Salmonella typhimurium* acid-inducible promoters by differential fluorescence induction. *Molecular Microbiology*. 1996;22:367-78.
- Zubkov MV, Fuchs BM, Eilers H, Burkill PH, Amann R. Determination of total protein content of bacterial cells by SYPRO staining and flow cytometry. *Applied and Environmental Microbiology*. 1999;65:3251-3257.

Clinical Microbiology

- Arrese JF, Goffin V, Avila-Camacho M, Greimers R, Piérard GE. A pilot study on bacterial viability in acne. Assessment using dual flow cytometry on microbials present in follicular casts and comedones. *International Journal of Dermatology*. 1998;37:461-464.
- Durodie J, Coleman K, Simpson IN, Loughborough SH and Winstanley DW. Rapid detection of antimicrobial activity using flow cytometry. *Cytometry* 1995;21:374-377.
- Gottfredsson M, Erlendsdóttir H, Sigfússon Á, Gudmundsson S. Characteristics and dynamics of bacterial populations during postantibiotic effect determined by flow cytometry. *Antimicrobial Agents and Chemotherapy*. 1998;42:1005-1011.

Environmental and Aquatic Microbiology

- Ferrari BC, Vesey G, Davis KA, Gauci M and Veal D. A Novel Two-Color Flow Cytometric Assay for the Detection of *Cryptosporidium* in Environmental Water Samples. *Cytometry* 2000; 41:216-222.
- Gasol JM, Zweifel UL, Peters F, Fuhrman JA, Hagström Å. Significance of size and nucleic acid content heterogeneity as measured by flow cytometry in natural planktonic bacteria. *Applied and Environmental Microbiology*. 1999;65:4475-4483.

Lebaron P, Parthuisot N, Catala P. Comparison of blue nucleic acid dyes for flow cytometric enumeration of bacteria in aquatic systems. *Applied and Environmental Microbiology*. 1998;64:1725-1730.

Marie D, Brussaard CPD, Thyrrhaug R, Bratbak G, Vaulot D. Enumeration of marine viruses in culture and natural samples by flow cytometry. *Applied and Environmental Microbiology*. 1999;65:45-52.

Porter J, Edwards C, Morgan JA, Pickup RW. Rapid, automated separation of specific bacteria from lake water and sewage by flow cytometry and cell sorting. *Applied and Environmental Microbiology*. 1993;59:3327-3333.

Servais P, Courties C, Lebaron P, Troussellier M. Coupling bacterial activity measurements with cell sorting by flow cytometry. *Microbial Ecology*. 1999;38:180-189.

Troussellier M, Courties, Lebaron P, Servais P. Flow cytometric discrimination of bacterial populations in seawater based on SYTO 13 staining of nucleic acids. *FEMS Microbiology Ecology*. 1999;29:319-330.

Vives-Rego J, Lebaron P, Nebe-von Caron G. Current and future applications of flow cytometry in aquatic microbiology. *FEMS Microbiology Reviews*. 2000;24:429-48.

Industrial, Food and Fermentation Microbiology

Arroyo G, Sanz PD, Préstamo G. Response to high-pressure, low-temperature treatment in vegetables: determination of survival rates of microbial populations using flow cytometry and detection of peroxidase activity using confocal microscopy. *Journal of Applied Microbiology*. 1999;86:544-556.

Borazjani RN, May LL, Noble JA, Avery SV, Ahearn DG. Flow cytometry for determination of the efficacy of contact lens disinfectant solutions against *Acanthamoeba* spp. *Applied and Environmental Microbiology*. 2000;66:1057-1061.

Gunasekera TS, Attfield PV, Veal DA. A flow cytometry method for rapid detection and enumeration of total bacteria in milk. *Applied and Environmental Microbiology*. 2000;66:1228-1232.

Wang X, Slavik MF. Rapid detection of *Salmonella* in chicken washes by immunomagnetic separation and flow cytometry. *Journal of Food Protection*. 1999;62:717-723.

General Yeast Biology

Bell PJJ, Deere D, Shen J, et al. A flow cytometric method for rapid selection of novel industrial yeast hybrids. *Applied and Environmental Microbiology*. 1998;64:1669-1672.

Deere D, Shen J, Vesey G, Bell P, Bissinger P, Veal D. Flow Cytometry and cell sorting for yeast viability assessment and cell selection. *Yeast*. 1998;14:147-160.

Virus

Marie D, Brussaard CPD, Thyrrhaug R, Bratbak G, Vaulot D. Enumeration of marine viruses in culture and natural samples by flow cytometry. *Applied and Environmental Microbiology*. 1999;65:45-52.



recycled



FM 32438

For Research Use Only. Not for use in diagnostic procedures.

© 2002 Becton, Dickinson and Company

23-6290-00 4/2002

BD Biosciences www.bdbiosciences.com

Clontech	Discovery Labware	Immunocytometry Systems	PharMingen			
Asia Pacific Tel (65) 6-860-1437 Fax (65) 6-860-1593	Brazil Tel (55) 11 5185-9995 Fax (55) 11 5185-9895	Canada Tel (888) 259-0187 Fax (905) 542-9391	Europe Tel (32) 53-720-211 Fax (32) 53-720-450	Japan Tel 0120-8555-90	Mexico Tel (52) 5-284-8281 Fax (52) 5-284-8288	United States Tel (877) 232-8995 Fax (408) 954-2347

BD Biosciences

Clontech
Discovery Labware
Immunocytometry Systems
PharMingen

