

Biological Applications for Multisizer. List of available articles.

Aspirin Analogues and Flow of Erythrocytes through Narrow Capillaries, Rao P.R., Jones S., Sun Z., Hosmane R.S., Abugo O. and Rifkind J., Dept. Mech. Eng. and Dept. Chem. & Biochem., Univ. Maryland, Baltimore, MD, and Gerontol. Res. Center, Nat. Inst. on Aging, NIH, Baltimore, MD, *Clin.Haemorheol.* , 15, No. 6, 877-887, 1995

Erythroid maturation and globin gene expression in mice with combined deficiency of NF-E2 and nrf-2, Martin F., van Deursen J.M., Shivdasani R.A., Jackson C.W., Troutman A.G. and Ney P.A., Department of Biochemistry, Saint Jude Children's Research Hospital, Memphis, TN, USA, *Blood* , 91, No. 9, 3459-3466, 1998.

Solvent/Detergent-Treated Plasma Suppresses Shear-Induced Platelet Aggregation and Prevents Episodes of Thrombotic Thrombocytopenic Purpura, Moake J., Chintagumpala M., Turner N., McPherson P., Nolasco L., Steuber C., Santiago-Borrero P., Horowitz M. and Pehta J., Baylor Coll. Med., and Rice Univ., Houston TX; Univ. Puerto Rico Med. Sch., San Juan, Puerto Rico; and New York Blood Center, *Blood* , 84, No. 2, 490-497, 1995

Photodynamic Treatment of Yeast Cells with the Dye Toluidine Blue : All-or-None Loss of Plasma Membrane Barrier Properties, Paardekooper M., Ven den Broek P.V.A., De Bruijne A.W., Elferink J.G.R., Dubbelman T.M.A.R. and Van Steveninck J., Dept. of Med. Biochem., State Univ. Leiden, The Netherlands, *Biochim.Biophys.Acta* , 1108, 86-90, 1992

Particle Release from Haemodialysers, Hoenich N.A., Thompson J., McCabe J. and Appleton D.R., Dept. of Med., School of Clin. Med. Sci., The Med. School; Dept. of Prosthodontics, The Dental School; and Div. of Med. Stat., School of Health Care Sci., The Med. School; The Univ., Newcastle upon Tyne, *Intl.J.Artif.Organs* , 13, No. 12, 803-808, 1990

Novel Method for Assessing Postantibiotic Effect by Using the Coulter Counter, Li R.C., Lee S.W. and Lam J.S., Depts. Pharm. and Ped., Fac. Med., Chinese Univ. Hong Kong, *Antimicrobial Agents and Chemotherapy* , 40, No. 7, 1751-1753, 1996

Assessment of the Cytocompatibility of Biomaterials by Analysis of Cellular Viability in a Coulter Multisizer Multichannel Analyser, Roux H., Duval J.L., Sigot-Luizard M.F. and Sigot M., Univ. de Tech. de Compiègne - Lab. de Biol. Cell. Exp., Compiègne, France, *Advances in Biomaterials* , 10, 81-87, 1992.