

MoFlo™ Chromosome Sorting

Introduction

Flow cytometric chromosome classification^{1,2} (flow karyotyping) has been routinely used in the analysis of mammalian karyotypes as well as to assess chromosomal abnormalities.^{3,4} Various chromosome isolation methods⁵⁻⁷ have been developed in the preparation of isolated chromosomes from mitotic cells for flow karyotyping or flow sorting. To generate a good chromosome preparation with minimal debris formation and large numbers of individual chromosomes for a well-resolved flow karyotype, protocols have made use of stabilizing agents such as divalent cations⁷ (magnesium ions) and cationic polyamines⁵ (spermine, spermidine) in the isolation buffer. Univariate⁸⁻¹⁰ or bivariate^{11,12} flow karyotype analysis on the isolated chromosomes can be performed by labeling with one (propidium iodide, ethidium bromide) or two DNA specific fluorescent dyes (Hoechst 33258, Chromomycin A3). Highly purified fractions of chromosomes can then be isolated rapidly from the labeled chromosomes using the MoFlo high performance cell sorter.

Materials and Methods

Chromosome preparation and staining. Block cells at metaphase using 0.1 µg/ml demecolcine for an optimized amount of time to maximize the percentage of mitotic cells (e.g. ~6 hours for a lymphoblastoid cell line). Pellet (289 g, 5 min) and resuspend cells in hypotonic KCl buffer for 10 mins at room temperature. Pellet (289 g, 5 min) and resuspend cells in ice cold polyamine isolation buffer⁵ and vortex for 20 seconds. Spin the chromosome suspension at 200 g for 2 mins, filter the supernatant through 20 µm mesh filter and stain overnight with 5 µg/mL Hoechst 33258, 40 µg/mL Chromomycin A3 and 10 mM MgSO₄. Add 10 mM of sodium citrate and 25 mM of sodium sulphite to the stained preparation one hour before flow analysis.

Instrument set up. Analyze the stained chromosome suspension on a MoFlo equipped with two water-cooled argon ion lasers spatially

separated at the flow chamber. Tune the first laser to emit multi-line UV (330-360 nm) to excite Hoechst 33258 and the second laser to emit light at 457 nm to excite Chromomycin A3. Maintain the power of both lasers at 300 mW. Trigger the signal based on Hoechst 33258 fluorescence and collect Forward Scatter (FSC) using a 351/10 nm band pass filter, Hoechst 33258 fluorescence using a 400 nm long pass filter combined with a 480 nm short pass filter and Chromomycin A3 fluorescence using a 490 nm long pass filter. Prealign the MoFlo for minimum peak coefficient of variance for both fluorescence parameters and forward scatter using 3 µm fluorescent beads. Configure the MoFlo for high speed sorting with optimal settings of the sheath pressure to ~60 psi and the drop drive frequency to ~95 KHz using a 70 µm Cytonozzle tip. Select the high purity sort option of single mode per single drop envelope on a 4WAY™ Sort.

Flow analysis and sorting. For data analysis, acquire a total of 100,000 events at a data rate of 1,000 events per second and analyze on Summit Software. Display data as a bivariate flow karyogram of Hoechst 33258 versus Chromomycin A3 fluorescence (Figure 1) after gating on low forward scatter and high Hoechst 33258 fluorescence to exclude some debris and clumps.

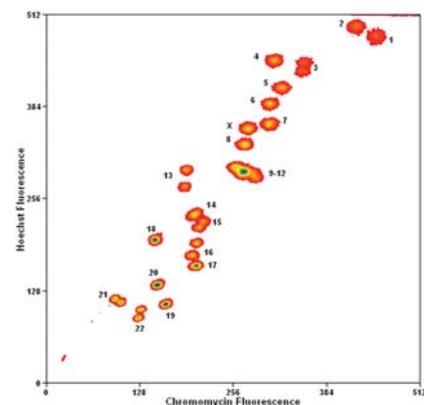


Figure 1

Bivariate flow karyotype of a normal human cell line. The chromosomes are stained with Hoechst 33258 and Chromomycin A3. The position of each chromosome type is indicated. In Summit Software, for analysis, enable smoothing. Smoothing method: Gaussian, # of Passes = 2.

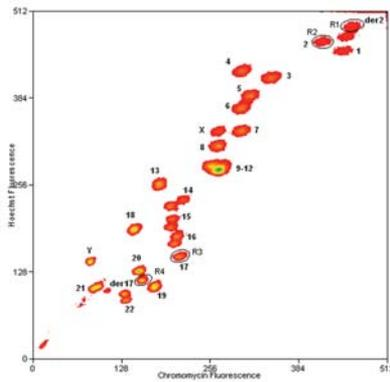


Figure 2

Bivariate plot of Hoechst 33258 vs. Chromomycin A3 fluorescence of a human cell line with translocation between chromosome 2 and chromosome 17. Sort gates are drawn on the translocated chromosomes, derivative 2 (R1) and 17 (R4) and its normal counterpart, chromosome 2 (R2) and 17 (R3). In Summit Software, for analysis, enable smoothing. Smoothing method: Gaussian, # of Passes = 2.

For chromosome sorting, set up the sort decision using the Sort Logic Editor by creating sort gates on the selected chromosome clusters on a bivariate flow karyogram (Figure 2) and isolate the required number of chromosomes into the collection vials.

Discussion

Flow karyotyping and purification of chromosomes isolated from mammalian cell lines can be achieved using the MoFlo high performance cell sorter. The sorting set up of the MoFlo can be customized for rapid isolation of chromosomes with high purity. With the MoFlo, chromosomes can be flow sorted at a data rate of approximately 8,000-12,000 chromosomes per second while still maintaining a good resolution of chromosome peaks. Following purification, DNA as well as proteins can be extracted from the flow sorted chromosomes for genomics¹³⁻¹⁵ and proteomics¹⁶ studies.

References

- Gray JW et al. High-speed quantitative karyotyping by flow microfluorometry. *Clin Chem* (21) 1975; 1258-62.
- Stubblefield E, Cram, S. & Deaven, L. Flow microfluorometric analysis of isolated Chinese hamster chromosomes. *Exp Cell Res* (94) 1975; 464-8.
- Gray JW et al. Application of flow karyotyping in prenatal detection of chromosome aberrations. *Am J Hum Genet* (42) 1988; 49-59 (1988).
- Gray JW and Langlois RG. Chromosome classification and purification using flow cytometry and sorting. *Annu Rev Biophys Chem* (15) 1986; 195-235.
- Sillar R and Young BD. A new method for the preparation of metaphase chromosomes for flow analysis. *J Histochem Cytochem* (29) 1981; 74-8.
- Bijman JT. Optimization of mammalian chromosome suspension preparations employed in a flow cytometric analysis. *Cytometry* (3) 1983; 354-8.
- Van den Engh G, Trask B, Cram S and Bartholdi M. Preparation of chromosome suspensions for flow cytometry. *Cytometry* (5) 1984; 108-17.
- Cram LS, Bartholdi MF, Ray FA, Travis GL and Kraemer PM. Spontaneous neoplastic evolution of Chinese hamster cells in culture: multistep progression of karyotype. *Cancer Res* (43) 1983; 4828-37.
- Green DK et al. Karyotyping and identification of human chromosome polymorphisms by single fluorochrome flow cytometry. *Hum Genet* (66) 1984; 143-6.
- Young BD, Ferguson-Smith MA, Sillar R and Boyd E. High-resolution analysis of human peripheral lymphocyte chromosomes by flow cytometry. *Proc Natl Acad Sci USA* (78) 1981; 7727-31.
- Van den Engh, GJ, Trask BJ, Gray JW, Langlois RG and Yu LC. Preparation and bivariate analysis of suspensions of human chromosomes. *Cytometry* (6) 1985; 92-100.
- Langlois RG, Yu LC, Gray JW and Carrano AV. Quantitative karyotyping of human chromosomes by dual beam flow cytometry. *Proc Natl Acad Sci USA* (79) 1982; 7876-80.
- Fiegler H et al. Array painting: a method for the rapid analysis of aberrant chromosomes using DNA microarrays. *J Med Genet* (40) 2003; 664-70.
- Gribble SM et al. Applications of combined DNA microarray and chromosome sorting technologies. *Chromosome Res* (12) 2004; 35-43.
- Gribble S, Ng BL, Prigmore E, Burford, DC and Carter NP. Chromosome paints from single copies of chromosomes. *Chromosome Res* (12) 2004; 143-51.
- Wray W and Wray VP. Proteins from metaphase chromosomes treated with fluorochromes. *Cytometry* (1) 1980; 18-20.

For Research Use Only. Not for use in Diagnostic Procedures.

Australia, Gladesville (61) 2 9844-6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028
 Czech Republic, Prague (420) 272 01 73 32 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707
 France, Villepinte (33) 1 49 90 90 00 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431
 India, Mumbai (91) 22 3080 5101 Italy, Cassina de' Pecchi, Milan (39) 02 953921 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146
 Latin America (1) (305) 380 4709 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Mijdrecht (31) 297 230630 Puerto Rico (787) 747 3335
 Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 564 3203 Spain, Madrid (34) 91 3836080
 Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2378 3456 Turkey, Istanbul (90) 216 309 1900
 UK, High Wycombe (44) 01494 441181 USA, Fullerton, CA (1) 800 742 2345