

CyAn™ ADP Detection of Soluble Cytokines Using Multiplexed Beads Assays

Introduction

Multiple soluble molecules in body fluids or culture supernatants can be measured by Flow Cytometry using multiplex bead-based assays. This method requires a smaller sample volume (about 25 μ L) and has a higher sensitivity and speed compared to other techniques, such as ELISA.¹

This application note illustrates the use of the CyAn ADP multicolor flow cytometer to acquire 11-plex samples prepared with the Bender MedSystems FlowCytomix Kit used on the culture supernatant of a CD8+ T cell line, HLA-A*0201-specific, obtained with repeated NV9 peptide (part of CMV pp65 antigen, NLVPMVATV) pulsing and IL-2 stimulation (see Figures 2 and 3).

This kit uses two sets of size-coded beads (5,5 microns shown in R1, and 4,4 microns, shown in R2) composed of 5 / 6 different beads subsets, fluorescence-coded by different intensities of the Starfire Red** dye (excitation from UV to red, emission max at 675nm). Each bead is coated with a different antibody, specific for a single analyte.

The analyte concentration is measured staining the cytokine + bead complex with a secondary, biotin-conjugated antibody, followed by streptavidin-PE staining.

Different PE intensities represent different analyte concentrations.

The kit includes a set of standards for each analyte. Thus, the measured intensities can easily be translated into concentration units.

Materials and Methods

A Th1/Th2 11-plex Kit was used to provide the concentration of IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, TNF- α , TNF- β , IFN- γ .

The procedure method in summary is:

- Step 1: Incubate beads, sample and biotin conjugates
- Step 2: Wash
- Step 3: Incubate with PE-streptavidin
- Step 4: Wash
- Step 5: Acquire sample on CyAn ADP 9 Color with Summit software v4.3.

Results

FSC and SSC were used to identify the two bead populations (Figure 1: R1 and R2). Two plots with FL2 (PE) vs. FL8 (Starfire Red) were created and gated on the two beads sets to separate the different Starfire Red levels, each representing an analyte, and to monitor the PE mean fluorescences

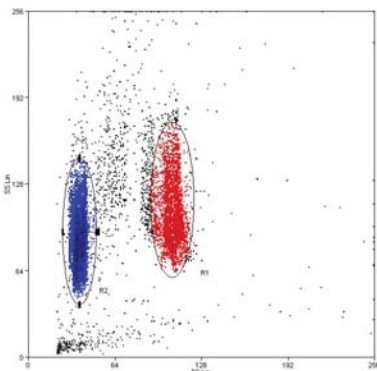


Figure 1

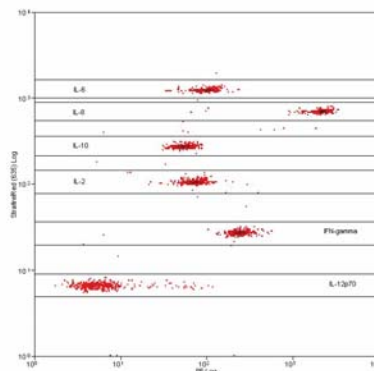


Figure 2

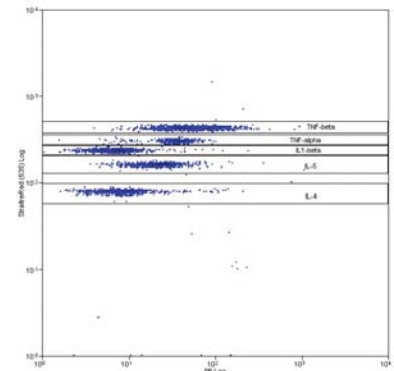


Figure 3

(Figures 2 and 3) representing the analyte concentration.

The FL8 channel gave the necessary resolution and compensation was not required between Starfire Red and PE. For each analyte the required minimum of 300 beads was acquired, 5,000 beads were collected in total.

Discussion

As shown, CyAn ADP can easily resolve the two different sized beads and analytes. A dedicated software (Bender FCMix Pro 2.1, included in the Kit) automatically calculates the cytokine concentration (pg/mL) by defining and applying 11 reference curves generated with corresponding 11 Standards set (e.g. for IL-4, see Figure 4).

Note

Please refer to the manuals and website provided by Bender MedSystems for detailed sample preparation and analysis procedures.

References

1. Paul WE and Seder RA. Lymphocyte responses and cytokines. *Cell* 76: 241-51 (1994).
2. Rooney JW et al. A common factor regulates both Th1- and Th2-specific cytokine gene expression. *EMBO Journal* 13:625-33 (1994).
3. Lederer JA et al. Regulation of cytokine gene expression in T helper cell subsets. *Journal of Immunology* 152: 77-86 (1994).

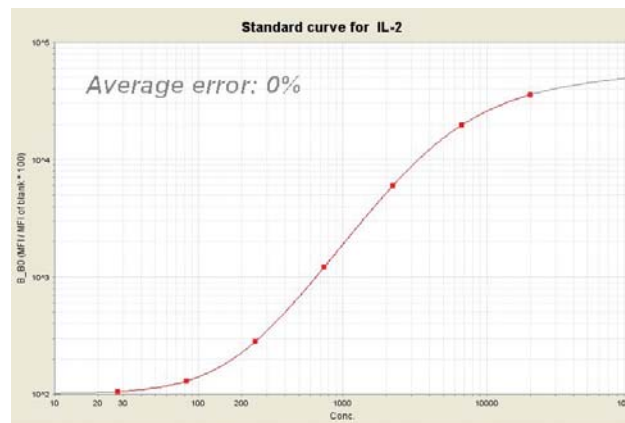


Figure 4

Contributors

Dr. Federico Ivaldi, D.ssa Giuseppina Li Pira – Unità di Immunologia Virale, Centro di Biotecnologie Avanzate – Genova, Italy.

Technical Tips

- Beads agglutination may occur in lipemic samples. Centrifugation of lipemic samples (5 minutes, 16,000 rcf) before analysis may help creating the needed sample fraction located between the pellet and the lipemic layer at the top.

**Starfire Red is a registered trademark of Bangs Laboratories, Inc.

For Research Use Only. Not to be used 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