

MoFlo ASTRIOS^{EQ}™ FORWARD SCATTER: CELL SORTING OF NANO AND LARGE PHYTOPLANKTON

SIMULTANEOUSLY WITH HIGH PURITY

CARLEY D. ROSS PH.D., AND ROBIN MORRIS BECKMAN COULTER LIFE SCIENCES, FLOW CYTOMETRY BUSINESS CENTER, FORT COLLINS, CO



B158

ABSTRACT

BACKGROUND: Phytoplankton conversion of light on the upper limits of the ocean consists of half of the photosynthesis on the Earth. The population densities indicate the health of not only the phytoplankton, but the entire aquatic ecosystem. The isolation and sorting of aquatic samples using flow cytometry allows for quick and effective population analysis. The forward scatter on the MoFlo Astrios allows for differentiation of small and large particles from 0.2 to 30 µm on FSC. This design provides researchers greater flexibility to isolate and sort specific phytoplankton of different sizes while utilizing the 7 laser, 42 parameter MoFlo Astrios.

METHODS: Plankton species; *Chlorella*, *Phormidium inunatum*, *Phormidium persicinum*, *Cryptomonas*, *Rhodospira*, *Synechococcus*, *Skeletonema*, *Fremyella*, were acquired from the UTEX: The Culture of Algae and grown in photobioreactors and cultured in specialized salt and fresh water media. Plankton species *Prochlorococcus marinus* and *Emiliana huxleyi* were grown in sterile 2 L containers supplied with 0.2 µm filtered air in salt water with fertilizer. Instant Ocean, 1/2 cup per gallon of deionized water, was added to the culture with Microalgae Grow Mass Pack with Silicate. Cells were harvested by gentle centrifugation, roughly 300 x g for 5-10 minutes. The supernatant was decanted/aspirated and the pellet resuspended in a sterile saline solution to achieve 1x10⁶ cells/mL. Plankton were then stained with SYTOX Green (Invitrogen, S7020), at a maximum concentration of 5 µM for 20 minutes after vortexing. Samples were filtered with a 70 µm Partec filter and kept on ice before flow cytometric analysis.

FLOW CYTOMETRY: The MoFlo Astrios^{EQ} was configured with 7 lasers and setup with a 100 µm tip to accommodate the larger phytoplankton (*Cryptomonas*). Cells were selected on their "live" status by being highly fluorescent in the red channels (chlorophyll) and low in the green channels (Sytox -). For small particle analysis, *Prochlorococcus* and *Synechococcus* were simultaneously analysed on FSC-Log parameters. Populations were sorted based on fluorescence and size as a 6-way sort into 5 mL tubes. The plankton were sorted at 25K eps to collect at least 100,000 events per each population using sort mode Purify 1-2.

RESULTS: Plankton populations were distinguishable using fluorescence and scatter patterns on both log and linear scales simultaneously. With the Astrios optical flexibility, the plankton fluorescence spectra were optimized for signal to noise. Isolation of the *Prochlorococcus*, *Synechococcus* and other plankton species using cell sorting achieved 99% purity for all populations.

CONCLUSIONS: The forward scatter and optical collection design of the MoFlo Astrios^{EQ} provide flexibility to detect large and small populations and sort them with high purity.

INTRODUCTION

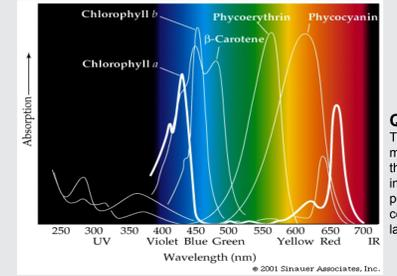
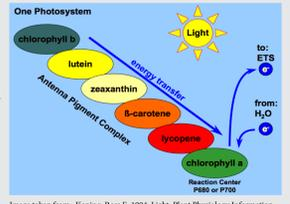
Marine biology, specifically photosynthetic plankton, is another sorting application for the new MoFloTM Astrios^{EQ} forward scatter design. Flow cytometry has previously been used with phytoplankton for rapid and objective evaluation of water quality (Trask), to analyze grazing (Cucco), cell viability (Dorsey), rapid strain identification (Simon), diversity assessment (Marie D), and to characterize plankton populations through molecular (Shi, Marie D) and genetic approaches (Shi).

Photosynthetic Pigments Found in Phytoplankton

Using the flow cytometer to analyze plankton is useful in the ability to identify plankton populations as well as sort the populations for further population identification. The high speed sort capabilities and small particle detection of the MoFlo Astrios^{EQ} expand the researchers capabilities to identify and isolate phytoplankton populations.

PLANT PIGMENTS ABSORPTION SPECTRA

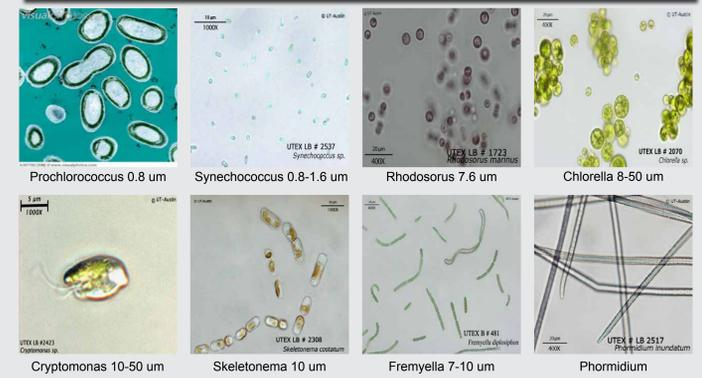
In a general photosystem, as seen in the figure to the right, the chlorophyll (b) transfers energy through the antenna pigment complex to chlorophyll (a) reaction center. The accessory pigments, such as Phycocyanin, Allophycocyanin, Allophycocyanin B, and Phycocerythrin, provide a composite absorption spectrum such that a wider range of visible and infrared radiation is absorbed by plants and algae.



QUANTITY OF PIGMENTS IN PHYTOPLANKTON

The quantity and concentration of the plant pigments found in the phytoplankton vary depending on the organism size and oceanic environment (including the habitat depth from ocean surface). Plankton populations may be distinguished by their fluorescence spectra using flow cytometry using multiple laser lines and emission filters.

PLANKTON

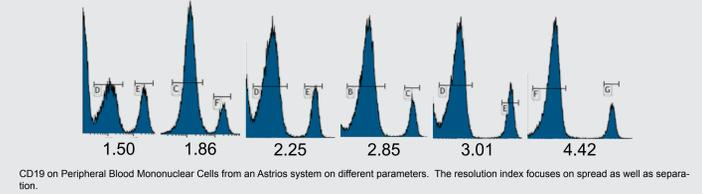


Fisher Distance

Quantifies the separation between a stained and unstained population in terms of the population widths

$$\text{Fisher Distance} = \frac{|Mean_1 - Mean_2|}{SD_1 + SD_2}$$

Requirements: Median of auto-fluorescent population needs to be above the baseline

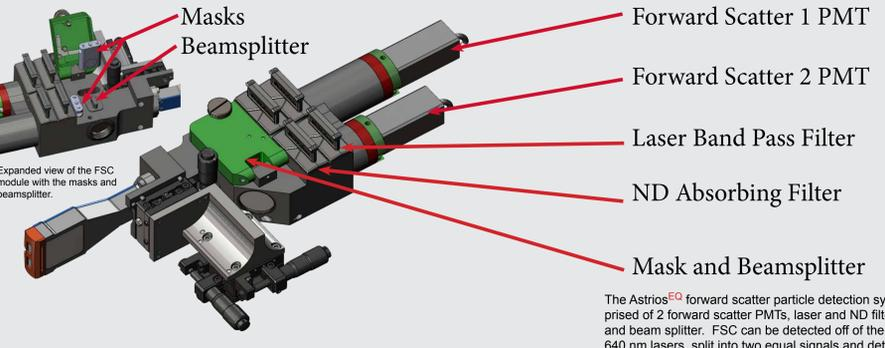


FLUORESCENCE

The Fisher Distance was used to determine the optimum excitation laser and emission bandpass using Kaluzi 1.2 Software. The Kaluzi software was used to find the initial optimum median channel and then used to calculate the Fisher Distance in the analysis package for quick data analysis. Below is an example of the fluorescence of the *Cryptomonas* plankton (left) and the resulting Fisher Distance of all of the plankton analyzed in Kaluzi (right). The red indicates parameters of poor resolution and blue indicates parameters of high separation.

Fluorescence Parameter	Chlorella	Cryptomonas	E. huxleyi	Fremyella	Phormidium inunatum	Phormidium persicinum	Phormidium unguiculatum	Prochlorococcus	Rhodospira	Skeletonema	Synechococcus
488-576/21	0.01	1.96	0.01	1.18	0.01	0.01	0.01	2.64	0.09		
488-620/29	0.01	1.96	0.01	1.18	0.02	0.01	0.01	2.89	0.09		
488-664/22	2.53	2.07	3.23	1.32	3.83	0.08	3.11	2.97	2.01		
488-710/45	2.12	1.80	3.18	0.68	1.69	0.40	1.01	2.66	2.64	1.69	
488-750/70	1.84	1.77	3.07	0.97	1.71	0.13	2.60	2.42	2.42	1.79	
512-576/21	0.01	1.73	0.02	0.02	0.05	0.01	0.01	3.24	0.01		
512-622/22	0.01	1.56	0.01	1.63	0.01	0.03	0.01	2.96	0.01		
512-664/21	2.90	1.98	3.03	1.04	2.02	0.53	0.69	3.07	3.14	2.99	
512-692/18	2.29	1.41	3.11	1.07	1.95	1.00	1.06	1.34	2.99	2.72	1.50
512-736/47	1.87	1.35	2.88	1.00	1.94	0.56	0.89	1.37	2.71	2.55	1.52
565-576/16	0.01	0.09	0.05	0.58	1.75	0.01	0.03	2.98	0.01		
565-614/20	0.01	0.09	0.05	0.58	1.66	0.13	0.39	1.23	2.83	3.15	
565-692/75	3.25	1.53	2.90	1.07	1.98	0.54	0.64	1.92	3.10	2.89	2.36
592-620/29	0.02	2.02	3.35	1.63	1.72	3.18	1.71	2.74	2.99	0.01	
592-671/47	4.42	1.72	2.95	1.17	1.94	0.54	0.74	2.36	3.14	2.99	0.01
592-722/44	4.42	1.59	2.81	1.20	2.06	1.11	1.18	2.20	2.84	2.63	0.15
592-795/70	2.58	1.88	3.13	1.46	2.01	0.45	0.72	2.26	3.17	2.96	0.26
640-671/30	5.21	1.88	3.03	0.99	1.89	0.25	0.59	2.26	3.51	3.09	2.61
640-722/44	4.42	1.73	2.86	1.10	1.84	0.47	0.95	1.99	3.13	2.77	0.15
640-795/70	2.76	1.95	2.98	1.18	1.81	0.93	1.30	2.17	2.97	2.59	0.26

FSC ENHANCED PARTICLE DETECTION SYSTEM



Two key components to the small particle detection are the air and sheath filters which eliminate the small particle debris. It is important to be aware of the sample media, tubes and other plastic components that may cause submicron background and blur the small particle resolution.

The new Astrios^{EQ} has a new FSC small particle detection system that allows users to customize the mask and laser FSC onto two FSC PMTs. The dual FSC parameter analysis gives the unique opportunity to view populations and isolate populations based on those scatter patterns as seen on the right with the *Synechococcus* sort.

OPTIMIZING FSC

MASKS

M - Material

- Better relation to particle size over large ranges of particles
- Designed to pull apart particles of different material types

S - Separation

- Reduce the sensitivity to different particle types
- Better for pulling out small differences in similar material types

P - All-Purpose

- Overall general performance
- Best when matched against "M" or "S" masks to pull out details

Counts (Histograms for M1, S1, P1)

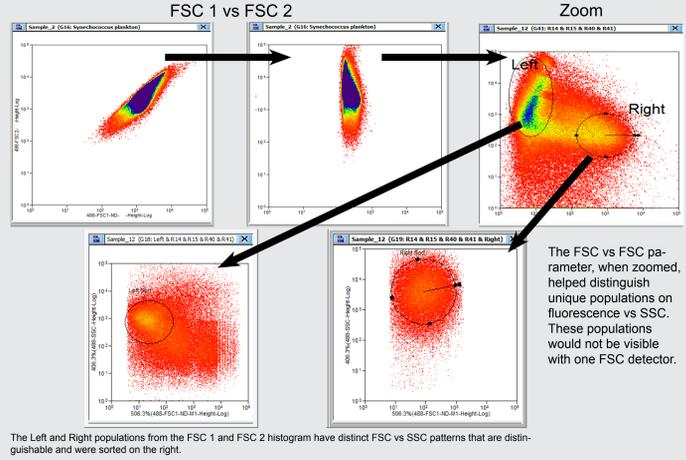
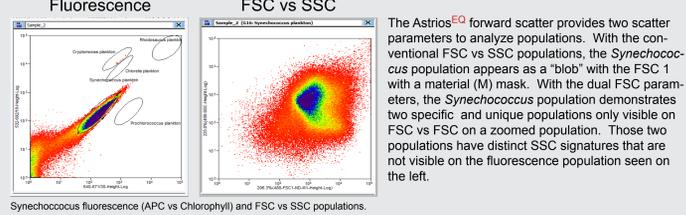
FSC Height (Histograms for S1, S2)

Above: Live/Dead separation of Mouse Spleen using S Masks.

Left: Using M masks and ND filters to visualize 0.2 to 30 µm polystyrene beads.

Forward scatter with whole blood on both FSC 1 and FSC 2 with P masks.

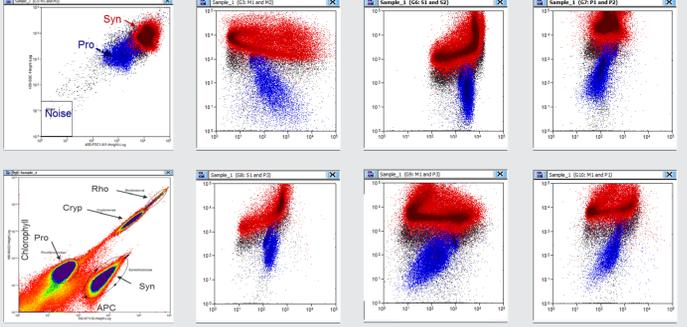
DUAL FSC USED TO FIND UNIQUE POPULATIONS



The Left and Right populations from the FSC 1 and FSC 2 histogram have distinct FSC vs SSC patterns that are distinguishable and were sorted on the right.

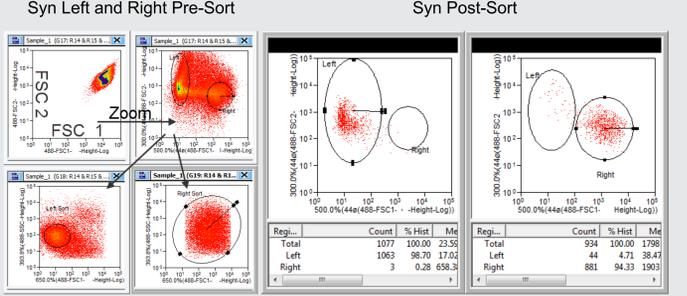
RESULTS

MASK SELECTION OF PRO AND SYN



The *Synechococcus* and *Prochlorococcus* cyanobacteria have distinct fluorescence characteristics and they were used to color gate to find unique populations of each plankton. Note the "Noise" gate was very far from the submicron plankton populations.

5-WAY HIGH SPEED SORT



The sort contained populations of *Rhodospira*, *Cryptomonas*, *Prochlorococcus* and the two populations found in the *Synechococcus* (as seen above). The sort was run at 50K EPS on Purify 1-2 into 5 mL tubes. The purity was: Syn Right - 94.33%, Syn Left - 98.70%, *Prochlorococcus* 98.56% (left), *Rhodospira* 98.59% (not pictured) and *Cryptomonas* 99.54%.

CONCLUSIONS

- The MoFlo Astrios^{EQ} new FSC module provided the ability to sort unique FSC populations for the *Synechococcus* population with high purity and high speed.
- The FSC module was able to distinguish from noise two submicron populations (*Prochlorococcus* and *Synechococcus*).
- The plankton populations were sorted with high purity and speed, indicating the robust nature of the MoFlo Astrios^{EQ} sorting system.

REFERENCES

Trask BJ, et al. Analysis of phytoplankton by flow cytometry. *Cytometry*. 1982 Jan;2(4):258-64.

Cucco TL, et al. Using phytoplankton and flow cytometry to analyze grazing by marine organisms. *Cytometry*. 1989 Sep;10(5):659-69.

Dorsey L, et al. Rapid analytical technique for the assessment of cell metabolic activity in marine microalgae. *Cytometry*. 1989 Sep;10(5):622-8.

Simon N, et al. Fluorescent in situ hybridization with rRNA-targeted oligonucleotide probes to identify small phytoplankton by flow cytometry. *Appl Environ Microbiol*. 1995 Jul;61(7):2509-13.

Marie D, et al. Use of flow cytometric sort to better assess the diversity of small photosynthetic eukaryotes in the English Channel. *FEMS Microbiol Ecol*. 2010 May;72(2):165-78.

Shi XL, et al. Groups without cultured representatives dominate eukaryotic picoplankton in the oligotrophic South East Pacific Ocean. *PLoS One*. 2009 Oct;29(4):10:e7657.

Marie D, et al. Eukaryotic picoplankton communities of the Mediterranean Sea in summer assessed by molecular approaches (DGGE, TTGE, QPCR). *FEMS Microbiol Ecol*. 2006 Mar; 35(3):403-15.

Shi XL, et al. Plastid 16S rRNA gene diversity among eukaryotic picoplankton sorted by flow cytometry from the South Pacific Ocean. *PLoS One*. 2011 Apr; 28; 6(4):e18979.

*** FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

