

# BD Accuri™ C6 Flow Cytometer

Nil Emre, PhD

BD Biosciences

R&D Manager, Stem Cells and Cell  
Function



# The BD Accuri C6 Flow Cytometer System

- An affordable, full-featured, easy-to-use flow cytometer
- Two lasers and six detectors



# The BD Accuri C6 Flow Cytometer

---

## Innovations in all the major components of a flow cytometer

- **Fluidics:** Peristaltic pumps and pulse dampeners allow miniaturization and direct-volume measurement
- **Optics:** Locked-down alignment
- **Signal detection:** Broad dynamic range obviates voltage adjustments
- **Software:** Intuitive



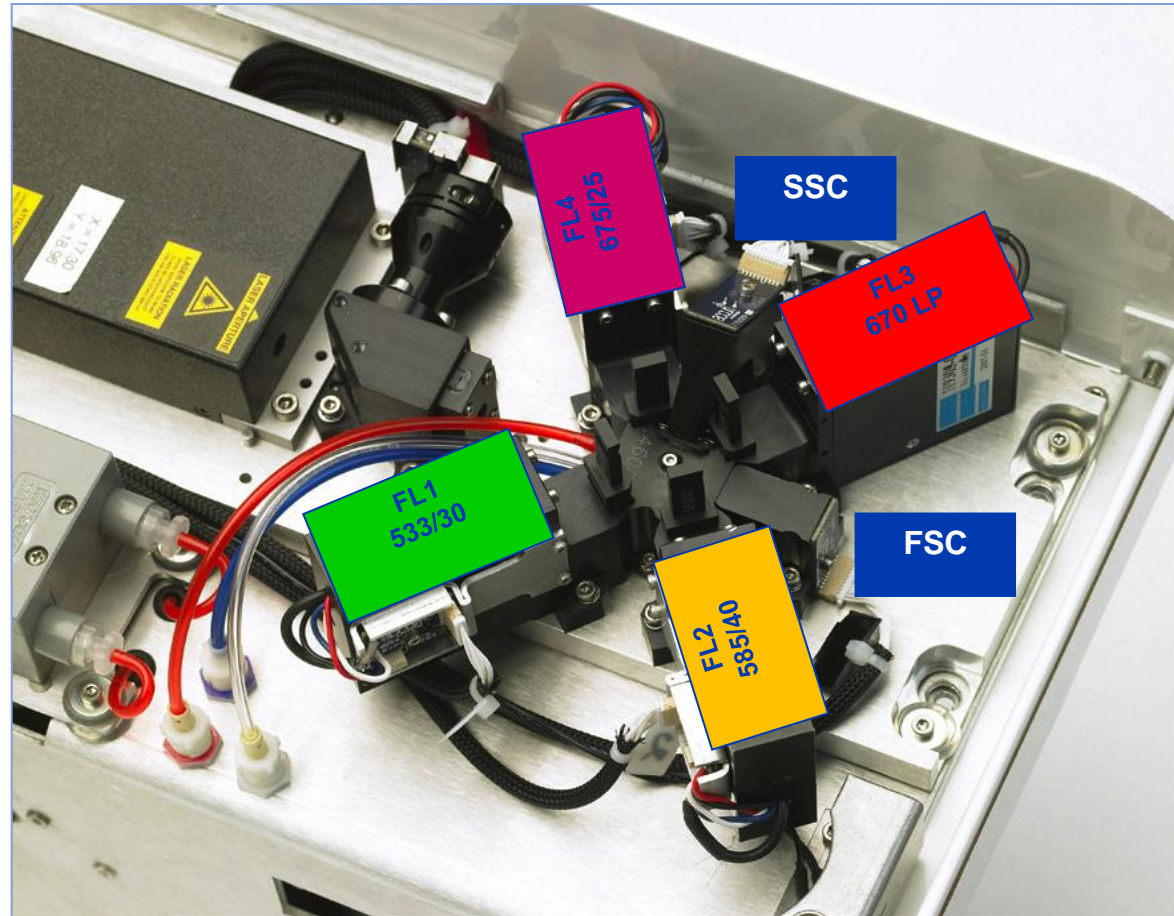
# Alignment and Signal Detection are Optimized and Locked Down

488-nm  
solid state laser

640-nm diode  
laser

PMTs for  
fluorescence  
detection

Diodes for light  
scatter detection



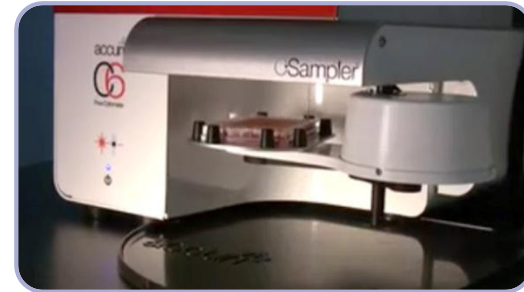
# Advantages of Pre-optimized Detector Settings

---

- Greatly reduces the risk of lost data due to improper setup
  - Saves time and sample
- No specialist training or dedicated operator required
- Predictable, reproducible analysis relative to the sample type and application
- Predictable fluorescence spillover

# Enhanced Sample Handling

- Direct volume measurement
- Many types of sample tubes may be used.
  - Flow cytometry tubes
  - Microcentrifuge tubes
  - Ninety-six-well plates with the BD CSampler™ accessory
- Open system conducive to kinetic studies
- BD CSampler™ accessory for automated sample introduction



# Intuitive Software

Sample Grid

Cytometer Status

Fluidics Controls

Run Criteria

Real-Time Updates

Collect
Analyze
Statistics

**A4** HPB CD3-F CD4, CD45, CD8

	1	2	3	4	5	6	7	8	9	10	11	12
A	A1	A2	A3	A4	A5	A6	A7	A8	A9	A10	A11	A12
B	B1	B2	B3	B4	B5	B6	B7	B8	B9	B10	B11	B12
C	C1	C2	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D	D1	D2	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E	E1	E2	E3	E4	E5	E6	E7	E8	E9	E10	E11	E12
F	F1	F2	F3	F4	F5	F6	F7	F8	F9	F10	F11	F12
G	G1	G2	G3	G4	G5	G6	G7	G8	G9	G10	G11	G12
H	H1	H2	H3	H4	H5	H6	H7	H8	H9	H10	H11	H12

C6 is connected and ready.

**Run Limits**

Run Unlimited

125000 events

in **Ungated Sample**

0 Min  0 Sec

0.0  $\mu$ L

**Backflush** **Unlog**

**Fluidics**

Slow  Medium  Fast

Flow Rate 14  $\mu$ L/min

Core Size 10  $\mu$ m

Custom

Flow Rate 11  $\mu$ L/min

**Set Core Size**

Core Size 5  $\mu$ m

**Threshold**

**Set Threshold**

None

80,000 on FSC-H

ADD to A4

Set Color Compensation

**Last Run**

0 Events

0.00 Time

0.0 Microliters

0 Events / Sec

0.0 Events /  $\mu$ L

**Cumulative**

229,303

0.00

0.0

0

0.0

**Delete Sample Data**

Warn before deleting

Data Capacity Used

8% of 10,000,000 Events

Histogram, Dot Plot, and Density Plot Display Area

Analysis and Gating Tools

Plot Statistics

# Kits and Templates on the BD Accuri C6

---

- “As Easy as Cell Analysis Is Going to Get” by combining Cell Biology Kits with acquisition and analysis templates.
- Free downloadable BD Accuri C6 Software Templates are matched to each kit at:
  - [www.bdbiosciences/go/templates](http://www.bdbiosciences/go/templates)
- Kits and Templates lead to quick and easy set-up and analysis of cell populations.



# Kits and Templates on the BD Accuri C6

## Available Templates:

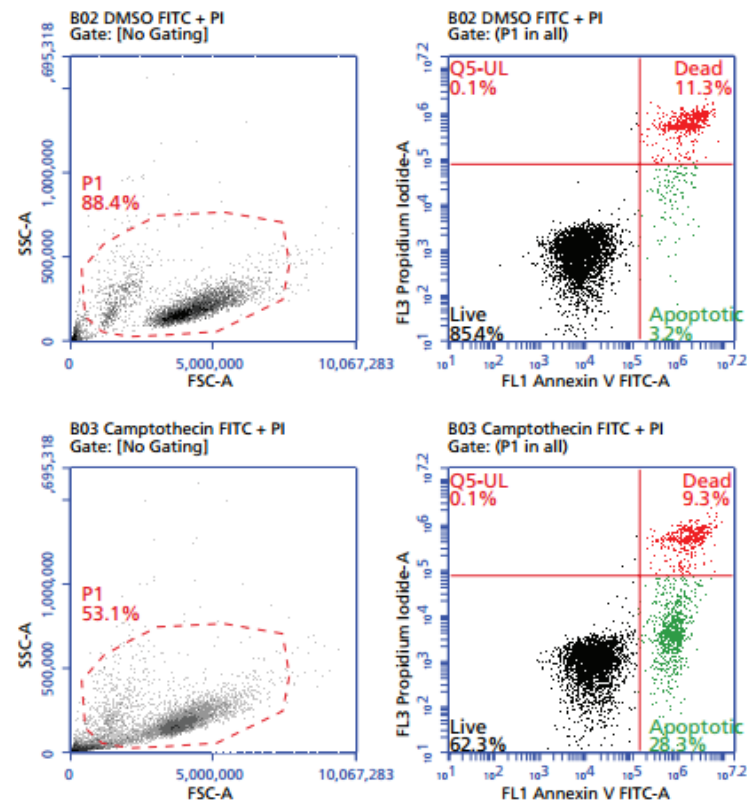
Category	Product Information Sheet	Brand	Kit	Cat. No.	Template
Cell Biology	<a href="#">BD Apoptosis Kits and Templates</a>	BD Pharmingen™	Annexin V FITC Apoptosis Detection Kit II	556570	Download
		BD Pharmingen™	Annexin V PE Apoptosis Detection Kit I	559763	Download
		BD™	MitoScreen (JC-1) Kit	551302	Download
		BD Pharmingen™	Caspase-3 PE Assay Kit	550914	Download
		BD Pharmingen™	Caspase-3 FITC Assay Kit	550480	Download
	<a href="#">BD Cell Cycle and DNA Kits and Templates</a>	BD Cycletest™ Plus	DNA Reagent Kit	340242	Download
		BD Pharmingen™	FITC BrdU Flow Kit	559619	Download
		BD Pharmingen™	APC BrdU Flow Kit	552598	Download

# Kits and Templates on the BD Accuri C6

## Available Templates:

Category	Product Information Sheet	Brand	Kit	Cat. No.	Template
Cell Biology	<a href="#">BD Apoptosis Kits and Templates</a>	BD Pharmingen™	Annexin V FITC Apoptosis Detection Kit II	556570	Download
		BD Pharmingen™	Annexin V PE Apoptosis Detection Kit I	559763	Download
		BD™	MitoScreen (JC-1) Kit	551302	Download
		BD Pharmingen™	Caspase-3 PE Assay Kit	550914	Download
	<a href="#">BD Cell Cycle and DNA Kits and Templates</a>	BD Pharmingen™	Caspase-3 FITC Assay Kit	550480	Download
		BD Cycletest™ Plus	DNA Reagent Kit	340242	Download
		BD Pharmingen™	FITC BrdU Flow Kit	559619	Download
		BD Pharmingen™	APC BrdU Flow Kit	552598	Download

## Sample Data:



# BD Accuri C6 Promotion



The graphic features a dark blue background. On the left, the text reads "BEST IN CLASS meets BEST TIME TO BUY" in white and orange. In the center, two overlapping circles are connected by a white plus sign. The left circle contains a white BD Accuri C6 flow cytometer with a red top and is labeled "10% off" and "BD Accuri™ C6". The right circle contains a box of BD reagents, a large white bottle, and several smaller vials, labeled "40% off" and "BD Reagents".

Note:

US Region Only

Promotion Period: Oct 1, 2014 – Dec 31, 2014



# For Additional Information...

---

If you have further questions:

Contact Technical Support (US) at:

877-232-8995, Prompt 3, 2

or email: [ResearchApplications@bd.com](mailto:ResearchApplications@bd.com)

Please visit our BD Accuri C6 resources site at:

[www.bdbiosciences.com/resources/accuri](http://www.bdbiosciences.com/resources/accuri).





*Dr. Alfonso Blanco Fernández*

Scientific Director of Conway Core Facilities  
Director of Flow Cytometry  
UCD - Conway Institute  
University College Dublin  
Belfield, Dublin 4. IRELAND

e-mail: [alfonso.blanco@ucd.ie](mailto:alfonso.blanco@ucd.ie)

Tel: 00353(0)1 716 6966/6947

# *REAL TIME* Flow Cytometry



International Society for Advancement of Cytometry

ISAC Scholars Program



# CORE

## TECHNOLOGIES AT UCD CONWAY INSTITUTE



# Flow Cytometric Measurements of Cytoplasmic Calcium Changes in Human Platelets<sup>1</sup>

Theresa A. Davies, Daniel Drotts, Gary J. Weil, and Elizabeth R. Simons<sup>2</sup>

Department of Biochemistry, Boston University School of Medicine, Boston, Massachusetts 02118

Received for publication June 26, 1987; accepted September 30, 1987

**Measurements of cytosolic  $Ca^{++}$  in individual cells.** Platelets loaded with Indo-1 were diluted to  $10 \times 10^6$  per ml in HEPES buffer, pH 7.4. Fluorescence measurements were obtained by using a FACS 440 dual laser system (Becton Dickinson) equipped with a Consort 40/PDP-11/23 microcomputer (Digital Corp.). A HEPES

of probe. Thrombin (0.001–0.05 U/ml) was then added directly to the sample and data were collected at the rate of 4,000 cells per time point after stimulation, at approximately 5-s intervals, with the first time point at 10 s. All data were collected in the linear mode to allow

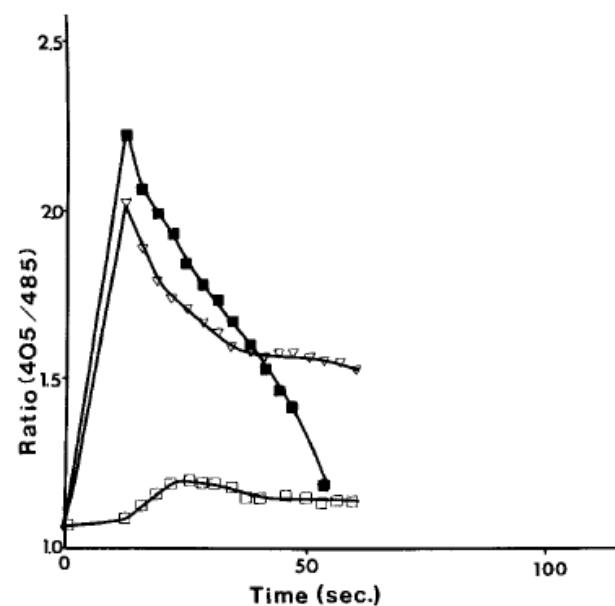


FIG. 2. Time course of the thrombin-induced change in  $[Ca^{++}]_{in}$  in responding and nonresponding populations of human platelets. The 405-nm/485-nm ratio of Indo-1 loaded platelets after thrombin stimulation by a 0.005 U/ml (subsaturating) thrombin dose vs. time. The responding population (■) and the nonresponding population (□) were determined by gating above baseline on the FACS. Also plotted is the total platelet response, which illustrates an average of responding and nonresponding populations (▽).



# Stop-flow method



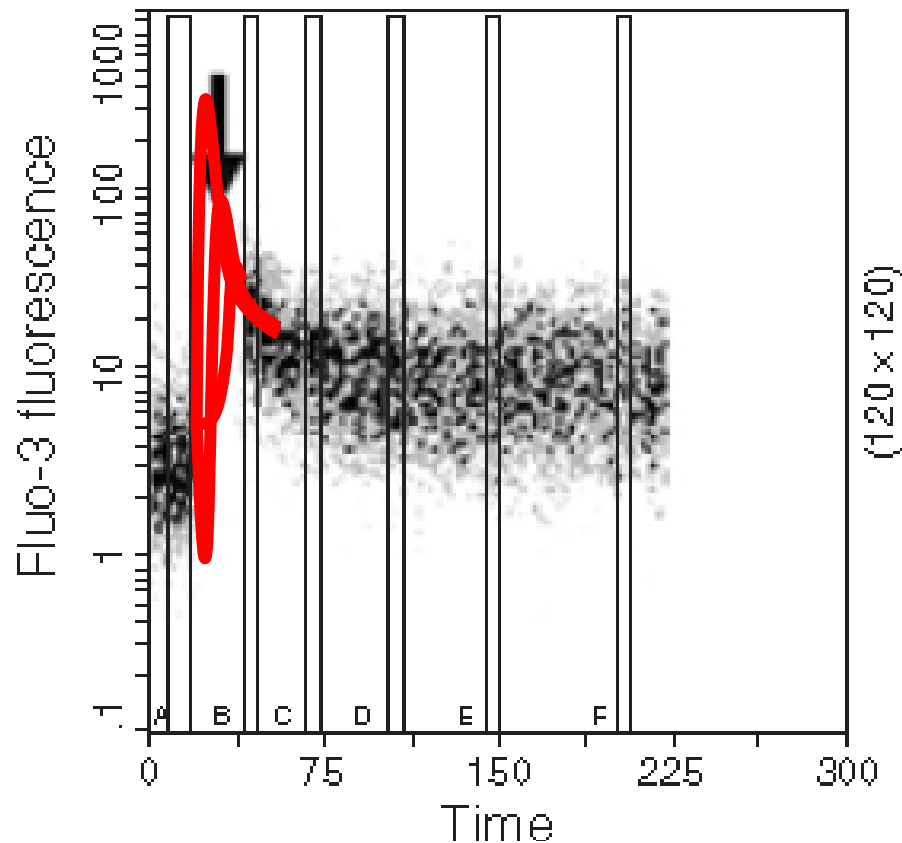
## Flow Cytometric Analysis of Calcium Mobilization in Whole-Blood Platelets

Contributed by Maria-do-Céu Monteiro, Maria-José Gonçalves, Filipe Sansonetty, and José-Enrique O'Connor

*Current Protocols in Cytometry* (2003) 9.20.1-9.20.8

Copyright © 2003 by John Wiley & Sons, Inc.

D



Panel D shows the procedure and time course of a typical kinetic assay of intracellular  $\text{Ca}^{2+}$  mobilization

# Continuous measurement



2009



TECHNICAL NOTE

## Cytometry

PART A  
Journal of the  
International Society for  
Advancement of Cytometry

### A Flow-Cytometric Method for Continuous Measurement of Intracellular $\text{Ca}^{2+}$ Concentration

Alice Vines,<sup>1</sup> Gethin J. McBean,<sup>1</sup> Alfonso Blanco-Fernández<sup>2\*</sup>

<sup>1</sup>UCD School of Biomolecular and Biomedical Science, UCD-Conway Institute, University College Dublin, Belfield, Dublin 4, Ireland

<sup>2</sup>Flow Cytometry Core Facilities, UCD-Conway Institute of Biomolecular and Biomedical Research, University College Dublin, Belfield, Dublin 4, Ireland

Received 26 November 2009; Revision Received 21 July 2010; Accepted 25 August 2010

Grant sponsor: Health Research Board of Ireland.

\*Correspondence to: Alfonso Blanco-Fernández, Flow Cytometry Core Facilities, UCD-Conway Institute of Biomolecular and Biomedical Research, University College Dublin, Belfield, Dublin 4, Ireland

Email: alfonso.blanco@ucd.ie

Published online 24 September 2010 in Wiley Online Library

#### • Abstract

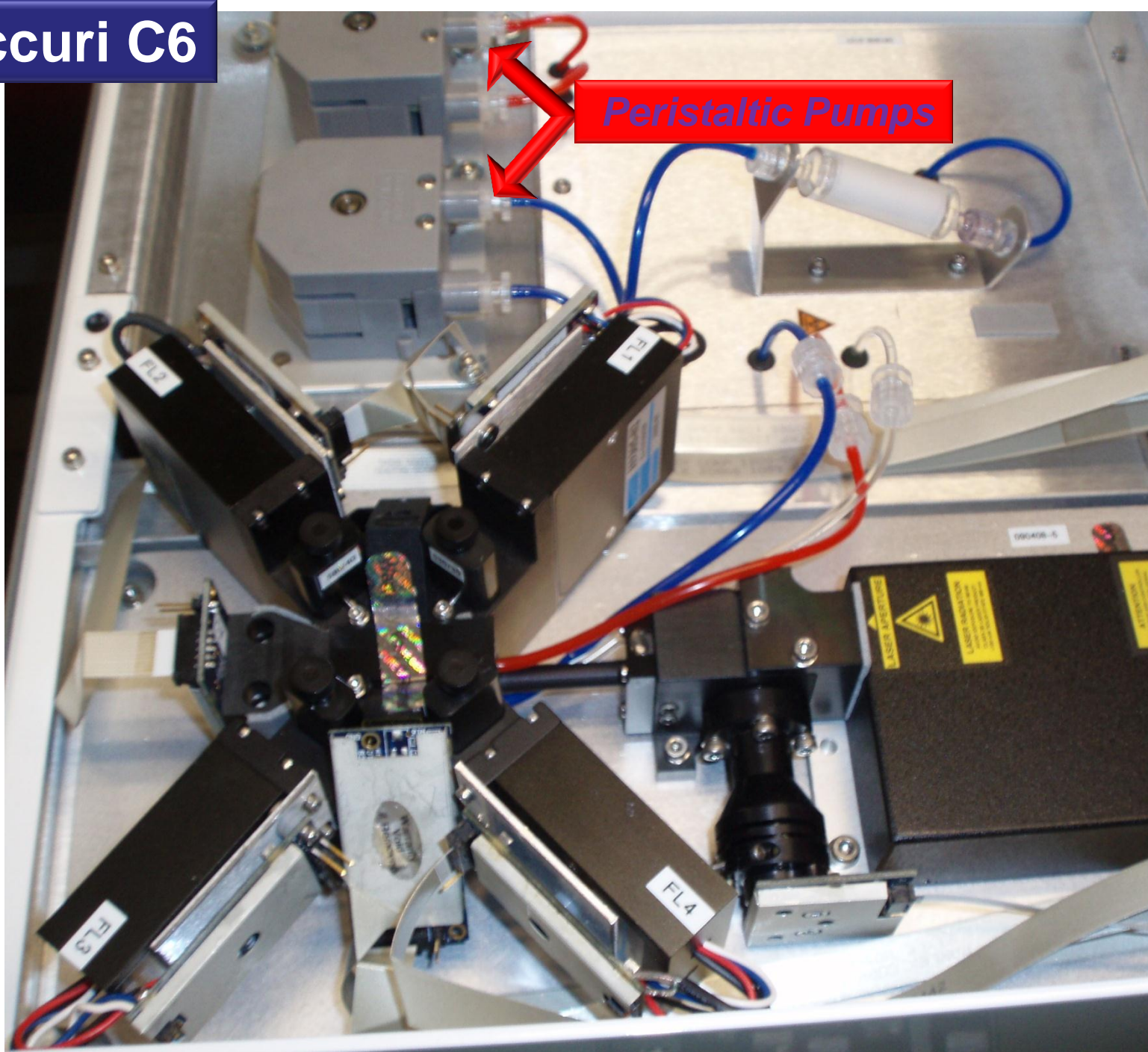
Alterations in intracellular  $\text{Ca}^{2+}$  concentration are amongst the most rapid responses to a variety of stimuli in mammalian cells. In the nervous system in particular, responses occur within nanoseconds. A major challenge in intracellular  $\text{Ca}^{2+}$  analysis is to achieve measurements within this very fast time frame. To date, the dynamic intracellular  $\text{Ca}^{2+}$  concentration has been monitored by confocal microscopy, plate-based assays, spectrofluorometry, and flow cytometry, although there are issues with the number of cells analyzed or gaps in recording due to the addition of compounds, with significant loss of detail of a rapid  $\text{Ca}^{2+}$  response. The new generation of flow cytometers (such as Accuri C6) resolves this problem by allowing the addition of test compounds with continuous monitoring of thousands of cells, providing a method for dynamic  $\text{Ca}^{2+}$  measurements. This system was tested with commonly used  $\text{Ca}^{2+}$  modulating agents in C6 glioma cells. Thapsigargin (TG), a blocker of  $\text{Ca}^{2+}$  uptake into the endoplasmic reticulum (ER), causes a significant increase in the intracellular calcium concentration via ER emptying followed by  $\text{Ca}^{2+}$  entry via store-operated  $\text{Ca}^{2+}$  channels (SOCC). This well-established pathway can be partially inhibited by 2-aminoethoxydiphenyl borate (2-APB), a blocker of SOCC. Both the increase with TG alone and the partial increase when coincubated with 2-APB were observed with continuous recording along with calibration curves using an Accuri C6 flow cytometer. With these new cytometers, dynamic  $\text{Ca}^{2+}$  concentration measurement becomes extremely accessible and accurate, while also providing extensive and valuable data regarding population health and responsiveness. © 2010 International Society for Advancement of Cytometry

#### • Key terms

continuous calcium measurement; calcium flux;  $\text{Ca}^{2+}$ ; live cells; kinetic flow cytometry



# BD Accuri C6

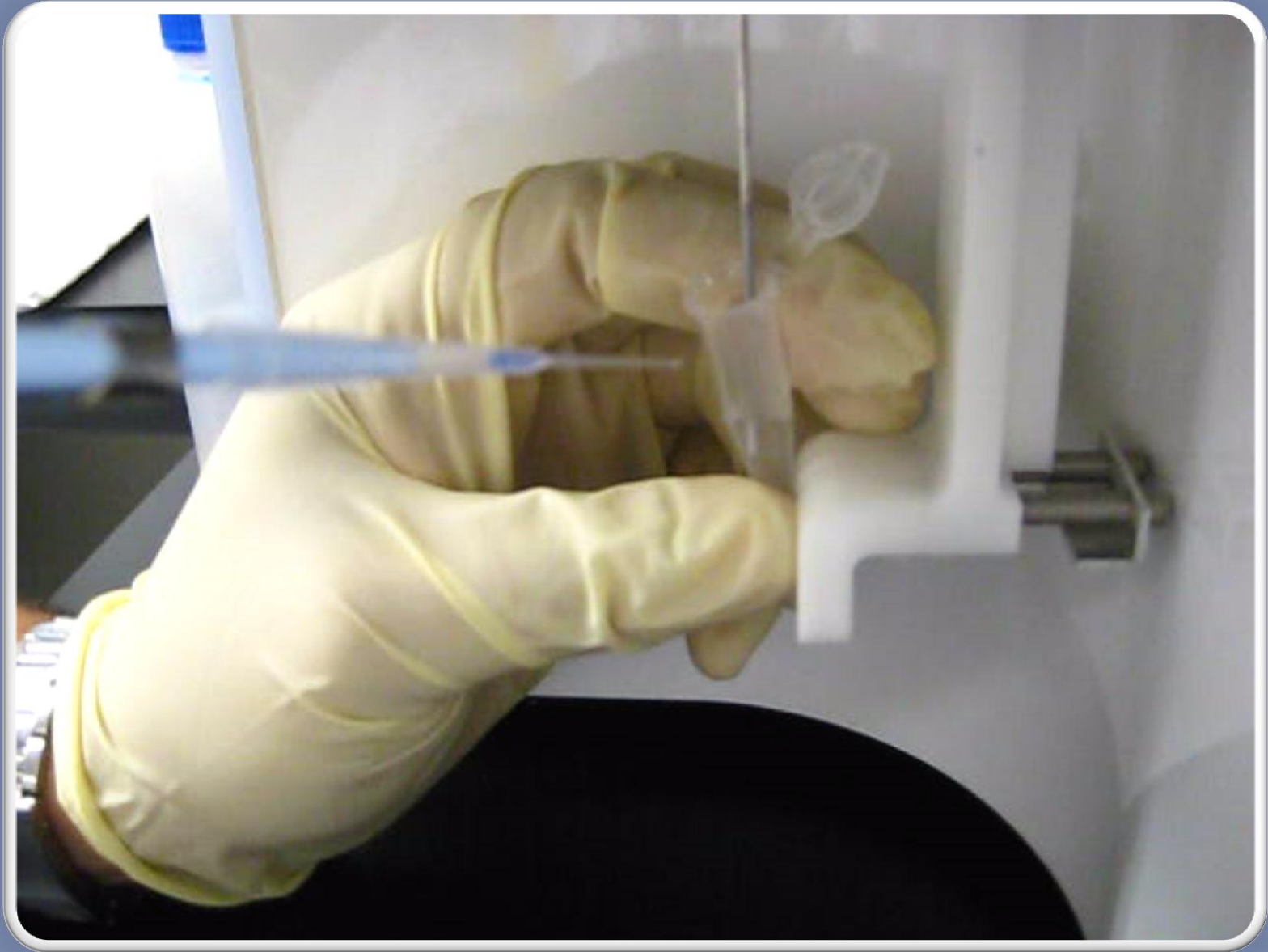


*Peristaltic Pumps*

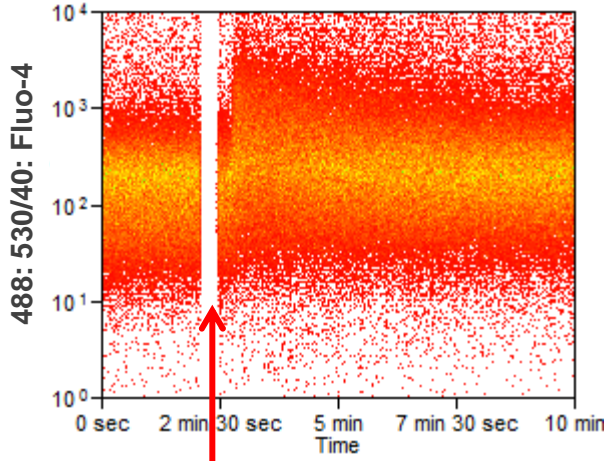
# BD Accuri C6



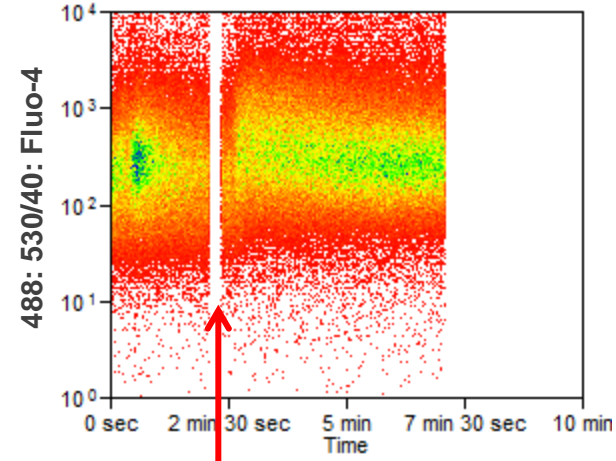
# BD Accuri C6



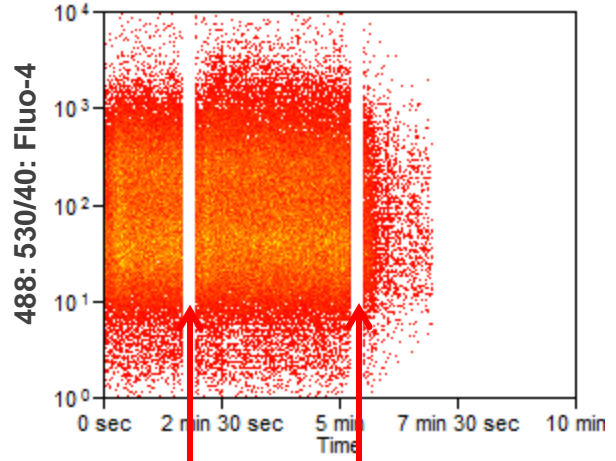
**Beckman Coulter Cyan ADP:**



**A23187**

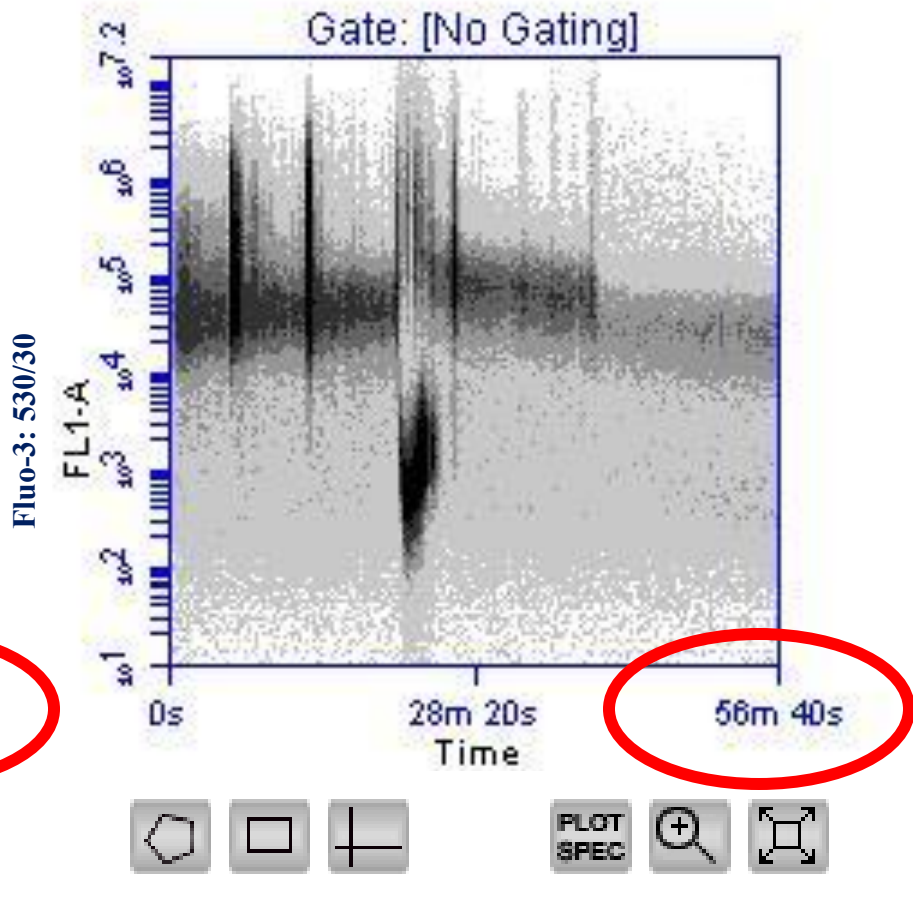
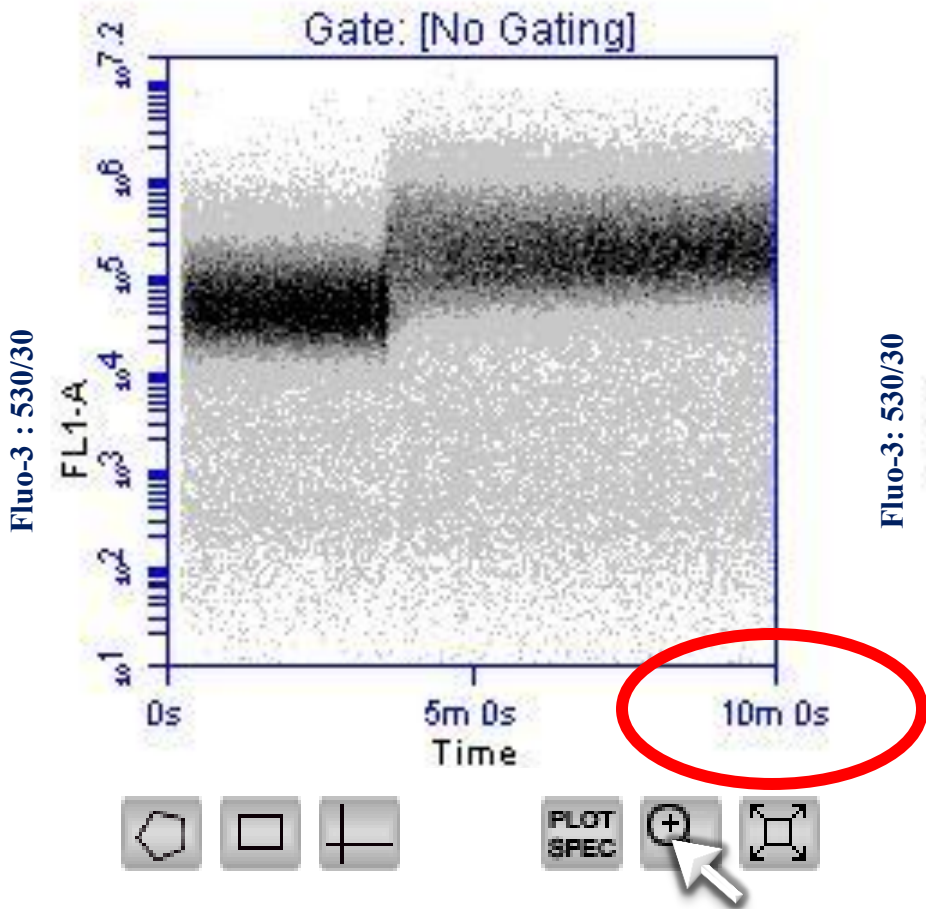


**A23187**



**Thapsigargin A23187**

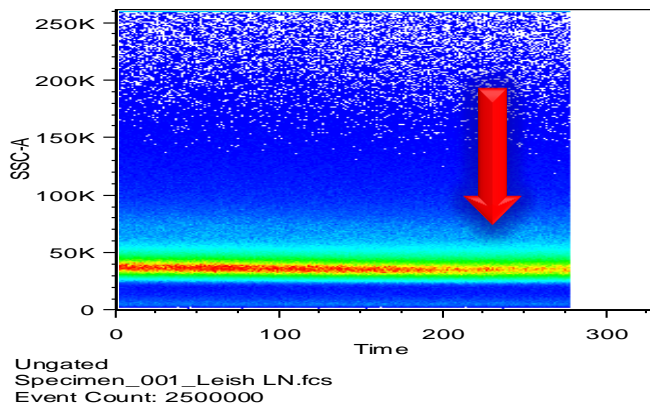
# BD Accuri C6. Time limitations?



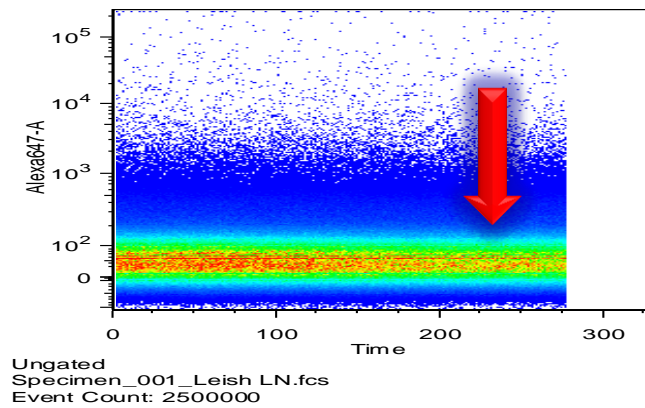


**Attention!!**

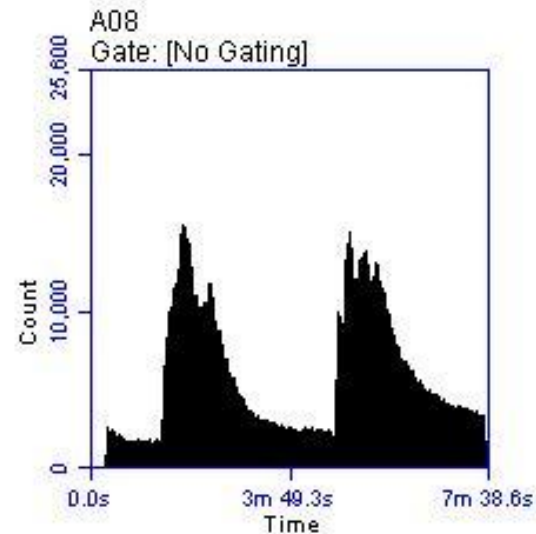
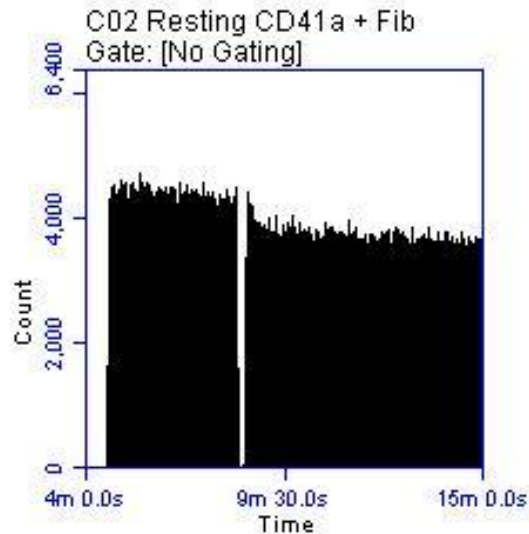
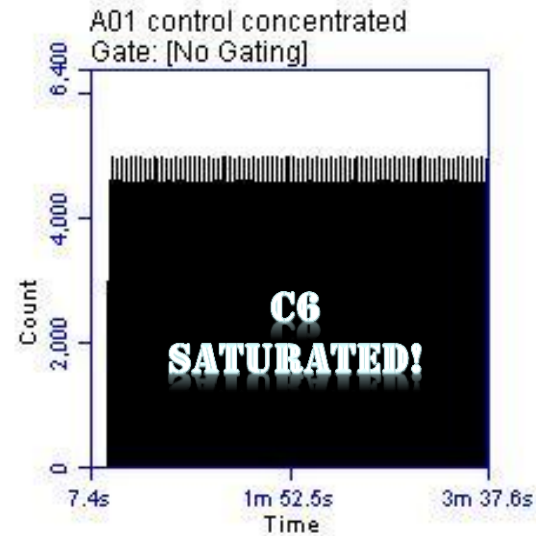
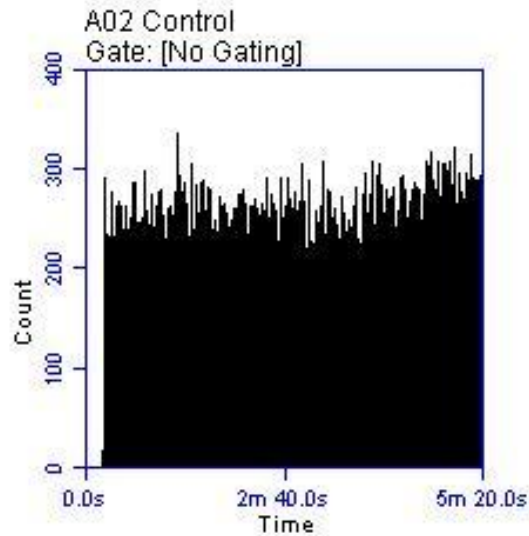
*Scatter*



*Fluorescence*



# Attention!!

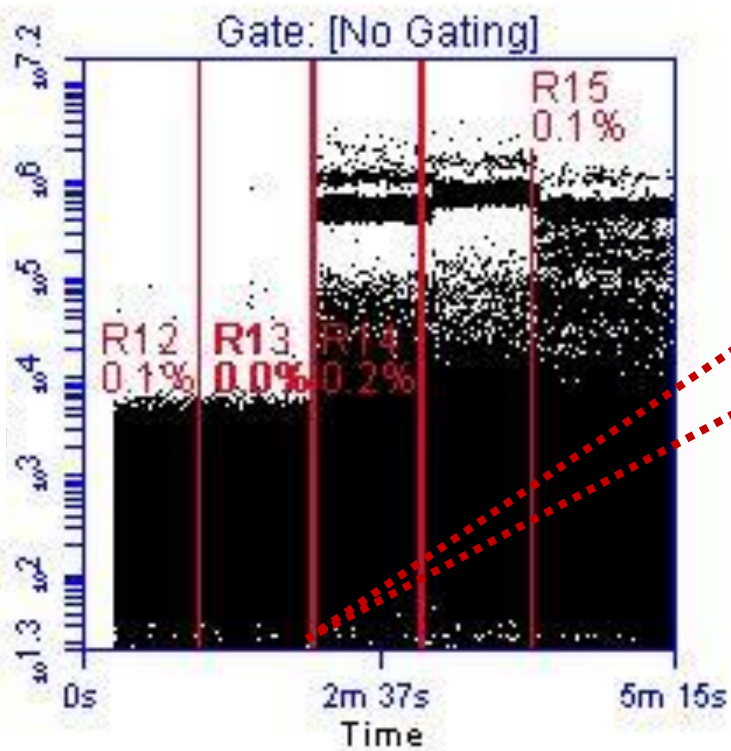


**Remember! Maximum Number of Events/Well =  $10^6$**   
**Optimise Number of Events/ sec**  
**based on the time of your experiment**

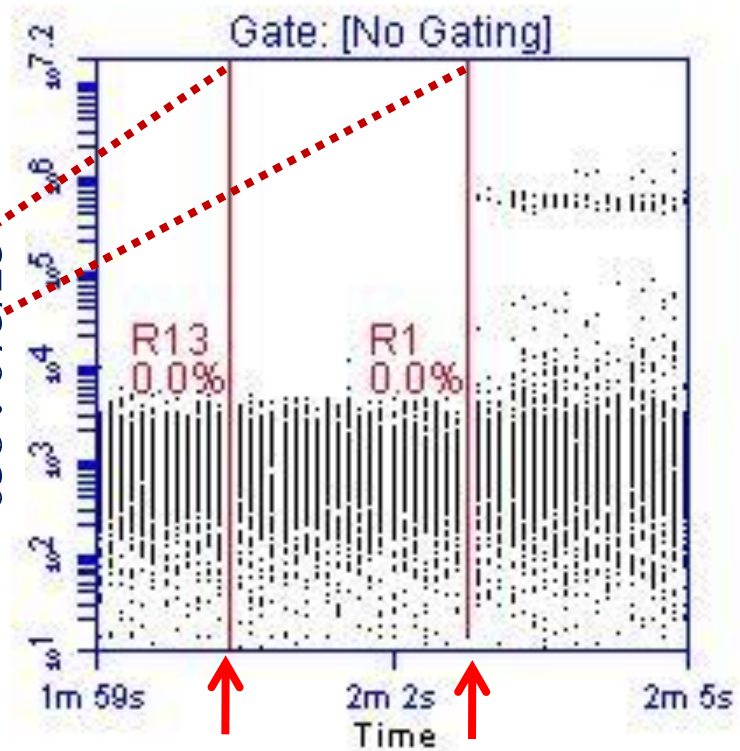


# How much time between addition and analysis?

638 : 675/25



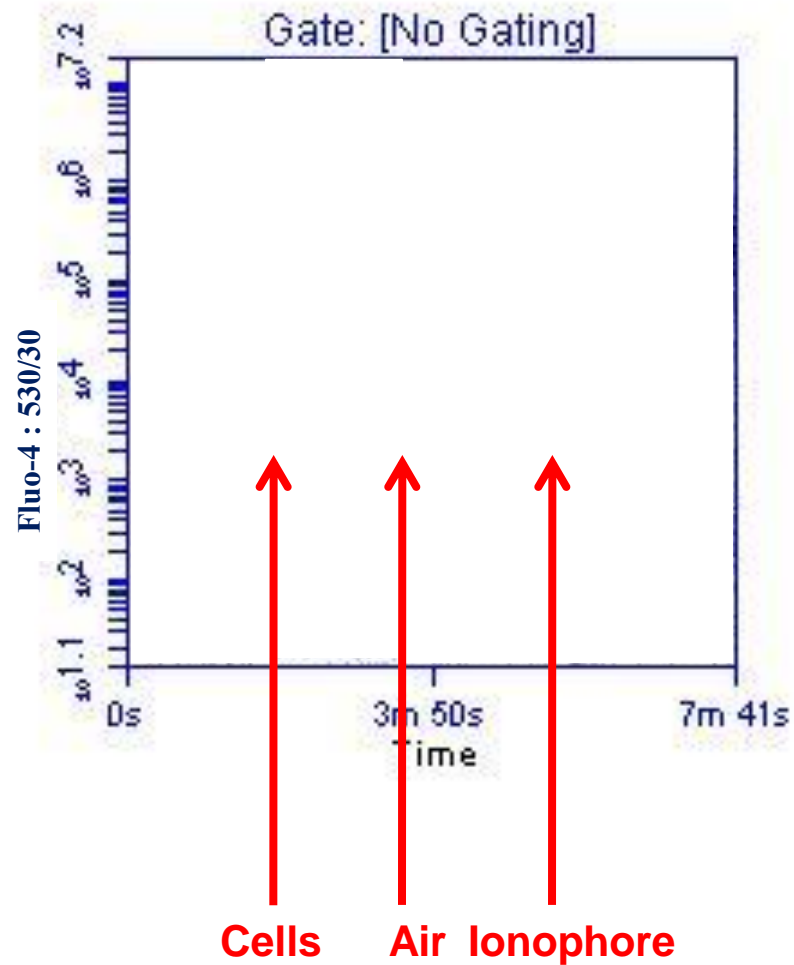
638 : 675/25



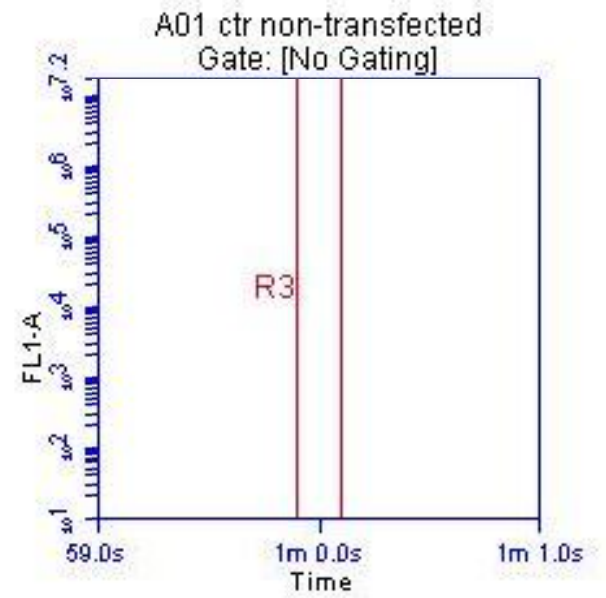
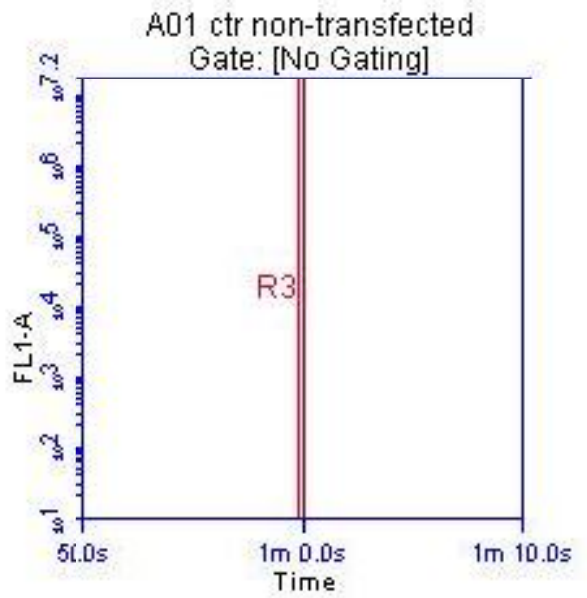
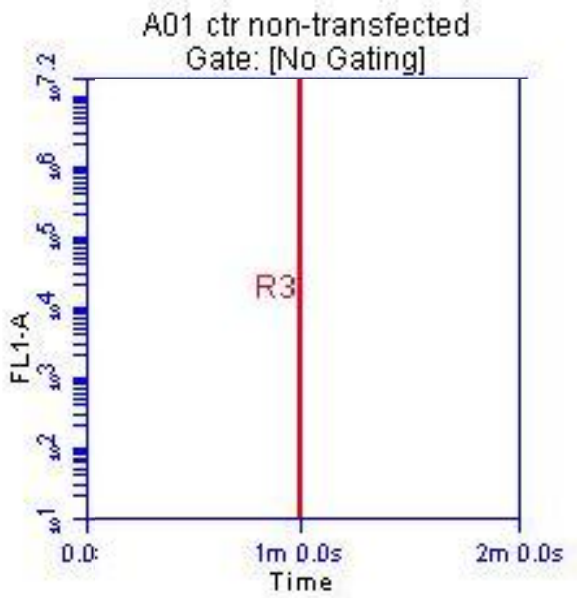
**Addition Time**      **Analysis Time**



# Do air bubbles affect the flow rate?



# Can you control the addition time?





Apoptosis

Calcium/  
other ions

GFP

Applications

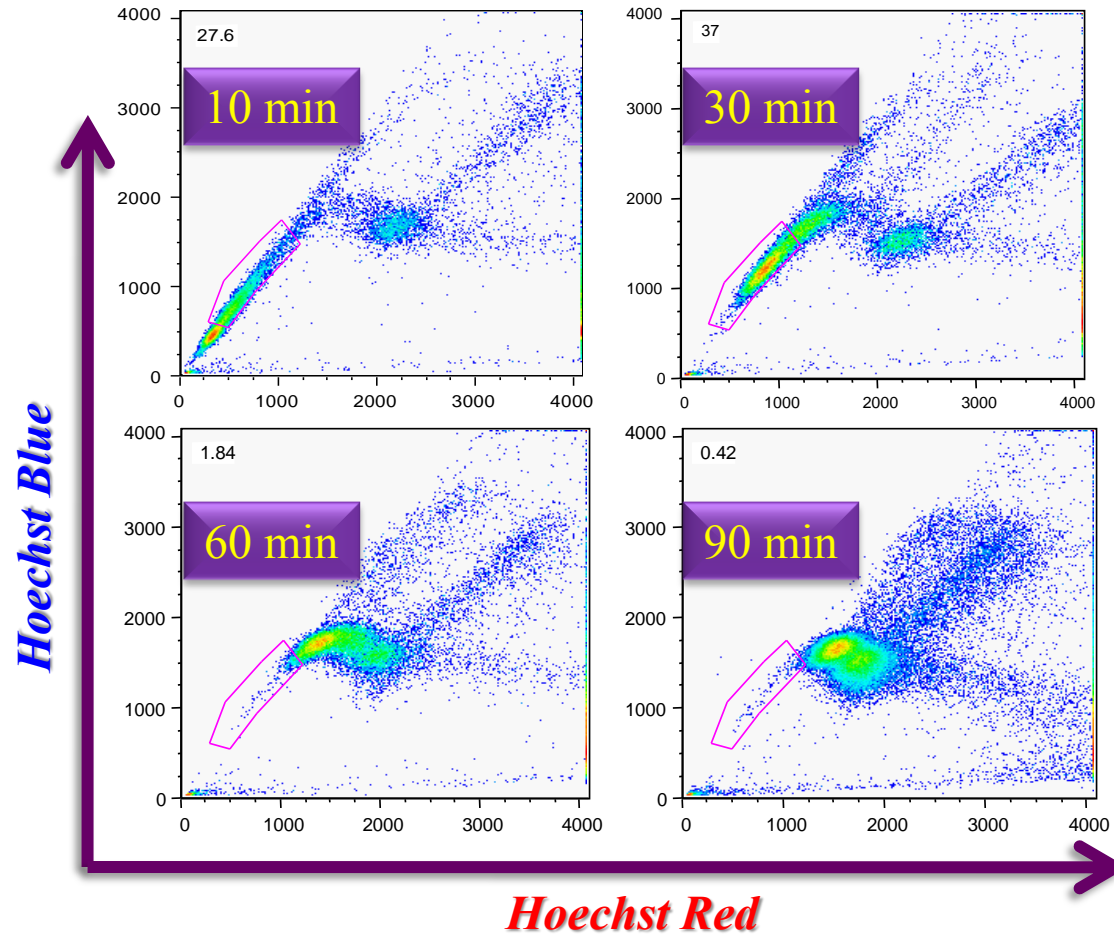
Staining  
optimization

Platelet  
activation

Phagocytosis/  
Nanoparticle uptake



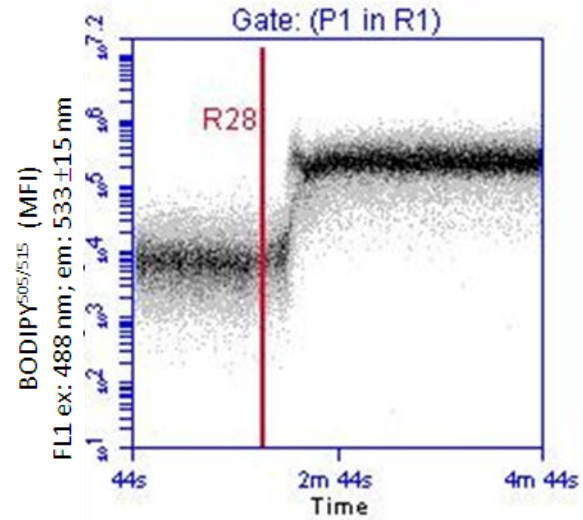
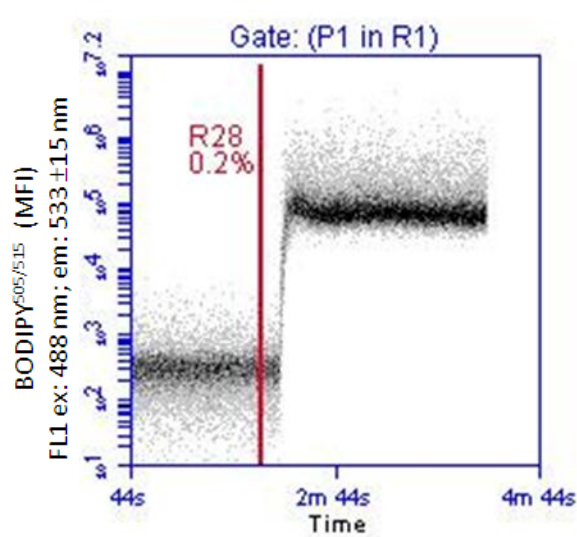
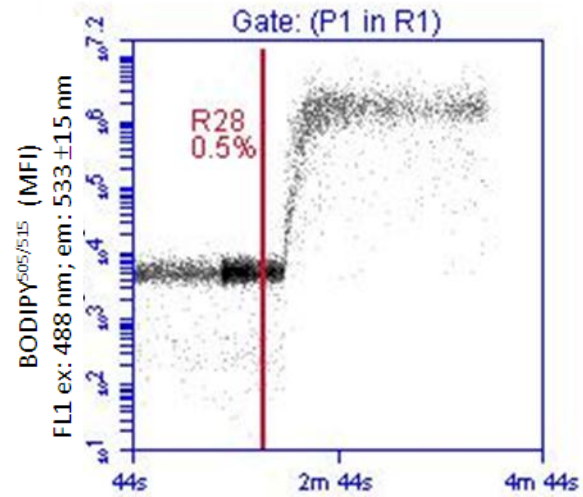
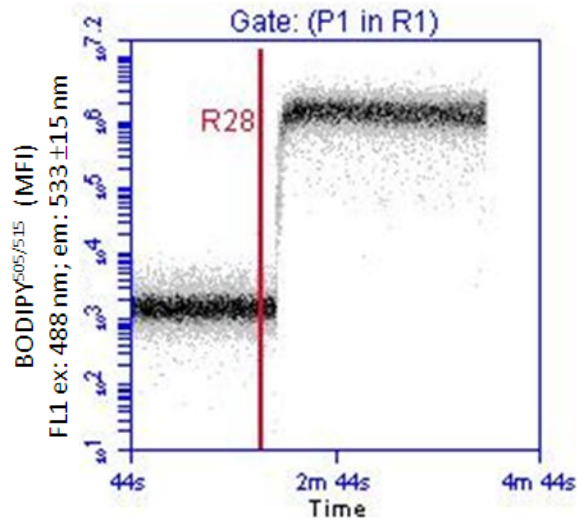
## Murine Bone Marrow - 5 $\mu$ g/ml



# Analysis of Algae physiology



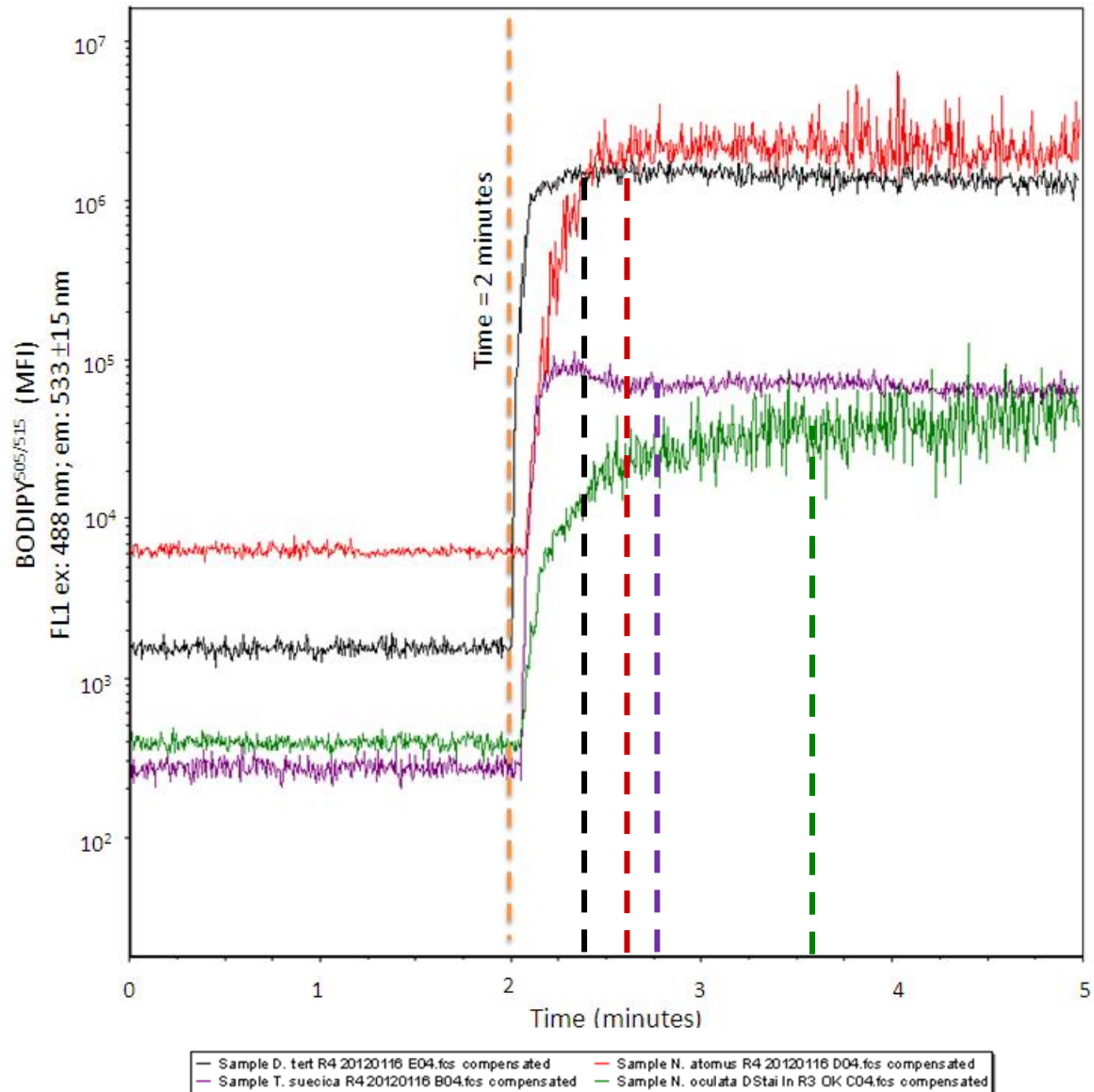
## Kinetics of 4 different algae



# Analysis of Algae physiology

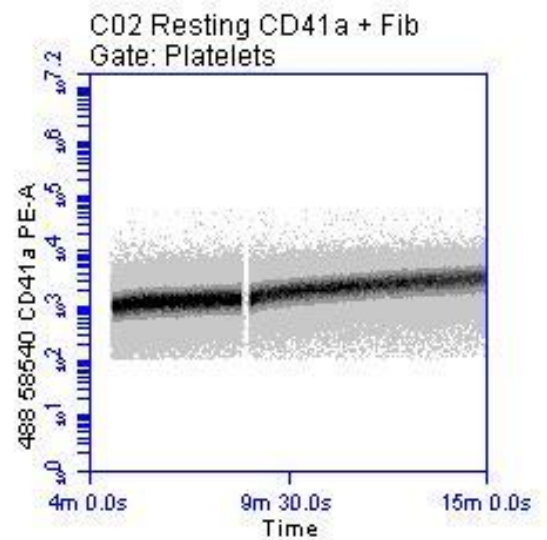
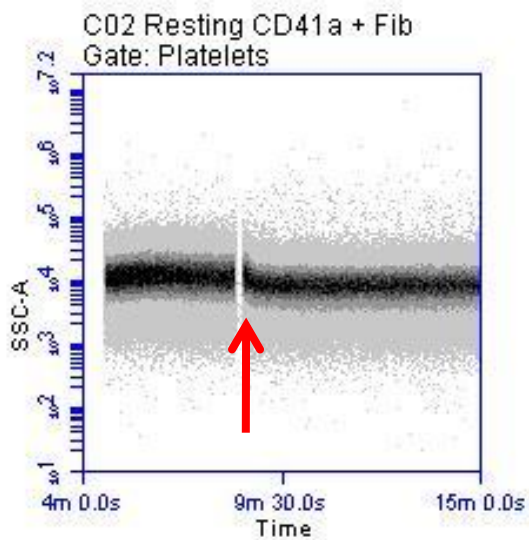
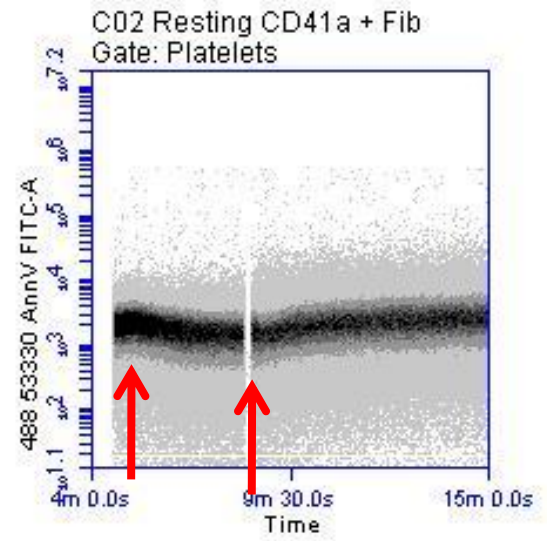
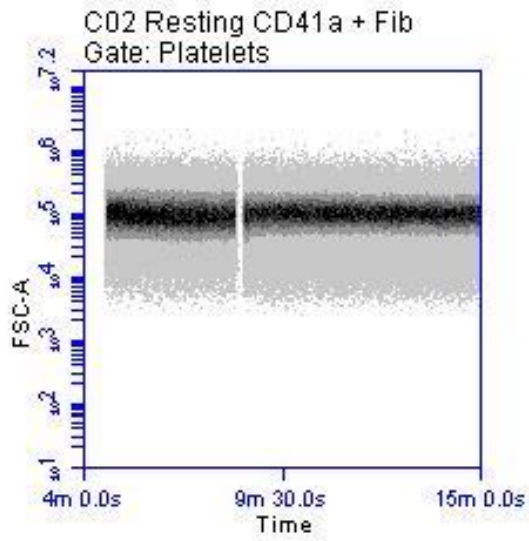
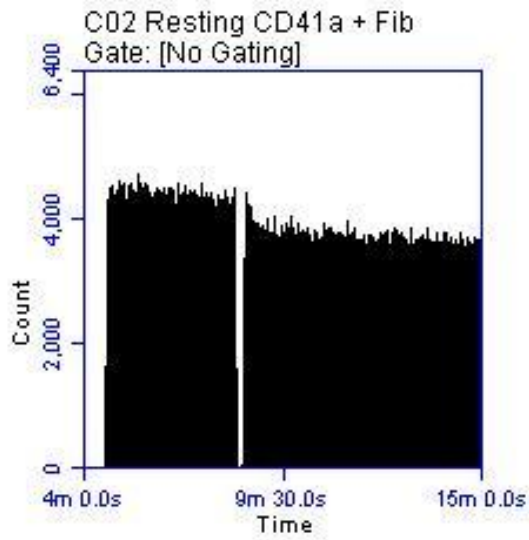


## Kinetics of 4 different algae (FCS Express v.4)

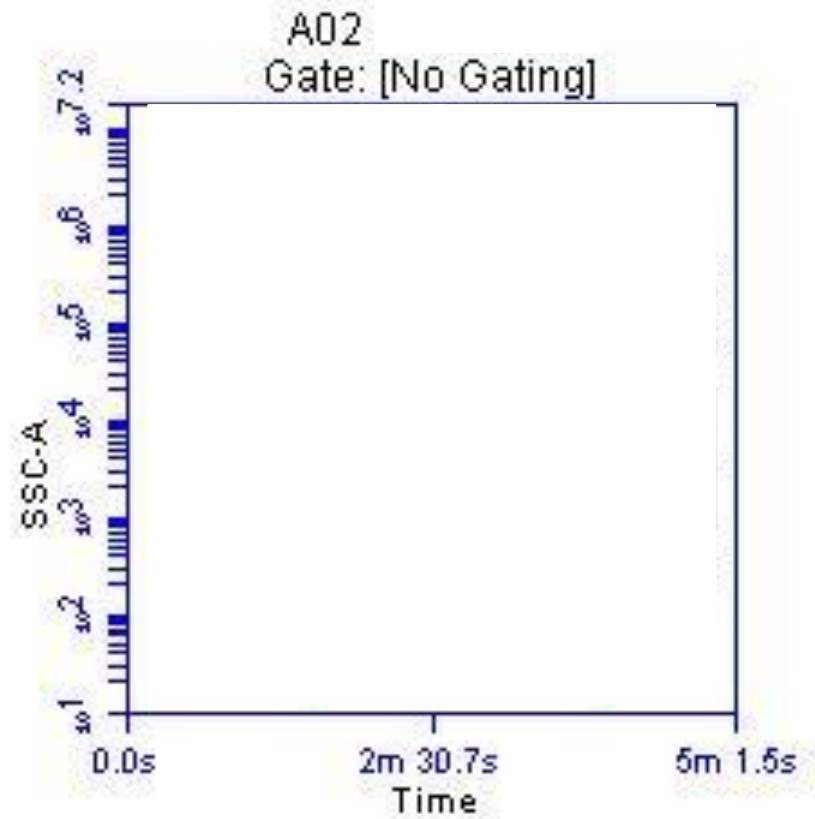
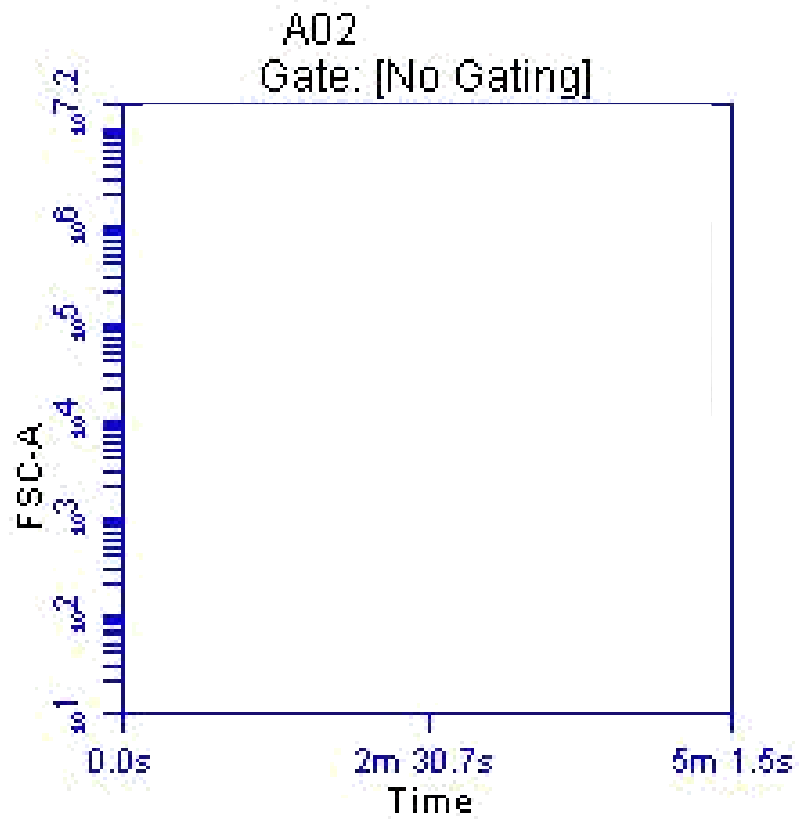




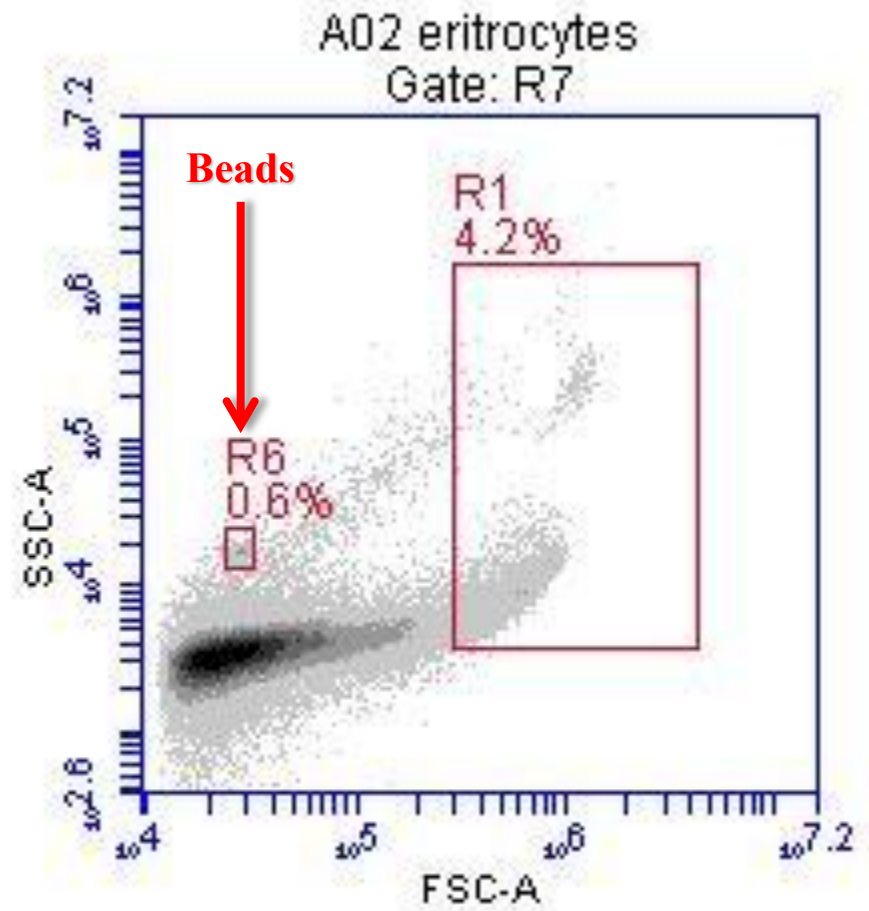
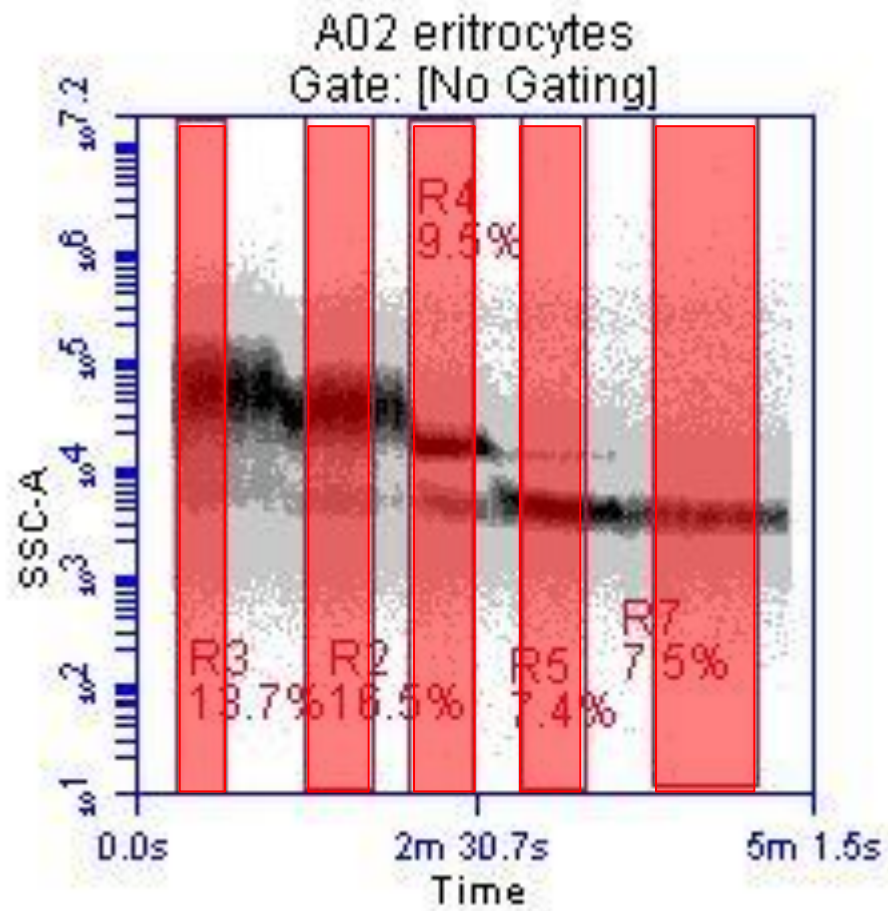
# Platelets



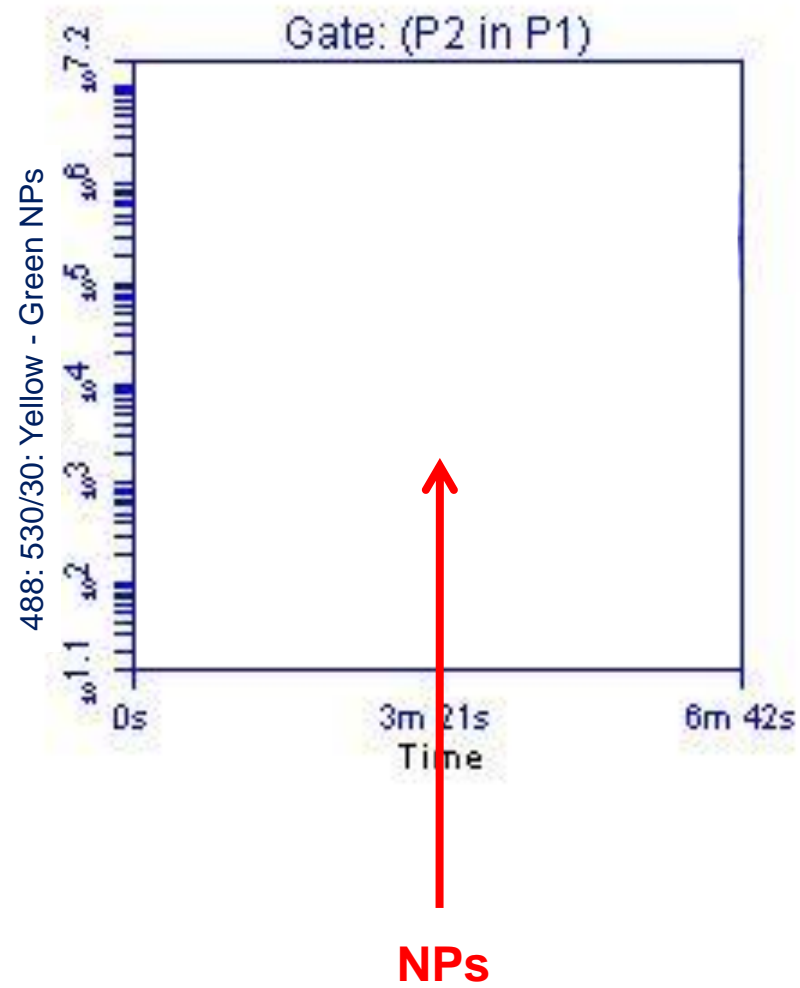
# Erythrocytes Lysing



# Erythrocytes Lysing



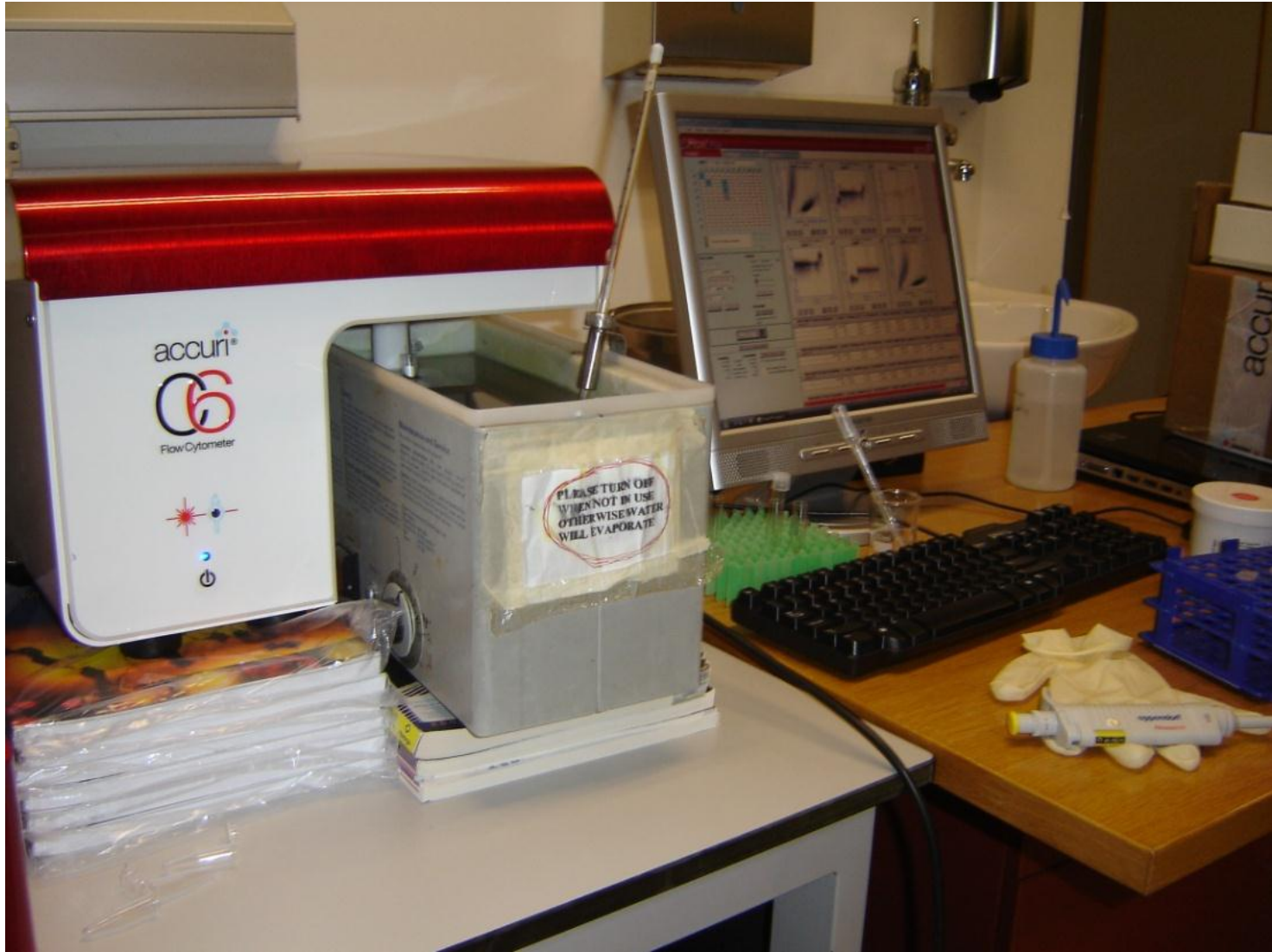
# Nanoparticles uptake



***Nanoparticles uptake***

A photograph of a baby sitting in a blue plastic tub filled with water. The baby is wearing a white headband and has their arms resting on the edges of the tub. The background shows a tiled floor and a white wall.

***Temperature control?***



***What if my sample has to be at 4°C?***







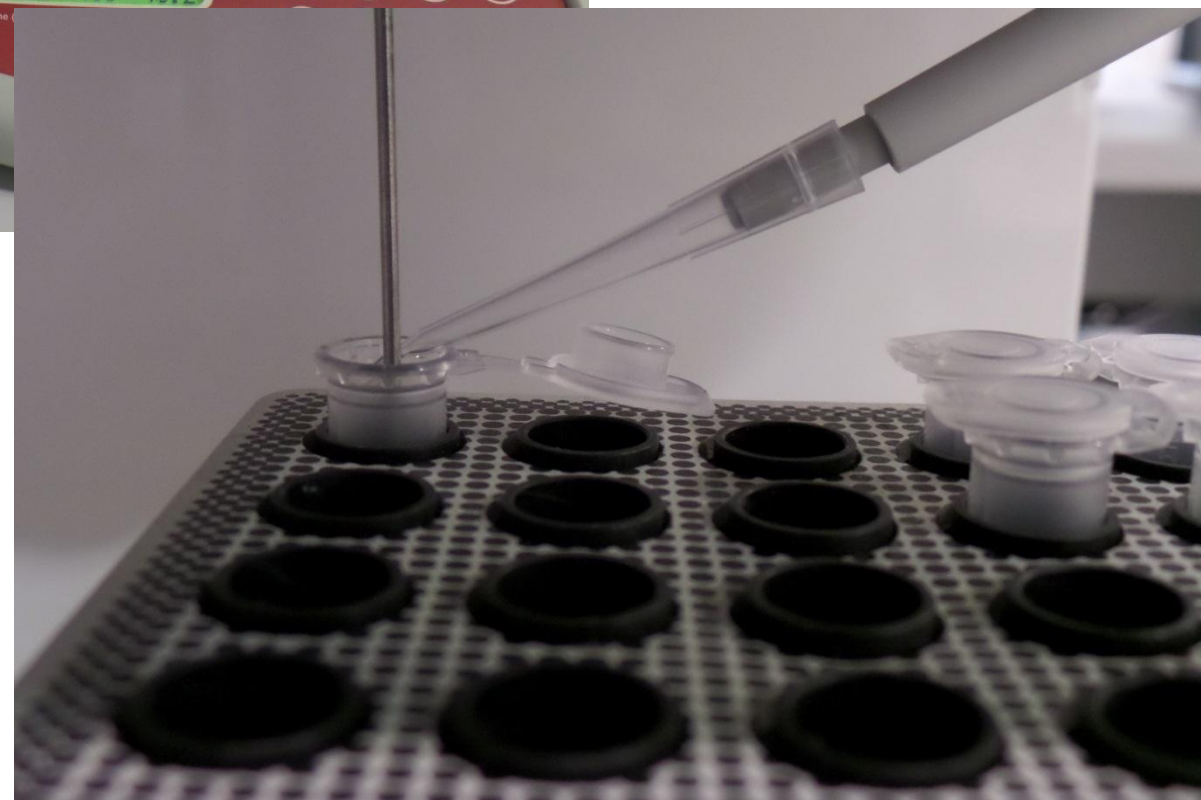
BD Accuri™



SB! biometra  
a part of Analytik Jena

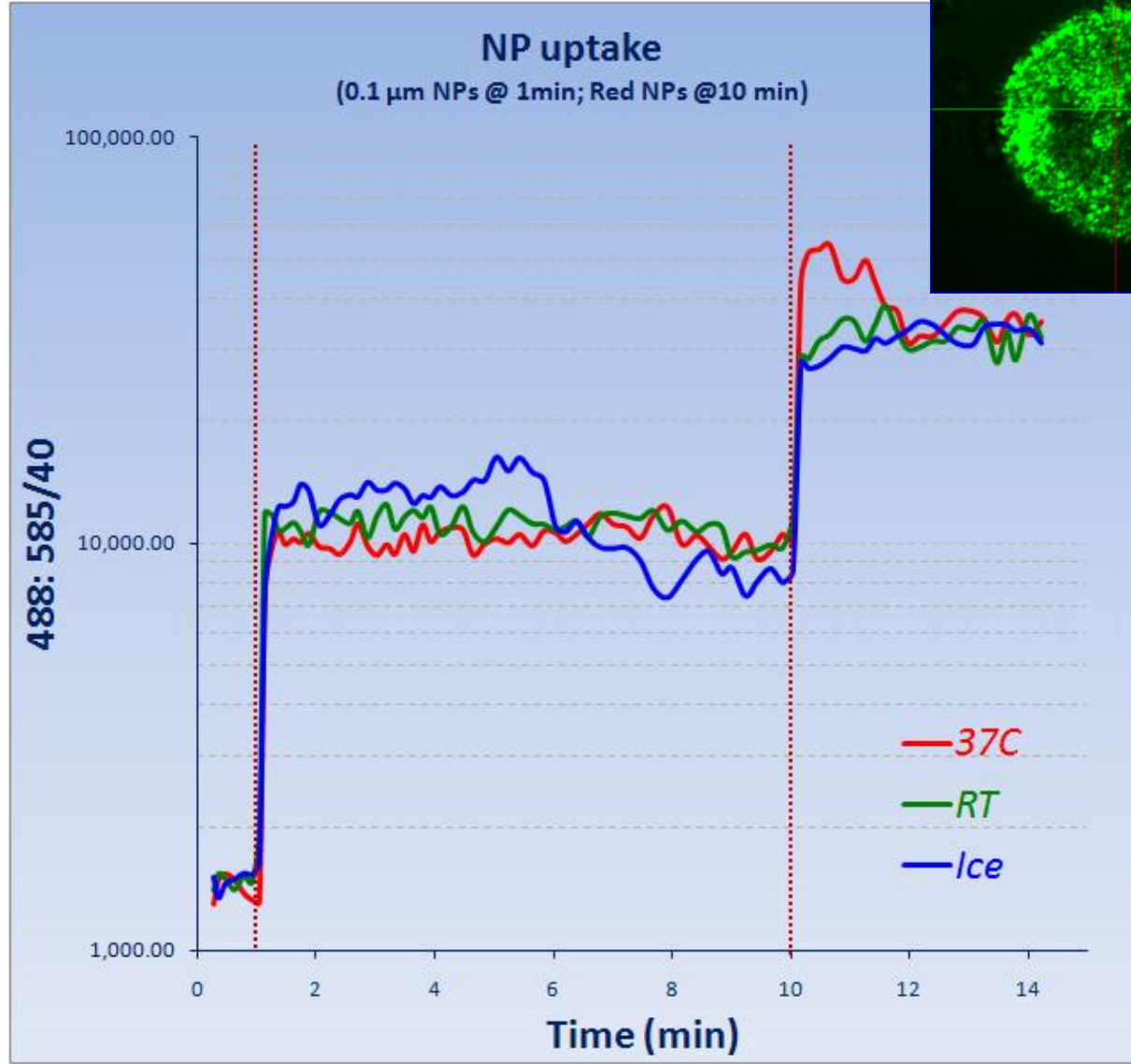
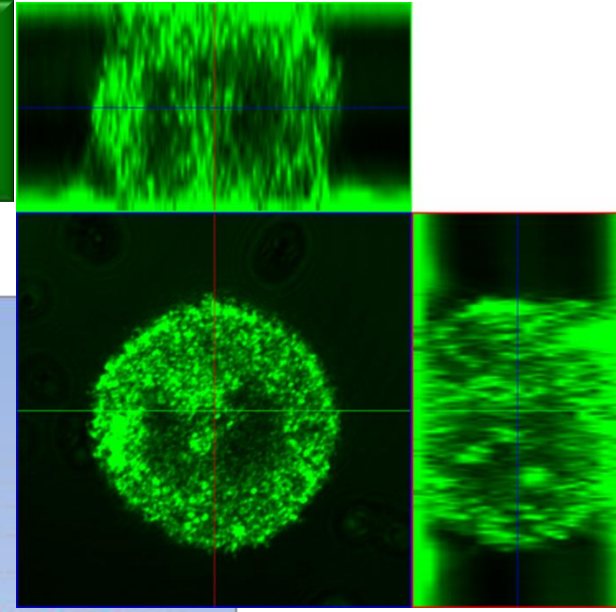
Tsi ThermoShaker

00:02	420	40.0	Set
00:00	418	40.2	Actual



# Nanoparticles uptake

Fixed cell in suspension covered by nanoparticles, after incubation at 4°C



## Measurement of Intracellular Calcium Ions by Flow Cytometry

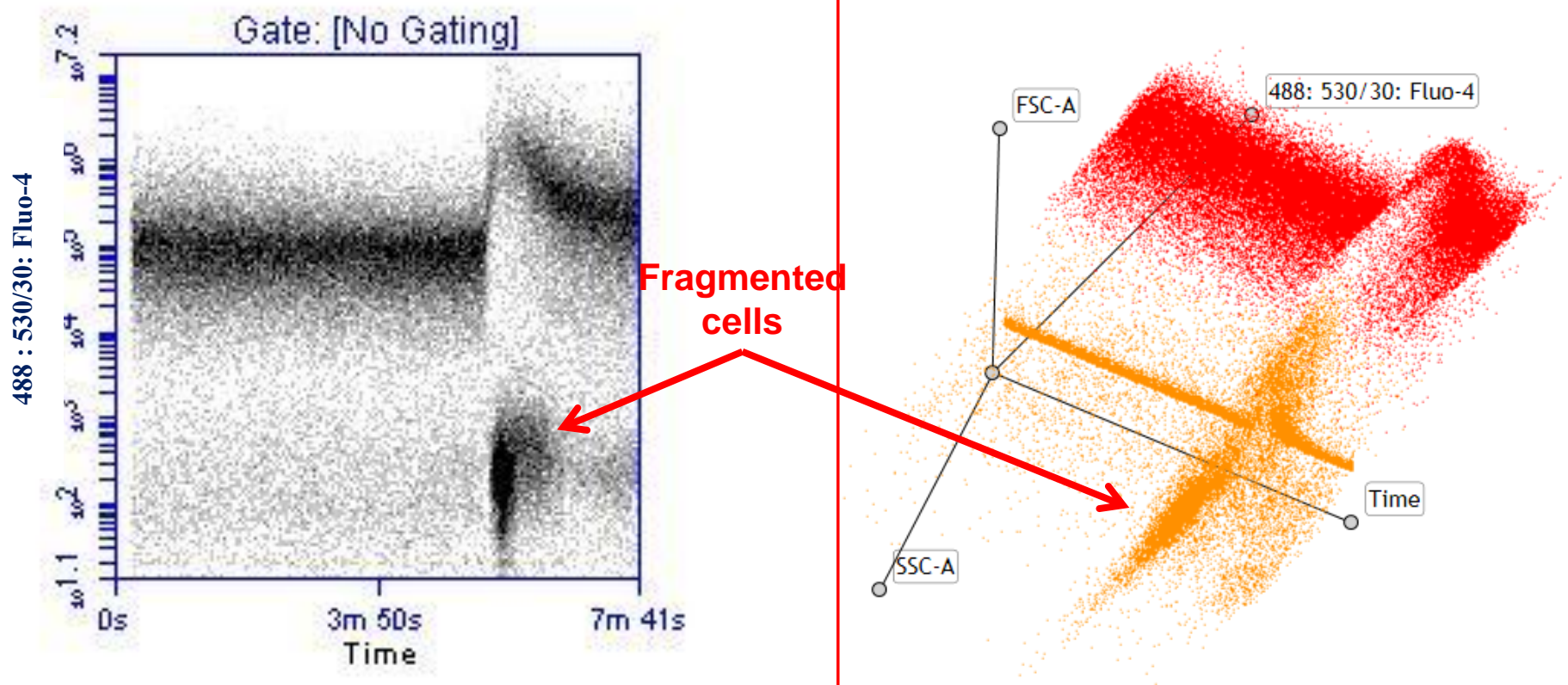
Contributed by Carl H. June, Ryo Abe, and Peter S. Rabinovitch

*Current Protocols in Cytometry* (1997) 9.8.1–9.8.19

Copyright © 1997 by John Wiley & Sons, Inc.

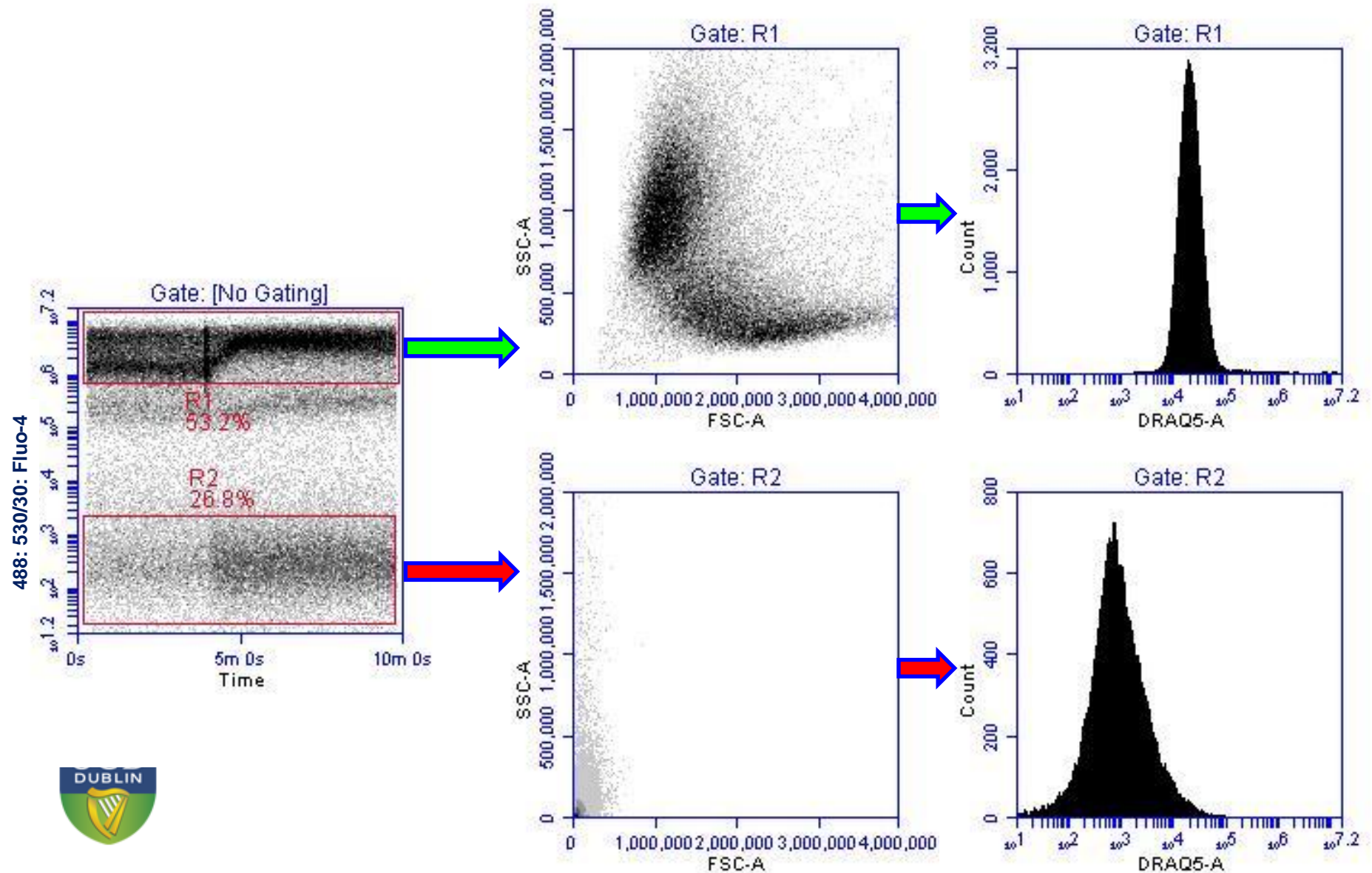
### Analyze fluo-3-loaded cells

6. Analyze cells by gating on forward and right-angle light scatter. Exclude dead cells by gating out cells without fluo-3 fluorescence.

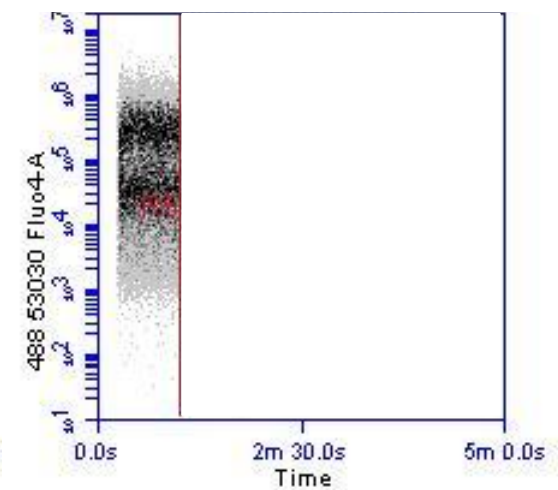
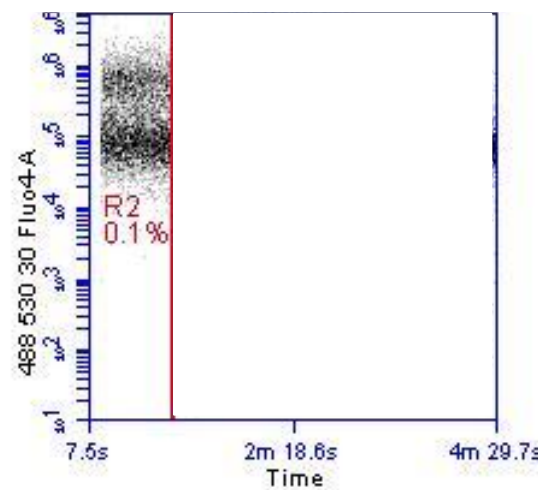
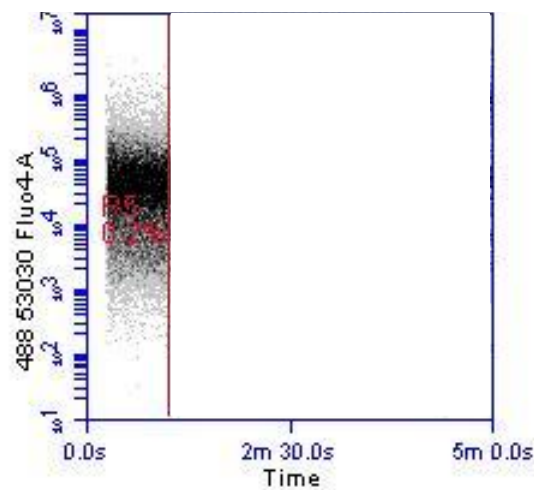
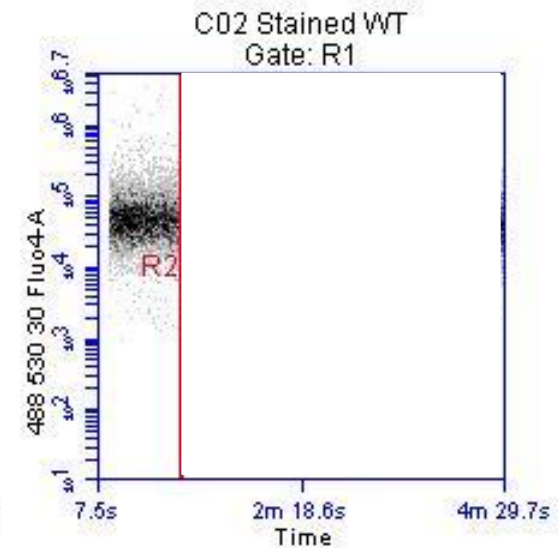
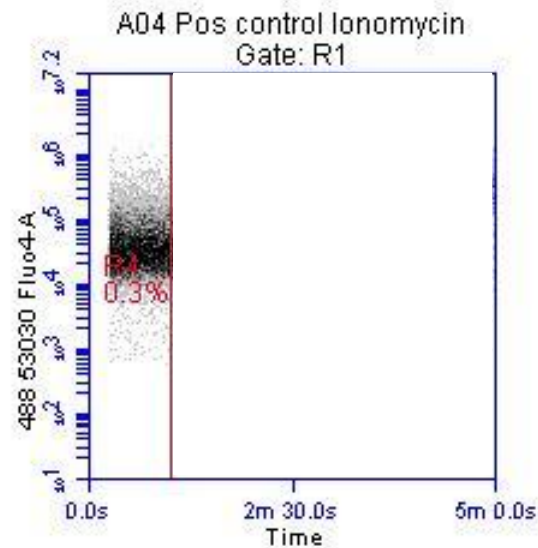
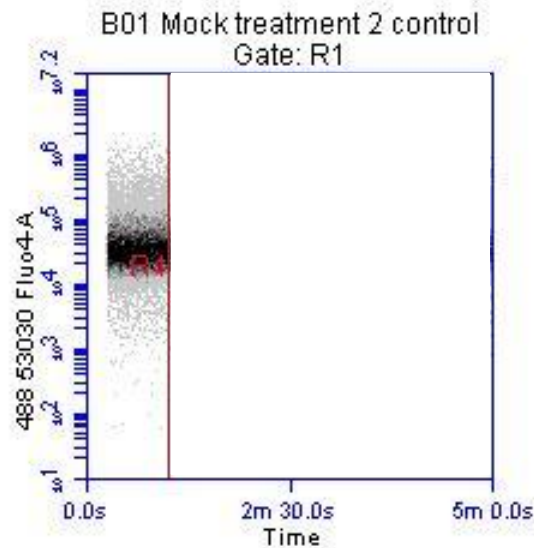


## Analyze fluo-3-loaded cells

- Analyze cells by gating on forward and right-angle light scatter. Exclude dead cells by gating out cells without fluo-3 fluorescence.

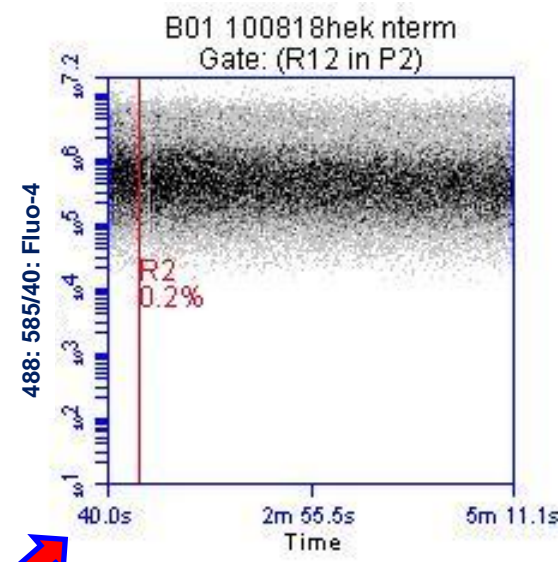
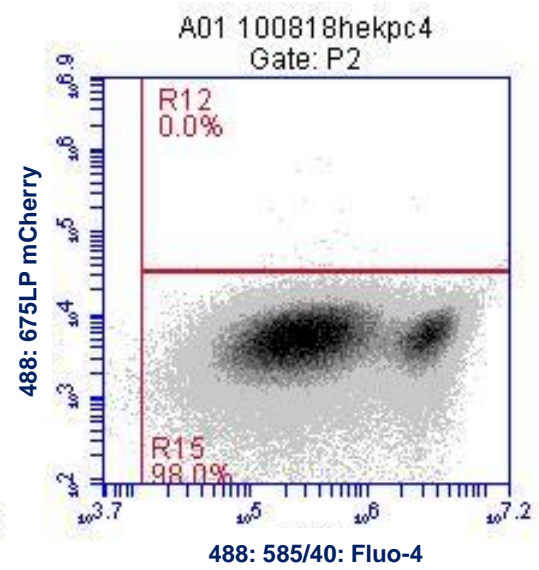
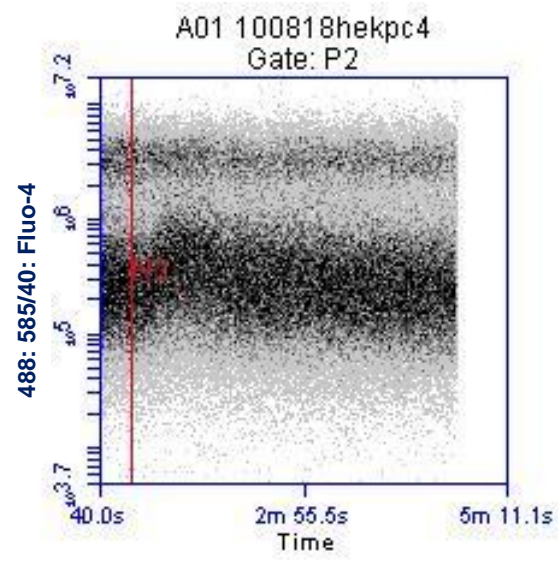


# Some Calcium profiles

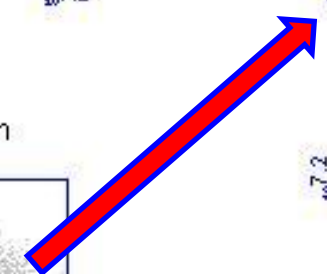
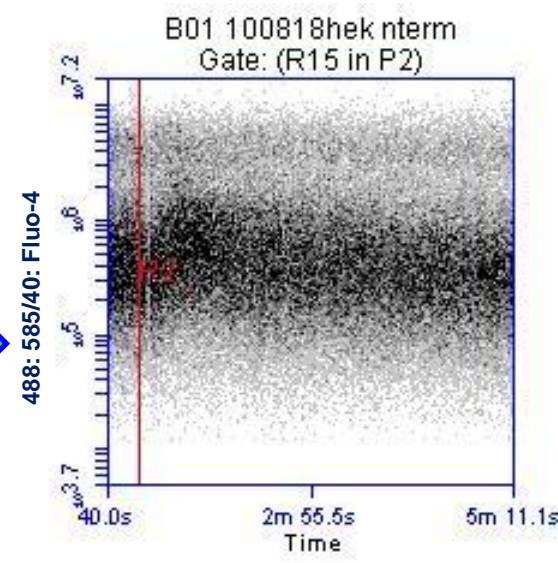
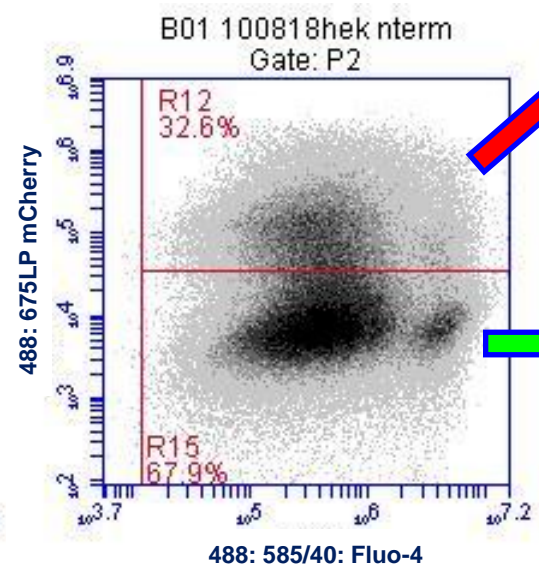
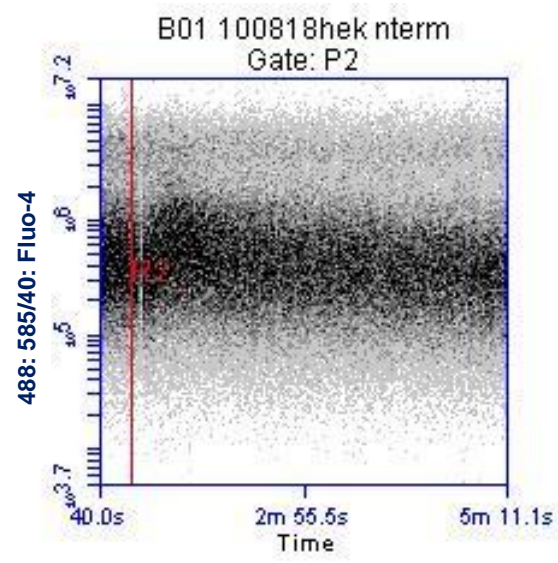


# Calcium flux in some mCherry transfected cell lines

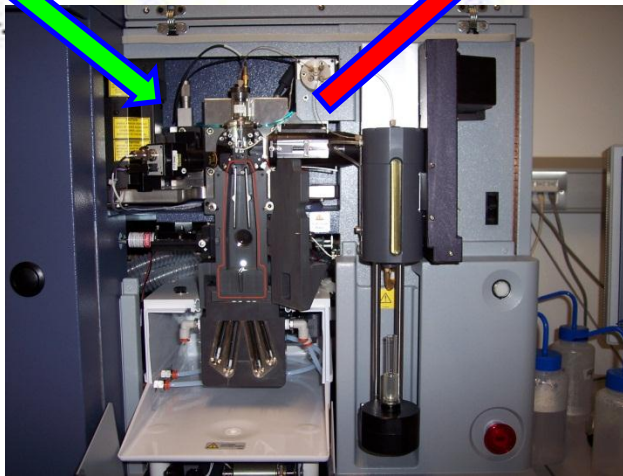
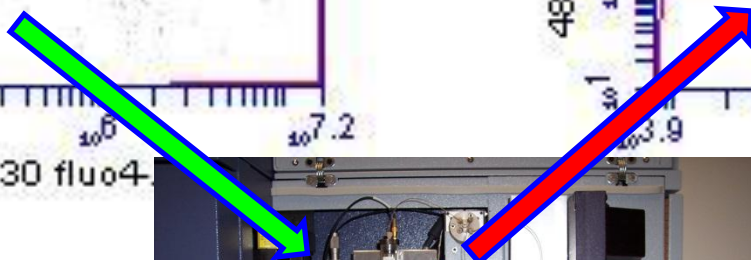
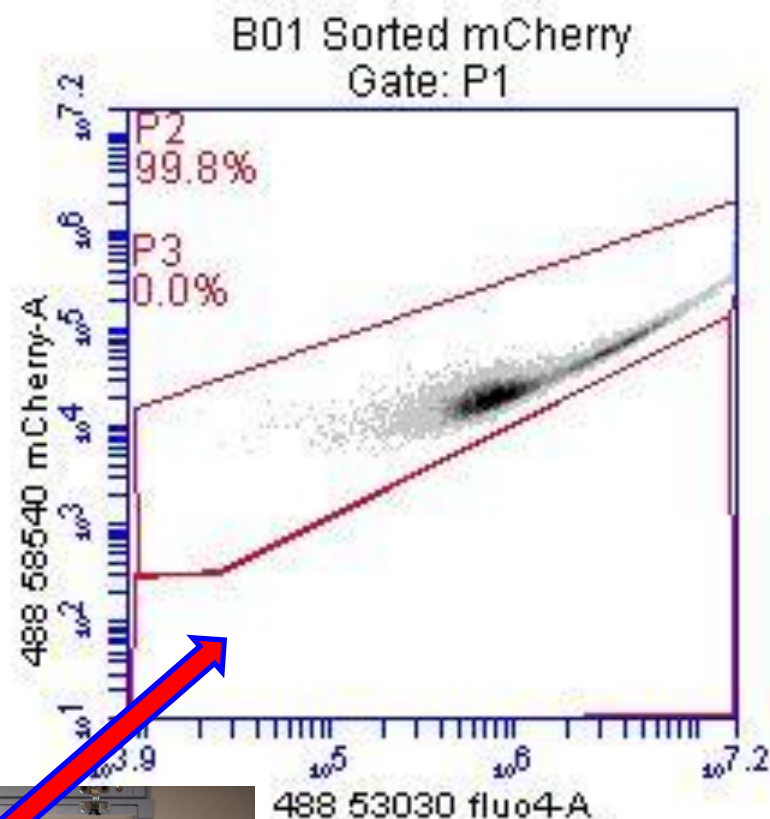
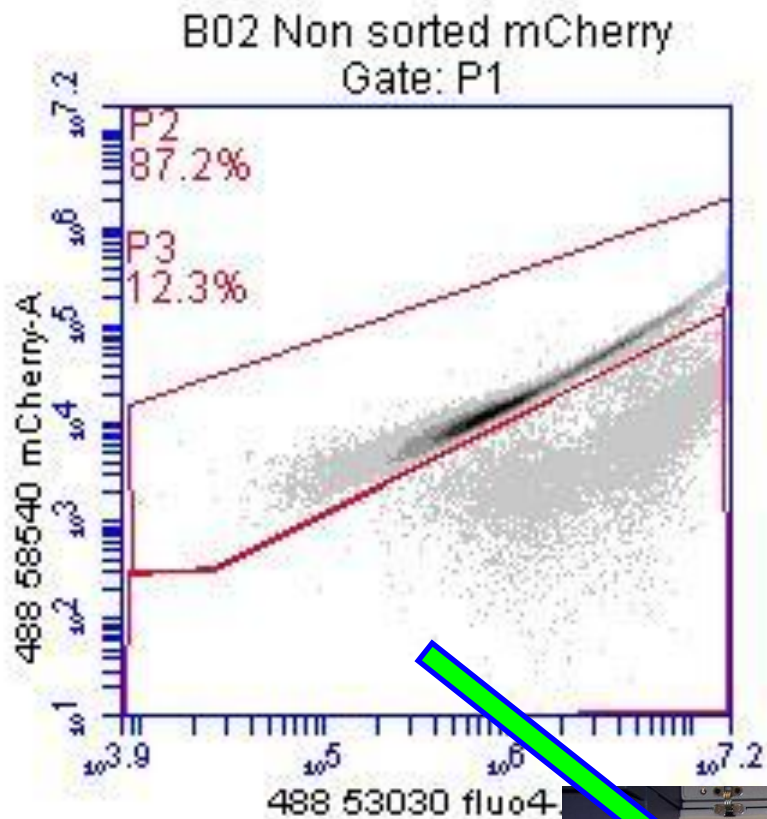
WT



mCherry



# Calcium flux in some mCherry transfected cell lines



# Sharing files and Templates

The screenshot shows the Microsoft Outlook interface. The title bar indicates the current mailbox is 'Inbox - alfonso.blanco@ucd.ie - Microsoft Outlook'. The ribbon includes 'File', 'Home', 'Send / Receive', 'Folder', 'View', and 'Search Tools'. The 'Search Tools' ribbon is active, showing options like 'Ignore', 'Clean Up', 'Junk', 'Delete', 'Reply', 'Reply All', 'Forward', 'More', 'Meeting', 'Viaje a Faro 201...', 'To Manager', 'Team E-mail', and 'Create New'. The left sidebar shows the folder hierarchy, with 'alfonso.blanco@ucd.ie' selected and 'Inbox (9421)' highlighted. The main pane displays an email titled 'Calcium Queries' from Cherise Dunn <DNNCHE002@myuct.ac.za> to Alfonso Blanco, dated Monday, 4/14/2014 12:14 PM. The email content discusses a calcium experiment and asks for advice on cell contamination and fluorescence quantification. A '7 KB' attachment icon is visible at the bottom right of the email content area.

**Calcium Queries**  
Cherise Dunn <DNNCHE002@myuct.ac.za>  
You replied to this message on 4/29/2014 8:42 AM.  
Sent: Mon 4/14/2014 12:14 PM  
To: Alfonso Blanco

Hi Alfonso,

I hope you've been well :) I need some advice please. While my calcium experiment seems to be working now, I found that with one of my cell lines that I was using, I was picking up 3 populations in the FSC vs. SSC plot. Do you think this is cell contamination, debris or a mixture of these? Could there be any other explanation?

Also, I find that the calcium response does not last as long as your examples in the BD paper. I'm thinking this may be cell specific.. have you ever used any other cells apart from glioma cell lines?

My last question was whether you know if there is a way of quantifying the change in Fluo-4 fluorescence before and after drug treatment? I thought you'd be the best source of advice for this. Thank you so much for all your help.

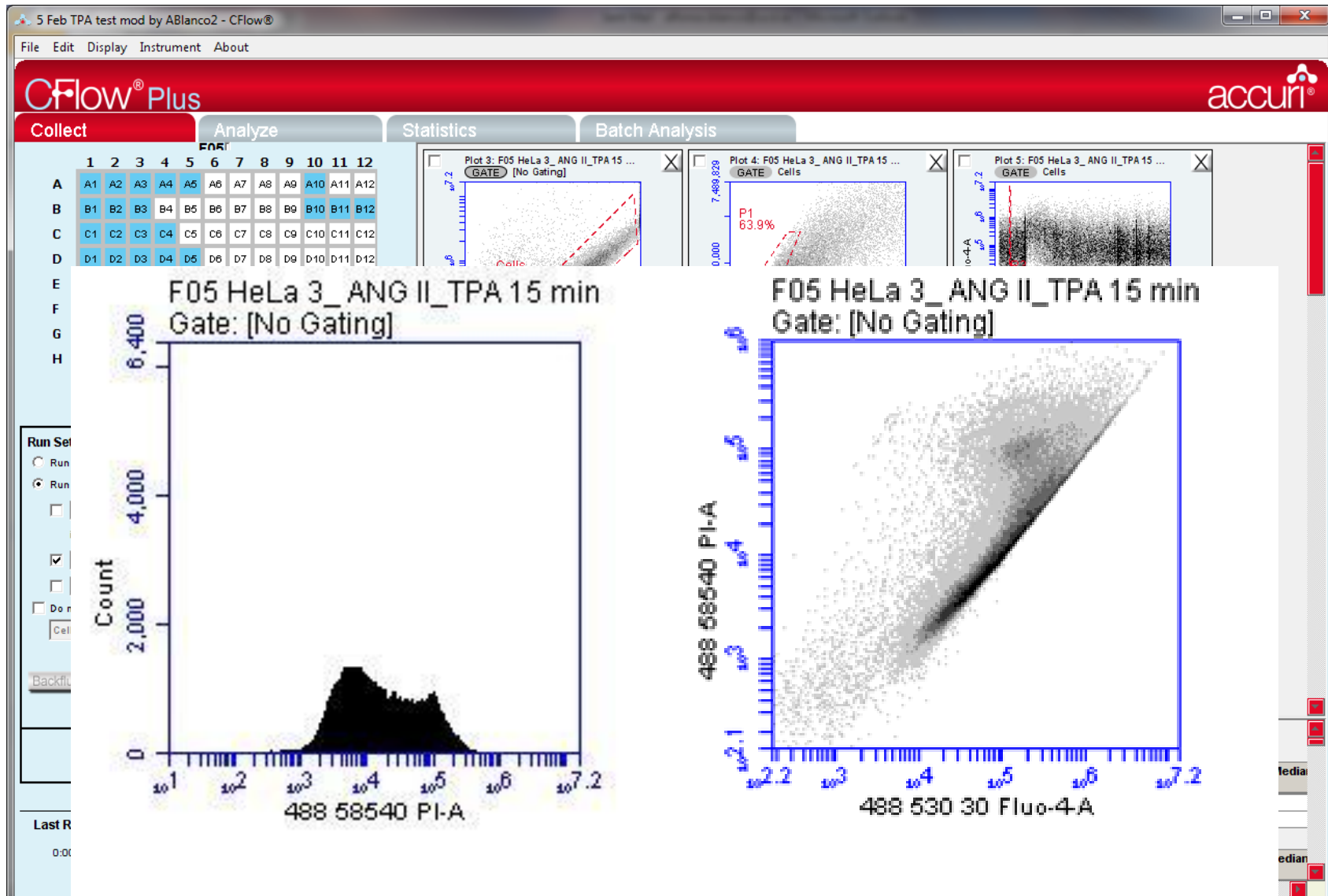
Warm Regards,  
Cherise  
BOOTCAMP (C:)

7 KB

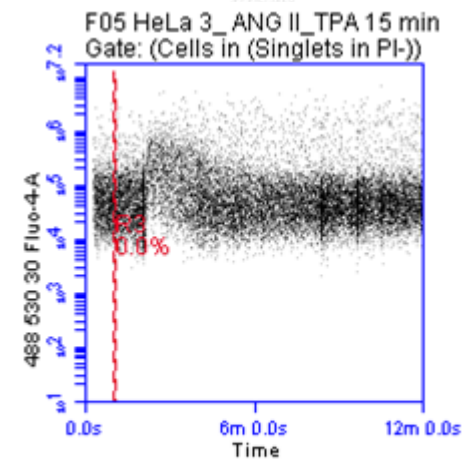
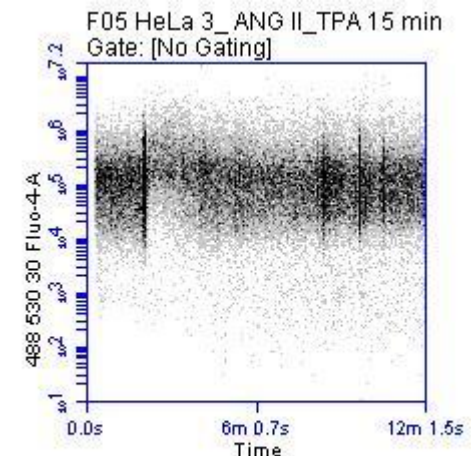
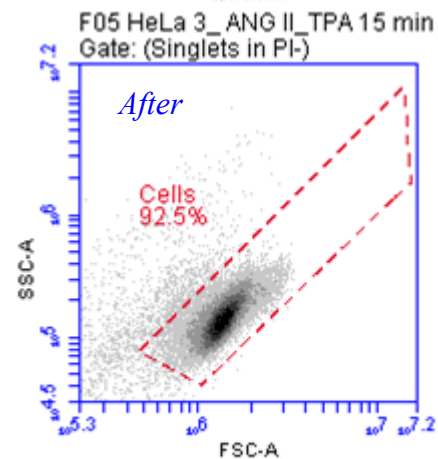
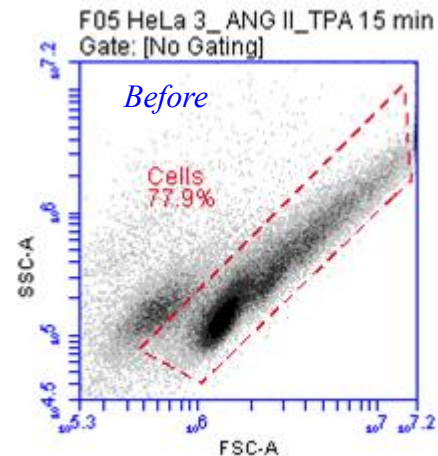
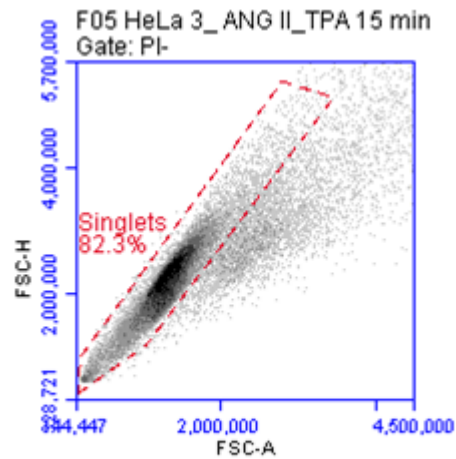
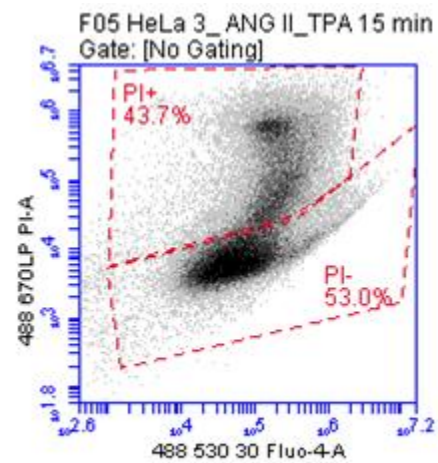


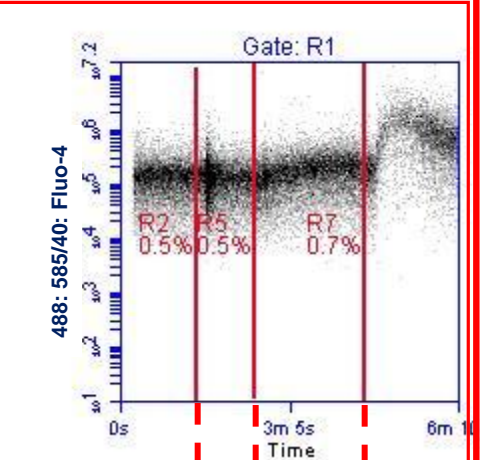
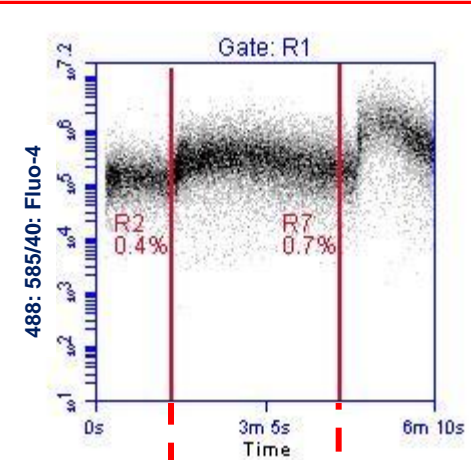
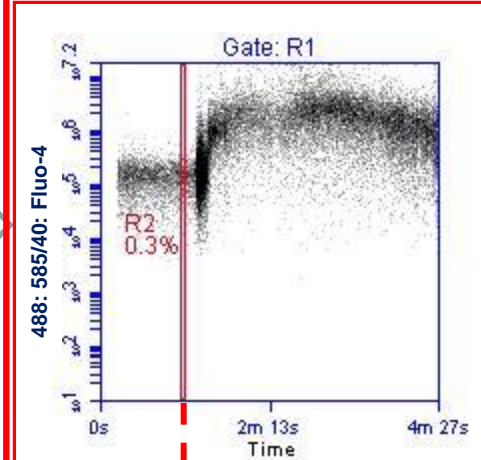
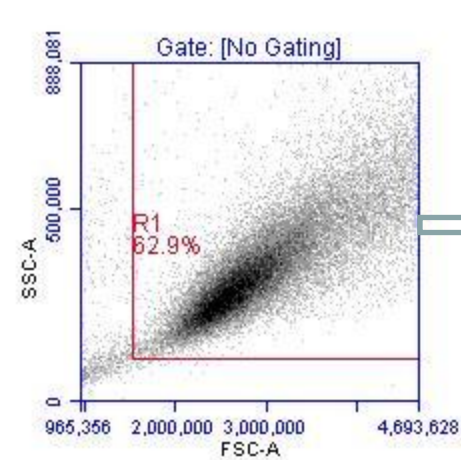


# Sharing files and Templates

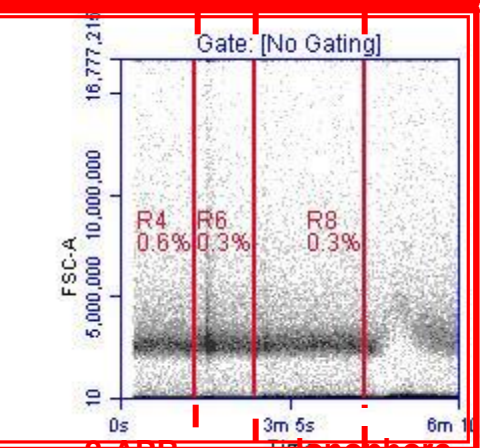
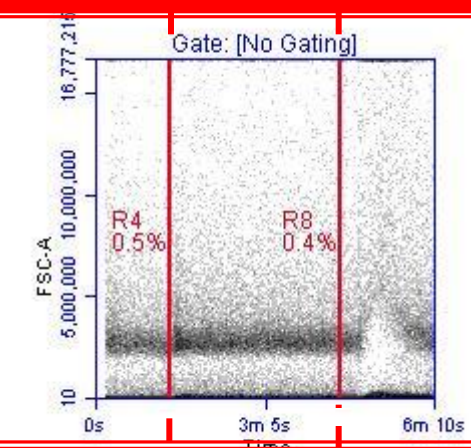
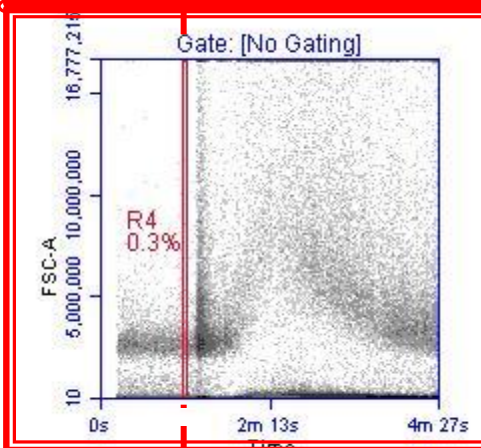
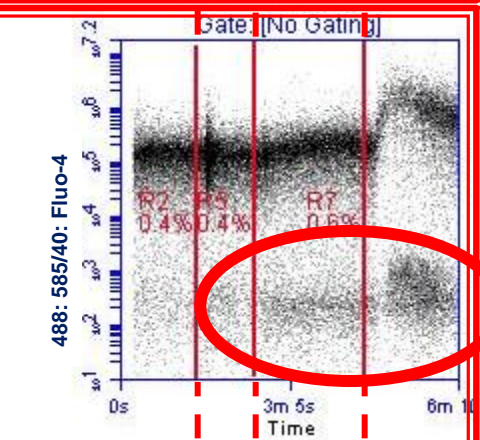
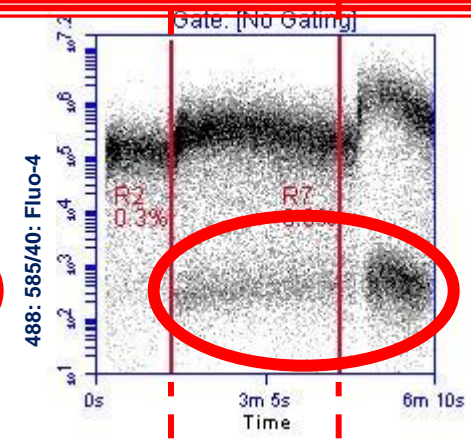
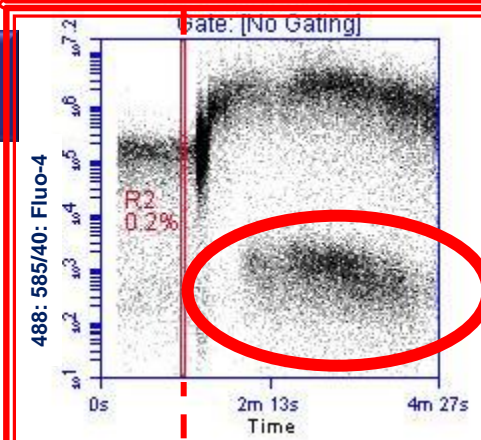


# Sharing files and Templates





# Calcium flux



**Ionophore**

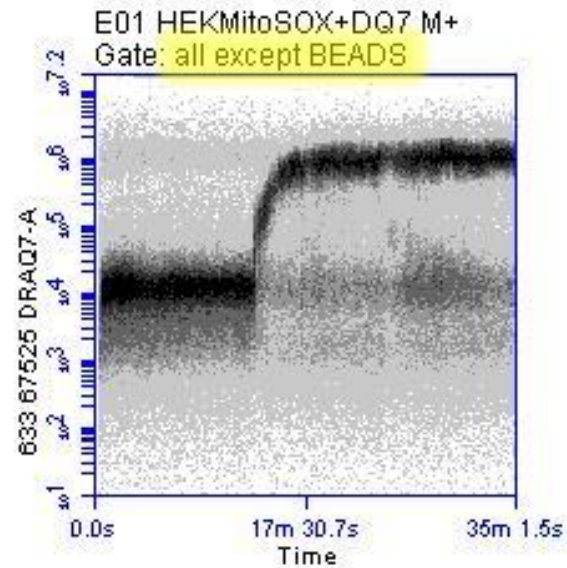
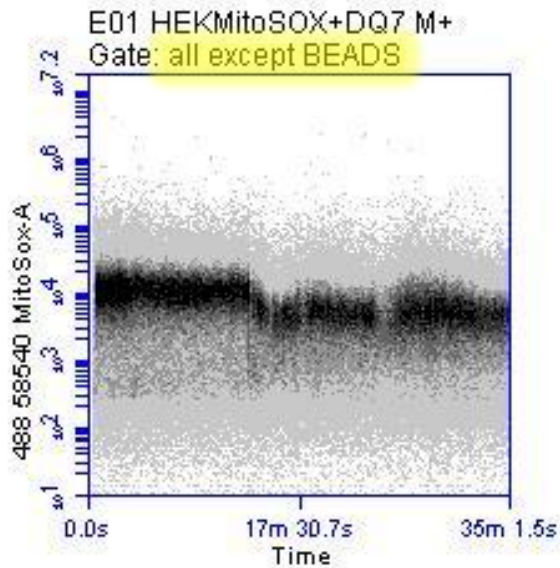
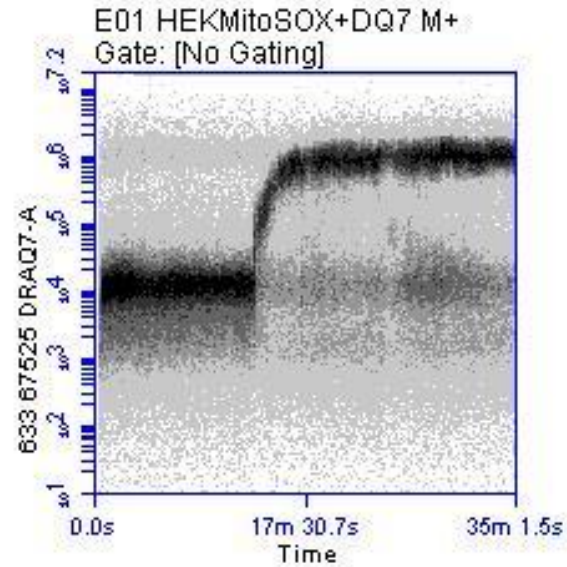
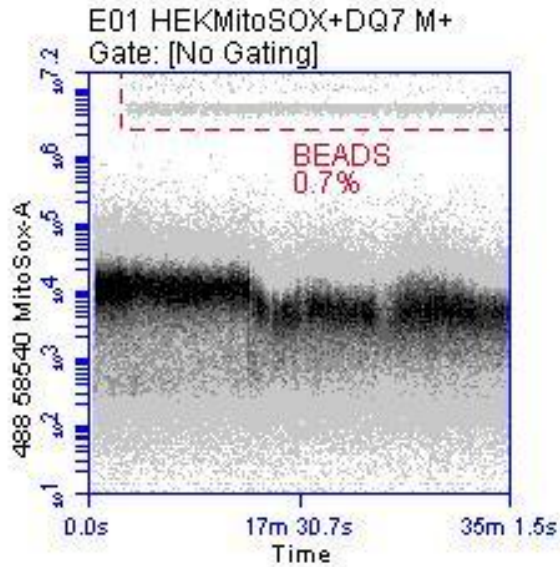
**Thapsigargin Ionophore**

**2-APB**

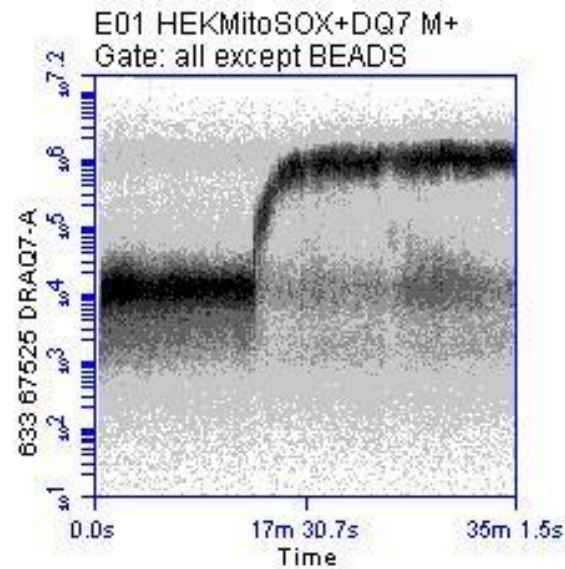
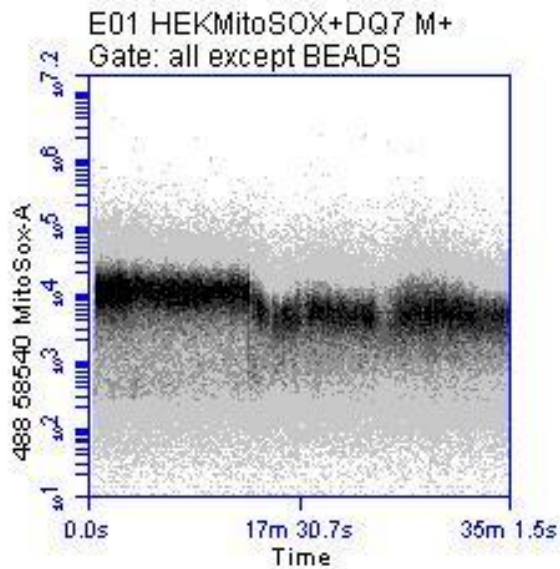
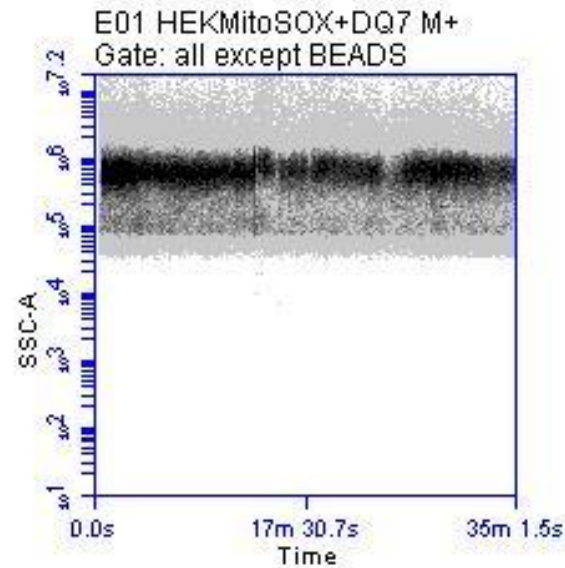
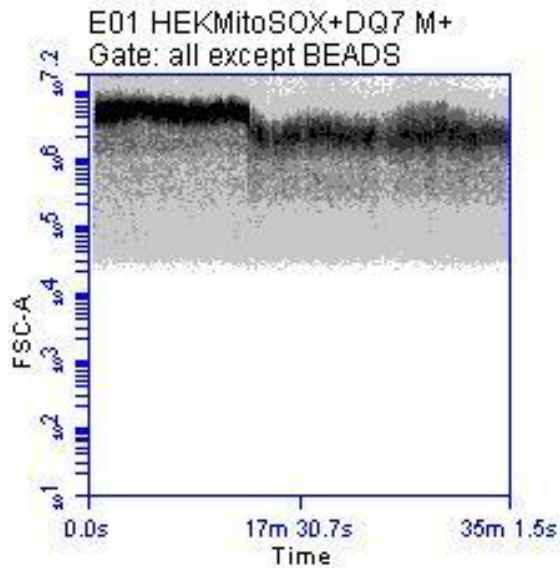
**Thapsigargin**

**Ionophore**

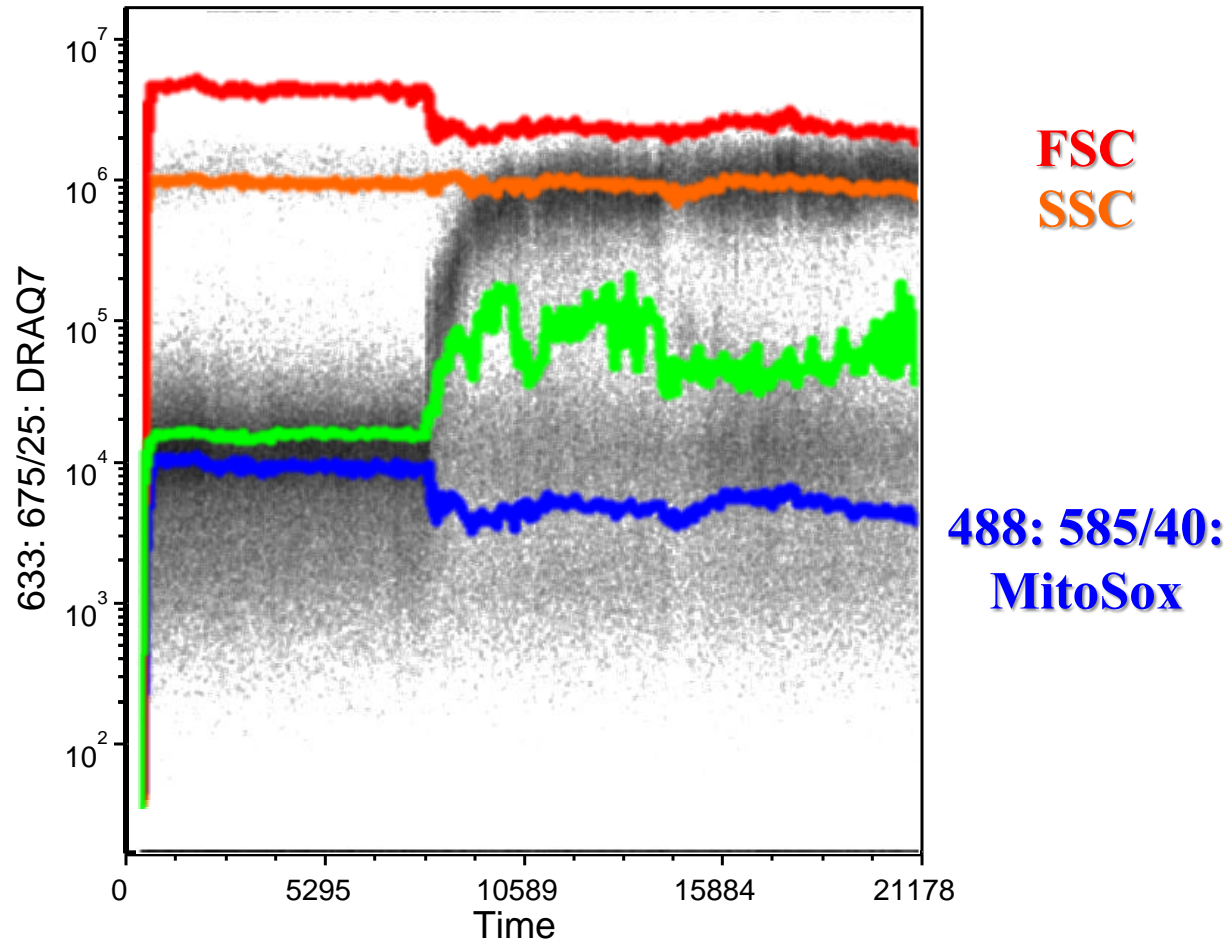
# Apoptosis / Viability



# Apoptosis / Viability



# Apoptosis / Viability



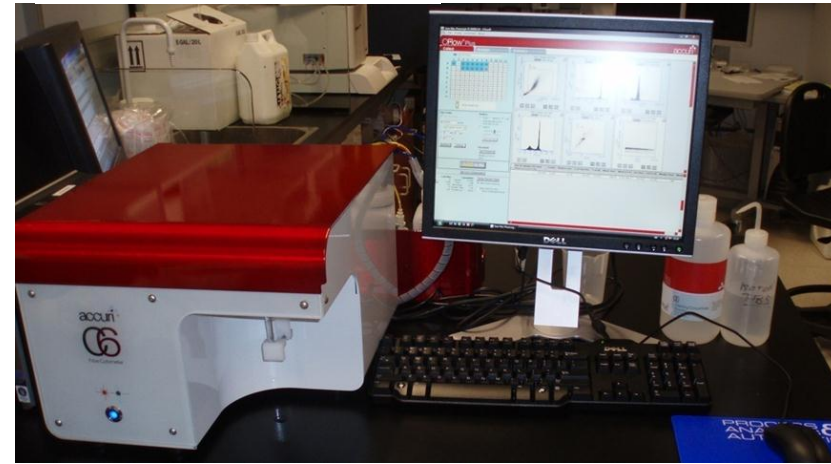
# Collaborations

# VABREMA biosystems MITOPLICATOR

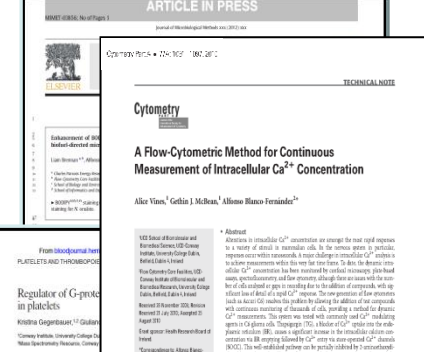
May 2012



## Conference & Tutorials



## Paper per review



## White Paper



## BD Brochure



## Video in Select Science





**VABREMA**  
biosystems

**Department of  
Electrical Engineering**







# VABREMA

biosystems



## MITOPLICATOR

### Mitoplicator – Hybridoma



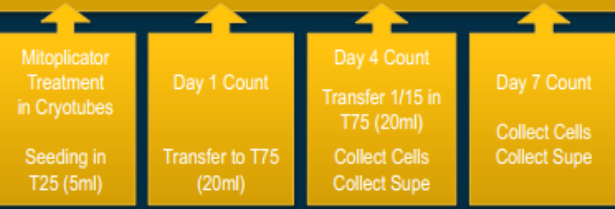
HYBRIDOMA GROWTH

TOTAL AB

AB / CELL

*(Data table & experimental settings were taken from the Hycult report, boxes indicate the median, based on which percentages were calculated)*

Condition:	Treated:	Total # cells (... * 10 <sup>6</sup> alive / dead)								Total µg Ab (IgG2b)		µg Ab / living cell	
		11-11		12-11		15-11		18-11		15-11	18-11	15-11	18-11
1	yes	1.4	0.15	9.0	4.2	12.6	10.6	33.0	36.0	32.6	393.9	2.6	11.9
2	yes	1.4	0.15	10.4	3.8	11.8	14.4	24.0	48.0	33.0	397.4	2.8	16.6
3	yes	1.4	0.15	8.0	0.4	10.6	12.6	24.0	33.0	31.4	397.3	3.0	16.6
4	no	1.4	0.15	9.4	0.6	11.2	12.4	21.0	39.0	31.7	258.1	2.8	12.3
5	no	1.4	0.15	12.0	0.4	8.4	13.8	15.0	30.0	31.2	230.8	3.7	15.4
6	no	1.4	0.15	11.0	2.4	10.4	17.2	9.0	33.0	31.8	256.4	3.1	28.5



**TREATED CELLS VS CONTROLS** → Slight increase in dead cells | 13% more cells at day 4 | 60% more cells at day 7 | 55% more Ab at day 7 | Ab Production / Cell similar at day 7

... treatment of this hybridoma seems to have a positive effect on the speed of cell division, no effect on protein production per cell and the protein produced is functional. "

*(Results & quote: courtesy of Helma Rutjes, PhD, Head of R&D, HyCult)*





## Stimulation of Capacitative Calcium Entry in HL-60 Cells by Nanosecond Pulsed Electric Fields\*

Received for publication, October 9, 2003, and in revised form, March 2, 2004  
Published, JBC Papers in Press, March 16, 2004, DOI 10.1074/jbc.M311135200

Jody A. Winter<sup>1</sup>, Peter T. Brackmores<sup>1</sup>, Karl H. Schoenbach<sup>1</sup>, and Stephen G. Beebe<sup>2</sup>

## Lipid nanopores can form a stable, ion channel-like conduction pathway in cell membrane

Andrei G. Pakhomov<sup>a,\*</sup>, Angela M. Bowman<sup>a</sup>, Bennett L. Ibey<sup>b</sup>, Franck M. Andre<sup>a</sup>,  
Olga N. Pakhomova<sup>a</sup>, Karl H. Schoenbach<sup>a</sup>

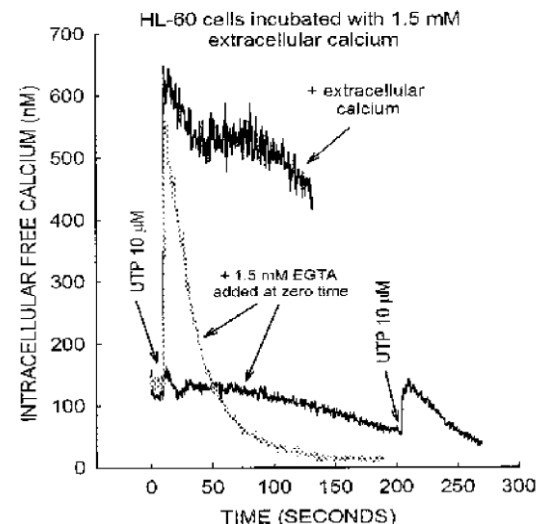
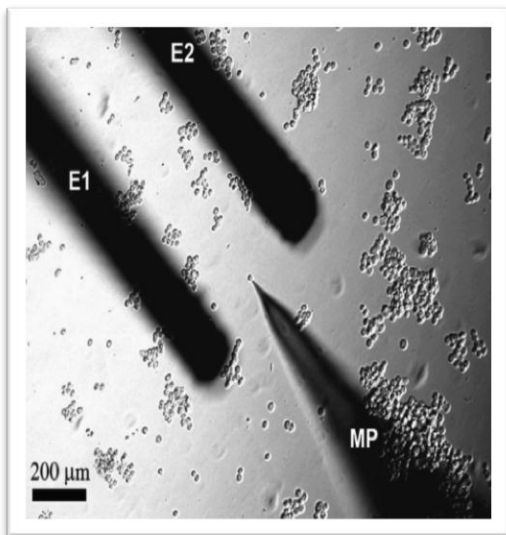
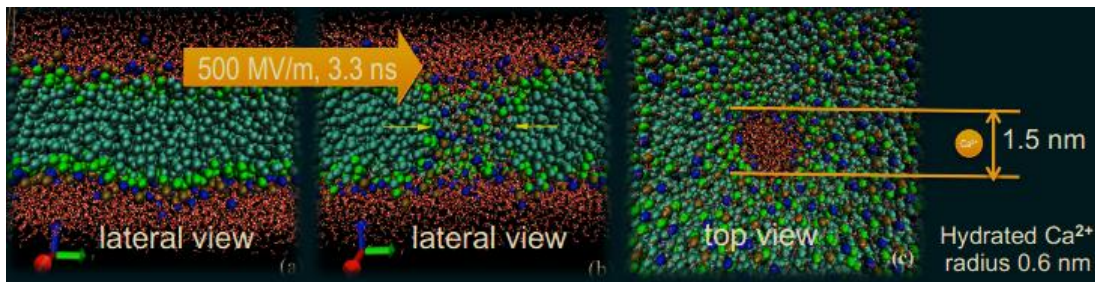
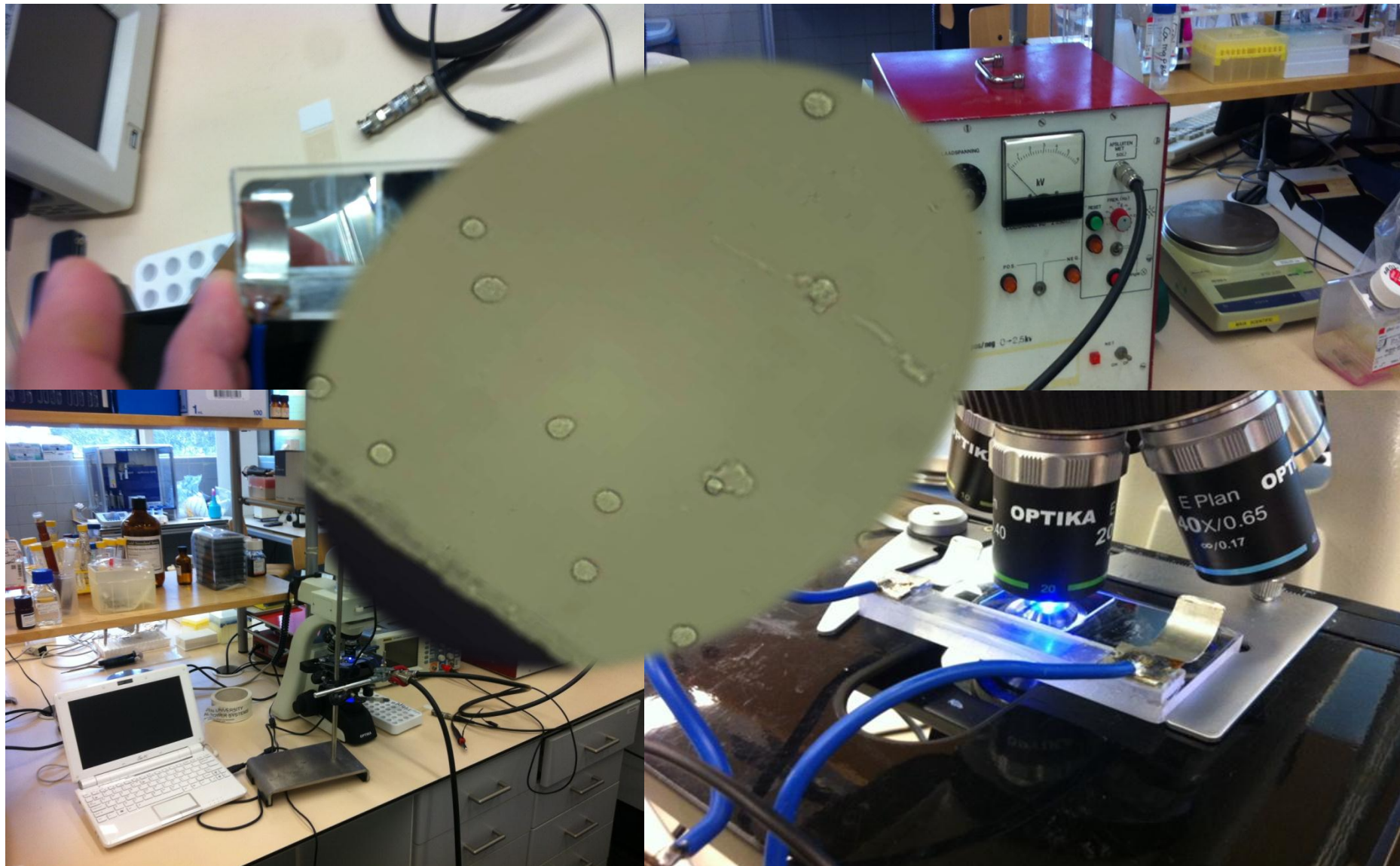


FIG. 2. HL-60 cells, stimulated by nsPEF, respond by releasing internal calcium stores. HL-60 cells incubated with extracellular calcium (1.5 mM) respond to nsPEF with a substantial and relatively prolonged release of  $[Ca^{2+}]_i$ . Cells incubated in calcium-deficient media with added 1.5 mM EGTA respond with similar magnitude; however, the response is quite transient (gray trace). This response represents release of internal stores of calcium. HL-60 cells incubated in the same calcium-deficient conditions were allowed to lose some of their calcium before treating with UTP. The gradual loss of calcium before treatment strongly suggests that calcium influx through the plasma membrane is not occurring and that the response seen to UTP is because of internal store release.



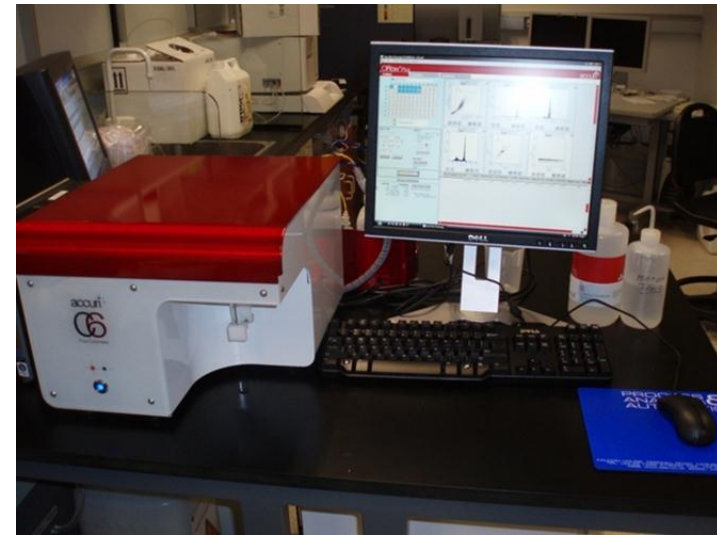


VABREMA  
biosystems

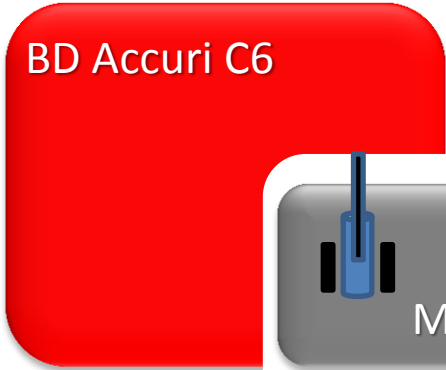




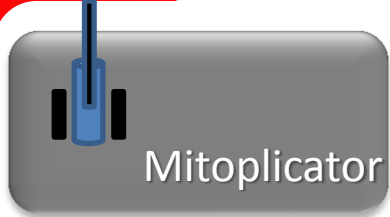
# MITOPLICATOR



## Initial Idea



BD Accuri C6



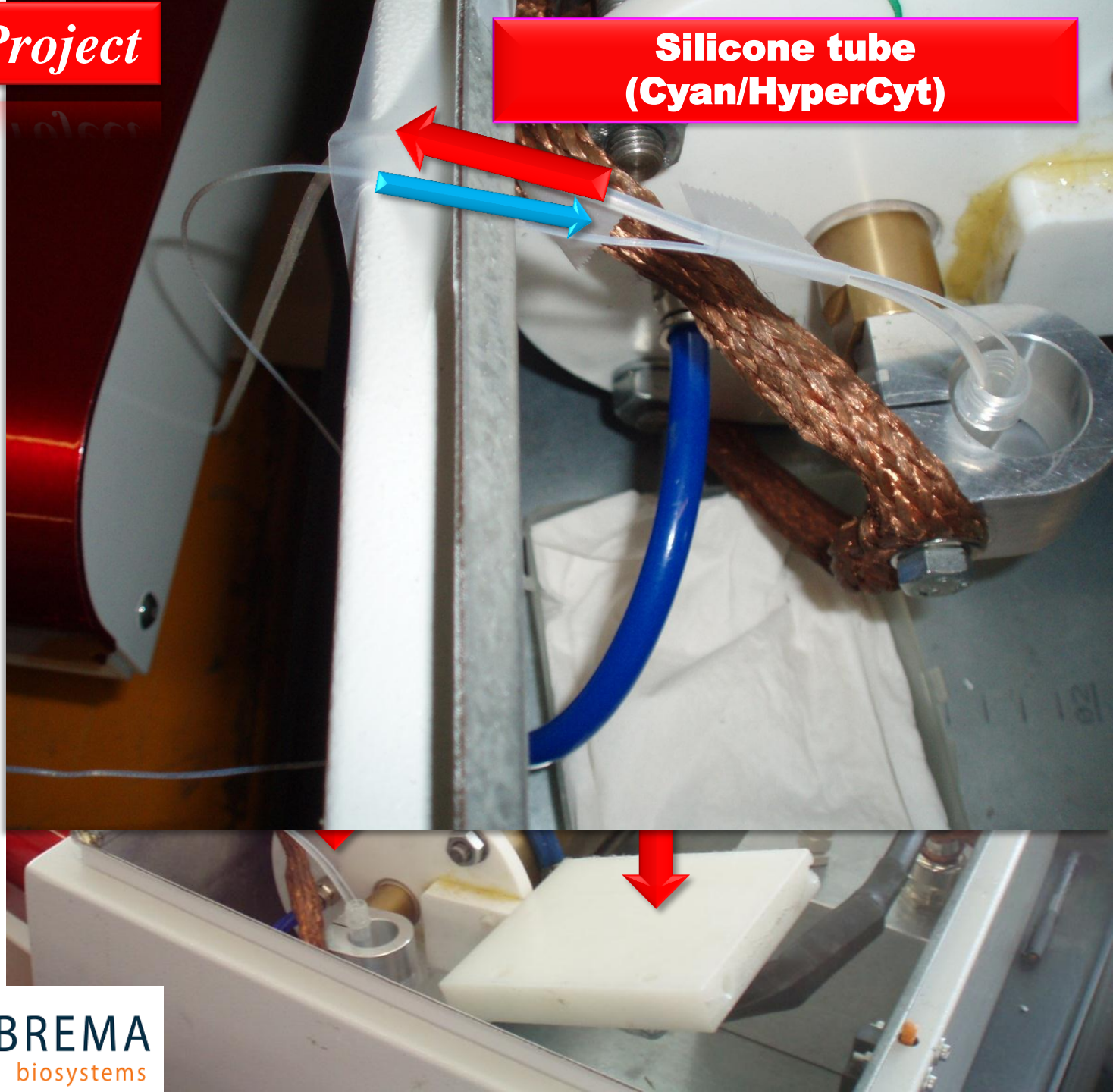
Mitoplicator



VABREMA  
biosystems

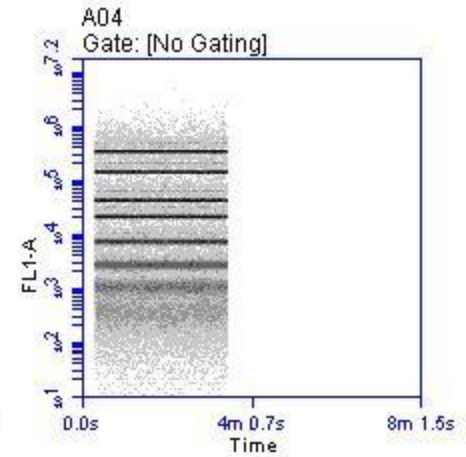
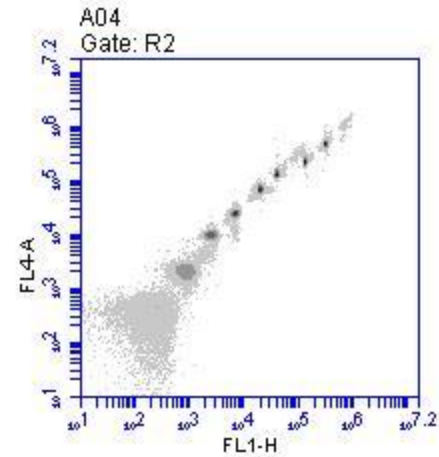
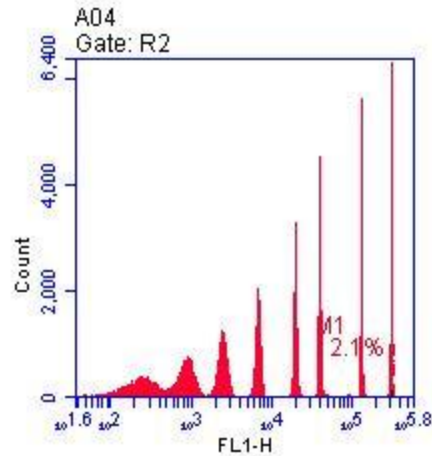
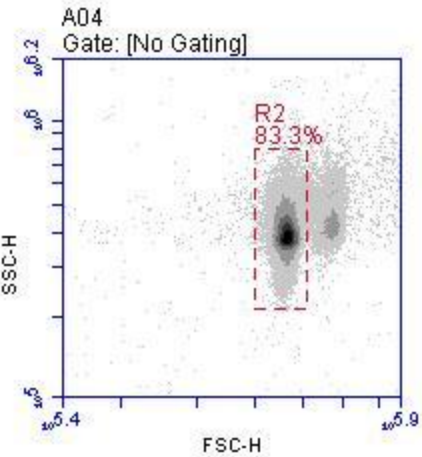
# Mitoplicator Project

**Silicone tube  
(Cyan/HyperCyt)**

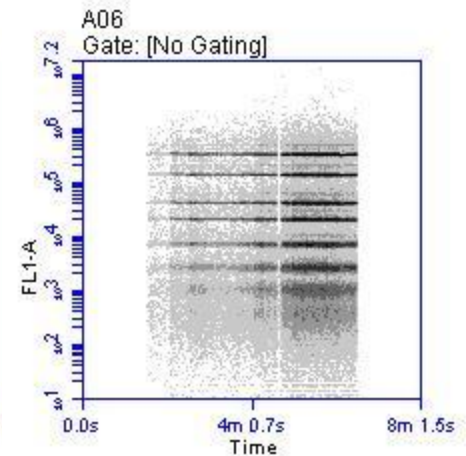
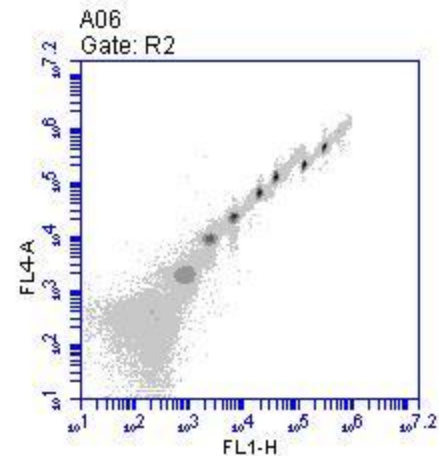
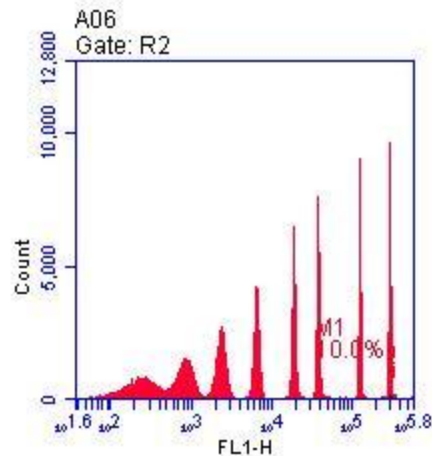
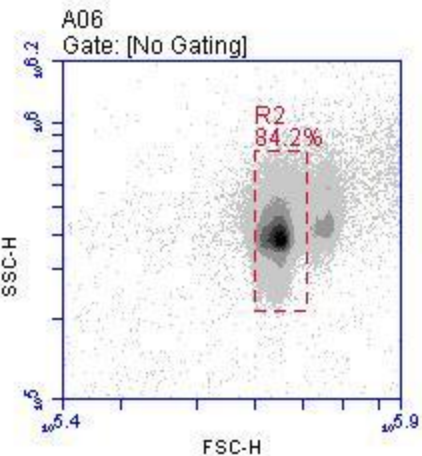


**VABREMA**  
biosystems

# One of the 1<sup>st</sup> experiments: Stability...



## 8 Peak beads in tube holder of the Accuri C6



## 8 Peak beads changing speeds in Mitoplicator



VABREMA  
biosystems

# Mitoplicator Project



**VABREMA**  
biosystems

# Mitoplicator Project

- CFSE – Proliferation
- Fluo4 – Calcium
- DRAQ7 – Viability
- Mitosox – ROS
- BODIPY – Lipids
- GFP –
- ...

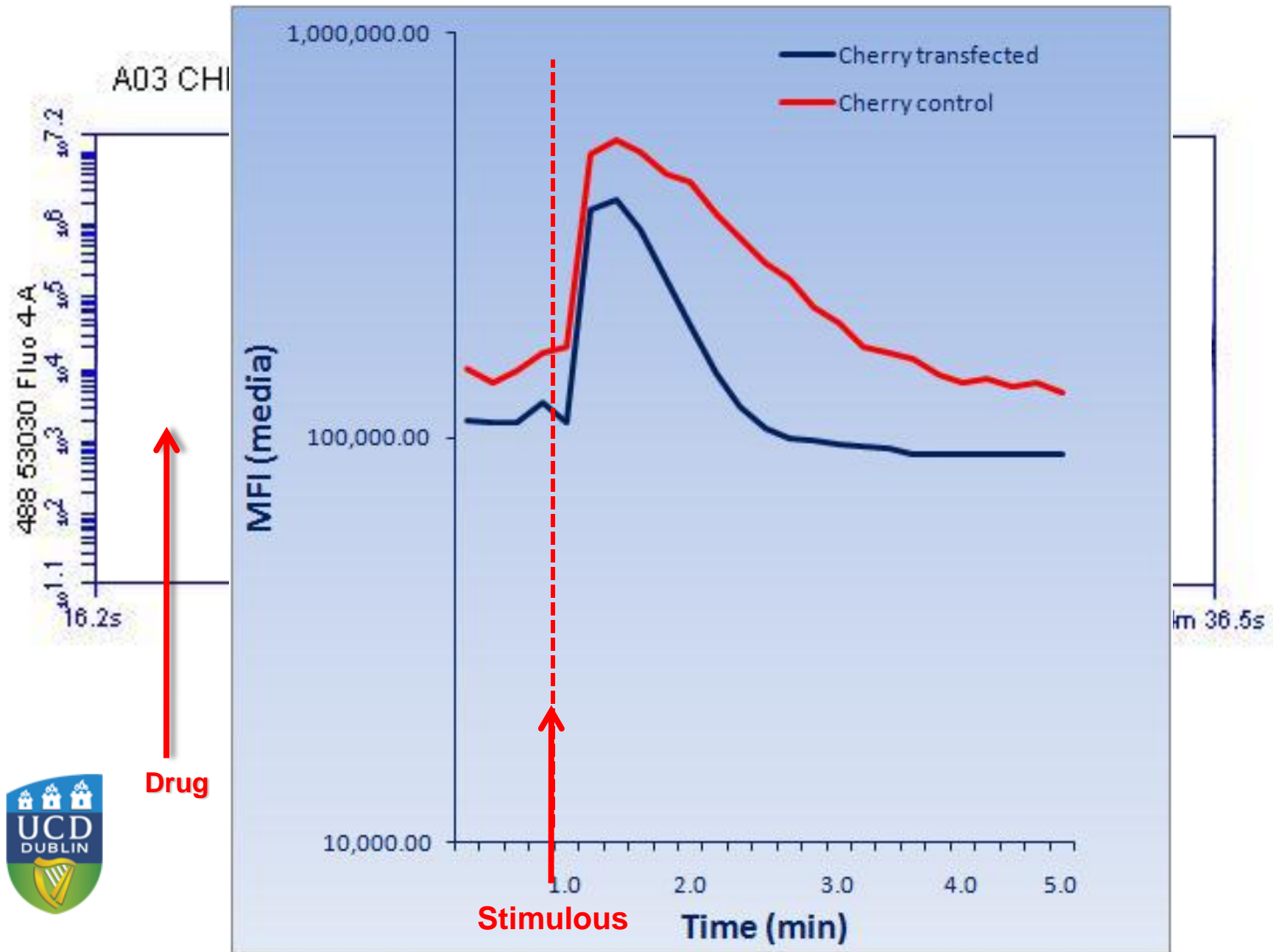
- C6 glial
- HEK
- CHO
- Jurkats
- Platelets
- ...



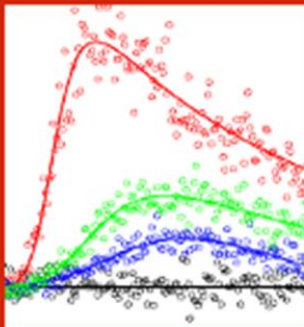
VABREMA  
biosystems



# Calcium flux in some mCherry transfected cell lines



# Kinetic analysis using FacsKin software



FacsKin

*Kinetic Analysis of Flow Cytometry Data*



**ISAC**

International Society for Advancement of Cytometry

ISAC Scholars Program

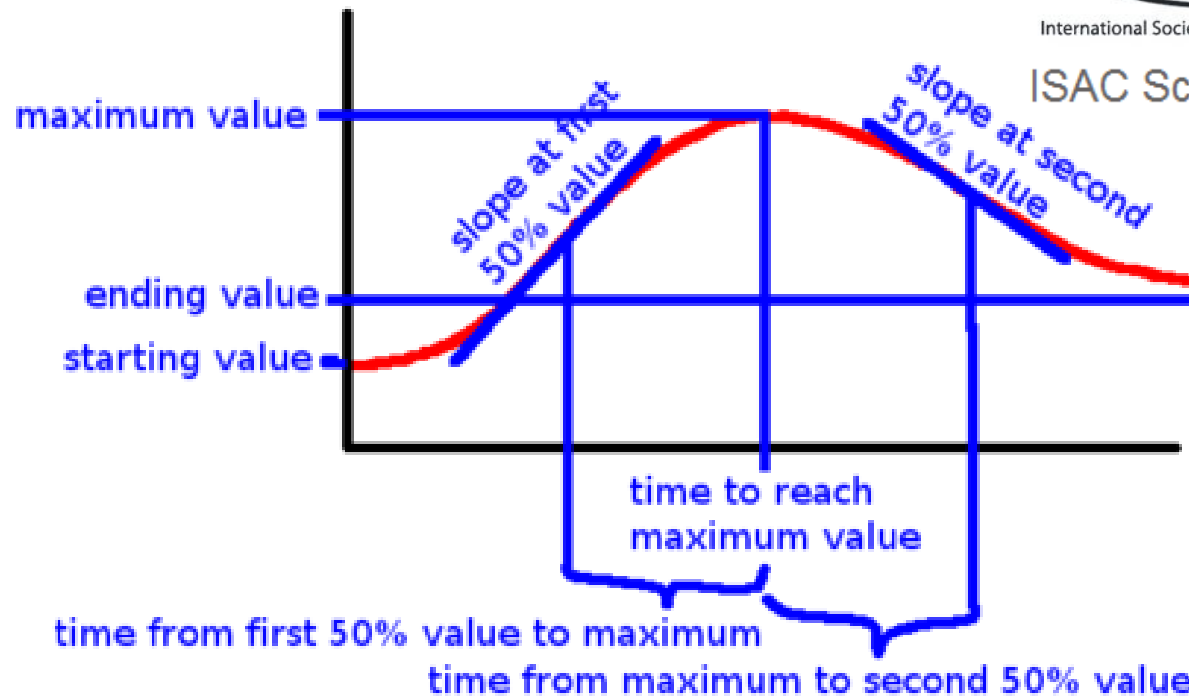
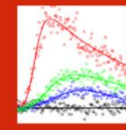
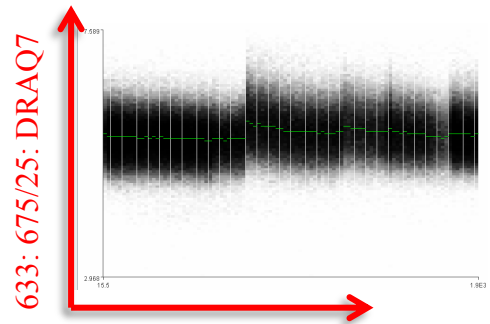
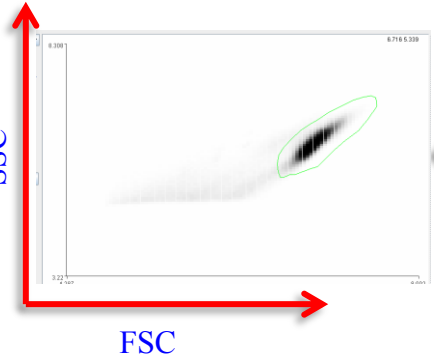


Figure 12. Parameters of dlogist+ function

# Kinetic analysis using FacsKin software

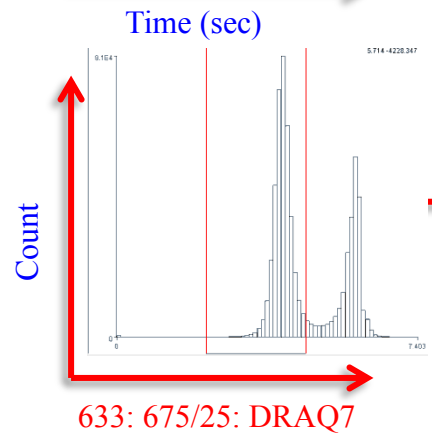


FacsKin  
Kinetic Analysis of Flow Cytometry Data

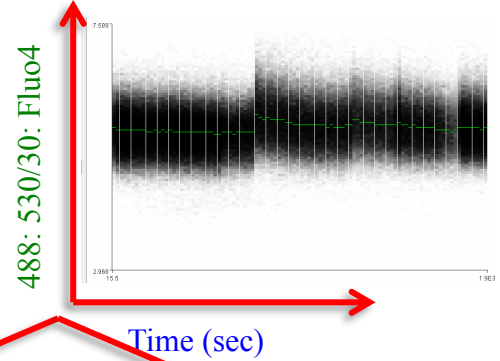


**Gate one:** Contains 120s baseline and 600s treatment/control (0-720s) (Mitoplicator on/off)

**Analysed parameters:**  
DRAQ7

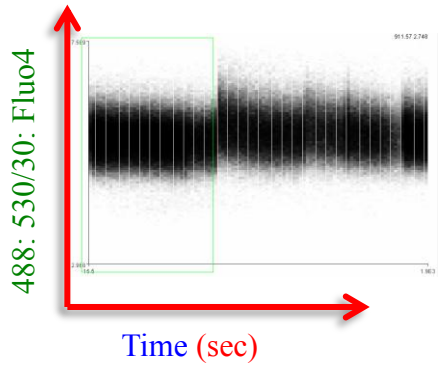


**Gate out:**  
DRAQ7+



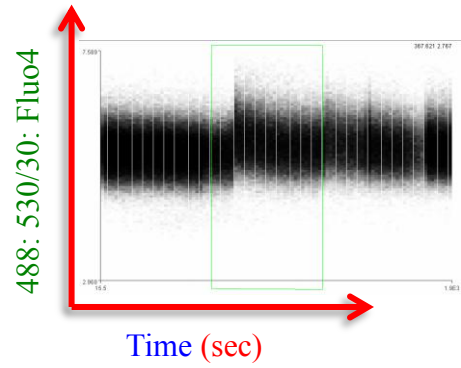
**Gate two:** Contains 120s baseline and 600s treatment/control (0-720s) (Mitoplicator on/off)

**Analysed parameters:**  
FLUO4, FSC SSC

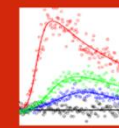


**Gate three:** Contains 100s baseline and 300s quenching time after the addition of 6ul 1mg/ml ionomycin (600 -1020s)

**Analysed parameters:**  
FLUO4, FSC SSC

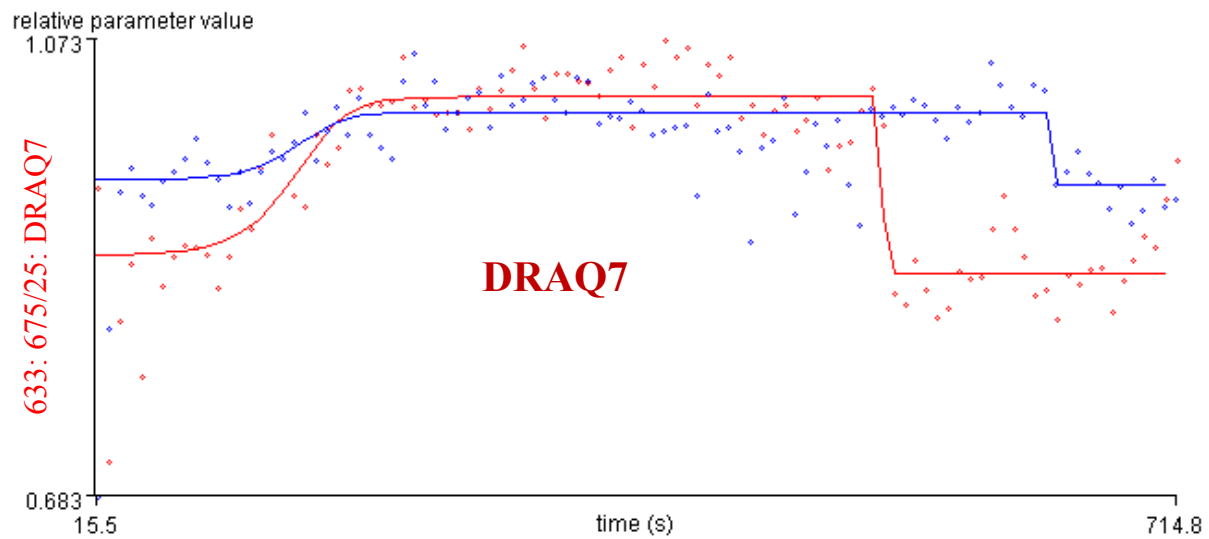
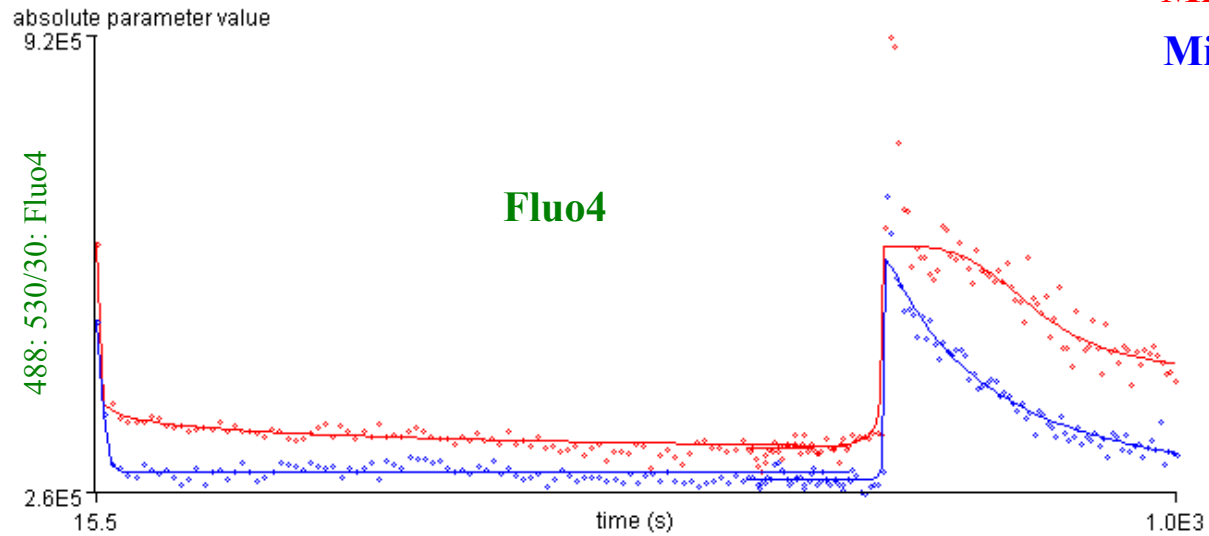


# Kinetic analysis using FacsKin software

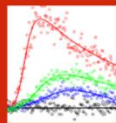


FacsKin  
Kinetic Analysis of Flow Cytometry Data

**Mitoplicator OFF**  
**Mitoplicator ON**



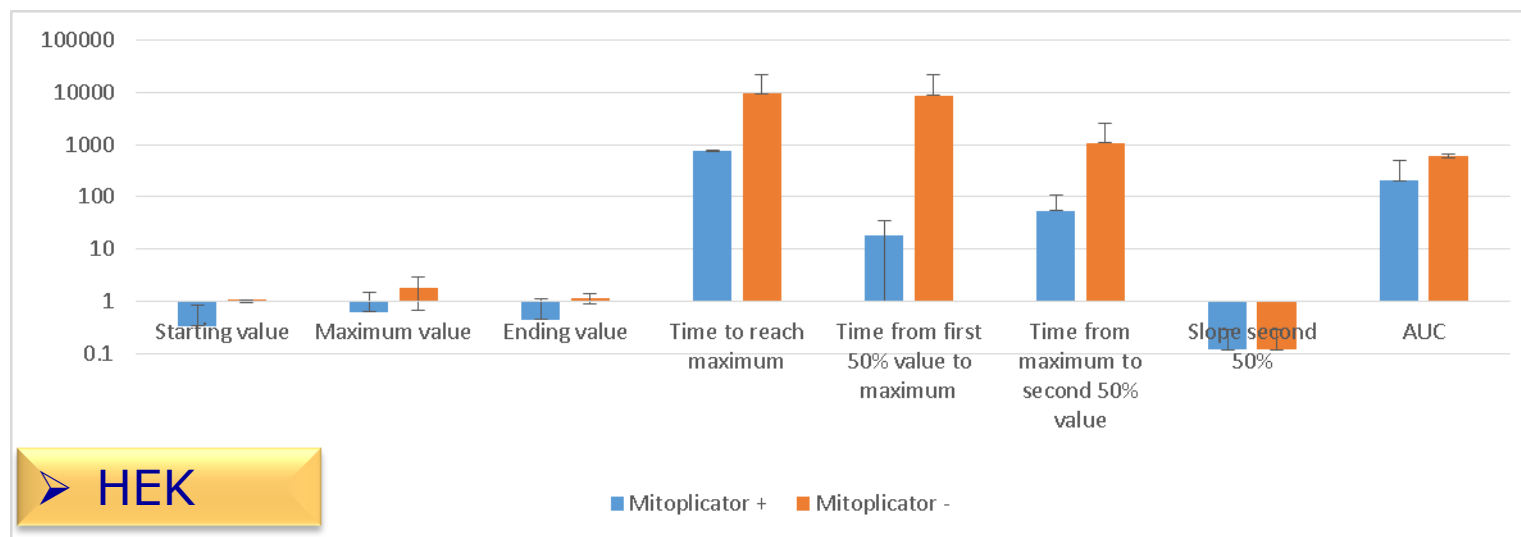
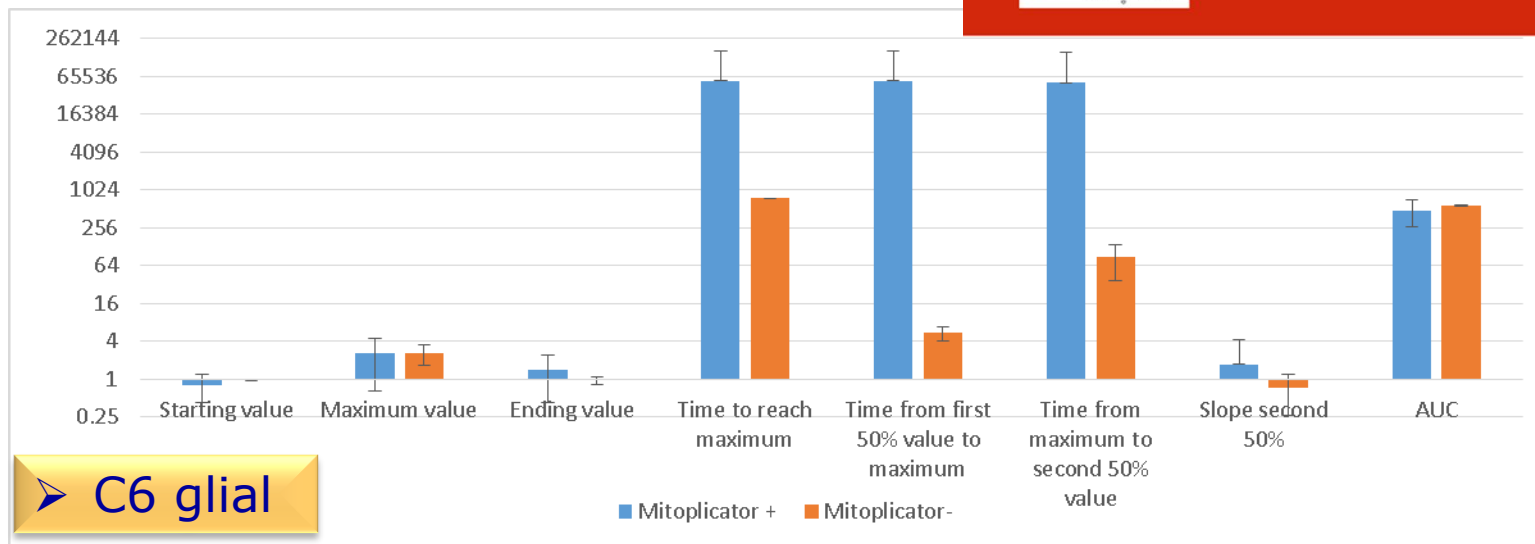
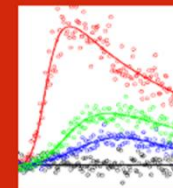
# Kinetic analysis using FacsKin software



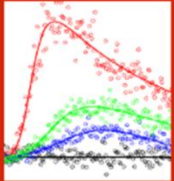
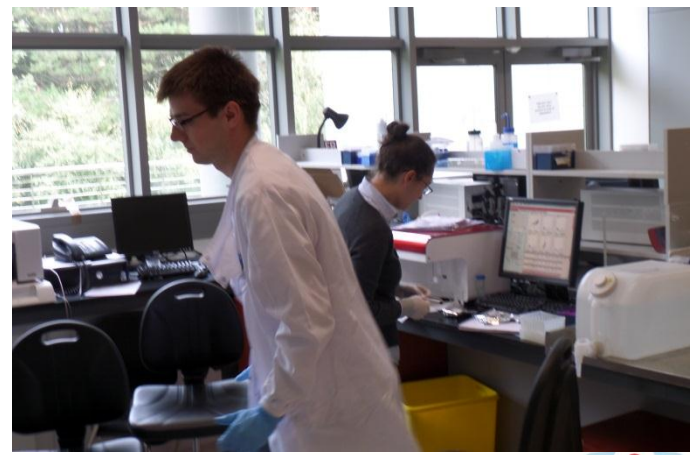
FacsKin  
Kinetic Analysis of Flow Cytometry Data

Cellular parameter	Mitoplicator treatment	Starting value	Maximum value	Ending value	Time to reach maximum	Time from maximum to second 50% value	Slope at first 50% value	AUC
I.C. calcium (0-720 s)	+	0,99 [0,49-1,07]	1,42 [1,02-1,81]	0,49 [0,00-0,96]	138,64 [0,01-41774,00]	82,24 [17,34-2527,98]	0,11 [0,01-0,69]	608,47 [327,72-613,18]
	-	1,03 [1,02-1,10]	1,37 [1,04-1,76]	0,78 [0,36-1,00]	110,34 [12,08-255,49]	330,91 [9,53-500,67]	0,11 [0,02-0,44]	611,61 [601,92-623,05]
	P=	0,69	0,89	0,49	0,89	0,20	1,00	0,49
I.C. calcium (620-1020 s)	+	1,00 [0,92-1,03]	2,35 [1,98-5,02]	1,72 [0,58-2,68]	751,69 [748,49-771,29]	67,12 [38,06-140,71]	0,67 [0,19-7,49]	599,74 [549,39-621,62]
	-	0,99 [0,97-1,01]	1,89 [1,03-3,78]	0,94 [0,86-41646,00]	752,00 [749,92-754,08]	70,97 [29,35-123,40]	0,74 [0,52-1,46]	591,73 [578,91-607,59]
	P=	0,89	0,49	0,89	1,00	0,89	1,00	0,89
SSC (0-720s)	+	1,01 [0,94-1,03]	1,03 [1,01-1,05]	0,96 [0,88-0,98]	318,24 [0,00-467,47]	166,51 [66,07-265,25]	0,01 [0,00-0,04]	597,72 [592,20-603,35]
	-	1,02 [0,95-1,04]	1,03 [1,02-1,05]	0,91 [0,36-0,95]	203,21 [15,50-451,66]	208,85 [127,30-1895,53]	0,15 [0,00-0,26]	600,47 [599,39-604,29]
	P=	0,69	1,00	0,34	0,89	0,34	0,34	0,49
SSC (620-1020s)	+	1,00 [0,99-1,01]	1,02 [1,01-41646,00]	0,99 [0,92-1,06]	825,72 [710,93-948,71]	145,97 [1,60-17905,04]	0,09 [0,00-0,21]	597,01 [592,13-603,02]
	-	0,99 [0,98-1,00]	1,06 [1,04-1,10]	1,02 [1,00-1,07]	967,99 [841,52-26407,23]	29,84 [1,29-23846,09]	0,02 [0,00-0,08]	593,56 [588,79-598,02]
	P=	0,34	0,34	0,20	0,11	0,69	0,34	0,34
FSC (0-720s)	+	1,01 [0,90-1,02]	1,04 [1,02-41640,00]	0,95 [0,90-41642,00]	340,39 [250,42-552946,52]	129,13 [0,00-19262,83]	0,06 [0,00-0,21]	603,04 [583,39-611,28]
	-	1,01 [0,93-1,02]	1,02 [1,01-1,04]	0,95 [0,94-0,97]	206,13 [0,00-255,22]	325,43 [0,00-487,41]	0,00 [0,00-0,08]	599,99 [599,38-600,77]
	P=	0,69	0,25	1,00	0,06	0,69	0,49	1,00
FSC (620-1020s)	+	0,99 [0,99-1,00]	1,04 [0,99-1641,00]	0,95 [0,92-1,00]	718,60 [0,01-865,07]	562,66 [20,33-41833,00]	0,17 [0,00-0,52]	595,18 [592,25-599,96]
	-	0,99 [0,97-1,00]	1,05 [1,00-41645,00]	0,99 [0,98-1,04]	842,28 [746,04-955,50]	102,79 [37,45-137,67]	0,13 [0,06-42]	595,46 [580,96-01,10]
	P=	0,89	0,77	0,20	0,34	0,34	0,89	0,89
Draq 7 (0-720s)	+	0,95 [0,89-0,99]	1,02 [1,01-1,03]	0,98 [0,87-1,02]	371,82 [258,03-435,34]	167,58 [126,82-11311,25]	0,00 [0,00-0,00]	597,86 [584,82-602,63]
	-	0,95 [0,86-0,99]	1,02 [1,00-1,04]	0,94 [0,79-0,97]	409,59 [332,14-737,63]	533,43 [170,16-1528,91]	0,00 [0,00-0,00]	595,58 [581,24-601,18]
	P=	0,88	0,89	0,34	0,49	0,49	0,66	0,69

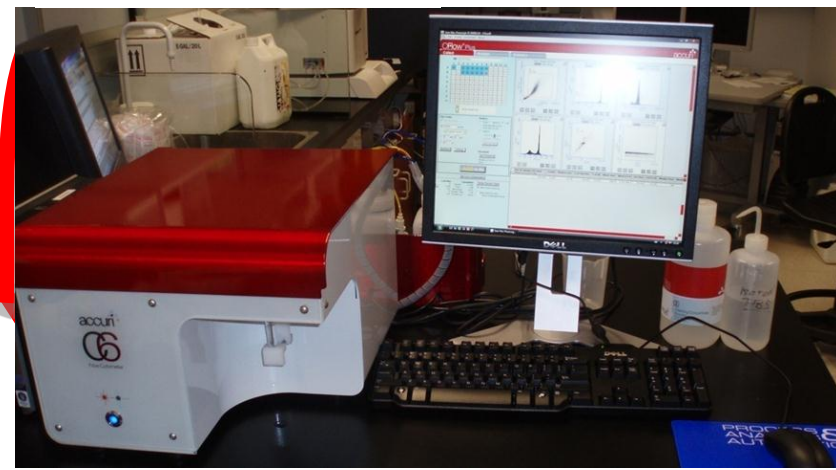




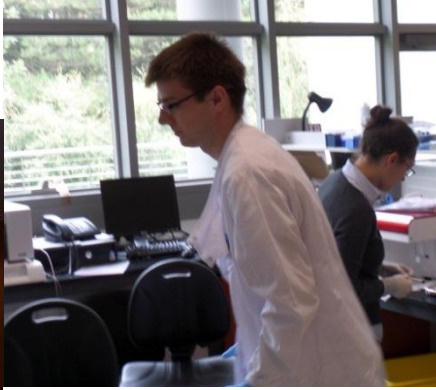
# Collaborations



**FacKin**  
*Kinetic Analysis of Flow Cytometry Data*



# Acknowledge





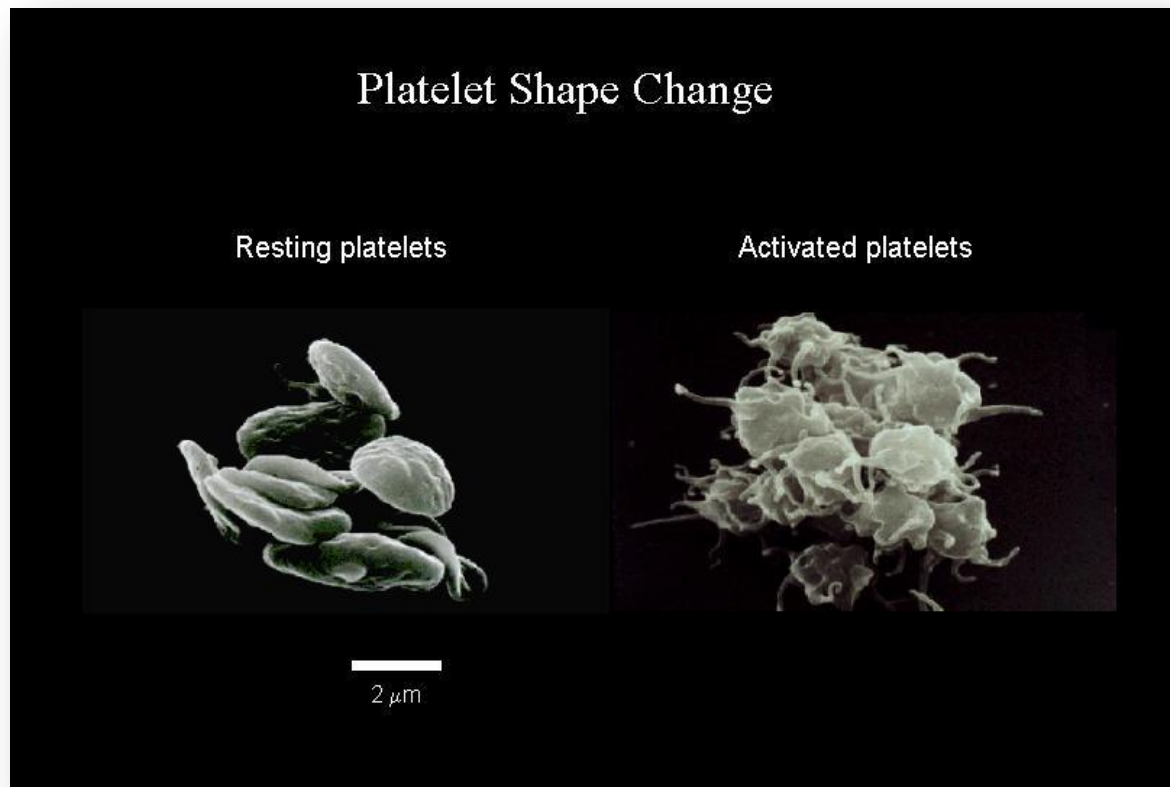
# Cell Biology Applications of Real- Time Flow Cytometry Platelets

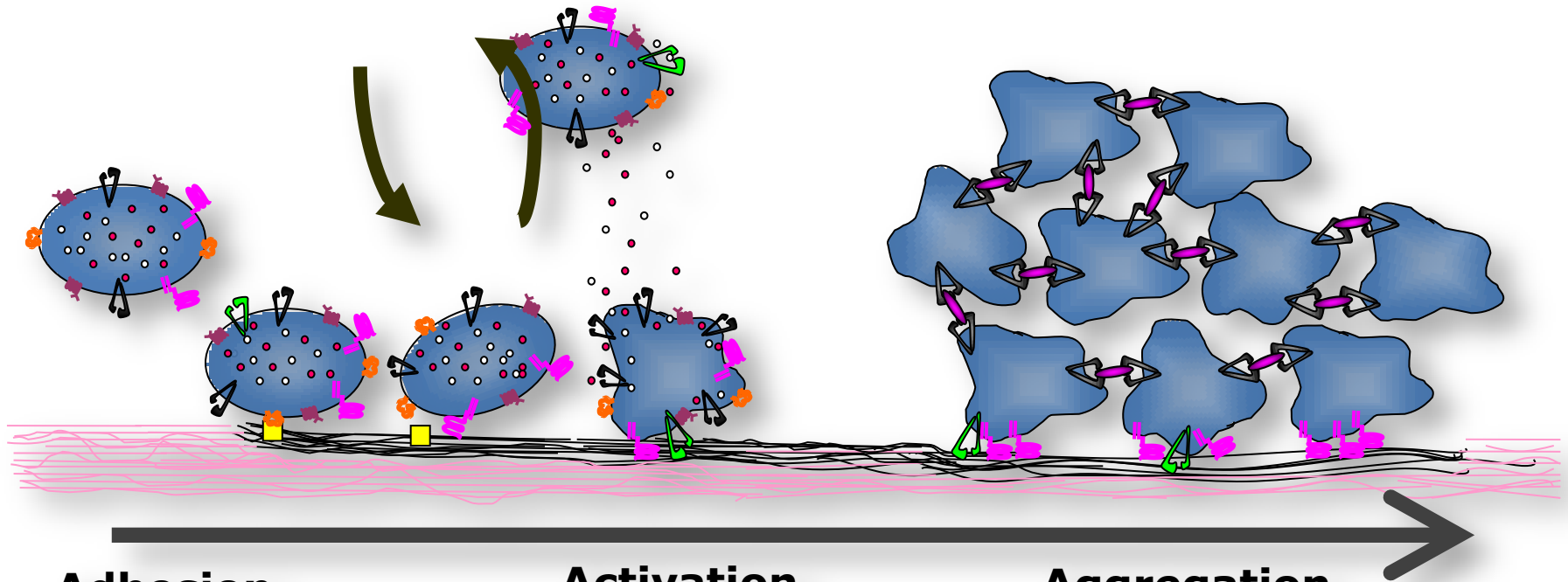
Chris Jones

ICMR

University of Reading

## Platelet function





**Adhesion**  
(transient tethering)

**Activation**  
(secretion)

**Aggregation**

 GPIb complex

  $\alpha$ -granule

  $\alpha$ IIb $\beta$ 3


 GPIIb/IIIa

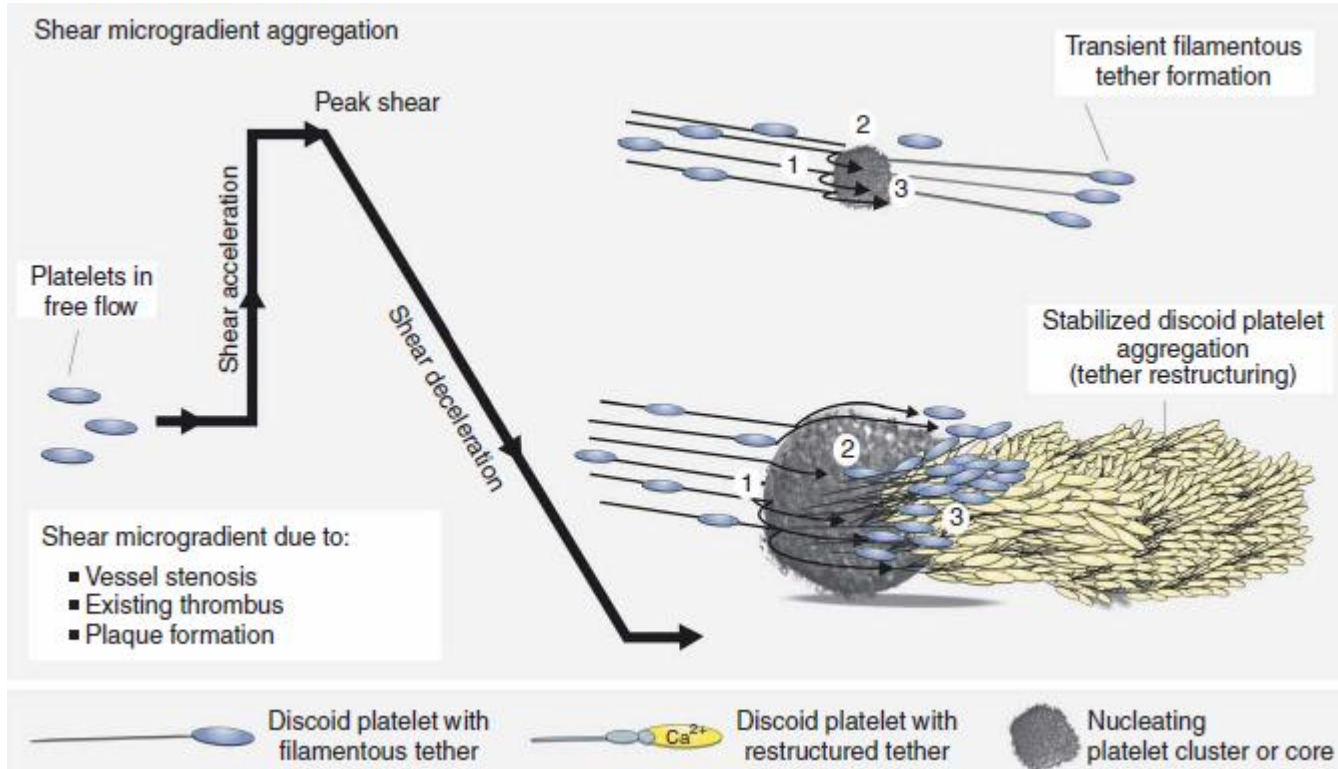
 Dense granule

 Fibrinogen

 GPCR

 VWF

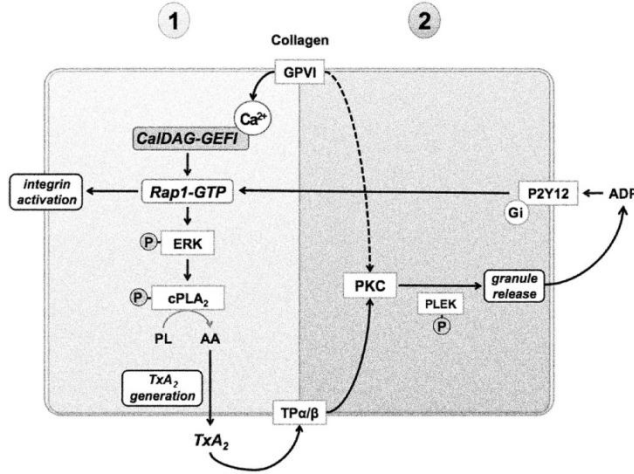
  $\alpha$ 2 $\beta$ 1



## A shear gradient–dependent platelet aggregation mechanism drives thrombus formation

Warwick S Nesbitt<sup>1,5</sup>, Erik Westein<sup>1,5</sup>, Francisco Javier Tovar-Lopez<sup>2</sup>, Elham Tolouei<sup>3</sup>, Arnan Mitchell<sup>2</sup>, Jia Fu<sup>1</sup>, Josie Carberry<sup>3</sup>, Andreas Fouras<sup>4</sup> & Shaun P Jackson<sup>1</sup>

**NATURE MEDICINE** VOLUME 15 | NUMBER 6 | JUNE 2009



## CalDAG-GEFI and protein kinase C represent alternative pathways leading to activation of integrin $\alpha$ IIb $\beta$ 3 in platelets

Stephen M. Cifuni, Denisa D. Wagner and Wolfgang Bergmeier  
Blood. 2008 Sep 1;112(5):1696-703.

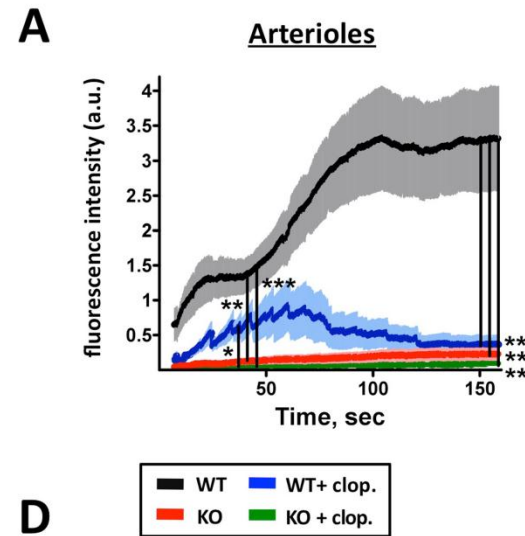
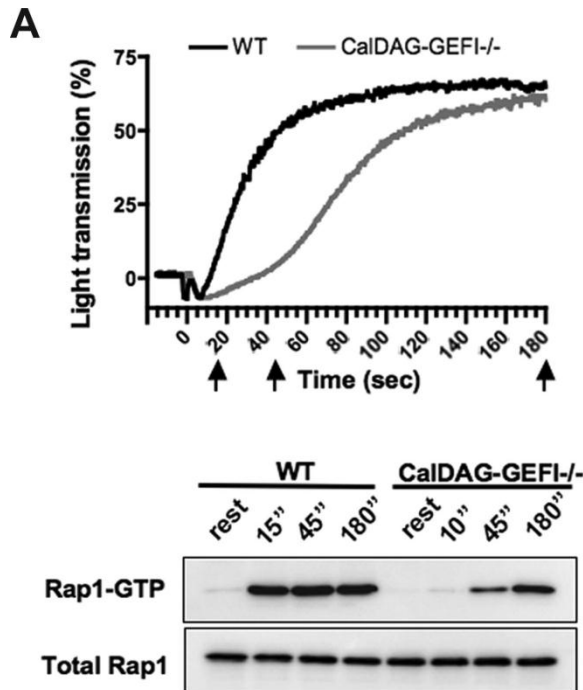
## CalDAG-GEFI is at the nexus of calcium-dependent platelet activation

Lucia Stefanini,<sup>1</sup> R. Claire Roden,<sup>1</sup> and Wolfgang Bergmeier<sup>1</sup>  
Blood 2009 114:2506-2514

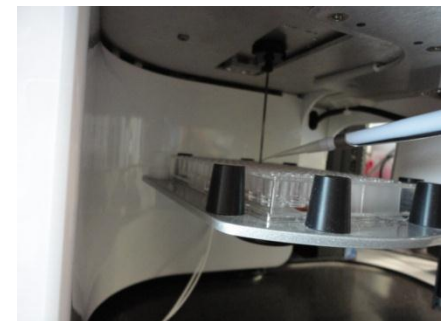
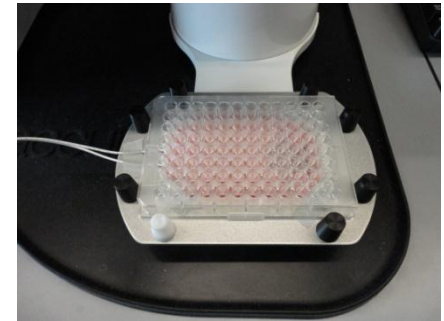
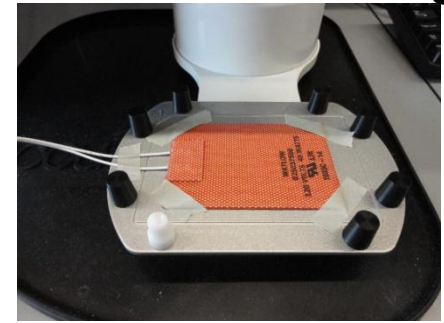
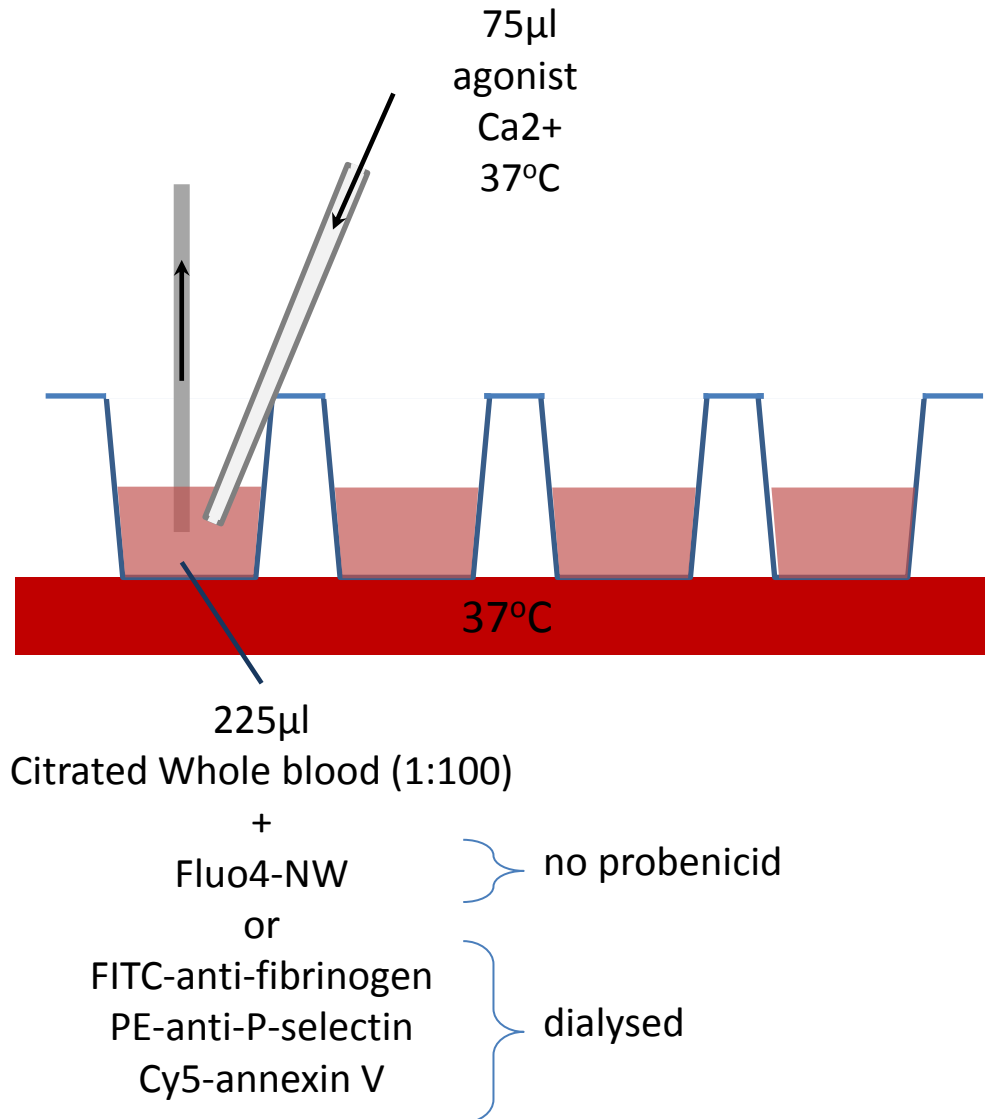
## The kinetics of $\alpha$ IIb $\beta$ 3 activation determines the size and stability of thrombi in mice: implications for antiplatelet therapy

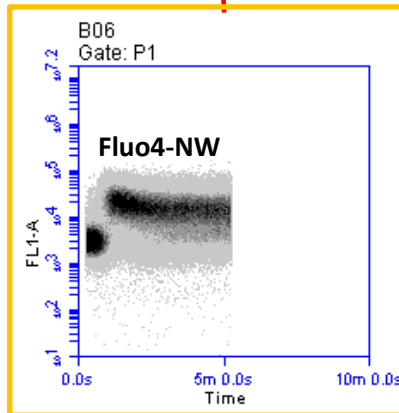
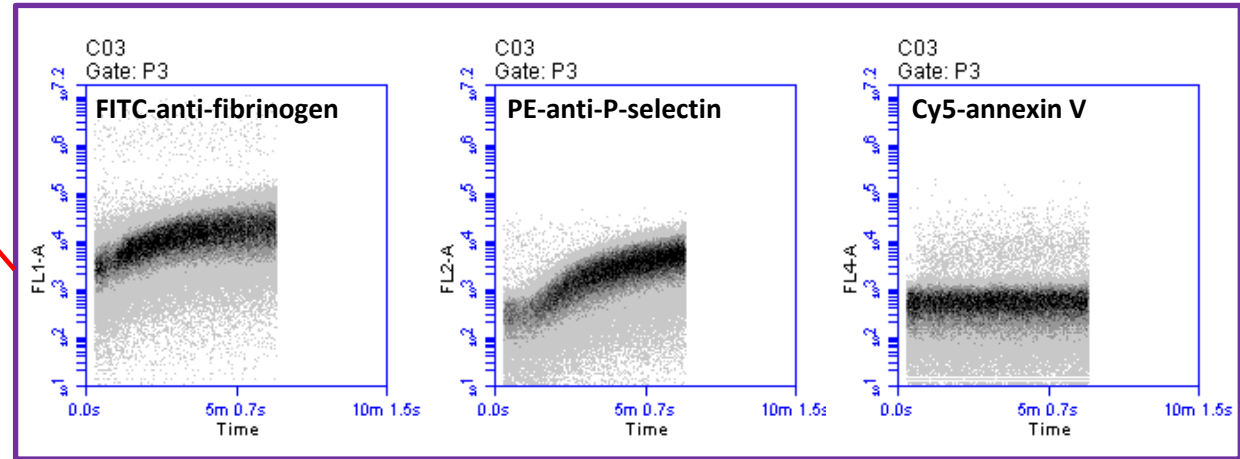
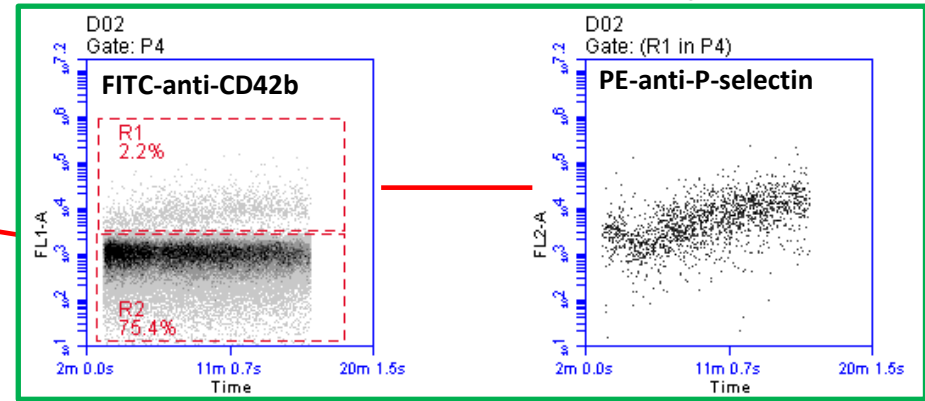
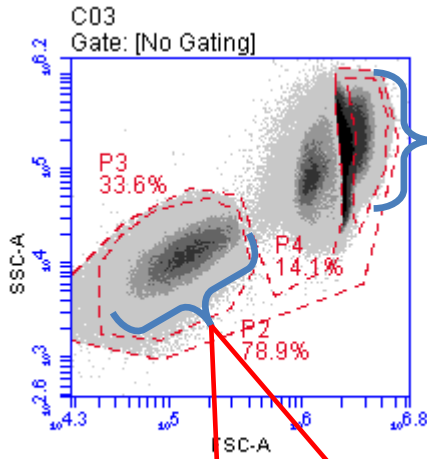
Moritz Stolla, Lucia Stefanini, R. Claire Roden, Massiel Chavez, Jessica Hirsch, Teshell Greene, Timothy D. Ouellette, Sean F. Maloney, Scott L. Diamond, Mortimer Poncz, Donna S. Woulfe and Wolfgang Bergmeier

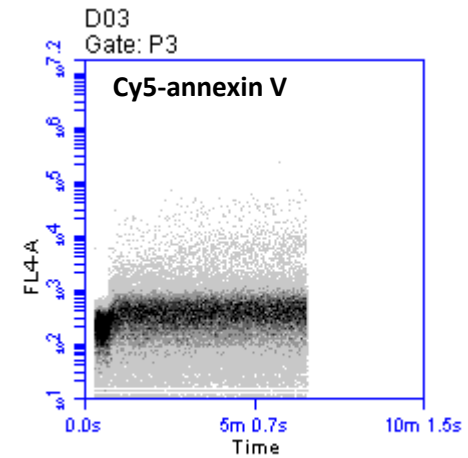
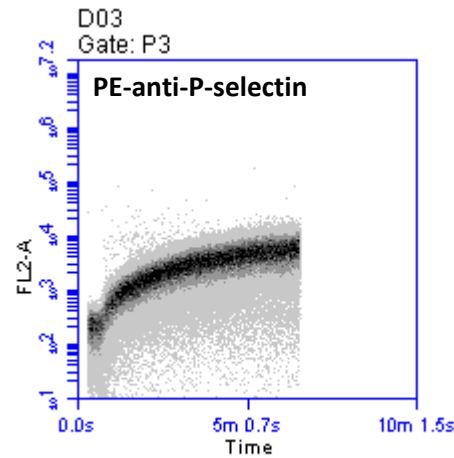
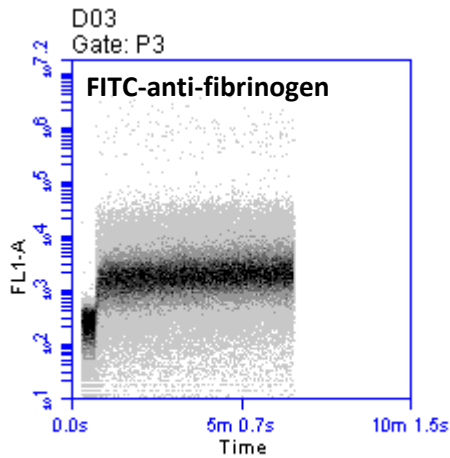
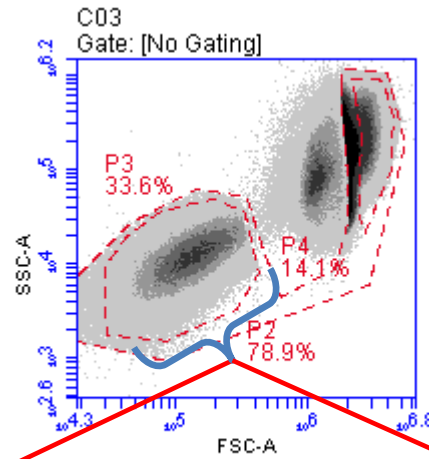
Blood. 2011 Jan 20;117(3):1005-13



# The assay

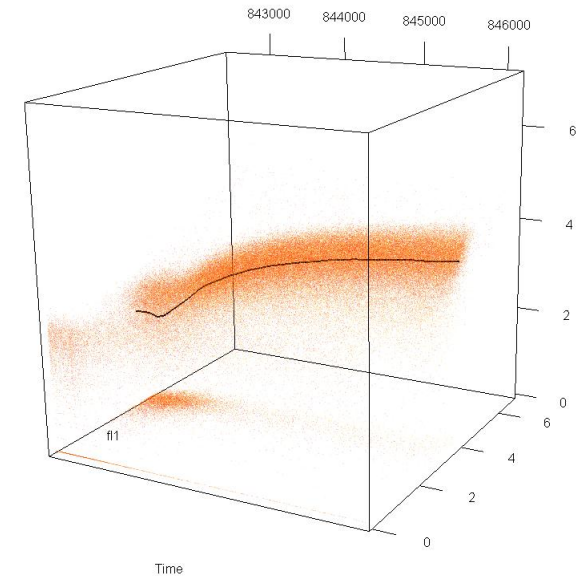
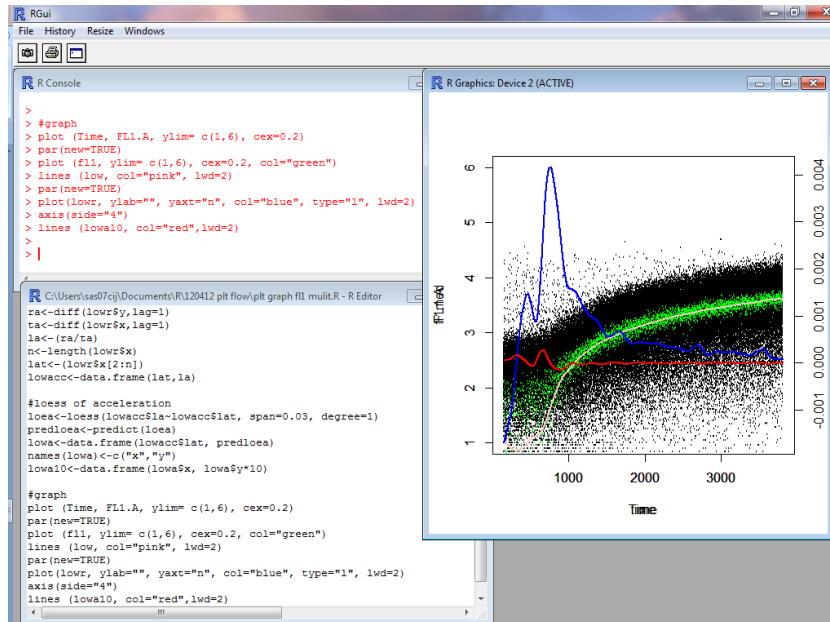
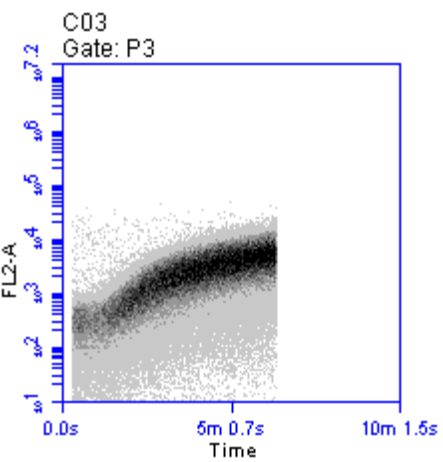




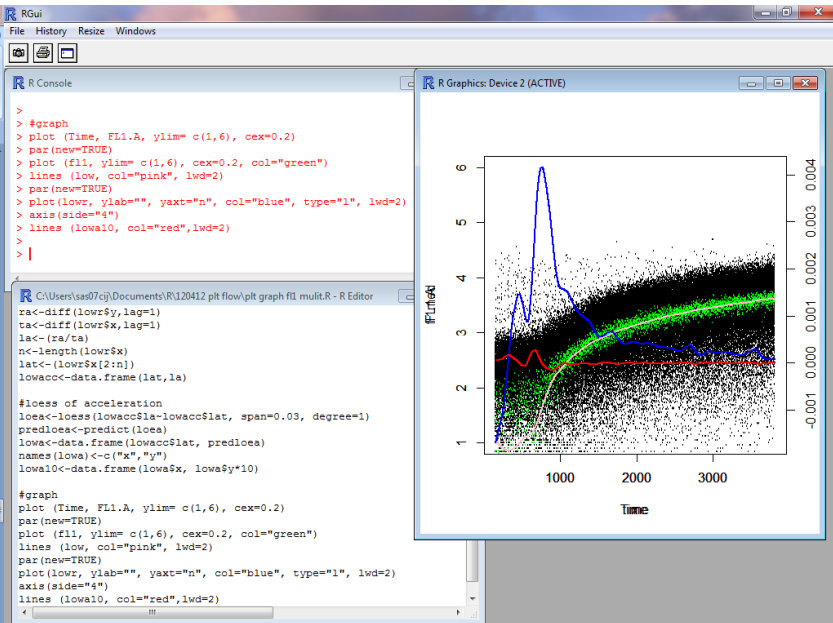




FSC-A	SSC-A	FL1-A	FL2-A	FL3-A	FL4-A	FSC-H	SSC-H	FL1-H	FL2-H	FL3-H	FL4-H	Width	Time
130381	21588	2439	379	386	58	262130	37432	2724	215	263	89	38	142
95998	6941	1046	338	105	307	200804	12846	1084	202	163	324	36	142
138125	7572	3538	472	80	235	259279	11645	3494	335	293	275	39	142
34556	8110	740	355	202	120	82669	16567	787	238	409	49	22	142
205857	14330	2784	397	437	153	396674	25294	3136	270	537	209	42	142
33899	2495	750	164	110	0	76473	5316	605	178	152	62	24	142
195063	13104	1419	340	293	319	392536	21766	1616	246	658	211	40	142
131735	8282	3198	504	394	0	269096	14464	3421	423	417	217	38	142
33757	2425	961	210	31	296	87982	4849	945	178	3	264	23	142
114836	7388	1307	173	0	152	244855	12029	1577	201	111	194	37	142
65290	4829	1608	263	0	0	146061	9801	1709	209	0	45	27	142
186328	22116	1646	138	0	0	388523	39867	1850	110	0	52	41	142
90824	13633	1400	138	110	65	198368	24023	1401	211	278	118	32	142
150220	6475	1388	607	380	11	320760	10668	1507	408	664	285	38	142



[www.r-project.org](http://www.r-project.org)

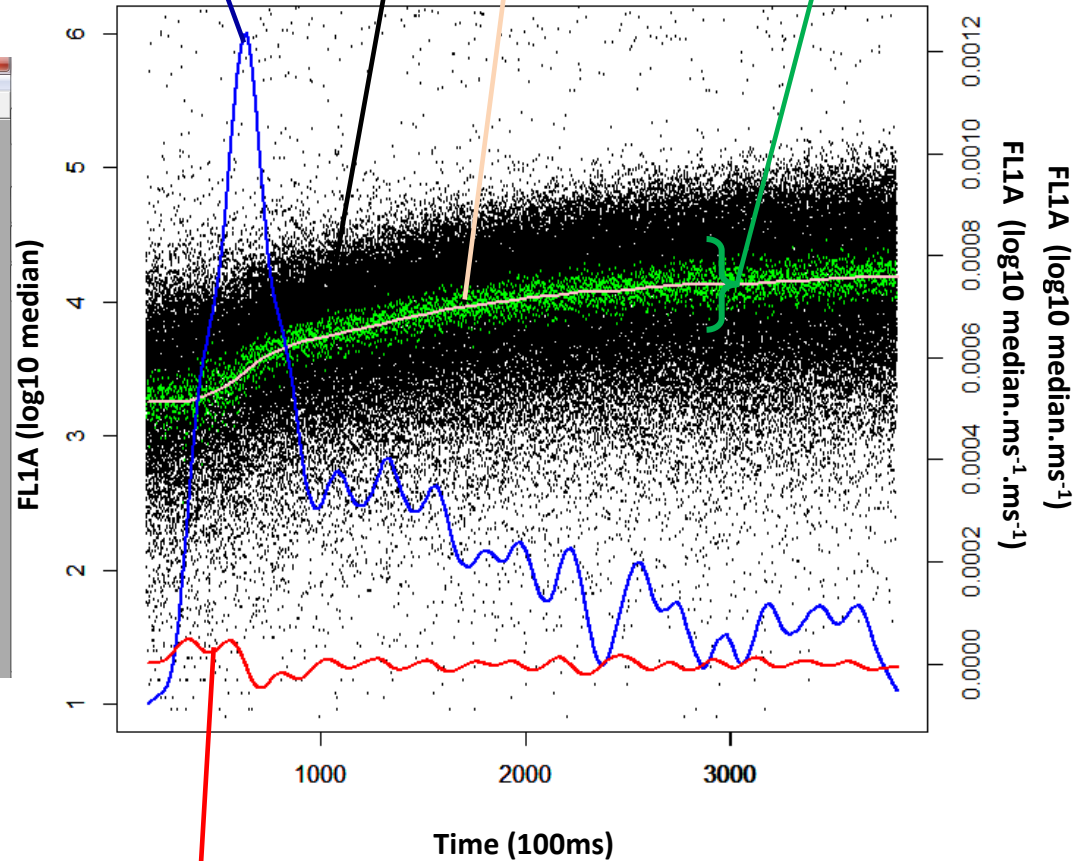


Rate of change of LOESS curve

Loess curve of median data

Raw data

Median fluorescence per 100ms



Acceleration of LOESS curve

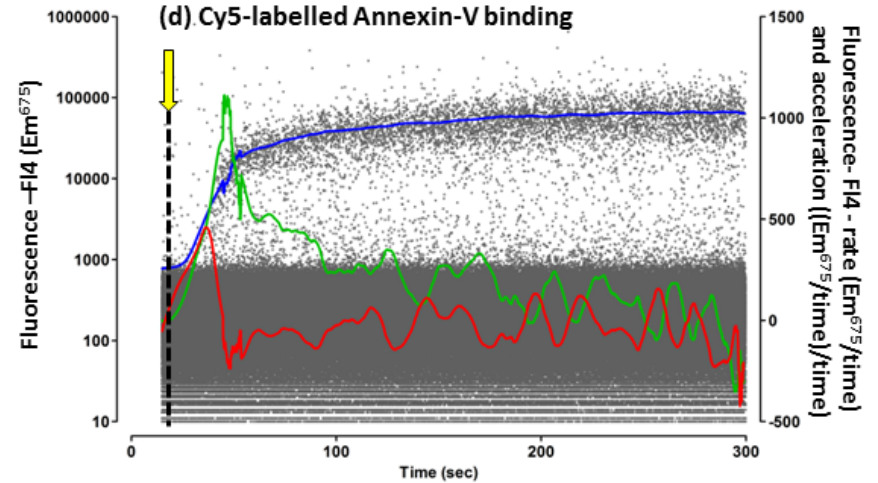
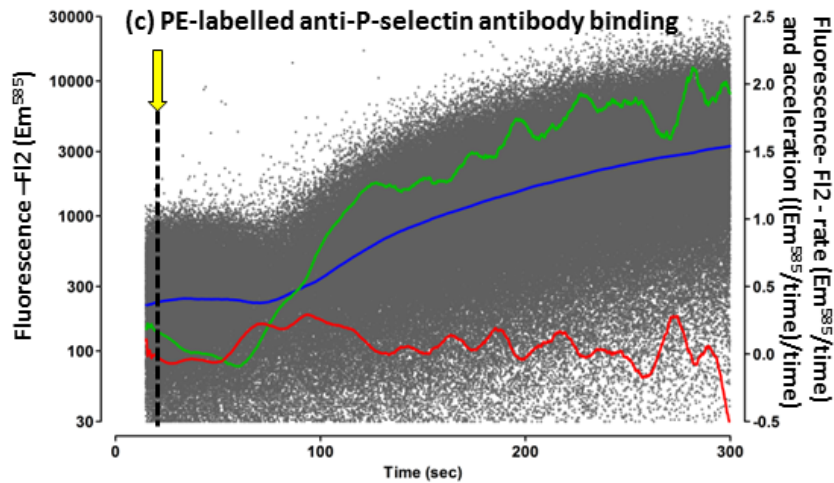
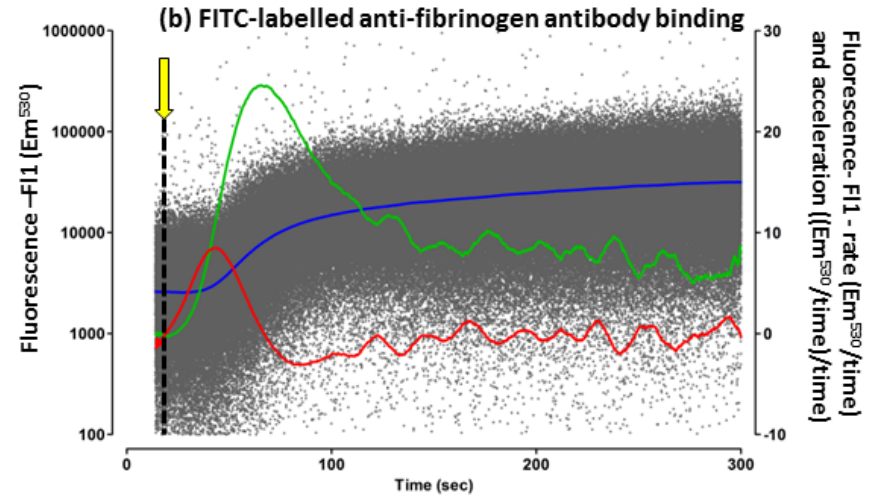
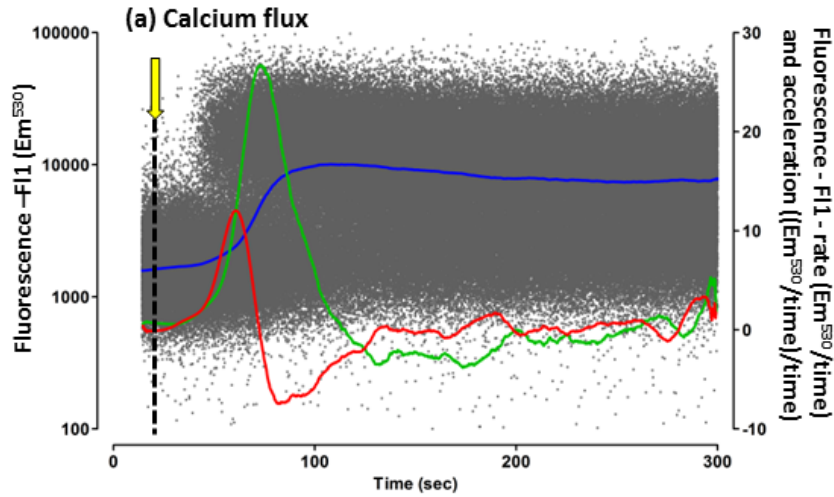
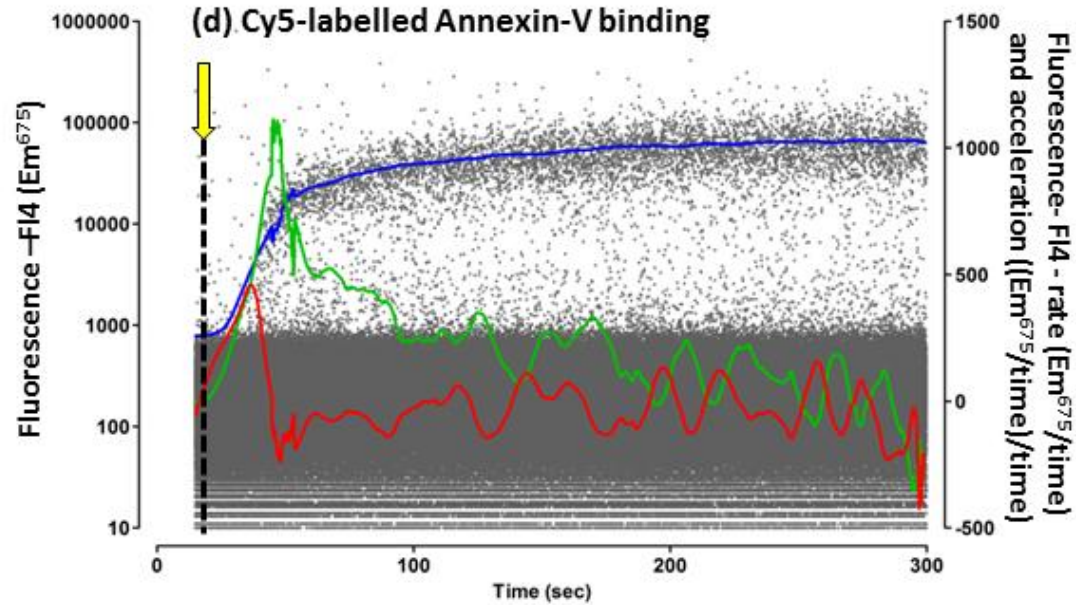


Table 1		CV
Ca <sup>2+</sup> flux	Max.	0.03
	Max. rate	0.03
	Max. acceleration	0.06
Fibrinogen binding	Max.	0.08
	Max. rate	0.06
	Max. acceleration	0.06
P-selectin exposure	Max.	0.08
	Max. rate	0.06
	Max. acceleration	0.16
Annexin-V binding	Max.	0.06
	Max. rate	0.07
	Max. acceleration	0.09

Table 1		Correlation with maximum	
		r <sup>2</sup>	p
Ca <sup>2+</sup> flux	Max.	.....	.....
	Max. rate	0.004	0.92
	Max. acceleration	0.594	0.13
Fibrinogen binding	Max.	.....	.....
	Max. rate	0.458	0.21
	Max. acceleration	0.260	0.38
P-selectin exposure	Max.	.....	.....
	Max. rate	0.001	0.97
	Max. acceleration	0.009	0.27
Annexin-V binding	Max.	.....	.....
	Max. rate	0.001	0.98
	Max. acceleration	0.434	0.02



**(e) Forward scatter, fibrinogen binding and P-selectin exposure on platelets binding Annexin-V**

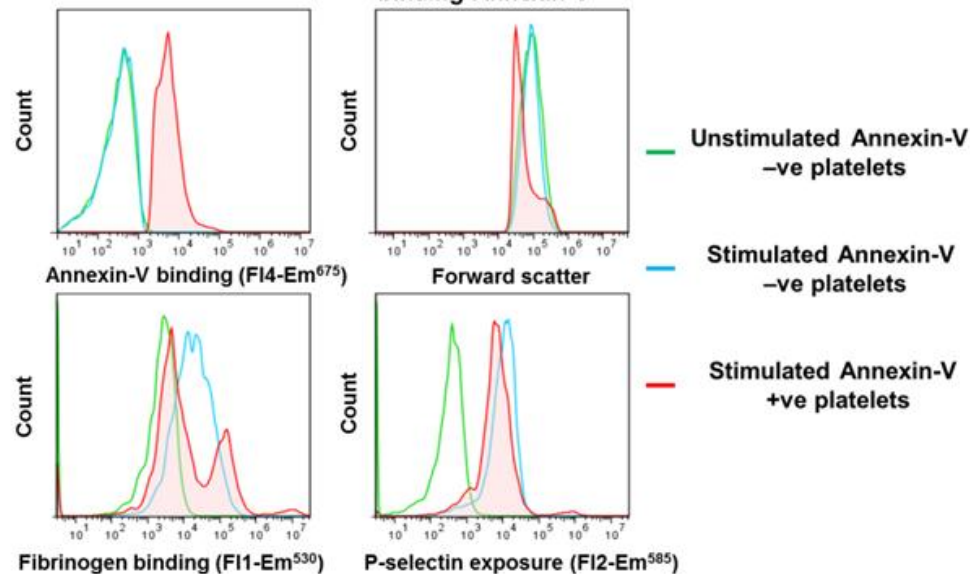
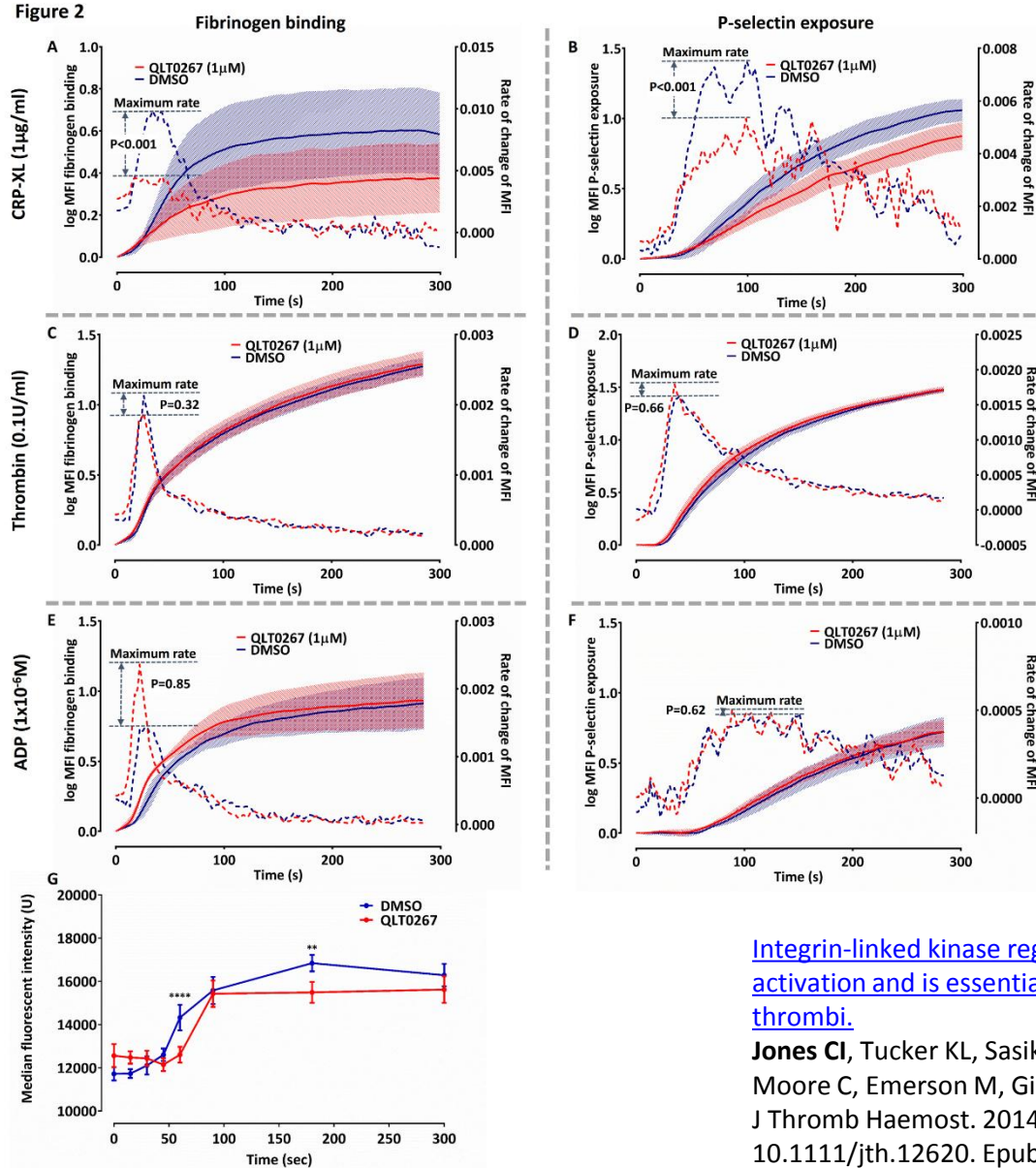


Figure 2



[Integrin-linked kinase regulates the rate of platelet activation and is essential for the formation of stable thrombi.](#)

Jones CI, Tucker KL, Sasikumar P, Sage T, Kaiser WJ, Moore C, Emerson M, Gibbins JM.

J Thromb Haemost. 2014 Aug;12(8):1342-52. doi: 10.1111/jth.12620. Epub 2014 Jul 31.

- Real time assay for the measurement the calcium flux or fibrinogen binding , P-selectin expression and Annexin V binding of populations of individual platelets
- Analysis methodology capable of efficiently extracting appropriate data for large studies or identifying interesting subpopulations



# Acknowledgements



**Prof. Jon Gibbins**



**Steve Garner**



This work was funded by the British Heart Foundation