



## European CytoFLEX Platform Meeting

Beckman Coulter Life Sciences invites you to join us at our two-day event for CytoFLEX users and those that would like to be one, to hear from leading scientists and experts about their research findings and insights as they share their technical expertise. Each day at the end of each session we will host our roundtable discussions where you can pose your questions to the speakers.

**VENUE: Park Plaza Utrecht, Westplein 50, 3531 BL Utrecht, The Netherlands**

The meeting is free of charge to attend but places are limited so please secure your seat without delay by registering.

We hope you will find our programme engaging and look forward to seeing you all in Utrecht.

### Day 1 Tuesday 14<sup>th</sup> June 2022

12.00	<b>Registration opens</b>
13.30	<b>Welcome Address</b> <i>Markus Kaymer and Alison Bell</i>
13.35	<b>Myeloid derived suppressors cells: new dampers in immune cell based therapy</b> <i>Nicola Tumino</i> <i>Ospedale Pediatrico Bambino Gesù, Rome, Italy</i>
14.15	<b>Comprehensive phenotyping of immune cell subsets from whole blood in people living with HIV</b> <i>Adriana Navas, MSc</i> <i>Radboud UMC, Nijmegen, The Netherlands</i>
14.55	<b>Refreshment break</b>
15.40	<b>To the power of 5! Five countries, Five LX's...</b> <i>Karen Hogg, University of York, UK</i>
16.20	<b>Roundtable Discussion</b>
17.10	<b>Closing comments Day 1</b> <b>Connecting the CytoFLEX !</b>
17.25	<i>Uwe Killemann - Particle</i>
17.40	<i>Maria Savino - Liquid Handling and Genomics</i>
17.55	<i>Lutz Ehrhardt - Centrifugation</i>
18.10	<b>Networking Event</b>
21.00	View the Virtual Reality experience for 'The CytoFLEX that Sorts' - the CytoFLEX SRT Cell Sorter.
	An opportunity to exchange information with your peers and other users of CytoFLEX.
	Chat with experts from Beckman Coulter Life Sciences.

### Day 2 Wednesday 15<sup>th</sup> June 2022

9.00	<b>CytoFLEX – from new optical technology to an entirely new product portfolio</b> <i>Mario Koksch</i> <i>Vice President and General Manager, Flow Cytometry Business Unit</i>
9.30	<b>Hematopoietic stem cell panel transfer from analyzer to CytoFLEX SRT cell sorter</b> <i>Juan Rubio-Lara</i> <i>University of York, UK</i>
10.10	<b>Refreshment Break</b>
10.40	<b>The CytoFLEX SRT Benchtop Cell Sorter: A Core Facility's Perspective</b> <i>Marjolijn Hameetman, BSc and</i> <i>IJsbrand Reyneveld, MSc</i> <i>Leiden University Medical Centre, Leiden, The Netherlands</i>
11.20	<b>Roundtable Discussion</b>
12.00	<b>Lunch</b>
13.00	<b>Microbiological 'Gold Standards' vs Flow Cytometry Method in Food and Beverage Industry</b> <i>Katarzyna Sielatycka and Joanna Śliwa-Dominiak</i> <i>Sanprobi Sp. z.o.o. Sp. K, Szczecin, Poland</i>
13.50	<b>New Generation Flow-Cytometry to Measure Circulating Microvesicles In Cancer</b> <i>Claudia Maria Radu</i> <i>University of Padua, Padua, Italy</i>
14.30	<b>Roundtable Discussion</b>
15.00	<b>Closing Comments</b>