



sartorius

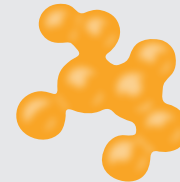
Research Xchange Forum 2018

Insights | Translation | Solutions

Trends and Challenges in Regenerative Medicine & Cell Therapy



March 20 – 21, 2018
Sartorius Goettingen,
Germany



The Sartorius & Science
Prize for Regenerative
Medicine & Cell Therapy

turning science **into solutions**



What to expect

Learn about the latest Insights in regenerative medicine and cell therapy and exchange ideas as to their Translation into viable Solutions.

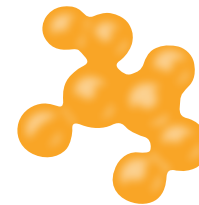
Our Research Xchange Forum brings together academic and industry experts to foster exchange and to stimulate dialog about the next generation of interdisciplinary research. Establish new contacts and explore the potential for collaborations with like-minded companies and academic professionals to gain a better understanding of their processes, workflows and challenges.

Deepen your discussions during a joint gala dinner and be there when the winner and the 3 finalists of "The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy" receive their awards.
Winner: U.S. \$25,000 + publication in *Science* print and online;
3 finalists: U.S. \$5,000 per winner + publication in *Science* online.

We look forward to seeing you at our Research Xchange Forum.

Highlights

- Keynote lectures by the winner and lectures by the 3 finalists of "The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy", including a speech by *Science* editor Dr. Priscilla N. Kelly
- Keynote lectures on cellular immunotherapy by industry experts
- Inspirational speech by the Nobel Prize laureate of 2014 in chemistry Prof. Dr. Stefan W. Hell
- Panel discussion with stakeholders from industry and academia including *Science* editor Dr. Priscilla N. Kelly
- Gala dinner and award ceremony for the exceptional scientists receiving "The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy"
- Thought-provoking dialog and networking opportunities with academic and industry experts
- Exclusive inside view into the work processes of Sartorius during an optional plant tour



The Sartorius & *Science*
Prize for Regenerative
Medicine & Cell Therapy

Awarded by



sartorius

Science



Speaker line-up*

Guest Contributors



Dr. Priscilla N. Kelly,
Washington DC, USA

AAAS/*Science*
Editor

Topic: "How to publish in *Science*"



Prof. Dr. Stefan W. Hell,
Goettingen, Germany

Max Planck Institute for Biophysical
Chemistry

Managing Director and Nobel Prize laureate
of 2014 in chemistry

Topic: "How to win a Nobel Prize"

© Peter Badge/
Typos1 in cop. with
Foundation Lindau
Nobel Laureate
Meetings – all
rights reserved
2017.

Academic Speakers

Keynote lecture by the winner and lectures by the 3 finalists of
"The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy"



Prof. Dr. Stefan Wölfl
Heidelberg, Germany

University of Heidelberg, Institute for
Pharmacy and molecular Biotechnology

Topic: Small molecules to enable, facilitate
and improve iPSC generation and cell
reprogramming

Overview: The generation of patient
specific specialized cell-types is a major
task in regenerative medicine. Despite the
success of genetic reprogramming protocols,
it is still a challenge to generate, cultivate
and expand patient specific induced plurip-
otent stem cells (iPSCs). The generation of
iPSCs requires the expression of specific

transcription factors, expression of which
is maintained to conserve the pluripotent
state. For reprogramming these factors are
inhibited and tissue specific sets of gene
expression are induced. We focus on the
development of chemical compounds to
replace genetic reprogramming and e.g.
have identified Oct4 inducing small mole-
cules. Although we cannot yet fully replace
genetic reprogramming factors, our small
molecules improve reprogramming
efficiency, overcome reprogramming
resistance and improve cultivation of iPSCs
for reprogramming.

Speaker line-up*

Industry Speakers



Dr. Ulf Geumann,
Munich, Germany

apceth Biopharma GmbH, Group Leader
Preclinical Development

Topic: Mesenchymal stem cells in
immunotherapy

Overview: apceth is harnessing the potential of mesenchymal stem cells (MSC), in which the mechanism of therapeutic effects are very different from that of CAR-T cells. The aim is to provide innovative cell-based gene therapies for the treatment of major indications such as cancer, chronic lung and kidney diseases and for immunomodulation.

apceth has created its allogenic therapeutic platform based entirely on the genetic engineering of MSC isolated from healthy donors.



Francesca Bellintani,
Milan, Italy

MolMed, Vector Downstream Process
Development Manager

Topic: Retroviral and lentiviral production:
development and scaling up of upstream
and downstream process steps

Overview: MolMed has implemented a large-scale process to meet the increasing demand for clinical and commercial grade vector production. There is a special focus on identification of challenges in the downstream | purification strategy to ensure efficient contaminant removal and preservation of viral infectivity.



Dr. Thierry Wurch
Suresnes, France

Servier, Director of External Studies in
Immuno-oncology

Topic: Allogeneic approach developed by
Servier and the preclinical data of UCART19

Overview: Servier is applying the Collectis' TALEN gene editing technology to overcome current limitations and provide an allogeneic, frozen, "off-the-shelf" T cell based medicinal product. UCART19 is an allogeneic CAR T-cell product candidate being developed for treatment of CD19-expressing hematological malignancies. It is currently being evaluated in a phase 1 clinical trials in relapsed | refractory acute lymphoblastic leukemia (ALL) of adult and pediatric patients. The selection of UCART19 as a candidate was based on the impressive clinical benefits leading to the marketing of the two first autologous CAR T-cell products Kymriah (Novartis) and Yescarta (Kite Gilead).



Dr. Nina G. Bauer
Basel, Switzerland

Lonza Assoc. Director of Commercial
Development, Autologous Cell Therapies

Topic: Manufacturing strategies for
patient scale cell therapies

Overview: The "Future Technologies" business area under the Pharma & Biotech division of Lonza has a focus on personalized medicine & cell therapy. One of the major challenges in these fields is the production of the patient-specific and patient-scale CAR-T cells or stem cells for treatment. Lonza has partnered with Octane on the Octane Cocoon™ platform, which offers a ground-breaking innovation to address this challenge. Unlike current, labor-intensive production methods, this platform has the potential to enable the automated manufacturing of patient-specific cell therapies for maximum efficiency, traceability and scale-up.

Agenda

Tuesday, March 20

Chair: Ruth McDermott

09:40	Transfer to Sartorius College (from hotel)	
10:00 – 11:00	Registration	
11:00 – 11:30	Welcome & Introduction	Gerry MacKay, Sartorius, Germany Member of the Group Executive Committee and Executive Vice President LPS Marketing Sales Services
11:30 – 12:00	Inspirational speech: How to win a Nobel Prize	Prof. Dr. Stefan Hell, Max Planck Institute for Biophysical Chemistry, Germany Managing Director and Nobel Prize laureate of 2014 in chemistry
12:00 – 12:30	How to publish in Science	Dr. Priscilla N. Kelly, AAAS Science, USA Editor
12:30 – 13:15	Lunch break Application Session	
13:15 – 14:00	Keynote presentation	Sartorius & Science Prize Winner
14:00 – 14:30	Lecture	Sartorius & Science Prize Finalist 1
14:30 – 15:00	Coffee break	
15:00 – 15:30	Lecture	Sartorius & Science Prize Finalist 2
15:30 – 16:00	Lecture	Sartorius & Science Prize Finalist 3

16:00 – 17:00	Panel Discussion In search of excellence –Scientific publications in the 21 st century	Moderation: Dr. Reinhard Baumfalk, Sartorius, Germany Vice President R&D Instrumentation & Control Participants: Prof. Dr. Stefan Hell, Dr. Priscilla N. Kelly, Dr. Ulf Geumann
17:00 – 17:15	Sum up & Closing	
17:20	Transfer to hotel (from Sartorius College)	
18:15	Transfer to award ceremony (from hotel)	
18:30 – 23:30	Award ceremony	
23:30	Transfer to hotel	

Agenda

Wednesday, March 21

Chair: Del Trezise

08:15	Transfer to Sartorius College Sartorius Trainingscenter (from hotel)	
08:30 – 10:00	Optional Plant Tour (Sartorius College) or Upscaling Lab Demonstration Tour (Sartorius Trainingscenter)	
09:40	Transfer to Sartorius College (from Sartorius Trainingscenter)	
10:00 – 10:15	Opening Day 2	
10:15 – 11:00	Manufacturing strategies for patient scale cell therapies	Dr. Nina G. Bauer, Lonza, Switzerland Assoc. Director of Commercial Development, Autologous Cell Therapies
11:00 – 11:30	Allogenic approach development by Servier and the preclinical data of UCART19	Dr. Thierry Wurch, Servier, France Director of External Studies in Immuno-oncology
11:30 – 12:30	Lunch break Application & Poster Session	
12:30 – 13:00	Small molecules to enable, facilitate and improve iPSC generation and cell reprogramming	Prof. Dr. Stefan Wöfl University Heidelberg, Germany Institute for Pharmacy and molecular Biotechnology
13:00 – 13:30	Mesenchymal stem cells in immunotherapy	Dr. Ulf Geumann, apceth Biopharma GmbH, Germany Group Leader Preclinical Development
13:30 – 14:00	Retroviral and lentiviral production: development and scaling up of upstream and downstream process steps	Dr. Francesca Bellintani, MolMed, Italy Vector Downstream Process Development Manager
14:00 – 14:10	Sum up & Closing	

Registration

March 20 – 21, 2018

To register and to obtain further information, please visit our website:
promotions.sartorius.com/rxf2018

The registration fee includes admission, gala dinner, all coffee breaks, lunches and the optional plant tour. Advance registration is required.

We offer an early bird discount until January 15, 2018

	Industry Business	Academia	PhD Students Students
Early Bird	395 EUR	295 EUR	105 EUR
Regular	595 EUR	495 EUR	195 EUR
Group Ticket	Pay the 1st ticket and get 50% off for every additional registration		

Venue:

Sartorius College
Otto Brenner Strasse 20
37079 Goettingen, Germany
www.sartorius.com

Your Contact:

Romi Jeromis
Sartorius Lab Instruments GmbH & Co. KG
Phone: +49.551.308.2659
forum@Sartorius.com

If you have any questions,
just contact us at any time – we would
be happy to help you.

Please note that photos and videos will be
taken during the event. If you do not wish
to be photographed, please let us know
upon registration.

Follow us on Twitter #rxf2018

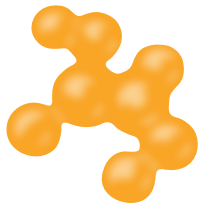


Research Xchange Forum 2018

Get insights into the latest developments from internationally renowned academic and industry speakers.

Special Section:

Lectures from "The Sartorius & *Science* Prize for Regenerative Medicine & Cell Therapy" Winner and Finalists.



The Sartorius & *Science*
Prize for Regenerative
Medicine & Cell Therapy

Awarded by



sartorius

Science

March 20 – 21, 2018
Register now!

To register and to obtain further information, please visit our website:

promotions.sartorius.com/rxf2018



Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen