



sartorius stedim
biotech

Upstream & Downstream Technology Forum 2018
June 5 – 7, 2018
Sartorius College, Goettingen, Germany



turning science into solutions

Invitation

Upstream and Downstream Processing: One Common Goal

Successful process optimization in upstream has led to increased cell densities and product titers. But what does this mean for downstream processing and purification of the final drug product? These modern, intensified production processes are pushing the bottleneck more and more towards downstream. The need to address this has become stronger than ever before.

Following the success of 2016 combined event also this year we bring together our established European Upstream Technology Forum and European Downstream Technology Forum to one joint event. Gain knowledge and share experience with experts from the industry on new trends and bottlenecks for both upstream and downstream. Take a look at the agenda and join us for this exciting event.

You will find up-to-date information posted on our website: www.sartorius.com/upstream-downstream-forum-2018

Broaden your process understanding and update your knowledge in upstream and downstream processing while meeting experts from the industry.

Agenda

Chairs

Jörg Weyand, Director Marketing Fermentation Technologies, Europe & LA
Dr. Grant Cameron, Director Marketing Purification Technologies, Europe
Sartorius Stedim Biotech, Germany

Tuesday, June 5, Upstream

11:30 – 12:00

Registration

12:00 – 13:00

Lunch | Exhibition

13:00 – 13:15

Welcome

Dr. Oscar-Werner Reif,
Chief Technology Officer,
Sartorius Stedim Biotech, Germany

13:15 – 14:00

Continuous and intensified processing: Current state of the art and future outlook

Dr. Gerben Zijlstra,
Platform Marketing Manager
for Continuous BioManufacturing,
Sartorius Stedim Biotech,
Netherlands

14:00 – 14:45

Joint development project of ambr[®] 250 ht perfusion

Heather Nunn,
Scientist, Pivotal Cell Sciences
and Technologies, Amgen, USA

14:45 – 15:30

Coffee Break | Exhibition

15:30 – 16:00

Evaluation of body feed harvest for cell culture fluids with high host cell protein and DNA load

Dr.-Ing. Detlef Eisenkrätzer,
Director Fermentation
Development Roche Pharma,
Roche Diagnostics GmbH, Germany

16:00 – 16:30

The next frontier in biomanufacturing: Novel therapeutic modalities and 'old challenges' in the 21st century

Dr. Jesús Zurdo, Senior Director
Strategic Innovation,
Lonza Pharma & Biotech, UK

16:30 – 17:00

Q&A | Discussion

17:15

Transfer to Welcome Dinner

21:00

Transfer back to Hotel

Agenda

Wednesday, June 6, Upstream

08:30 – 09:00
Coffee | Exhibition

09:00 – 09:30

Scale-up of lentiviral vectors for use in gene and cell-based therapies
Dr. Carol Knevelman,
Head of the Process R&D department, Oxford Biomedica, UK

09:30 – 10:00

Process development goes "viral"
Dr. Eugenio Macchiarulo,
Vector Development Associate Director, Adaptimmune Ltd., UK

10:00 – 10:45
Coffee Break | Exhibition

10:45 – 11:15

They are what they eat –
A compendium about the power of cell culture media
Dr. Aziz Cayli,
Founder and CEO of Florabio, Florabio AS, Turkey

11:15 – 11:45

Evaluation of single-use bioreactors to optimize cell growth and scale-up of viral veterinary vaccine production
Dr. Lidia Garcia,
Principal Scientist, Zoetis Manufacturing Research S.L., Spain

11:45 – 12:00

Q&A | Discussion

12:00 – 13:00
Lunch | Exhibition

► **Joint Session:**
Where Upstream Meets Downstream

13:00 – 13:30

USP & DSP challenges and technologies for a production of a freeze-dried vaccine
Jordi Ruano Bou,
Production Director of HIPRA S.A., Spain

13:30 – 14:00

The role of data analytics in improving biomanufacturing processes
Robert Zuban,
Senior Data Scientist, Sartorius Stedim Data Analytics AB, Sweden

14:00 – 14:30
Coffee Break | Exhibition

14:30 – 15:00

Strategic decision making progress for the sizing and implementation of a new single-use 2000 L biosimilar plant
Luisa Amoedo,
Site Director, mAbxience, Spain

15:00 – 15:30

Implementation of membrane chromatography for adenovirus downstream process intensification
Dr. Helge Abrecht,
Expert Scientist, GSK Vaccines, Belgium

15:30 – 16:00

Q&A | Discussion

16:15
Transfer to Hotel

18:15
Transfer to Networking Dinner

21:45
Transfer back to Hotel

Agenda

Thursday, June 7, Downstream

08:30 – 09:00
Coffee | Exhibition

09:00 – 09:30

ambr[®] crossflow: First impressions of a novel high throughput process development system

Dan Gibson,
Biopharm Process Research,
GSK, UK

09:30 – 10:00

Effective viral clearance – more than high reduction factors

Dr. Horst Ruppach,
Scientific Director Viral Safety,
Charles River, Germany

10:00 – 10:45
Coffee Break | Exhibition

10:45 – 11:15

Virus removal capacity in the downstream process of new format antibodies

Dr. Moritz Bennecke,
Downstream Processing and
Recovery Manager,
Roche Diagnostics GmbH, Germany

11:15 – 11:45

Implementation and validation of a single-use mixing system for virus inactivation with solvent | detergent

David Balbuena,
Head of Production,
LFB Biomanufacturing, France

11:45 – 12:15

Scale down capability and benefits for tangential flow filtration

William Man, Senior Supervisor,
Genentech, USA

12:15 – 13:00

Q&A | Discussion
Meeting Closure

13:00 – 14:00
Lunch | Exhibition
End of Conference

We look forward to welcoming you in Goettingen, Germany.

Conference Address

Sartorius College
Otto-Brenner-Strasse 20
37079 Goettingen, Germany
Phone +49.551.308.2101

Welcome Dinner | Tuesday

Restaurant Bullerjahn

Markt 9
37073 Goettingen
Phone: +49.551.3070100

Networking Dinner | Wednesday

Restaurant Graf Isang

Seestrasse 37
37136 Seeburg
Phone: +49.5551.919880

We encourage you to forward this invitation to interested colleagues.



Sartorius College, Goettingen

Venue

Sartorius College
Otto-Brenner-Strasse 20
37079 Goettingen, Germany
www.sartorius.com
Phone +49.551.308.2101

Transportation

The nearest airport is Hanover, about 1½ hours by train or car to Goettingen. The train station is 10 min from the College by taxi.

Please find a detailed map on the next page.

Directions

On motorway A7 (Hannover-Kassel) take the exit for Goettingen Nord (72). Take first exit for Industriegebiet, take first exit in roundabout. Then at the next light, take a left. The College will appear on the left-hand side.

Your Contacts

Antje Schneider
Sartorius Stedim Biotech,
Goettingen
Phone +49.551.308.1397
antje.schneider@sartorius-stedim.com

Vanessa Henze
Sartorius Stedim Biotech,
Goettingen
Phone +49.551.308.2032
vanessa.henze@sartorius-stedim.com

Nadine Engel
Sartorius Stedim Biotech,
Goettingen
Phone +49.551.308.4129
nadine.engel@sartorius-stedim.com

Registration

Early bird rate (until April 27):

€ 395 + 19 % VAT
€ 150 + 19 % VAT for academia

Standard price (from April 28):

€ 550 + 19 % VAT
€ 170 + 19 % VAT for academia

This includes the welcome dinner, networking dinner, all breaks and lunches.

To register, please either
– register on the web:
www.sartorius.com/upstream-downstream-forum-2018
– call or e-mail listed contacts

Accommodation

We have arranged limited allotments in the following hotels.

Please mention "Upstream & Downstream Technology Forum 2018".

Hotel Rennschuh

Kasseler Landstrasse 93
37081 Goettingen
Phone: +49.551.9009.0

www.rennschuh.de
€ 55 incl. breakfast

Novostar Hotel Goettingen

Kasseler Landstrasse 25d
37081 Goettingen
Phone +49.551.99770

www.novostar.de
€ 57 – 77 incl. breakfast

Park Inn by Radisson Goettingen

Kasseler Landstrasse 25c
37081 Goettingen
Phone +49.551.270707.520

www.parkinn-goettingen.de
€ 78 – 113 incl. breakfast

Further hotel information:
www.sartorius.com/
upstream-downstream-forum-2018

1 Sartorius Campus Headquarters
Otto-Brenner-Strasse 20
37079 Goettingen

2 Sartorius Stedim Plastics GmbH
Karl-Arnold-Strasse 21
37079 Goettingen

3 Sartorius Application Center
Marie-Curie-Strasse 7
37079 Goettingen

4 Sartorius Service Center CE (Distribo)
Groner Siekanger 1
37081 Goettingen

5 Sartorius Off-Campus Site
Telekom Building
Philipp-Reis-Strasse 2a
37075 Goettingen

P Parking Building
Access via Willi-Eichler-Strasse 28
Parking for visitors and employees

Where to stay
1. Hotel Freizeit In
2. Hotel Rennschuh
3. Novostar Hotel
4. Park Inn by Radisson

Contact:
Phone +49.551.308.0
E-Mail
empfang.campus@sartorius.com

For more information about Sartorius, please visit
www.sartorius.com

Shuttle Service

We arrange a shuttle service from the Sartorius College to the hotels above and vice versa. You will find detailed information in your conference information and at the hotel receptions.

We advise you that photographs | videos will be taken at the event which may be published later. In case you do not consent to such publication, we ask you to inform us.