



sartorius stedim
biotech

European Upstream & Downstream Technology Forum 2016 June 07 – 09, 2016

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

Agenda

Chairs

Dr. Christel Fenge, Davy De Wilde Fermentation Technologies,
Dr. Fritjof Linz Purification Technologies
Sartorius Stedim Biotech, Germany

Tuesday, June 07, Upstream

11:30 – 12:00

Registration

12:00 – 13:00

Lunch | Exhibition

13:00 – 13:15

Welcome

Dr. Oscar-Werner Reif,
Executive Member of the GEC,
Research and Development,
Bioprocess Solutions, Sartorius
Stedim Biotech, Germany

► **Smart upstream
development & scale up**

13:15 – 13:45

Advanced process development using automated micro bioreactors shortens timelines and provides process solutions for a rapid scale up

Dr. Jincal Li,
Executive Director, Biologics
Process Development, Wuxi AppTec
Biopharmaceuticals Co., Ltd., China

13:45 – 14:15

SAM – Sedimentation in an automated microbioreactor: Using the ambr[®] 15 as a scale down tool for perfusion cultures

Steffen Kreye,
USP Development Scientist,
Glycotope, Germany

14:15 – 14:45

ambr[®] 250: from process screening up to production scale in cell culture development

Mira Drolshagen,
USD project engineer, Boehringer
Ingelheim, Germany

14:45 – 15:15

Coffee Break | Exhibition

15:15 – 15:45

Continuous processing in biopharmaceutical manufacturing

Dr. Tina Lütke-Eversloh,
Senior process manager,
Rentschler Biotechnologie,
Germany

15:45 – 16:15

Overcoming the challenge of data deluge from high throughput bioreactors. Gaining insight from gigabyte datasets using an informatics approach.

Doug Marsh,
Senior scientist, Biopharm
Process Research, GSK Pharma, UK

16:15 – 16:45

Media optimization for enhanced processes

Michael Gillmeister,
Senior scientist, Lonza Biologics, US

16:45 – 17:00

Q&A | Discussion

17:15

End of Day

17:30

Transfer to Welcome Dinner

Agenda

Wednesday, June 08, Upstream

► Optimizing large scale manufacturing

08:30 – 09:00
Coffee | Exhibition

09:00 – 09:30

Technical & process considerations when establishing a 2000 L production SU facility

Roel de Waard,
Process Engineer Upstream Process Development, Synthron Biopharmaceuticals BV, Netherlands

09:30 – 10:00

Virus production utilizing Single Use Bioreactors

Jordi Ruano,
Biologics Production Director, Hipra, Spain

10:00 – 10:30
Coffee Break | Exhibition

10:30 – 11:00

SDD – Standard Disposables Design™: Application to Media and Buffer preparation

John Levesley,
Senior Process Engineer PM Group, Ireland

11:00 – 11:30

MFCS/win based automation for XD high cell density cultivation processes

Jurjen de Jong,
Bioprocess scientist, Patheon Biologics, Netherlands

11:30 – 12:00

Scale-up in the single use age: Does geometry matter?

Colin Jaques,
Senior Principal Scientist, Lonza Biologics, UK

12:00 – 13:00

Lunch | Exhibition

► **Joint Session:**
Where upstream meets downstream

13:00 – 13:45

Design and Implementation of Novel USP and DSP Continuous Biomanufacturing Concepts and Technologies: Experiences, practical examples and future outlook

Gerben Zijlstra,
Senior consultant, Xendo, Netherlands

13:45 – 14:15

Risk Assessment Strategies for Lifecycle Management using Multivariate Data Analysis

Prof. José Menezes,
Founder & Managing Director, 4Tune Engineering GmbH, Germany

14:15 – 15:40

Coffee Break | Exhibition

15:40 – 16:10

Process Analytical Technology (PAT) for Bioprocessing
Cyrus Agarabi,
Sr. Regulatory Research | Review Officer, FDA, US

16:10 – 16:40

Continuous Manufacturing, Hybrid Options Transitioning from Batch to Continuous Economics in Downstream Processing

Andrew Sinclair,
President & Founder, Biopharm Services, UK

16:40 – 17:00

Q&A | Discussion

17:00
Transfer to Hotel

18:15
Transfer to Networking Dinner

21:45
Transfer back to Hotel

Agenda

Thursday, June 09, Downstream

08:30 – 09:00
Coffee | Exhibition

09:00 – 09:30

Single Use Equipment as enabler of ADC conjugation in a standard protein plant using closed systems
Dr. Berthold Boedeker,
Chief Scientist, Global Biological Development WUP, Bayer, Germany

09:30 – 10:00

Smarter DSP design based on early-stage development data
Dr. Michael Box,
Purification Development, Lonza Biologics, UK

10:00 – 10:30
Coffee Break | Exhibition

► **Virus Removal Trends and Challenges**

10:30 – 11:00

The viral load and the use of pre-filter in viral clearance studies of virus filtration
Dr. Horst Ruppach,
Global Director, Viral Clearance and Virology, Charles River, Germany

11:00 – 11:30

Improving DSP operations of emerging biopharmaceuticals: reinventing old school processes
Cristina Peixoto,
Head Downstream Process Development, Animal Cell Technology Unit, IBET, Portugal

11:30 – 12:00

Virus Safety Aspects in the Development of a Downstream Process for Recombinant Proteins
Franz Nothelfer,
Associate Director Protein Science, Boehringer Ingelheim Pharma, Germany

12:00 – 13:00
Lunch | Exhibition

13:00 – 13:30

Optimizing sulfated cellulose membrane adsorbers to purify influenza virus
Raquel Fortuna, PhD student,
Bioprocess Engineering group, Max-Planck Institute – Magdeburg, Germany

13:30 – 13:50

Q&A | Discussion

13:50 – 14:00

Meeting closure

14:00 – 15:30

Plant tour

We look forward to welcoming you in Goettingen, Germany.

Conference address

Sartorius College
Otto Brenner Str. 20
37079 Goettingen, Germany
Phone +49.551.308.2101

Welcome dinner – Tuesday

Restaurant Orient Lounge (Freizeit In)

Dransfelder Str. 3
37079 Goettingen
Phone: +49.551.90010

Networking dinner – Wednesday

Hotel FREIgeist

Am Gesundbrunnen
37154 Northeim
Phone: +49.5551.607.0

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

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

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

Registration

Early bird rate (until April 29th):

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

Standard price (from April 30th):

€ 550 + 19 % VAT
€ 150 + 19 % VAT for academia

This includes the networking dinner, all breaks and lunches.

To register, please either

- register on the Web:
www.sartorius.com/upstream-downstream-forum
- call or e-mail listed contacts

Accommodation

We have arranged limited allotments in the following hotels.

Please mention "Upstream & Downstream Technology Forum".

Hotel Rennschuh

Kasseler Landstrasse 93
37081 Goettingen
Phone: +49.551.9009.0

www.rennschuh.de
€ 57 incl. breakfast

Park Inn by Radisson Goettingen

Kasseler Landstrasse 25c
37081 Goettingen
Phone +49.551.270707.520

www.parkinn-goettingen.de
€ 108 – 128 incl. breakfast

BEST WESTERN Parkhotel Ropeter

Kasseler Landstraße 45
37081 Göttingen
Phone +49.551.9020

www.bestwestern.de/hotels/Goettingen/
BEST-WESTERN-Parkhotel-Ropeter
€ 81 incl. breakfast

Best Western Hotel am Papenberg

Hermann-Rein-Strasse 2
370759 Goettingen
Phone +49.551.30550

www.hotel-am-papenberg.bestwestern.de
€ 95 incl. breakfast



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.

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