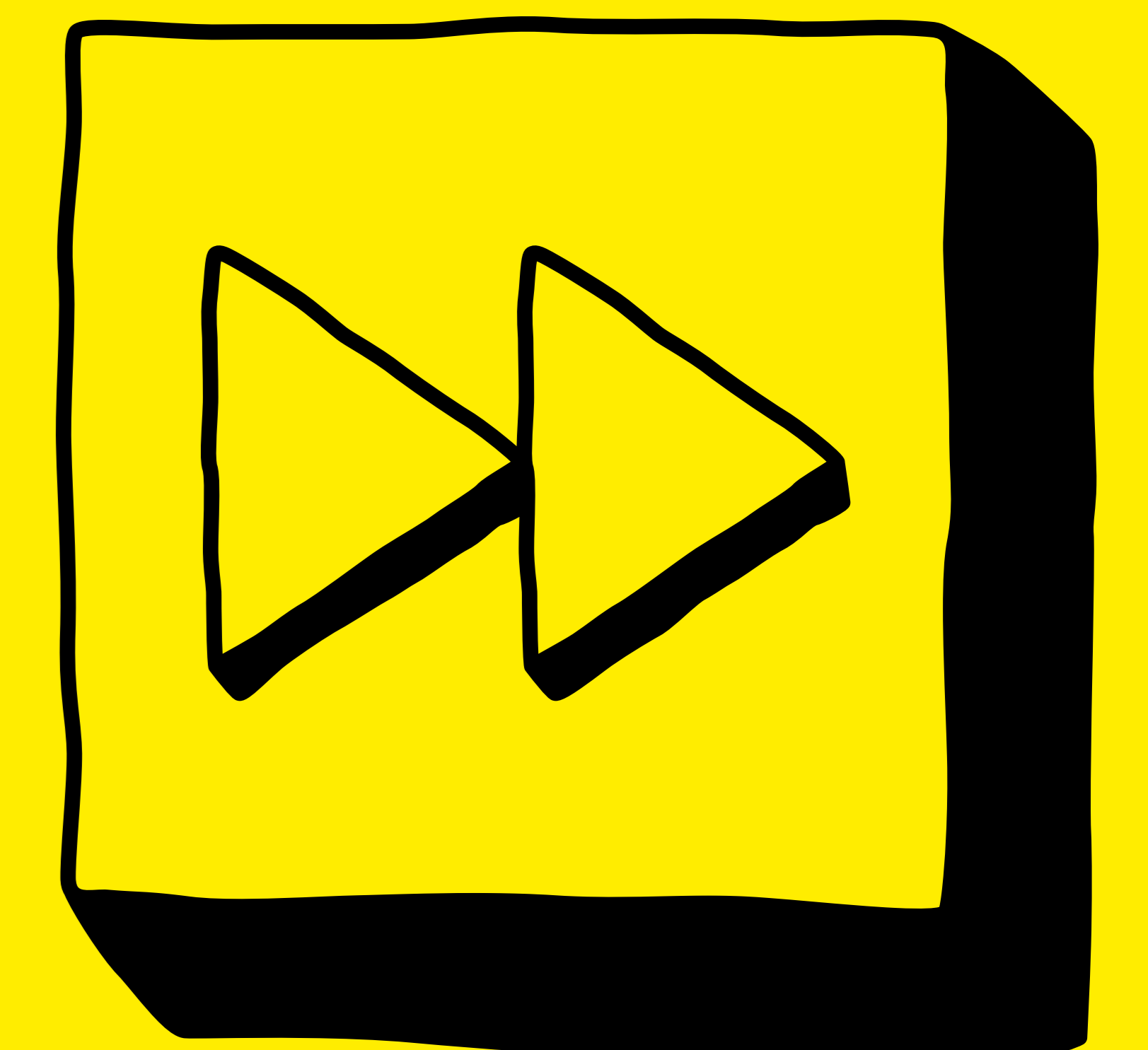
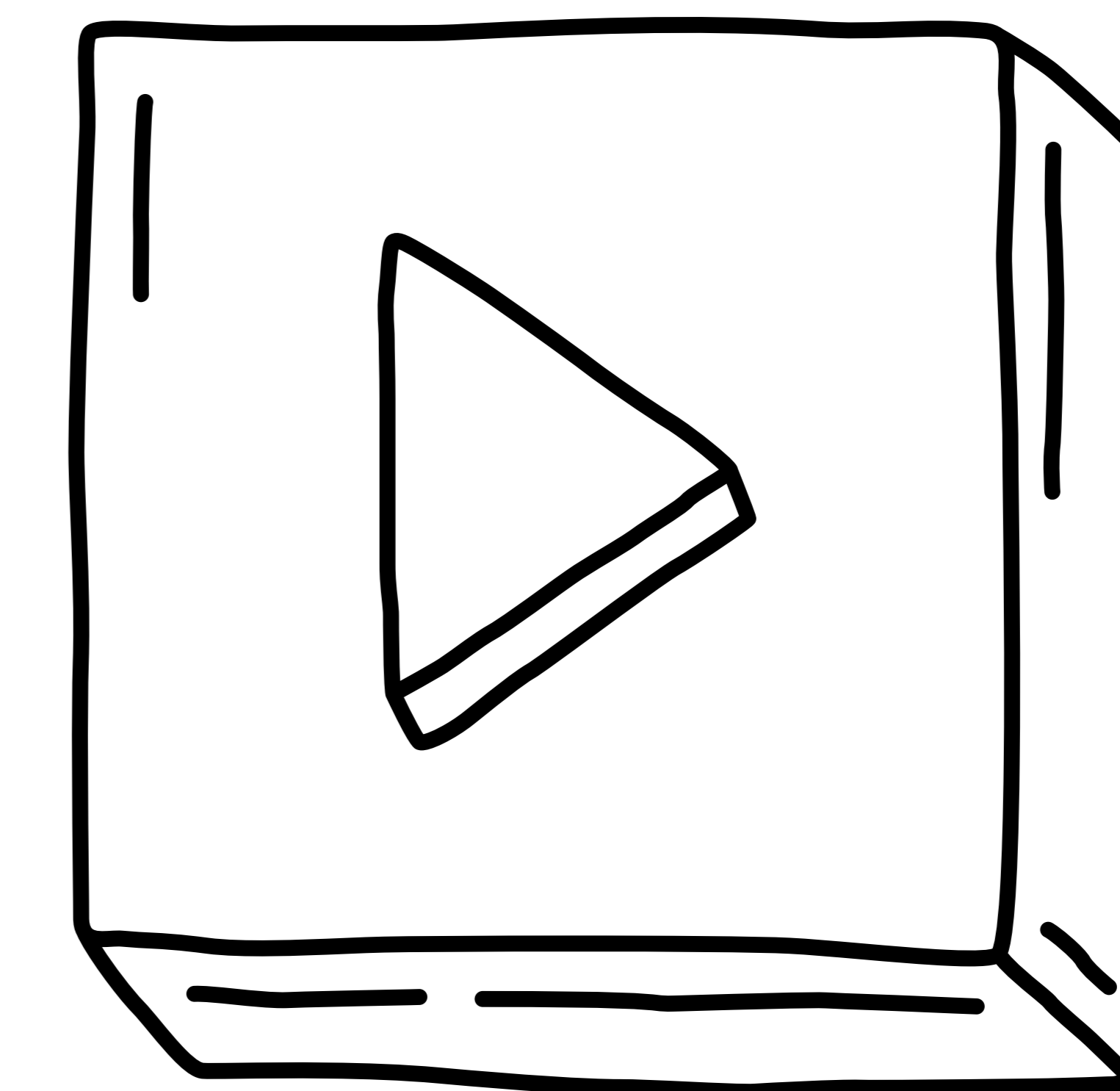


# SARTORIUS

September 15-17 | Göttingen, Germany

Sartorius PI Forum 2026 —  
It's time to intensify



**Sartorius PI Forum Germany 2026**  
**Day 1 | Tuesday, September 15 | Morning**

Time	Main plenary room	Speaker	Virtual experience (w/ Meta glasses)	Poster Session
8:00 a.m.	Registration & coffee			
9:00 a.m.	Keynote lectures			
9:00 a.m.	Welcome   Opening   It's time to intensify!	<b>Miriam Monge</b> , Head of Customer & Industry Advocacy Strategy BPS, Sartorius		
9:20 a.m.	Title of talk: tbc	<b>Konstantin Konstantinov</b> , PhD, CTO (Ring Therapeutics) & Operating Partner (Flagship Pioneering)		
9:50 a.m.	Integrated continuous bioprocessing to enable significant productivity gains via process intensifications	<b>Weichang Zhou</b> , PhD, Chief Technology Officer, MediLink Therapeutics		
10:20 a.m.	Coffee break			
10:50 a.m.	Redefining biologics manufacturing: Process intensification from development to the clinic and launch	<b>Nihal Tugcu</b> , PhD, Global Head, Mammalian Platform, Global CMC Development, Sanofi		
11:20 a.m.	Continuous precipitation and flocculation: An underutilized unit operation in integrated continuous biomanufacturing	<b>Alois Jungbauer</b> , PhD, Professor, BOKU University		
11:50 a.m.	Panel discussion with keynote speakers			
12:15 p.m.	Networking lunch			
1:15 p.m.	The boardroom playbook – Economic insight for next-gen manufacturing	<b>Suzanne Farid</b> , PhD, Head of Department & Professor of Bioprocess Systems Engineering, UCL Biochemical Engineering <b>Nitin Chopra</b> , Senior Expert Business Intelligence, Sartorius <b>John Erickson</b> , Senior Fellow, NIIMBL	Facility of the future with full virtual PI experience	Poster session
2:20 p.m.	Transfer to break-out rooms			

**Event homepage:**  
[Sartorius PI Forum Germany 2026](#)

Sartorius PI Forum Germany 2026  
Day 1 | Tuesday, September 15 | Afternoon

Time	Main plenary room	Speaker	Virtual experience (w/ Meta glasses)	Poster Session	Workshops (Visitor Center)	Workshops (Visitor Center)	Demo (R&D)	Demo (Application & Training Center)	Demo (Application & Training Center)	General Application Center Tours
DSP intensification: From theory to practice			Facility of the future with full virtual PI experience	Poster session	Workshop data analytics AI on the BioProcess shop floor: Real data, real processes, real results	Workshop USP Engineering the future of CHO: Smarter cells, smarter media, higher yields	R&D Demo PI in action: Scaling mAbs from development to early GMP			
2:30 p.m.	Advantages of protein A membrane chromatography: High productivity, improved resolution, and more	<b>Li Yifeng, PhD</b> , Executive Director, Head of Downstream Process Development, WuXi Biologics								
3:00 p.m.	Affinity chromatography modelling to leverage multicolumn chromatography development	<b>Alan Deduyer</b> , Downstream Process Scientist, Sanofi								
3:30 p.m.	Coffee break									
USP   Advanced process control			Facility of the future with full virtual PI experience			Process modelling workshop Using ExCIT to identify the best PI technology for your process	R&D Demo PI in action: Scaling mAbs from development to early GMP	Pionic® Demo Our platform makes DSP intensification easier using a turnkey, flexible solution	Resolute® BioSMB Demo See how easy it is to develop MCC and scale it up to cut costs and buffer consumption	
4:00 p.m.	TFF perfusion for process intensification: Lessons learned and future directions	<b>Patrick Romann</b> , Bioprocess Innovation Manager, Levitronix GmbH								
4:30 p.m.	Intensified upstream processing: Fed-batch and perfusion options for robust scale-up	<b>Nuno Buxo Carinhas</b> , Senior Expert in Science & Technology, Novartis Pharma AG								
5:00 p.m.	Collaborative workshop   Ensuring a successful implementation of PI technologies	<b>Geoffrey Pressac</b> , Senior Expert Sales Development Chromatography, Sartorius								
5:45 p.m.	End of sessions									
6:00 p.m.	Bus transfer to dinner									
7:00 p.m.	PI Forum gala dinner									

Sartorius PI Forum Germany 2026  
Day 2 | Wednesday, September 16 | Morning

Time	Main plenary room	Speaker	Virtual experience (w/ Meta glasses)	Poster Session	Workshops (Visitor Center)	Workshops (Visitor Center)	Demo (R&D)	Demo (Application & Training Center)	Demo (Application & Training Center)	General Application Center Tours
8:00 a.m.	Check-in & coffee									
	Keynote presentations									
8:30 a.m.	Combining new generation membrane adsorber with connected DSP systems and advanced data monitoring	<b>Benoît Mothes</b> , Head of Purification Process Development, Sanofi, Vitry								
9:00 a.m.	tbc									
9:30 a.m.	Transfer to breakout rooms									
	USP   Advanced process control									
9:45 a.m.	Title of talk: tbc	<b>Charles Heise</b> , PhD, Associate Director (BPSD, Process Development), FUJIFILM Biotechnologies	Facility of the future with full virtual PI experience	Poster session	Workshop data analytics AI on the BioProcess shop floor: Real data, real processes, real results	Process modelling workshop Using ExCIT to identify the best PI technology for your process	R&D Demo PI in action: Scaling mAbs from development to early GMP	Pionic® Demo Our platform makes DSP intensification easier using a turnkey, flexible solution	Resolute® BioSMB Demo See how easy it is to develop MCC and scale it up to cut costs and buffer consumption	
10:15 a.m.	Title of talk: tbc	<b>Sadia Ahmed Sugulle</b> , USP Process Expert, Merck								
10:45 a.m.	Coffee break									
	DSP intensification: From theory to practice									
11:15 a.m.	Title of talk: tbc	<b>Benjamin Perruchoud</b> , DSP Scientist, Merck	Facility of the future with full virtual PI experience		Workshop data analytics AI on the BioProcess shop floor: Real data, real processes, real results	Workshop USP Unlocking ultra-high cell density: Bioreactor innovations driving process intensification		USP Demo Latest PAT, advanced automation, data analytics and integrated hardware platforms for PI		Application Center Tour to experience true end-to-end solutions
11:45 a.m.	Title of talk: tbc	<b>Innocent Bekard</b> , Director, Technical Operations, CSL Limited								
12:15 p.m.	Networking lunch			Poster session						

Sartorius PI Forum Germany 2026  
Day 2 | Wednesday, September 16 | Afternoon

Time	Main plenary room	Speaker	Virtual experience (w/ Meta glasses)	Poster Session	Workshops (Visitor Center)	Workshops (Visitor Center)	Demo (R&D)	Demo (Application & Training Center)	Demo (Application & Training Center)	General Application Center Tours
DSP intensification: From theory to practice			Facility of the future with full virtual PI experience	Poster session		Workshop USP Advancing continuous bio-manufacturing: Innovations in mAb development, scale-up, and early GMP for the Pionic® Platform		USP Demo Latest PAT, advanced automation, data analytics and integrated hardware platforms for PI		Application Center Tour to experience true end-to-end solutions
1:15 p.m.	Facilitating process intensification through mechanistic modelling: Application to the purification of a plasma protein via aptamer affinity chromatography	<b>Damien Bataille</b> , Head of Process Development Dept., LFB Biotechnologies <b>Edouard Nicoud</b> , Co-CEO, Ypso-Facto								
1:45 p.m.	Design of experiments (DoE)-driven optimization of sequential multi-column chromatography for IEX capture and flowthrough polishing of plasma-derived IgG	<b>Daniel Strohmer</b> , Scientist at the Plasma Innovation Lab, Takeda Manufacturing Austria AG								
2:15 p.m.	Title of talk: tbc	Asymbio								
2:45 p.m.	Coffee break									
USP   Advanced process control			Facility of the future with full virtual PI experience	Poster session		Process modelling workshop Using ExCIT to identify the best PI technology for your process	R&D Demo PI in action: Scaling mAbs from development to early GMP	Pionic® Demo Our platform makes DSP intensification easier using a turnkey, flexible solution	Resolute® BioSMB Demo See how easy it is to develop MCC and scale it up to cut costs and buffer consumption	
3:15 p.m.	Automating nutrient control in bench-scale continuous bioreactors using capacitance	<b>Lorraine Peters</b> , Associate Scientist, AstraZeneca								
3:45 p.m.	The McMaster-Sartorius partnership: Empowering innovation and future leaders in bioprocess automation	<b>David LaTulippe</b> , Professor, McMaster University <b>Brandon Corbett</b> , Principal Scientist, Advanced Data Analytics, Sartorius								
4:15 p.m.	Scalability of Lonza's intensified fed batch process with automation	<b>Anthony Beaney</b> , Principal Scientist, Lonza Biologics <b>Anita Dabek</b> , Associate Principal Scientist, Lonza Biologics								
4:45 p.m.	Workshop feedback session & wrap-up									
5:30 p.m.	Bus transfer									
6:30 p.m.	PI Forum networking celebration									