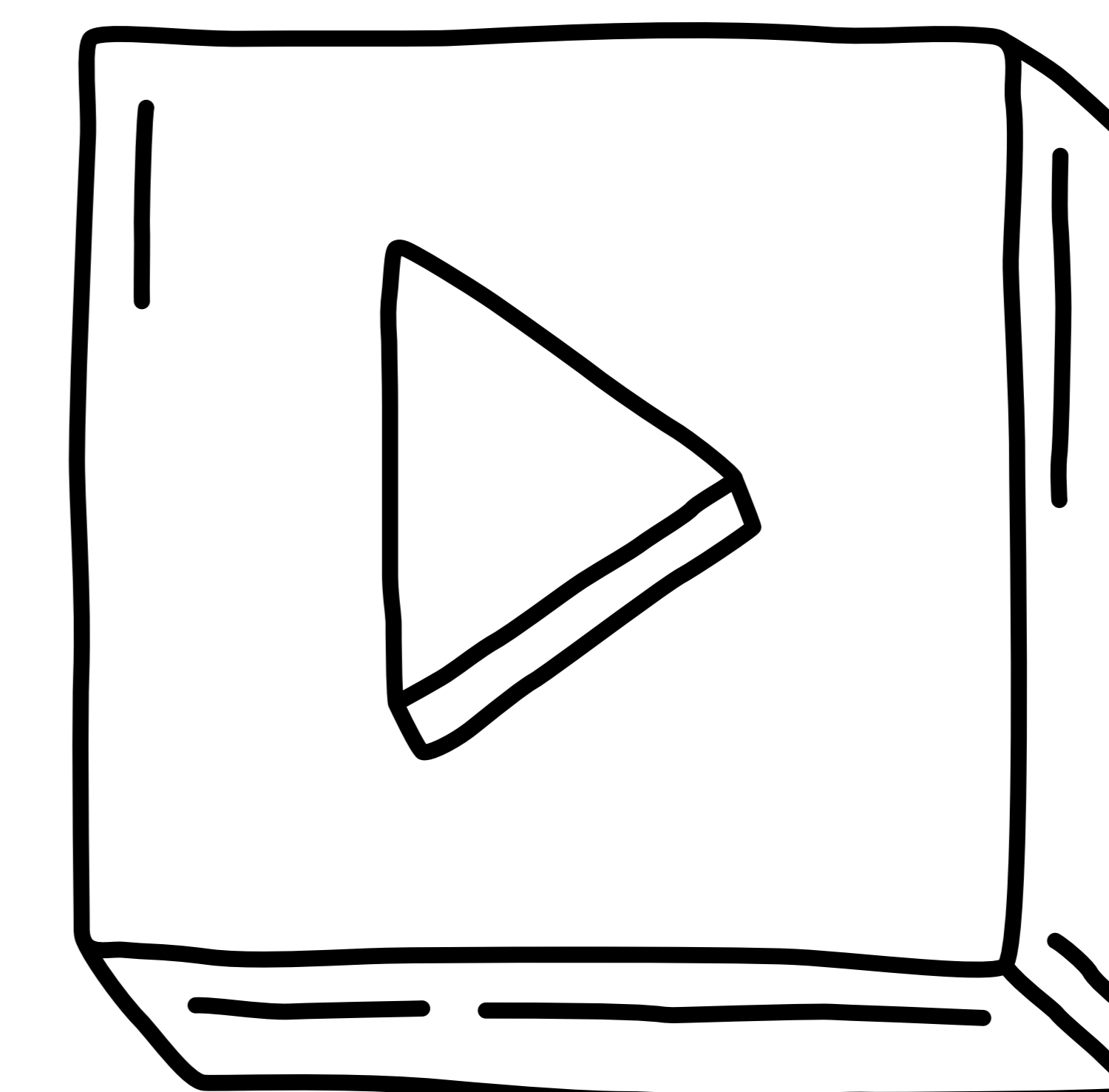


SARTORIUS



June 3-4 | Hyderabad, India

Sartorius PI Forum India 2026 —
It's time to intensify

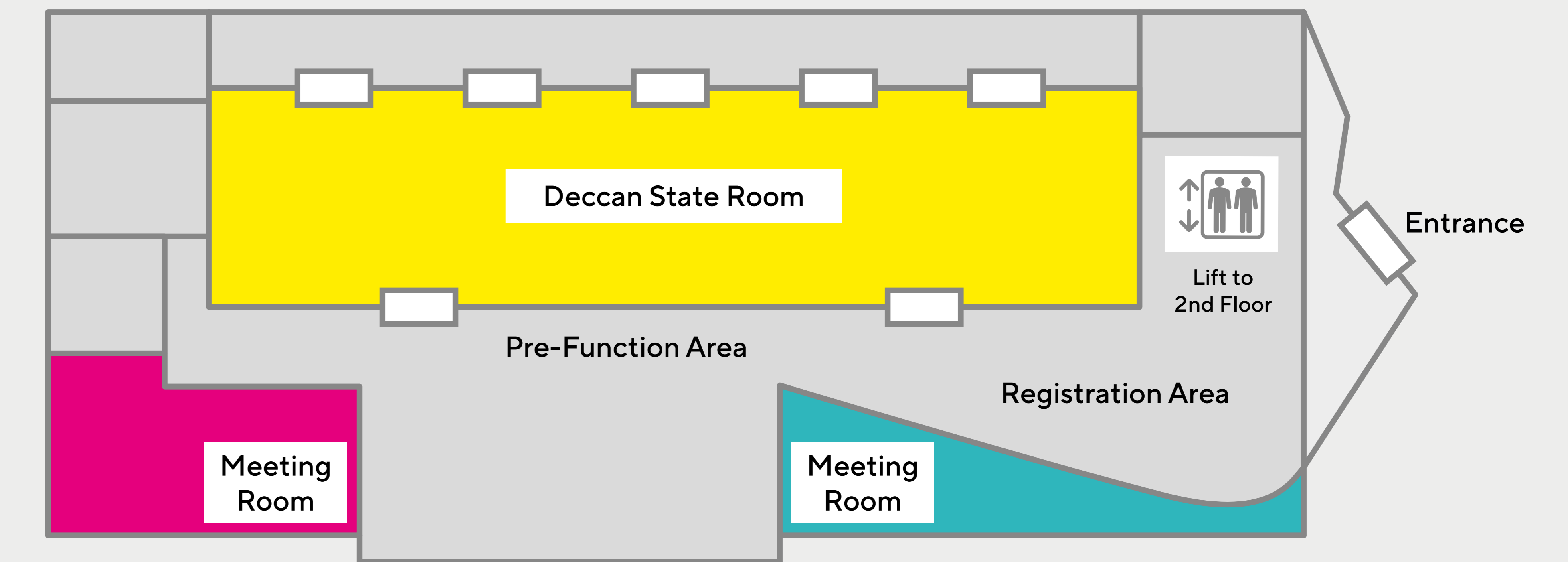


Sartorius PI Forum India 2026
Day 1 | Wednesday, June 3

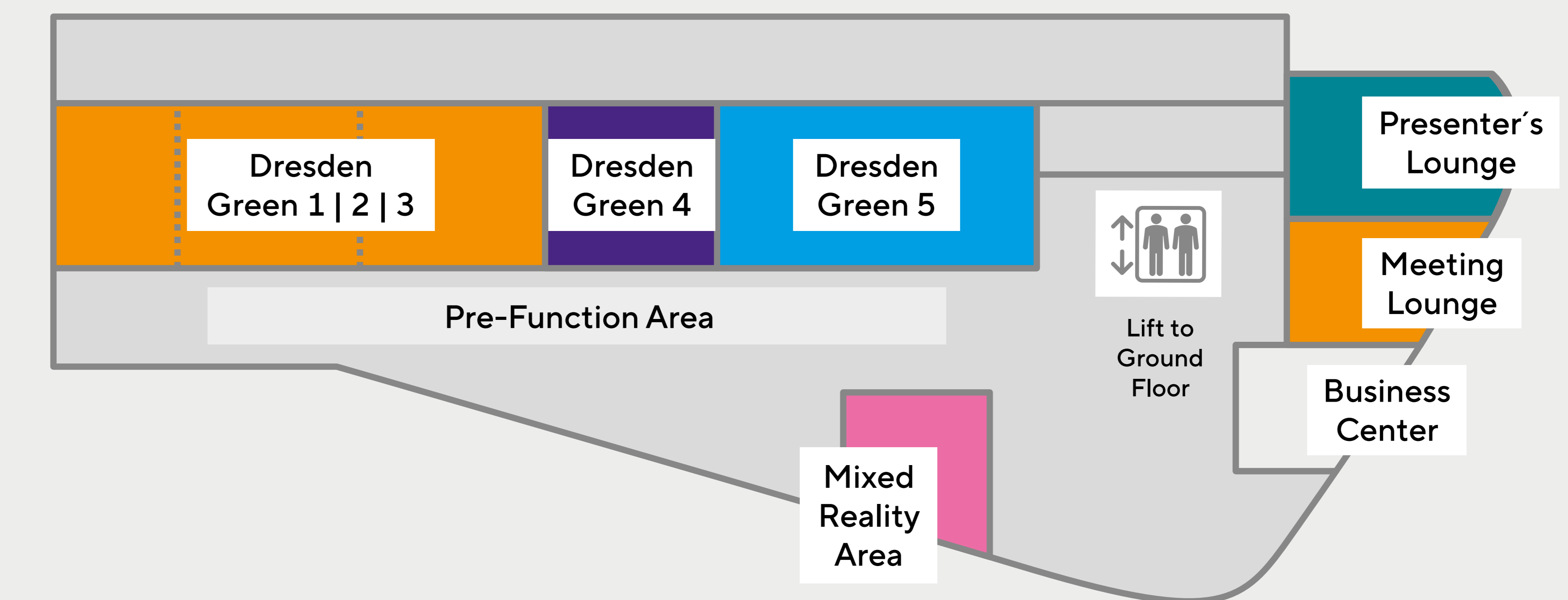
Time	Track 1 Deccan State Room	Speaker	Virtual experience (with Meta glasses)
9:00 a.m.	Registration & coffee		
	Plenary talks	Moderator: Miriam Monge, Sartorius	
10:00 a.m.	Welcome Opening It's time to intensify!	Miriam Monge , Head of Customer & Industry Advocacy Strategy BPS, Sartorius	
10:30 a.m.	Accelerating intensified process development: Evolving from custom workflows to platform approach	Karthik P. Jayapal , Director, Biologics Process Development, Merck, Sharp & Dohme	
11:00 a.m.	Title of talk: tbc	Speaker: Shilpa Gadgil , VP, Head of Bioprocess Development, Head of CDMO Development, Enzene Biosciences Ltd	
11:30 a.m.	Coffee break		
12:00 p.m.	Modernizing bioprocessing: Continuous manufacturing adoption, COGs optimization, and workforce challenges	Kandarp N. Raval , Associate Vice President Manufacturing, Kashiv Biosciences LLC	
12:30 p.m.	Process intensification at Lonza: A platform journey from seed train to DSP	Hemanth Kaligotla , Director of External Innovation, Lonza Biopharma	
1:00 p.m.	Networking lunch		Facility of the future with full virtual PI experience
	Intelligent bioprocessing	Moderator: Priyanka Gupta	
2:00 p.m.	The boardroom playbook – Economic insight for next-gen manufacturing including panel discussion: Hemanth Kaligotla, Kandarp N. Raval, Karthik P. Jayapal, Shilpa Gadgil, Sandeep Majumdar and Miriam Monge	Nitin Chopra , Senior Expert Business Intelligence, Sartorius Priyanka Gupta , Senior Expert Bioprocess Solutions, Sartorius	
3:30 p.m.	Coffee break		
4:00 p.m.	AI-driven virtual cell platforms for biologics development and manufacturing	Ling Wu , President & Co-Founder, ChemT Biotechnology	
4:30 p.m.	Transforming biosimilars scale-up through process intensification	Sandeep Majumdar , Site Head & Associate Executive Vice President, Intas Pharmaceuticals	
5:00 p.m.	Regulation of biologics in India and provisions for adoption of innovative technologies	Arvind Kukrety , Deputy Drug Controller, CDSCO, HQ, New Delhi	
5:30 p.m.	Workshop Overcoming the bottlenecks to implement PI in commercial manufacturing		
6:15 p.m.	End of sessions		Facility of the future with full virtual PI experience
7:00 p.m.	Networking dinner		

Floorplan:
ITC Kohenur, a Luxury Collection Hotel, Hyderabad, India

Ground Floor: Deccan State Room



Second Floor: Breakout Rooms



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

Sartorius PI Forum India 2026
Day 2 | Thursday, June 4 | Morning

Time	Track 1 Deccan State Room	Speaker	Virtual experience (with Meta glasses)	Workshops (breakout rooms)	Workshops (breakout rooms)	Workshops (breakout rooms)
8:30 a.m.	Check-in & coffee	Moderator: Priyanka Gupta				
	Plenary talks					
9:30 a.m.	Rethinking biopharma manufacturing: From scale to agility through process intensification	Dr. Rhonda Duffy, COO, Biocon Ltd				
10:00 a.m.	Bio-pharmaceuticals process intensification: Hypes or hopes	Dr. Sanjeev Gupta, Sr. Vice President & Head Biosimilar Advanced Biotech Lab, IPCA Laboratories				
10:30 a.m.	Coffee break & transfer to breakout rooms					
	USP Advanced process control	Moderator: Gerben Zijlstra	Facility of the future with full virtual PI experience	Downstream breakout rooms		
11:00 a.m.	Making intensified upstream processes work at scale: Strategies, challenges and practical solutions	Ankur Bhatnagar, Vice President & Head of R&D CMC, Biocon Biologics Limited	Facility of the future with full virtual PI experience	DSP Pionic® workshop	DSP workshop: Virus clearance strategies	DSP workshop: Membrane process
11:30 a.m.	Biomass specific perfusion rate as a control lever for the continuous manufacturing of biosimilar antibodies	Eunice Leong, Senior Scientist, A*STAR Bioprocessing Technology Institute		Breaking the downstream bottleneck: Enabling cost effective mAb manufacturing through integrated continuous bioprocessing	Overcome challenges in virus clearance validations in connected and continuous downstream processing	Membrane chromatography for mAbs: Intensifying processes, cutting time and cost
12:00 p.m.	Enhancing AAV titers in HEK293 cells: The power of perfusion and addressing high cell density transfection barriers	Dr. Dennis Karthaus, Manager of Media and Processes, Media and Process Development Xell, Sartorius				
12:30 p.m.	Networking lunch					
	Fitting DSP to higher titers and stronger cost pressure	Moderator: Jean-Luc Beulay	Facility of the future with full virtual PI experience	Upstream breakout rooms		
1:30 p.m.	Our perspective on process intensification: Enabling FCCMTM an intensified process platform for cost effective production of biologics	Abijar Bhuri, Associate Vice President (AVP), Enzene Biosciences Ltd. Pune	Facility of the future with full virtual PI experience	USP workshop	USP PAT – digital demo	
2:00 p.m.	Accelerating downstream efficiency through capture process intensified multi-column capture chromatography	Vijay Sudalaimuthu, Vice President & Head of DS Biosimilars manufacturing, Biocon Biologics Limited		Interactive technology overview: Where upstream PI can be implemented and what value it will bring	Experience Sartorius' latest upstream PI technologies up-close and hands-on	
2:30 p.m.	Transitioning from AEX resin to membrane chromatography: A fit-for-purpose downstream strategy	Arpit Jain, Vice President, Bio Bulk Manufacturing, Intas Pharmaceutical-Biopharma Division				
3:00 p.m.	Coffee break					

Sartorius PI Forum India 2026
Day 2 | Thursday, June 4 | Afternoon

Time	Track 1 Deccan State Room	Speaker	Virtual experience (with Meta glasses)	Workshops (breakout rooms)	Workshops (breakout rooms)	Workshops (breakout rooms)
	USP Advanced process control	Moderator: Henry Weichert	Facility of the future with full virtual PI experience	Downstream breakout rooms		
3:30 p.m.	Engineering immune tolerance with regulatory T cells: Design of experiments as a decision tool for cell therapy manufacturing	Esmond Lee Yan Zhao , Senior Scientist, Bioprocessing Technology Institute, A*Star Singapore		DSP Pionic® workshop	DSP workshop: Virus clearance strategies	DSP workshop: Membrane process
4:00 p.m.	Process intensification and analytics: A duo to harness complexity in bioprocessing	Kumaran S V , Fuctional Leader (Upstream), Biocon Biologics Limited	Facility of the future with full virtual PI experience	Breaking the downstream bottleneck: Enabling cost effective mAb manufacturing through integrated continuous bioprocessing	Overcome challenges in virus clearance validations in connected and continuous downstream processing	Membrane chromatography for mAbs: Intensifying processes, cutting time and cost
4:30 p.m.	Coffee break					
	Fitting DSP to higher titers and stronger cost pressure	Moderator: Jean-Luc Beulay	Facility of the future with full virtual PI experience	Upstream breakout rooms		
5:00 p.m.	Title of talk: tbc	Dr. Rahul Bhambure , Principal Scientist, National Chemical Laboratory		Digital Solutions Workshop	USP PAT – digital demo	
5:30 p.m.	tbc	tbc		Showing the capabilities of PAT, automation and data analytics	Experience Sartorius' latest upstream PI technologies up-close and hands-on	
6:00 p.m.	Panel discussion Chair – Miriam Monge	Esmond Lee Yan Zhao , Rhonda Duffy , Dr. Sanjeev Gupta , Yatindra Tirunagari , Dr. Gerben Zijlstra & Priyanka Gupta				
6:45 p.m.	End of sessions					
7:00 p.m.	Networking dinner					