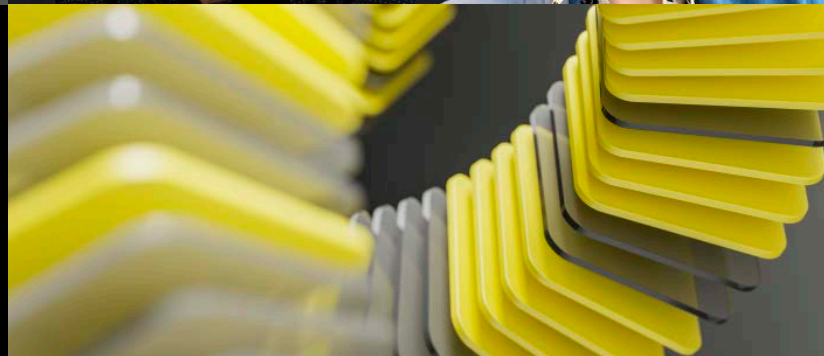
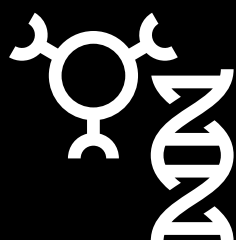
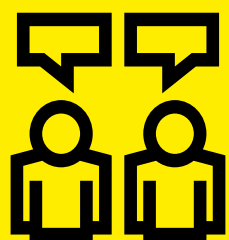
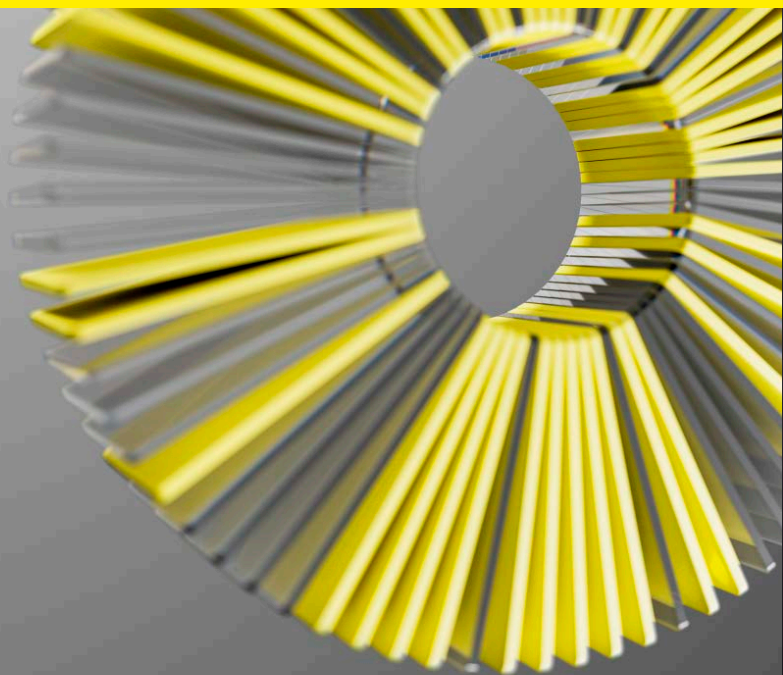


SARTORIUS

Sartorius Cell & Gene Therapy Forum

Driving the Future of Innovation
in Cell & Gene Therapy, Together

October 23, 2025
Boston, MA, USA



Sartorius Cell & Gene Therapy Forum

We are pleased to invite you to this year’s Sartorius Cell & Gene Therapy Forum, in Boston, MA, on October 23. The day will begin with a plenary session and a dynamic panel discussion featuring leaders from across the cell and gene therapy landscape.

In the afternoon, attendees may choose between dedicated cell therapy and gene therapy tracks, followed by an opportunity to engage with peers and thought leaders during our cocktail reception.

As a pre-event activity, you have the option to sign up for a guided tour of the new Sartorius Center for Bioprocess Innovation in Marlborough, MA, on the afternoon of October 22.

Please see our detailed agenda and speaker information below.

Venue:
MassBio
700 Technology Square
Cambridge, MA 02139, USA



Event homepage
www.sartorius.com/en/company/exhibition-conferences/local-events/cgt-forum-2025-boston

Sartorius CGT Forum – At a Glance

Wed, Oct 22		Thursday, Oct 23	
		Registration Meet & Greet	
		8:30 a.m. • Welcome & Opening – Priyanka Gupta , Sartorius Maya Fürstenau-Sharp , Sartorius Maurice Phelan , Sartorius • Carl Schoellhammer , DeciBio Consulting • Regulatory Perspectives in ATMP Development and Manufacturing – Basak Clements , BioMatria Doli Patel , Sartorius • Fireside Chat: Navigating Investor Expectations in Cell & Gene Therapy – Phil Vanek , Redline Bioadvisors • Reinventing Cell & Gene Therapy Development and Manufacturing with Agentic AI – Sanjay Srivastava , Accenture • Strategies for CGT Start-Ups – Tony Ting , Kijji Therapeutics • Panel Discussion with Keynotes	
		12:15 p.m. Networking Lunch	
1:00 p.m. - 5:45 p.m.		1:15 p.m.	Gene Therapy Track Cell Therapy Track
Sartorius Center for Bioprocess Innovation Marlborough Tour		• Case Study in AAV8 End-to-End Bioprocessing – Trent Lyman, Matica Biotechnology • Empowering 5.0 BioSolutions Through Realtime PAT and Computational Controls Reducing Waste, While Saving Time and Cost – Edita Botonjic-Sehic & Aaron Cowley, RecBioPharm • Kinetically Controlled Assembly and Machine-Learning Optimization of Gene-Delivery Nanoparticles – Hai-Quan Mao, Johns Hopkins University • Case Study: Leveraging Platform Technologies for Plasmid Production – Dennis Powers, G-CON Miriam Monge, Sartorius • The Power of Data – Combining Analytics and Model-Driven Approaches to Optimize Media and Processes in Viral Vector Production – Sandra Klausning, Sartorius • Scaling Cardiomyocyte Cell Production Towards 3D Bioprinting of Organ-Scale Tissues – Mark Skylar-Scott, Stanford University • Driving Cell Therapy Innovation: Integrated Automation and Non-Viral Gene Editing – Chesney Michels, Elevate Bio • Advances in Scale-up and Optimization of Operating Conditions for iPSC Bioprocess Development – Sarah Gilpin, Sartorius • Removing Barriers to Sustainable Autologous Therapies – Pierre Springuel, UCL • Academic Medical Cell Therapy Development, Manufacturing, and Delivery at Dana-Farber Cancer Institute – Julie Porter Sarah Nikiforow, Dana-Farber Cancer Institute	
6:00 p.m. - 9:00 p.m.		Closing Remarks & Feedback Session	
Dinner in Marlborough and bus back to Boston		5:30 p.m. - 7:30 p.m. Networking Cocktail Reception	

Agenda

Wednesday, October 22

Join us for an exclusive tour of the Sartorius Innovation Center in Marlborough, MA.

Designed to foster collaboration and real-world application, this state-of-the-art facility showcases the latest Sartorius technologies driving cell and gene therapy forward. Experience firsthand our next-gen automated cell therapy platforms, LNP generation tools, and Process Development & Clinical GMP Services. Discover how our experts, workflows, and the innovative technologies developed by Corporate Research are shaping the future of therapeutic development.

Start	Activity
1:00 p.m.	Check-in at Residence Inn Boston, 120 Broadway, Cambridge, MA 02142
1:30 p.m.	Bus Departs from Hotel
2:45 p.m.	Arrival at Destination
2:50 p.m.	Welcome and Introductions
3:00 p.m.	Guided Tour Begins
5:30 p.m.	Tour Concludes
5:45 p.m.	Board Bus for Dinner
6:00 p.m.	Group Dinner in Marlborough
9:00 p.m.	Return to Residence Inn Boston

Thursday, October 23

Start	End	Topic	Speaker Chair: Maya Fuerstenau-Sharp
8:30 a.m.	8:30 a.m.	Registration	
8:30 a.m.	8:50 a.m.	Welcome to the CGT Forum Opening Remarks	Priyanka Gupta , Project Lead CGT Forum, Advocacy Team, Sartorius Maya Fürstenau-Sharp , Head of Marketing ATS, Sartorius Maurice Phelan , President of Sartorius North America Inc.
8:50 a.m.	9:15 a.m.	Vectors, Cells, and Valuations: Mapping the CGT Market Moment	Carl Schoellhammer , PhD, Partner, DeciBio Consulting
9:15 a.m.	9:55 a.m.	Regulatory Perspectives in ATMP Development and Manufacturing	Basak Clements , Founder and Senior Advisor, Biomatria Doli Patel , Head of Quality and Regulatory Affairs BPS, Sartorius Moderator: Mark Hites , Regulatory Affairs Professional, Sartorius
9:55 a.m.	10:25 a.m.	Fireside Chat Funding the Future: Navigating Investor Expectations in Cell & Gene Therapy	Philip Vanek , Founder & CEO, Redline Bioadvisors Chief Commercialization Officer, ISCT Moderator: Tiffany Pogue , Business Development Professional, Advanced BioProcessing, Sartorius
10:25 a.m.	10:45 a.m.	Coffee Break	
10:45 a.m.	11:15 a.m.	Reinventing Cell & Gene Therapy Development and Manufacturing with Agentic AI	Sanjay Srivastava , Managing Director, Cell & Gene Therapy CoE Lead, Accenture
11:15 a.m.	11:45 a.m.	Strategies for CGT Startups	Tony Ting , Chief Scientific Officer, Kiji Therapeutics
11:45 a.m.	12:15 p.m.	Panel Discussion with Keynotes	Carl Schoellhammer , Philip Vanek , Sanjay Srivastava , Tony Ting Moderator: Steve Binnering , Head of Collaborative Technology Development, Sartorius
12:15 p.m.	1:15 p.m.	Lunch	

Agenda

Thursday, October 23

Start	End	Topic	Speaker Chair: Paul Cashen
Gene Therapy & NA Track			
1:15 p.m.	2:00 p.m.	Case Study in AAV8 End-to-End Bioprocessing	Trent Lyman , Head of Process Development, Matica Biotechnology
2:00 p.m.	2:45 p.m.	Empowering 5.0 BioSolutions Through Realtime PAT and Computational Controls Reducing Waste, While Saving Time and Cost	Edita Botonjic-Sehic , Head of Process Analytics and Data Science, RecBioPharm Aaron Cowley , Chief Scientific Officer, RecBioPharm
2:45 p.m.	3:15 p.m.	Coffee Break	
3:15 p.m.	3:45 p.m.	Kinetically Controlled Assembly and Machine-Learning Optimization of Gene-Delivery Nanoparticles	Hai-Quan Mao , Director, Institute for NanoBioTechnology; Professor, Materials Sci & Eng, and Biomed Eng, Johns Hopkins University
3:45 p.m.	4:15 p.m.	Case Study: Leveraging Platform Technologies for Plasmid Production	Dennis Powers , Senior Vice President of Product & Strategy, G-CON Miriam Monge , Head of Customer & Industry Advocacy Strategy BPS, Sartorius
4:15 p.m.	4:45 p.m.	The Power of Data - Combining Analytics and Model-Driven Approaches to Optimize Media and Processes in Viral Vector Production	Sandra Klausing , Head of Product Development Cell Line and Media Solutions (CMTS), Sartorius
4:45 p.m.	5:00 p.m.	Short Break	
5:00 p.m.	5:30 p.m.	Closing Remarks & Feedback Session	Priyanka Gupta , Project Lead CGT Forum, Advocacy Team, Sartorius Maya Fürstenau-Sharp , Head of Marketing ATS, Sartorius
5:30 p.m.	7:30 p.m.	Networking Cocktail Reception	

Start	End	Topic	Speaker Chair: Shanya Jiang
Cell Therapy Track			
1:15 p.m.	1:45 p.m.	Scaling Cardiomyocyte Cell Production Towards 3D Bioprinting of Organ-Scale Tissues	Mark A. Skylar-Scott , Assistant Professor, Stanford Department of Bioengineering
1:45 p.m.	2:15 p.m.	Driving Cell Therapy Innovation: Integrated Automation and Non-Viral Gene Editing	Chesney Michels , Director of Innovation and Platform for Cell & Gene Therapy, Elevate Bio
2:15 p.m.	2:45 p.m.	Advances in Scale-up and Optimization of Operating Conditions for iPSC Bioprocess Development	Sarah Gilpin , Principal Scientist, Process Development Services for Advanced Therapies, Sartorius
2:45 p.m.	3:15 p.m.	Coffee Break	
3:15 p.m.	4:00 p.m.	Autologous Immune Cell Therapy Manufacturing	Pierre Springuel , PhD, University College London
4:00 p.m.	4:45 p.m.	Academic Medical Cell Therapy Development, Manufacturing, and Delivery at Dana-Farber Cancer Institute	Julie Porter , Vice President, Cellular Therapies Operations, Dana-Farber Cancer Institute Sarah Nikiforow , Technical Director, DFCI Immune Effector Cell Therapy Program, Dana-Farber Cancer Institute
4:45 p.m.	5:00 p.m.	Short Break	
5:00 p.m.	5:30 p.m.	Closing Remarks & Feedback Session	Priyanka Gupta , Project Lead CGT Forum, Advocacy Team, Sartorius Maya Fürstenau-Sharp , Head of Marketing ATS, Sartorius
5:30 p.m.	7:30 p.m.	Networking Cocktail Reception	

SARTORIUS