

SARTORIUS

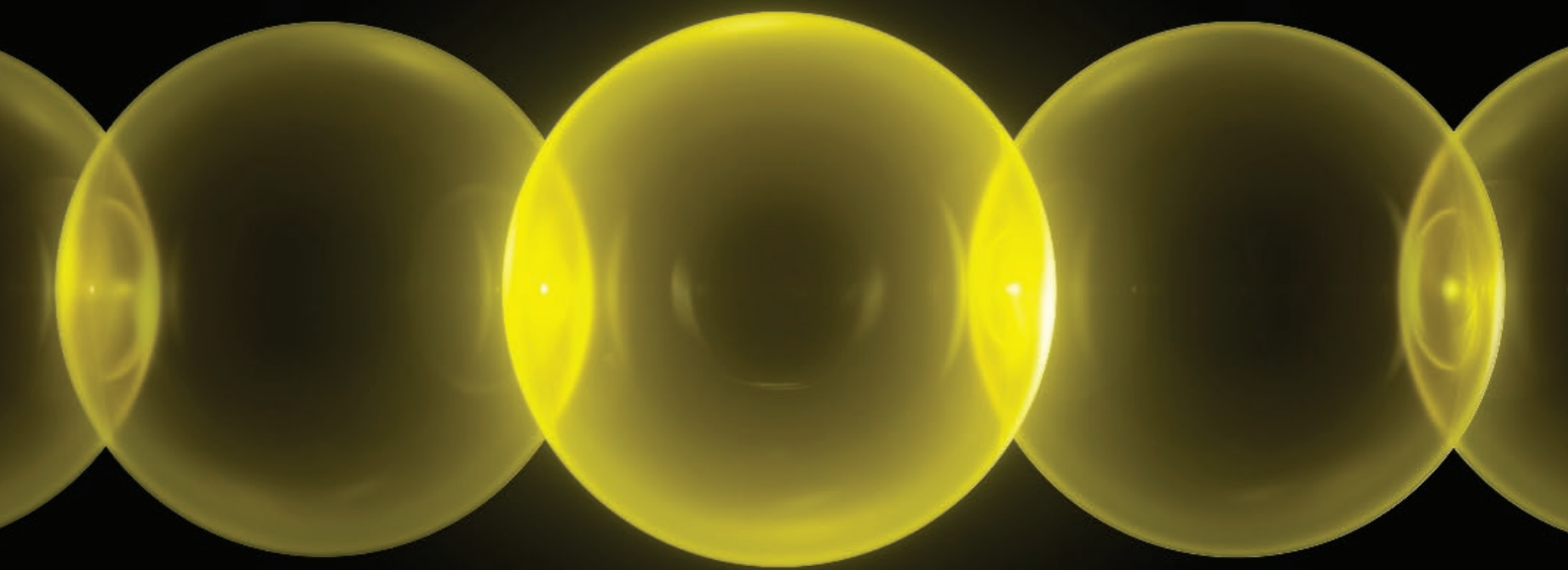
2025

Pharma Xchange Forum

Quality Testing and Characterization of Biologics

July 15

Vinpearl Landmark 81,
Ho Chi Minh City, Vietnam



Date:
July 15, 2025

Venue:
Vinpearl Landmark 81,
Ho Chi Minh City, Vietnam



Agenda | AM

Main Forum					
From	Until	Topic	Location	Presenter	
07:30	08:25	Registration / Check-in			
08:25	08:30	Welcome Address by Regional Leaders	Main Stage		
08:30	08:50	Opening Address Rise of Biologics	Main Stage	Eirik Pettersen	Head of Sales, LPS APAC
08:50	09:15	From Molecules to Biologics: The Principles of Octet	Main Stage	Paul Choi	Application Scientist, BioAnalytics
09:15	09:45	Assessing Critical Quality Attributes and Protein Quantitation	Main Stage	Ishan Subudhi Susheelendra Vaidya	Business Manager, BioAnalytics Technical Expert, BioAnalytics
09:45	10:00	Coffee Break			
10:00	10:40	Concept Of Minimum Sample Weight And Cost Efficiency	Main Stage	Eszter Paldy, PhD	Manager of Product Management Premium Weighing
10:40	11:20	Data Integrity for Sterility Testing	Main Stage	Glenn Manning, PhD	Regional Microbiology QC Expert
11:20	12:00	The Importance of 21 CFR in Final Lot Release Testing	Main Stage	Ms. Nguyễn Thị Thu Hường	Deputy Head, National Institute for Drug Quality Control
12:00	13:00	Networking Lunch			

Agenda | PM

Hands-on Workshop					
From	Until	Topic	Location	Trainer	
13:00	13:20	QC Workshop 1: Weighing Accuracy in Sample Preparation	Hands-on Workshop	Dylan Chan	Application Scientist, Lab Weighing & Lab Water
13:24	13:44	QC Workshop 2: Pre-Analytic Sample Preparation QC	Hands-on Workshop	Shuyun Lim	Application Scientist, Lab Filtration & Liquid Handling
13:47	14:07	QC Workshop 3: Microbiology QC and Sterility Testing	Hands-on Workshop	Glenn M., Randall C.	Application Scientist, Microbiology
14:10	14:30	QC Workshop 4: Bioanalytical QC – Assessing Critical Quality Attributes	Hands-on Workshop	Joanna Goh	Application Scientist, BioAnalytics
14:30	14:45	Coffee Break & Feedback			
Panel Discussion					
From	Until	Topic	Location	Presenter	
14:45	15:25	Expert Q&A: Ensuring Excellence: Innovations and Challenges in Pharma and Biologics QC	Main Stage		
15:25	15:30	Closing remark & Lucky draw			
15:30	16:00	Gift Redemption & End			

In partnership with:

