

SARTORIUS

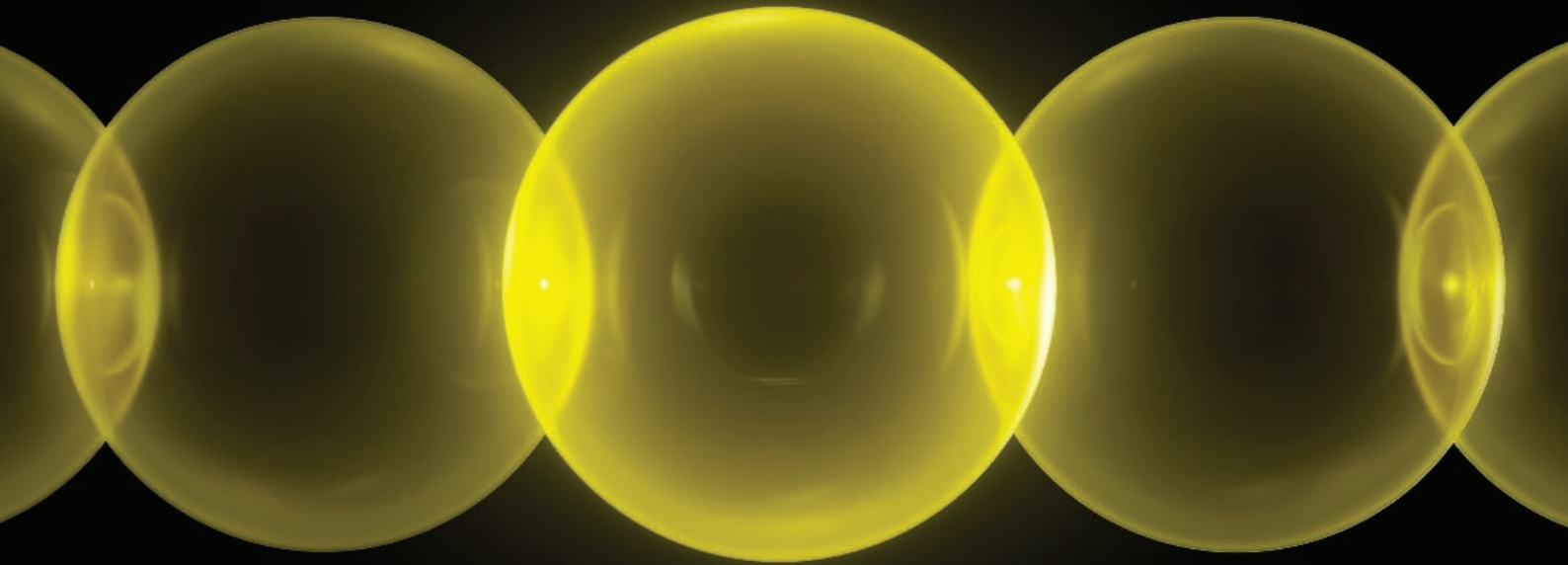
2025

Pharma Xchange Forum

Quality Testing and Characterization of Biologics

July 17

InterContinental Jakarta,
Indonesia



Date:
July 17, 2025

Venue:
InterContinental Jakarta,
Indonesia



Agenda | AM

Main Forum					
From	Until	Topic	Location	Presenter	
08:00	08:55	Registration / Check-in			
08:55	09:00	Welcome Address by Regional Leaders	Main Stage		
09:00	09:20	Opening Address Rise of Biologics	Main Stage	Eirik Pettersen	Head of Sales, LPS APAC
09:20	09:45	From Molecules to Biologics: The Principles of Octet	Main Stage	Paul Choi	Application Scientist, BioAnalytics
09:45	10:15	Assessing Critical Quality Attributes and Protein Quantitation	Main Stage	Ishan Subudhi Susheelendra Vaidya	Business Manager, BioAnalytics Technical Expert, BioAnalytics
10:20	10:30	Coffee Break			
10:30	11:10	Concept Of Minimum Sample Weight And Cost Efficiency	Main Stage	Eszter Paldy, PhD	Manager of Product Management Premium Weighing
11:10	11:50	Data Integrity for Sterility Testing	Main Stage	Glenn Manning, PhD	Regional Microbiology QC Expert
11:50	12:30	Practical Implementations of Data Integrity in the Pharmaceutical Industry	Main Stage	Apt. Arthur Henry, S.Si, MM	PT Elpharma Multi Solution
12:30	13:30	Networking Lunch			

Agenda | PM

Hands-on Workshop					
From	Until	Topic	Location	Trainer	
13:30	13:50	QC Workshop 1: Weighing Accuracy in Sample Preparation	Hands-on Workshop	Dylan Chan	Application Scientist, Lab Weighing & Lab Water
13:54	14:14	QC Workshop 2: Pre-Analytic Sample Preparation QC	Hands-on Workshop	Shuyun Lim	Application Scientist, Lab Filtration & Liquid Handling
14:17	14:37	QC Workshop 3: Microbiology QC and Sterility Testing	Hands-on Workshop	Glenn M., Randall C.	Application Scientist, Microbiology
14:40	15:00	QC Workshop 4: Bioanalytical QC - Assessing Critical Quality Attributes	Hands-on Workshop	Joanna Goh	Application Scientist, BioAnalytics
15:00	15:15	Coffee Break & Feedback			
Panel Discussion					
From	Until	Topic	Location	Presenter	
15:15	15:55	Expert Q&A: Ensuring Excellence: Innovations and Challenges in Pharma and Biologics QC	Main Stage		
15:55	16:00	Closing remark & Lucky Draw			
16:00	16:30	Gift Redemption & End			

In partnership with:

