

# Beyond the Barrier: Integrating Aseptic Containment Technology with Microbial Excellence

**Date:** 12 March 2026 (Thu)

**Time:** 09:30 - 15:30 SGT

**Venue:** Sartorius Singapore Pte. Ltd. 30 Pasir Panjang Road, #06-32, Mapletree Business City, Singapore 117440

**Featuring Experts from:** Sartorius | SKAN (supported by ITS Science & Medical Pte Ltd)



# Beyond the Barrier: Integrating Aseptic Containment Technology with Microbial Excellence

**Date:** 12 March 2026 (Thu)

**Time:** 09:30 - 15:30 SGT

**Venue:** Sartorius Singapore Pte. Ltd. 30 Pasir Panjang Road, #06-32,  
Mapletree Business City, Singapore 117440



In an era of tightening regulatory scrutiny, your facility's weakest link often lies where barrier systems meet monitoring methods. True contamination control requires more than a physical barrier—it demands a seamlessly validated ecosystem.

Join us for an exclusive deep dive into building a robust Contamination Control Strategy (CCS). This session bridges advanced barrier technology with applied microbiology to address real-world challenges in aseptic processing, environmental monitoring, and sterility testing.

## Program

- 09:30 Welcome & Check-in
- 10:00 Introduction and Application Center Tour
- 10:15 In Focus: *Continuous Microbial Air Monitoring*  
*Overview of regulatory requirements for continuous microbial air monitoring under EU GMP Annex 1 and EN 17141 including practical implementation using active air samplers and gelatine membrane filters.*
- 10:45 Break
- 11:00 In Focus: *SKAN Solutions*  
*How to achieve the most reliable sterility test results with Spectra, the latest version of SKAN isolator technology, especially developed for aseptic and aseptic-toxic applications.*
- 12:00 Lunch
- 13:00 In Focus: *Modernizing Sterility Testing: Innovation and Compliance*  
*Key aspects of sterility testing focusing on regulatory compliance, including 21 CFR Part 11 and EU GMP Annex 11, and best practices using sterility testing pumps and closed canister systems.*
- 13:45 Break
- 14:00 Hands-on Sessions by Sartorius & Skan
- 15:00 Panel Q&A
- 15:15 Closing Remark

Please be informed that we may take pictures and record videos during this event. Please notify us if you do not agree to have any images of you published in such pictures taken or video recordings made.

## Featuring Experts from: Sartorius & SKAN



**Eric Clement**  
Manager of Product Management, Reagent & Consumable Solutions | Sartorius



**Lim Shu Yun**  
Field Application Specialist, SEA & Taiwan | Sartorius



**Dr. Roland Madoerin**  
Head Sales Asia & Latin America | SKAN



**Frank Homberger**  
Regional Sales Director Asia | SKAN



**[Host] Randall Chan**  
Customer Advocacy Manager LE, Marketing Ops APAC | Sartorius

Save the Date:  
March 12, 2026 (Thu)