

## Unlock New Skills in Live-Cell Analysis

September 19 - 20, 2024 | Adelaide | 4 North Terrace | AHMS, Level 6

### Incucyte® Workshop Learn, Apply, & Excel!

Are you tired of having to take your cells out of the incubator to check on them?

Are you tired of missing those critical timepoints time and time again?



Ask new questions



Protect your cells



Improve productivity



Get new answers

Guest speakers include:



**Dr Agatha Labrindis**

AHMS Facility  
Manager  
The University of  
Adelaide



**Dr Nick Eyre**

Senior Research Fellow  
in Microscopy and  
Group Leader  
Flinders University



**Jenna Shean**

Product Manager,  
Incucyte  
Sartorius



Register now

This event is free of charge for you.  
Food and beverages will be provided.

Scan the QR code  
to reserve your spot  
in person or virtual



Questions? Contact us

Nabiha Elias, PhD  
Field Application Specialist, Incucyte  
Sartorius  
Phone: +61 487 051 263  
Email: [nabiha.elias@sartorius.com](mailto:nabiha.elias@sartorius.com)

Anthony Boghdadi, PhD  
Field Account Manager, BioAnalytics  
Sartorius  
Phone: +61 455 339 551  
Email: [anthony.boghdadi@sartorius.com](mailto:anthony.boghdadi@sartorius.com)