

ImageXpress High-Content Imaging User Group Meeting Agenda



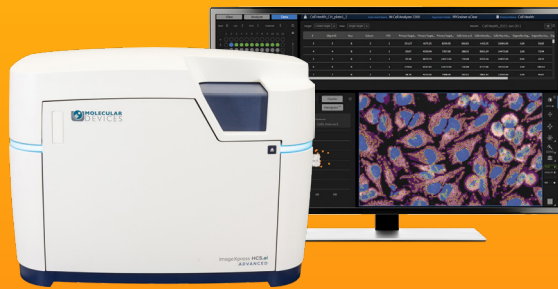
Thursday, October 23, 2025 | 9:30 am – 4:00 pm PDT

AC Hotel San Francisco Airport, Bayview Room, Floor 2,
1333 Veterans Blvd, South San Francisco, CA 94080

Lunch and light refreshments will be served

Time	Session	Presenters
9:30 – 10:00	Registration	
10:00 – 10:15	Introduction & Agenda	Lars Kristiansen Vice President, Product, Strategy & Innovation, Molecular Devices
10:15 – 10:45	Overcoming the challenges of live cell imaging of hiPSC-derived neurons using IN Carta Image Analysis Software	Mie Yamanaka, MD Postdoctoral Fellow, Gladstone Institute – Shinya Yamanaka Lab
10:45 – 11:15	Next-Gen Solutions: ImageXpress HCS.ai and Portfolio Innovations	Kayla Hill Sales Specialist, Molecular Devices
11:15 – 11:45	Advancing 3D in vitro CNS models: imaging assay development for high-throughput screening	Sneha Rao, PhD Senior Scientist, Arc Institute
11:45 – 12:30	Lunch	
12:30 – 1:30	Breakout Session	
1:30 – 1:45	High-content imaging to elucidate killing kinetics for patients with familial hemophagocytic lymphohistiocytosis (fHLH)	Allison Bien Research Technician, UC Berkeley at the Innovative Genomics Building
1:45 – 2:15	Multiplexed analysis to explore and refine human pluripotent stem cells differentiation trajectories in vitro	Yves Maury, PhD Director of Discovery, Neurona Therapeutics
2:15 – 3:00	High-Content Imaging Simplified: Introducing the new HCS.ai platform with AI-enabled IN Carta Image Analysis	Angeline Lim, PhD and Misha Bashkurov, PhD Product Managers, Cellular Imaging, Molecular Devices
3:00 – 3:30	Screening organoid and epithelial cells to read out maturity and functional states for drug development	Sarah Gierke, PhD and Max Tauc-Adrian, PhD Scientists, Genentech, Center for Advanced Light Microscopy and Electron Microscopy (CALM-EM)
3:30 – 4:00	Closing Remarks	Serena Cervantes Senior Field Application Scientist, Cellular Imaging, Molecular Devices

ImageXpress® HCS.ai
High-Content
Screening System



AI-powered **IN Carta®**
Analysis Software

Our speakers



Mie Yamanaka, MD
Postdoctoral Fellow
Gladstone Institute –
Shinya Yamanaka Lab



Sneha Rao, PhD
Senior Scientist
Arc institute



Allison Bien
Research Technician
UC Berkeley at the Innovative
Genomics Building



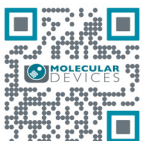
Yves Maury, PhD
Director of Discovery
Neurona Therapeutics



Sarah Gierke, PhD
Scientist
Genentech – the Center for
Advanced Light Microscopy and
Electron Microscopy (CALM-EM)



Max Tauc-Adrian, PhD
Scientist
Genentech – the Center for
Advanced Light Microscopy and
Electron Microscopy (CALM-EM)



For additional information or to register, please scan the QR code or visit:
<https://info.moleculardevices.com/dd/img/imaging-user-meeting-south-san-francisco-2025>

Contact Us

Phone: +1.800.635.5577
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current
listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+1.800.266.5338

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

