

SLAS 2024. TECAN'S AGENDA.



5TH FEBRUARY.

Spotlight presentation, by Mimetas (20 min.)

12:00-12:20 pm

Solutions Spotlight Theater (Exhibition 1290)

Membrane-free high-throughput gut-on-chip models for phenotypic evaluation of intestinal epithelial responses. Workflow powered by Tecan.

Poster presentation, by Mimetas

12:00-1:00 pm

Exhibit Halls AB Poster Session A

A high-throughput gut-on-a-chip model integrating adult stem cell-derived organoids for disease modeling and drug discovery. Workflow powered by Tecan.

Poster presentation, Dr. Lynda O'Leary

2:00-3:00 pm

Poster Assignment Number: 1405-B

Session: Poster Session B

Multiplex analysis of Predictive Oncology's 3D tumor models using Tecan's Spark® Cyto i maging multimode reader.

A standardized, high-throughput platform for automated, rapid, and extensive plasma proteome characterization (30 Min.)

4:30-5:00 pm

Room 210C

Dr. Manuel Bauer., Associate Director Mass Spectrometry, Tecan.



6TH FEBRUARY.

Standardization and automation for liquid biopsy solutions in oncology (90 Min.)

8:30-10:00 am

Room 104C

Lauren Leimann, Executive Director, BloodPAC; Alanna Church from Boston Children's Hospital/Harvard U and Alexandra Sommer, Senior Product Manager Clinical Diagnostics, Tecan.

SLAS 2024. TECAN'S AGENDA.



6TH FEBRUARY.

Innovative automation solutions for 3D cell biology, single-cell insights and proteomics workflows (90 Min.)

8:30-10:00 am

Room 151A

Rene Overmeer, Head of Assay development, and Automation at HUB Organoid, and Bogdan Budnik from Wyss Institute Boston, Alexander Schmidt, Head of Proteomics Core Facility, Biozentrum Basel and Lesley Schultz, Director Enterprise Cellomics-Proteomics, Tecan.

Lab of Today - Get the digital advantage now (90 Min.)

8:30-10:00 am

Room 153B

Marco Ravot-Licheri, Head of Digital Life Sciences Business, Tecan, Jason Meredith, Team Leader Product Manager Software, Tecan and Hal Wehrenberg, Vice President Digital Transformation, Tecan.

Poster presentation, Dr. Michael Brilhante

12:00 - 1:00 pm

Poster Assignment Number: 1126-C

Session: Poster Session C

The pioneering NETSseq platform leverages an automated RNA sequencing workflow to drive innovation in precision medicine against neurodegenerative diseases.

Poster presentation Dr. Michael Brilhante

2:00 - 3:00 pm

Poster Assignment Number: 1131-D

Session: Poster Session D

Automation Revolution: Streamlined Sample Collection and DNA Extraction for High-Throughput Microbiome Analysis.