

Automated Workflows: Connected with Confidence

SLAS2026 | Agilent Poster Presentations



Learn more about the many exciting ways Agilent can confidently connect you to automation solutions.

Poster Presentations:



Date: Tuesday February 10, 2026

Time: 12:00 PM - 1:00 PM

Poster #: 1317-C

High-Throughput Automation of mRNA Library Preparation for Reliable RNA-seq Workflows

Author: Lauren Alcoser, Field Application Scientist, Agilent Technologies

Co-authors: Cole Rohrbaugh, Jimmy Perrott, Thomas Wu, Eleanor Matrullo



Date: Tuesday February 10, 2026

Time: 2:00 PM - 3:00 PM

Poster #: 1318-D

High-Throughput Automation of Standard and Customized Cell Painting Assays: Integrated Liquid Handling and Imaging Workflow

Author: Joe Clayton, Technical Marketing Director, Agilent Technologies