

Expert-Led Training Calendar

Welcome to our Expert-Led Training Calendar, designed to elevate your skills and knowledge. Explore the upcoming sessions led by industry experts. **Training Calendar**

Product Line	Training title	Location	Date/Time	Details & Registration	Details
In Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, May 7, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
High Content Screening	HCS Advanced – Live Training	Hamburg Germany	Tues, May 13-15, 2025	Register Here	This class is full
Image Cytometry	Celigo™ Image Cytometer overview and operation training	Webinar via MS Teams	Tues, May 20, 2025 1-2 pm ET	Register Here 	The Celigo image cytometer is a plate-based high-throughput screening tool for whole-well imaging and advanced cell-based assays. Discover the diverse applications you can perform directly in a plate and gain hands-on experience with the instrument's operation. Additionally, we will host a Q&A session to address any queries related to our cell counting instruments
In Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, June 4, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
In Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, July 2, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
In Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Aug 6, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Sept 3, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Oct 1, 2025 8-10 am PT	Register Here	Click link to register

Expert-Led Training Calendar

Welcome to our Expert-Led Training Calendar, designed to elevate your skills and knowledge. Explore the upcoming sessions led by industry experts. **Training Calendar**

Product Line	Training title	Location	Date/Time	Details & Registration	Details
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Nov 5, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Dec 3, 2025 8-10 am PT	Register Here	Overview of hardware, software and imaging best practices
Maestro Advanced Training	Live Training	Live Training, Hamburg, Germany	Tues, Oct 14-16, 2025	Register Here	For users who want to independently set-up new (NGS) applications, alter existing (NGS) applications, optimize pipetting performance and perform basic maintenance tasks on the Sciclone G3 and Zephyr G3 (NGS) workstations.

How to Register

To explore detailed information and to register for any session, click on the respective link in the table above. This will direct you to our platform.

Note:

Training locations and dates are subject to change. Please refer to the schedule for the latest updates. For inquiries, contact us at customertraining@revvity.com