

# 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, April 2, 2025, 8-10 am PT	<a href="#">Register Here</a>	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, May 7, 2025, 8-10 am PT	<a href="#">Register Here</a>	Overview of hardware, software and imaging best practices
High Content Screening	HCS Advanced	Hamburg Germany	May 13-15	<a href="#">Register Here</a>	<b>This class is full</b>
In Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, June 4, 2025 8-10 am PT	<a href="#">Register Here</a>	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	<a href="#">Register Here</a>	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	<a href="#">Register Here</a>	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	<a href="#">Register Here</a>	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	<a href="#">Register Here</a>	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	<a href="#">Register Here</a>	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
High Content Screening	Maestro Advanced	Hamburg Germany	October 14-16	To be added	Class dates being finalized
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Nov 5, 2025 8-10 am PT	<a href="#">Register Here</a>	Click link to register
Vivo University – IVIS	Online training for Revvity's IVIS™ instrument users	Webinar via MS Teams	Wed, Dec 3, 2025 8-10 am PT	<a href="#">Register Here</a>	Click link to register

## 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](mailto:customertraining@revvity.com)