

Global Training Calendar 2024

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

To explore detailed information, scroll to training details section. To register for any session, click on the respective link in the table, this will direct you to our platform.

Training Calendar

Product Line	Training Title	Location	Dates - 2024	Registration	Details
In Vivo Imaging	In Vivo University	Hopkinton, MA - USA	March 19-21 June 25-27 Aug 20-22 Nov 19-21	Click to request	For complete details of the trainings, scroll to the section Training Details, below. If the class is full, contact us to add you to the list for the next session.
Next Generation Sequencing (NGS)	Maestro Advanced Training	Hamburg, Germany - Europe	March 19-21	Registration	
Automated Workstations	Janus G3 User Training	Hamburg, Germany - Europe	March 19-21	Registration	
High Content Screening	Revvity High Content Basic Training	Hamburg, Germany - Europe	April 9-11	Registration	
High Content Screening	High Content Advanced Training	Hamburg, Germany - Europe	May 14-16	Registration	

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

Global Training Calendar - 2024

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

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.

Training Details

In Vivo University

Description: This live classroom training is conducted by our In Vivo Imaging Scientist and will include guest speakers from Center's Applied Biology, R&D and Reagents teams. The topics will range from practical aspects of in Vivo imaging to applied techniques, specific methodologies and optimization of instruments, reagents and software.

Location: Revvity Hopkinton, MA - USA

Dates: March 19-21, June 25-27, August 20-22, Nov. 19-21

[Link to register](#)

Maestro Advanced Training

Description: 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.

Dates: March 19-21

Location: Revvity Cellular Technologies GmbH – Hamburg, Germany

[Link to register](#)

Janus G3 User Training

Description: Users who want to independently set up new automated applications, optimize pipetting performance, and perform basic maintenance tasks on the JANUS G3 Liquid Handler.

Location: Revvity Cellular Technologies GmbH – Hamburg, Germany

Dates: March 19-21

[Link to register](#)

Revvity High Content Basic Training

Description: You will learn the principles of High Content Screening. After attending the training, you will be able to run the system and create new image acquisition protocols. For the image analysis, you will learn the key principles behind the Harmony™ software, how to optimize existing algorithms and design your own algorithm using the Harmony™ building blocks.

Location: Revvity Cellular Technologies GmbH – Hamburg, Germany

Dates: April 9-14

[Link to register](#)

Revvity High Content Advanced Training

Description: You will learn the principles of High Content Screening. After attending the training, you will be able to run the system and create new image acquisition protocols. For the image analysis, you will learn the key principles behind the Harmony™ software, how to optimize existing algorithms and design your own algorithm using the Harmony™ building blocks.

Location: Revvity Cellular Technologies GmbH – Hamburg, Germany

Dates: May 14-16

[Link to register](#)

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