



# Flexible training options for increased productivity.

## Revvity radiometric detection training

A respected leader in the global scientific community, Revvity has been delivering top-quality products and services for more than 70 years. We continue that legacy today with a range of radiometric detection training offerings that deliver the necessary level of training to ensure your success through familiarization with your instrument and beyond.



Tri-Carb/Quantulus GCT



Wizard<sup>2</sup>



MicroBeta<sup>2</sup>

## Basic familiarization training

Basic instrument training occurs at the time of installation and includes performance testing and customer familiarization of key features and functions as well as safety measures. The training is based on Standard Operating Procedures and is performed by one of Revvity's certified customer service engineers.

## Web-based user training (*online*)

Links to our comprehensive, web-based training courses are provided by a radiometric detection specialist following installation of your radiometric detection instrument. Revvity's courses include interactive content for the following products:

Tri-Carb®, Wizard<sup>2</sup>®, Quantulus™, MicroBeta<sup>2</sup>®.

Users can register with their sales representative by providing their contact details, instrument model and serial number.

**Fee listed on Instrument Quote PN-7009002**

## Rad detection remote support (*fee per hour*)

Revvity offers remote customer training for those who prefer to have a product specialist deliver training remotely online. This support can include basic detection theory, instrument operation, protocol setup instruction, or \*remote protocol programming.

*\*(Instrument with direct internet access required.)*

**Fee for Service PN-N0236514**

## 1-Day on-site new instrument operational training

Revvity offers on-site customer training for those customers who prefer to have a product specialist come to their site. This advanced customer training program includes basic detection theory, instrument operation and protocol setup instruction. On-site training is individually scheduled following purchase.

**Fee for Service PN-7009003**

The Revvity logo is displayed in a lowercase, sans-serif font. It is positioned in the bottom right corner of the page, above a large yellow wavy graphic that spans the width of the page.