



# Transform your research with advanced radiometric detection.

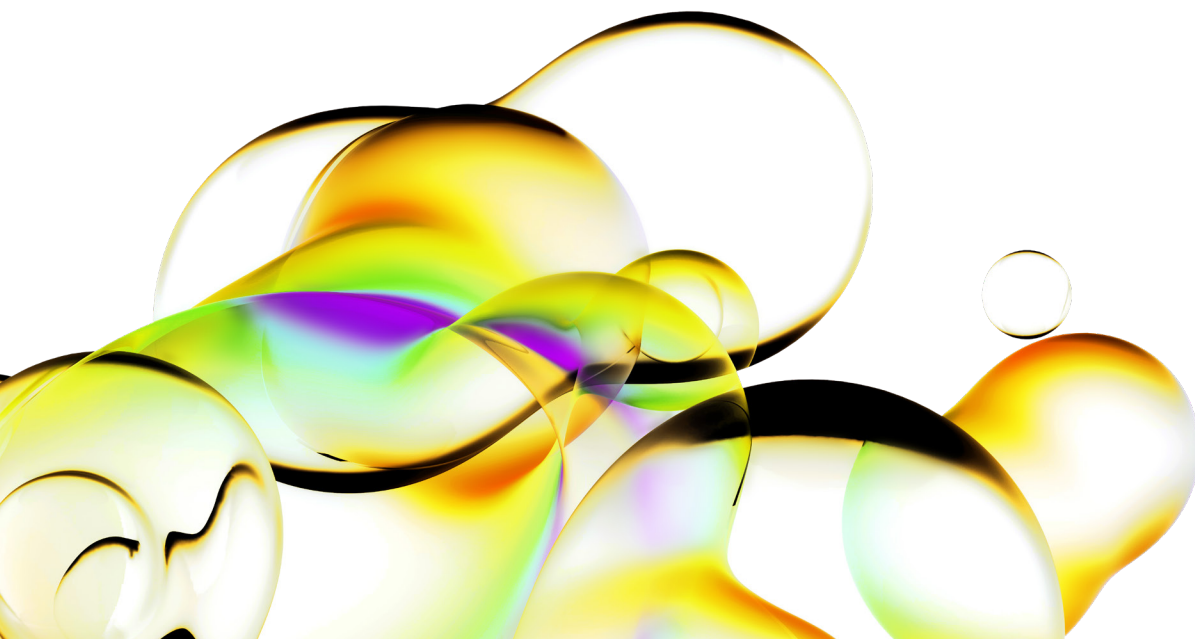
## Instrument trade-in program

Whether you're working in environmental research, industrial applications, radio-carbon, or drug discovery and development, now is the perfect time to upgrade your lab.

Enhance your precision, efficiency, and data quality with the latest technology. Trade in your old instruments and step up to Revvity's Tri-Carb™ and Quantulus™ liquid scintillation counters, MicroBeta<sup>2</sup>™ microplate counters, sample oxidizers, or Wizard<sup>2</sup>™ gamma counters.



[Learn more](#)



Contact your sales representative to learn more about Revvity's radiometric research solutions.

### Tri-Carb™ and Quantulus™ liquid scintillation counters

Ideal for everything from basic research to advanced, low-level environmental analysis. Able to exceed requirements for your most demanding applications, backed by our 60 years of experience and innovation.

[Learn more](#)



### MicroBeta2™ microplate counter

Provides advanced radiometric and luminescent detection abilities for GPCRs, kinases, and reporter gene assays. Delivers superior sensitivity and performance along with flexibility.

[Learn more](#)



### Sample Oxidizer

Utilizes the patented "open/close" flame oxidation technology to expand your experimental designs in liquid scintillation counting.

[Learn more](#)



### Wizard2™ gamma counter

Offers outstanding sensitivity together with high-level security, connectivity, and ease of use, giving you greater command over your gamma counting process.

[Learn more](#)



revvity