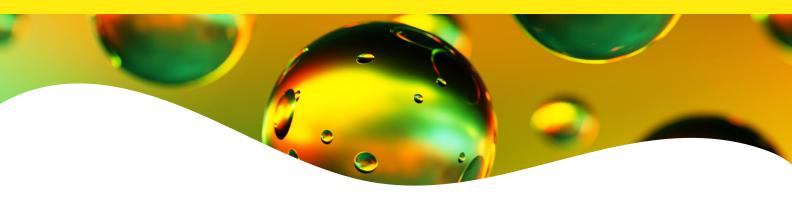
# revvity

# Radiometric detection systems and reagents.



For highest sensitivity and accurate, quantitative results, radiometric detection is the gold standard. Our radiometric detection systems are designed around your application needs related to gamma counting, beta counting, and other types of radiochemical detection and measurement.

## Tri-Carb liquid scintillation counters

The Tri-Carb® benchtop vial-based liquid scintillation analyzers have a strong reputation for high performance liquid scintillation counting with sensitive detectors for measuring small amounts of beta, gamma, and alpha radioactivity. The Tri-Carb's superior reliability meets the requirements of even the most demanding applications, including basic academic and pharmaceutical research, ADME/Tox research, as well as environmental, nuclear power applications, biofuel, and other bio-based material analyses.



# Quantulus GCT liquid scintillation spectrometer

The Quantulus GCT™ 6220 Liquid Scintillation Counter is a fully loaded benchtop instrument with unsurpassed performance for measuring man-made, cosmic, and other natural radionuclides. Patented GCT technology combined with BGO guard background reduction accurately measures to near-background levels. Proprietary Dual PSA Discriminator and PSA Histogram improves accuracy for mixed and unknown samples.



# WIZARD<sup>2</sup> automatic gamma counter

The WIZARD<sup>2®</sup> automatic gamma counters are designed for superior counting performance with all types of samples for every gamma counting application. The newest generation of gamma counters offer significant hardware and software enhancements. Unique well-type detectors, sample changer system advanced robotics, and highly effective lead shielding result in high counting efficient with constant background and minimal crosstalk.



# MicroBeta<sup>2</sup> plate based liquid scintillation and luminescence counters

The MicroBeta² is designed for researchers engaged in all major radiometric and luminescence applications. By combining LSC reliability with a plate reader's simplicity, it allows vast savings in time, consumables, and waste. The MicroBeta² is ideal for assays using ³H ¹⁴C, ³³P, ³⁵S, ⁵¹Cr, ¹²⁵I, and glow-type luminescence labels as well as for ScintiPlate, FlashPlate, and other scintillation proximity assays. It counts filters and tubes in addition to microtitration plates.



# Model A307 sample oxidizer

The 307 Sample Oxidizer is an automatic preparation and oxidization system for both single and dual radiolabeled samples containing 3H and/or 14C for use in liquid scintillation counting. The patented "open/close" flame oxidation technology will expand your experimental designs in liquid scintillation counting. Reliable combustion of biological, environmental, and industrial sample material increases accuracy in liquid scintillation counting due to no self-absorption, no chemiluminescence, no color quench, no spectral overlap, and no variable chemical quench.



#### Cell harvesters

The FilterMate™ Universal Harvester allows for the simple, simultaneous harvesting of samples from 96-and 24-well microplates. This compact, versatile system harvests 2 or 96 samples simultaneously onto solid filter supports or into microplates. Speed and simplicity are the primary strengths of the Filtermate Harvester.

#### Radiochemicals

Revvity has been a leading supplier of radiochemicals for over 60 years. We're on the cutting edge when it comes to safe and appropriate packaging and handling to ensure confidence and consistent performance in your radiometric applications. Our radioactivity portfolio includes radioactive nucleotides, radiolabeled ligands, radiotherapeutic nuclides, radionuclides, a variety of steroids, carbohydrates, cyclic nucleotides, sugars and amino acids, CAT assay reagents, lipids, and microspheres, along with custom services for specialized requirements.

## Liquid scintillation cocktails

Only the right cocktail choice, paired with a correct sample preparation procedure and instrument, will result in accurate and reproducible counting results. We offer high performance cocktails with high counting efficiency, high quench resistance, very high flash point for simple transportation, no special storage requirements, very low vapor pressure, with low toxicity.

# Liquid scintillation counting vials

Revvity offers high quality glass and plastic vials. The best vial to choose is dependent on the type and volume of sample to be counted and the cocktail that will be used. Glass vials are manufactured from low potassium glass tubing. The uniform wall thickness contributes to excellent counting reproducibility, and plastic vials are injection (blow) molded to exacting specifications from virgin high-density (linear) polyethylene (HDPE) with lower background level and higher counting efficiency.

www.revvity.com 2

### Microplates

A microplate is not just a microplate. It's important that researchers choose the correct microplate for the correct application. The correct selection of a microplate can help improve signal, reduce background, provide higher image clarity, and provide for more relevant biological models in a higher throughput format. Revvity microplates can help you achieve optimal results in your radiometric assays.

# Scintillation proximity assay (SPA)

Combine the sensitivity of radioactive assays with the convenience of a homogeneous (no wash steps) assay using SPA technology. Scintillation proximity assay (SPA) is a homogeneous and versatile technology for the rapid and sensitive assay of a wide range of biological processes including use with enzyme targets and receptor targets, radioimmunoassays, and molecular interactions. The assay format requires no separation steps and is amenable to automation.



