



revvity

# Scintillating results at your lowest concentrations.

## Newly redesigned to make the best even better

For over 60 years Revvity has been providing unmatched versatility and sensitivity in radiometric detection. That's why research and environmental labs around the world recognize the top brands for their liquid scintillation application needs — our Tri-Carb® and Quantulus™ GCT families of liquid scintillation counters. Tri-Carb LSC systems are newly redesigned with enhanced features and optional PSA (Pulse Shape Analysis) Alpha and Beta discrimination counting mode technology. And the new Quantulus GCT, with patent-pending technology, delivers the ultralow-level sensitivity you trust, but in a lighter benchtop footprint that fits any lab.

The Quantulus GCT 6220 has extremely high sensitivity, and unmatched detection of low-level Alpha and Beta radioactivity. The Bismuth Germanium Oxide guard works in conjunction with patent-pending Guard Compensation Technology (GCT) to further lower the instrument background, enhancing sensitivity to accurately measure near-background sample activity. Now 62% smaller and 76% lighter, the Quantulus GCT allows flexibility and mobility in the lab.



### New instrument features:

- Modern easy-to-use Windows® 8.1 based software
- Redesigned ergonomic arm - enter data either sitting or standing, or fold away saving space
- Instrument operational status indicator LED for clear assay status at a glance

- The replay assay reprocessing allows you to change parameters without re-running the assay
- Greater throughput with 408 standard 20 mL vials, or 720 small 4 mL or 7 mL vials capability
- Three spectrum memories for storage of alpha, beta, and cosmic or environmental gamma events simultaneously for optimized resolution



Quantulus GCT 6220 Liquid Scintillation Counter on portable bench top cart

The new Tri-Carb portfolio of liquid scintillation counters have been redesigned with optional PSA (Pulse Shape Analysis) to provide better particle discrimination for the detection of Alpha, Beta.

- The Tri-Carb live spectral display and plotting and chemiluminescence detection with optional correction allows for flexibility in your applications
- Single photon counting for bioluminescence assays
- Custom printing and electronic report output for versatility in reporting
- Color-corrected single- and dual-labeled DPM; and optional 60 user protocols with unlimited assays to meet your throughput needs



Tri-Carb Liquid Scintillation Counter

## Tri-Carb and Quantulus GCT system configurations at a glance

Feature	Tri-Carb 4810TR	Tri-Carb 4910TR	Tri-Carb 5110TR	Quantulus GCT 6220
Built-in computer with Windows® 8.1 operating system, 4 GB RAM (minimum), and 250 GB hard disk (minimum), three high-speed USB ports, and Dual Gigabit Ethernet support	S	S	S	S
19-inch-wide LCD monitor with DVI-D video output, keyboard, and ergonomic arm	S	S	S	S
QuantaSmart Software for Windows® 8.1 operating system	S	S	S	S
LED System Status Indicator	S	S	S	S
TR-LSC background electronics	S	S	S	S
EasyView Spectrum Analysis	O	O	O	S
Spectraworks™ Software	O	O	O	S
Enhanced Security (21 CFR Part 11 Compatibility)	NA	O	O	O
Patent-pending Guard Compensation Technology	NA	NA	NA	S
BGO detector guard	NA	NA	NA	S
High-Sensitivity Count Mode (HSCM)	O	O	S	NA
Ultralow-level Count Mode (ULLCM)	NA	O	O	S
Alpha/Beta discrimination (PSA)	NA	O	O	S
Pulse amplitude comparison (PAC) standard when purchased with LLCM	NA	O	O	S
Sample numeric data screening	S	S	S	S
Direct DPM	S	S	S	S
Single/dual color-corrected DPM	O	S	S	S
Triple Label DPM	O	O	S	S
Enhanced IPA (Instrument Performance Assessment)	NA	O	S	S
Live spectral display and plotting	S	S	S	S
Group PrioStat automatic priority interrupt	S	S	S	S
Sample PrioStat manual special function interrupt	NA	O	S	S
Luminescence correction	O	O	S	S
Single photon counting for bioluminescence assays	S	S	S	S
Enhanced Replay sample recall and reprocessing	O	S	S	S
Varisette sample changer	O	S	S	S
Number of user acquisition protocols with unlimited assays	15, 60, optional	30, 60, optional	60	60
2D Barcode reader	O	O	O	O
Worklist	O	O	S	S
Temperature Control	O	O	O	S

S = Standard O = Optional NA = Not Available

Contact your local representative at 1-888-723-7946 for a quote or visit [www.revivity.com](http://www.revivity.com) for more information.

