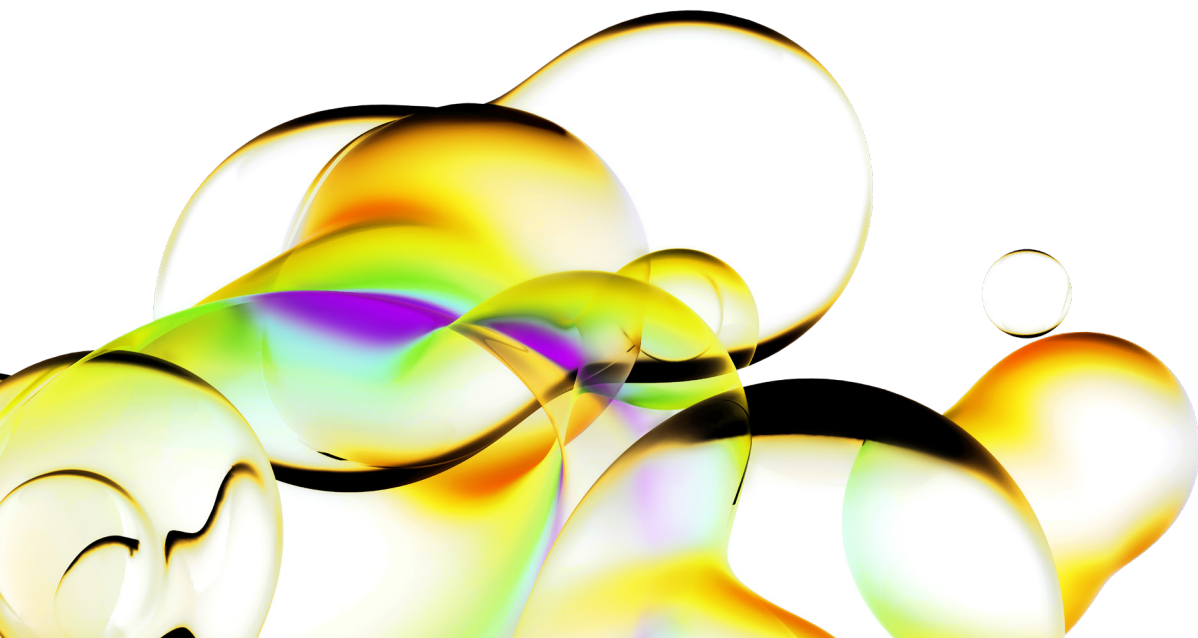




Delivering superior low-level detection.

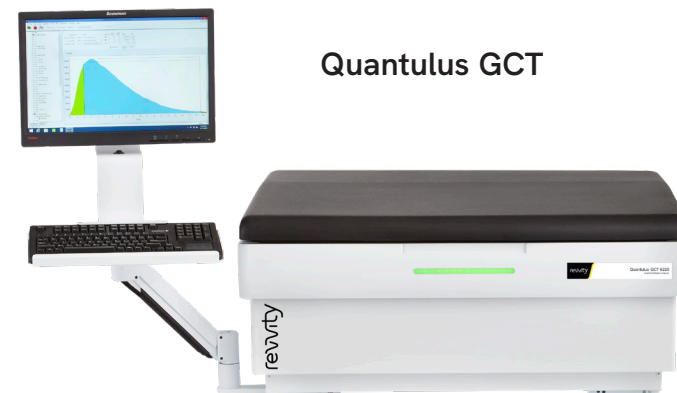


The Quantulus™ GCT is an advanced liquid scintillation counter for multiuser labs that perform environmental and routine counting for research applications. Especially ideal for environmental applications when you need to detect extremely low-level Alpha and Beta radioactivity. It offers unsurpassed performance for measuring man made, cosmic and other natural radionuclides.

While the Quantulus GCT is particularly well suited for the detection of extremely low-level Alpha and Beta applications, its high sample capacity allows a greater throughput which means more flexibility to address an increased number of demanding liquid scintillation counting applications.

Key features:

- **Patented bismuth germanium oxide (bgo) and guard compensation technology (GCT)** enhances sensitivity to accurately measure near-background sample activity.
- **QuantaSmart™ software** offers robust multitasking, easy networking environment with unlimited assays in a secure multiuser setting.
- **SpectraWorks 2** provides spectral analysis that determines optimal counting regions and calculates figure of merit automatically.



Applications

Environmental



TENORM-Technologically enhanced naturally occurring radioactive material identification & mapping

Industrial



From workers urine routine tests to radiation waste screening to district contamination monitoring

Drinking water



Determine gross alpha / gross beta, tritium, radon as well as other radionuclides in water for human use

Food



Strontium (^{90}Sr) in milk, synthetic materials (^{14}C), radio-caesium (^{134}Cs , ^{137}Cs)

Radio-carbon



Quantifying bio fractions in fossil mixes (biofuels, plastics, food, spirits, waste)

Drug development



The use of radiochemicals in drug development is still considered as the gold standard screening tool

For more information, visit www.revivity.com

