

TriCarb[®] and Quantulus[™] GCT liquid scintillation analyzers.

TriCarb & Quantulus [™] GCT system configurations	TriCarb 4810TR	TriCarb 4910TR	TriCarb 5110TR	Quantulus [™] GCT 6220
Feature				
Built-in computer with Windows 8.1 operating system 4 GB (minimum) RAM and 250 GB hard disk (minimum) 3 high speed USB ports and Dual Gigabit Ethernet support NEW	S	S	S	S
19 inch Wide LCD monitor with DVI-D video output and ergonomic keyboard NEW	S	S	S	S
QuantaSmart Software for Windows [®] 8.1 operating system	S	S	S	S New
LED System Status Indicator NEW	S	S	S	S
TR-LSC background electronics	S	S	S	S
EasyView Spectrum Analysis	O	O	O	S NEW
Spectraworks Software	O	O	O	S
Enhanced Security (21 CFR Part 11 Compatibility)	NA	O	O	O
Patent Pending Guard Compensation Technology NEW	NA	NA	NA	S
BGO detector guard	NA	NA	NA	S
HSCM (High Sensitivity Count Mode)	O	O	S	NA
ULLCM (Ultra Low Level Count Mode)	NA	O	O	S
Alpha/Beta discrimination (PSA),	NA	O New	O New	S
Pulse amplitude comparison (PAC) Std. when purchased with LLCM	NA	O New	O New	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

TriCarb & Quantulus™ GCT system configurations	TriCarb 4810TR	TriCarb 4910TR	TriCarb 5110TR	Quantulus™ GCT 6220
Feature				
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

