

Guava® easyCyte specifications

Guava® easyCyte HT Systems



Guava® easyCyte HT System	5HT	5HT HPL	6HT-2L	8HT	11HT	12HT	BG HT	BGR HT	BGV HT
Catalog No.	0500-4005	0500-4009	0500-4007	0500-4008	0500-4020	0500-4012	0500-4015	0500-4025	0500-4030
Violet (405 nm) Laser					•	•			•
Blue (488 nm) 50 mW Laser	•		•						
Blue (488 nm) 150 mW Laser		•		•	•	•	•	•	•
Green (532 nm) Laser							•	•	•
Red (642 nm) Laser			•	•		•		•	
FSC	•	•	•	•	•	•	•	•	•
SSC	•	•	•	•	•	•	•	•	•
Blue-V (450/45 nm)					•	•			•
Green-V (512/18 nm)									•
Green-V (525/30 nm)					•	•			
Yellow-V (575/25 nm)									•
Yellow-V (583/26 nm)					•	•			
Orange-V (620/52 nm)									•
Red-V (695/50 nm)					•	•			
NIR-V (785/70 nm)					•	•			
Green-B (512/18 nm)							•	•	•
Green-B (525/30 nm)	•	•	•	•	•	•			
Yellow-B (575/25 nm)							•	•	•
Yellow-B (583/26 nm)	•	•	•	•	•	•			
Red-B (695/50 nm)	•	•	•	•	•	•	•	•	•
NIR-B (785/70 nm)				•	•	•	•	•	•
Yellow-G (575/25 nm)							•	•	•
Orange-G (609/30 nm)								•	
Orange-G (620/52 nm)							•		•
Red-G (695/50 nm)							•	•	•
NIR-G (785/70 nm)							•	•	•
Red-R (664/20 nm)			•	•		•		•	
NIR-R (785/70 nm)				•		•		•	
Microcapillary Fluidics	•	•	•	•	•	•	•	•	•
Direct, Absolute Cell Counts	•	•	•	•	•	•	•	•	•
Automation-plate and tubes	•	•	•	•	•	•	•	•	•
Mixing	•	•	•	•	•	•	•	•	•
Dell® Laptop	•	•	•	•	•	•	•	•	•
InCyte™ Software	•	•	•	•	•	•	•	•	•
Digital Signal Processing	•	•	•	•	•	•	•	•	•

Guava® easyCyte Single Sample Systems



Guava® easyCyte Single Sample System	5	5 HPL	6-2L	8
Catalog No.	0500-5005	0500-5009	0500-5007	0500-5008
Violet (405 nm) Laser				
Blue (488 nm) 50 mW Laser	•		•	
Blue (488 nm) 150 mW Laser		•		•
Green (532 nm) Laser				
Red (642 nm) Laser			•	•
FSC	•	•	•	•
SSC	•	•	•	•
Blue-V (450/45 nm)				
Green-V (512/18 nm)				
Green-V (525/30 nm)				
Yellow-V (575/25 nm)				
Yellow-V (583/26 nm)				
Orange-V (620/52 nm)				
Red-V (695/50 nm)				
NIR-V (785/70 nm)				
Green-B (512/18 nm)				
Green-B (525/30 nm)	•	•	•	•
Yellow-B (575/25 nm)				
Yellow-B (583/26 nm)	•	•	•	•
Red-B (695/50 nm)	•	•	•	•
NIR-B (785/70 nm)				•
Yellow-G (575/25 nm)				
Orange-G (609/30 nm)				
Orange-G (620/52 nm)				
Red-G (695/50 nm)				
NIR-G (785/70 nm)				
Red-R (664/20 nm)			•	•
NIR-R (785/70 nm)				•
Microcapillary Fluidics	•	•	•	•
Direct, Absolute Cell Counts	•	•	•	•
Automation-plate and tubes				
Mixing				
Dell® Laptop	•	•	•	•
InCyte™ Software	•	•	•	•
Digital Signal Processing	•	•	•	•

Feature	Description
Dynamic Range	5 decades
Instrument Size	51 cm x 25 cm x 59 cm
Weight	67 lbs., 31 Kg
Absolute Cell Count Capability	Yes, without reference beads
Particle Size	0.2- 60 μ m
Fluidics System	Positive displacement pump
Sample Format	0.5 mL and 1.5 mL tubes and 96-well plate
Minimum Sample Volume	100 μ L for wells, 150 μ L for 0.5 mL tubes and 900 μ L for 1.5 mL tubes
Fluorescence Precision	< 4% CV for CEN
Acquisition Speed	10-1,000 events/second
Operating Temperature	16°-30°C; 80% humidity
Computer Specifications	Dell™ laptop running Windows® 7 Ultimate (32-bit only), and including Microsoft Excel. Minimum configuration: Intel Core i5-3210M processor (2.5 GHz, 3M cache); 4 GB, DDR3-1600 MHz SDRAM; 320 GB hard drive; 2 USB ports
Operating System	Windows® 7 or Mac OSX (analysis only)
Software	guavaSoft™ 3.0
Light Scatter Detection	Forward (0°, +/-9) Side (90°, +/-30)
Flow Rate	7-70 μ L/min
Sheath Fluid	None
Waste	50 mL at 8 hours of continuous use
Warranty	1 year