

SX-8G IP-Star Automated System General Specifications

IP-Star Automated System Specifications	
Magnetic Separation Device	<ul style="list-style-type: none"> • Dispenser instrument with 8 nozzles • Nozzle Pitch: 9 mm • Magnet Strength: 3000 Gauss
Plate locations	<ul style="list-style-type: none"> • 6 modules for 96-well plates
Pipette tip locations	<ul style="list-style-type: none"> • 4x 96 tip modules and 3x 24 tip modules
Compatible Consumables	<ul style="list-style-type: none"> • 96-well plates with round bottom. (Diagenode) • Polypropylene filter tips (Diagenode) • 12-strip tubes/ 8-strip tubes and caps (Diagenode)
Validated Kits	<ul style="list-style-type: none"> • Automated ChIP • Automated MeDIP • Automated MethylCap
Pipetting Volumes	<ul style="list-style-type: none"> • 5-200 μl
Pipetting Precision (I)	<ul style="list-style-type: none"> • 5-9 μl: \leftarrow 20% • 10-20μl: \leftarrow 5% • 21-200 μl: \leftarrow 2%
Temperature Blocks	<ul style="list-style-type: none"> • Two 96 well Peltier blocks for heating and cooling
Block Ramp Rate	<ul style="list-style-type: none"> • Heating (4°C-95°C): +0.12°C/sec • Cooling (95°C-4°C): -0.08°C/sec
Block Temperature Range	<ul style="list-style-type: none"> • 4°C-95°C
Temperature Accuracy	<ul style="list-style-type: none"> • +/- 2°C (above this values a system error occurs)
Temperature Uniformity Across Block	<ul style="list-style-type: none"> • At 4°C (+/- 0.6) • At 95°C (+/- 1.9)
Peltier's longevity	<ul style="list-style-type: none"> • 10000 to 12000 runs.
Cooling lower housing parts (boards and drives)	<ul style="list-style-type: none"> • Air inlet on both sides through dust filters
Air Flow on stage	<ul style="list-style-type: none"> • Air outlet by two fans
Decontamination Capability	<ul style="list-style-type: none"> • UV lamp controlled via PC-software • Safety switch off when front door is open
Drop catcher	<ul style="list-style-type: none"> • Liquid drop catcher for all 8 nozzles • Located below tips during travel of nozzle-head
Tip waste slide	<ul style="list-style-type: none"> • Automated discard location (waste chute) • Stainless steel, autoclavable and cleaning with ethanol
Tip waste bag	<ul style="list-style-type: none"> • Plastic bag, transparent, replaced in customers responsibility
Electrical Requirements	<ul style="list-style-type: none"> • Option 1: Voltage: 200 - 240 VAC \pm 10% • Option 2: Voltage: 100 - 120 VAC \pm 10% • Frequency: 50-60 Hz \pm 1% • Maximum Current Draw: xA
Overvoltage category	<ul style="list-style-type: none"> • II

Computer System	<ul style="list-style-type: none"> • Processor: AMD ATHLON 5400B (2.8GHz, 512KBX2) [Included in Price] • Operating System: Genuine Windows Vista Business Bonus-Windows XP Professional downgrade • Memory: 1.0GB DDR2 Non-ECC SDRAM, 800MHz, (1DIMM) [Included in Price] • Hard Drive: 80GB SATA, 7200 RPM Hard Drive with Data Burst Cache. [Included in Price] • Monitor: 1024 x 768 monitor (65535) (Size – 14") • Slot: optional riser-card cage Slot
Software	<ul style="list-style-type: none"> • SX8-8GV52
Instrument Physical Specifications	
Width	<ul style="list-style-type: none"> • 1070 mm (42.5 inches)
Height	<ul style="list-style-type: none"> • 780 mm (31 inches)
Depth	<ul style="list-style-type: none"> • 650 mm (26 inches)
Weight	<ul style="list-style-type: none"> • Approximately 130 kg (300 lbs)
General Instrument Design	<ul style="list-style-type: none"> • Benchtop style for standard lab tables • Front door transparent for a good view inside the instrument • Fully covered housing, impervious to UV-light from inside instrument and from outside lab • Good access when front door is opened for sample loading and instrument cleaning
Instrument Environmental Requirements	
Room Temperature	<ul style="list-style-type: none"> • 18-22°C (64.4°F to 71.6°F)
Recommended Humidity	<ul style="list-style-type: none"> • 30 - 50%
Maximum Humidity	<ul style="list-style-type: none"> • 80% humidity at 30°C
Temperature Storage	<ul style="list-style-type: none"> • -40°C to 65°C (-40°F to 149°F)
Storage Place	<ul style="list-style-type: none"> • For indoor use only
Maximum Altitude	<ul style="list-style-type: none"> • 2000 m
Pollution Level	<ul style="list-style-type: none"> • 2 (Normally only non-conductive pollution occurs)