



Maxwell[®] CSC System for Clinical Labs

Gain Efficiency & Confidence

Maxwell CSC



cGMP-Compliant Clinical Sample Concentrator

Worry-Free Nucleic Acid Extraction



Manufactured Under cGMP: Current Good Manufacturing Practices

The new Maxwell® Clinical Sample Concentrator automated nucleic acid extraction system is specifically designed for clinical laboratories needing to process multiple samples for critical downstream assays.

Increase Confidence in Results

Manufactured to a quality standard that assures consistent performance for every run.

Rely on Quality

Each system is backed by the Promega tradition of quality, service and support.

Maintain Electronic Records Easily

Meets 21 CFR Part 11 guidelines for electronic records.

Extraction Report

Protocol: Blood v0.9 Status: Completed Started at: 7/23/2012 2:14:41 PM Duration: 00:27:38 Operator: Promega-Maxwell/Promega	Software: Maxwell® CSC Software Version: 1.8 Instrument Name: PromegaInstrument Serial Number: beta unit 12 Firmware Version: 3.29.0 Maintenance Due: Yes	Catalog Number: n1121 Use Confirmations: Sample pre-processing is complete as per the Blood Technical Manual Deck tray has been prepared as follows and placed in the instrument: • Cartridges with seal completely removed placed at positions 1-16 • Elution tubes placed at positions 1-16 • Appropriate volume of elution reagent has been added to each elution • Plunger has been placed in well 8 of each cartridge Plungers are not present on instrument plunger bar
--	--	---

#	Sample ID	Lot Number	Exp. Month	Cartridge ID	Elution Tube ID	Custom1	Custom2
1	prototype 1	123456	12/2012				
2	production 106	123456	12/2012				
3	ps1 114	123456	12/2012				
4	ps1 116	123456	12/2012				
5	prototype 2	123456	12/2012				
6	production 63	123456	12/2012				
7	ps1 115	123456	12/2012				
8	ps1 106	123456	12/2012				
9	prototype 3	123456	12/2012				
10	production 93	123456	12/2012				
11	ps1 164	123456	12/2012				
12	ps1 101	123456	12/2012				
13	prototype 4	123456	12/2012				
14	production 55	123456	12/2012				
15	ps1 109	123456	12/2012				
16	ps1 111	123456	12/2012				

System Report

Protocol: SelfTest v0.9 Status: Completed Started at: 7/23/2012 8:11:01 AM Duration: 00:00:18 Operator: PromegaService	Software: Maxwell® CSC Software Version: 1.3 Instrument Name: PromegaInstrument Serial Number: beta unit 12 Firmware Version: 3.29.0 Maintenance Due: Yes
InitializationTest: Pass MotionTest: Pass HeaterTest: Pass	

Set Up a Run Quickly & Easily

Enhance Laboratory Workflow



Adapt to Your Workflow Needs

Maxwell® CSC provides higher throughput and automation that saves time and money, while the agile platform allows you to respond to changing laboratory demands on a daily or hourly basis.

Set up a Run Quickly

The intuitive touch-screen interface and preprogrammed protocols make it easy to set up a run, minimize variability and save time.

Reduce Processing Time

Extract nucleic acids from 1-16 samples in 38 minutes, with minimal pre-processing.

Maximize Bench Space

Easily fits into your laboratory, even if space is limiting. The compact footprint (330.2 × 457.2 × 330.2mm) allows laboratories to increase throughput and automate sample processing, while at the same time conserving critical laboratory bench space.

Increase Efficiency

Automating nucleic acid extraction allows higher throughput sample processing with minimal hands-on time, saving labor and decreasing costs.



Ensure Success in Downstream Assays

Quality Nucleic Acid Redefined



Increase Confidence in Your Results

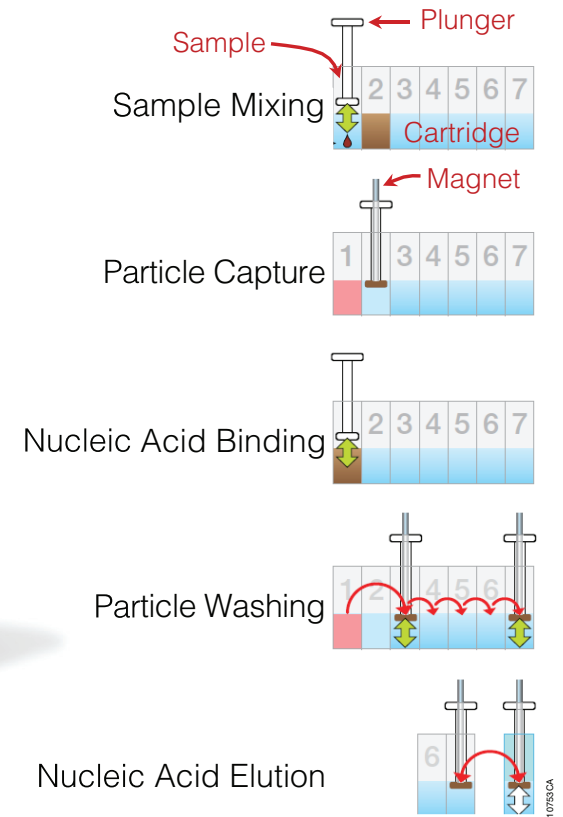
Because the Maxwell® CSC is manufactured to strict specifications, you can count on consistent, contamination-free sample preparation.

Magnetic Particle Mover, Not Liquid Handler

No liquid handling or splashing during processing means minimal risk of cross-contamination.

Performance-Tested Chemistries

Performance-tested chemistries means higher quality sample preparation and less variability between runs.



Magnetic particles are moved through the Maxwell cartridge wells.

Keep Your Lab Up & Running

Premier Service Package & Support

Plan with Assurance

The Maxwell® CSC Instrument includes the premier service package during the warranty period, routine preventive maintenance, field service and access to a loaner instrument, if required.

Improve Your Budget Planning

Service packages provide predictable support costs.

Enjoy Consistent and Reliable Performance

Preventive maintenance, replacement parts, and trained service personnel ensure your instrument is functioning properly.

Maintain Your Workflow

Field support and instrument loaner program ensure minimal instrument downtime for your laboratory.

All of this gives you confidence that your laboratory will experience minimal instrument downtime, allowing you to maintain throughput and productivity.

Available Service Options

SERVICE AGREEMENT BENEFIT	STANDARD	PREMIER
Period	Renewable annually	Renewable annually
Parts and Labor	Yes	Yes
Shipping	Yes	Yes
Loaner Instrument	Yes, 5-7 business days	Yes, 24 hours
Field Support	No	Yes, 48 hours
Annual Preventive Maintenance	No	Yes, 1 per 12 months

Our standard and premier service packages allow you to choose your level of support after the warranty period expires.

Ordering Information

Product	Size	Catalog Number
Maxwell® CSC Instrument	1 each	AS4000 ^{1,3}

Related Products	Size	Catalog Number
Maxwell® CSC Blood DNA Kit	48 preps	AS1321 ^{2,3}
Standard Service Agreement		SA1110 ⁴
Premier Service Agreement		SA1120 ⁴
Preventive Maintenance		SA1130 ⁴
Maxwell® CSC Installation Qualification		SA1140
Maxwell® CSC Operational Qualification		SA1150
Maxwell® CSC Installation and Operational Qualification		SA1160

¹ For In Vitro Diagnostic Use. Professional use only.

² Sufficient for 48 automated isolations from 300µl of whole blood samples. The Maxwell® CSC Cartridges are single use only.

³ Products are only available in the US.

⁴ Service offerings may differ by geography. Contact your local Promega representative for details on product availability.