

Maxwell[®] CSC Instruments

Automated Nucleic Acid Extraction for
Diagnostic and Clinical Research Applications





Maxwell[®] CSC Instruments

Maxwell[®] CSC Instruments easily integrate into any laboratory workflow. Their compact size minimizes the impact on precious laboratory space and maximizes sample capacity. Choose an instrument to suit your laboratory throughput needs. Automatically process 1–16 samples with the Maxwell[®] CSC System, or increase your capacity with the Maxwell[®] CSC 48 System to process 1–48 samples.

Reliable

The Maxwell[®] CSC Instruments use prefilled cartridges and preprogrammed methods to reduce the potential for user error. The Maxwell[®] CSC 48 includes an integrated vision system to verify correct deck tray setup prior to a run, further ensuring error-free operation. The purified nucleic acids are ready for use in diagnostic assays and are compatible with a wide range of downstream amplification-based applications.

Flexible

The Maxwell[®] CSC Instruments are a high-value laboratory addition, capable of both clinical research (RUO) and diagnostic (IVD) applications with dual-mode software. The instruments and chemistries work with a variety of sample types, such as human blood, plasma, swabs, tissues, and formalin-fixed, paraffin-embedded (FFPE) samples. No need to wait to process full runs. The Maxwell 1:1 Sample to Reagent Cartridge system allows you to process anywhere from 1 to 16 samples (Maxwell[®] CSC System) or 1 to 48 samples (Maxwell[®] CSC 48 System) simultaneously without wasting reagents.

Intuitive

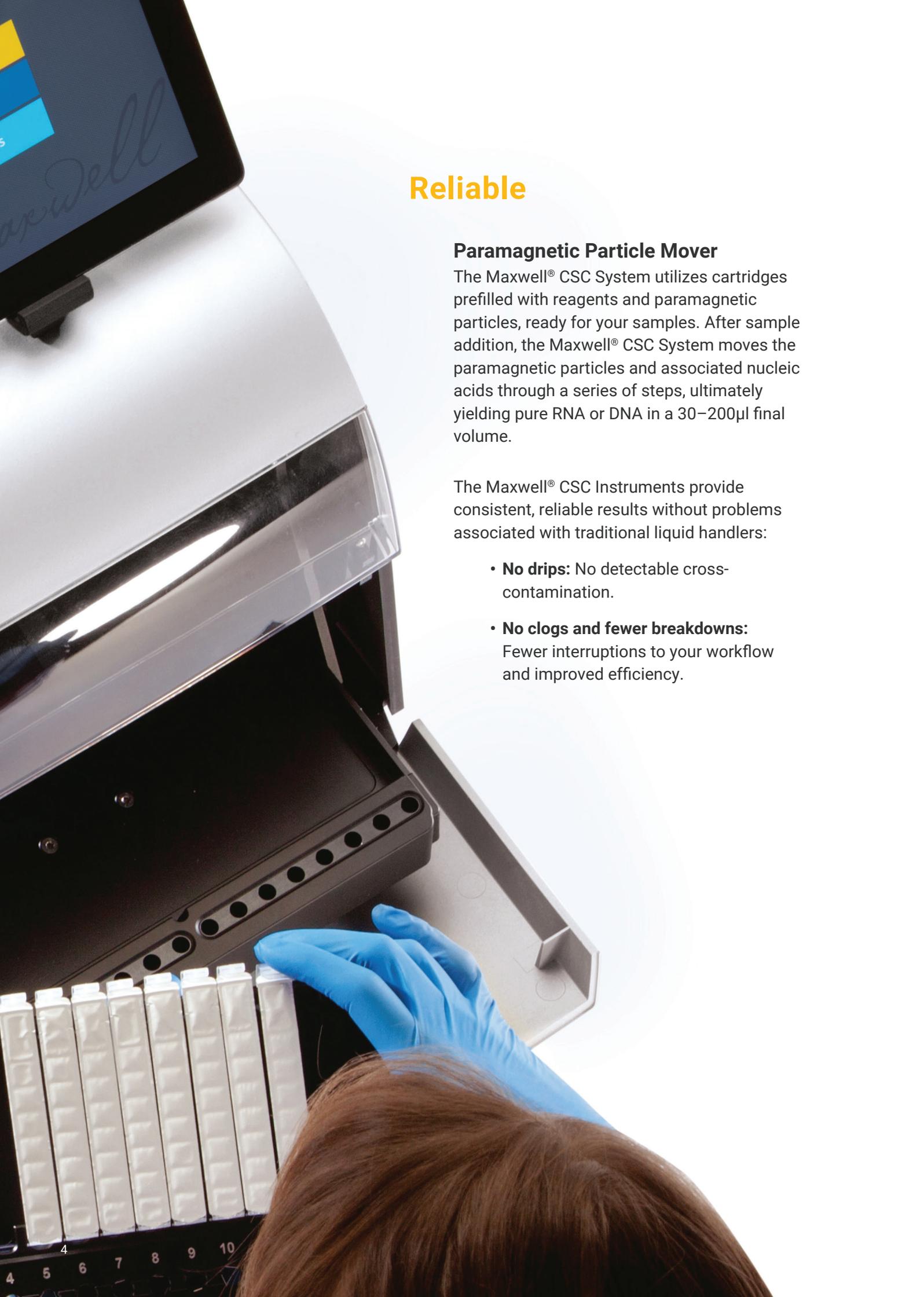
Maxwell[®] CSC software uses an easy to use and intuitive graphical user interface. You'll be able to set up and perform your first run in a matter of minutes after initial instrument setup.

Efficient

Automating nucleic acid extraction frees up your time. The integrated bar code reader eliminates the need to manually enter sample IDs and creates a streamlined, simple sample tracking report that easily exports to your laboratory notebook or is accessible to your laboratory informatics system (LIMS).

Supported

The Maxwell[®] CSC Instruments are backed by a comprehensive one-year warranty, our expert service team and a full line of additional service products (see page 11).



Reliable

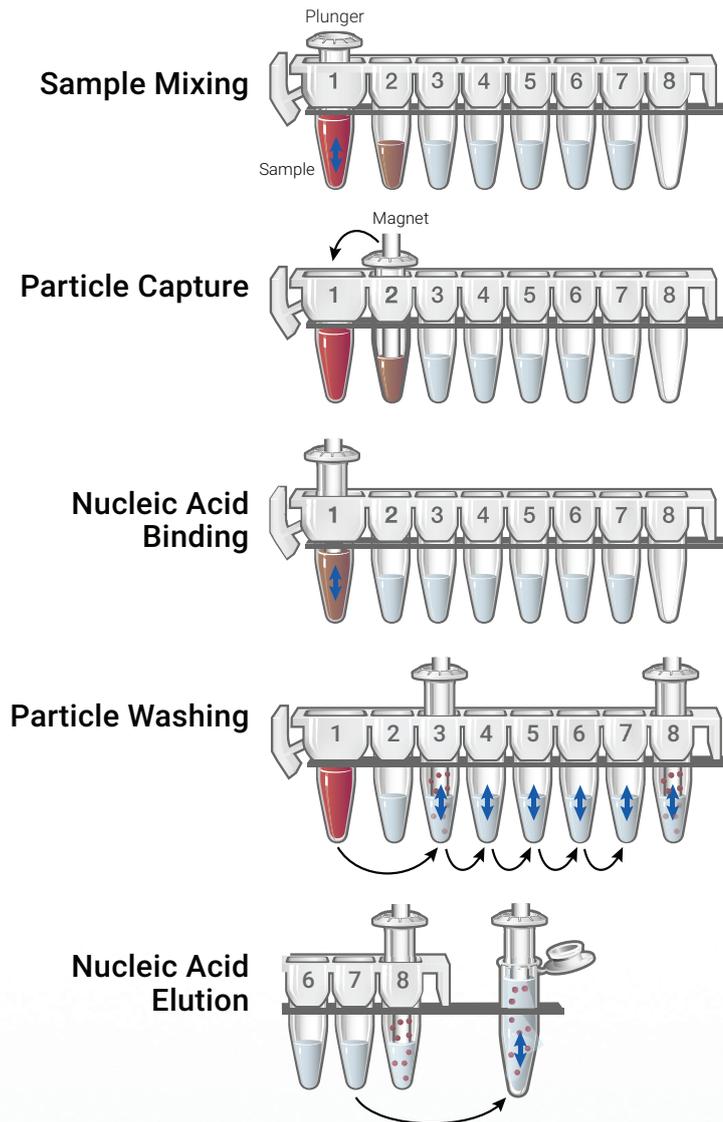
Paramagnetic Particle Mover

The Maxwell® CSC System utilizes cartridges prefilled with reagents and paramagnetic particles, ready for your samples. After sample addition, the Maxwell® CSC System moves the paramagnetic particles and associated nucleic acids through a series of steps, ultimately yielding pure RNA or DNA in a 30–200µl final volume.

The Maxwell® CSC Instruments provide consistent, reliable results without problems associated with traditional liquid handlers:

- **No drips:** No detectable cross-contamination.
- **No clogs and fewer breakdowns:** Fewer interruptions to your workflow and improved efficiency.

Paramagnetic Particle Workflow



A magnetic bar assembly, covered by a plastic sheath (plunger), moves the paramagnetic particles from well to well, eliminating clogs, breakdowns and cross-contamination.

Flexible

A single instrument platform capable of both clinical research (RUO) and diagnostic applications (IVD).

CE-MARKED, IVD-LABELED KITS			
AS1321	Maxwell® CSC Blood DNA Kit	AS1580	Maxwell® CSC Rapid ccfDNA Kit
AS1350	Maxwell® CSC DNA FFPE Kit	AS1780	Maxwell® CSC Viral Total Nucleic Acid Purification Kit
AS1360	Maxwell® CSC RNA FFPE Kit	AS1820	Maxwell® CSC Whole Blood DNA Kit
AS1410	Maxwell® CSC RNA Blood Kit	AS1850	Maxwell® CSC Genomic DNA Kit
AS1560	Maxwell® CSC XtractAll FFPE DNA/ RNA Kit	AS1860	Maxwell® CSC Pathogen Total Nucleic Acid Kit

RESEARCH USE ONLY KITS			
AS1330	Maxwell® RSC Viral Total Nucleic Acid Purification Kit	AS1590	Maxwell® RSC Rapid ccfDNA Kit
AS1340	Maxwell® RSC simplyRNA Tissue Kit	AS1600	Maxwell® RSC PureFood GMO and Authentication Kit
AS1370	Maxwell® RSC Cell DNA Purification Kit	AS1610	Maxwell® RSC Tissue DNA Kit
AS1380	Maxwell® RSC simplyRNA Blood Kit	AS1620	Maxwell® RSC Cultured Cells DNA Kit
AS1390	Maxwell® RSC simplyRNA Cells Kit		
AS1400	Maxwell® RSC Blood DNA Kit	AS1630	Maxwell® RSC Stabilized Saliva DNA Kit
AS1440	Maxwell® RSC RNA FFPE Kit	AS1640	Maxwell® RSC Buccal Swab DNA Kit
AS1450	Maxwell® RSC DNA FFPE Kit	AS1660	Maxwell® RSC PureFood Pathogen Kit
AS1460	Maxwell® RSC miRNA Tissue Kit	AS1680	Maxwell® RSC miRNA Plasma and Serum Kit
AS1480	Maxwell® RSC ccfDNA Plasma Kit	AS1700	Maxwell® RSC Fecal Microbiome DNA Kit
AS1490	Maxwell® RSC Plant DNA Kit	AS1720	Maxwell® RSC FFPE Plus DNA Kit
AS1500	Maxwell® RSC Plant RNA Kit	AS1831	Maxwell® RSC Enviro TNA Kit
AS1510	Maxwell® RSC Xcelerate DNA FFPE Kit	AS1840	Maxwell® RSC ccfDNA LV Plasma Kit
AS1520	Maxwell® RSC Whole Blood DNA Kit	AS1880	Maxwell® RSC Genomic DNA Kit
AS1540	Maxwell® RSC Buffy Coat DNA Kit	AS1890	Maxwell® RSC Pathogen Total Nucleic Acid Kit
AS1550	Maxwell® FSC DNA IQ™ Casework Kit		
AS1570	Maxwell® RSC XtractAll FFPE DNA/ RNA Kit	AS2110	Maxwell® RSC Purewater Kit

If you don't see the kit you want, we could be developing it right now. Contact your local Promega Representative for the most up-to-date information.

Intuitive

Intuitive Software Interface

Easy-to-use touch screen interface for walk-away, automated nucleic acid extraction at your fingertips.

Extraction in Only a Few Short Steps

1. Prepare samples and add to cartridge.
2. Scan kit to select appropriate protocol.
3. Select cartridge positions and assign sample IDs.
4. Start run.



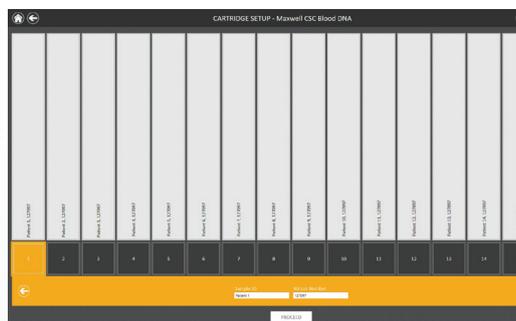
Prepare your sample, add it to a cartridge and load it into the Maxwell® CSC.



Select Start.



Scan your kit to select your method.



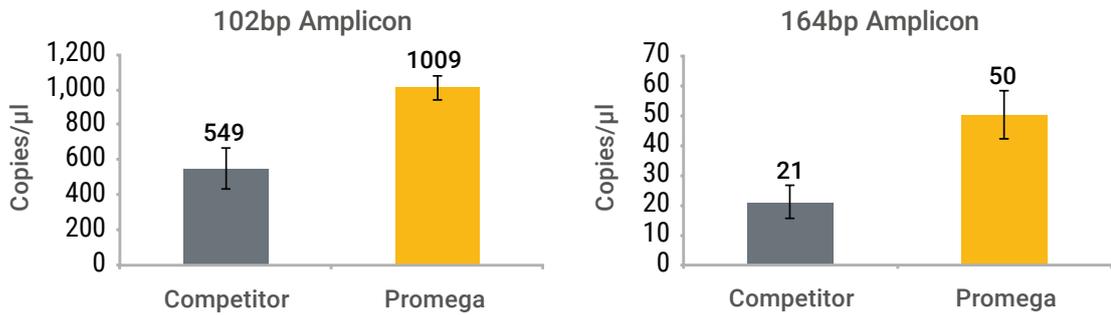
Assign sample IDs to their corresponding cartridge positions and run.

Versatile Instrument with an Easy-to-Use Interface

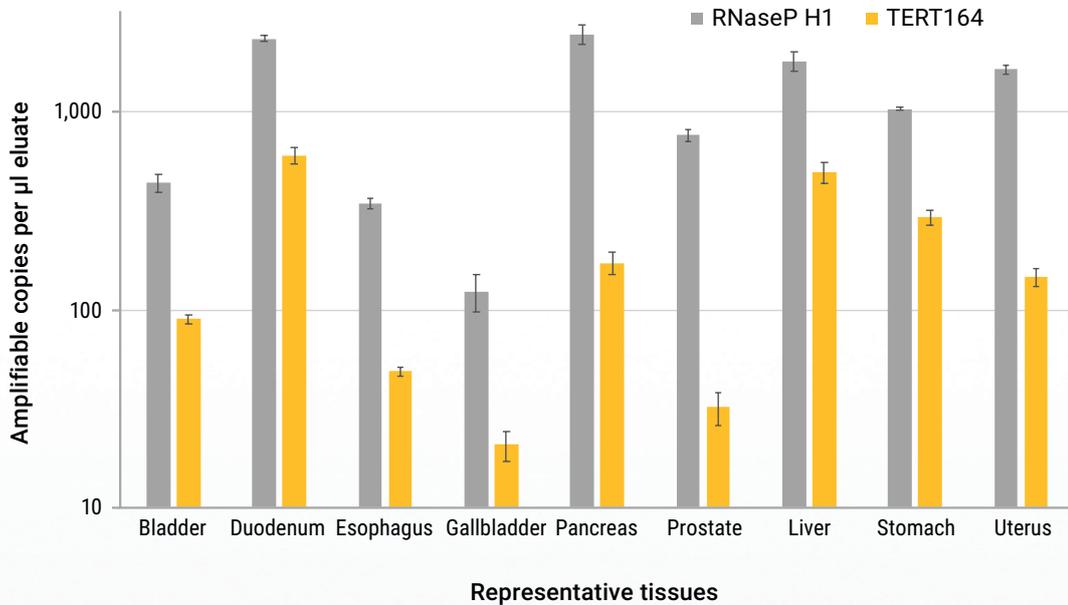
Quality and Performance

The Maxwell® CSC Instruments and chemistries provide high-quality nucleic acid for use in your downstream assays.

Obtain more Amplifiable DNA

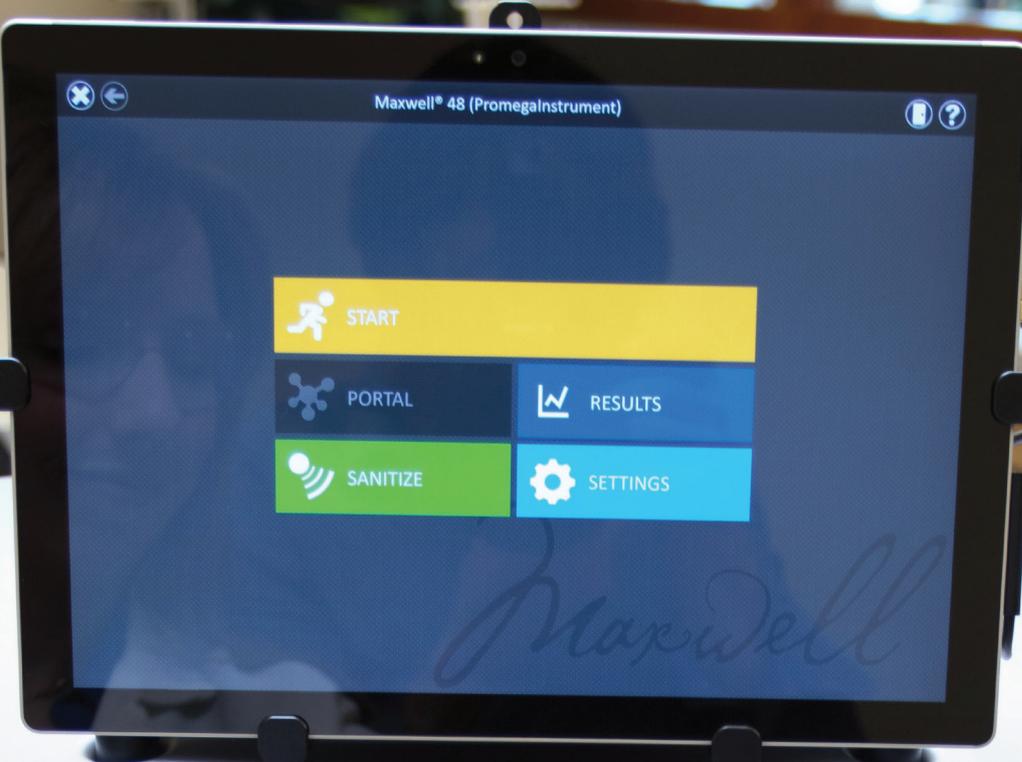


Broad Tissue Type Extraction of Amplifiable DNA from FFPE



Purification of amplifiable DNA from diverse tissue types using the Maxwell® CSC DNA FFPE Kit. Single FFPE sections were processed in triplicate using the Maxwell® CSC DNA FFPE Kit workflow on the Maxwell® CSC Instrument. Eluates were assessed by qPCR using a short amplicon (RNase P H1; 102bp) and with a longer amplicon (TERT; 164bp) to assess size distribution.





Supported

The Maxwell® CSC Instruments are backed by a comprehensive standard warranty, expert service and a full line of additional service products.

With a comprehensive menu of service and support options, you can be confident that your equipment will be working properly and ready when you need it.

Installation Qualification (IQ) and Operational Qualification (OQ) services ensure that your instrument is fully functional and provide documentation that the instrument is operating to defined specifications.

Available Service Options

SERVICE AGREEMENT	MAXWELL® CSC CAT.#	MAXWELL® CSC 48 CAT.#
Premier Warranty Upgrade	SA1111	SA1450
Standard Service Agreement	SA1110	SA1451
Premier Service Agreement	SA1120	SA1452
Preventive Maintenance	SA1130	SA1456
Installation Qualification	SA1140	SA1357
Operational Qualification	SA1150	SA1358
IQ/OQ Combination Package	SA1160	SA1359

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, 48 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.

*Where available

Ordering Information

Product	Cat.#
Elution Magnet, 16-Position	AS4017
Elution Magnet, 24-Position	AS4018
Maxwell® CSC Instrument	AS6000
Maxwell® CSC 48 Instrument	AS8000
Maxprep™ Liquid Handler*	AS9105/AS9205

Products may be covered by pending or issued patents or may have certain limitations.

Please visit our web site for more information.

Maxwell® and Maxprep™ are a trademark of Promega Corporation.

Note: The Maxwell® CSC products are not available in all regions.

*Maxwell® CSC instruments operated in RUO mode are compatible with the Maxprep™ Liquid Handler and Portal Software.

For more information, visit:

www.promega.com/maxwell-csc

www.promega.com/maxwell-csc-48



Promega GmbH
 Gutenbergring 10
 69190 Walldorf, Germany
www.promega.com

Germany/Austria
 Phone +49 6227 6906-290
de_techserv@promega.com

Poland
 Phone +48 22 531 06 68
pl_techserv@promega.com

Central/Eastern Europe
 Phone +49 6227 6906-0
de_ceesales@promega.com

