



Flexible NUCLEIC ACID *Solutions*

for your lab



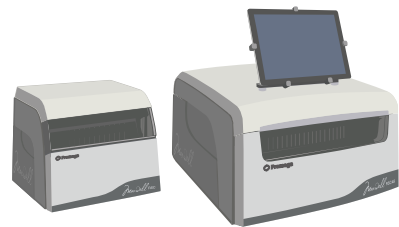
Nucleic Acid Preparation

made easy

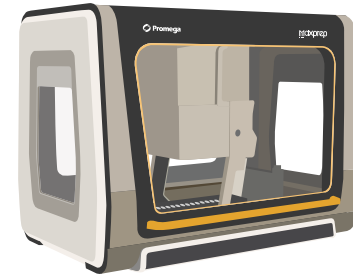
Our modular nucleic acid preparation solutions let you adapt your workflow as your laboratory needs change.

Sample Prep

- › Reagent Additions
- › Incubation
- › Elution Buffer Addition
- › Plunger Placement



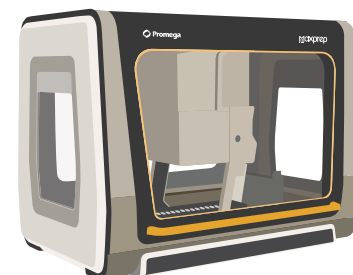
With prefilled cartridges and preprogrammed methods, the Maxwell® RSC Instruments offer reliable, consistent nucleic acid purification with the flexibility to process only the sample number you need.



Preprogrammed to work with the Maxwell® RSC Nucleic Acid Purification Kits, the Maxprep™ Liquid Handler simplifies and automates sample and Maxwell® cartridge preparation.

Extraction

Maxwell® RSC (1–16 Samples)
Maxwell® RSC 48 (1–48 samples)



With post-extraction methods for nucleic acid quantitation, normalization, eluate transfers and amplification setup, the Maxprep™ Liquid Handler simplifies your downstream assay preparation.

Post-Extraction

- › Quantitation Dye Addition
- › Sample Normalization, eluate transfers
- › PCR, qPCR, RT-qPCR Prep



Streamline Your Sample Handling *with one instrument*

Increased Consistency

Whether you are processing samples for nucleic acid purification with your Maxwell® RSC or RSC 48 Instrument or preparing purified samples for downstream amplification reactions, the Maxprep™ Liquid Handler ensures that every sample is handled the same way every time regardless of the operator. The easy-to-use software interface means anyone in your lab will be able to quickly run samples using the preprogrammed protocols. With no methods to create or protocols to write, the Maxprep™ Liquid Handler is ready to do the tedious work for you.

With preloaded methods for sample preprocessing and post-purification applications, the Maxprep™ Liquid Handler easily integrates wherever you need it in your laboratory workflow.

Preloaded Post-Purification Protocols

ELUATE TRANSFER

Transfer eluted nucleic acid from Maxwell® elution tubes to more permanent storage containers. Reformat and combine multiple Maxwell® runs from elution tubes to microplate formats.

NORMALIZATION

Import quantitation data into the Maxprep™ software and normalize your samples:

- › Normalize on mass or concentration
- › Normalize entire eluate volumes
- › Normalize small aliquots of eluates, while preserving the rest of the sample for later use

QUANTITATION

Add quantitation dyes such as Promega Quantifluor® Dyes

PCR SET UP

- › qPCR
- › RT-qPCR
- › PCR
- › RT-PCR



Flexible, Modular *nucleic acid purification*



Increased Flexibility

As your personal purification assistant, the Maxwell® RSC or RSC 48 Instrument easily integrates into your laboratory workflow, while their compact size minimizes the impact on precious laboratory space. Use the Maxwell® RSC for smaller sample batches up to 16 samples, or use the Maxwell® RSC 48 Instrument for sample batches up to 48 samples.

New for Maxwell® RSC 48! A large LED Indicator and an integrated vision system for detecting proper cartridge setup further reduces the potential for user error.

FLEXIBLE

The Maxwell® RSC Instruments work with a variety of sample types, including human blood, cells and tissues. The protocols are fast and easy. You can process anywhere from 1 to 16 samples simultaneously with the Maxwell® RSC and 1 to 48 samples with the Maxwell® RSC 48.

RELIABLE

Using prefilled cartridges and preprogrammed methods reduces the potential for user error. Purified nucleic acids are compatible with a wide range of downstream applications such as PCR, qPCR, Sanger and next generation sequencing (NGS), micro arrays and digital PCR.

EFFICIENT

Automating nucleic acid purification frees up your time. The integrated quantitation eliminates the need to re-enter sample IDs and creates a streamlined, simple sample tracking report that easily exports to your laboratory notebook.

INTUITIVE

Maxwell® RSC software is so easy to use and intuitive that you'll be able to set up and perform your first run in a matter of minutes after initial instrument setup.



Your Complete *nucleic acid purification system*

Continuous Purification

Create a virtual purification assembly line, preparing Maxwell® cartridges on the Maxprep™ Liquid Handler while your Maxwell® RSC Instrument extracts cartridges that have already been prepared. Multiple modules can be placed anywhere in the laboratory where space is available and communicate sample tracking information from one device to the next. Simply transfer the Maxwell RSC deck tray from one module onto the next and all the sample information will be there ready for you to process.



Preprocess samples
with the Maxprep™
Liquid Handler

Purify nucleic acids
with the Maxwell®
RSC or RSC 48

Keep working.
Preprocessing and
purification can happen
simultaneously.

Streamline your post-
purification process
(normalize, quantitate,
PCR, RT-qPCR, qPCR)
with Maxprep™

Increased Traceability

The Promega Portal software tracks samples as they move from one module to the next. The Portal software automatically shares sample IDs and recommends protocols for next steps as you move from one module to the next.

Sample Tracking

The Maxprep™ Liquid Handler automatically scans the sample bar codes as they are placed into the instrument. Once you scan the desired Maxwell® RSC Kit and the bar code on the Maxwell® Deck Trays, the Promega Portal software does the rest. Portal monitors where each sample is placed in the Deck Tray, what Maxwell® kit was used, and passes this information directly to the Maxwell® RSC or RSC 48 Instrument for walk-up extraction.

Adaptable

The Maxprep™ Liquid Handler can accommodate up to two Maxwell® RSC Deck Trays for sample and cartridge preparation.



World Class *service & support*

Our 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), Operational Qualification (OQ) and Preventive Maintenance (PM) services ensure that your instrument is fully functional and provide documentation that the instrument is operating to defined specifications.

Maxprep™ Liquid Handler

The Maxprep™ Liquid Handler is installed by qualified Promega personnel who will test the instrument, document installation actions and train lab personnel to successfully operate the instrument.

Ordering *information*

Expand and adapt your laboratory automation based on your throughput needs. Our modular workflow allows you to incorporate existing Maxwell® RSC Instruments into the Maxprep™ Liquid Handler workflow and add additional units as your throughput needs grow.

Maxprep™ Configurations	Cat.#
Maxprep™ Liquid Handler, RSC Carriers	AS9100
Maxprep™ Liquid Handler, RSC 48 Carriers	AS9200
Maxprep™ Liquid Handler, RSC Carriers w/ UV Light	AS9101
Maxprep™ Liquid Handler, RSC 48 Carriers w/ UV Light	AS9201

Maxprep™ Accessories and Consumables	Cat.#
Maxprep™ Carrier, Maxwell® RSC	AS9402
Maxprep™ Carrier, Maxwell® RSC 48 Front	AS9403
Maxprep™ Carrier, Maxwell® RSC 48 Back	AS9404
Maxprep™ 50µl Conductive Filter Tips (60 Tip Boxes)	AS9301
Maxprep™ 300µl Conductive Filter Tips (60 Tip Boxes)	AS9302
Maxprep™ 1000µl Conductive Filter Tips (40 Tip Boxes)	AS9303

Maxwell® Instruments	Cat.#
Maxwell® RSC 48 Instrument	AS8500
Maxwell® RSC Instrument	AS4500

Maxwell® Accessories	Cat.#
Maxwell® RSC Bar Code Reader	AS3200
Maxwell® RSC Deck Tray	SP6019
Maxwell® RSC 48 Front Deck Tray	AS8401
Maxwell® RSC 48 Back Deck Tray	AS8402



Instrument *specifications*

Maxwell® RSC

Sample Capacity: 1–16 Samples

Intended Use:
Research Use Only

Dimensions (W × D × H):
13 × 13.6 × 11.8 inches
(330.2 × 345.2 × 299.7 mm)

Weight: 24.2 lbs (11 kg)

Main Power Voltage:
100–240V AC, 50/60Hz;
autosensing; 2.5 Amps

Operating Environment:
15–25°C

Service and Warranty: 1-year
limited warranty on parts and labor.
Extended service contracts

UV-Bulb: Average lifetime
approximately 6,000 hours

Maxwell® RSC 48

Sample Capacity: 1–48 Samples; (2)
24 position, removable trays

Intended Use:
Research Use Only

Dimensions (W × D × H):
21 × 20 × 12.5 inches
(533 × 508 × 317.5 mm)

Weight: 70 lbs (31.75 kg) est.

Main Power Voltage:
100–240V AC, 50/60Hz;
autosensing; 4 Amps

Operating Environment:
15–25°C

Service and Warranty: 1-year
limited warranty on parts and labor.
Extended service contracts

UV-Bulb: Average lifetime
approximately 6,000 hours

MaxPrep™ Liquid Handler

Sample Capacity: 1–48 Samples; (2)
24 position Maxwell RSC 48 or (2) 16
position Maxwell RSC removable trays

Intended Use:
Research Use Only

Dimensions (W × D × H):
42.1 × 27.8 × 33 inches
(106.9 × 70.6 × 83.3 cm)

Note: the total instrument height with the
door open is 48.4inches (123cm)

Weight: 220 lbs (98.6 kg) est.

Main Power Voltage:
Input: 100–240VAC, 50/60Hz, 5A

Operating Environment:
15–25°C

Service and Warranty: 1-year
limited warranty on parts and labor.
Extended service contracts

UV-Bulb: Average lifetime
approximately 10,000 hours