

TECHNICAL MANUAL

Maxwell[®] 16 LEV Blood DNA Kit and Maxwell[®] 16 Buccal Swab LEV DNA Purification Kit

Instructions for Use of Products
AS1290 and AS1295

Caution: Handle cartridges with care; seal edges may be sharp.



Maxwell[®] 16 LEV Blood DNA Kit and Maxwell[®] 16 Buccal Swab LEV DNA Purification Kit

All technical literature is available at: www.promega.com/protocols/
Visit the web site to verify that you are using the most current version of this Technical Manual.
E-mail Promega Technical Services if you have questions on use of this system: techserv@promega.com

1.	Description.....	1
2.	Product Components and Storage Conditions	2
3.	Before You Begin.....	3
3.A.	Preparation of Whole Blood Samples (Cat.# AS1290)	3
3.B.	Preparation of Buccal Swab Samples (Cat.# AS1295)	4
3.C.	Maxwell [®] 16 Cartridge Preparation	4
4.	Instrument Run: AS2000 and AS3000 Instruments	6
5.	Reference	8
6.	Troubleshooting.....	9
7.	Related Products.....	9
8.	Summary of Changes	10

1. Description

The Maxwell[®] 16 LEV Blood DNA Kit and Maxwell[®] 16 Buccal Swab LEV DNA Purification Kit are used with the Maxwell[®] 16 Instrument to provide an easy method for efficient, automated purification that concentrates genomic DNA (gDNA) from whole blood and buccal swab samples, respectively. The Maxwell[®] 16 Instrument is supplied with preprogrammed purification procedures and is designed for use with the predispensed reagent cartridges, maximizing simplicity and convenience. The instrument can process up to 16 samples in 40 minutes. The purified DNA can be used directly in a variety of downstream applications, including PCR and agarose gel electrophoresis.

The Maxwell[®] 16 Instrument purifies samples using a novel paramagnetic particle, called the MagnaCel[™] particle, which provides a mobile solid phase that optimizes sample capture, washing and purification of gDNA. This particle utilizes cellulose-based binding of nucleic acids and provides a higher bind capacity and cleaner eluate than traditional silica-based DNA purification. The Maxwell[®] 16 Instrument is a magnetic particle-handling instrument that efficiently binds gDNA to the paramagnetic particle in the first well of a prefilled cartridge and mixes during processing. The approach to magnetic capture avoids common problems such as clogged tips or partial reagent transfers that result in suboptimal purification processing by other commonly used automated systems.



2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
Maxwell® 16 LEV Blood DNA Kit	48 preps	AS1290

For Laboratory Use. Sufficient for 48 automated isolations from 300µl of whole blood samples. Includes:

- 2 × 1ml Proteinase K (PK) Solution
- 20ml Lysis Buffer
- 48 Maxwell® 16 LEV Cartridges (MCD)
- 50 LEV Plungers
- 50 Elution Tubes (0.5ml)
- 20ml Elution Buffer

Storage Conditions: Store the Maxwell® 16 LEV Blood DNA Kit at 15–30°C.

PRODUCT	SIZE	CAT.#
Maxwell® 16 Buccal Swab LEV DNA Purification Kit	48 preps	AS1295

For Laboratory Use. Sufficient for 48 automated isolations from buccal swab samples. Includes:

- 2 × 1ml Proteinase K (PK) Solution
- 20ml Lysis Buffer
- 48 Maxwell® 16 LEV Cartridges (MCD)
- 50 LEV Plungers
- 50 Elution Tubes (0.5ml)
- 20ml Elution Buffer
- 50 Clearing Columns

Storage Conditions: Store the Maxwell® 16 Buccal Swab LEV DNA Purification Kit at 15–30°C.

Available Separately (recommended for sample extraction)

PRODUCT	SIZE	CAT.#
ClickFit Microtube, 1.5ml	1,000/pack	V4741

Safety Information: The reagent cartridges contain ethanol and isopropanol. These substances should be considered flammable, harmful and irritants.



The Maxwell® 16 reagent cartridges are designed to be used with potentially infectious substances. Users should wear appropriate protection (e.g., gloves and goggles) when handling infectious substances. Users should adhere to their institutional guidelines for the handling and disposal of all infectious substances when used with this system.

3. Before You Begin

Maxwell® 16 Instrument Hardware and Firmware Setup

To use the Maxwell® 16 LEV Blood DNA Kit or Buccal Swab LEV DNA Purification Kit, the Maxwell® 16 Instrument must be configured with LEV hardware. If your Maxwell® 16 Instrument contains standard elution volume (SEV) hardware, it will need to be reconfigured using the Maxwell® 16 LEV Hardware Kit (Cat.# AS1250). Reconfiguring the instrument is simple and easy. Refer to the *Maxwell® 16 Instrument Technical Manual* specific for your instrument for directions.

Materials to Be Supplied by the User

- optional, rotating tube mixer for liquid blood samples
- benchtop vortex mixer
- pipettors and pipette tips for sample transfer into pre-filled reagent cartridges
- 1.5–2.0ml tubes for incubation of samples (e.g., ClickFit Microtube, 1.5ml [Cat.# V4741]; recommended to prevent the cap from opening during heating)
- heating block set at 56°C
- microcentrifuge (buccal swab protocol)
- buccal swabs (e.g., Puritan Medical Products Cat.# 25-806 1PD or 25-806 1PC)

3.A. Preparation of Whole Blood Samples (Cat.# AS1290)

Whole Blood Sample Processing Capacity

The total yield of genomic DNA from whole blood samples depends on the sample volume and number of white blood cells/ml. Each cartridge supplied in the Maxwell® 16 LEV Blood DNA Purification Kit is designed to purify genomic DNA from up to 300µl of whole blood, assuming an average number of white blood cells in the range of 4×10^6 to 1.1×10^7 /ml whole blood (values for a normal healthy adult; 1).

Note: Whole blood samples collected in EDTA, ACD or heparin tubes can be used. These samples may be either fresh or frozen. Frozen samples should be thawed before processing. We recommend mixing all blood samples before use. EDTA blood collection tubes are preferred if the purified DNA will be used in downstream amplification assays.

1. Mix all blood samples for at least 5 minutes at room temperature.
2. Prepare and label incubation tubes compatible with heating block.
3. Add 30µl of Proteinase K (PK) Solution to each incubation tube.
4. Add liquid blood (up to 300µl) to each incubation tube.
5. Add 300µl of Lysis Buffer to each incubation tube.
6. Vortex each tube for 10 seconds.
7. Incubate each tube in the heating block (set to 56°C) for 20 minutes. During this incubation, prepare cartridges as described in Section 3.C.
8. Transfer each blood lysate sample from the incubation tube to well #1 of each cartridge. (Well #1 is the well closest to the cartridge label and furthest from the user.)

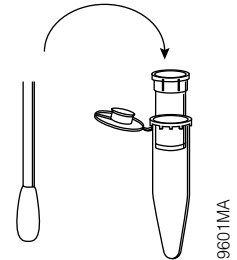
3.B. Preparation of Buccal Swab Samples (Cat.# AS1295)

Buccal Swab Sample Processing Capacity

The total yield of genomic DNA from buccal swab samples depends on the cellular material contained on the swabs. Each cartridge supplied in the Maxwell® 16 Buccal Swab LEV DNA Purification Kit is designed to purify genomic DNA from 1 or 2 buccal swabs.

Note: Customers have performed this protocol successfully with samples collected using Pur-Wraps Sterile Polyester Tipped Applicators, Puritan Medical Products Cat.# 25-806 1PD or Pur-Wraps Sterile Cotton Tipped Applicators, Puritan Medical Products Cat.# 25-806 1PC. We also recommend using ClickFit Microtube, 1.5ml (Cat.# V4741) for sample preparation.

1. Collect samples with a standard buccal swab collection procedure.
2. Assemble a Clearing Column/microtube for each sample.
3. Cut the head off the applicator stick. Add dried swab head to the Clearing Column/microtube assembly.
4. In a separate tube, mix 300µl Lysis Buffer + 30µl Proteinase K for each sample.
5. Add 330µl of Lysis Buffer/Proteinase K to swab head in the Clearing Column/microtube assembly.



6. Close tube over the Clearing Column and vortex for 10 seconds.

Note: If using tubes other than the recommended Cat.# V4741, the tube may not close.

7. Incubate for 20 minutes at 56°C.

Note: Some flowthrough from the column may be observed in the microtube after incubation. This is normal.

8. Centrifuge the Clearing Column/microtube assembly with swab for 2 minutes at maximum speed.
9. Remove the Clearing Column with swab head and discard.
10. Add flowthrough to well #1 of the Maxwell® 16 LEV cartridge (see Section 3.C for cartridge preparation).
11. Process using Maxwell® 16 in LEV Research Mode using the LEV Blood method (see Section 4).
12. Once the extraction is complete, remove and cap each elution tube. Store appropriately until use.

3.C. Maxwell® 16 Cartridge Preparation

1. Change gloves before handling cartridges, LEV Plungers and Elution Tubes. Place the cartridges to be used in the Maxwell® 16 LEV Cartridge Rack (Cat.# AS1251). Place each cartridge in the rack with the label side facing away from the Elution Tubes. Press down on the cartridge to snap it into position. Carefully peel back the seal so that all plastic comes off the top of the cartridge. Ensure that all sealing tape and any residual adhesive are removed before placing cartridges in the instrument.
2. Place one plunger into well #8 of each cartridge.

- Place an empty Elution Tube into the Elution Tube position for each cartridge in the Maxwell® 16 LEV Cartridge Rack. Add 50µl of Elution Buffer to the bottom of each Elution Tube.

Notes:

- If you are processing fewer than 16 samples, center the cartridges on the platform.
- Specimen or reagent spills on any part of the Maxwell® 16 LEV Cartridge Rack should be cleaned with a detergent-water solution, followed by a bacteriocidal spray or wipe, then water. Do not use bleach on any instrument parts.



Figure 1. Maxwell® 16 LEV DNA Purification Cartridge. This figure shows the contents of a cartridge. In all cases, lysate sample is added to well #1.



Figure 2. Setup and configuration in the Maxwell® 16 LEV Cartridge Rack. Elution Buffer is added to the Elution Tubes as indicated.



4. Instrument Run: AS2000 and AS3000 Instruments

Setup for AS2000 Maxwell® 16 Instruments

Refer to the *Maxwell® 16 Instrument Operating Manual #TM295* for more detailed information.

To run the “Blood” protocol, you must have Maxwell® 16 firmware version 4.71 or higher installed on your instrument.

1. Turn on the Maxwell® 16 Instrument. The instrument will power up, display the firmware version number, proceed through a self-check and home all moving parts.
2. Verify that the instrument settings indicate an “LEV” hardware configuration and “Rsch” operational mode setting.
3. Select “Run” on the Menu screen, and press the Run/Stop button to start the method.
4. Select “DNA” on the menu screen, then select “OK” at the Verification screen.
5. Select “Blood” on the Menu screen, then select “OK” at the Verification screen. “Blood” Method is used for both kits (Cat.# AS1290 and AS1295).
6. Open the door when prompted to do so on the screen. Press the Run/Stop button to extend the platform.



Warning: Pinch point hazard.

7. Transfer the Maxwell® 16 LEV Cartridge Rack containing the prepared cartridges on the Maxwell® 16 Instrument platform. Ensure that the rack is placed in the Maxwell® 16 Instrument with the Elution Tubes closest to the door. The rack will only fit in the instrument in this orientation. If you have difficulty fitting the rack on the platform, check that the rack is in the correct orientation. Ensure that the cartridge rack is level on the instrument platform.

Note: Hold the Maxwell® 16 LEV Cartridge Rack by the sides to avoid dislodging cartridges from the rack.

8. Verify that samples were added to well #1 of the cartridges, cartridges are loaded on the instrument, Elution Tubes are present with 50µl of Elution Buffer and LEV Plungers are in well #8.
9. Press the Run/Stop button. The platform will retract. Close the door.



Warning: Pinch point hazard.

10. The Maxwell® 16 Instrument will immediately begin the purification run. The screen will display the steps performed and the approximate time remaining in the run.

Notes:

1. Pressing the Run/Stop button or opening the door will pause the run.
 2. If the run is abandoned before completion, the instrument will wash the particles off the plungers and eject the plungers into well #8 of the cartridge. The sample will be lost.
11. When the automated purification run is complete, the LCD screen will display a message that the method has ended.

End of Run

12. Follow on-screen instructions at the end of the method to open door. Verify that plungers are located in well #8 of the cartridge at the end of the run. If plungers are not removed from the magnetic plunger bar, push them down gently by hand to remove them.
13. Press the Run/Stop button to extend the platform out of the instrument.
14. Remove the Maxwell® 16 LEV Cartridge Rack from the instrument. Remove Elution Tubes containing DNA, and close the tubes.



Note: Following the automated purification procedure, the LEV Cartridge Rack will be warm. It will not be too hot to touch. To remove the rack from the instrument platform, hold onto the sides of the rack.

15. Remove the cartridges and plungers from the Maxwell® 16 LEV Cartridge Rack, and discard as hazardous waste.



Do not reuse reagent cartridges, LEV Plungers or Elution Tubes.

Setup for AS3000 Maxwell® 16 MDx Instruments

Refer to the *Maxwell® 16 MDx Instrument Technical Manual #TM320* for detailed information. To run the “Blood” protocol, you must have the Maxwell® 16 Firmware version 1.10 or higher installed on your instrument

1. Turn on the Maxwell® 16 MDx Instrument. The instrument will power up, display the firmware version number, proceed through a self-check and home all moving parts.
2. Verify that the Home screen indicates “LEV” and the LEV hardware is present. Press “Run” to continue.
3. Enter user and PIN, if this option is enabled.
4. At the Protocols screen, select “DNA”.
5. At the Method screen, select “Blood”.
6. On the next screen, verify that the correct user was chosen. The protocol should read “DNA”. Select “Run/Stop” to continue.
7. Open the door when prompted on the screen, then select “Run/Stop”.



Warning: Pinch point hazard.

8. Follow on-screen instructions for bar code reader input if this option is enabled.
9. Transfer the Maxwell® 16 LEV Cartridge Rack containing the prepared cartridges on the Maxwell® 16 Instrument platform. Ensure that the rack is placed in the Maxwell® 16 Instrument with the Elution Tubes closest to the door. The rack will only fit in the instrument in this orientation. If you have difficulty fitting the rack on the platform, check that the rack is in the correct orientation. Ensure the rack is level on the instrument platform.

Note: Hold the Maxwell® 16 LEV Cartridge Rack by the sides to avoid dislodging cartridges from the rack.

Setup for AS3000 Maxwell® 16 MDx Instruments (continued)

10. Verify that samples were added to well #1 of the cartridges, cartridges are loaded on the instrument, Elution Tubes are present with 50µl of Elution Buffer and LEV Plungers are in well #8.
11. Press the Run/Stop button. The platform will retract. Close the door.



Warning: Pinch point hazard.

The Maxwell® 16 Instrument will immediately begin the purification run. The screen will display the approximate time remaining in the run.

Notes:

1. Pressing the Run/Stop button or opening the door will pause the run.
 2. If the run is abandoned before completion, the instrument will wash the particles off the plungers and eject the plungers into well #8 of the cartridge. The samples will be lost.
12. When the automated purification run is complete, follow instructions on the screen for data transfer. For detailed instructions, refer to the *Maxwell® 16 MDx Instrument Technical Manual #TM320* and *Maxwell® Sample Track Software Technical Manual #TM314*.

End of Run

13. Follow on-screen instructions at the end of the method to open door. Verify that plungers are located in well #8 of the cartridge at the end of the run. If plungers are not removed from the magnetic plunger bar, push them down gently by hand to remove them.
14. Press the Run/Stop button to extend the platform out of the instrument.
15. Remove the Maxwell® 16 LEV Cartridge Rack from the instrument. Remove Elution Tubes containing DNA, and cap the tubes.



Note: Following the automated purification procedure, the LEV Cartridge Rack will be warm. It will not be too hot to touch. To remove the rack from the instrument platform, hold onto the sides of the rack.

16. Remove the cartridges and plungers from the Maxwell® 16 LEV Cartridge Rack, and discard as hazardous waste.



Do not reuse reagent cartridges, LEV Plungers or Elution Tubes.

For the Maxwell® 16 MDx Instrument, ensure samples are removed before the UV light treatment to avoid damage to the nucleic acid.

5. Reference

1. Henry, J.B. (2001) *Clinical Diagnosis and Management by Laboratory Methods*, 20th ed., W.B. Saunders Company, 509.

6. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com. E-mail: techserv@promega.com

Symptoms

Lower than expected A_{260}
(lower than expected yield)

Causes and Comments

Blood that has undergone multiple freeze-thaw cycles may have degraded DNA. Use fresh samples whenever possible, or avoid multiple freeze-thaw cycles.

Proteinase K Solution was not added. The lysis and yield are dependent upon complete extraction with proteinase K. If proteinase K was not added in Section 3.A, Step 3, the resulting blood sample will be red. Proteinase K-treated samples turn greenish brown. This can be used as a quick diagnostic method of determining whether or not the proteinase K was added.

Whole blood sample contained low white blood cell count. The yield of genomic DNA from blood samples depends on the number of white blood cells present in the sample

Whole blood sample was not mixed before processing. Be sure to mix whole blood samples before processing to ensure that the white blood cells are in suspension.

In some cases, total RNA can be copurified with the genomic DNA. To remove copurified RNA, an RNase treatment can be performed. Add 5 μ l of RNase A (Cat.# A7973) per milliliter of Elution Buffer.

If yield is low for the buccal swab protocol, your swab may not contain enough cellular material. Examine the buccal swab collection method.

7. Related Products

Product	Size	Cat.#
Maxwell® 16 LEV Hardware Kit	1 each	AS1250
RNase A Solution, 4mg/ml	1ml	A7973
ClickFit Microtube, 1.5ml	1,000/pack	V4741



8. Summary of Changes

The following change was made to the 8/17 revision of this document:

1. Added “whole blood and buccal swab“ to the text near the end of the first sentence of Section 1.

© 2010–2017 Promega Corporation. All Rights Reserved.

Maxwell is a registered trademark of Promega Corporation. MagnaCel is a trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.