Maxwell® RSC Stabilized Saliva DNA Kit

Instructions for Use of Product AS1630.



Preparing Stabilized Saliva Samples for DNA Purification

This Quick Protocol provides instructions for use of the Maxwell® RSC Stabilized Saliva DNA Kit to purify DNA from stabilized saliva samples. Purification has been tested with the DNA Genotek Oragene® DISCOVER DNA Collection Kit, the IsoHelix GeneFiXTM Saliva DNA Collection Device and the Biomatrica DNAgard® Saliva HT Collection Device. For detailed instructions, including information on instrument setup and troubleshooting, please refer to the *Maxwell® RSC Stabilized Saliva DNA Kit Technical Manual* #TM480 available at: www.promega.com/protocols/

Materials to be Supplied by the User

- benchtop vortex mixer
- pipettors and pipette tips for sample transfer
- Stabilized Saliva Collection Device

Collecting Saliva Samples

1. Collect samples with a standard saliva collection procedure.

Note: Storing the collectors for more than two days after collection may improve extraction performance.

- 2. Vortex the stabilized saliva until the sample is homogenous (>10 seconds).
- 3. Add up to 1ml of stabilized saliva to well #1 (the largest well) of the Maxwell® RSC cartridge.

Maxwell® Automated DNA Purification

Cartridge Preparation

- 1. Place the cartridges to be used in the deck tray with well #1 (the largest well) facing away from the elution tube.
- 2. 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 the cartridge in the instrument.
- 3. Place one plunger in well #8 of each cartridge. Well #8 is the well closest to the elution tube.
- 4. Place an empty elution tube into the elution tube position for each cartridge. Add 150µl of Elution Buffer to the bottom of each elution tube.

Note: Use only the Elution Tubes (0.5ml) provided with the kit; other tubes may be incompatible with supported Maxwell[®] Instruments.

Instrument Run on the Maxwell® Instruments

Follow the instrument run instructions in the Maxwell® RSC Stabilized Saliva DNA Kit Technical Manual #TM480.

Maxwell is a registered trademark of Promega Corporation.

DNAgard is a registered trademark of Biomatrica, Inc. GeneFix is a trademark of Cell Projects Ltd. Oragene is a registered trademark of DNA Genotek, a subsidiary of OraSure Technologies, Inc.

Additional protocol information is in Technical Manual #TM480, available online at: www.promega.com

