

# Maxwell® RSC Whole Blood DNA Kit

Instructions for Use of Product AS1520.



Quick Protocol

## Preparing Whole Blood Samples for DNA Purification

### Materials to Be Supplied by the User

- **optional:** rotating tube mixer for liquid blood samples
- pipettors and pipette tips for sample transfer into prefilled Maxwell® reagent cartridges

The total yield of genomic DNA from whole blood samples depends on the sample volume and number of white blood cells/ml.

**Note:** Fresh or frozen whole blood samples collected in EDTA, ACD or heparin tubes can be used.

1. Mix all blood samples for at least 5 minutes at room temperature.
2. Transfer 50–500µl of each blood sample from the starting blood tube to well #1 (the largest well) of each cartridge.

## Maxwell® Automated Purification

### Cartridge Preparation

1. Place the cartridge to be used in the Deck Tray(s) 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 cartridges in the instrument.
3. Place a plunger in well #8 of each cartridge. Well #8 is the well closest to the elution tube.
4. Place an empty elution tube in the elution tube position for each cartridge. Add at least 60µl of Elution Buffer to the bottom of each elution tube.

**Notes:** Use only the Elution Tubes (0.5ml) provided in 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 Whole Blood DNA Kit Technical Manual* #TM455.

Maxwell is a registered trademark of Promega Corporation.

Additional protocol information is in Technical Manual #TM455, available online at: [www.promega.com](http://www.promega.com)

