

Maxwell® RSC Cell DNA Purification Kit

INSTRUCTIONS FOR USE OF PRODUCT AS1370.

Preparation of Cells for DNA Purification

Materials to Be Supplied by the User

- Microtubes, 1.5ml (Cat.# V1231) or larger
- PBS (10.1mM Na₂HPO₄, 1.8mM KH₂PO₄, 140mM NaCl, 2.7mM KCl [pH 7.3])

Preprocessing Samples

1. Pellet cells from the sample by centrifugation.
2. Remove the supernatant, being careful not to disturb the cell pellet.
3. Resuspend the pellet in up to 100µl of PBS.

Maxwell® RSC Automated DNA Purification

Cartridge Preparation

1. Place the cartridge to be used in the Deck Tray with the printed side 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.

Note: If you are processing fewer than 16 samples, center the cartridges on the Deck Tray.

3. Place a plunger in well #8 of each cartridge. Well #8 is the well closest to the elution tube.
4. Place elution tubes in the front of the Deck Tray. Add 50µl of Elution Buffer to the bottom of each elution tube.

Notes:

1. If Elution Buffer is on the side of the tube, the elution may be suboptimal.
2. Use only the Elution Tubes (0.5ml) provided with the kit; other tubes may be incompatible with the Maxwell® RSC Instrument.
5. Transfer the cell suspension to well #1 of the cartridge. Well #1 is the well closest to the printed side and furthest from the elution tube.

Note: The total volume of the sample should not exceed 100µl.

Instrument Run on the Maxwell® RSC Instrument

1. Refer to the *Maxwell® RSC Instrument Operating Manual* #TM411 for detailed information.
2. Follow the instrument run instructions in the *Maxwell® RSC Cell DNA Purification Kit Technical Manual* #TM418.

Ordering and Technical Information

www.promega.com • Phone 608-274-4330

Toll free in USA 800-356-9526 • Fax 608-277-2601