

DNA from Tissue using Maxwell® RSC Blood DNA Kit

Purify high quality DNA from Tissue using the Maxwell® RSC Blood DNA Kit on the Maxwell® RSC Instrument.

Kit: Maxwell® RSC Blood DNA Kit (Cat.# AS1400)

Analyses: TapeStation

Sample Type(s): Tissue

Input: Up to 200mg

Materials Required:

- Maxwell® RSC Blood DNA Kit (Cat.# AS1400)
- Maxwell® RSC Instrument (Cat.# AS4500)
- Tissue Lysis Buffer (Cat.# A5091)
- Heat block

This protocol was developed by Promega Applications Scientists and is intended for research use only.

Users are responsible for determining suitability of the protocol for their application.

For further information, see Technical Manual TM419, available at:

www.promega.com/protocols

or contact Technical Services at: techserv@promega.com

Protocol:

1. Add 200–1000µl of Tissue Lysis Buffer to the tissue sample. Add sufficient buffer to submerge tissue. Higher amounts of tissue will require more Tissue Lysis Buffer for complete homogenization.
2. Add 20µl Proteinase K (PK) Solution.
3. Incubate at 56°C for 1 hour. Vortex 2-3 times during the incubation. This should result in complete homogenization of the tissue.
4. *Optional* if debris remains, centrifuge homogenate to pellet any remaining solids.
5. Transfer 200µl to well #1 of a Maxwell® RSC Blood DNA Cartridge.
6. Place a plunger in well #8 of the cartridge.
7. Add 100µl of Elution Buffer to Elution Tubes.
8. Select the Maxwell® RSC Blood DNA method and run.

Results:

High-quality DNA was purified from 200mg human breast tissue using 1ml of Tissue Lysis Buffer, without the optional centrifugation step, using the Maxwell® RSC Blood DNA Kit.

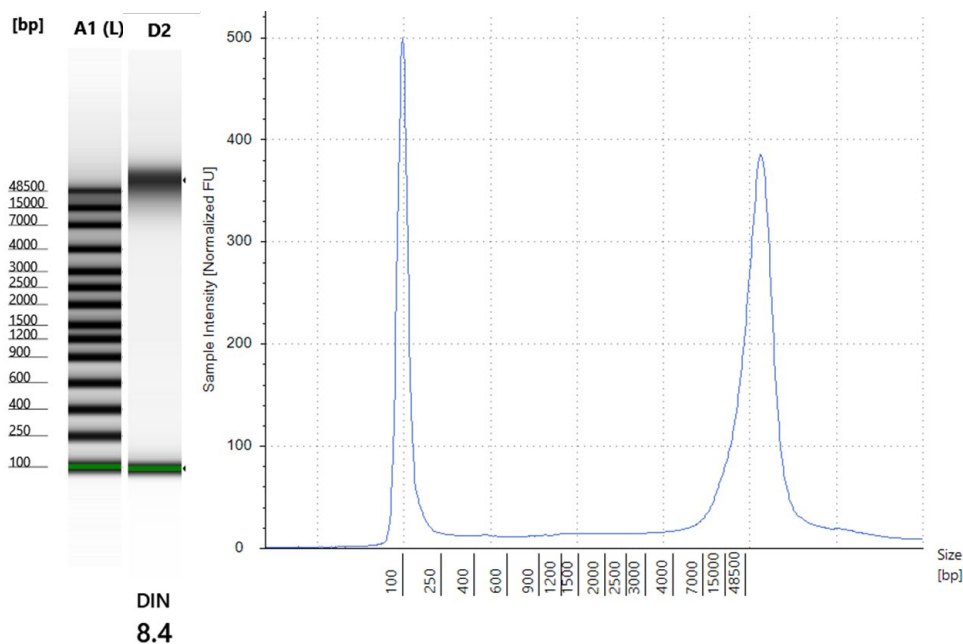


Figure 1. Genomic DNA TapeStation electrophoresis of DNA purified from breast tissue. DNA was purified from 200mg of human breast tissue using the Maxwell® RSC Blood DNA Kit (Cat.# AS1400) and Maxwell® RSC Instrument (Cat.# AS4500). DNA was then analyzed by electrophoresis on an Agilent TapeStation 4200 using the Genomic DNA ScreenTape and Reagents.