

DNA Purification from Oil Palm Leaf using the Maxwell® RSC System

Isolate high-quality, amplifiable DNA from oil palm leaf tissue using the Maxwell® RSC System.

Kit: Maxwell® RSC Plant DNA Kit (Cat.# AS1490)

Analyses: NanoDrop, QuantiFluor® quantitation, agarose gel electrophoresis, qPCR

Sample Type(s): Oil Palm, *Elaeis guineensis*

Input: 1–5 punches (~1.7mg/punch)

Materials Required:

- Maxwell® RSC Instrument (Cat.# AS4500)
- Maxwell® RSC Plant DNA Kit (Cat.# AS1490)
- 2.0ml screw-top tubes
- homogenization steel bead
- bead-beating device (MP Biomedicals Fast-Prep®-24 Instrument)
- microcentrifuge

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 TM458, available at:

www.promega.com/protocols

or contact Technical Services at:
techserv@promega.com

Protocol:

1. Using a 4mm punch, place up to 5 punches into a 2ml screw-top tube.
2. Add 300µl of Tissue Lysis Buffer and 10µl of RNase A solution to each sample tube.
3. Using the bead-beating device, homogenize samples for desired time (e.g., FastPrep®-24 Instrument at 4M/S, 20 seconds × 4, with 20-second delay between each time).
4. Centrifuge the samples in a microcentrifuge at maximum speed for 2 minutes.
5. Add 300µl of Nuclease-Free Water to Well #1 of the Maxwell® cartridge.
6. Transfer the entire volume of supernatant to Well #1 of the Maxwell® cartridge.
7. Place one of the supplied elution tubes into the sample rack, and add 50µl of the supplied Elution Buffer for each sample.
8. Place the plunger in Well #8 of the cartridge.
9. On the Maxwell® RSC Instrument, select the Plant DNA Kit protocol. Start Run.

Results:

DNA was purified from oil palm leaves with the Maxwell® RSC Plant DNA Kit, using 1–5 punches of input (approximately 0.0017–0.0085g).

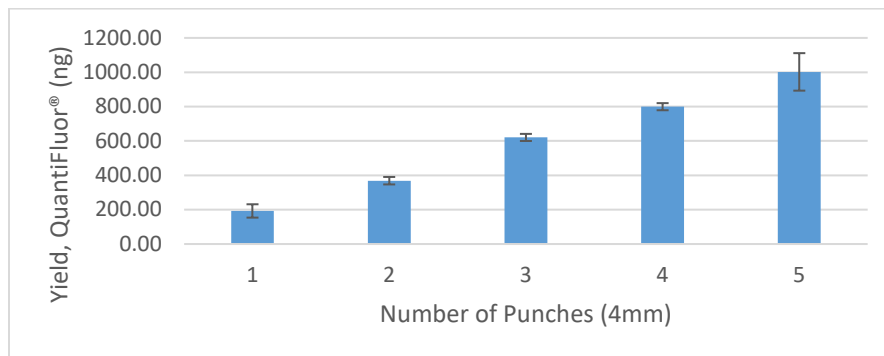


Figure 1. DNA yield purified from oil palm leaf and measured by QuantiFluor® ONE dsDNA System. Example of the linearity of DNA yield from from 1–5 punches of oil palm leaf input.

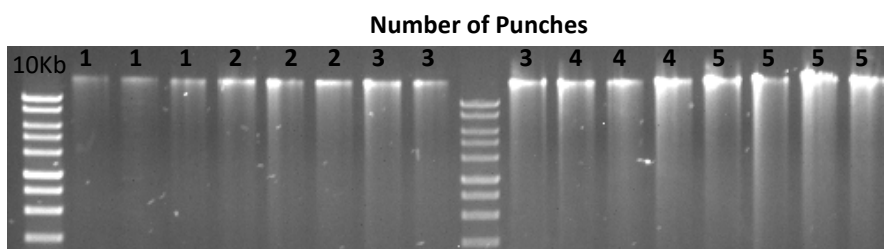


Figure 2. DNA integrity for purified oil palm DNA. Examples of high molecular weight DNA purified from indicated number of punches and electrophoresed on a 1% agarose gel.

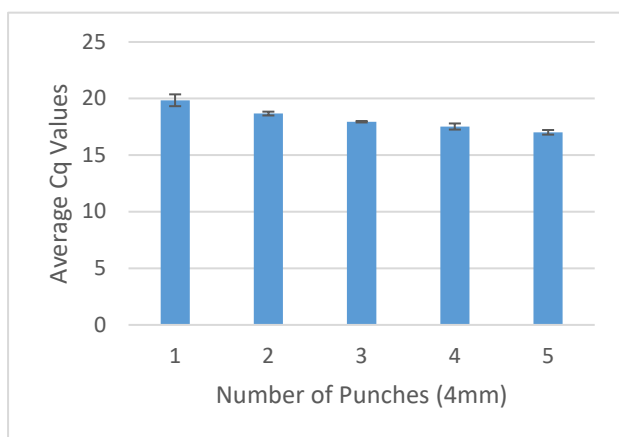


Figure 3. qPCR amplification of purified oil palm DNA. GoTaq® qPCR amplification with Universal Plant Primers (ATP-1) using 2µl of undiluted DNA eluate per reaction.

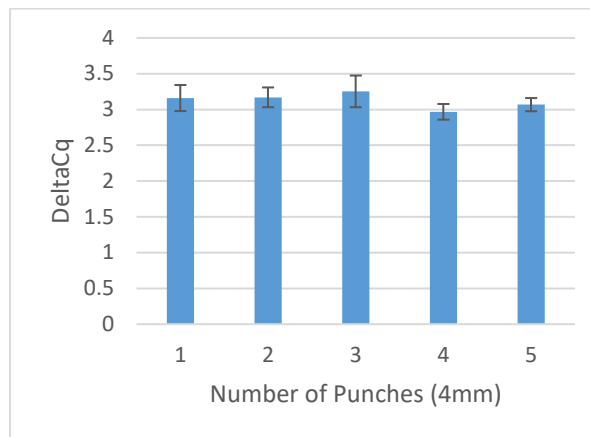


Figure 4. qPCR amplification inhibition measured by ΔCq values for DNA purified from oil palm. Samples were amplified undiluted, and diluted 1/10. A ΔCq of ~3.3 indicates no qPCR inhibition.