

DNA Isolation from Cotton Leaf Tissue using the ReliaPrep™ gDNA Tissue Miniprep System

Isolate high-quality, amplifiable DNA from cotton leaf tissue using the ReliaPrep™ gDNA Tissue Miniprep System.

Kit: ReliaPrep™ gDNA Tissue Miniprep System (Cat.# A2051)

Sample Type: Cotton leaf tissue

Input: Up to 25mg

Materials Required:

- ReliaPrep™ gDNA Tissue Miniprep System (Cat.# A2051)
- 2.0ml screw-top tubes
- homogenization steel bead
- bead-beating device (e.g., MP Biomedicals FastPrep®-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 TM345, available at:

www.promega.com/protocols

or contact Technical Services at:
techserv@promega.com

Protocol:

1. Using a 5mm punch, place desired number of punches (up to 25mg) into a 2ml screw-top tube.
2. To each sample add:
 - 100µl of Tail Lysis Buffer (TLA)
 - 300µl of Cell Lysis Buffer (CLD)
 - 20µl of RNase A Solution
 - 20µl of Proteinase K
3. Using the bead-beating device, homogenize samples for desired time (e.g., FastPrep®-24 Instrument at 6m/s, 40 seconds).
4. Centrifuge samples in a microcentrifuge at max speed for 1 minute.
5. Incubate at room temperature for 10 minutes.
6. Centrifuge samples at max speed for 1 minute to reduce foaming.
7. Add 250µl of Binding Buffer (BBA) to each sample and vortex for 10 seconds.
8. Centrifuge samples at max speed for 2 minutes, and transfer liquid supernatant to a ReliaPrep™ Binding Column inside a collection tube.
9. Centrifuge samples at max speed for 1 minute. Transfer column to a new collection tube; discard the flowthrough and used collection tube.
10. Add 500µl of Column Wash Solution (CWD) to the sample and centrifuge at max speed for 2 minutes. Repeat this wash step for a total of 3 times, discarding liquid and collection tubes after every wash.
11. Place the column in a clean 1.5ml microcentrifuge tube.
12. Add 50µl Nuclease-Free Water to the column. Centrifuge for 1 minute at max speed. Eluates are ready for use in downstream applications.

Results:

Table 1. Cotton leaf DNA concentration, yield and purity based on quantitation using the QuantiFluor® ONE dsDNA System (Cat.# E4871) and the NanoDrop®-1000. DNA of high purity was recovered with purity ratios for samples >1.7. N=3.

Sample	NanoDrop			QuantiFluor® ONE	
	ng/μl	A ₂₆₀ /A ₂₈₀	A ₂₆₀ /A ₂₃₀	ng/μl	Yield (μg)
Cotton Leaf	81.62	2.02	1.75	33.67	1.52

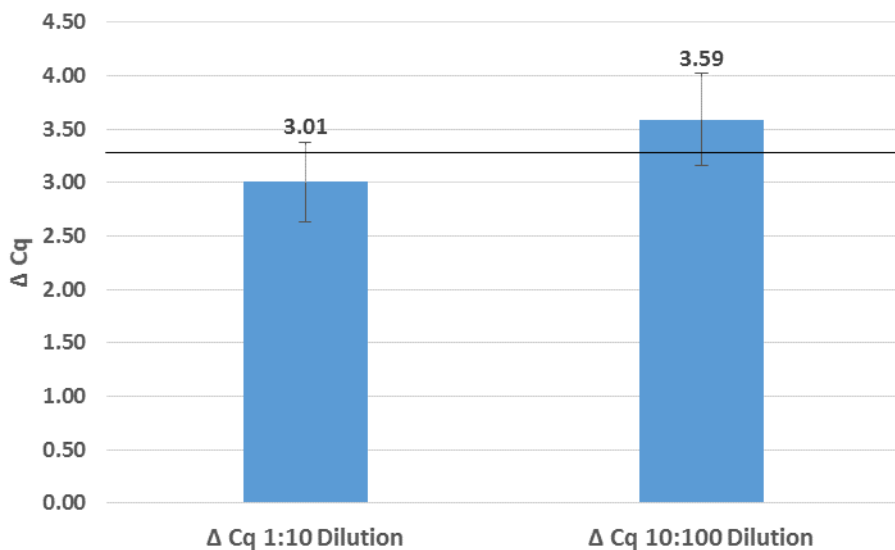


Figure 1. Inhibition analysis of purified cotton leaf DNA. DNA samples were serially diluted 1:10 and 10:100. For a sample diluted tenfold, ΔCq values are expected to be 3.3. ΔCq values significantly less than 3.3 may indicate the presence of inhibitors. ΔCq values of DNA from cotton leaf samples indicate little to no inhibition of the serially diluted eluates. N=3.