Instructions for Use of Products A2351 and A2352.



DNA Isolation with Deparaffinization Using Mineral Oil

Materials to Be Supplied By the User

- 95-100% ethanol
- 80°C heat block
- 56°C heat block
- 1.5 or 2ml microcentrifuge tubes
- equivalent of ≤100µm tissue sections (see Technical Manual #TM352)

Notes:

- Add Blue Dye to the Lysis Buffer before starting the procedure (see Technical Manual #TM352).
- Add 95–100% ethanol to the 1X Wash Solution before starting the procedure.
- Perform all centrifugations at room temperature.

Deparaffinization Using Mineral Oil

- 1. Add mineral oil to the sample:
 - For sections \leq 50 microns, add 300µl of mineral oil.
 - For sections > 50 microns, add 500µl of mineral oil.
- 2. Incubate at 80°C for 1 minute.
- 3. Vortex to mix.

Sample Lysis

- 1. Add 200ul of Lysis Buffer (with Blue Dye added) to the sample.
- 2. Centrifuge at $10,000 \times g$ for 15 seconds. Two phases will be formed, a lower blue (aqueous) phase and an upper (oil) phase.
- 3. Add 20µl of Proteinase K directly to the lower phase and mix by pipetting.
- 4. Incubate at 56°C for 1 hour.
- 5. Incubate at 80°C for 4 hours.

Note: Optimal recovery of amplifiable DNA is obtained with a 4-hour incubation at 80°C. Incubation may be reduced to 1 hour, but it will result in lower DNA yields.

6. Allow the sample to cool to room temperature. Centrifuge briefly to collect any condensation.

Optional storage: After incubating at 80°C, samples may be stored overnight at 2–10°C. If samples are stored this way, allow them to warm to room temperature prior to adding RNase and proceeding with the protocol.

RNase Treatment

- 1. Add 10µl of RNase A directly to the lysed sample in the lower phase. Mix the lower phase by pipetting.
- 2. Incubate at room temperature (20–25°C) for 5 minutes.



DNA Isolation with Deparaffinization Using Mineral Oil (continued)

Nucleic Acid Binding

- 1. Add 220µl of BL Buffer to the lysed sample.
- 2. Add 240 μ l of ethanol (95–100%). Vortex briefly to mix.
- 3. Centrifuge at $10,000 \times g$ for 15 seconds. Two phases will be formed, a lower blue (aqueous) phase and an upper (oil) phase.
- 4. For each sample to be processed, place a Binding Column into one of the Collection Tubes provided. **Note:** Wear gloves when handling the columns and tubes.

 Transfer the entire lower blue (aqueous) phase of the sample, including any precipitate that may have formed, to the Binding Column/Collection Tube assembly, and cap the column. Discard the remaining mineral oil.

Note: The mineral oil is inert and will not interfere with the extraction procedure if some of the oil phase is carried over to the Binding Column.

- 6. Centrifuge the assembly at $10,000 \times g$ for 30 seconds.
- 7. Discard the flowthrough, and reinsert the Binding Column into the Collection Tube.
- 8. Proceed immediately to column Washing and Elution.

Column Washing and Elution

- 1. Add 500µl of 1X Wash Solution (with ethanol added) to the Binding Column. Cap the column.
- 2. Centrifuge at $10,000 \times g$ for 30 seconds.
- 3. Discard the flowthrough, and reinsert the Binding Column into the same Collection Tube.
- 4. Add 500µl of 1X Wash Solution to the Binding Column. Cap the column.
- 5. Centrifuge at 10,000 \times *g* for 30 seconds.
- 6. Discard the flowthrough, and reinsert the Binding Column into the same Collection Tube.
- 7. Open the cap on the Binding Column, and centrifuge the Binding Column/Collection Tube assembly at $16,000 \times g$ for 3 minutes to dry the column.

Note: Centrifuging with the cap open ensures thorough drying of the column. It is important to dry the column to prevent carryover of ethanol to the eluate.

- 8. Transfer the Binding Column to a clean 1.5ml microcentrifuge tube (not provided), and discard the Collection Tube.
- 9. Add 30–50µl of Elution Buffer to the column, and cap the column.
- 10. Centrifuge at 16,000 \times *g* for 1 minute. Remove and discard the Binding Column.
- 11. Cap the microcentrifuge tube, and store the eluted DNA at -30 to -10° C.

Additional protocol information in Technical Manual #TM352, available online at: www.promega.com