

ReliaPrep™ FFPE Miniprep Systems

The ReliaPrep™ FFPE Miniprep Systems provide complete, streamlined methods for manual purification of genomic DNA (gDNA) and total RNA from formalin-fixed paraffin-embedded (FFPE) tissue.

FFPE samples are precious and the fixation process can be damaging to DNA and RNA embedded in these samples. The optimized lysis, binding, and elution reagents in the ReliaPrep™ FFPE Miniprep Systems result in increased functional yield from FFPE samples without using harsh organic solvents or overnight digestions.

Isolate Quality, Intact Nucleic Acids: Optimized lysis and binding conditions reverse modifications introduced by the fixation process, resulting in intact, amplifiable nucleic acids.

Safely Deparaffinize Your Sample: Deparaffinization step occurs without harsh organic solvents.

Save Time: Purify gDNA or total RNA from FFPE tissue in less than two and one-half hours with minimal hands-on time. No overnight digestion required.

Easy to Use: Minimal preparation time; simply add ethanol and go!

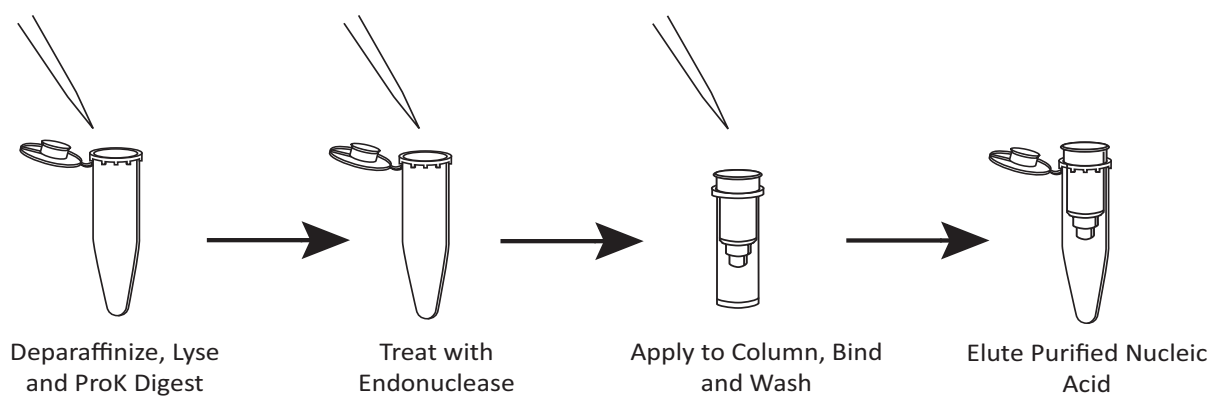


Figure 1. The streamlined workflow allows for rapid and effective deparaffinization, endonuclease treatment (in lysis buffer), binding and elution. Total time—including reagent preparation and deparaffinization is approximately 2½ hours for gDNA and 2 hours for total RNA.

Assessing Fragmentation of Purified Nucleic Acids

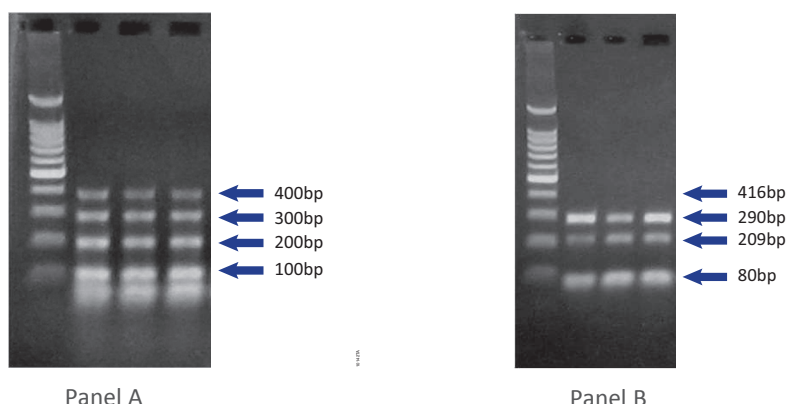


Figure 2. Multiplex PCR of variable length, end-point PCR fragments of GAPDH gene from nucleic acid purified from 10µm slices of FFPE human liver. Panel A. 10µl purified gDNA used as template in multiplex PCR (GoTaq® Hot Start). Fragments ranging from 100-400bp detected. Panel B. 5µl purified RNA used as template in GoScript™ Reverse Transcription System. 5µl of cDNA product used as template for the multiplex PCR (GoTaq® Hot Start). Fragments from 80bp-290bp detected.

All reactions performed in triplicate from sequential slices. 10µl of PCR product were run on a 2% agarose gel.

Downstream Applications Data

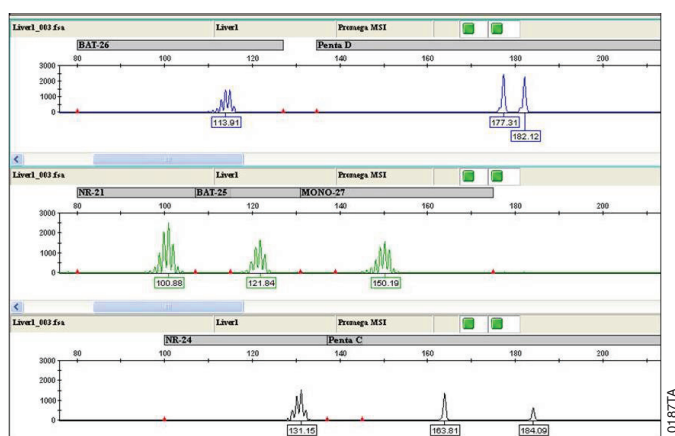


Figure 3. Genomic DNA was purified from 10µm human liver FFPE sections using the ReliaPrep™ FFPE gDNA Miniprep System. STR analysis was performed using Promega's Microsatellite Instability (MSI) Analysis System v. 1.2 (Cat.# MD1641). This system is compatible with STR amplification and MSI detection.

Ordering Information

PRODUCT	SIZE	CAT.#
ReliaPrep™ FFPE gDNA Miniprep System	10 reactions	A2351
	100 reactions	A2352
ReliaPrep™ FFPE Total RNA Miniprep System	10 reactions	Z1001
	100 reactions	Z1002