

FEATURES
AND BENEFITS

NUCLEIC
ACID INTEGRITY

FUNCTIONAL YIELD

WORKFLOW

RELATED
PRODUCTS

ORDERING
INFORMATION

RESOURCES



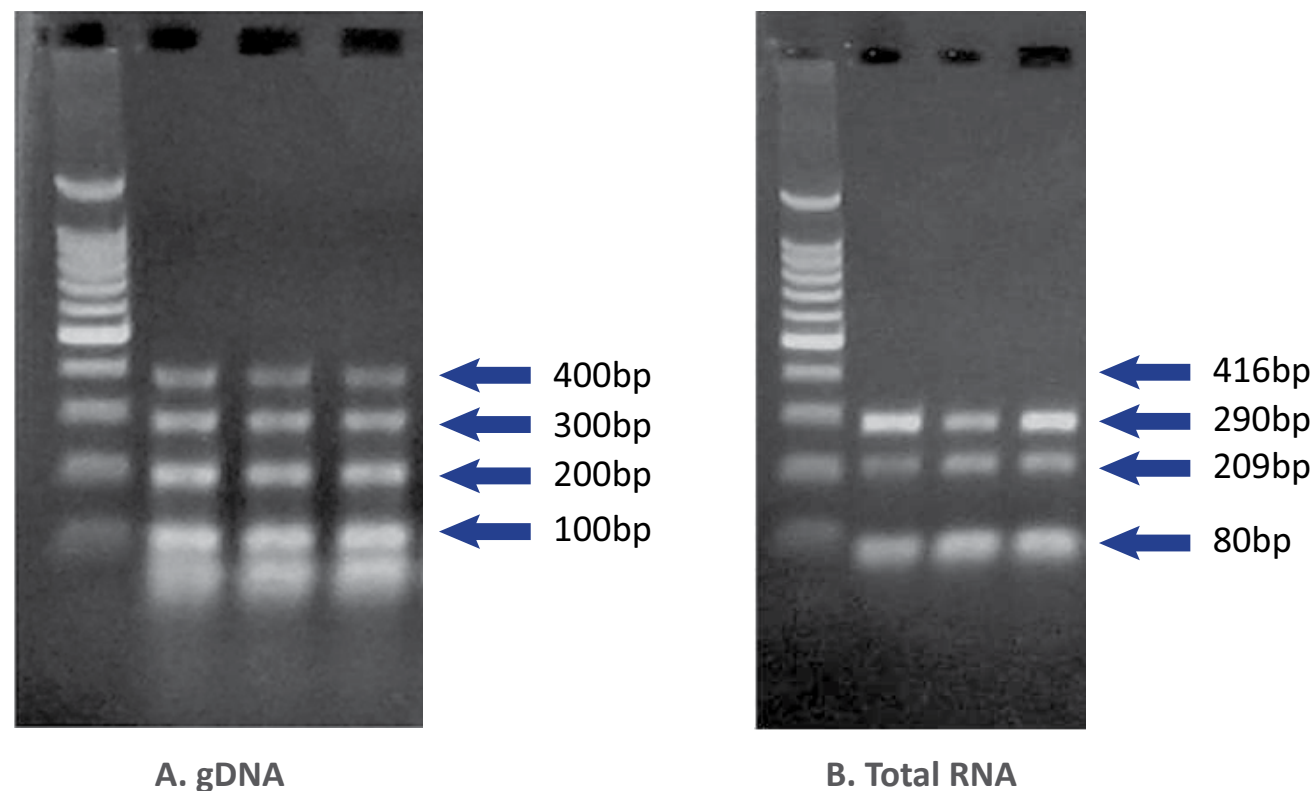
ReliaPrepTM FFPE Miniprep Systems

*Purify High-Quality, Intact Nucleic Acids from FFPE
Tissue Without Harsh 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 and total RNA from FFPE tissue in less than two and half hours with minimal hands on time. No overnight digestion required.
- ✓ **Easy to use:** Minimal preparation time; simply add ethanol and go!

Nucleic Acid Integrity

Nucleic acids from FFPE tissue are characteristically fragmented. ReliaPrep™ FFPE Miniprep Systems help preserve the integrity of your nucleic acids.



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 PCR product run on a 2% agarose gel.

Multiplex PCR of variable length, end-point PCR fragments of GAPDH gene from nucleic acid purified from 10µm slices of FFPE human liver.

FEATURES
AND BENEFITS

NUCLEIC
ACID INTEGRITY

FUNCTIONAL YIELD

WORKFLOW

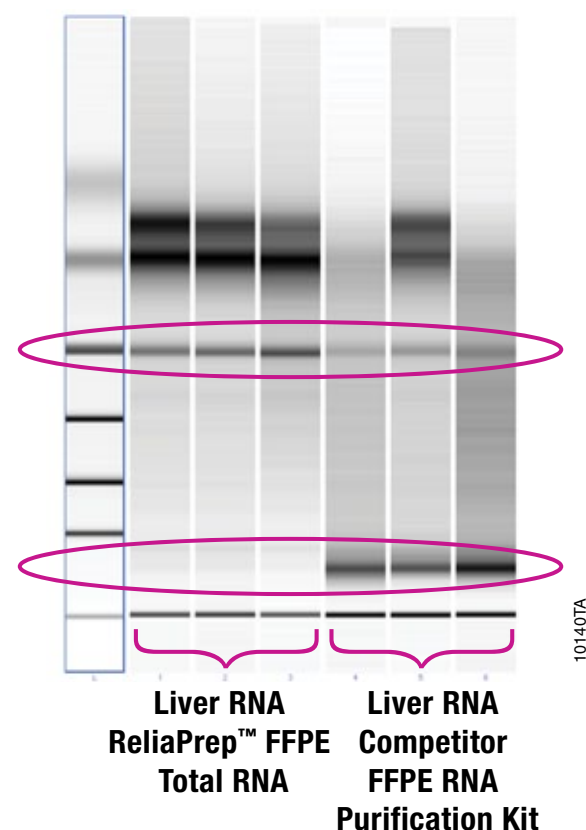
RELATED
PRODUCTS

ORDERING
INFORMATION

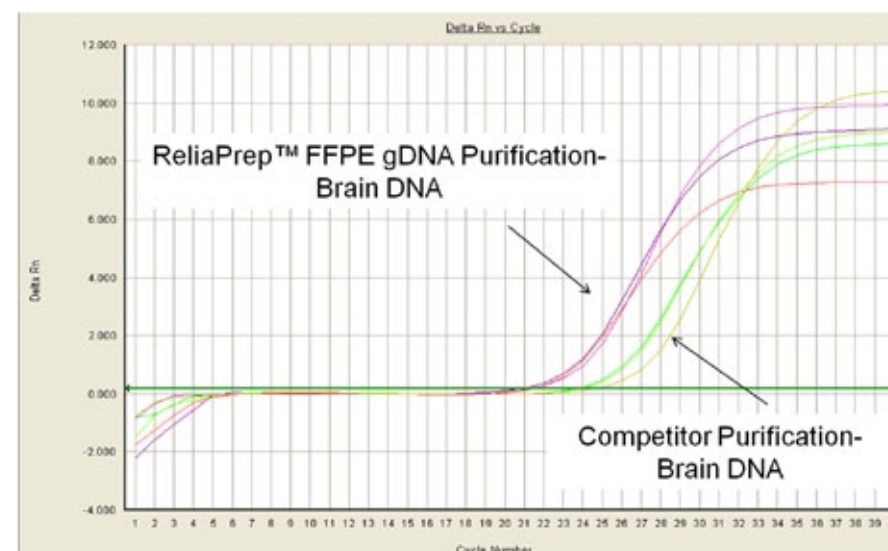
RESOURCES

Functional Yield

Yield is important, especially with precious samples. Gain more functional yield from your FFPE samples and see the difference in your downstream applications.



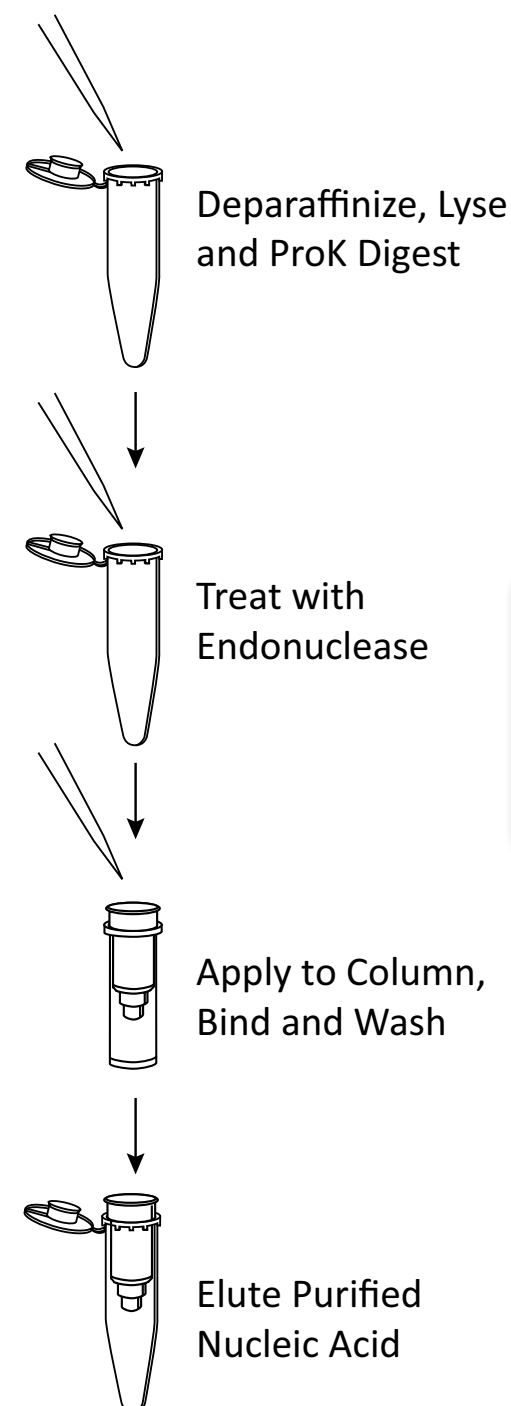
Total RNA purified from sequential 10µm mouse liver FFPE sections analyzed on an Agilent Bioanalyzer. More large-fragment RNA was purified with the ReliaPrep™ FFPE Total RNA Miniprep System than with competitor kits.



Genomic DNA purified from a single 10µm mouse brain section using either the ReliaPrep™ FFPE gDNA Miniprep System or competitor kit were analyzed using GoTaq® qPCR System. Earlier C_q values were obtained with the ReliaPrep™ FFPE System.

Workflow

- Safely deparaffinize, lyse, and ProK digest your sample without overnight digestions or harsh organic solvents
- DNase or RNase treatment occurs in the lysis buffer, minimizing unwanted nucleic acid carryover
- Quickly and efficiently bind and wash your sample
- Elute purified DNA or RNA and begin your nucleic acid analysis



Total time:
DNA: ~2 ½ hr
RNA: ~1 ½ hr

PROMEGA ISLIDES

FEATURES
AND BENEFITS

NUCLEIC
ACID INTEGRITY

FUNCTIONAL YIELD

WORKFLOW

RELATED
PRODUCTS

ORDERING
INFORMATION

RESOURCES



RELIAPREP™ FFPE MINIPREP SYSTEMS

Related Products

PRODUCT	SIZE	CATALOG NUMBER
Maxwell® 16 FFPE Plus LEV DNA Purification Kit	48 preps	AS1135
Maxwell® 16 LEV simplyRNA Cells Kit	48 preps	AS1270
Maxwell® 16 LEV simplyRNA Tissue Kit	48 preps	AS1280
Maxwell® 16 Instruments	Various	See online catalog
ReliaPrep™ Blood gDNA Miniprep System	100 preps	A5081
	250 preps	A5082
ReliaPrep™ gDNA Tissue Miniprep System	100 preps	A2051
	250 preps	A2052
GoTaq® Amplification Family	Various	See online catalog
GoTaq® MDx Hot Start Polymerase	Various	See online catalog
GoScript™ Reverse Transcription System	50 reactions	A5000
	100 reactions	A5001
Wizard® SV Gel and PCR Clean-Up System	Various	See online catalog
SV Total RNA Isolation System	Various	See online catalog
Recombinant RNasin® Ribonuclease Inhibitor	2,500u	N2511
	10,000u	N2515

PROMEGA ISLIDES

FEATURES
AND BENEFITS

NUCLEIC
ACID INTEGRITY

FUNCTIONAL YIELD

WORKFLOW

RELATED
PRODUCTS

ORDERING
INFORMATION

RESOURCES



RELIAPREP™ FFPE MINIPREP SYSTEMS

Ordering information

PRODUCT	SIZE	CATALOG NUMBER
ReliaPrep™ FFPE gDNA Miniprep System	10 reactions	A2351
	100 reactions	A2352
ReliaPrep™ FFPE Total RNA Miniprep System	10 reactions	Z1001
	100 reactions	Z1002

FEATURES
AND BENEFITS

NUCLEIC
ACID INTEGRITY

FUNCTIONAL YIELD

WORKFLOW

RELATED
PRODUCTS

ORDERING
INFORMATION

RESOURCES



Resources

The Technical Manuals provide detailed information to guide you through the purification process. The Quick Protocols give concise, step-by-step instructions for purification.

ReliaPrep™ FFPE gDNA Miniprep System
Technical Manual and Quick Protocol

ReliaPrep™ FFPE Total RNA Miniprep System
Technical Manual and Quick Protocol

Inquiry about ReliaPrep™ FFPE
Miniprep Systems:



TechServ@promega.com

GoTaq, Maxwell, RNasin and Wizard are registered trademarks of Promega Corporation. GoScript and ReliaPrep are trademarks of Promega Corporation. Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

