



FormaPure DNA FFPE Extraction Kit

SAMPLE TYPE (FORMALIN-FIXED, PARAFFIN-EMBEDDED TISSUE)

Increasingly, Formalin-Fixed, Paraffin-Embedded (FFPE) samples are an essential part of many next-generation sequencing (NGS) sample prep workflows, specifically in the field of cancer research. High quality extraction protocols focused on yield and purity are essential to achieving reliable results. FormaPure DNA now provides an automation-ready, SPRI-based extraction reagent kit to support your evolving research needs.

Most notably, it provides significant time savings in the form of reduced turnaround times and less than 15 minutes of hands-on time when automated on a Biomek liquid handler.

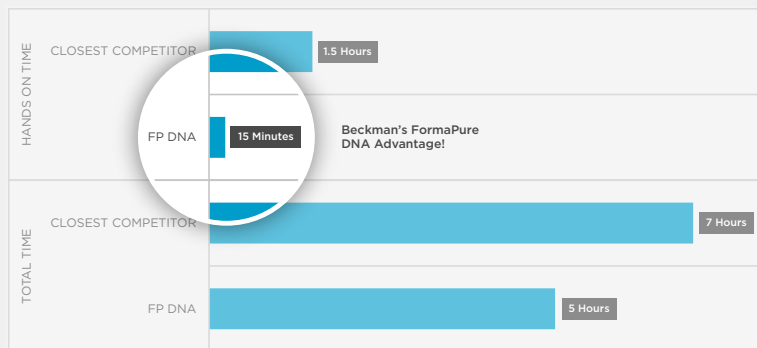
- Minimal hands-on time
- Dramatically increase productivity through automation - low to high throughput
- Proven compatibility with NGS and other downstream applications
- Consistent extraction performance
- Enhanced lab safety



Kit Components

- Mineral Oil
- Wash
- Lysis
- RNase A
- Bind
- Proteinase K

TURBO CHARGE THE FFPE WORKFLOW: 48 SAMPLES ON BIOMEK 4000



Our turbo charged workflow offers six times the productivity of the closest competitor while providing the purity, yield and downstream application compatibility needed to meet today's FFPE workflow challenges.

Figure 1. Comparison of Hands-On Time and Total Time for FormaPure DNA and the closest competitor FFPE DNA extraction kit.

NEXT-GENERATION SEQUENCING COMPATIBILITY

POSITION	GENE	FORMAPURE DNA SAMPLES WITH VARIANT CALL (MAX 6)	FORMAPURE DNA VARIANT FREQ (%)	EXPECTED ALLELIC FREQ (%)
140453136	BRAF	6	12.2	10.5
55599321	KIT	6	9.2	10
55242464	EGFR Δ	6	1.8	2
55259515	EGFR	6	3.1	3
55241707	EGFR	6	25	24.5
25398281	KRAS	6	13.7	15
25398284	KRAS	6	5.3	6
115256530	NRAS	6	17.2	12.5
178952085	PIK3CA	6	17.5	17.5
178936091	PIK3CA	6	9.1	9

Table 1. Extraction with FormaPure DNA supports detection of low frequency variants.

FormaPure DNA maintains detectability of low frequency variants without introducing extraction based experimental bias. DNA extracted using FormaPure DNA is suitable for the most stringent downstream applications including Next-Generation Sequencing.

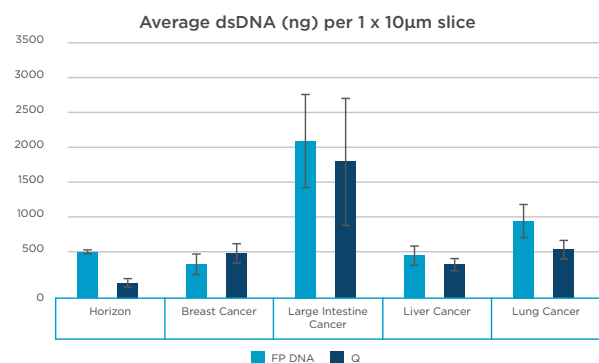


Figure 2. Average (n=8) yield per 10 µm slice. FormaPure DNA (FP DNA) compared to competitor (Q) FFPE DNA extraction kit. Multiple formalin fixed paraffin embedded (FFPE) cancer tissues were compared, including the Horizon Discovery- HD200 FFPE reference tissue.

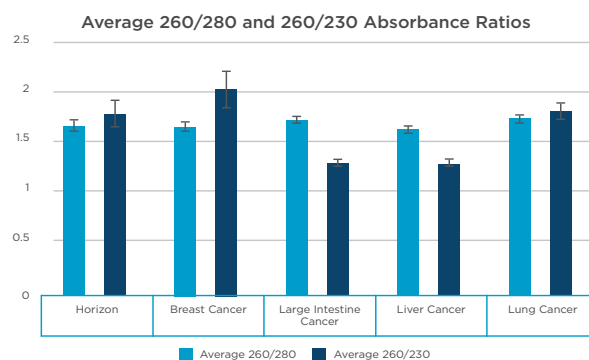


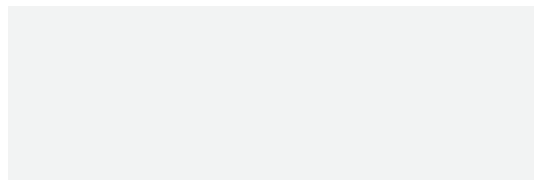
Figure 3. Average (n=8) 260/280 absorbance ratios per 10 µm slice from FormaPure DNA (FP DNA) extraction method. Multiple formalin fixed paraffin embedded (FFPE) cancer tissues were evaluated, including the Horizon Discovery- HD200 FFPE reference tissue.

FormaPure DNA is available in three kit sizes based on your throughput needs. Contact your local sales representative or visit beckman.com to request a quote.

PRODUCT ORDERING INFORMATION

B89230	Formapure DNA-50 Preps
B89231	Formapure DNA-96 Preps
B89232	Formapure DNA-384 Preps

For more information, please contact:



© 2016 Beckman Coulter, Inc. All rights reserved.
Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com

AAG-2041DS10.16