

High Recovery of DNA for Downstream Applications

Agencourt Genfind v2 System Genomic DNA Isolation Kit

The Agencourt Genfind v2 DNA purification system isolates and purifies genomic DNA (gDNA) from whole blood and serum. Additional protocols are available for extracting genomic DNA from cultured eukaryotic cells, bacteria, FTA cards, and fresh or frozen tissue¹. The kit is powered by cutting-edge SPRI (Solid Phase Reversible Immobilization) paramagnetic bead-based technology to effectively produce a high recovery of DNA for downstream applications such as PCR[†] and genotyping. SPRI allows for fast separation, easy manipulation and simple automation compared to traditional filtration technologies. The method can be run manually in a 2 mL tube format or 96-well format, or automated in 96-well format on the Beckman Coulter Biomek NX^P or FX^P workstations.

Key Features:

- Whole blood processing from 50 to 400 μ L volumes
- High gDNA recovery up to 6 μ g from 200 μ L of whole blood
- No centrifugation or vacuum filtration required
- Biomek Span-8 and 96-Multi-channel methods are available for whole blood
- Automated extraction of 96 samples in approximately 2.5 hours

Consistent and High Recovery of gDNA

Figure 1 demonstrates that Agencourt Genfind v2 delivers high quality and consistent yields in comparison to competitor filtration and bead-based methods. The unique properties of the SPRI technology allows for the effective capture of gDNA in the presence of various anticoagulants.

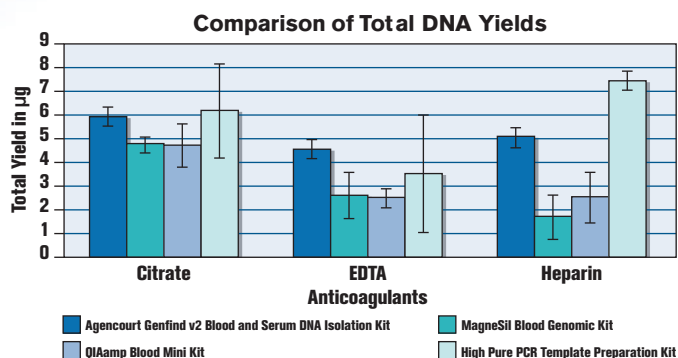


Figure 1. Replicates of eight 200 μ L samples of frozen human blood were manually extracted and eluted in 200 μ L H₂O (High Pure was eluted in 10 mM tris, pH 8.5); for all kits analyzed, aliquots were taken from the same patient sample for each anticoagulant used. DNA concentrations and yields were determined using the Beckman Coulter DTX plate reader to measure the A260 absorbances.

DS-17164A

Table 1 - Ct Values

Sample	Mean Ct
Agencourt Genfind v2 10 ng	21.50 \pm 0.37
Agencourt Genfind v2 1 ng	24.19 \pm 0.17

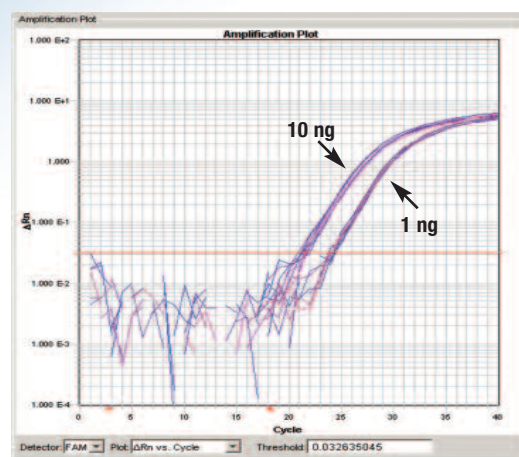


Figure 2. Ct values for nucleic acid purified using the Agencourt Genfind v2 system.

Ct values obtained in qPCR reactions indicate that samples purified by the Agencourt Genfind v2 system produce amplifiable gDNA with a high sample-to-sample consistency (Table 1, Figure 2).

High Quality gDNA

The Agencourt Genfind v2 DNA isolation method isolates and purifies gDNA that is intact and suitable for downstream PCR and genotyping. Genomic DNA from human whole blood samples was isolated with Agencourt Genfind v2 and analyzed by gel electrophoresis (Figure 3A).

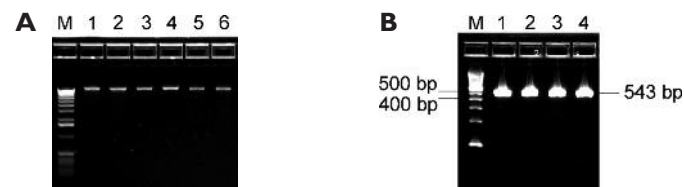


Figure 3A. Gel electrophoresis of 1 μ L genomic DNA through a 0.8% agarose gel. DNA was isolated from 200 μ L blood with the Agencourt Genfind v2 purification method and eluted in 200 μ L H₂O. Key: M = 1 kb ladder, lanes 1 & 2 = Citrate, lanes 3 & 4 = EDTA, and lanes 5 & 6 = Heparin.

Figure 3B. PCR for the human ADP ribosylation factor 1 (ARF1) gene was performed using 2 μ L of extracted gDNA in a 20 μ L reaction and 10 μ L of each reaction was resolved through a 4% agarose gel. Key: M = 100 bp ladder, lanes 1 – 2 = EDTA, and lanes 3 – 4 = Heparin.

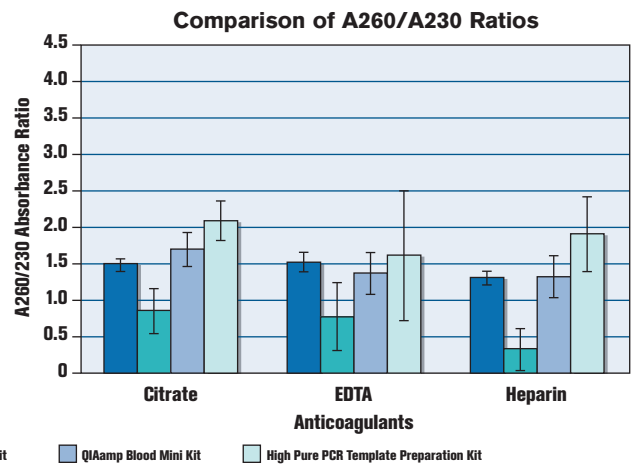
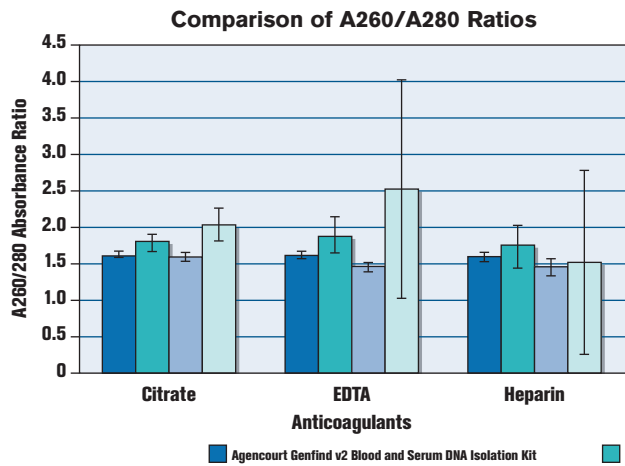


Figure 4. Absorbances were determined using 2 μ L for Nanodrop* analysis. Replicates of 8 were performed for each anticoagulant sample.

No Inhibition from Anticoagulants

Heparin can be a powerful PCR inhibitor that can make DNA amplification a challenge. Agencourt Genfind v2 efficiently removes common anticoagulants from blood such as Citrate, EDTA and Heparin. Figure 3B shows that the DNA from a blood sample containing Heparin amplifies as easily as the DNA from blood containing EDTA.

High Purity

Contaminants, such as salt and protein, can negatively affect downstream application of genomic DNA. As seen in Figure 4 Agencourt Genfind v2 produces gDNA with consistent A260/A280 and A260/A230 ratios improving success in downstream applications.

Summary

The Agencourt Genfind v2 system is efficient and automation-friendly. By automating the process using Agencourt software methods on either the Biomek Span-8 or 96 Multi-channel workstation, gDNA can be isolated and purified in approximately 2.5 hours. With the power of SPRI chemistry, the Agencourt Genfind v2 system produces highly pure DNA and consistent data thus allowing researchers to minimize retesting of precious samples.

Kit Components

- Lysis Buffer
- Wash I Buffer
- Wash II Buffer
- Binding Buffer
- Proteinase K
- Proteinase K Buffer



Ordering Information

For more information, please visit our website at www.beckmancoulter.com or contact your local sales representative.

Product	Size	Product #
Agencourt Genfind v2 DNA Isolation 2 mL Tube Kit	50 ² preps, 400 μ L	A41499
Agencourt Genfind v2 DNA Isolation 96-well Plate Kit	384 ² preps (4 x 96), 200 μ L	A41497
Agencourt Genfind v2 96 Batch Software Method, v3.x		A42568
Agencourt Genfind v2 Span-8 96 Batch Software Method, v3.x		A42569

Related Products	Size	Product #
Agencourt AMPure PCR Purification 60 mL Kit	1333 preps (25 μ L PCR reaction volume)	A29152

1. Please inquire with your local sales representative for these additional protocols.
2. Based on whole blood sample.

For Research Use Only. Not for use in diagnostic procedures.

† The PCR process is covered by patents owned by Roche Molecular Systems, Inc., and F. Hoffman-La Roche, Ltd.

* All trademarks are property of their respective owners.

Beckman Coulter, the stylized logo, Agencourt, AMPure, Biomek, Genfind, and SPRI are trademarks of Beckman Coulter, Inc. and are registered with the USPTO.

