

NucleoMag[®] Blood 200 μ L

Automated purification of genomic DNA from 200 μ L blood samples on the platform KingFisher[®] Flex



Introduction

The isolation of genomic DNA from whole blood samples is an initial and crucial step for various diagnostic workflows. Clinical applications such as genotyping, HLA typing, biomarker discovery, newborn screening, and pharmacogenetics are widely performed in laboratories worldwide. To enable a state of the art molecular diagnostic, the sensitivity and performance of biomolecular detection methods like qPCRs, next-generation sequencing, and microarray analysis, is constantly improved. A main aspect to face these workflow requirements, is to facilitate the extraction of highly pure DNA in substantial amounts. To provide a fast and consistent sample processing, MACHEREY-NAGEL designed the NucleoMag[®] Blood 200 μ L kit, specialized for automated purification of genomic DNA from 200 μ L whole blood samples (fresh or frozen, EDTA or citrate treated) in a 96-well format. The optimized protocol allows the processing of 96 samples within 45 minutes including a sample lysis on the KingFisher[®] platform.

Product at a glance

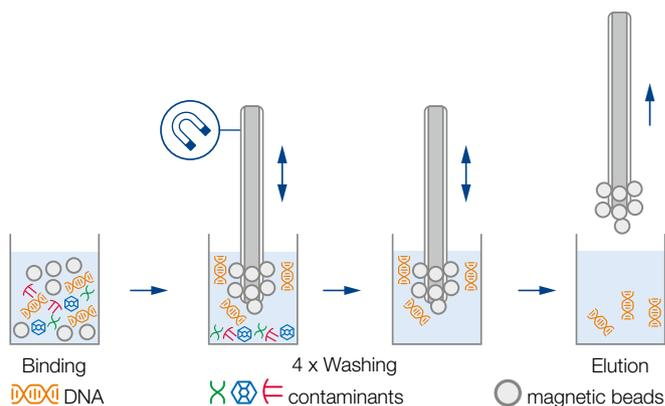
NucleoMag [®] Blood 200 μ L	
Technology	Magnetic bead technology
Sample material	\leq 200 μ L whole blood (fresh or frozen, EDTA or citrate treated)
Preparation time	Approx. 45 min on KingFisher [®] Flex for 96 samples incl. sample lysis
Typical yield	2–8 μ g (200 μ L blood; depending on sample quality)
Elution volume	50–100 μ L
Binding capacity	0.4 μ g/ μ L beads
King Fisher [®] Flex	
Sample volume	20–5000 μ L
Capacity	24 / 96 samples (8 plates per deck)
Heating / cooling	4–96 °C
Size / weight	60 x 38 x 68 cm / 28 kg

Material and Methods

Whole blood (fresh, frozen, treated with EDTA or citrate) is lysed at room temperature with Lysis Buffer MBL1 and Proteinase K. Following lysis incubation, binding of DNA to the NucleoMag[®] B-Beads is achieved by the provision of Binding Buffer MBL2.



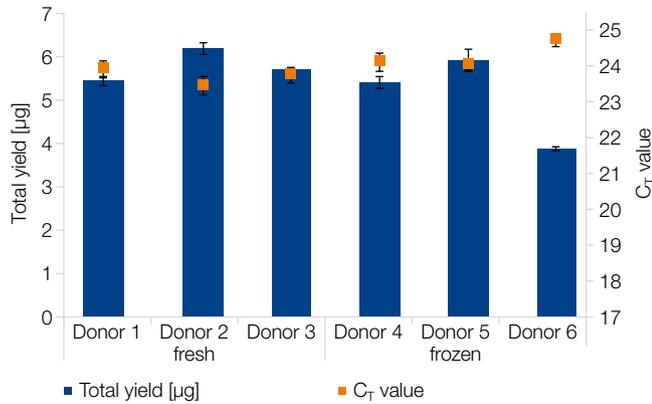
Subsequent DNA isolation is performed on the automation platform KingFisher[®] Flex. The isolation principle is based on reversible adsorption of nucleic acids to paramagnetic NucleoMag[®] B-Beads under appropriate buffer conditions.



Workflow on automation platform

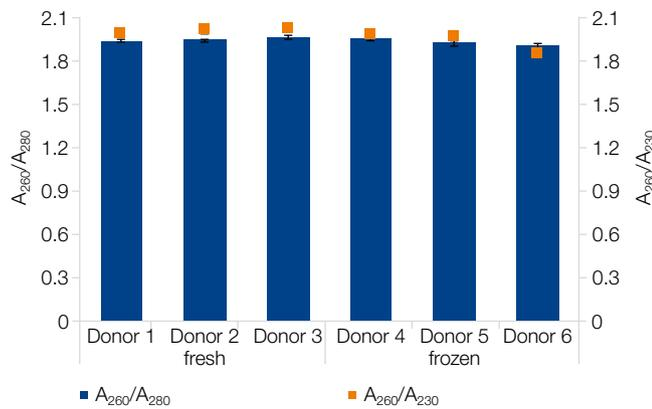
After magnetic separation, the NucleoMag[®] B-Beads are washed four times to remove contaminants and salts using three different wash buffers (MBL3, 80 % ethanol and MBL4). Highly pure genomic DNA is finally eluted under low ionic strength conditions in a slightly alkaline elution buffer (MBL5).

Application data



Isolation of genomic DNA from fresh and frozen human blood samples

DNA was isolated from fresh and frozen 200 µL human blood samples (n = 8) using the NucleoMag® Blood 200 µL kit on a KingFisher® Flex platform. The total yield was determined by UV-spectrometry (dark blue bars). A subsequent qPCR analysis (orange squares) was performed with a Taqman® Probe for a 250 bp β-Actin amplicon using the SensiFast™ Probe Lo-ROX kit from Bioline on an Applied Biosystems® 7500 Real-Time PCR System.



Purity of isolated nucleic acids from fresh and frozen human blood samples

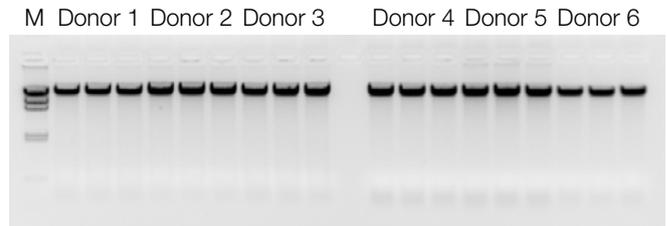
DNA was isolated from fresh and frozen 200 µL human blood samples (n = 8) using the NucleoMag® Blood 200 µL kit on a KingFisher® Flex platform. The purity was determined by UV-spectrometry resulting into an average A₂₆₀/A₂₈₀ value of 1.93 ± 0.2 (dark blue bars) and into an average A₂₆₀/A₂₃₀ value (orange squares) of 1.96 ± 0.05.

Ordering information

Product	Specifications	Preps	REF
NucleoMag® Blood 200 µL	Kit based on magnetic bead technology for the isolation of genomic DNA from 200 µL blood samples including NucleoMag® B-Beads, buffers, and Proteinase K	1 x 96	744501.1
		4 x 96	744501.4
KingFisher® Accessory Kit B*	Deep-well Plates, Deep-well Tip Combs, and Elution Plates for 4 x 96 NucleoMag® Blood 200 µL preps using the KingFisher® Flex platform	4 x 96	744951

*For use on KingFisher® Flex

NucleoMag® is a registered trademarks of MACHEREY-NAGEL; KingFisher® and Applied Biosystems® are registered trademarks of Thermo Scientific Inc.; SensiFast™ is a trademark of Bioline Reagents Ltd.



Integrity of isolated nucleic acids

The integrity of the isolated nucleic acids from fresh (donor 1-3) and frozen (donor 4-6) blood samples was analyzed by gel electrophoresis (10 µl per eluate; 1 % TAE-gel; M: Lambda DNA/Hind III – Thermo Scientific).

Automate your DNA extraction from blood samples

MACHEREY-NAGEL delivers a ready to go solution for your high throughput DNA extraction from various blood samples, such as fresh, frozen, treated with EDTA or citrate. We adapted the NucleoMag® Blood 200 µL kit on instruments of the KingFisher® series to speed up your nucleic acid purification workflow.

- Reliable performance and excellent yields from various blood sample material
- Speed up your DNA extraction by processing of 96 samples in less than 45 minutes (including sample lysis)