Automated DNA/RNA Isolation

Automated Nucleic Acid Isolation Fulfilling Biobanking Needs



Biobanking applications, biospecimen repositories and Mircrobiome research projects are facing demanding and diverse sample materials like whole blood, FFPE-tissue to feces-material in all kind of sample volumes and conditions. To fulfill the needs of sensitive genetic research applications such as whole genome sequencing, real time/digital PCR, and methylation analysis etc., the extraction of high quality and high yield nucleic acids is an essential initial step, especially when it comes to long-term storage of the extracted DNA/RNA samples as required in Biobanking.

PerkinElmer offers the perfect solutions for this challenge by providing

- Automated Nucleic Acid Isolation, highly flexible in sample volume (10 µl 10 ml), sample material and throughput
- Huge variety of processable sample materials (blood, buffy coat, saliva, tissue, feces, etc.)
- Maximum yields and best purities (up to 50 µg DNA/ml blood; OD_{260/280} 1.9; OD_{260/230} >2.0)
- Long term stable nucleic acids

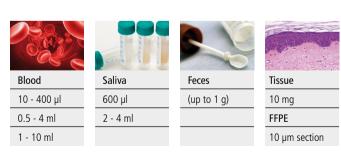


Fig. 3: Protocols optimized for Biobanking

Further chemagic Kits are available on request.

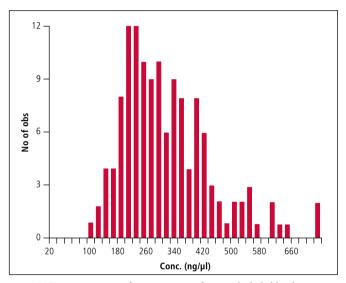


Fig. 1: DNA concentration of 133 extractions from 4 ml whole blood, elution vol. 500 μ l. Mean conc.: 314 ng/ μ l, mean yield: 157 μ g

Genomic DNA isolation from \leq 7 ml human blood (with chemagic DNA Blood Kit special CMG-715), elution volume = 500 μ l

Achieving high average yields: $63 \mu g/ml$ blood, DNA concentration: $346.6 ng/\mu l \pm 163.01$ High DNA quality: $A_{260}/A_{260} = 1.98 \pm 0.26$

- > 2,500 DNA extractions per year
- > 200 runs per year



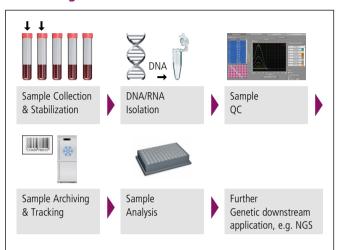
Fig. 2: Customer results of Biobank HCB-IDIBAPS, Barcelona, Spain, Blood and Fluid Bank, Veronica Fernandez Pascual, PhD



Your Tool for DNA/RNA Isolation Workflows in Biobanking

Based on your automation needs PerkinElmer offers full workflow solutions covering high throughput DNA/RNA isolation, liquid handling robotics, Nucleic Acid quality control and analysis, optimized for Biobanking and Biorepository demands.

Biobanking Workflow



All from a single-source PerkinElmer delivers Biobanking application support from DNA to Data.

Next Generation Sequencing Workflow



For more information please visit www.perkinelmer.de/category/genomic-analysis

For Research Use Only. Not for use in Diagnostic Procedures. US: For Laboratory Use Only. Not Intended For Use in Diagnostic Procedures.

PerkinElmer, Inc.

940 Winter Street Waltham, MA 02451 USA Phone: (+1) 800-762-4000 or (+1) 203-925-4602

www.perkinelmer.com

PerkinElmer chemagen Technologie GmbH

Arnold-Sommerfeld-Ring 2 52499 Baesweiler, Germany Phone: (+49) 2401 805-501 Fax: (+49) 2401 805-519

support.chemagen@perkinelmer.com, www.chemagen.com



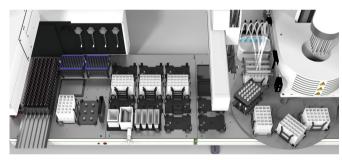
chemagicTM 360 Nucleic Acid Extractor

Key features

- Sample volumes from 10 μl 10 ml
- High throughput
- LIMS compatible log files, barcode reading

Benefits

- Ready to use DNA/RNA up to 50 μg DNA/ml blood
- Compact benchtop design
- Integrates with Liquid Handling Systems



chemagic 360
i Nucleic Acid Extractor integrated in Perkin Elmer JANUS
° Liquid Handling Workstation

Automation solutions

- High throughput primary sample transfer and DNA/RNA Isolation
- Optional: normalization, PCR setup applications, eluate transfer into storage formats
- Numerous sample tube and plate formats for primary sample input and eluate output
- Barcode sample and eluate tracking for seamless integration to LIMS systems, sample security and storage solutions

