

# Automated nucleic acid purification.



revvity

chemagic™ Technology

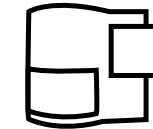
For research use only. Not for use in diagnostic procedures.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Established chemagic technology

## Reliable automated nucleic acid purification

As one of the pioneers of magnetic bead-based nucleic acid purification, chemagic technology has been meeting the rigorous demands of laboratories world-wide for over 25 years. Combining proprietary M-PVA Magnetic Beads<sup>1</sup> with advanced automation, chemagic technology provides reliable nucleic acid integrity with proven success in challenging downstream assays. The flexible, robust, and high-end automation caters to diverse workflow needs while ensuring sample integrity, reliable performance and streamlined lab efficiency. As a trusted partner in diagnostics, biobanking and scientific research, benefit from our personalized technical support and extensive experience in handling diverse human sample matrices. For any enquiries, please contact us at [support.chemagen@revvity.com](mailto:support.chemagen@revvity.com).



> 2000 chemagic Instruments



Installed Over 84 Countries



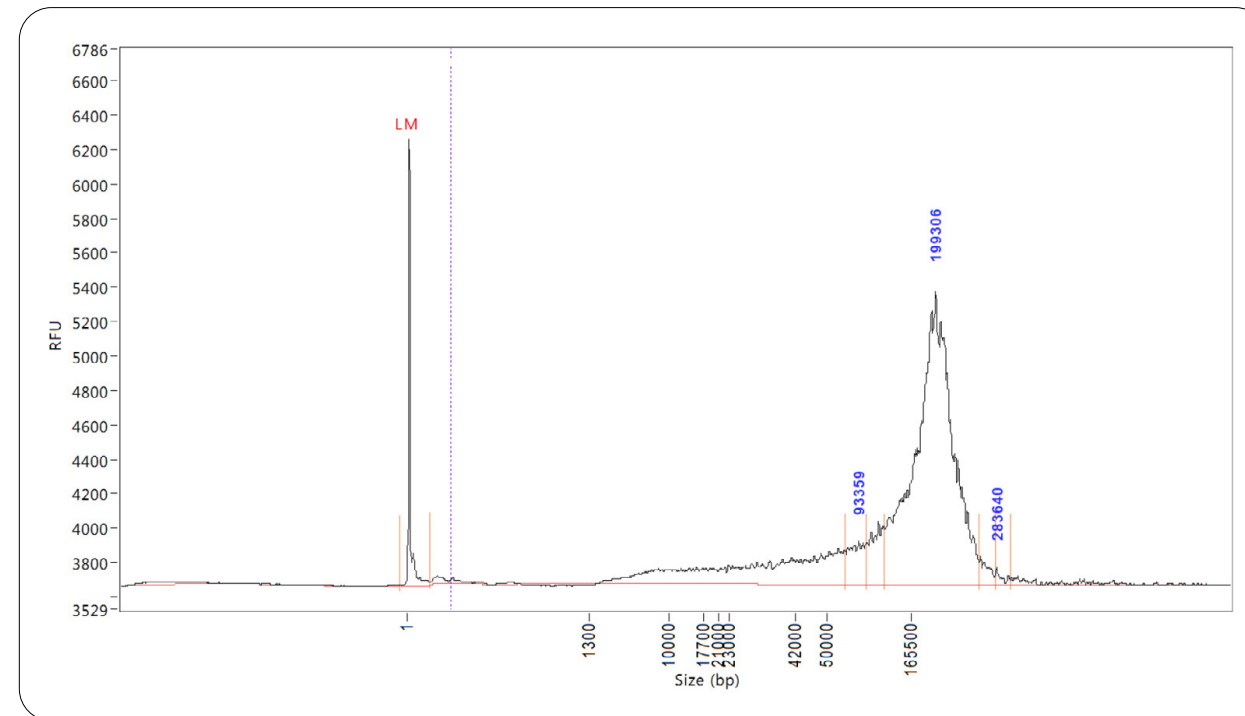
With > 2500 Citations



# Advantages of our technology

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

High nucleic acid integrity (sizes up to ~ 200 kb)  
ideal for long-term storage and challenging downstream assays



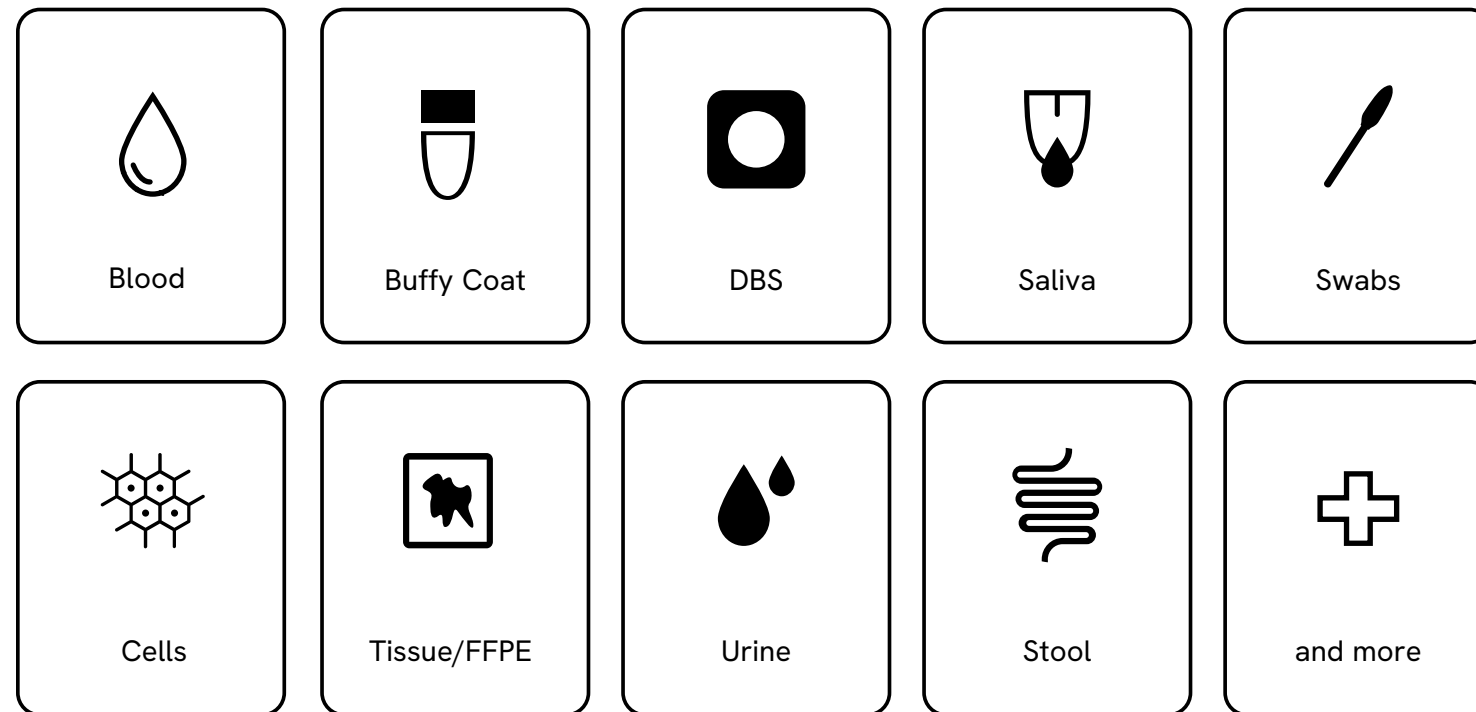
DNA is not degraded during extraction. DNA was extracted from 4 mL samples of whole blood using the chemagic 360 instrument and diluted 1:1000 for analysis of molecular weight using an Agilent® FEMTO Pulse™ system.



# Advantages of our technology

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Flexible sample volume inputs (50  $\mu$ l to 18 ml) and ability to process diverse sample matrices.





# Advantages of our technology

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

No detectable cross-contamination and carry-over of inhibitors from efficient M-PVA Magnetic Bead handling.

	1	2	3	4	5	6	7	8	9	10	11	12
A	16.92	undet.	16.73	undet.	17.44	undet.	16.58	undet.	16.65	undet.	16.63	undet.
B	undet.	16.49	undet.	16.83	undet.	16.99	undet.	16.78	undet.	16.93	undet.	16.74
C	16.57	undet.	16.96	undet.	16.85	undet.	16.94	undet.	16.74	undet.	17.13	undet.
D	undet.	16.87	undet.	16.64	undet.	17.45	undet.	17.59	undet.	17.18	undet.	16.98
E	16.68	undet.	16.92	undet.	17.65	undet.	17.53	undet.	16.85	undet.	16.94	undet.
F	undet.	17.39	undet.	17.01	undet.	17.04	undet.	17.16	undet.	16.89	undet.	16.71
G	16.92	undet.	16.73	undet.	17.44	undet.	16.58	undet.	16.65	undet.	16.63	undet.

- milliQ water
  - 5 µl HSV positiv direct control



# Advantages of our technology

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Reduced hands-on time and running costs from optimal buffer dispensing system.



Technical support and consultation available for verification of sample types and workflow requirements.





- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Solutions to fulfill your laboratory needs

Whether you are processing one or multiple sample types in a single workflow, have low or high sample volumes, or require specific extraction conditions, there is a chemagic solution that will meet your laboratory demands. Benefit from universal kit options for diverse sample types, single and multi-sample type specific kits to optimize for runtimes and cost, or custom kits to fulfil specific requirements.

The chemagic technology has been applied successfully in diverse applications ranging from biobanking, molecular genetic analysis, donor screening, oncology and infectious disease research.





- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Genetic analysis and biobanking

## Genomic DNA from diverse sample types for genetic workflows and biobanking

The high, archival integrity of nucleic acids obtained with chemagic technology is ideal for biobanks and research laboratories. The chemagic separation technology automates the efficient extraction of high molecular weight (HMW) DNA, enabling successful application in diverse downstream assays for genetic analysis such as HLA typing<sup>2</sup> for donor screening, SNP arrays for pharmacogenetics<sup>3</sup>, MLPA for copy number detection<sup>4</sup>, NGS from dried blood spots for newborn research<sup>5</sup> and long read sequencing in genetic research<sup>6</sup>

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Diverse sample material	Up to 400 µl blood, 500 µl buffy coat, 600 µl saliva, DBS, stool, and FFPE samples	chemagic Universal H96 Kit	CMG-718
gDNA	Blood, buffy coat, saliva, bone marrow cells	Up to 4 ml	chemagic DNA Blood 4K Kit H24	CMG-1074

[Read more](#)

"More than 10 years ago, we chose chemagen instruments and reagents for isolating DNA from blood or buccal swabs. After performing HLA genotyping on over 10 million samples, we remain extremely satisfied with this decision and the consistent performance."

- Dr. Vinzenz Lange, CTO, DKMS Life Science Lab, Germany





- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Precision oncology

## cfDNA from liquid biopsies

The unique large volume processing capability of chemagic instruments and high specificity of M-PVA Magnetic Beads combine to boost extraction yields of cfDNA from liquid biopsies, increasing chances of detection downstream. Easily scale from low to high throughputs and manual to automated kit options. chemagic cfDNA extraction has been used to support the clinical research of colorectal<sup>7</sup>, lung<sup>8</sup>, prostate and anal cancer<sup>9</sup> with downstream methods including digital PCR and pan-cancer NGS panels.

Target	Sample type	Sample volume	Product name	Product no.
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	3 - 5 ml	chemagic cfDNA 5k Kit H24	CMG-1304
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	6 - 10 ml	chemagic cfDNA 10k Kit H24	CMG-1310
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	1 - 5 ml	chemagic cfDNA 5k Kit	CMG-134

[Read more](#)

“We use cfDNA extracted with chemagic™ technology in digital droplet PCR (ddPCR) assays to support research of colorectal cancer, lung cancer, and anal cancer (based on HPV cfDNA detection) with a sample to result turnaround time down to 24 h.”

- Professor Niels Pallisgaard, Dept of Pathology, Zealand University Hospital, Denmark



- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

## DNA from frozen/FFPE tissue, blood, or bone marrow

To support the growing demand, nucleic acid extraction from blood, bone marrow cells, or tumor tissue that is either fresh, frozen, or formalin-fixed, paraffin-embedded (FFPE) is often performed. The chemagic kits consistently produce high-quality, NGS grade DNA and RNA from these sample types and more. Commonly cited downstream applications include RT-PCR, methylation arrays, MLPA and pan-cancer NGS panels<sup>10,11</sup>.

Target	Sample type	Sample volume	Product name	Product no.
gDNA	FFPE tissue	10 µm section	chemagic FFPE DNA Kit H96	CMG-1099
gDNA	Blood, buffy coat, saliva, bone marrow	Up to 400 µl	chemagic DNA Blood 400 Kit H96	CMG-1091
gDNA	Diverse sample material	Up to 400 µl blood, 500 µl buffy coat, 600 µl saliva, stool and FFPE samples	chemagic Universal H96 Kit	CMG-718

[Read more](#)




- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Translational research and biomarker development

## RNA from cells, tissue, and blood

The chemagic RNA isolation technology has been applied to various sample types including blood, tissues, and cells and provides researchers consistent high integrity RNA suitable for a variety of downstream applications including RNA sequencing<sup>12</sup>.

Target	Sample type	Sample volume	Product name	Product no.
RNA	Tissue, cells from cell cultures, blood	Up to 2.5 ml blood or 10 mg tissue	chemagic RNA Kit H96	CMG-1212*
RNA	Blood from PAXGene Tubes	Up to 2.5 ml blood	chemagic RNA PGBlood Kit H96	CMG-1219

\*CMG-1212 is a universal kit solution for RNA while CMG-1219 is an optimized kit solution for PAXGene blood tubes to maximize for yield and performance.

[Read more](#)



- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

## miRNA from plasma, serum, or exosomes

Despite the small size and low concentrations of cell free circulating microRNA (miRNA), the chemagic miRNA kits obtain high, consistent yields comparable to manual methods with only 10 minutes of hands-on time. Suitable for plasma, serum, extracellular vesicles and saliva, miRNAs obtained have been analyzed by rtqPCR and NGS.

Target	Sample type	Sample volume	Product name	Product no.
miRNA (cfRNA)	Plasma, serum, exosomes, saliva	Up to 200 $\mu$ l	chemagic miRNA 200 Kit H96	CMG-1224
miRNA (cfRNA)	Plasma, serum, exosomes, saliva	1 ml	chemagic miRNA 1K Kit H24	CMG-1223

[Read more](#)

"We have started the testing of cfRNA isolations with our chemagic 360 instrument - the RNA yields appear fine from different plasma and serum samples; thus we look forward to use the kit for high throughput applications."

- Dr. Maija Puhka, Head of HiPrep and EV Core, Institute for Molecular Medicine



- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Infectious disease research

## Pathogen nucleic acids from body fluids

Fast turn-around times, diverse sample type processing and high, consistent yields make chemagic nucleic acid kits with chemagic instruments the choice for many research labs working on infectious diseases. Automated extraction workflows from gram positive and negative bacteria, intra and extracellular viruses, yeast, fungi, and protozoa have been performed from body fluids like saliva, sputum, serum, plasma, lavage, urine, semen, amniotic fluid as well as swabs in various transport media. Challenging sample types such as stool, ticks, cultures, and wastewater have also been processed. Applications cited have been wide-ranging from blood-borne parasite screening to wastewater surveillance of infectious agents<sup>13,14,15</sup>.

Target	Sample type	Sample volume	Product name	Product no.
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 300 µl	chemagic Viral DNA/RNA 300 Kit H96 *	CMG-1033-S
Pathogen NA	Body fluids, blood, ticks, cultures, supernatant of stool samples	Up to 200 µl	chemagic Viral DNA/RNA 200 Kit H96	CMG-1049-A
Pathogen NA	Body fluids, blood, ticks, cultures, supernatant of stool samples	Up to 1 ml	chemagic Body Fluid 1K Kit H96	CMG-1131

\* The kit "chemagic Viral DNA/RNA 300 Kit H96" has also been used for pathogen detection from wastewater.

[Read more](#)

"The chemagic 360 instrument has been functioning well, satisfying the requirements of our COVID-19 testing (4000 samples/day) and meeting our expectations for quality and reliability."

- Department of Microbiology, Government of Rajasthan, S.M.S. Medical School, Jaipur

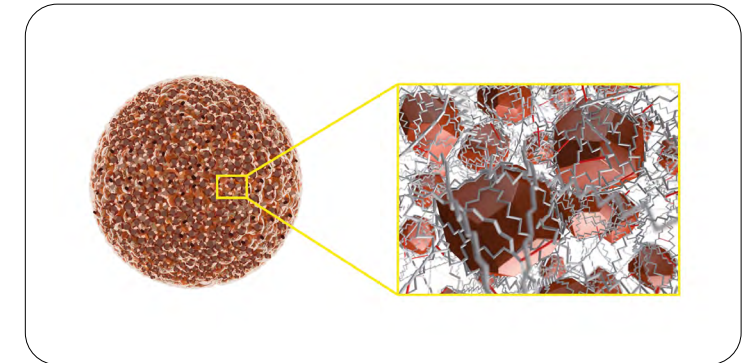
# M-PVA Magnetic Beads

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

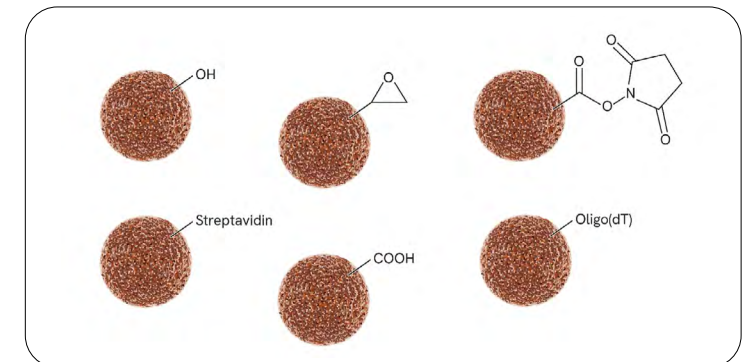
chemagic M-PVA Magnetic Beads consist of superparamagnetic nanometer sized magnetite particles encapsulated in a matrix of crosslinked polyvinyl alcohol.

Unique benefits:

- High magnetite content, enabling fast separation even from large volumes.
- Mechanically robust, facilitating use in automation, stable at RT and even autoclavable.
- Hydrophilic PVA surface sees no measurable protein binding, resulting in highly pure nucleic acids.
- Chemical functionality of the matrix enables different modifications with high binding capacity.



| Schematic image of a M-PVA Magnetic Bead



| M-PVA Magnetic Beads with functionalized groups



- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# Maximize quality with chemagic automation

## Where flexibility meets high performance

To achieve high nucleic acid integrity, the chemagic instruments employ the use of rotating magnetizable steel rods that automate resuspension of M-PVA Magnetic beads for gentle yet efficient extraction of nucleic acids. Utmost flexibility in sample volume range (50  $\mu$ l – 18 ml) and throughput, together with fast run-times, are achieved with the easily interchangeable 12-, 24- or 96-format chemagic Rod Heads of the chemagic 360 instrument and chemagic Prime™ instrument.

For low to medium throughput processing, the chemagic Prepito™ and chemagic Prime Jr instruments enable up to 12 or 48 samples per run respectively with sample volumes of up to 1 ml.



- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

chemagic automated nucleic acid purification

### Fully automated workflows

### Sample throughput

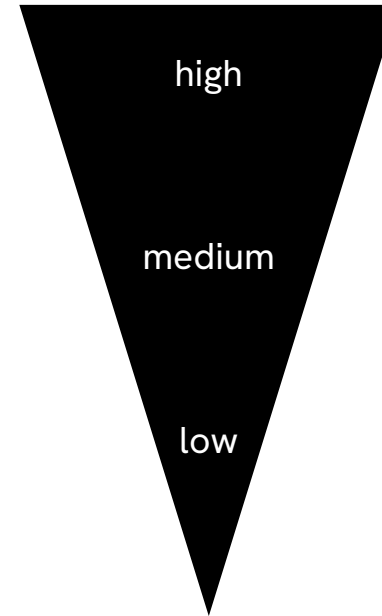
### For combined workflows



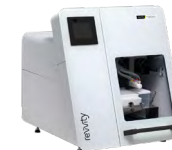
chemagic Prime Instrument



chemagic Prime Jr Instrument



chemagic 360 Instrument



chemagic Prepito Instrument

Instrument	Sample throughput	Samples per run	Product no.
chemagic Prime Instrument	High	1-192	2039-0020
chemagic 360 Instrument chemagic 360-D Instrument*	High	1-96	2024-0020 2024-0010
chemagic Prime Jr Instrument chemagic Prime Jr-D Instrument*	Medium	1-48	CMG-2029-0020 2029-0010
chemagic Prepito Instrument chemagic Prepito-D Instrument*	Low	1-12	2022-0030 2022-0020

\* Products may not be licensed in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability.





- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

# chemagic Kits

## Ready-to-use solutions for diverse needs

Catering to a wide variety of extraction targets and sample types, chemagic nucleic acid kits are stable at room temperature and produced in an ISO environment. The chemagic Kits come ready-to-use with plastic consumables included. Being a full value chain owner, benefit from a dependable supply and the ability to customize your kit specifications.

- Ability to configure kit to custom specifications.
- Stable at room temperature with a long shelf life of up to 24 months.  
In-use stability of up to 3 months allows for partial kit use.
- Ready-to-use with plastic consumables included.



## chemagic Kits on third party instruments

To benefit from M-PVA Magnetic bead technology and optimized buffer chemistries, the chemagic kits can be used with automated liquid handlers from Revvity or other providers.

[Available kits](#)

## chemagic Kits for manual workflows

For labs with smaller throughputs or just beginning to implement automation, manual chemagic Kits can be used together with magnetic stand accessories in a manual workflow.

[Manual kits](#)

[Magnetic stands](#)



# Product list

## chemagic 360 Kits for chemagic 96 Rod Head

Preps/Kit: 960 (unless otherwise stated\*) | Reaction vessel: 96 deep-well plate | All plastic consumables included.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Diverse sample material	Up to 400 µl blood, 500 µl buffy coat, 600 µl saliva, stool and FFPE samples	chemagic Universal H96 Kit	CMG-718
gDNA	Blood, buffy coat, saliva, bone marrow	Up to 400 µl	chemagic DNA Blood 400 Kit H96	CMG-1091
gDNA	Blood, buffy coat, saliva, bone marrow	Up to 800 µl	chemagic DNA Blood 800 Kit H96*	CMG-1218
gDNA	Buffy coat	Up to 200 µl	chemagic DNA Buffy Coat 200 Kit H96	CMG-713-MK
gDNA	Saliva	Up to 600 µl	chemagic DNA Saliva 600 Kit H96	CMG-1037
gDNA	Epithelial cells from the inner cheek lining	1 swab	chemagic DNA Buccal Swab Kit H96**	CMG-748
gDNA	Dried blood spot (DBS)	3 - 13 mm punch out	chemagic DNA Blood Spot 13 Kit H96	CMG-779
gDNA	FFPE tissue	10 µm section	chemagic FFPE DNA Kit H96	CMG-1099
gDNA	Tissue, cells	Up to 10 mg	chemagic DNA Tissue 10mg Kit H96	CMG-723
RNA	Tissue, cells from cell cultures, blood	Up to 2.5 ml blood or 10 mg tissue	chemagic RNA Kit H96***	CMG-1212
RNA	Blood from PAXGene Tubes	Up to 2.5 ml blood	chemagic RNA PGBlood Kit H96	CMG-1219
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	Up to 1.5 ml	chemagic cfDNA 1.5k Kit H96	CMG-1396
miRNA (cfRNA)	Plasma, serum, exosomes, saliva	Up to 200 µl	chemagic miRNA 200 Kit H96	CMG-1224

\* The "chemagic DNA Blood 800 Kit H96" is applied with the 96-Rod-Head but in 48-deep-well-plates. The amount of preps/Kit is 480.

\*\* Sample collection can be done directly or with swabs.

\*\*\* This is a universal kit for RNA extraction. An optimized kit solution for PAXGene blood tubes is also available (CMG-1219).



## chemagic 360 Kits for chemagic 96 Rod Head

Preps/Kit: 960 | Reaction vessel: 96 deep-well plate | All plastic consumables included.

Target	Sample type	Sample volume	Product name	Product no.
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 300 $\mu$ l	chemagic Viral DNA/RNA 300 Kit H96*, **	CMG-1033-S
Pathogen NA	Body fluids, blood, ticks, cultures, supernatant of stool samples	Up to 200 $\mu$ l	chemagic Viral DNA/RNA 200 Kit H96*	CMG-1049-A
Pathogen NA	Body fluids, blood, ticks, cultures, supernatant of stool samples	Up to 1 ml	chemagic Body Fluid 1K Kit H96*	CMG-1131
Pathogen NA, gDNA	Stool	Up to 1 g	chemagic DNA Stool Kit H96	CMG-1076

\* Sample collection can be done directly or with swabs.

\*\* The kits "chemagic Viral DNA/RNA 300 Kit H96" and "chemagic Viral DNA/RNA 9.6K Kit H12" can be also used for pathogen detection from wastewater (COVID).

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact



## chemagic 360 Kits for chemagic 24 Rod Head

Preps/Kit: 240 | Reaction vessel: 24 deep-well plate | All plastic consumables included.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Blood, buffy coat, saliva, bone marrow cells	Up to 2 ml	chemagic DNA Blood 2K Kit H24	CMG-1097
gDNA	Blood, buffy coat, saliva, bone marrow cells	Up to 4 ml	chemagic DNA Blood 4K Kit H24	CMG-1074
gDNA	Buffy coat	Up to 500 µl	chemagic DNA Buffy Coat 500 Kit H24	CMG-1080
gDNA	Epithelial cells from the inner cheek lining	Up to 3 swabs	chemagic DNA Buccal Swab Kit H24*	CMG-1048
gDNA	Saliva	Up to 4 ml	chemagic DNA Saliva 4K Kit H24	CMG-1081
gDNA	Tissue, cells	Up to 100 mg	chemagic DNA Tissue 100mg Kit H24	CMG-1207
RNA	Tissue, cells from cell cultures, blood	Up to 2.5 ml blood or 10 mg tissue	chemagic RNA Kit H24**	CMG-1084
RNA	Blood from PAXGene Tubes	Up to 2.5 ml blood	chemagic RNA PGBlood Kit H24	CMG-1100
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	1 - 2 ml	chemagic cfDNA 2k Kit H24	CMG-1302
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	3 - 5 ml	chemagic cfDNA 5k Kit H24	CMG-1304
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	6 - 10 ml	chemagic cfDNA 10k Kit H24	CMG-1310
miRNA (cfRNA)	Plasma, serum, exosomes, saliva	1 ml	chemagic miRNA 1K Kit H24	CMG-1223
Pathogen NA	Body fluids, blood, ticks, cultures, supernatant of stool samples	Up to 1 ml	chemagic Body Fluid 1K Kit H24***	CMG-1103

\* Kits require swabs for sample collection.

\*\* This is a universal kit for RNA extraction. An optimized kit solution for PAXGene blood tubes is also available (CMG-1100).

\*\*\* Sample collection can be done directly or with swabs.



## chemagic 360 Kits for chemagic 12 Rod Head

Preps/Kit: 250 | Reaction vessel: 50 ml Falcon Tubes | Some plastics (tip sleeves) included. | 50 ml Falcon tubes not provided with kit.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Blood, buffy coat, saliva	Up to 10 ml	chemagic DNA Blood 10K Kit H12	CMG-704
gDNA	Blood, buffy coat, saliva	Up to 7 ml	chemagic DNA Blood 7K Kit H12	CMG-715
gDNA	Buffy coat	Up to 2 ml	chemagic DNA Buffy Coat 2K Kit H12	CMG-728
gDNA	Saliva	Up to 4 ml	chemagic DNA Saliva 4K Kit H12	CMG-1035
RNA	Tissue, cells from cell cultures, white blood cells	Up to 2.5 ml blood or 10 mg tissue	chemagic Total RNA Kit H12	CMG-1083
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	10, 15 or 18 ml	chemagic cfDNA 18k Kit H12	CMG-1318
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 9.6 ml	chemagic Viral DNA/RNA 9.6K Kit H12*, **	CMG-749

\* Sample collection for Kits can be done directly or with swabs.

\*\* The kits "chemagic Viral DNA / RNA 300 Kit H96" and "chemagic Viral DNA/RNA 9.6K Kit H12" can be also used for pathogen detection from wastewater (COVID).



## chemagic Prime Kits for chemagic 96 Rod Head

Preps/Kit: 960 | Reaction vessel: 96 deep-well plate | All plastic consumables included.

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Blood, buffy coat, saliva	Up to 400 µl	chemagic Prime DNA Blood 400 Kit H96	CMG-1491
gDNA	Buffy coat	Up to 200 µl	chemagic Prime DNA Buffy Coat 200 Kit H96	CMG-1413
gDNA	Saliva	Up to 600 µl	chemagic Prime DNA Saliva 600 Kit H96	CMG-1437
gDNA	FFPE tissue	10 µm section	chemagic Prime FFPE DNA Kit H96	CMG-1499
RNA	Tissue, cells from cell cultures, blood	Up to 2.5 ml blood or 10 mg tissue	chemagic Prime RNA Kit H96	CMG-1412
Pathogen NA	Blood, body fluids	Up to 200 µl	chemagic Prime Viral NA/gDNA 200 Kit H96	CMG-1449
Pathogen NA	Blood, body fluids	Up to 1 ml	chemagic Prime Viral NA/gDNA 1K Kit H96	CMG-1431
Pathogen NA	Stool	Up to 1 g	chemagic Prime DNA Stool Kit H96	CMG-1476

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact



## chemagic Prime Kits for chemagic 24 Rod Head

Preps/Kit: 240 | Reaction vessel: 24 deep-well plate | All plastic consumables included.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Blood, buffy coat, saliva	Up to 2 ml	chemagic Prime DNA Blood 2K Kit H24	CMG-1497
gDNA	Blood	Up to 4 ml	chemagic Prime DNA Blood 4K Kit H24	CMG-1474
gDNA	Saliva	Up to 4 ml	chemagic Prime DNA Saliva 4K Kit H24	CMG-1481
RNA	Tissue, cells from cell cultures, white blood cells	Up to 2.5 ml blood or 10 mg tissue	chemagic Prime Total RNA 2.5K Kit H24	CMG-1484
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	Up to 5 ml	chemagic cfDNA 5k Kit H24	CMG-1444
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 1 ml	chemagic Prime Body Fluid 1K Kit H24*	CMG-1406

\* Sample collection can be done directly or with swabs.



## chemagic Prime Jr Kits

Preps/Kit: 480 | Reaction vessel: 96 deep-well plate | All plastic consumables included.

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Diverse sample types, blood, buffy coat, saliva, plasma, serum, lavage, tissue, cells, DBS, bacterial culture	200 - 800 $\mu$ l	Junior gDNA Kit*	CMG-1534
Pathogen NA	Diverse sample types, blood, buffy coat, saliva, plasma, serum, lavage, tissue, cells, DBS, bacterial culture	200 - 1,000 $\mu$ l	Junior Pathogen NA Kit*	CMG-1521
RNA	Tissue, cells from cell cultures, white blood cells	Up to 2.5 ml blood or 10 mg tissue	Junior Total RNA Kit	CMG-1512

\* Sample collection can be done directly or with swabs.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact





## chemagic Prepito Kits

Preps/Kit: 180 | Reaction vessel: Micronic Tube / 96 deep-well plate | All plastic consumables included.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Blood, buffy coat, saliva	250 µl	Prepito DNA Blood250 Kit	CMG-2002
gDNA	Blood, buffy coat, saliva	400 µl	Prepito DNA Blood400 Kit	CMG-2003
gDNA	Blood, buffy coat, saliva	600 µl	Prepito DNA Blood600 Kit	CMG-2004
gDNA	Diverse sample types, blood, buffy coat, saliva, plasma, serum, lavage, tissue, cells, DBS, bacterial culture	250 µl, 10 mg tissue, 3 - 5 ml amniotic fluid	Prepito DNA Universal Kit	CMG-2034
gDNA	FFPE tissue	10 µm sections	Prepito FFPE Kit	CMG-2027
gDNA	Tissue, cells	10 mg tissue	Prepito DNA Tissue10 Kit	CMG-2010
RNA	Tissue, cells from cell cultures, white blood cells	250 µl, 10 mg tissue	Prepito RNA Kit**	CMG-2039
RNA	Blood from PAXGene Tubes	Up to 2.5 ml blood	Prepito RNA PGBlood Kit	CMG-2041
cfDNA (ctDNA)	plasma, serum, exosomes, urine	Up to 1.5 ml	Prepito cfDNA 1.5K Kit	CMG-2025
miRNA (cfRNA)	plasma, serum, exosomes, saliva	Up to 200 µl	Prepito miRNA 200 Kit	CMG-2042
Pathogen NA	Blood, body fluids	Up to 200 µl	Prepito NA Body Fluid Kit	CMG-2021
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 300 µl	Prepito Viral DNA/RNA300 Kit*	CMG-2017
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 1 ml	Prepito Viral DNA/RNA1k Kit*	CMG-2018

For research use only.  
Not for use in diagnostic procedures.

\* Sample collection can be done directly or with swabs. | \*\* This is a universal kit for RNA extraction. An optimized kit solution for PAXGene blood tubes is also available (CMG-2041).



## chemagic Kits for third party instruments

Preps/Kit: 960 (unless otherwise stated\*) | Reagents only.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Target	Sample type	Sample volume	Product name	Product no.
PCR products	Reaction mixes	20 µl	chemagic PCR Pure Kit LH	CMG-450
Sequencing products	Reaction mixes	20 µl	chemagic SEQ Pure Kit LH	CMG-458
gDNA	Dried blood spot (DBS)	3 - 13 mm punch out	chemagic DNA Blood Spot Kit LH	CMG-1202
gDNA	Blood	Up to 200 µl	chemagic DNA blood 200 kit LH	CMG-460
gDNA	Tissue	10 mg tissue	chemagic NA Tissue Kit LH	CMG-476-HP
gDNA	Blood, buffy coat, saliva	400 µl	chemagic NA Universal Kit*	CMG-437
Pathogen NA	Body fluids, saliva, sputum, serum, plasma, lavage, semen, amniotic fluid	Up to 300 µl	chemagic Viral DNA/RNA Kit	CMG-1033-P

\* The "chemagic NA Universal Kit" is delivered with 480 preps/Kit.



## chemagic Kits for manual workflows

Preps/Kit: 40-50 | Reaction vessel: Tubes | Reagents only.

Target	Sample type	Sample volume	Product name	Product no.
gDNA	Diverse sample types, blood, buffy coat, saliva, plasma, serum, lavage, tissue, cells, DBS, bacterial culture	100 µl / 250 µl blood, 10 mg /40 mg tissue material, cell pellets from 1 - 5 ml amniotic fluid, DBS punches and bacteria pellets	chemagic DNA Universal Kit (50 preps/kit)*	CMG-196
cfDNA (ctDNA)	Plasma, serum, exosomes, urine	1 - 5 ml	chemagic cfDNA 5k Kit (40 preps/kit)	CMG-134

\* Sample collection can be done directly or with swabs.

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact



## Magnetic Stand Accessories

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

Compatible tubes/plates	Number of positions	Protocols applied with	Product name	Product no.
1.5 ml or 2 ml tubes	12 magnetic and 12 non-magnetic positions	Manual workflows for up to 2 ml preparation volumes	chemagic Magnetic Stand 2x12	CMG-300
96 Deep-well plate	96 magnetic positions	Automated 96-well format workflows	chemagic Magnetic Stand 96	CMG-301
50 ml tubes	2 magnetic positions	Manual workflows for up to 50 ml preparation volumes	chemagic Magnetic Stand 50k Type A	CMG-305
50 ml tubes	2 magnetic positions	Manual workflows for up to 15 ml preparation volumes	chemagic Magnetic Stand 50k Type B	CMG-303
50 ml tubes	1 magnetic position (Position secures tube for discard of supernatant by pouring)	Manual workflows for up to 50 ml preparation volumes	chemagic Magnetic Stand 50k Type E	CMG-304
50 ml tubes	1 magnetic position (Position secures tube for discard of supernatant by pouring)	Manual workflows for up to 15 ml preparation volumes	chemagic Magnetic Stand 50k Type F	CMG-302
50 ml tubes	12 magnetic positions	Manual workflows for up to 15 ml preparation volumes	chemagic Magnetic cfDNA Stand	CMG-306
PCR plates	96 magnetic positions	For automated PCR clean-up	chemagic Stand 96 PCR	CMG-307



# References

- chemagic technology
- Advantages of our technology
- Solutions for your needs
- Genetic analysis / biobanking
- Precision oncology
- Translational research
- Infectious disease research
- M-PVA Magnetic Beads
- chemagic instruments
- chemagic kits
- Product list
- Contact

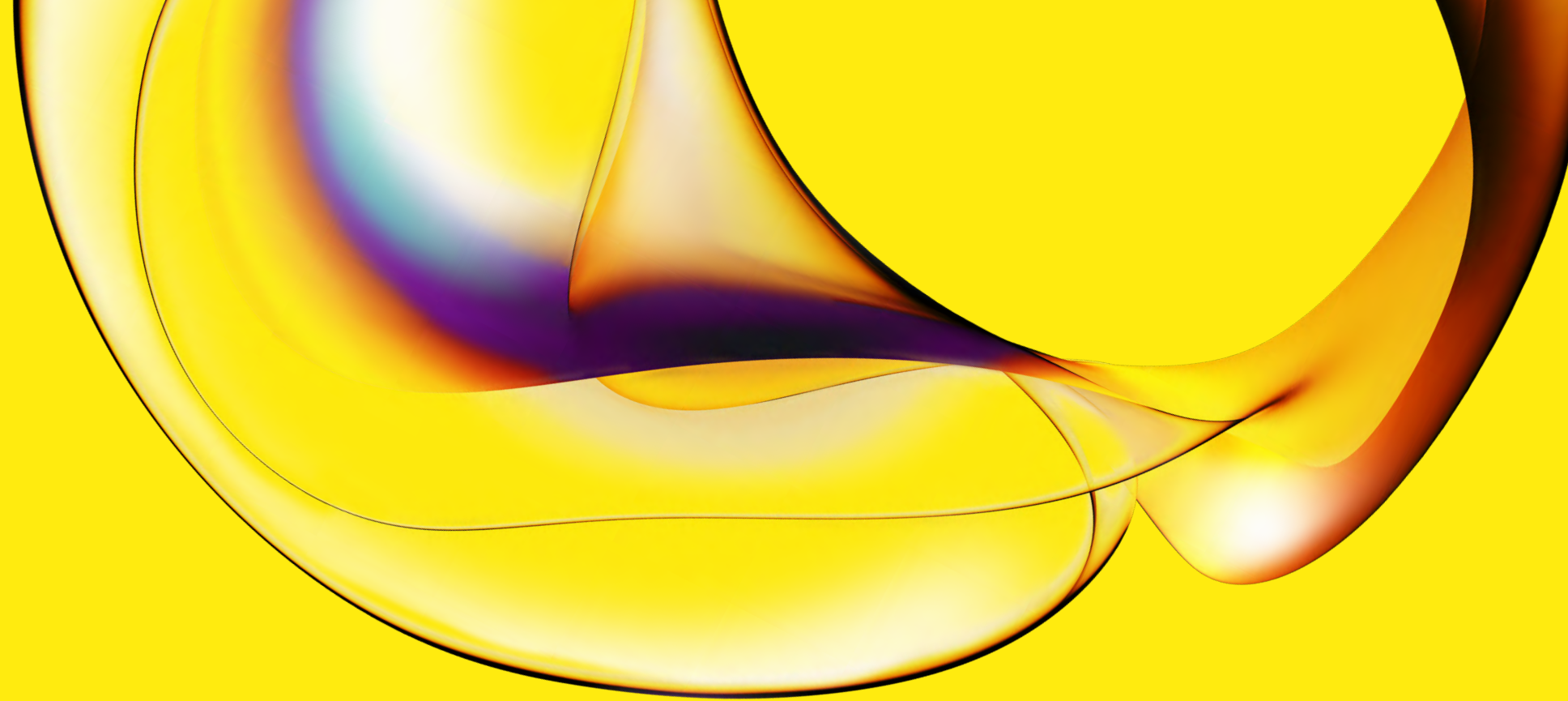
1. Oster J, Parker J, Brassard LÅ. Polyvinyl-alcohol-based magnetic beads for rapid and efficient separation of specific or unspecific nucleic acid sequences. *J Magn Magn Mater.* 2001;225:145-150. doi: 10.1016/S0304-8853(00)01243
2. Schöfl G, Lang K, Quenzel P, et al. (2017). 2.7 million samples genotyped for HLA by next generation sequencing: lessons learned. *BMC Genomics* 18:1, 18(1), 1-16. <https://doi.org/10.1186/S12864-017-3575-Z3>.
3. Macha S, Koenen R, Sennewald R, et al. (2014). Effect of Gemfibrozil, Rifampicin, or Probenecid on the Pharmacokinetics of the SGLT2 Inhibitor Empagliflozin in Healthy Volunteers. *Clinical Therapeutics*, 36(2), 280-290.e1. <https://doi.org/10.1016/j.clinthera.2014.01.003>
4. Evans DG, Lalloo F, Ryan NA, Bowers N, Green K, Woodward ER et al. (2022). Advances in genetic technologies result in improved diagnosis of mismatch repair deficiency in colorectal and endometrial cancers. *Journal of Medical Genetics*, 59(4), 328-334. <https://doi.org/10.1136/JMEDGENET-2020-107542>
5. Balciuniene, Jorune et al. (2023). At-Risk Genomic Findings for Pediatric-Onset Disorders From Genome Sequencing vs Medically Actionable Gene Panel in Proactive Screening of Newborns and Children. *JAMA Network Open*, 6(7), e2326445-e2326445. <https://doi.org/10.1001/JAMANETWORKOPEN.2023.26445>
6. Beyter D, Ingimundardottir H, Oddsson A, et al. (2021). Long-read sequencing of 3,622 Icelanders provides insight into the role of structural variants in human diseases and other traits. *Nature Genetics*, 53(6), 779-786. <https://doi.org/10.1038/s41588-021-00865-4>
7. Callesen LB, Hansen TF, Andersen RF, et al. (2023). ctDNA-guided adjuvant treatment after radical-intent treatment of metastatic spread from colorectal cancer-the first interim results from the OPTIMISE study. *Acta Oncologica (Stockholm, Sweden)*, 62(12), 1742-1748.
8. Frank MS, Andersen CSA, Ahlborn LB, Pallisgaard N, Bodtger U, Gehl J. (2022). Circulating Tumor DNA Monitoring Reveals Molecular Progression before Radiologic Progression in a Real-life Cohort of Patients with Advanced Non-small Cell Lung Cancer. *Cancer Research Communications*, 2(10), 1174-1187. <https://doi.org/10.1158/2767-9764.CRC-22-0258>
9. Lefèvre AC, Pallisgaard N, Kronborg C, Wind KL, Krag SRP, Spindler KG. (2021). The clinical value of measuring circulating hpv dna during chemo-radiotherapy in squamous cell carcinoma of the anus. *Cancers*, 13(10). <https://doi.org/10.3390/cancers13102451>
10. Benhamida JK, Hechtman JF, Nafa K, et al. (2020). Reliable Clinical MLH1 Promoter Hypermethylation Assessment Using a High-Throughput Genome-Wide Methylation Array Platform. *The Journal of Molecular Diagnostics*, 22(3), 368-375. <https://doi.org/10.1016/J.JMOLDX.2019.11.005>
11. Sorrells S, McKinnon KE, McBratney A, Sumey C. (2021). Longitudinal and multi-tissue molecular diagnostics track somatic BRCA2 reversion mutations that correct the open reading frame of germline alteration upon clinical relapse. *Npj Genomic Medicine* 2021 6:1, 6(1), 1-6. <https://doi.org/10.1038/s41525-021-00181-0>
12. Schieck, Maximilian et al. (2020) Implementation of RNA sequencing and array CGH in the diagnostic workflow of the AIEOP-BFM ALL 2017 trial on acute lymphoblastic leukemia. *Annals of hematology* vol. 99(4) 809-818. <https://doi.org/10.1007/s00277-020-03953-3>
13. Costa E, Rocha D, Lopes JIF, et al. (2024). Detection of Plasmodium spp. in asymptomatic blood donors by the new Brazilian NAT PLUS HIV/HBV/HCV/Malaria Bio-Manguinhos kit. *Transfusion*, 64(3), 501-509. <https://doi.org/10.1111/TRF.17726>



- **chemagic technology**
- **Advantages of our technology**
- **Solutions for your needs**
- **Genetic analysis / biobanking**
- **Precision oncology**
- **Translational research**
- **Infectious disease research**
- **M-PVA Magnetic Beads**
- **chemagic instruments**
- **chemagic kits**
- **Product list**
- **Contact**

## References

14. LaTurner, Zachary W et al. (2021) Evaluating recovery, cost, and throughput of different concentration methods for SARS-CoV-2 wastewater-based epidemiology" Water research vol. 197: 117043. doi:10.1016/j.watres.2021.117043 (2021). Evaluating recovery, cost, and throughput of different concentration methods for SARS-CoV-2 wastewater-based epidemiology. Water Research, 197. (2021).
15. West NW, Vasquez AA, Bahmani A, et al. (2022) Sensitive detection of SARS-CoV-2 molecular markers in urban community sewersheds using automated viral RNA purification and digital droplet PCR. Sci Total Environ. 847:157547. doi:10.1016/j.scitotenv.2022.157547



[www.revivity.com](http://www.revivity.com)

revvity

Copyright ©2024, Revvity, Inc. All rights reserved.

1426889