

Streamlined
walk-away
automation.



revvity

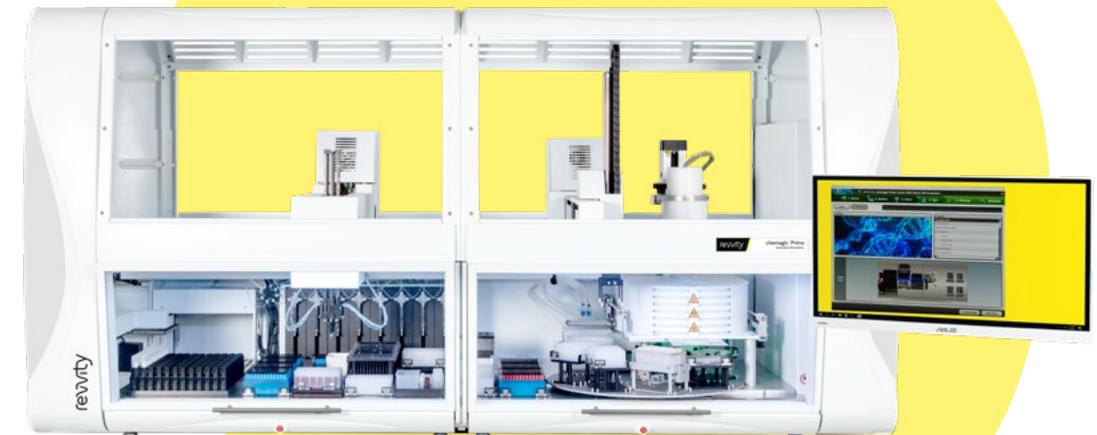
chemagic™ Prime™ instrument

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



It's prime time.

chemagic Prime:
For all human
sample materials.



The chemagic Prime
Dimensions: L 2.3 m, W 0.9 m, H 1.9 m (incl. table)

Integrated automation of nucleic acid isolation to assay setup

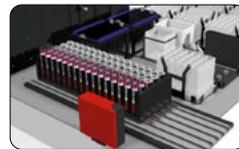
The chemagic Prime instrument is a fully automated research solution for primary sample transfer, DNA and RNA isolation, optional normalization, and PCR setup applications, that offers streamlined, walk-away sample processing. It delivers contaminant-free, extracted DNA and RNA in high concentrations, offering researchers a validated, single supplier workflow.

Nucleic acid isolation is a critical prerequisite for many downstream applications. The chemagic Prime instrument uses patented magnetic bead technology (M-PVA-Magnetic Beads) with fully automated liquid handling to provide high quality and high yield isolation of nucleic

acids from a variety of sample types, suitable for NGS, MLPA, genotyping and PCR. Unlike other automated solutions, the chemagic Prime instrument uses magnetized rods instead of magnetic plates to separate nucleic acids from solutions. By transferring the beads instead of the process solutions, the contamination risk is minimized, and higher purity, more intact DNA and RNA can be isolated. Reagent kits are available for the isolation of DNA and RNA from various human samples including whole blood, saliva, plasma, tissues, FFPE samples, and feces.



Make your lab a prime site



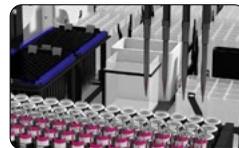
Full traceability

- Barcode reading of sample tubes
- Barcode reading of output tubes/ plates racks
- Reagent lotnumber tracking
- Bidirectional LIMS communication



Integrated chemagic dispenser

- Flexible reagent volumes
- Economic bulk reagents
- Fast processing times



Integrated liquid handling

- Handling of primary samples
- Dispensing of magnetic beads, elution buffer and required enzymes
- Final eluate transfer with integrated removal of magnetic bead traces
- Optional addition of internal controls
- Tracking of all liquid operations
- Safe travel path



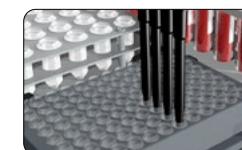
Superior chemagic kit and technology performance

- Transfer of magnetic beads instead of liquids
- High yields, recoveries and purities
- Outstanding performance for human sample materials
- Delivering high molecular weight DNA



Double sample set capacity

- Walkaway automation for 2 x 24 and 2 x 96 samples
- Minimizes hands on time
- Gripper hardware and cradles/racks optimized for transports



Optional add ons

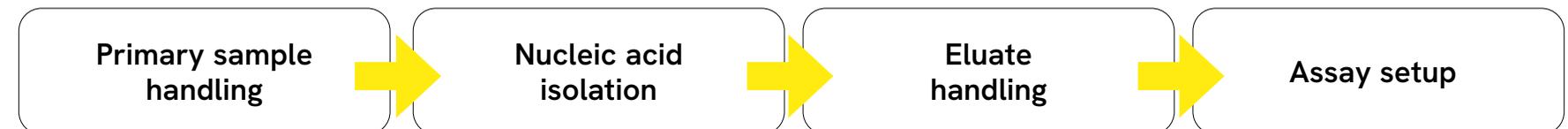
- PCR setup
- Eluate aliquoting
- Normalization

Supported applications

- Genomic DNA and RNA isolation from blood, cells, saliva and tissue for biobanking
- Isolation of viral/bacterial nucleic acids from serum/plasma pools
- Processing human body fluids for research applications
- Extraction of high molecular weight DNA for long read sequencing
- Isolation of DNA from various samples for microbiome analysis



Workflow



Meet another chemagic Prime family member

The chemagic Prime Junior instrument - designed with flexibility in mind. Streamlined workflow for up to 48 (2 x 24) low volume samples (up to 1ml).





Kits information

Selection of chemagic kits to be applied on the chemagic Prime instrument using the chemagic 96 Rod Head

Target	Sample type	Sample volume	Product name	Product number
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	Blood, buffy coat, saliva	up to 2 ml	chemagic Prime DNA Blood 2K Kit H24	CMG-1497
gDNA	Saliva	up to 600 µl	chemagic Prime DNA Saliva 600 Kit H96	CMG-1437
gDNA	Saliva	up to 4 ml	chemagic Prime DNA Saliva 4K Kit H24	CMG-1481
gDNA	FFPE tissue	10 µm section	chemagic Prime FFPE DNA Kit H96	CMG-1499
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
RNA	Tissue, cells from cell cultures, white blood cells	up to 2.5 ml blood or 10 mg tissue	chemagic Prime RNA Kit H96	CMG-1412

Kits information

Selection of chemagic Kits to be applied on the chemagic Prime instrument using the chemagic 24 Rod Head

Target	Sample type	Sample volume	Product name	Product number
Pathogen NA	Stool	up to 1 g	chemagic Prime DNA Stool Kit H96	CMG-1476
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
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



Ordering information

Product	Product number
chemagic Prime 8 instrument (8 pipetting channels) Streamlined walkaway instrument for primary sample transfer, DNA/RNA isolation and assay setup	CMG-404
Mandatory items	
chemagic Prime IQ/OQ	CMG-13005663
chemagic Prime ID Sync QA	CMG-13005665
Required accessories (at least one to be ordered with the instrument)	
chemagic Prime 24 Rod Head Set	CMG-387
chemagic Prime 96 Rod Head Set	CMG-388
Optional accessories	
chemagic Prime UV-light option	CMG-CLS145057
2D Flatbed Tube Reader Upgrde Kit (Ziath DataPaq ZTS-EXT)	CMG-13009695

Please refer to the chemagic Prime order guide for more optional accessories.



www.revivity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2024, Revvity, Inc. All rights reserved.

1320685