



# Fast, compact, fully automated: chemagic Prepito.

The chemagic™ Prepito™ nucleic acid extraction unit is based on chemagic proven technology for magnetic particle separation and represents the top quality sample preparation system in a compact benchtop instrument.

It utilizes many years of experience in automated nucleic acid isolation gained with the high throughput chemagic 360 instrument. In combination with the Prepito Kits, it delivers high integrity, yield, and purity of DNA/RNA, and ensures the success of your downstream application.

This innovative instrument realizes cost effectiveness through automated dispensing of buffers into standard plastic devices instead of using expensive prefilled cartridges. The chemagic Prepito instrument includes bar code reading and a USB port data transfer to support a high quality assurance. Hands-on time is minimized. Processing time is as short as 40 minutes for 1 - 12 samples per batch (for e.g. 250 µl blood).



| chemagic Prepito nucleic acid extraction unit

## Features at a glance

- 1 - 12 samples per run
- Small benchtop solution
- Integrated buffer dispensing
- Revolutionary resuspension technology
- Standard plastic devices (included in the Prepito Kits)
- Prepito Kits and pre-installed protocols for a huge variety of sample materials
- High integrity, yield, and purity of RNA and DNA (move to top of this list)
- Input volumes up to 1 mL
- Barcode reading
- LIMS-compatible log files
- USB port

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

Product name	Product number	No. of samples	Sample volume	Barcode reading	Dispensing
chemagic Prepito instrument*	2022-0030	1 - 12	100 µL - 1 mL	included	integrated

## Prepito Kits - Ready-to-use for diverse needs

Product name	Sample type	Sample size	Product no.
Prepito DNA Blood250 Kit	Whole Blood	250 µL	CMG-2002
Prepito DNA Blood400 Kit Blood	Blood, buffy coat, saliva	400 µl	CMG-2003
Prepito DNA Blood600 Kit	Whole Blood	600 µL	CMG-2004*
Prepito DNA Tissue10 Kit	Tissue	10 mg liver, 2 - 4 mm mouse tail	CMG-2010
Prepito viral DNA/RNA200 Kit	Serum/plasma	200 µL	CMG-2015
Prepito viral DNA/RNA300 Kit	Serum/plasma	300 µL	CMG-2017
Prepito viral DNA/RNA 1 k Kit	Serum/plasma	1000 µL	CMG-2018
Prepito miRNA 200 Kit	Serum/plasma	200 µl	CMG-1242
Prepito NA Body Fluid Kit	Blood, serum/plasma, swabs, feces supernatant urine, bacterial o/n culture	200 µL	CMG-2021
Prepito cfDNA 1.5 K Kit	Plasma/serum	1500 µL	CMG-2025
Prepito FFPE Kit	FFPE tissue	up to 5 mg FFPE tissue material	CMG-2027
Prepito DNA Cyto Pure Kit	Various sample materials	e.g. 250 µL blood, 10 mg tissue, 3 - 5 ml pelleted amniotic fluid	CMG-2034
Prepito RNA Kit	Tissue, blood from EDTA tubes	Up to 2.5 ml blood, up to 10 mg tissues	CMG-2039
Prepito RNA PGBlood Kit	Blood from PAXGene tubes	Up to 2.5 ml	CMG-2041

### Prepito Kits are for Research Use Only. Not for use in Diagnostic Procedures.

All Prepito Kits\*\* contain ready-to-use reagents and all required plastic consumables. Standard kit size: 180 preps./kit for 96 well plates, for the processing of 1-12 samples in parallel.

\*90 preps./kit for 48 well plates, for the processing of 1-6 samples in parallel.

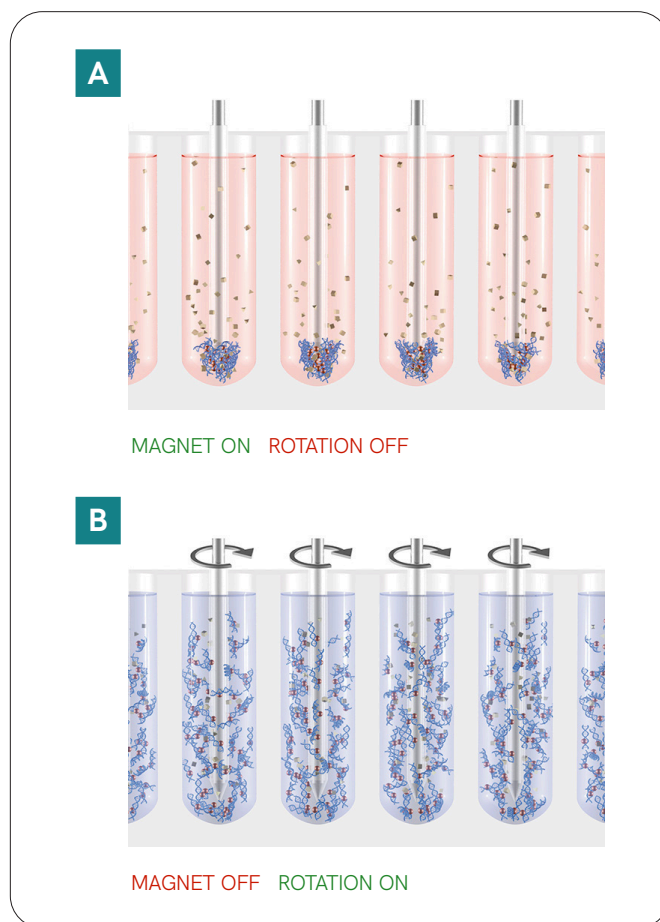
## Magnetic separation

The magnetic separation is based on the use of metal rods that are lowered into a process solution (A). To collect beads from the solution, the rods are magnetized. Pellets form at the tips of the rods, and the rods are withdrawn from the solution with the pelleted beads attached.

Resuspension into the next process solution, for example, wash or elution buffer, is achieved by switching off the magnetism while rotating the rods (B). This gentle but efficient resuspension technology ensures ultrapure, highly intact and high yields of nucleic acids obtained with minimal manual intervention.

### Benefits of chemagic automation

- Pure, intact and high yields of DNA/RNA
- Scaleable sample volumes from 100  $\mu$ L - 1 mL
- No detectable cross-contamination and carry-over of inhibitors
- Different sample materials in one run possible
- Reliable and reproducible results, even old, compromised, fresh or frozen, EDTA or citrate stabilized blood samples can be processed without any restriction
- Outstanding robustness of the instrument
- Cost-effective sample preparation
- Enhanced success rates for your downstream applications



| Figure 1: chemagic Separation Technology

revvity