



Nucleic acid extraction automation

Leveraging the chemagic™ 360 instrument and explorer™ G3 workstation to scale throughput and minimize manual intervention.

Why automate?

- **Increase throughput:** integrate multiple chemagic™ 360 instruments
- **Increase hands-off time:** scalable walk-away capacity
- **Reduce errors:** by decreasing human touchpoints
- **Increase reproducibility:** with equal sample treatment through automating critical steps
- **Add additional functions:** concentration measurements, normalization, NGS library preparation, (q)PCR, transfection

Why automate with Revvity?

- **Application expertise:** with our broad product portfolio we support your application from sample extraction to endpoint analysis
- **One point-of-contact:** we offer complete solutions including 3rd party devices and serve as your first contact for all requests
- **Dedicated team:** built over 25 years to support labs across the world

Where to start?

Identify your goals:

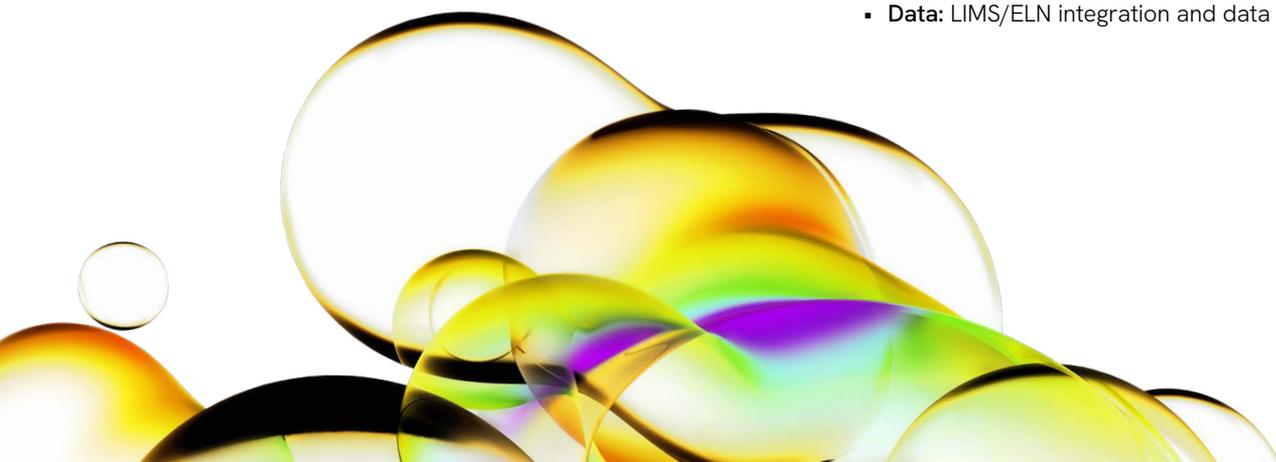
- increase throughput
- decrease hands-on time
- add additional functionalities

Identify workflows that will benefit most from automation:

- Which workflow creates the biggest bottleneck?
- Which workflow puts the largest strain on staff?
- Where do you have the highest risk for human error?
- Where do you need to reduce contamination risks?

Identify your requirements:

- **Functionality:** pre- and post-extraction sample processing requirements
- **Samples:** type, volume, number, container
- **Throughput:** requirements for throughput and hands-off time
- **Containment and contamination:** sample and/or operator protection
- **Data:** LIMS/ELN integration and data analysis



Application examples

Extraction from DBS, NGS library prep and biobanking

- **DNA extraction:** extract DNA from dried blood spots
- **Normalization and quality:** measure DNA concentration and fragment size, normalize samples
- **NGS:** automate NGS library preparation on liquid handler
- **Biobanking:** store DNA in sealed plates or tubes



Liquid biopsy

- **Tube handling:** sort, decap and scan sample tubes
- **Sample preparation:** tube-to-plate transfer and reagent addition
- **cfDNA extraction:** extract from 0.5 - 8 ml of plasma, urine or cerebral spinal fluid (CFS)
- **Sample analysis:** automate concentration measurement, fragment analysis, dPCR/qPCR, and NGS library preparation



Pathogen detection

- **Sample preparation:** tube-to-plate transfer and reagent addition
- **Nucleic acid extraction:** extract from a variety of microorganisms and sample materials
- **Sample analysis:** automate concentration measurement, fragment analysis, dPCR/qPCR, and NGS library preparation



Demos available!

We also support remote demos.
Contact us to schedule an appointment!



Nucleic acid
purification

Scan or click
for more information



Contact us

Scan or click
for more information



Integrated lab
automation

Scan or click
for more information

