



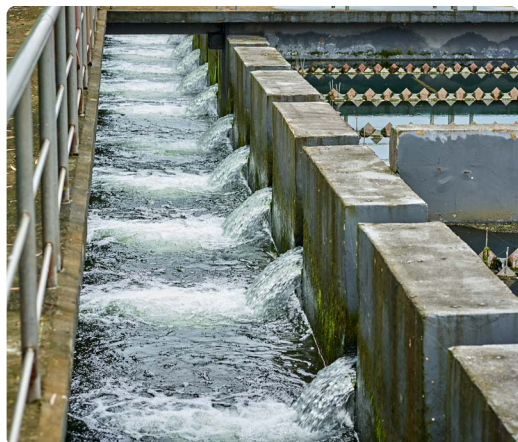
Scale up, effortlessly.

Automated nucleic acid purification for diverse applications

Gain an edge with established chemagic™ automated nucleic acid purification, trusted by small to large laboratories worldwide in challenging downstream applications.

Large sample volume processing

- Detect low abundance targets!
- From rare ctDNA mutations to low viral / bacterial copies in wastewater.

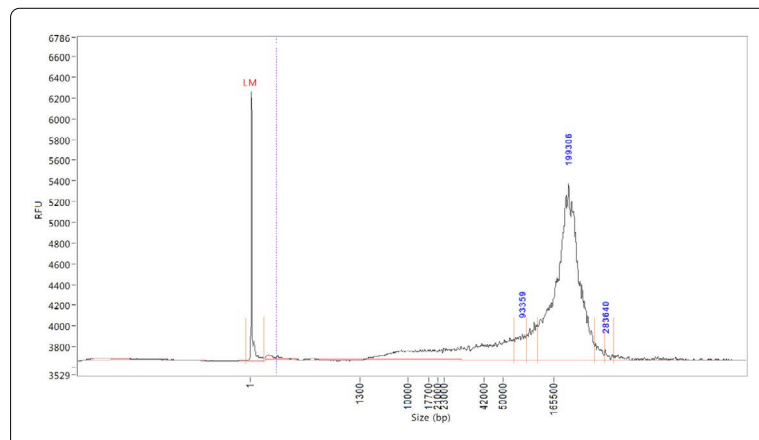


"More than 10 years ago, we chose chemagen instruments and reagents for isolating DNA from blood or buccal swabs. After performing HLA genotyping on over 10 million samples, we remain extremely satisfied with this decision and the consistent performance."

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

Efficient high molecular weight DNA extraction

- Obtain stable HMW DNA > 100 kb in ~1 hour!
- Proven in large-scale population studies involving long read sequencing.





Reliable automation with excellent support

- Increase instrument uptime and maximize efficiency!
- Single supplier solution for faster response times and customized reagent solutions.

Fully automated workflows

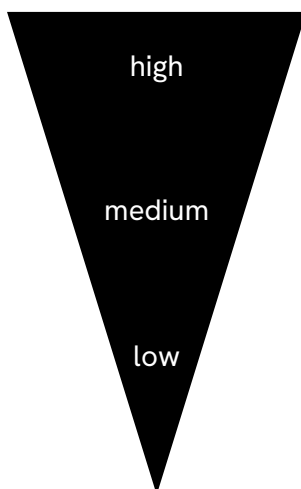


chemagic Prime

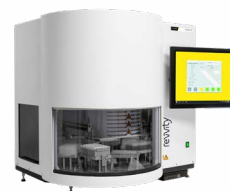


chemagic Prime Jr

Sample throughput



For combined workflows



chemagic 360



chemagic Prepito



Achieve similar extraction performance across chemagic platforms - less time optimizing, more time delivering results!



Begin your
free trial*

* Terms and conditions apply

revvity