



revvity

COMING THIS SUMMER

The new generation of automated nucleic acid purification.

chemagic™ Neo Instrument

Scale without limits.

From small studies to high-throughput, one instrument independent from batch size, or sample type.

Results you can count on.

Consistent, reliable performance, run per run.

High yield. Less compromise.

M-PVA magnetic bead technology delivers optimal purity for your downstream applications.

More samples. Less time.

Accelerate your workflow without sacrificing quality.

Complexity handled.

Smart automation takes care of the hard part, so you can focus on your research.



www.revvity.com

Copyright ©2026, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc.

All other trademarks are the property of their respective owners.

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

2734324