

Customize your high-throughput nucleic acid purification workflows with our expertise

Breadth of NAP Expertise

Promega has a breadth of nucleic acid purification (NAP) expertise that can address virtually all application and extraction needs. We offer chemistry and instrument capabilities at all scales and a range of nucleic acid extraction products that generate high yields from many sample types.

Breadth of Portfolio

Our unique chemistries are optimized for purification from a variety of samples, including:

- Saliva
- Plasma
- Fresh cells
- Bacteria
- Plant
- Serum
- Whole blood
- Frozen cells
- Formalin-fixed, paraffin-embedded (FFPE) samples

Build Your High-Throughput Solution with Our Experts

1



Automation Consultation

We consult with you to deeply understand your unique needs.

2



Proposal

Our Field Support Scientist team will draft a Statement of Work (SOW) to ensure we are aligned with your needs.

3



On-site Implementation

We schedule our onsite work so that it best fits your schedule.

Adapt Product Format for Efficient, High-Throughput Workflow

4



Custom packaging and kitting as required.

5



Safe and on-time delivery of your custom solution.

6



Ongoing Technical Support.

Breadth of Scale

From single-prep manual solutions, to cartridge-based Maxwell® Systems, to plate-based processing compatible with high-throughput liquid handling systems, Promega chemistries cover the range of your purification needs.

Flexible Formats

One size does not fit all. Rarely are high-throughput NAP workflows the same between laboratories. We can adapt NAP product formats to optimize your laboratory's efficiency.

Custom High-Throughput Nucleic Acid Purification

Primary Manufacturer

We have a team dedicated to DNA/RNA purification resin and buffer manufacturing, with over 45 years of combined NAP experience. We have dedicated laboratory manufacturing space and equipment capable of handling buffer batch sizes up to several thousand liters and multi-kilogram resin batches. We also have a team dedicated to physical and functional characterization and advanced QC methodology of our resins and buffers. Since we are the primary manufacturer of purification reagents and resins, you will collaborate directly with the scientists who have designed, manufactured and tested the products that are delivered to you.



High-Throughput Purification with Experts Included

Our team of automation experts will help you choose an optimized high-throughput solution that meets your laboratory's needs. We can implement Maxwell® HT chemistries on a variety of sample types and instrumentation platforms. Maxwell® HT Systems allow purification of DNA or RNA at scale on any laboratory liquid handler in 24-, 96- and 384-well SLAS format. Maxwell® purification chemistries use novel magnetic particle-based solutions that naturally decrease carryover contamination.

We can purify high-quality DNA and RNA at any scale. Our customized chemistries and easy-to-use protocols are designed for your specific sample type, ensuring consistent, reliable results in downstream assays.

Expert Support

In addition to trusted chemistry, you will gain expert support to get started with automation or optimize your current high-throughput workflow. Our team of automation experts can offer assistance with most of the leading laboratory automation providers in the world and help you develop and implement an automated NAP solution customized to the needs of your laboratory.

Promega has created specialized teams of Field Support Scientists and Field Application Scientists with a combined scientific and programmatic expertise of more than 125 years. We believe this service sets us apart from other providers who may be unable to work with third-party platforms.

Variety of Platforms

In addition, we have a variety of high-throughput platforms in our Applications and Training centers. Applications Scientists partner with our field-based experts to develop and verify new methods, provide proof-of-principle test customer samples, and push the limits of methods and chemistries. Most of our applications scientists have advanced scientific degrees,

and collectively we have over 90 years of experience adapting, developing and supporting Promega technologies. We have three Applications and Training Centers globally (Madison, Wisconsin; Charbonnières-les-Bains, France; and Singapore) to meet regional needs. Each center has permanent applications scientists and all are equipped to do applications testing with Promega products.

From manufacturing, QC, shipment and support, Promega has been at the leading edge of purification technology and automation for over 35 years. Our unique expertise sets us apart as a perfect partner for your high-throughput chemistry needs.

Let's **TALK**
CUSTOM

Contact us today!

One of our scientists will be in touch to discuss your needs, address your challenges and serve as your champion in finding a custom solution.

www.promega.com/CustomSolutions

custom@promega.com

