

Product format and packaging

Finding the right custom manufacturing partner can be challenging. There are many options from which to choose and many decisions to be made in order to match your particular needs to the capabilities of any potential supplier. It is important to find a manufacturer who is equipped to meet both your future and current needs, who can consistently deliver a quality product on time to your specifications and who understands and applies quality and regulatory requirements in accordance with internationally recognized standards. When evaluating a manufacturing partner, it is important to think beyond the product itself. Bottles/tubes, labeling and packaging are critical components of a complete custom solution. Promega has multiple custom packaging options to meet your needs, from bulk to private label to packaging designed to match your workflow.

Support

As part of our Custom process, Promega offers a personal service to ensure that our experts are accessible to support customers during project planning and implementation. Promega scientists provide technical consultation on custom product design to help you during the planning and manufacturing process. Our Packaging Development Department, staffed by Packaging Engineers, can help accelerate your product development by proving out specifications, providing expertise to assist with design decisions and offering advice based on previous manufacturing experience. Our team is ready to consult on primary, secondary and shipping containers, including labels.

Filling

Promega's four bottle-filling lines and three tube-filling lines provide for flexibility in meeting your needs today and as your business grows. In addition, we have cartridge-dispensing lines and glass-filling lines for products to be lyophilized.

Our filling lines have the capability to dispense product ranging from 10 microliters to 3.5 liters. In our newest Optima line, product quantities of 40,000 bottles can be dispensed per shift. We also maintain manual dispensing capabilities for smaller volume runs or unique container types.



Primary Packaging

Promega provides a choice of bottles and tubes. Our packaging team will work with you to ensure you are selecting the best primary container that not only meets your needs but maintains product integrity. When selecting a container, the following considerations are critical: Product compatibility, distribution environment, light sensitivity, end customer needs, and fit with your existing packaging brand strategy.

We have numerous containers in a variety of sizes:

- HDPE round bottles (natural, white)
- Polypropylene round bottles (natural)
- HDPE round bottles (amber)
- PETG square bottles (amber)
- Polypropylene skirted tubes (natural, amber)
- Polypropylene conical tubes (natural, amber)
- Glass vials (amber, clear)



Lyophilization

Promega has multiple lyophilizers offering a variety of capacity options for our customers. We use glass-filling lines that integrate rubber stoppers and aluminum seals for a complete lyophilization solution.



Custom Packaging

Labeling

Promega offers a variety of labeling choices, dependent on your intended use of the product. Our Custom team will outline your choices and work with you in label selection for both primary and secondary packaging. Label availability:

- Promega standard labels
- Unbranded blank stock
- Customer-specific labels

We have capabilities for manual and automated label application. We also have the capabilities to do 100% vision inspection on some of our lines. Labels can be pre-printed, partially printed (variable data) or full in-house printed.



Secondary Packaging

Promega has the ability to design and produce paperboard cartons, corrugated shippers, foam and labels. We can also utilize our extensive supplier network to meet unique customer needs. We now have the in-house capabilities to perform ship testing, as well as ASTM 4169 ISTA and custom distribution protocols.



Quality

We dispense and package thousands of products. We often get asked how we guarantee that there is no cross-contamination. This is controlled by:

- Line clearance and cleaning procedures
- Filling in Environmentally Controlled Rooms
- Following our Quality System procedures that are governed by ISO 9001 and ISO 13485
- Validation Standard Operating Procedures (SOPs)

Promega Sustainable Packaging

Our new packaging is part of our commitment to sustainability and corporate responsibility efforts that benefit our planet, people, communities and products. Now Promega uses Sustainable Forestry Initiative (SFI) Chain of Custody level certified fiber in our printed package material. The SFI label indicates a commitment to tracking fiber, from forest content to end product, using a third-party internationally recognized sustainability standard. In addition, the ink, coating and glue are sourced from renewable products containing no controversial materials.



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

