

CUSTOM PRODUCTS AND SERVICES - EXACTLY WHAT YOU WANT

Kathryn D. Hebble and Kristina A. Pearson

Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, USA

Promega service and support surround you; from design to delivery. Promega offers custom service in lot reservation, global delivery and scheduled ordering. Most Promega standard catalog products can be customized to meet a customer's specific needs. Here we provide examples of customized formulations, dispensed sizes and kitting options. Envision the exact formulation that you want – a custom solution is mixed to your specification every time. Dispensing flexibility ensures that Promega can provide product in the format that you need. This includes your choice of product volume in a variety of vials and bottles. With a complete range of dispensing and packing options, you choose: format, size container and labeling. Look around your lab. What would help make your work more efficient and effective?