

# Freedom EVO<sup>®</sup> / REMP SSS Factory Package

## Safer sample storage and processing

Reference 10001BO100397



Fully automated sample management solution for closed-loop automation

The Freedom EVO / REMP SSS Factory offers researchers fully automated sample storage and retrieval based on the unique REMP Tube Technology<sup>™</sup>. This technology enables individual sample access and manipulation by simply pushing the tube through the source rack to the destination rack, increasing the efficiency over other tube rack systems. REMP Tube Technology also offers the benefits of minimizing freeze/thaw cycles, library cleaning and compression, and avoiding oxidation and water uptake.

REMP storage systems complement the flexibility of the Freedom EVO. The combination of these two systems allows automation of a wide range of sample preparation processes. Laboratories can automate the storage of dry and liquid samples, and complete the downstream sample processing to offer closed-loop automation. Sophisticated software tracks your samples at all stages and allows them to be chosen for retrieval from imported sample lists or queries. After samples are picked in the -20 °C environment they are delivered in delivery racks to the Freedom EVO. The Freedom EVO then handles thawing, de-capping and aliquoting of samples to assay ready plates. The delivery rack is then returned to the REMP Small-Size Store and the software automatically updates the volume in each tube, ensuring subsequent tube deliveries have sufficient volume for your next experiment.

Scientific instrumentation; not for use in human, clinical or diagnostic procedures. All systems or products are tailor-made one-offs that were exclusively developed on customer demand. All displayed pictures and images originate from development, test or installation phases and do not claim to represent the final installation at customer site.

Tecan and Freedom EVO are in major countries registered trademarks and Reatrix is a trademark of Tecan Group Ltd., Männedorf, Switzerland. REMP is a registered trademark and REMP Tube Technology, REMP Tube Punching Module and REMP Small-Size Store are trademarks of REMP AG, Oberdiessbach, Switzerland.

© 2009, Tecan Trading AG, Switzerland, all rights reserved.