

## REMP Small-Size Store™ (SSS)

### Safer sample storage

Reference 10001BO100399

---



REMP Small-Size Store™ (SSS) for laboratory sample management

The REMP Small-Size Store (SSS) offers researchers fully automated sample storage and retrieval based on the unique REMP Tube Technology™.

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.

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 for further processing. The delivery rack is then returned to the SSS and the software automatically updates the location 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 is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland. REMP is a registered trademark, and REMP Small-Size Store and REMP Tube Technology are trademarks of REMP AG, Oberdiessbach, Switzerland.  
© 2009, Tecan Trading AG, Switzerland, all rights reserved