

Freedom EVO[®] and REMP Small-Size Store[™] (SSS) integration

Tecan has seamlessly integrated the REMP Small-Size Store[™] (SSS) with the Freedom EVO[®] liquid handling workstation, creating a fully automated liquid handling and sample management platform. The REMP SSS is a fully automated storage and retrieval system that fits perfectly into a research laboratory.

It can handle samples in a variety of formats, including microplates and the REMP Tube Technology[™] Consumables. Integration of the SSS with the Freedom EVO and Freedom EVOware[®] enables walkaway operation for sample retrieval and processing steps. REMP Tube Technology enables desired samples to be cherry picked within the store's controlled environment. The samples are easily transferred to Tecan's Freedom EVO workstation using an integrated plate shuttle. Samples are decapped using the REMP ACD96[™] and processed as required on the Freedom EVO workstation. Upon completion, the samples are subsequently capped and returned to storage when they are no longer needed. The combination of the sample storage and processing platform minimizes the time that samples are exposed to the environment and ensures they are returned to the SSS as soon as processing is completed.



Fully integrated Freedom EVO and REMP Small-Size Store (SSS)

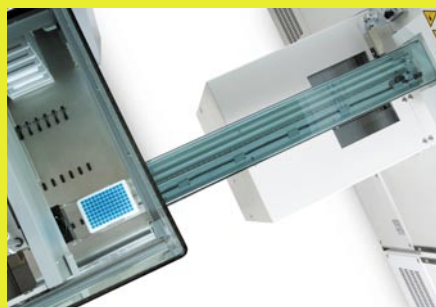
Integration, furthermore, enables increased process security. The REMP Reatrix 2D Scanner can be used on-board the Freedom EVO for tube and

plate identification. Additionally, sample volumes, concentrations and identification numbers can be fully tracked within the software. Data

handling is perfectly integrated, making this platform ideal for compound dissolution and reformatting factories, or genomic sample factories.



The plate carrier transports samples between the Small-Size Store (SSS) and Freedom EVO



The REMP Reatrix 2D Scanner can be used on-board the Freedom EVO