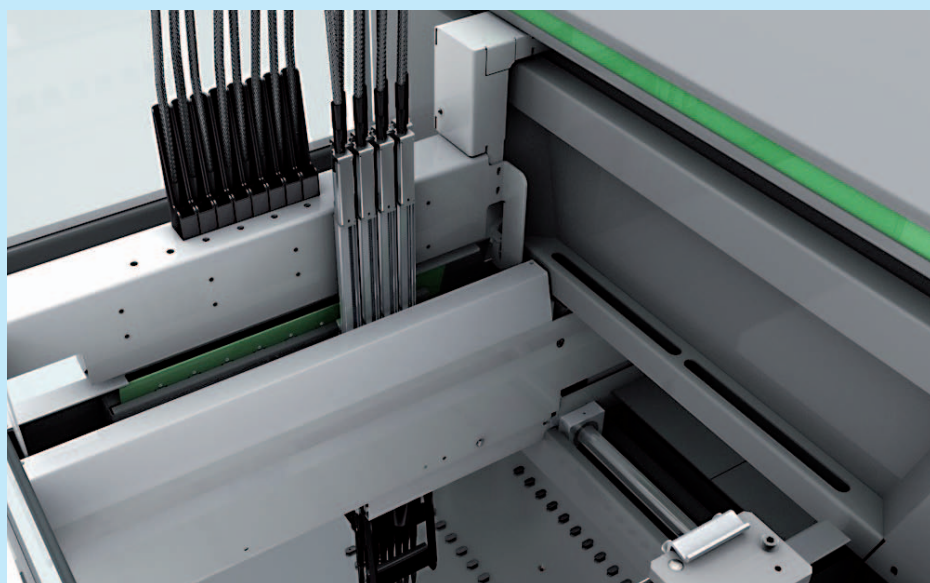
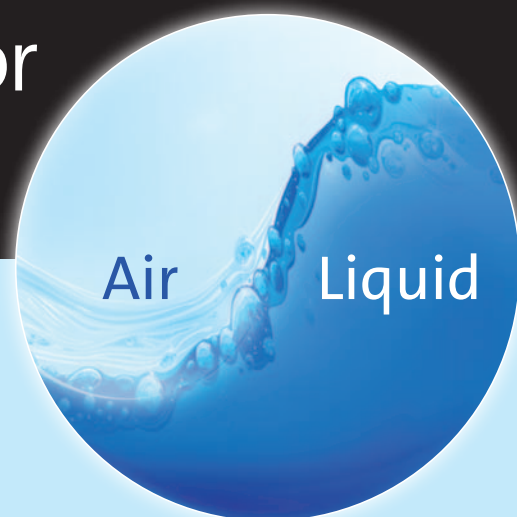


Tecan to launch Air LiHa for Freedom EVO[®] platform

With over 30 years of experience in advanced liquid handling for life sciences, Tecan understands that flexibility is crucial to making laboratory automation work for you. Our FlowThru™ Technology-based liquid displacement systems and automation expertise have established Tecan as a market leader in laboratory liquid handling solutions,

and we will soon be offering even greater choice by expanding our portfolio to include an air displacement pipetting arm for our Freedom EVO workstations.

Air and liquid displacement pipetting systems each have their own advantages. Liquid displacement offers a choice of



The new Air LiHa will provide a choice of pipetting technologies for Freedom EVO workstations

fixed or disposable tips, a large dynamic range and ultimate system cleanliness, whereas air displacement eliminates the need for system liquids and associated daily maintenance operations. The new Air Liquid Handling Arm™ (Air LiHa) for the Freedom EVO platform will provide our customers with greater versatility than ever before. By uniquely offering a choice of pipetting technologies on the same platform, or even combining both air and liquid displacement on a single workstation, Tecan will be able to provide unrivalled flexibility to suit the needs of your application and laboratory workflow and increase productivity.

To find out more on Tecan's Air LiHa, visit www.tecan.com/airliha

5 ml disposable tips now available for LiHa Arm

Tecan has further extended the flexibility of the Freedom EVO Liquid Handling (LiHa) Arm with the launch of a 5 ml disposable tip. Previously only available to Tecan customers working with fixed tips, the unrivalled pipetting volume range offered by the LiHa Arm can now be exploited by customers who require the highest levels of sample security.

The 5 ml LiHa disposable tips are available in either filtered or non-filtered options in Tecan Pure format, and are ideally suited to a wide range of applications in the pharmaceutical, academic and clinical sectors. The new tips can be used with

existing LiHa Arms, requiring only minor hardware modifications, and allow customers to take advantage of the complete range of LiHa disposable tips – from 10 µl to 5 ml sizes – using a single pipetting arm. This offers maximum throughput and flexibility for applications such as biobanking, large volume extractions and veterinary, food and clinical testing, allowing bulk liquid transfers to be rapidly performed without additional pipetting hardware.

To find out more on Tecan's 5 ml LiHa disposable tips, visit

www.tecan.com/consumables



Rack of 5 ml disposable tips for LiHa Arm