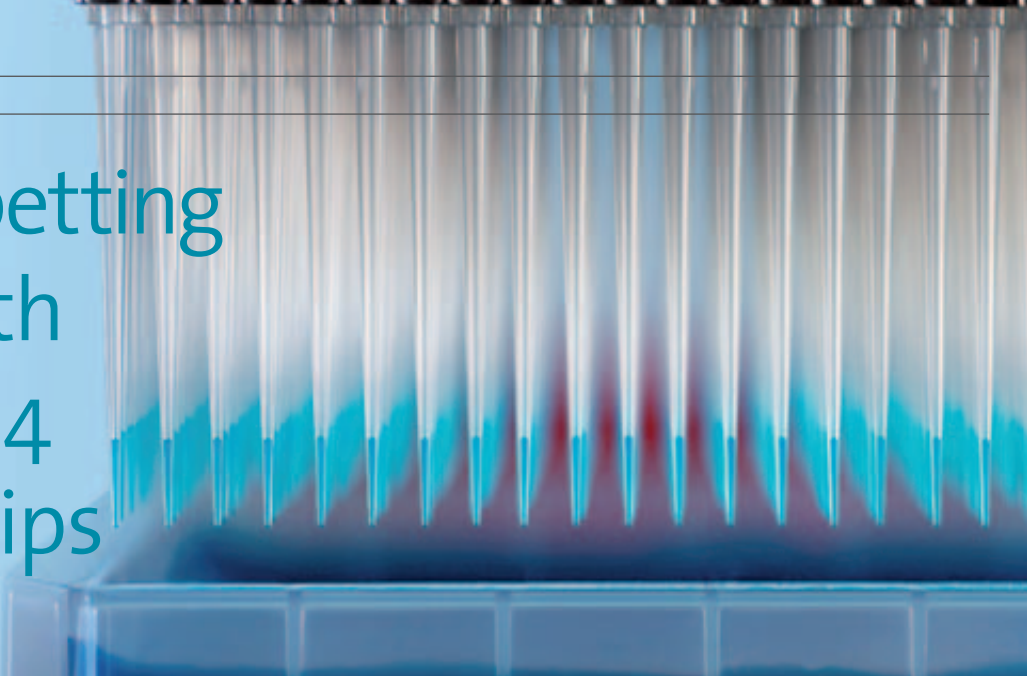


Enhanced pipetting flexibility with new MCA 384 Disposable Tips

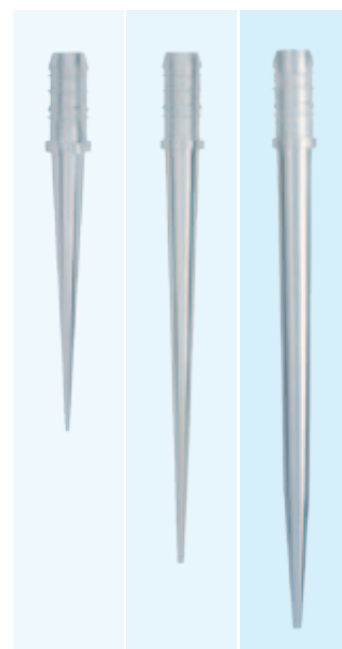


New 15 µl MCA 384 Disposable Tips with filter

Tecan has launched a new generation of high quality MCA 384 Disposable Tips designed to optimize the performance and reliability of the versatile MultiChannel Arm™ (MCA) 384 pipetting option on Freedom EVO® workstations. The all new range of MCA 384 Disposable Tips has been extended to include a 15 µl tip size alongside the existing 50 µl and 125 µl sizes, providing high quality, reliable pipetting across a broad volume range.

All MCA 384 Disposable Tips are now available in our enhanced Tecan Pure and Tecan Sterile purity levels, with filtered versions of Tecan Sterile tips also available in all sizes. Tips are precision manufactured to deliver reliable 96- or 384-channel pipetting, as well as single row or column pipetting, and have a patented design which ensures high accuracy and low CVs even at low volumes, helping to save valuable reagents and reduce maintenance requirements. The new range offers maximum format flexibility, including the use of 1,536-well plates with 15 µl tips, full depth access to 96 deep-well plates with 125 µl tips and use of typical 0.7 ml and 1.4 ml 96 storage tube formats, providing a comprehensive solution to help you meet your automation needs. In addition, each product is supplied with a Certificate of Conformity (CoC) to ensure full traceability and guarantee that the product meets the stated specifications, helping you to meet regulatory and quality documentation requirements.

To find out more on Tecan's MCA 384 Disposable Tips, visit www.tecan.com/consumables



15 µl, 50 µl and 125 µl MCA 384 Disposable Tips