

Application-focused consumables for life sciences

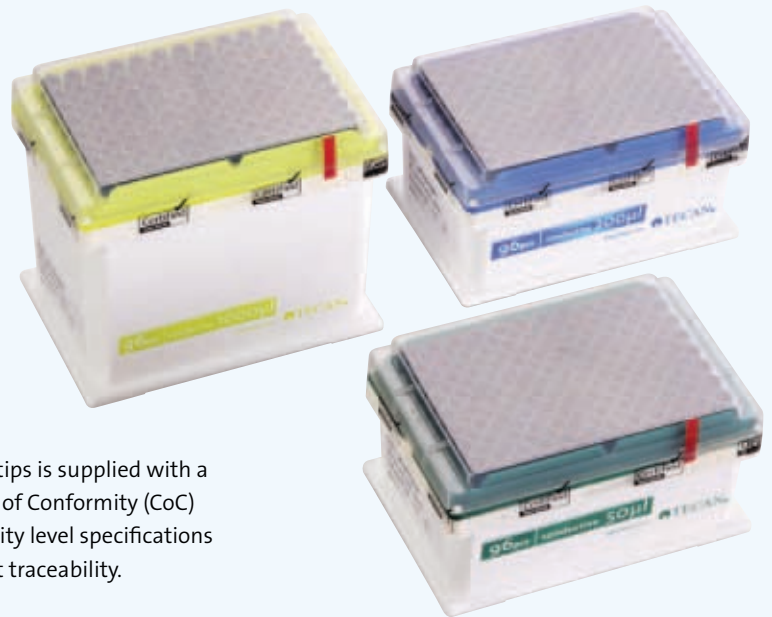
Tecan continuously strives to provide quality consumables which meet the changing needs of life sciences research. The recent trend towards high content, cell-based screening has seen an increase in demand for automation of this type of assay, and Tecan now offers disposable tips for Freedom EVO liquid handling (LiHa) arms in Tecan Sterile purity.

Previously only available in Tecan Standard or Tecan Pure formats, LiHa disposable tips are now available in Tecan's highest purity level, and are ideally suited to cell-based assays and screens, as well as microbiology applications. Available in 50, 200 and 1,000 µl volume sizes, either as filtered or non-filtered options, tip racks are individually packaged in new SBS-format boxes with an automation-friendly lid to minimize the risk of contamination. The SBS-format increases flexibility on the worktable, as well as offering higher throughput and longer walkaway times, by ensuring full compatibility with the Freedom EVO's storage and automation devices. In accordance with Tecan's high quality standards, all Tecan Sterile LiHa disposable tips are lot tested by an external, accredited laboratory to ensure tips are sterile and free from human DNA, RNase, DNase, PCR inhibitors and pyrogens/endotoxins. In

addition, each box of tips is supplied with a lot specific Certificate of Conformity (CoC) which guarantees purity level specifications and offers full product traceability.

Tecan has also recently launched a new 25 ml trough designed to minimize waste of reagents when using the LiHa arm. The conical base design of this trough reduces the dead volume for conductive liquid level detection (cLLD) by a third compared to the standard shaped Tecan trough, making it ideally suited for use with the valuable reagents now commonplace in many assays. Currently available in Tecan Pure format, this innovative yet simple product helps to reduce reagent costs while being fully compatible with existing hardware.

To find out more on Tecan's high quality consumables, visit www.tecan.com/consumables



Tecan Sterile LiHa disposable tips



New Tecan trough design minimizes dead volume