

New products and purity levels for MCA 96 disposable tips

Tecan has introduced new purity levels for the MCA 96 disposable tips, to satisfy a broader range of applications. All Tecan Pure MCA 96 disposable tips are free of human DNA, RNase, DNase and PCR inhibitors, making them suitable for genomics research and forensics use. Tecan Sterile disposable tips offer the same high purity standards, with the additional benefits of being sterile and free from pyrogens/endotoxins, and are ideally suited to cell biology, microbiology and cell-based screening applications. Nested tips for the MCA 96 are now also available in Tecan Sterile purity, extending the benefits of Tecan nested tips into these fields.

Tecan Pure and Tecan Sterile disposable tips are manufactured in a controlled environment, and subject to periodic purity testing according to USP31 and ISO 11137 guidelines. Tecan's new Certificate of Conformity provides access to detailed information on purity level specifications, lot number and expiry date, offering full product traceability.



Sterile nested tips



Tecan is also introducing wide bore tips for the MultiChannel Arm™ 96. The wider orifice of these tips minimizes shearing and fragmentation, allowing effective pipetting of fragile cell suspensions or genomic DNA. Handling of viscous fluids is also improved, because of the decreased flow resistance. The narrower aperture of conventional disposable tips can get clogged, making wide bore tips a superior alternative in areas such as crop research, where suspensions of plant material – such as leaves and seeds – are handled. They are available as non-filtered Tecan Pure or Tecan Sterile with filters, and have already completed successful external testing at Bayer CropScience in Belgium, where the filtered version is now in routine laboratory use.

Tecan Pure and Tecan Sterile disposable tips are manufactured in a controlled environment, and subject to periodic purity testing according to USP31 and ISO 11137 guidelines.

To find out more, visit www.tecan.com/purity and www.tecan.com/mca96diti

All products: Not for use in human clinical or diagnostic procedures.