

# New 50 µl disposable tip for liquid handling platforms



50 µl tips pipetting into microplate

Our family of high quality, conductive disposable tips has now been completed with the introduction of the 50 µl tip size, complementing our existing size range of 10, 200 and 1,000 µl. These disposable tips have been custom-developed for our automated liquid handling workstations, enhancing performance and minimizing cross-contamination in a broad range of genomics, forensics and ELISA-based applications.

The 50 µl tip's precision and volume range (1.0 to 50 µl) is optimal for PCR set-up and DNA quantification and normalization, as well as for other small volume pipetting applications such as ELISAs. The unique geometry of the 50 µl tip makes it compatible with many different types of labware, including microplates, deep-well plates and sample tubes; it is also able to pierce septa. Its universal nose cone is designed for our existing Freedom EVO® series of liquid handling instruments and, in particular, for the recently launched 8 Plus 1 Access™ liquid handling arm for biopharmaceutical research and the HID EVOLution™ platform for forensic applications.

Our disposable tips are produced to meet critical technical specifications, maintaining high standards of quality and reproducibility. Each tip is individually inspected and tested during manufacturing. The validated production process meets the requirements of ISO 9001 and ISO 13485.

All Tecan conductive disposable tips

- minimize risk of carryover
- reduce the chance of biological samples being contaminated (eg. RNase)
- increase throughput by minimizing tip washing steps
- enhance pipetting performance and results through optimized design
- are available with or without filter.

For more information on all our consumables products, visit our new consumables pages at [www.tecan.com/consumables](http://www.tecan.com/consumables)

●●● and don't forget

## PowerScanner™ Tecan's new flagship microarray scanner

Who wouldn't want the best possible microarray scanner in their lab?

The much anticipated, compact high resolution microarray scanning system will soon become reality: Tecan's PowerScanner™, a new generation, automated high resolution microarray scanner, is due for launch in August 2008. As the new flagship of Tecan's 'microarray suite' range, the PowerScanner features an ozone-protected environment for your valuable microarrays, and offers state-of-the-art resolution, sensitivity, accuracy and reliability for automated processing of up to 48 slides in one run.

For more information on the PowerScanner, visit

[www.tecan.com/PowerScanner](http://www.tecan.com/PowerScanner), where you can also request updates and enter a competition to win a NanoQuant Plate™ (for Infinite® 200), LaserCheck™ (for LS Reloaded™), hybridization chamber set (for HS 400™ Pro) or an Apple iPod touch 16 GB.

*Apple and iPod are trademarks of Apple Inc., registered in the US and other countries.*

