

Quality tips you can count on

New purity level certification ensures reliable results with LiHa disposable tips.

Tecan's range of high quality disposable tips is constantly evolving to help you achieve the highest performance levels on your Tecan liquid handling instruments. As well as developing new products and improving tip design, we are focused on stringent quality standards and tailoring purity levels towards specific applications.

As ever, Tecan's Liquid Handling Disposable Tips are manufactured to the highest production standards; each tip is optically inspected during production and, to improve process security and minimize the risk of user errors, tips are now supplied in trays color-coded for different tip volumes. The purity of

blister-packaged Liquid Handling Disposable Tips is now tested by an external laboratory, guaranteeing tips are free from human DNA, PCR inhibitors, RNase and DNase.

In addition, each case of tips is shipped with a Tecan Certificate of Conformity (CoC), using Tecan's validated quality assurance processes to provide full product traceability.

Free from	Specification	Method
Human DNA	<2 pg (less than in a human cell)	Polymerase chain reaction
PCR inhibitors	<10 amplifiable targets (equivalent to control)	Polymerase chain reaction
RNase	<1.0 $\times 10^{-9}$ Kunitz units	Not detectable by RNA digestion
DNase	<1.0 $\times 10^{-6}$ Kunitz units	Not detectable by DNA digestion

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

Hands-on training for microRNA profiling

Tecan has teamed up with Exiqon, manufacturers of the high specificity Locked Nucleic Acid (LNA™) detection technology, to offer customers practical training in automation of high throughput microRNA profiling. Using Exiqon's miRCURY™ LNA microRNA arrays and Tecan's HS 400™ Pro hybridization station, the course includes hands-on laboratory work to help users achieve robust and reliable results, as well as tips, tricks and presentations from experienced microRNA specialists.

To find out more about these workshops, visit www.tecan.com/microRNA

LNA and miRCURY are trademarks of Exiqon A/S, Vedbaek, Denmark.



Practical training for high throughput microRNA profiling