

# Transfer Tool for

# Nested Disposable Tips.

**INCREASE THROUGHPUT AND WALKWAY TIME WITHOUT  
REQUIRING A GRIPPER**

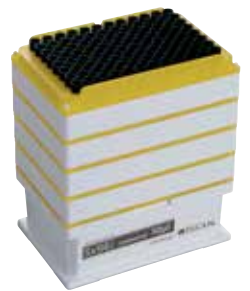
Increasing sample numbers drive many laboratories to look for opportunities to automate and optimize their liquid handling processes. One of the dilemmas that laboratories face when configuring their automated liquid handling systems is the need to balance capacity and functionality against space and budget constraints.

Tecan has introduced a cost-effective solution that makes it possible to benefit from the space-saving advantages of Nested Liquid Handling (LiHa) disposable tips without requiring a gripper to remove empty tip trays. This patent pending solution is based on the pick-up and disposal of empty tip trays using a unique consumable – the Disposable Transfer Tool – and is compatible with the LiHa and Air LiHa (Freedom EVO®) and FCA and Air FCA (FLuent®) pipetting arms.

## **Main features and benefits**

- Fully automated Nested LiHa disposable tip handling without a gripper – save equipment costs
- Five-fold increase in tip capacity per deck position – save instrument and laboratory space
- Quick set up and selection in the software\* – reduce installation time

\*Disposable Transfer Tool handling is supported by Freedom EVOware® 2.6 SP1 and FluentControl® 1.2 onwards.



## Increased workstation capacity and longer walkaway times

The Disposable Transfer Tool allows users to benefit from the extra capacity of the Nested LiHa disposable tips without needing a gripper module. When a tray of tips is empty, the pipetting arm fetches a Disposable Transfer Tool from the bespoke holder\* and uses this to transfer the empty tray to the waste.

## Straightforward implementation

Tip and empty tray handling with the Disposable Transfer Tool is conveniently automated and easily implemented through Freedom EVOware and FluentControl. Systems currently set up for non-SLAS disposable tips can be easily converted with the addition of a specific carrier. For further space-saving, carriers are also available with an integrated waste slide for tip and tray disposal.

\*The Transfer Tool Holder can be quickly and easily filled by hand prior to beginning a series of runs.

## Quality and purity

Tecan's high quality Nested LiHa disposable tips are designed to optimize pipetting performance and instrument reliability, with a patented self-centering spacer between layers for enhanced process security. The tips are available in both Tecan Pure and Tecan Sterile purity levels to match the requirements of your application.

## Nested LiHa disposable product portfolio

### Tecan Pure

Certified to be free from human DNA, DNase, RNase and PCR inhibitors.



**350 µl, Non-filtered**

300 83 400

**10 µl, Non-filtered**

301 04 977

**10 µl, filtered**

301 04 978

### Tecan Sterile

Certified to be sterile and free from endotoxins, pyrogens, human DNA, DNase and RNase.



**350 µl, Non-filtered**

300 83 401

**10 µl, Non-filtered**

301 04 979

### Disposable Transfer Tool

Disposable Transfer Tool for Nested LiHa disposable tips, compatible with LiHa or Air LiHa pipetting options.



300 83 403

### Transfer Tool Holder

Holder for 16 Disposable Transfer Tools, which fits into a three-position trough carrier.



300 92 506

[www.tecan.com](http://www.tecan.com)

**TECAN.**

**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **Denmark** +45 70 23 44 50 **France** +33 4 72 76 04 80  
**Germany** +49 79 51 94 170 **Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74  
**Sweden** +46 8 750 39 40 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33

© 2016 Tecan trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)