

Nested Liquid Handling (LiHa)

Disposable Tips.

BOOST EFFICIENCY, FLEXIBILITY AND WALKAWAY TIME WITH TECAN'S NESTED DISPOSABLE TIPS

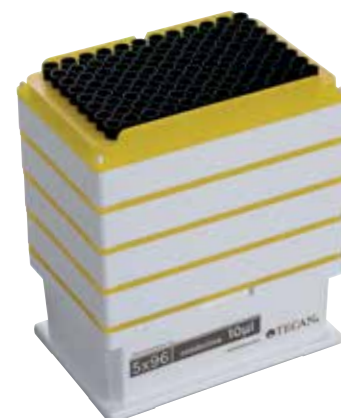


The demand for higher throughput triggers many laboratories to look for methods to increase the capacity of their automated systems. Tecan has now addressed this need by creating Nested LiHa Disposable Tips, allowing five trays of tips to be stacked on a single SLAS carrier. This solution allows optimized use of deck space through increasing tip capacity per carrier position and minimizes the need for manual intervention for Freedom EVO® and Fluent® liquid handling workstations. Flexibility of the system is enhanced as the 350 µL tip volume offers greater versatility with the same pipetting precision as the 200 µL LiHa Disposable Tip, whereas the 10 µL Tip allows pipetting in in PCR and 384 well plates.

Main features and benefits

- Five-fold increase in tip capacity on the workdeck – longer walkaway times
- Increased tip volume range & labware compatibility – higher flexibility without compromising on performance
- Quick set-up and selection in the software* – reduced installation time

*From Freedom EVOware® V2.6 and FluentControl® V1.2 onwards



Increased workstation efficiency

The nested design allows tips to be stored in stacks of five racks of 96 tips (480 tips) per SLAS position on the deck of the workstation. This increases tip capacity and/or allows integration of additional analytical modules on the instrument.

Increased tip volume range and improved labware compatibility

A broad range of 3 to 350 µl – with CVs of ≤5 % and ≤0.5 % respectively* – makes the new 350 µl nested tip the disposable tip of choice for a wide variety of applications. The 10 µl nested tip is designed in such way, that you can easily reach the bottom of a PCR or 384 well plate.

*Using the Air LiHa Arm with water or serum; methods upon request.

Straightforward implementation

The tip and rack handling for the nested tips is conveniently fully automated and easily implemented through the Freedom EVOware® and FluentControl® software. Empty tip racks can be removed and disposed of using a Robotic Manipulator (RoMa) Arm, Robotic Gripper Arm (RGA) or MCA Gripper option.

Systems currently set up for non-SLAS disposable tips can be easily converted by adding a specific carrier. For further space-saving, carriers are also available with an integrated waste slide for tip and rack disposal.

Quality and purity

Tecan's high quality Nested LiHa Disposable Tips are designed to optimize pipetting performance and instrument reliability. Two purity levels are available to match the requirements of your applications:

Nested LiHa disposable product portfolio

Tecan Pure

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



Tecan Sterile

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



350 µl, Non-filtered

300 83 400

10 µl, Non-filtered

301 04 977

10 µl, filtered

301 04 978

350 µl, Non-filtered

300 83 401

10 µl, Non-filtered

301 04 979

www.tecan.com



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

