



Disposable Tips for MultiChannel Arm™ 384

Reliable and flexible 384- and 96-channel pipetting with Tecan's quality disposable tips

The MultiChannel Arm (MCA) 384 option for Freedom EVO® workstations provides automated liquid handling with unrivalled throughput and flexibility for life science applications.

Tecan's high quality MCA 384 Disposable Tips, with certified purity levels and filtered options, have been specifically designed to support the flexibility and optimize the performance of the MultiChannel Arm 384. MCA 384 Disposable Tips have a patented design which ensures accurate, reproducible low volume pipetting results and a broad volume range, offering a reliable solution to maximize productivity.

Precision pipetting over broad volume range

The MCA 384 option is designed to maximize format and application flexibility, even allowing switching between 96- and

384-channel dispensing during a run, and the MCA 384 Disposable Tips range complements this versatile module.

MCA 384 Disposable Tips are precision manufactured to offer high accuracy (the ability to dispense the desired volume) and low CVs (the variation of pipetting results), even at low volumes. Three tip sizes – 15 µl, 50 µl and 125 µl – are available to cover a volume range of 0.5 µl to 125 µl for 96- or 384-channel pipetting, as well as single row or column pipetting.

This volume range can be further extended using the Extended Volume Adaptor (EVA), allowing rapid and automated switching to high volume 96-channel pipetting, even during a run. This offers unrivalled flexibility by allowing use of the complete range of Tecan MCA 96 Disposable Tips, including nested tips and 500 µl tips.



Precision fit ensures reliability

The advanced design of the MCA 384 Disposable Tips incorporates multiple stabilization guides and flexible sealing rings to provide excellent seal integrity without the use of O-rings. This patented sealing system maintains high pipetting accuracy and low CVs over time, while reducing maintenance requirements.

Compatibility

The MCA 384 tip dimensions are compatible with regular 384 half-deep-well plates, and 15 µl tips can also be used with standard 1,536-well plate formats. In addition, 125 µl tips offer full depth access to 96 deep-well plates and typical 0.7 ml and 1.4 ml 96 storage tube formats, allowing rapid sample transfer and assay plate generation in 96- or 384-well formats.

Certified quality

Each MCA 384 Disposable Tip is manufactured and checked in a fully automated production line to the highest quality standards.

Each lot is tested and certified to guarantee reliable quality, helping you to deliver optimal results. To ensure regulatory compliance and meet your quality documentation requirements, Tecan provides a Certificate of Conformity (CoC) with each product.

Tecan purity level and filtered options

MCA 384 Disposable Tips are available in two purity levels to suit the needs of your application

Tecan Pure

- Free from human DNA, DNase, RNase and PCR inhibitors
- Targeted towards molecular diagnostics, forensics and genomics applications

Tecan Sterile

- Sterile
- Free from human DNA, DNase, RNase, pyrogens and endotoxins
- Intended for applications in cell biology, microbiology and cell-based screening

All Tecan Sterile MCA 384 Disposable Tip sizes are also available with a filter to prevent aerosol formation.

Disposable Tip ordering information

	Tecan Pure		Tecan Sterile	
	Non-filtered	Non-filtered	Non-filtered	Filtered*
15 µl	30051802	30051803	30051804	30051804
50 µl	30051805	30051806	30051807	30051807
125 µl	30051808	30051809	30051810	30051810

* Maximum pipetting volume 14.5 µl, 40 µl and 110 µl respectively

Main features

- Broad pipetting range from 0.5 to 125 µl (up to 500 µl in 96-channel format)
- Filtered options in all sizes
- All tips are certified Tecan Pure/Tecan Sterile
- Maintenance-friendly O-ring free sealing system

Specifications of disposable tips

Pipetting of aqueous solutions¹

With COMBO Adapter in 384 format

Tip size	Range	Performance ²
15 µl	0.5 – 15 µl	CV < 4 % at 0.5 µl ³
50 µl	1.0 – 50 µl	CV < 4 % at 1.0 µl
125 µl	2.0 – 125 µl	CV < 3 % at 2.0 µl

Pipetting of DMSO solutions¹

With COMBO Adapter in 384 format

Tip size	Range	Performance ²
15 µl	0.5 – 15 µl	CV < 4 % at 0.5 µl
50 µl	0.5 – 50 µl	CV < 4 % at 0.5 µl
125 µl	2.0 – 125 µl	CV < 3 % at 2.0 µl

¹ Photometry measurement, contact dispense, method and liquid handling parameters on request

² Accuracy = + - 5 % ³ Accuracy = + - 10 %

Austria +43 62 46 89 33 Belgium +32 15 42 13 19 China +86 21 28 96 63 33 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 48 174 Singapore +65 644 41 886 Spain +34 93 490 01 74 Sweden +46 31 75 44 000
Switzerland +41 44 922 89 22 UK +44 118 9300 300 USA +1 919 361 5200 Other countries +41 44 922 8125

The MultiChannel Arm option is a scientific instrumentation. Not for use in human clinical or diagnostic procedures.

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan and Freedom EVO are registered trademarks and MultiChannel Arm is a trademark of Tecan Group Ltd., Männedorf, Switzerland.

© 2011, Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com

