



## MultiChannel Arm™ 96 option

True parallel processing for the Freedom EVO® liquid handling workstation

Tecan's 96-channel pipetting head (MCA 96) brings higher productivity to almost any automaliquid handling process, from drug discovery and compound management to biotechnology or genomics applications. The MultiChannel Arm 96 module can be mounted onto the Freedom EVO 100, 150 or 200 platforms.

### True parallel processing

The combination of the MultiChannel Arm 96 with the Freedom EVO platform optimizes your automation, increasing the speed and efficiency of your pipetting processes. By separating different functionalities on up to three independent arms (e.g. liquid handling (LiHa), robotic manipulator (RoMa) and MultiChannel), the flexible Freedom EVO workstation is designed to speed up your laboratory workflow.

For example, the use of a MultiChannel Arm, combined with an independent robotic manipulator arm, allows the 96-channel head to perform pipetting functions, while the RoMa simultaneously positions labware in preparation for subsequent pipetting sequences.

### Switch between tip types, even during a run

Tecan's MultiChannel Arm has the flexibility to work with both disposable and washable tips, which can be automatically exchanged during a run. This means you can use washable tips when appropriate, for example while pipetting reagent, and then switch to disposable tips whenever your protocol demands a new tip.



Washable fixed tips reduce running costs.

### Pipette with just 8- or 12-channels when necessary

The 96-channel head is able to pick up just one column or row of disposable tips, subsequently working as either an 8- or 12-channel pipette, enabling pipetting into a single column or row of a 96-well microplate; it can even use just one disposable tip if required, for example in a hit picking application. For applications such as IC<sub>50</sub> assays, 8 or 12 channels can be used to perform a serial dilution over a 96-well

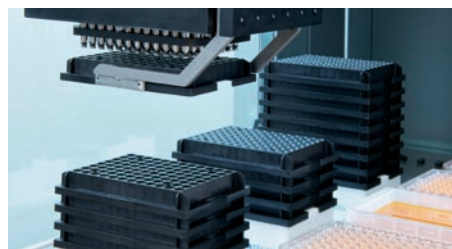


Pipetting with a single row of 12 disposable tips e.g. for serial dilutions.

plate. Used tips are discarded, and the MCA 96 can pick up further columns or rows of disposable tips to continue with the next steps of your protocol.

### Compatible with all Freedom EVO options

The MCA 96 is able to access up to 44 different SBS positions across the entire worktable, while still having plenty of space beneath the pipetting head for other Freedom EVO modules. It is compatible with a diverse range of Tecan equipment and



Optional gripper e.g. for moving disposable tip racks.

third party devices that are already integrated on the Freedom EVO. A timely addition of liquid with all 96 channels in parallel will optimize your assay results as the MCA 96 can directly pipette into labware placed on Tecan's Freedom EVO options, such as the Te-VacS™, Te-Shake™ orbital mixer or Te-MagS™ magnetic bead separator.

#### Main features:

- true parallel processing
- compatible with all Freedom EVO options
- 1-, 8- or 12-channel pipetting capability
- ability to switch between disposable and washable steel tips, and different tip sizes, during a run
- optional gripper module moves labware (microplates, tip boxes, plate transfers onto Te-VacS SPE manifold) on the worktable

## Specifications

<b>Pipetting head</b>	96-channel head					
<b>Volume range</b>	1 – 200 µl					
<b>Tips</b>	Disposable tips: 50 µl*, 100 µl*, 200 µl* and 200 µl wide bore					
	Capacity to work with:					
	1 to 12 columns of 8 disposable tips, 1 to 8 rows of 12 disposable tips, single disposable tip					
	Washable steel tips: Outer surface PTFE® coated or uncoated					
<b>Precision**</b>	Volume	50 µl disp.	100 µl disp.	200 µl disp.	50 µl uncoated wash.	200 µl uncoated wash.
	1 µl	< 6%	–	–	< 10%	–
	1,5 µl	–	< 6%	–	–	–
	2 µl	< 6%	< 6%	< 6%	< 6%	–
	5 µl	< 4%	< 4%	< 4%	< 4%	< 4%
	10 µl	< 3%	< 3%	< 3%	< 3%	< 3%
	> 10 µl	< 3%	< 3%	< 3%	< 3%	< 3%
<b>Speed</b>	Fetch tips, transfer 10 µl liquid and discard tips in 25 seconds					
<b>Options</b>	Active wash station with individual wash channels for each tip Disposable tip stacker with automated tip chute					
<b>Operating software</b>	Freedom EVOware® and Freedom EVOware Plus, version 1.4 or higher					
<b>Optional</b>	Gripper module for transporting labware on the worktable.					

\* also available with filter and sterile tips

\*\*photometry measurement of aqueous solutions, contact dispense; method available on request.

Austria +43 62 46 89 33 Belgium +32 15 42 13 19 China +86 21 289 86333 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 48 174 Portugal +351 21 000 82 16 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 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, Freedom EVO and Tecan EVOware are registered trademarks and MultiChannel Arm, Te-MagS, Te-Shake and Te-VacS are Trademarks of Tecan Group Ltd., Männedorf, Switzerland. PTFE is a registered trademark of E. I. Du Pont De Nemours, Wilmington, USA.

© 2010, Tecan Trading AG, Switzerland, all rights reserved.

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