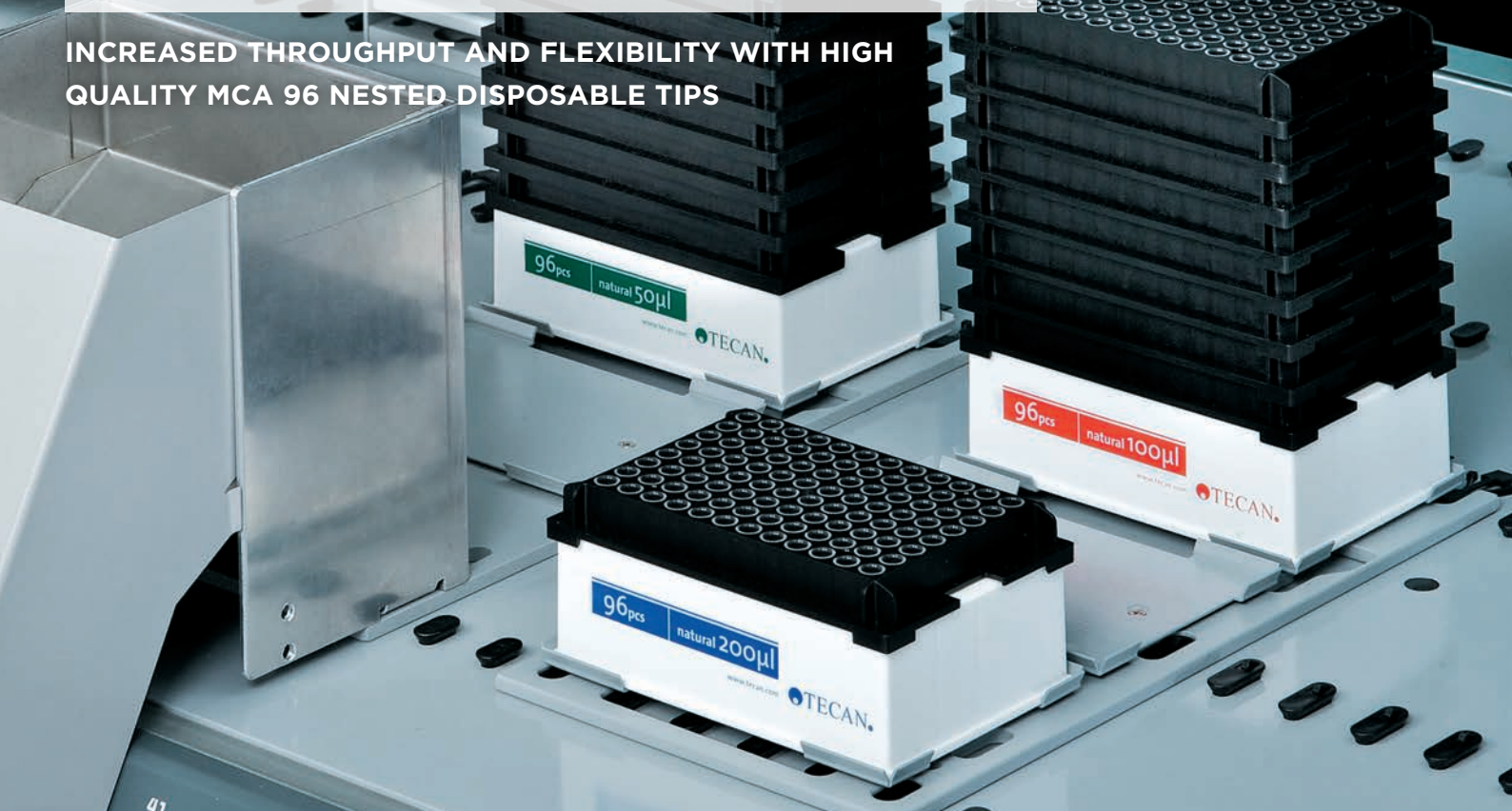


Nested MultiChannel Arm™ 96

(MCA 96) Disposable Tips.

INCREASED THROUGHPUT AND FLEXIBILITY WITH HIGH QUALITY MCA 96 NESTED DISPOSABLE TIPS



Tecan disposable tip portfolio offers a variety of pipetting solutions to complement the MultiChannel Arm 96 for Freedom EVO® and Fluent®* liquid handling workstations. With sterile, non-sterile and filtered options, the ANSI/SLAS-format product design has been optimized for increased throughput and reliability, providing a cost-effective solution for liquid handling applications.

Main features and benefits

- Improved throughput and walkaway times, thanks to nested format allowing more tips per carrier
- Increased flexibility on the workdeck, due to ANSI/SLAS format
- More environmentally friendly, through significant weight and waste reduction
- Cost effective



* Only possible with a MCA 384 arm and an Extended Volume Adapter (EVA).



Optimized workstation efficiency and higher flexibility on the deck

Non-filtered tips can be nested either on the deck in passive stacks of eight racks (768 tips) per ANSI/SLAS position or, for higher throughput requirements, externally in Te-Stack™ modules of up to 40 racks (3,840 tips) per column.

Switching to a fully ANSI/SLAS-compatible rack design increases flexibility and optimizes use of space on the workdeck. Specialized non-ANSI/SLAS labware carriers are no longer required, and systems fitted with Te-Stack modules can now use ANSI/SLAS-format columns for both tip racks and microplates.

Environmental and logistical benefits

The nested tip design simplifies storage, lowers transport costs and produces less waste – by reducing storage space requirements by 60 %, and using less packaging materials compared with non-nested formats.








Straightforward implementation

Systems currently set up for non-ANSI/SLAS disposable tips can be easily converted, and an optional waste slide for high throughput tip disposal is also available with the ANSI/SLAS-format carriers.

Quality and purity

All MCA 96 disposable tips are manufactured, tested and certified to guarantee reliable quality, helping you to acquire optimal results. Two purity levels are available to suit the needs of your application.

Nested MCA 96 disposable product portfolio

		Tecan Pure		Tecan Sterile		
		Certified to be free from human DNA, DNase, RNase and PCR inhibitors.		Certified to be sterile and free from endotoxins, pyrogens, human DNA, DNase und RNase.		
						
		Non-filtered	Non-filtered	Non-filtered	Filtered	Non-filtered
50 µl		30038606	30038609 30038610 ^b	30038607	30038608	30048822
100 µl		30038611	30038614	30038612	30038613	30048823
200 µl		30038616	30038619 30038620 ^b	30038617	30038618 ^c	30048824
200 µl Wide bore		30050348			30050349 ^c	
500 µl ^a		30046341			30046342 ^d	

a) Maximum volume only with MCA 384 with 4-to-1 Extended Volume Adapter (EVA); maximum volume on MCA 96 is 200 µl
 b) Nested for Te-Stack, c) maximum pipetting volume of 150 µl, d) maximum pipetting volume of 400 µl

www.tecan.com

Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **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 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11

© 2020 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

