

## Tecan Disposable Tips for Cavro® OEM Components

Maximized productivity and reliability with Tecan quality tips

### Quality matters

Tecan offers a wide range of disposable tips for use with the Cavro ADP and Cavro Omni Robot. Tecan Disposable Tips have been carefully designed and validated, and are produced to stringent standards of quality and reproducibility, ensuring reliable liquid handling performance. Manufacturing is fully automated in a controlled environment, using state-of-the-art production processes, integrated quality control procedures and premium raw

materials. During production, each tip undergoes in-line optical inspection using a CCD camera and image recognition software, checking all functional parts to ensure consistent pipetting accuracy and precision. The production process conforms to several ISO certification requirements, including ISO 9001 and ISO 13485. For full product traceability, each sales unit of Tecan Disposable Tips is supplied with a Certificate of Conformity (CoC), providing important product information such as part number, lot number, expiration date, and storage conditions. The CoC certifies that the product was made according to Tecan specifications, giving customers full confidence that all liquid handling steps will perform as expected to generate reliable assay results.



Tecan Disposable Tips are designed and produced to specific quality standards for highly precise and accurate pipetting with Cavro's Air Displacement Pipettor (ADP)

### Color-coded tip trays

All disposable tip trays are color-coded for easy visual determination of tip volume on the worktable, minimizing the risk of errors.

## Purity levels

Tecan Disposable Tips for Cavro OEM Components are available in two different purity levels, each tailored towards different applications:




- **Tecan Standard** purity level is targeted towards applications such as ELISA, sample handling and biochemical screening, and is supplied in space-saving cardboard packages.
- **Tecan Pure** purity level has been designed to match the needs of applications such as molecular diagnostics, forensics, and genomics research. In contrast to Tecan Standard, Tecan Pure disposable tips are packaged in blister packs, and are tested and certified to be free from human DNA, DNase, RNase, and PCR inhibitors by an external accredited laboratory using a standard PCR protocol.

### Tecan Disposable Tips purity levels

|  | Tecan Standard                                    | Tecan Pure  |
|--|---|---|
| Application examples/<br>customer segments | ELISA<br>Sample handling<br>Biochemical screening | Genomics research<br>Forensics<br>Molecular Diagnostics |
| Tecan Standard purity                      | ✓   | ✓   |
| Human DNA free                             | -   | ✓   |
| PCR inhibitor free                         | -   | ✓   |
| RNase/DNase free                           | -   | ✓   |
| Pyrogen/Endotoxin free                     | -   | -   |
| Sterile                                    | -   | -   |

## Disposable Tip product portfolio



|   | Tecan Standard                                   |            | Tecan Pure   |            |
|---|--|------------|--|------------|
| Application examples<br>customer segments   | ELISA · Sample handling<br>Biochemical screening |            | Genomics research<br>Forensics · Molecular Diagnostics |            |
| Features  | Non-filtered                                     | Filtered   | Non-filtered   | Filtered   |
| 50 µl    |  |            | 30 032 115   | 30 032 114 |
| 200 µl   | 300 644 15                                       | 300 644 16 | 30 064 848   | 30 064 849 |
| 1000 µl  | 300 644 17                                       | 300 644 18 | 30 064 860   | 30 064 861 |

.....

**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 2206 3206 **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 **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

.....

### Safety and regulatory compliance

Tecan Cavro OEM components are designed and manufactured to the highest quality standards, meeting ISO 9001 and ISO 13485 requirements, EU RoHS directives and the QSR compliant quality system. Most Tecan Cavro components are UL recognized components. Products shown are not to scale.

Tecan Group Ltd. makes every effort to include accurate and up-to-date information, 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. Changes 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 please contact your Tecan representative. This may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan and Cavro are registered trademarks of Tecan Group Ltd., Männedorf, Switzerland.

© 2013, Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

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