

FAQ

Changes to Enhance the LiHa Tip Supply Chain

Dear Tecan Valued Customers,

This FAQ addresses frequently asked questions regarding the changes we are making to enhance our products and manufacturing footprint for LiHa disposable tips. If you have further questions, please do not hesitate to contact us: <https://helpdesk.tecan.com/en/support/login>

Contents

1. General Supply Chain & Commercial	3
1.1. Why is Tecan enhancing its manufacturing footprint?	3
1.2. Are these changes permanent?	3
1.3. Will there be a change in article numbers?	3
1.4. Will hanging tip products still be available?	3
1.5. How long have the US and CH manufacturing sites been active for Tecan?	3
1.6. What are the key benefits of choosing Tecan's products?	3
1.7. Which product numbers should I use if I am a new customer or starting a new project?	4
1.8. Can I purchase transition stock?	4
1.9. How can I get product samples	4
2. Future Part Number List	5
3. Product Quality & Performance	7
3.1. Will there be any changes in product quality or performance?	7
3.2. Will there be any changes in specifications with this change?	7
3.3. What is the change in Z-position mount offset, and why does it matter?	8
3.4. Will Tecan share data demonstrating the equivalence of tip performance?	8
3.5. What do I have to do to adjust for this Z-position change?	8
3.6. Does this affect my pipetting performance or workflow?	8
3.7. What is the risk if I do not make this adaptation?	8
3.8. How can I assess the risk level of my project or workflow during this transition?	9
3.9. Will I need to revalidate assays or methods?	9
4. Logistics, Lead Times & Pricing	9
4.1. Will lead times or delivery reliability change?	9
4.2. Will pricing change because of these manufacturing changes?	9
5. Product Characteristics - Shelf Life, SKU sizes, Storage & Handling	9
5.1. Is there a change in shelf life or storage conditions for the tips?	9
5.2. Will the number of tips in a sales unit change due to the supply chain expansion?	9
5.3. Are packaging materials or labeling changing due to the site change?	9
5.4. When will the tips from the CH & US sites be available?	10
6. Sustainability	10
6.1. How does this change support Tecan's sustainability commitments?	10



FAQ

Changes to Enhance the LiHa Tip Supply Chain

7. Regulatory, Quality & Documentation	10
7.1. Do you have any IVD customers who already use tips from the expanded manufacturing sites?	10
7.2. How does Tecan ensure traceability of parts from multiple manufacturing sites?	10
7.3. Is there any change in the regulatory classification of the products?	10
7.4. Are there change control processes in place to manage this change?	11
8. Software & Labware Definition Compatibility	11
8.1. Will Tecan provide Labware Definitions for EVOware versions prior to 2.8?	11
8.2. For which versions of FluentControl will Labware Definitions be provided?	11
8.3. Will Tecan provide Labware definitions for EVOlyzer, Veya and Magni Flex?	11



FAQ

Changes to Enhance the LiHa Tip Supply Chain

1. General Supply Chain & Commercial

1.1. Why is Tecan enhancing its manufacturing footprint?

Tecan is increasing its production capacity at established sites in the United States (US) and Switzerland (CH) to strengthen supply security, increase scalability, and support stable pricing. By expanding locally, we reduce the risk of cross-border disruptions – such as tariffs and logistics interruptions – while bringing production closer to our customers and improving delivery reliability.

This initiative also supports our sustainability goals by manufacturing nearer to our customers, which helps minimize the environmental impact associated with long-distance transportation.

1.2. Are these changes permanent?

Yes, these changes are part of Tecan's long-term strategy to ensure a robust and reliable supply chain for our customers, now and in the future.

1.3. Will there be a change in article numbers?

Yes. With the exception of 10 µL and 350 µL tips, most disposable tips will transition to new article numbers to ensure clear traceability and alignment with internal and regulatory tracking. Some of the article number changes are associated with the introduction of the new ANSI SLAS format tray (Combi Tray Light™). In [Section 2](#) of this FAQ you can find a list of the new article numbers.

1.4. Will hanging tip products still be available?

Yes. Hanging tray products remain available. Please see [Section 2](#) for the product list

1.5. How long have the US and CH manufacturing sites been active for Tecan?

- United States (US): producing since 2023
- Switzerland (CH): producing since ~2003

Both manufacturing sites are qualified and validated to Tecan's standards, and all products are verified to meet Tecan's specifications and requirements.

1.6. What are the key benefits of choosing Tecan's products?

- **Reliable Manufacturing & Supply:** Tecan's global manufacturing network ensures long-term price stability, sustainability, and a reliable supply of products, minimizing the risk of disruptions and supporting consistent product availability
- **Product Enhancements:** Benefit from advanced features like the Combi Tray Light, barcoded trays, and more robust blisters for enhanced performance.
- **Regional Supply & Local Support:** With manufacturing and supply capabilities in multiple regions, Tecan offers faster delivery times and responsive local support to meet your needs.
- **Comprehensive Support:** Extensive documentation as well as product samples for testing during the transition period are available to assist you in making informed decisions.



FAQ

Changes to Enhance the LiHa Tip Supply Chain

1.7. Which product numbers should I use if I am a new customer or starting a new project?

Except for Nested and Sterile products, please use the new product numbers as outlined in [Section 2](#). For Combi Tray products, please consider we will start to introduce Combi Tray Light for SLAS format refill products in early 2026.

1.8. Can I purchase transition stock?

We would recommend that you first complete an impact assessment for your application. If there is minimal impact to your application, you should be able to transition as soon as you have completed any needed testing activities and depleted your inventory of products made in DE – for guidance on testing approaches, see 3.10. If you perceive a need for transition stock, we recommend you contact your Tecan Sales representative to discuss your needs.

1.9. How can I get product samples

Please contact your Tecan representative to organize product samples.



Changes to Enhance the LiHa Tip Supply Chain

Table 1 Part number and description changes for ANSI SLAS tray format products (blue highlighted part numbers change):

FAQ

Changes to Enhance the LiHa Tip Supply Chain

Table 2 Part number changes for Hanging tray format products (blue highlighted part numbers change):

CURRENT PART NUMBER		FUTURE PART NUMBER			
Part Number	Description	Part Number Change (Y/N)	New Part Number	Description	Implementation from
30000627	DITI LIHA 200µL CONDU. 2304 PCE.	Y	30064848	DITI LIHA 200µL CONDU.2304 PCE.	DEC 2025
30000629	DITI LIHA 200µL CONDU.FIL. 2304 PCE.	Y	30064849	DITI LIHA 200µL CONDU.FIL. 2304 PCE.	
30000630	DITI LIHA 1000µL CONDU.2304 PCE.	Y	30064860	DITI LIHA 1000µL CONDU.2304 PCE.	
30000631	DITI LIHA 1000UL CONDU.FIL.2304 PCE.	Y	30064861	DITI LIHA 1000µL CONDU.FIL.2304 PCE.	
30032114	DITI LIHA 50µL CONDU.FIL. 2304 PCE	Y	30200712	DITI LIHA 50µL CONDU.FIL. 2304 PCE	
30032115	DITI LIHA 50µL CONDU. 2304 PCE	Y	30200713	DITI LIHA 50µL CONDU. 2304 PCE	
30115239	DITI LIHA 1000UL COND.FIL. 960 PCE WIDE	Y	30201343	DITI LIHA 1000µL COND.FIL. 2304 PCE WIDE	
10612510	DITI LIHA 200µL CONDU.17280 PCE.	Y	10612552	DITI LIHA 200µL CONDU.17280 PCE.	
10612511	DITI LIHA 200µL CONDU.FIL.17280 PCE.	Y	10612553	DITI LIHA 200µL CONDU.FIL.17280 PCE.	DEC 2026
10612512	DITI LIHA 1000µL CONDU.9600 PCE.	Y	10612554	DITI LIHA 1000µL CONDU.9600 PCE.	NOV 2025
30104803	DITI LIHA 10µL CONDU. 2304 PCE.	N			N/A
30104804	DITI LIHA 10µL CONDU.FIL. 2304 PCE.	N			
10612555	DITI LIHA 1000µL CONDU.FIL.9600 PCE.	N			
10612513	DITI LIHA 1000µL CONDU.FL.9600 PCE.	N			



FAQ

Changes to Enhance the LiHa Tip Supply Chain

3. Product Quality & Performance

3.1. Will there be any changes in product quality or performance?

No. All disposable tips produced at all manufacturing sites have undergone the same thorough qualification and verification on Tecan platforms. They meet the same performance criteria for liquid-handling precision and pipetting accuracy.

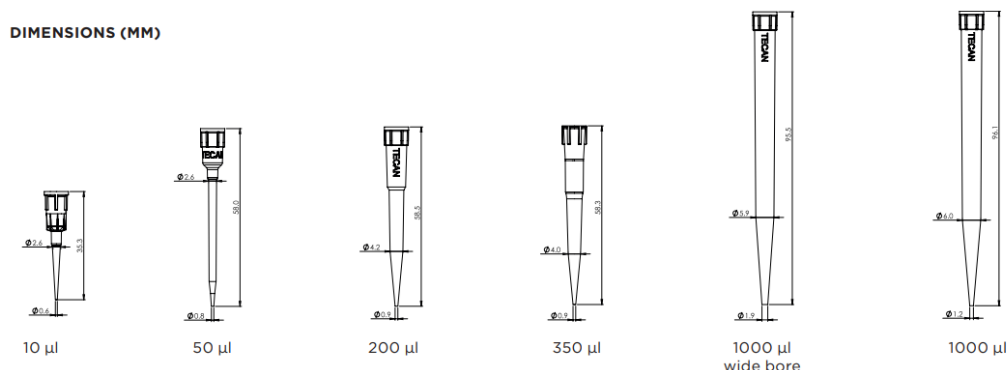
This means you can expect the same high-quality performance, reliability, and compatibility with your Tecan instruments.

All tips from CH and the US are manufactured in ISO 14644-certified cleanrooms – an upgrade from earlier controlled environments – to ensure consistent, contamination-controlled production.

3.2. Will there be any changes in specifications with this change?

Pipetting performance: Pipetting performance (CV & accuracy) of the US & CH disposable tips are within Tecan specification and equivalent to the legacy tips (country of origin DE).

Dimensions: The nominal dimensions of US & CH tips are within Tecan specifications and equivalent to the DE tips. The nominal tip dimensions are equivalent across all LiHa disposable tips:



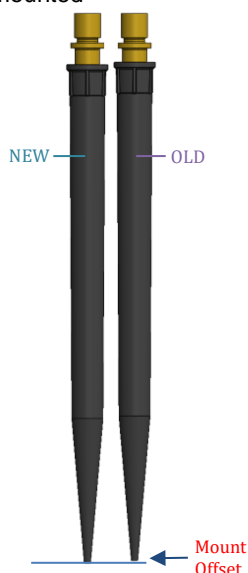
Tip mounting: The 50µL, 200µL, and 1000µL CH & US tips have a slightly thicker outer wall. This difference in the outer wall thickness means that when these tips mount onto the cone, there is a slight difference in the mounting compared to DE tips.

This results in a difference in the Z-position (see Figure 1), which is observed as a change in the nominal Z-axis, leading to a lower position of the mounted tips. The amount of this minimal offset in the Z-axis depends on the platform & tip size (Note: Z-position accuracy is not currently specified). Please refer to the LiHa data sheet 399504 for further information on product specifications. **Please be aware that we strongly recommend an update of system settings according to section 3.5 when using the disposable tips with the new part numbers.**

FAQ

Changes to Enhance the LiHa Tip Supply Chain

Figure 1 Mount offset: There is a 0.3 - 0.6 mm Z-position difference when products from the US & CH (NEW) are mounted



3.3. What is the change in Z-position mount offset, and why does it matter?

Tecan does not specify the absolute Z-position of the tip during pipetting, as this value depends on multiple system-specific factors. While the instrument manuals (e.g., *Fluent Operator Manual*, *Freedom EVO Manual*, and *Veya Labware Definition Guide*) provide Z-coordinates for the instrument only, they do not include tolerances for disposable tips.

For precision-sensitive applications, it is recommended to verify whether an adjustment of the z-positions is required. Tecan supports this process by providing preconfigured Labware Definitions reflecting the updated nominal Z-offset values. See below (3.5) for information about adjusting Z positioning.

3.4. Will Tecan share data demonstrating the equivalence of tip performance?

Yes, if required, a product comparison and equivalence statement can be provided.

3.5. What do I have to do to adjust for this Z-position change?

When making the switch to products manufactured in CH or the US, please follow the instructions on this web page to download and import labware definitions or make adjustments to your instrument: <https://tecan-link.com/LabDev>

3.6. Does this affect my pipetting performance or workflow?

If the Z-position is adjusted in the software, no impact is expected in system performance on Tecan platforms. The disposable tips continue to meet all defined performance specifications. Nonetheless, for critical applications or workflows, customer confirmation of workflow performance is recommended (e.g. assessment, testing, validation etc.).

3.7. What is the risk if I do not make this adaptation?



FAQ

Changes to Enhance the LiHa Tip Supply Chain

For applications or workflows that have steps going to Z-max with a 50µL, 200µL, or 1000µL tip from the US or CH using the incorrect labware definition, it is possible that the tip may reach the bottom of the labware, or you may have an incorrect aspiration.

3.8. How can I assess the risk level of my project or workflow during this transition?

To assess the risk, you need to review the details of your respective workflow:

If you use liquid level detection, the risk from the lower Z-position difference is low. In this case, you just need to update to the new labware definitions.

If you use Z-Max in your application with a legacy 50 µL, 200 µL, or 1000 µL tip, the risk is higher when switching to US or CH tips. In this case, updating to the new labware definitions and performing a functional verification to ensure proper operation is strongly recommended.

3.9. Will I need to revalidate assays or methods?

Tecan will provide documentation and support to assist with any required verification or requalification activities if needed. For highly regulated environments (e.g., IVD), you may need to perform a functional verification and/or validation, and Tecan is available to assist with documentation or qualification data as needed.

4. Logistics, Lead Times & Pricing

4.1. Will lead times or delivery reliability change?

Lead times are expected to improve or remain stable because of the expanded capacity, additional supply flexibility, and shorter delivery distances for many regions.

4.2. Will pricing change because of these manufacturing changes?

No, the manufacturing changes themselves will not directly impact pricing. By diversifying our manufacturing network and reducing reliance on cross-border shipments affected by tariffs, we are better positioned to absorb or mitigate cost increases. Regional production also helps Tecan reduce internal logistics costs. While these steps support our commitment to long-term price stability, standard price adjustments, such as those related to inflation (CPI), may still apply.

5. Product Characteristics - Shelf Life, SKU sizes, Storage & Handling

5.1. Is there a change in shelf life or storage conditions for the tips?

No. The shelf life and recommended storage conditions for Tecan disposable tips remain unchanged, regardless of manufacturing site. All packaging and product labeling continue to carry the same specifications.

5.2. Will the number of tips in a sales unit change due to the supply chain expansion?

No, the number of tips in a sales unit will remain the same.

5.3. Are packaging materials or labeling changing due to the site change?



FAQ

Changes to Enhance the LiHa Tip Supply Chain

While article numbers will change, packaging materials and labeling standards will continue to follow the same specifications as before. The label will solely include the country of origin (COO) and new article numbers (both where applicable). The COO may be either CH or the US and can be differentiated by the Lot number.

5.4. When will the tips from the CH & US sites be available?

The following product formats are already available:

CH site:

Standard Purity Products (cardboard envelope) in Hanging Tray, Filtered & Non-Filtered

US site:

Pure Purity Products (blister refill) in Hanging Tray, Filtered

Pure Purity Products (blister refill) in Combi Tray, Filtered – **Note:** will change to Combi Tray Light

For certain products, the transition and introduction timelines vary by part number and region. The planned transition dates, from which we anticipate having limited availability of legacy part numbers, are outlined in [Section 2](#).

6. Sustainability

6.1. How does this change support Tecan's sustainability commitments?

By manufacturing closer to customers, Tecan can significantly reduce the environmental impact associated with long-distance transportation and international shipping.

The distributed, regionalized production model also improves scalability and allows for faster production ramp-ups when needed, enhancing both sustainability and responsiveness to customers' orders.

7. Regulatory, Quality & Documentation

7.1. Do you have any IVD customers who already use tips from the expanded manufacturing sites?

Tecan has various customers already using tips from the expanded manufacturing sites.

7.2. How does Tecan ensure traceability of parts from multiple manufacturing sites?

Tecan ensures that all parts are fully traceable down to the specific production site, either through the part number or the Lot number format. This allows us to identify the origin of each part and respond quickly to any site-specific issues.

7.3. Is there any change in the regulatory classification of the products?

For US:

No. Tecan's catalogue consumable products themselves are not classified as medical devices, and most LiHa products are Class 1, 510(k) exempt.



FAQ

Changes to Enhance the LiHa Tip Supply Chain

For EU:

No. Tecan's catalogue consumable products themselves are not classified as medical devices and most LiHa products are either Class A according to (EU) 2017/746 or General Purpose.

7.4. Are there change control processes in place to manage this change?

Yes, these changes are managed through internal assessment and change management processes.

8. Software & Labware Definition Compatibility

8.1. Will Tecan provide Labware Definitions for EVOware versions prior to 2.8?

No. Labware definitions are only available for EVOware 2.8 and can be used with EVOware 2.8 SP7. For prior versions of EVOware software, manual adjustments will need to be made. Instructions for this are available here: <https://tecan-link.com/LabDev>

8.2. For which versions of FluentControl will Labware Definitions be provided?

We aim to support FluentControl versions back to 1.6.

8.3. Will Tecan provide Labware definitions for EVOlyzer, Veya and Magni Flex?

We will communicate solutions for these platforms or components at a later date.

