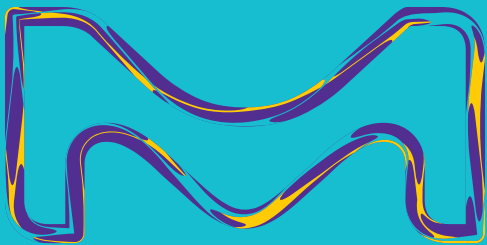
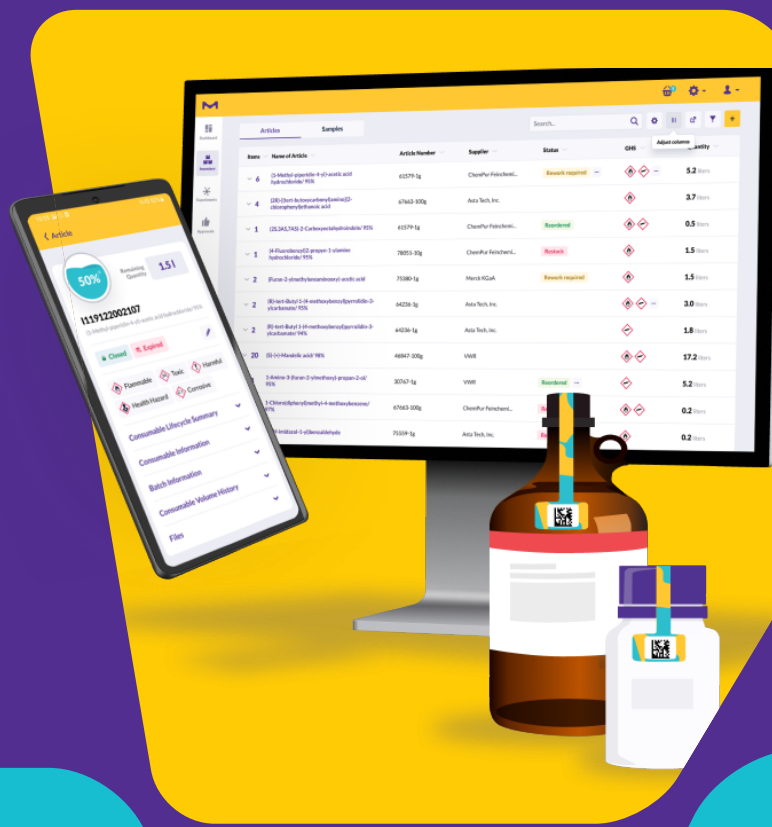


**All about the
LANEXO® System**

**YOU HAVE QUESTIONS.
WE HAVE ANSWERS.**



Want to tell your colleagues all about what the LANEXO® System can do for your lab? Got a question that occurred to you after the webinar or trade show? We've got you covered. Below, we answer the most common questions about the LANEXO® System.

1. Why do I need the LANEXO® System?

If you manage your laboratory inventory manually, you're throwing away money and time, and you may be putting your business at risk.

- Routine manual lab inventory upkeep takes up to 25% of scientists' time—time that could instead be devoted to scientific discovery.
- Transcription errors during manual registration cascade throughout the whole workflow, threatening data integrity and reproducibility and causing delays.
- Disorganized inventory records make predicting reagent usage nearly impossible, leading to overstocking and wasted materials.
- Lack of easy oversight also hinders root cause analysis.
- Poorly organized inventory puts your quality and safety compliance at risk.

The LANEXO® System **frees your scientists from the tedious, time-consuming administrative work of managing your chemical inventory**, so they can do science instead. Your staff can easily register and monitor consumables, their storage locations and their expiry dates using radiofrequency identification (RFID) labels that they place on bottles and vials and on cabinets, along with web and mobile applications.

2. What is the LANEXO® System?

The LANEXO® Lab Inventory, Safety and Compliance Management System is a Software-as-a-Service (SaaS) solution specifically designed for regulated QC laboratories: Pharma, Biopharma (incl. large, small/mid-sized and biotech), Industrial Testing and CROs/CDMOs.

The system consists of:

- An Android app that lets you register consumables simply by scanning them with a smartphone and that give you access to all the system's capabilities.
- A full range of radiofrequency identification (RFID) labels you can apply to consumables from any vendor, as well as your own in-house preparations:
 - Smart Seals (a) that detect when bottles are opened and automatically calculates their expiry dates based on opening date.
 - Identifier labels (b) for reagents, buffers and other consumables you prepare in-house—even samples.
 - Position labels (c) for tracking storage locations: cabinets, shelves, refrigerators, freezers, etc.
- A Web app that lets you browse consumables, set Rules for safety and restocking, set user permissions and search/report audit trails.
- User access cards (d) so that every user can log in to the app using a unique digital signature, for full 21 CFR Part 11 compliance.



a. Smart Seal



b. Identifier



c. Position label



d. User access card

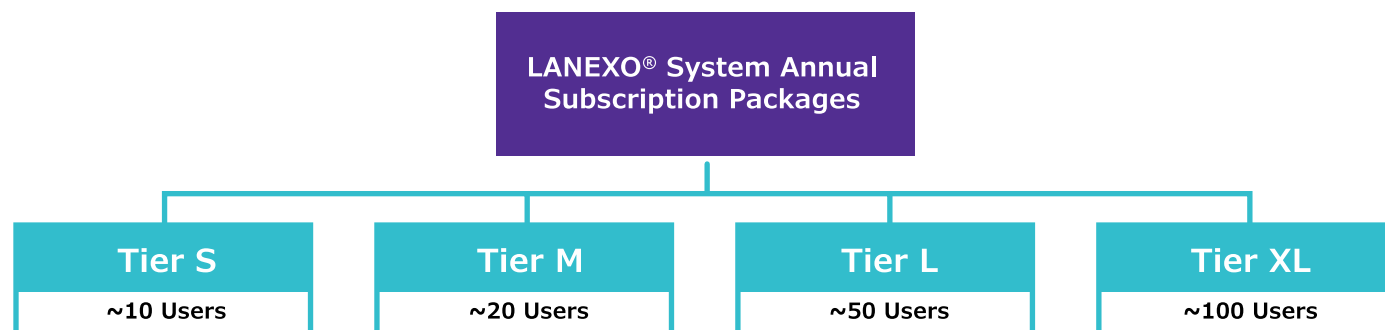
3. What subscription options do you offer for the LANEXO® System?

For annual LANEXO® System subscriptions, we offer package tiers from which you can choose based on your specific needs—that is, the number of users in your lab and thus the number of RFID labels you'll need.

Trials are available, letting you explore all LANEXO® System features and determine the exact yearly subscription that's right for your lab. Also included are:

- Approximately 6 hours of user training, depending on need.
- Weekly health checks on support status and product quality.
- Quarterly and annual business review meetings with all relevant stakeholders.

We recommend speaking with a **Lab Informatics Sales Specialist** for more details.



4. What are the specific benefits of the LANEXO® System?

Higher Lab Productivity

- **Get fast digital access to data**—such as stock levels, location, chain of custody, and documentation such as SDS and CoAs—for all reagents, including those prepared in-house.
- **Spend less time** on repetitive, error-prone operational and admin tasks like manual inventory control and expiration date monitoring.
- **Improve shelf life monitoring** with RFID “Smart Seals” that auto-register opening dates and calculate expiry dates, saving you the costs of buying new reagents and disposing of the expired ones.

Quality, Safety & Compliance Monitoring

- **Lower the risk of human error** and improve the reliability and traceability of your compliance documentation (reagent data, experimental workflows, etc.) with digital data capture.
- **Facilitate GxP compliance and be more audit ready:** Each reagent, including the ones you prepare in-house, can be instantly matched to a full audit report.
- **Avoid using expired reagents** thanks to easy identity checks and automated alerts along experimental workflows, minimizing errors and saving you from having to repeat experiments.
- **Improve safety and regulatory compliance** with real-time storage monitoring and automatic alerts in case of deviations (e.g., incompatible chemicals stored together) or low stock levels.
- **One centralized system for all data**, creating a single source of truth for your entire lab.

Easy to Set Up, Easy to Use

- **Rapid integration** into your existing workflows.
- **Easily—and accurately—capture data** with user-friendly Android and Web applications for adoption by all users, plus RFID labels.
- **Custom-developed RFID Smart Seals** developed with your lab's requirements in mind.
- **Achieve complete lab coverage** with a vendor-neutral system.
- **Harmonize access for all authorized users** through a secure web interface.

5. Do I need to make major investments in IT, or change anything else?

Nope! Implementing LANEXO® System in your lab is an easy four-step process:



1. Order your LANEXO® System trial and receive a package of RFID labels.



2. Schedule onboarding training for all authorized users.



3. Put labels on your consumables and storage locations.



4. You're good to go!

6. How can I get in touch with you?

You can reach out to us at:



Web:

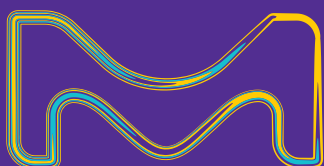
- www.lanexosystem.com
- www.SigmaAldrich.com/LANEXO



Email:

- North America: LANEXO@emdgroup.com

MilliporeSigma
400 Summit Drive
Burlington, MA 01803



MilliporeSigma is the U.S. and Canada Life Science business of Merck KGaA, Darmstadt, Germany.

© 2022 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M and LANEXO® System are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

MS_FL9876EN Ver. 1.0
43655 08/2022