



Global quality and regulatory affairs.

THE ROLE OF QUALITY MANAGEMENT IN SCALING HEALTHCARE
INNOVATION, FROM LIFE SCIENCE TO THE CLINIC.



Quality management

in the century of biology.



**Message from Günter Weisshaar,
Head of Quality and Regulatory Affairs, Tecan:**

Quality management continues to evolve. The healthcare sector needs quality and compliance systems that are adaptable and precise, capable of making swift decisions in real time.

Quality management and regulatory compliance systems are becoming integral to the digital ecosystems used to connect stakeholders, technologies and data sources, and ensure seamless collaboration, efficient information sharing and enhanced decision-making.

The trend for digitalization brings:

- **Advanced data analytics** enabling users to capture, analyze and share quality data in real time.
- **Automated quality processes** that use digitalized management modules to avoid human error in auditing, risk- and document management, etc.
- **Virtual quality work**, enabling remote audits, virtual-reality training and remote service calls.
- **Qualification and incentivization** of in-house experts to meet the ever-advancing needs of digital quality management within the broader digital ecosystem.
- **Connectivity** to broader digital ecosystems, for the ability to share information with customers, suppliers, and internal and external partners in a timely manner.

We are committed to continuous improvement and maintaining the highest standards of performance and regulatory compliance. We work to ensure the scalability and adaptability of our quality management systems to meet our customers' needs, wherever they are in the world.

Expertise and agility in quality and regulatory affairs.



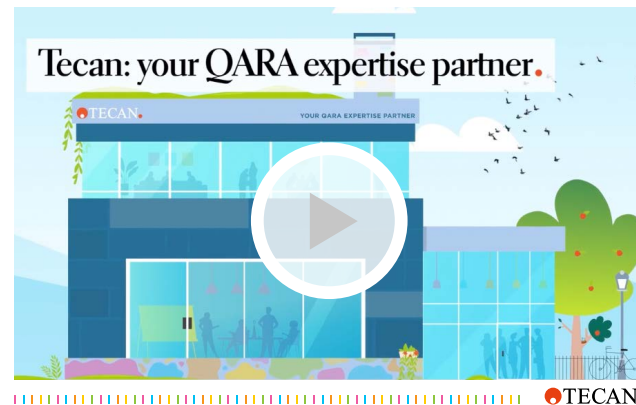
At Tecan, we offer unparalleled expertise in quality and regulatory affairs. Our skilled team is well-versed in the standards, regulations and frameworks that govern our industries. This knowledge is an essential component of our order fulfillment and quality management systems (QMS). It ensures that both we and our customers remain fully compliant.

And as the regulatory landscape evolves, we've fostered continuous learning. We're an active participant in regulatory and industry forums and, working with our partners, we've become an influential player, shaping new standards in IVD* and in medical devices.

These skills enabled us to become one of the first companies to gain IVDR** certification, well ahead of the EU's 2022 deadline. In just 12 months, we achieved certification for 21 products and lent our support to several partner companies. With our support, they completed their regulatory work and obtained IVDR certification on time.

*In Vitro Diagnostic Directive, **In Vitro Diagnostic Regulation

“
**Enhance your business with our
quality and regulatory expertise.**
”



WATCH NOW



Empowering customers.



As technology advances and regulations change, it's crucial to partner with an expert who understands the trends and stays ahead of requirements. Tecan is that partner. Knowledgeable and innovative, we have the agility to meet changing needs with minimal disruption to existing processes.

When COVID-19 took the world by surprise, Tecan was ready to help. The power of our OEM partnership was exemplified by Thermo Fisher Scientific's 'Amplitude™ Solution'. With Tecan's support, this high-throughput, automated PCR testing system for SARS-CoV-2 was developed and launched in record time, fully in line with all regulatory requirements.

We collaborate with our customers and support them on their quality and regulatory journey, and our products, services and solutions make lab and clinical procedures reproducible, scalable, and compliant.



To learn how we contribute to healthcare and the life sciences, visit the Tecan Journal and read our customer stories.

www.tecan.com/tecan-journal

Continuing advancement in quality and regulatory affairs.

The future of quality and regulatory affairs is digital. We are helping our customers to translate the constant and increasing flow of digital information into actionable insights, to support the long-term sustainability of their business.

We are committed to maintaining regulatory compliance through the course of this digital transformation. In line with our sustainability promise, we ensure that our business practices, like our products, add value to society.



For Tecan's sustainability policies, VISIT
www.tecan.com/annual-reports

> Tecan - Who we are

Tecan (www.tecan.com) improves people's lives and health by empowering customers to scale healthcare innovation globally from life science to the clinic. Tecan is a pioneer and global leader in laboratory automation. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments, components and medical devices that are then distributed by partner companies. Founded in Switzerland in 1980, the company has more than 3,000 employees, with manufacturing, research and development sites in Europe, North America and Asia, and maintains a sales and service network in over 70 countries. In 2022, Tecan generated sales of CHF 1,144 million (USD 1,192 million; EUR 1,144 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).

> Australia

+61 39 647 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

> 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 Trading AG, Switzerland, all rights reserved.

www.tecan.com

