

Tecan Group achieves ISO accreditation



Günter Weisshaar, Senior Vice President, Head of Quality Assurance and Regulatory Affairs

In recent years, Tecan has successfully embarked on a program to achieve ISO certification for all its global manufacturing and sales organizations, and now the Tecan Group parent company has been certified to ISO 13485 standard for its web-based management system.

The new, simpler and more efficient system ensures that Tecan employees have the time to concentrate their efforts on customer requirements and product developments. Acting at a global level, it provides a wealth of corporate information, document templates, SOPs and directives to all employees worldwide. The system is based around a comprehensive information repository that maintains and archives all documents, so that employees can be sure they have access to current and valid texts. The whole process is now paperless, completely automated and controlled with electronic signatures, from a request to change a management document to its release and upload onto the system. And, at every stage, the system saves time by notifying each employee of pending tasks.

The system also features a regularly updated organizational matrix chart of all Tecan employees and their roles, so that it is immediately clear who in the Company is best served to help in any situation, especially important in such a large multinational company. In addition, a series of iDashboards can be set up to monitor data, key performance indicators and analysis (PIAs), quality data, etc., and automatically upload the data in a standardized reporting format.

As a whole, the system has been a huge success and customers are undoubtedly benefiting from the increased efficiency and up-to-date information they now receive. It is a dynamic system with even more potential, and we at Tecan will constantly look for ways to improve.

Tecan locations



- Tecan sales office
- R&D and manufacturing site