

# Tecan Company Policy Statement

---

## Product Environmental Compliance

We are committed to providing our customers with safe, quality, and environmentally friendly products and to comply with all relevant product environmental legislations. We strive to continually identify and realize opportunities to reduce the environmental footprint of our products during product design, manufacture, use and disposal.

This statement specifically describes our commitment how we comply with the following laws and regulations:

- The European Union (EU) Directive on “**R**estriction of use of certain **H**azardous **S**ubstances in electrical and electronic equipment”, 2011/65/EU (RoHS2 Directive),
- The Management Methods for Restricted Use of Hazardous Substances in Electrical and Electronic Products (China RoHS),
- The European Union (EU) Regulation EC 1907/2006 on **R**egistration, **E**valuation, **A**uthorization and **R**estriction of **C**hemicals (REACH Regulation),
- The European Union (EU) Directive on **W**aste **E**lectrical and **E**lectronic **E**quipment Directive, 2012/19/EU (WEEE Directive),
- The U.S. Act – Section 1502 of the Dodd-Frank Wall Street Reform and Consumer Protection Act of 2010

We monitor changes and applicability of the above regulations to ensure continued compliance and we will update this statement as required accordingly.

## EU RoHS2 Directive

The EU enacted legislation to reduce the use of certain hazardous substances in electrical and electronic equipment being sold in the EU. Directive 2011/65/EU (RoHS2) bans the use of lead, mercury, hexavalent chromium, polybrominated biphenyls (PBB), and polybrominated diphenyl ethers (PBDE) above 0.1% and the use of cadmium above 0.01% in homogenous materials in electrical and electronic equipment.

Tecan products affected by EU RoHS2 fall within Category 8 (“medical devices”) or Category 9 (“monitoring and control instruments”) or are considered as electrical or electronic components.

Tecan complies with all requirements of EU RoHS2.

## China RoHS

China has published “The Management Methods for Restricted Use of Hazardous Substances in Electrical and Electronic Products” on January 6th, 2016 which went into effect on July 1st, 2016. The substances addressed and their thresholds match with those of EU RoHS2. The key difference is however that China RoHS does not ban these substances but mandates specific declaration and labelling if present above the threshold.

To comply with these requirements Tecan labels its equipment imported into China with the China RoHS logo specifying the “environment friendly use period” and secures that the table of hazardous substances in the format defined in standard SJ/T 11364-2014 accompanies the equipment.

## The EU REACH Regulation

With the EU REACH regulations (EC 1907/2006), the EU established on June 1st, 2007, an integrated system for the **R**egistration, **E**valuation, **A**uthorization and restriction of **C**hemicals (REACH).

REACH requires companies that manufacture or import chemicals in quantities exceeding 1t per substance to evaluate the risks to human health and environment resulting from the use of those chemicals and to take the necessary steps to manage any identified risk.

As Tecan does neither manufacture nor imports chemicals in quantities exceeding 1t it has no registration obligations under REACH. For its immunoassay business Tecan provides the safety data sheets or the necessary information to its customer as defined in Art. 31 & 32.

Tecan secures compliance with REACH substance restriction (Annex XVII) for all its products.

As a manufacturer of articles Tecan is committed to fulfil the duty to communicate information on the presence of substances of very high concern (SVHC) listed on the actual REACH candidate list as defined in Art. 33.

## The EU WEEE Directive

We are committed to and in support of compliance with the EU Directive on Waste Electrical and Electronic Equipment (WEEE – 2012/19/EU). We strive to ensure that our products are appropriately disposed of or recycled at the end of their lifespan. Tecan products that are subject to WEEE are marked with the “crossed-out wheel bin” WEEE symbol to indicate that the product should not be discarded along with other household waste but must be collected and treated separately.

## Conflict minerals

Tecan’s ultimate goal is to prevent the use of conflict minerals (gold, tin, tantalum, or tungsten) that originate from sources whose profits support armed groups and human rights abuses in the Democratic Republic of the Congo (DCR) and adjoining countries, yet support the use of materials from legitimate sources.

The US Dodd Frank Act (section 1502) from 2010, addressing US publicly listed companies, is the first legislation that tackles this problem. Even if not affected directly, Tecan is committed to support its customers to comply with this legislation and adheres to the applicable elements of the OECD Due Diligence Guidance for responsible supply chains of minerals from conflict affected and high-risk areas.

**Please do not hesitate to contact your Tecan sales representative, if you have any further questions or concerns on the above.**

DocuSigned by:

*Achim Von Leoprechting*

Signer Name: Achim Von Leoprechting  
Signing Reason: I have reviewed this document  
Signing Time: 2021-02-23 | 1:24:15 PM CET

B973F977F55044EE93C69557F875F539

---

Dr. Achim von Leoprechting, CEO

DocuSigned by:

*Günter Weisshaar*

Name des Unterzeichners: Günter Weisshaar  
Signiergrund: Ich habe dieses Dokument geprüft  
Signierzeit: 2021-02-23 | 12:08:20 PM CET

3249E468E1E442B78570095698CAE5B8

---

Günter Weisshaar, Senior VP QA/RA