

## **TAX PRINCIPLES 2023 - 2024**

### **Introduction**

Tecan is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics and clinical diagnostics. It is our mission to improve people's lives and health, by empowering our customers to scale healthcare innovation globally from life science to the clinic. We live our core business values "Ambition", "Highest standards" and "Trust".

Our behavior is governed by strict adherence to our ethical code, respect for our environment, and full compliance to applicable laws and regulations in all the jurisdictions where we operate. This is no different when it comes to taxation, which is an integral element of our overall corporate social responsibility.

Our Tax Principles are in line with our core business values and are designed to support Tecan in delivering its strategic ambitions.

### **Principle 1 Sustainability & Governance**

Tax is a core part of Corporate Sustainability and Governance and is overseen by the Board of Directors (the "Board"). The Group's Tax Principles are in line with the goals of the OECD/G20 Base Erosion and Profit Shifting project and with its core principles, coherence, substance and transparency.

The Board of Tecan regularly reviews the effectiveness of the Tax Principles and maintains a sound system of risk management and internal control. The Board is updated annually on tax risks and adherence to the Tax Principles.

Our Tax Principles are mandatory and apply to all the entities and employees of the Group.

### **Principle 2 Compliance**

We comply with the tax legislation of the jurisdictions in which we operate, adhering to both, its letter and spirit, and pay the right amount of tax at the right time.

All tax returns, claims, elections, disclosures, and payments shall be made accurately and on time.

The Tecan Group Transfer Pricing Policy is defined and implemented based on the internationally accepted arm's length principle, as described by the international tax conventions and by the OECD Guidelines, and as implemented in local rules and regulations. Transfer pricing methods follow a thorough analysis of the functions, risks and assets of the parties to the transaction.

To ensure that the Group complies with local tax laws in the jurisdictions in which it operates, and that solid and responsible tax planning is undertaken, we seek for adequate personnel, resources, up-to-date expertise, training and systems, and develop tax awareness across Tecan functions and businesses.

### **Principle 3 Business structure driven by commercial considerations**

Our business structure is driven by commercial considerations, is aligned with business activity and has genuine substance.

Our tax planning is based on reasonable, solid interpretations of applicable law and is aligned with the substance of the economic and commercial activity of our business. We do not use tax havens or opaque corporate structures to hide or reduce the transparency of our activities.

Our Tax Principles extend to our relationships with employees, customers and suppliers. We do not engage in arrangements whose sole purpose is to create a tax benefit.

#### **Principle 4 Constructive and Professional Relationships with Tax Authorities**

It is the Group's aim to maintain constructive and professional relationships with local tax authorities, based on mutual respect, transparency and trust.

We are open and transparent with tax authorities, responding to relevant tax authority enquiries in a straightforward and timely manner to assist in the evaluation of tax liability.

Where there are misunderstandings of fact or law, we will seek to work with tax authorities, where possible, to identify the issues and explore options to resolve those misunderstandings or disagreements.

We will not bribe or otherwise induce tax officials, government officials or ministers with the aim of obtaining more beneficial outcomes with respect to tax matters.

#### **Principle 5 Transparency**

We strive for a regular information to our stakeholders, including investors, employees, civil society and the general public, about our approach to tax and taxes paid. Disclosures are made in accordance with the relevant domestic regulations, as well as applicable reporting requirements and standards such as IFRS.

This document is produced in accordance with the requirements of Part 2 of Schedule 19 of the UK Finance Act 2016 and applies to our Tecan affiliate in the UK.