

SUSTAINABILITY AT TECAN

10 KEY POINTS



01. Sustainability Governance

Tecan has a Sustainability Committee, chaired by our CEO, and a global sustainability program. Our Sustainability Report, part of Tecan's Annual Report, is drafted to GRI (Global Reporting Initiative) standards and is available on [Tecan.com](https://www.tecan.com).

Tecan has a defined approach to managing our climate impact and the transition to a circular economy, including by implementing ecodesign principles in product research and development, and optimizing product lifespan.

02. Environment




03. Environment

Tecan has near-term and long-term scope 1, 2 and 3 greenhouse gas emissions reduction targets validated by the Science Based Targets initiative and has used 100% renewable electricity since 2025.

Tecan strives to create an innovative and positive workplace culture, enabling a beneficial social impact and strengthening business performance. Tecan has global programs and has been certified a Great Place to Work™ in seven countries, covering 92% of our global workforce.

04. Social Impact




05. Social Impact

We are building a culture where all individuals at Tecan can “stay unique and make it count.” We are committed to encouraging everyone to fully participate and drive an inclusive and psychologically safe work environment. We aim to remove discrimination, biases, and barriers to enable opportunities, career growth, and personal development.

Tecan maintains strong governance and ethical standards, as set out in the [Code of Conduct](#) and [supporting policies](#).

06. Governance




07. Governance

Executive compensation is linked to program achievements: The variable pay of Tecan's senior management is linked to Tecan's sustainability targets, as described in the company's [Governance Report](#).

Tecan's [Human Rights and Responsible Business Practices](#) policy sets out the company's commitments to stakeholders throughout its value chain.

08. Responsible Sourcing




09. Sustainability Reporting

Tecan is committed to transparency and reports on the above topics and more in its [Annual Report](#), as well as through a number of other platforms. Selected ratings are in the table below.

Survey	Current Rating
CDP - Climate Change	B
CDP - Water	C
CDP - SEA	A-
EcoVadis	Silver, 75/100
S&P CSA	54/100

10. Sustainability Ratings

