

Männedorf, 29 June 2020

Celebrating 40 years of innovation – Committed to diversity

This year we are celebrating Tecan's 40th anniversary. Since the beginning and to this very day, it is the passion and expertise of our employees that has made us the company we are – an innovating pioneer and global market leader in laboratory automation.

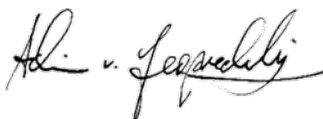
In 1980, it all started in a barn of a farm house on the banks of Lake Zurich in Switzerland. Today, we are a global company with manufacturing, research and development sites on three continents, a sales and service network spanning around the world and our exceptional colleagues from 50 different countries. We are proud of and value the diversity that has been built within Tecan as we approach 2,000 employees globally.

On the occasion of our 40th anniversary, we would therefore like to thank all former and current employees who made Tecan the company it is today. It is our quest to continuously seek new ways to empower both our customers and our colleagues to excel every day. We celebrate an open, diverse, compassionate and inviting culture where mutual respect, dignity and equality is a mindset and unwavering commitment to all of our colleagues and customers.

In our Tecan culture, we value diversity of ethnicity, origin, gender, religion or personal ideology, disability, age and/or sexual orientation. Further, we commit to continuously improving our diversity representation at all levels in our organization through the recruitment and development of our diverse colleagues. We see inclusion, diversity and equality as the foundation for our success and future progress since we draw energy and creativity from this multiplicity of experiences and viewpoints.

Our purpose is to positively impact healthcare and the life of all people through bridging the gap between life science research and clinical diagnostics. We believe the value we generate from our diverse perspectives will enable our success in delivering on our promise, We are always there for you.

Sincerely,



Achim von Leoprechting
CEO

