

Press Release

Tecan to present at upcoming Investor Conferences

Männedorf, Switzerland, January 8, 2024 - The Tecan Group (SIX Swiss Exchange: TECN) today announced that it will attend and present at the following investor conferences this week:

42nd Annual J.P. Morgan Healthcare Conference, San Francisco

Date: 10 January 2024 Presentation: 09:00 am PST (6:00 pm CET) Presenter: Dr. Achim von Leoprechting, CEO, Tecan Additional participant: Tania Micki, CFO Webcast live under the "Investor Relations" tab of Tecan's website www.tecan.com A replay of the webcast will be available after the presentation and will remain available for 30 days.

Baader Helvea Swiss Equities Conference, Bad Ragaz

Date: 11 January 2024 Participants: Klaus Lun, EVP, Head of the Life Sciences Business division Ralf Griebel, EVP, Head of the Partnering Business division Martin Brändle, SVP, Corp. Comm. & Investor Relations Not webcasted

Octavian Seminar 2024, Davos

Date: 12 January 2024 Participants: Klaus Lun, EVP, Head of the Life Sciences Business division Ralf Griebel, EVP, Head of the Partnering Business division Martin Brändle, SVP, Corp. Comm. & Investor Relations

Not webcasted

About Tecan

Tecan (www.tecan.com) improves people's lives and health by empowering customers to scale healthcare innovation globally from life science to the clinic. Tecan is a pioneer and global leader in laboratory automation. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments, components and medical devices that are then distributed by partner companies. Founded in Switzerland in 1980, the company has more than 3,500 employees, with manufacturing, research and development sites in Europe, North America and Asia, and maintains a sales and service network in over 70 countries. In 2022, Tecan generated sales of CHF 1,144 million (USD 1,192 million; EUR 1,144 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).

Tecan Group Ltd., Seestrasse 103, CH-8708 Männedorf, Switzerland T +41 44 922 88 88, F +41 44 922 88 89, info@tecan.com, www.tecan.com



2

For further information:

Tecan Group Martin Braendle Senior Vice President, Corporate Communications & IR Tel. +41 (0) 44 922 84 30 Fax +41 (0) 44 922 88 89 investor@tecan.com www.tecan.com