

Tecan schedules conference call on August 17th to discuss half year 2022 financial results

Männedorf, Switzerland, August 10, 2022 – The Tecan Group AG (SIX Swiss Exchange: TECN) will hold a conference call to discuss the financial results for the first half year 2022 on Wednesday, August 17th, 2022 at 09:00 CEST. The results will be published on August 17th, 2022 at 6:30 CEST.

The conference call will be audio-webcast with synchronized presentation slides. The webcast and a replay of the webcast will be made available in the "Investor Relations" section of Tecan's website www.tecan.com.

Interested parties can also listen to the conference call by phone.

The dial-in numbers for the call are as follows:

Participants from Europe: +41 (0) 58 310 50 00 or +44 (0) 207 107 0613 (UK)

Participants from the U.S.: +1 (1) 631 570 56 13

Participants should if possible dial in 15 minutes before the start of the event.

On August 17th, the full 2022 Interim Report will be available on the company website www.tecan.com under Investor Relations.

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 3,000 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 2021, Tecan generated sales of CHF 947 million (USD 1,041 million; EUR 877 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).

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