

Tecan to present at UBS Global Healthcare Virtual Conference

Männedorf, Switzerland, May 21, 2021 – The Tecan Group (SIX Swiss Exchange: TECN) today announced that it will present at the UBS Global Healthcare Virtual Conference next week.

UBS Global Healthcare Virtual Conference

Date: 25 May 2021

Presentation and Fireside Chat: 4:00 pm CET (10:00 am EST) with Dr. Achim von Leoprechting, CEO, Tecan

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.

About Tecan

Tecan (www.tecan.com) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics and clinical diagnostics. The company specializes in the development, production and distribution of automation solutions for laboratories in the life sciences sector. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries. In 2020, Tecan generated sales of CHF 731 million (USD 778 million; EUR 683 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