

Tecan to present at upcoming Investor Conferences

Männedorf, Switzerland, November 13, 2019 – The Tecan Group (SIX Swiss Exchange: TECN) today announced that it will present at the following investor conferences in the coming weeks:

CS Swiss EQ Mid Cap Conference, Zurich, Switzerland

Date: 15 November 2019, 10:30 am CET

Presenter: Dr. Achim von Leoprechting, CEO, Tecan

Not webcasted

Jefferies London Healthcare Conference, London, UK

Date: 20 November 2019, 1:20 pm GMT (2:20 pm CET)

Presenter: 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 90 days.

Berenberg European Conference 2019, Pennyhill Park, UK

Date: 3 and 4 December 2019

Presentation: 3 December 2019, 4:00 pm GMT (5:00 pm CET)

Presenter: Dr. Achim von Leoprechting, CEO, Tecan

Not webcasted

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 automated workflow 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 2018, Tecan generated sales of CHF 594 million (USD 606 million; EUR 516 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).



2

For further information:

Tecan Group

Martin Braendle

Vice President, Communications & Investor Relations

Tel. +41 (0) 44 922 84 30

Fax +41 (0) 44 922 88 89

investor@tecan.com

www.tecan.com