

Press Release

Tecan schedules Analyst and Media Conference on March 11 to discuss Full-Year 2013 Financial Results

Männedorf, Switzerland, March 3, 2014 – The Tecan Group AG (SIX Swiss Exchange: TECN) will hold an Analyst and Media Conference to discuss the full-year 2013 financial results on Tuesday, March 11, 2014 at 09:30 CET in Zurich. The results will be published on March 11, 2014 at 6:30 CET.

The conference will be audio-webcast with synchronized presentation slides. Interested parties may access the webcast via Tecan's homepage www.tecan.com. A link to the webcast will be provided immediately prior to the event.

Interested parties can also listen to the conference by phone.

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

Participants from Europe: +41 58 310 5000 or +44 203 059 58 62 (UK)

Participants from the U.S.: +1 866 291 4166

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

The recorded webcast will be available the same day on the website www.tecan.com together with the press release, the presentation and a preprint of the Financial Report 2013. An iPad app for Financial Reports of the Tecan Group is also available from the App Store.

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 2012, Tecan generated sales of CHF 391 million (USD 416 million; EUR 323 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).



2

For further information:

Tecan Group

Martin Braendle
Head of Corporate Communications & Investor Relations
Tel. +41 (0) 44 922 84 30
Fax +41 (0) 44 922 88 89
investor@tecan.com
www.tecan.com