

Tecan schedules conference call on August 14th to discuss half year 2013 financial results

Männedorf, Switzerland, August 7, 2013 – The Tecan Group AG (SIX Swiss Exchange: TECN) will hold a conference call to discuss the financial results for the first half year 2013 on Wednesday, August 14th, 2013 at 10:00 CET. The results will be published on August 14th, 2013 at 7:00 CET.

The conference call 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 call 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 (1) 866 291 4166

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

On August 14th, the full 2013 Interim Report will be available on the company website www.tecan.com under Investor Relations. 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).

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