

## **Tecan to present at J.P. Morgan Healthcare Conference**

**Männedorf, Switzerland, January 9, 2014** – The Tecan Group (SIX Swiss Exchange: TECN) today announced that it will present at the J.P. Morgan Healthcare Conference on January 16, 2014 in San Francisco, CA.

The investor presentation by David Martyr, CEO of Tecan, will begin at 1:30 p.m. Pacific time (10:30 p.m. CET) and will be webcast live. The webcast may be accessed under the “Investor Relations” tab of Tecan’s website [www.tecan.com](http://www.tecan.com).

A replay of the webcast will be available after the presentation and will remain available for 90 days.

### **About Tecan**

Tecan ([www.tecan.com](http://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](mailto:investor@tecan.com)

[www.tecan.com](http://www.tecan.com)