

## **Tecan announces Capital Markets Day**

**Maennedorf, Switzerland, June 9, 2016** - Tecan Group (SIX Swiss Exchange: TECN) will hold its Capital Markets Day on Thursday, June 16, 2016 close to its headquarters in Maennedorf/Zurich, Switzerland. The event is open for analysts, institutional investors and media representatives and will be hosted by CEO, Dr. David Martyr, along with other members of the management, who will provide an update on Tecan's strategy, priorities and new products in the Life Sciences Business as well as Partnering Business divisions.

Presentations will begin at 08:45 a.m. CET at an event location and will be webcast live. All presentations will be in English. The focus in the afternoon will be on product demonstrations and a site tour at the Tecan headquarters in Maennedorf.

The webcast and a replay of the webcast will be made available in the "Investor Relations" section of Tecan's website [www.tecan.com](http://www.tecan.com).

### **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 2015, Tecan generated sales of CHF 440 million (USD 459 million; EUR 411 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TECN; ISIN CH0012100191).

### **For further information:**

#### **Tecan Group**

Martin Brändle  
Vice President, 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)