

## **Tecan appoints Dr. Achim von Leoprechting to the Management Board**

**Männedorf, Switzerland, August 9, 2013** – The Tecan Group (SIX Swiss Exchange: TECN) today announced that Dr. Achim von Leoprechting has been appointed a member of the Management Board. As the new Head of the Partnering Business division, effective October 1, 2013, he will assume responsibility for Tecan's global OEM business, which has been led by Dr. David Martyr, CEO of the Tecan Group, on an interim basis since the beginning of the year.

Dr. David Martyr commented: "I am very pleased that Achim von Leoprechting will be strengthening our management team with his long-standing experience in the life science sector. This is a further important step in the process of reshaping our management team that I began at the beginning of the year. We now have the right divisional leadership in place for our Partnering Business and Life Sciences Business, with extensive domain knowledge, experience and capabilities to support our future accelerated growth."

Dr. Achim von Leoprechting has 14 years of experience in a wide range of management positions in the life science industry. Since 2002, he has worked in various areas and held diverse positions with increasing management responsibility at PerkinElmer Inc. (NYSE:PKI), most recently as Vice President and General Manager Europe, Middle East, Africa and India. In this role, Achim von Leoprechting was responsible for all Life Science and Technology business conducted by the U.S. group in these regions worldwide. Prior to this, he held positions with global responsibility for individual business divisions within the group of companies, for example as Vice President and General Manager In Vitro Solutions and Vice President and General Manager Imaging and Detection Technologies. In addition, as European Sales Director Molecular Medicine / Bio-discovery he was responsible for a regional sales organization for several years. Before working for PerkinElmer, Achim von Leoprechting held a number of positions in product management at Packard Bioscience, which today is also a part of PerkinElmer.

Achim von Leoprechting has a doctorate in biology from the University of Freiburg (Germany). He is 45 years old and a German national.

### **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



2

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 Brändle

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)