

Annual General Meeting 2017: Tecan shareholders endorse all motions proposed by the Board of Directors

- **Dividend of CHF 1.75 per share**
- **Rolf Classon, Heinrich Fischer, Oliver Fetzer, Lars Holmqvist, Karen Huebscher, Christa Kreuzburg and Gérard Vaillant re-elected as members of the Board of Directors**
- **Rolf Classon confirmed by shareholders as Chairman of the Board**
- **Christa Kreuzburg, Oliver Fetzer and Gérard Vaillant confirmed as members of the Compensation Committee**
- **All proposals relating to compensation approved**

Männedorf, Switzerland, April 11, 2017 – The Annual General Meeting of Tecan Group Ltd. (SIX Swiss Exchange: TECN) has today endorsed all motions proposed by the Board of Directors. Shareholders approved the Annual Report, the financial statements and the consolidated financial statements, and granted discharge to the Board of Directors and the Management Board. They also approved a dividend of CHF 1.75 per registered share, unchanged versus the previous year. The payout will take place on April 19, 2017.

All members of the Tecan Board of Directors were confirmed by the shareholders for a term of one year: Heinrich Fischer, Dr. Oliver Fetzer, Lars Holmqvist, Dr. Karen Huebscher, Dr. Christa Kreuzburg, Gérard Vaillant and Rolf Classon. Rolf Classon was re-elected as Chairman of the Board of Directors. The shareholders also confirmed Christa Kreuzburg, Oliver Fetzer and Gérard Vaillant as members of the Compensation Committee.

Ernst & Young AG, Zurich, was reappointed as the statutory auditors for fiscal year 2017. Shareholders approved the motion proposed by the Board of Directors to appoint Proxy Voting Services GmbH, Zurich, as the independent proxy for the period up to the close of the Tecan Group 32nd Annual General Meeting in 2018.

All proposals relating to compensation approved

Shareholders voted on the conditions governing the compensation of the Board of Directors and Management Board for the following fiscal year, according to the Ordinance Against Excessive Compensation in Listed Stock Companies of the Swiss government. The shareholders made a binding decision to endorse the proposed motion on the maximum total amount of compensation for 2018. The compensation report, which was submitted for an advisory vote for the sixth time, was also accepted by shareholders.



2

Key upcoming date

Tecan will publish the 2017 Interim Report on August 16, 2017.

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 2016, Tecan generated sales of CHF 506 million (USD 511 million; EUR 464 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

www.tecan.com