

Annual General Meeting 2011: Tecan shareholders endorse all motions proposed by the Board of Directors

- **Removal of the registration and voting limitations to further strengthen good Corporate Governance**
- **Oliver S. Fetzter and Erik Walldén elected as new members of the Board of Directors**
- **Rolf A. Classon, Heinrich Fischer, Lukas Braunschweiler, Dominique F. Baly and Gérard Vaillant re-elected for a term of one year; Rolf A. Classon confirmed as Chairman**
- **Dividend unchanged at CHF 1.00 per share**

Männedorf, Switzerland, April 19, 2011 – The Annual General Meeting of the Tecan Group (SIX Swiss Exchange: TECN) has today endorsed all motions proposed by the Board of Directors. The 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 an unchanged dividend on the previous year of CHF 1.00 per registered share. The dividend will be paid out from the available reserve for additional paid-in capital and is therefore not subject to withholding tax. The payout will take place on April 28, 2011.

Rolf A. Classon, Heinrich Fischer, Lukas Braunschweiler, Dominique F. Baly and Gérard Vaillant were re-elected as Members of the Board of Directors of Tecan for a term of one year. The shareholders elected Oliver S. Fetzter and Erik Walldén as new members of the Board. Rolf A. Classon, Chairman of the Board of Directors, bid farewell to members Jürg Meier and Peter Ryser, who are leaving, thanking them warmly for their commitment and valuable contributions to Tecan's development. Immediately after the Annual General Meeting, Rolf A. Classon was confirmed as Chairman by the Board.

Further improvements in Tecan's Corporate Governance

The shareholders voted in favor of removing the restrictions of registration and voting rights to 5% of the share capital. The Board of Directors of Tecan had proposed the corresponding amendment to the Articles of Incorporation in order to reinforce shareholders' participation rights, thereby further improving Corporate Governance. This shareholder-friendly measure also makes the Tecan share more attractive, particularly in an international environment and with institutional investors. The shareholders also approved the change proposed by the Board of Directors to the Articles of Incorporation to lift the ban on entries in the share register after invitations to the Annual General Meeting have been sent. This measure also creates better conditions for shareholders to exercise their rights.



2

KPMG AG, Zurich, was confirmed as statutory auditor for 2011.

Next date for the diary

The 2011 Interim Report will be published on Thursday, August 18, 2011.

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 2010, Tecan generated sales of CHF 371 million (USD 356 million; EUR 269 million). Registered shares of Tecan Group are traded on the SIX Swiss Exchange (TK: TECN/Reuters: TECZn.S/ ISIN: 12100191).

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

www.tecan.com