

Changes to the Tecan Development & Operations Division

- **Martin Brusdeilins will take over as Head of the Development & Operations Division with immediate effect**
- **Jürg Dübendorfer, up to now Head of the Development & Operations Division, is leaving Tecan at his own request by no later than the end of May 2012**

Männedorf, Switzerland, November 8, 2011 – The Tecan Group (SIX Swiss Exchange: TECN) announced today that Martin Brusdeilins will take over as Head of the Development & Operations Division with immediate effect. Jürg Dübendorfer, who has led the Division since January 2011, is to leave the company and the Group Management Board at his own request by no later than the end of May 2012.

Martin Brusdeilins has been a Member of the Group Management Board and Head of the Science & Technology organizational unit since 1 September 2011. He will continue in this role in addition to the new position. Martin Brusdeilins has extensive experience in a variety of management positions in the diagnostic and life science industry. Before joining Tecan, he worked at Ortho-Clinical Diagnostics Inc., New York, USA, for almost 8 years where he was responsible for global product development of all diagnostic instruments and reagents as well as the life cycle management of these products.

Jürg Dübendorfer joined Tecan in 2000 and has held various technical and management positions. As a Member of the Group Management Board, he led the newly created Services & Consumables business unit from 2006 until the end of 2010. Under his leadership, the business unit recorded significant growth rates with enhanced profitability.

Tecan's Board of Directors and CEO regret Mr Dübendorfer's decision. They would like to thank him warmly for his valuable contribution to the development of Tecan and wish him all the best for the future.

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



2

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