

Tecan successfully completes the acquisition of Sias

Männedorf, Switzerland, November 30, 2015 – The Tecan Group (SIX Swiss Exchange: TECN) announced today that it has successfully completed the acquisition of Sias to further expand its Partnering Business. On October 27, 2015, Tecan announced that it had reached agreement with the majority shareholders of Sias to acquire Sias AG, a leading OEM supplier of a wide range of modular and complete laboratory automation solutions. From December 1, 2015, Sias will be included in the consolidated financial statements of the Tecan Group as a part of the Partnering Business segment.

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 2014, Tecan generated sales of CHF 400 million (USD 437 million; EUR 331 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