

Invitation.

Analyst and Media Conference

Full-Year Financial Results 2018
Tecan Group AG
Zurich, 11 March 2019

Ladies and Gentlemen,

We cordially invite you to Tecan's 2018 Full-Year Financial Results Conference.

Date: **Wednesday, March 11, 2019**

Time: **9:00 am—10:00 am (CET)**

Venue: **Widder Hotel
Rennweg 7
8001 Zurich**

Speakers: **David Martyr, CEO
Achim von Leoprechting, EVP &
designated CEO
Rudolf Eugster, CFO**

The press release with the 2018 full-year financial results and a preprint of the Financial Report 2018 will be published on Monday, March 11, 2019 at 6:30 am (CET).

The conference will be audio-webcast with synchronized presentation slides. Interested parties may access the webcast via Tecan's homepage www.tecan.com. A link to the webcast together with the presentation will be provided immediately prior to the event.

Interested parties can also listen to the conference by phone. Please dial the following numbers 15 minutes before the meeting starts:

Europe: 9:00 am (CET) +41 58 310 50 00

UK: 8:00 am (GMT) +44 207 107 0613

USA: 3:00 am (EST) +1 631 570 5613

A replay of the webcast will be available the same day on our website www.tecan.com.

For organizational purposes, we kindly ask you to return the enclosed reply form before March 1, 2019.

We look forward to welcoming you in person or online.

Kind regards



Martin Braendle
VP, Communications & Investor Relations
T +41 44 922 8430, F +41 44 922 8889
investor@tecan.com
www.tecan.com

TECAN.



