



# Invitation.

## Analyst and Media Conference

**Full-Year Financial Results 2017**  
**Tecan Group AG**  
**Zurich, 14 March 2018**

Ladies and Gentlemen,

**We cordially invite you to Tecan's 2017  
Full-Year Financial Results Conference.**

Date: **Wednesday, March 14, 2018**  
Time: **9:00 am—10:00 am (CET)**  
Venue: **SIX Swiss Exchange  
Pfungstweidstrasse 110  
8005 Zurich**  
Room: **Auditorium**  
Speakers: **David Martyr, CEO  
Rudolf Eugster, CFO**

The press release with the 2017 full-year financial results and a preprint of the Financial Report 2017 will be published on Wednesday, March 14, 2018 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](http://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](http://www.tecan.com).

**For organizational purposes, we kindly ask you to return the enclosed reply form before March 1, 2018.**

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](mailto:investor@tecan.com)  
[www.tecan.com](http://www.tecan.com)



---

# Reply form.

## Analyst and Media Conference

---

Full-Year Financial Results 2017

Tecan Group AG

Zurich, 14 March 2018

---

Date & time:

**14 March 2018, 9:00 am**

Location:

**SIX Swiss Exchange Convention Point,  
Auditorium  
Pfingstweidstrasse 110, 8005 Zurich**

- Yes, I will attend the conference in Zurich
- I will participate in the webcast / conference call
- No, I am not able to participate

Please send your response by email or fax no later than  
1 March 2018 to:

Martin Braendle, [investor@tecan.com](mailto:investor@tecan.com) or  
Fax +41 (0)44 922 88 89

Last name:

---

First name:

---

Company:

---

Country:

---

Email:

---

Tecan Group Ltd.,  
Seestrasse 103,  
CH-8708 Männedorf,  
Switzerland  
T +41 44 922 88 88,  
F +41 44 922 88 89,  
[info@tecan.com](mailto:info@tecan.com),  
[www.tecan.com](http://www.tecan.com)

