

# Invitation.

## Analyst and Media Conference

Full-Year Financial Results 2019

Tecan Group AG

Zurich, 17 March 2020

Ladies and Gentlemen,

We cordially invite you to Tecan's 2019 Full-Year Financial Results Conference.

Date: **Tuesday, March 17, 2020**

Time: **9:00 am–10:00 am (CET)**

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

Speakers: **Achim von Leoprechting, CEO  
Rudolf Eugster, CFO (until 28.2.20)**

Other members of the management will also attend the event. The press release with the 2019 full-year financial results and the annual report will be published on Tuesday, March 17, 2020 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 5, 2020.**

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)

 **TECAN.**



---

# Reply form.

## Analyst and Media Conference

---

Full-Year Financial Results 2019

Tecan Group AG

Zurich, 17 March 2020

---

Date & time:

17 March 2020, 9:00 am

Location:

**Widder Hotel**  
**Rennweg 7**  
**8001 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  
5 March 2020 to:  
Martin Braendle, investor@tecan.com

Last name:

---

First name:

---

Company:

---

Country:

---

Email:

---

Tecan Group AG,  
Seestrasse 103,  
CH-8708 Männedorf,  
Switzerland  
T +41 44 922 88 88,  
F +41 44 922 88 89,  
info@tecan.com,  
www.tecan.com

The logo consists of a red circle with a white dot inside, followed by the word "TECAN." in a bold, black, sans-serif font.

