

Tecan Group Ltd, Männedorf

Buyback of own registered shares on the standard trading line on SIX Swiss Exchange

The Board of Directors of Tecan Group Ltd, Seestrasse 103, 8708 Männedorf («Tecan» or the «Company»), has decided on 11 August 2025 to buy back its own registered shares with a nominal value of CHF 0.10 each for a maximum acquisition value of CHF 120 million. The repurchased registered shares shall be used for general business purposes, including treasury purposes and the financing of potential acquisitions (the «buyback program»). Based on the closing price of the registered share on 8 August 2025 on the SIX Swiss Exchange, this corresponds to a maximum of 770,218 registered shares or a maximum of 6.01% of the share capital of the Company, which currently amounts to CHF 1,282,588.30 and is divided into 12,825,883 listed registered share with a par value of CHF 0.10 each. Due to future share price developments, the number of registered shares actually repurchased may deviate from the stated number of registered shares, however, under no circumstances will more than 10% of the share capital and voting rights entered in the Commercial Register be repurchased under the buyback program.

The buyback program is exempted from compliance with the provisions on public takeover bids provided for based on section 6.1 of Circular No. 1 of the Swiss Takeover Board dated 27 June 2013.

Duration of the buyback program

Shares may be purchased under the buyback program on the SIX Swiss Exchange between 13 August 2025 and 12 August 2027 inclusive. Tecan is under no obligation to repurchase registered shares at any time. Tecan will decide to make purchases based on the market conditions. Tecan reserves the right to terminate the buyback program early.

Delegation agreement

Tecan and Zürcher Kantonalbank have entered into a delegation agreement pursuant to Article 124 (2) a) and (3) FMIO. Under the agreement, Zürcher Kantonalbank can repurchase shares independently, subject to certain criteria. Tecan is however entitled to terminate the delegation agreement at any time without stating its reasons or modify the parameters in accordance with Article 124 (3) FMIO.

Maximum daily buyback amount

In accordance with Article 123 (1) c) FMIO, the maximum daily buyback amount is published on Tecan's website at: <https://www.tecan.com/sbp>

Publication of transaction details

Tecan will publish details of all transactions pursuant to Section 27 of the Swiss Takeover Board Circular No. 1 of 27 June 2013 on Tecan's website at: <https://www.tecan.com/sbp>

Non-public information

Tecan confirms that, as of the date of this advertisement, it does not have any non-public information which could have a considerable influence on the shareholders' ability to make a decision.

Own registered shares

As of 8 August 2025, Tecan held, directly or indirectly, 125,784 own registered shares (1.0% of the share capital and of the voting rights).

Shareholders with more than 3% of voting rights

According to the company's annual report and the disclosure notifications received until 8 August 2025, the following shareholders held more than 3% of the share capital and the voting rights of Tecan:

Chase Nominees Ltd., London, United Kingdom:
12.1% of the share capital and of the voting rights

UBS Fund Management (Switzerland) AG, Basel, Switzerland:
10.2% of the share capital and of the voting rights

BlackRock Inc., New York, NY, United States:
4.5% of the share capital and of the voting rights

Pictet Asset Management SA, Geneva, Switzerland:
5.01% of the share capital and of the voting rights

Israel Englander, New York, United States (direct/indirect holder:
Millennium Partners LP, George Town, KY Integrated Assets III LLC,
Wilmington, Delaware, United States):
3.003% of the share capital and of the voting rights

Credit Suisse Funds AG, Zurich, Switzerland:
3.0% of the share capital and of the voting rights

Swisscanto Fondsleitung AG, Zurich, Switzerland:
3.0% of the share capital and of the voting rights

Tecan has no knowledge whether the shareholders have any intentions to sell registered shares within the scope of the buyback program.

Applicable law and place of jurisdiction

Swiss law. The exclusive place of jurisdiction is Zurich.

Mandated bank

Tecan has mandated Zürcher Kantonalbank with the execution of the buyback program.

Swiss security number / ISIN / ticker symbol

Tecan Group Ltd registered share of CHF 0.10 nominal
1210019 / CH0012100191 / TECN

This notice does not constitute a prospectus for the purposes of Art. Art. 35 et seq. of FINSA.

This offer is not being and will not be made, directly or indirectly, in the United States of America and/or to US persons and may be accepted only by Non-US persons and outside the United States of America. Accordingly, copies of this document and any related materials are not being, and may not be, sent or otherwise distributed in or into or from the United States of America, and persons receiving any such documents (including custodians, nominees and trustees) may not distribute or send them in, into or from the United States of America.

12 August 2025