

SARTORIUS

Sartorius Aktiengesellschaft Göttingen

ISIN DE0007165607 and
ISIN DE0007165631

Dividend announcement

The Annual General Meeting of our Company held on March 26, 2026 resolved to appropriate the Company's retained profit totaling €338,851,995.67 for fiscal 2025 as follows:

Payment of a dividend of € 0.73 per dividend-bearing ordinary share	= €24,994,228.37
Payment of a dividend of € 0.74 per dividend-bearing preference share	= €25,762,079.54
Carried forward to new account	€288,095,687.76
Total:	€338,851,995.67

In accordance with Section 58(4)(2) of the German Stock Corporation Act (AktG), the entitlement to the dividend will be deemed payable on the third business day following the resolution of the Annual General Meeting, i.e. on April 1, 2025. The payment of the distribution amounts is generally made after deduction of 25% capital gains tax and 5.5% solidarity surcharge on the capital gains tax (26.375% in total) as well as church tax, if applicable, by the custodian banks.

The central paying agent is Commerzbank AG.

Domestic shareholders whose shares are held in a custody account at a domestic credit institution shall receive the dividend without deduction of capital gains tax and solidarity surcharge, provided that the credit institution is entitled to do so on the basis of a non-assessment certificate, a "permanent overpayer" certificate or an exemption instruction issued by the shareholder concerned.

In the case of foreign shareholders, the withheld capital gains tax including the solidarity surcharge may be reduced in accordance with European law or existing agreements for the avoidance of double taxation between the Federal Republic of Germany and the state concerned.

Göttingen, March 2025
The Executive Board