

Simplifying Progress

Diversity Fact Sheet

Sartorius Stedim Biotech 2023

SARTORIUS

“At Sartorius, diversity is not about statistics. It’s about culture. We foster diversity because it ultimately teaches us to be more open. And we consider openness as a key prerequisite for success.”

Joachim Kreuzburg, CEO



Welcome | 553 New People Joined Sartorius Stedim Biotech in 2023

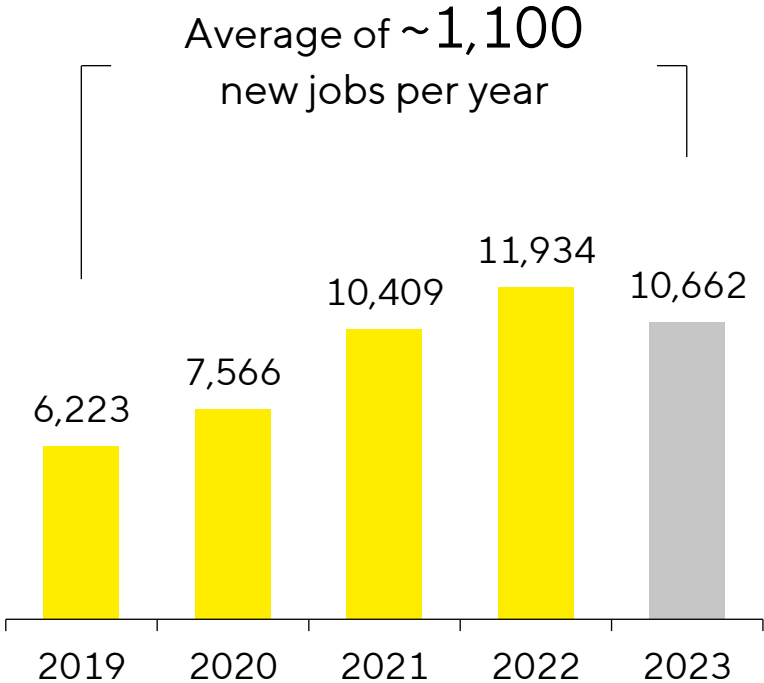
553 | New hires in 2023

264 | New employees from M&A in 2023

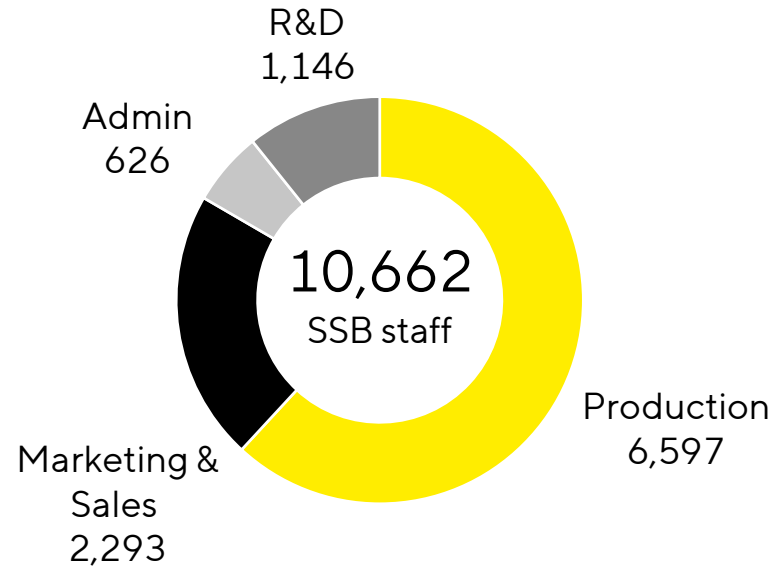
2,160 | Avg. hires annually 2019-2023

39% | Employees with higher education

Employee growth



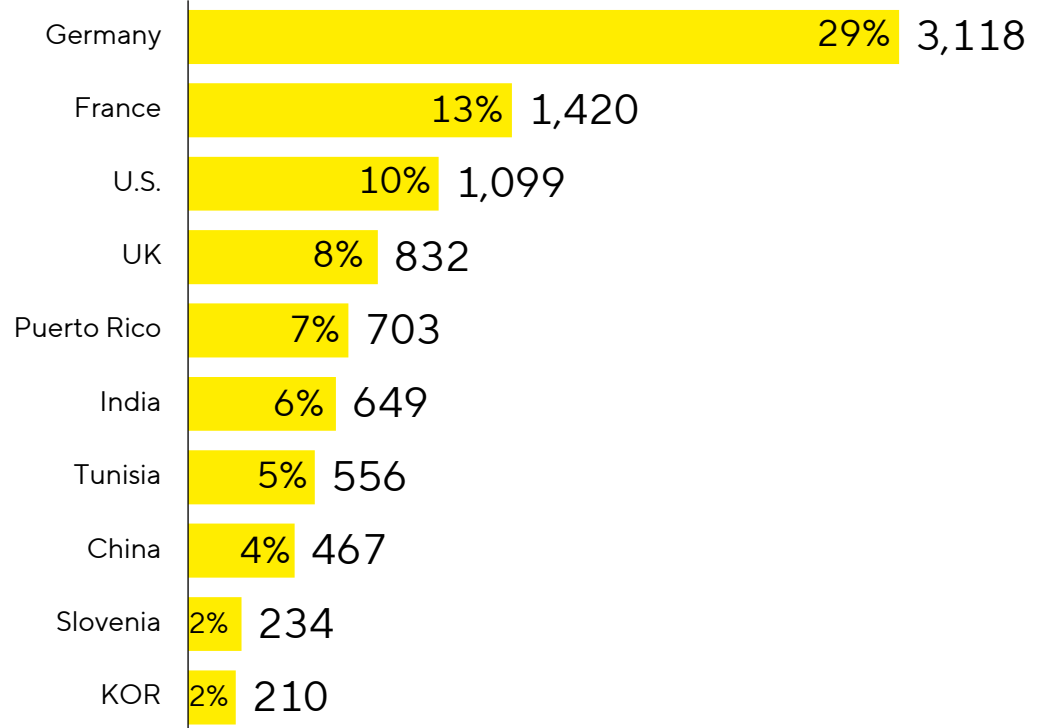
Employees by function



Nationality | SSB Unites People from 102 Nations

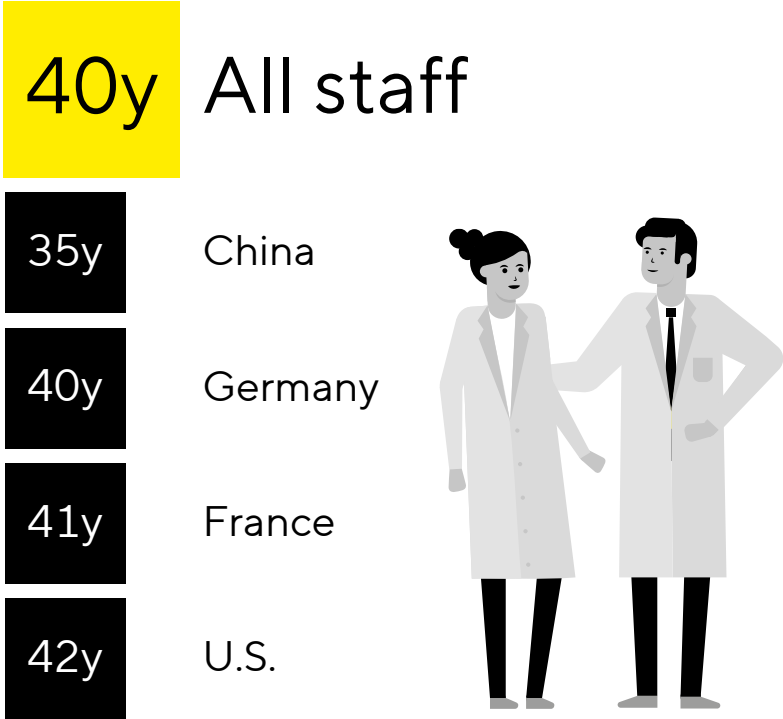


Top 10 nationalities – Share and number of staff

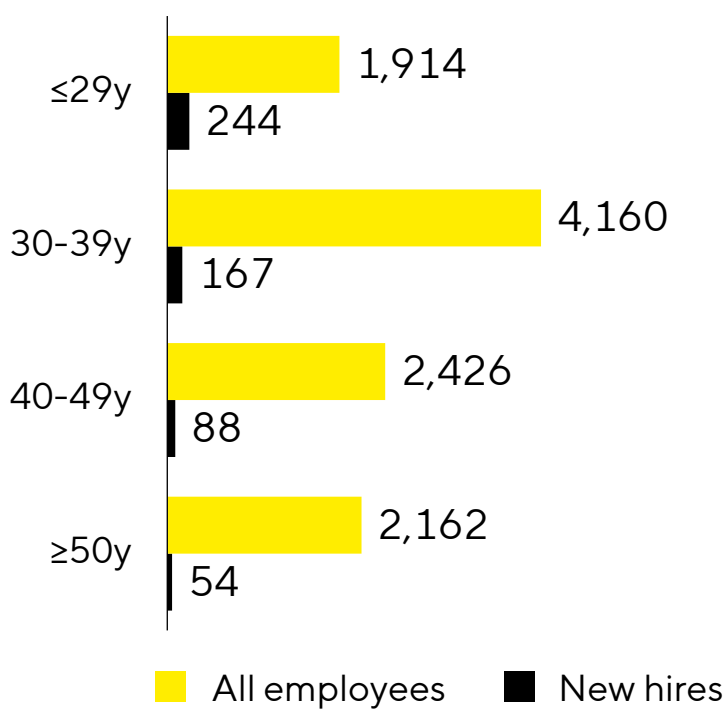


Age | Young Talent and Experienced Professionals Shape Our Workforce

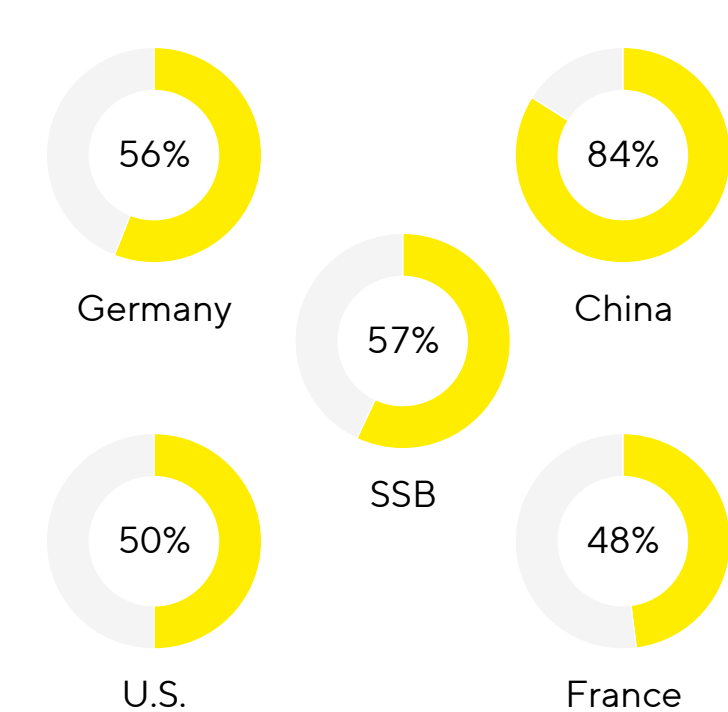
Average age



Number of staff by age

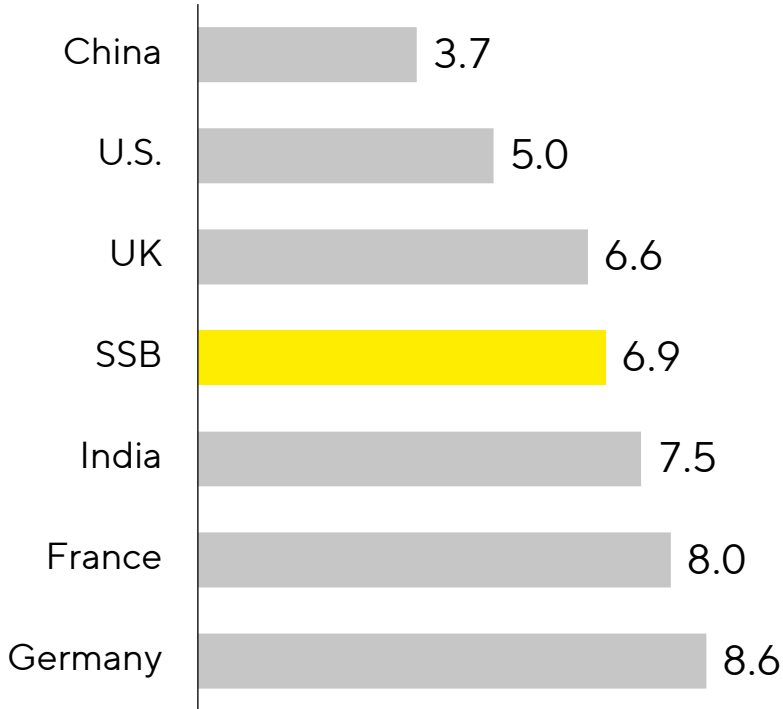


Share of staff younger than 40y

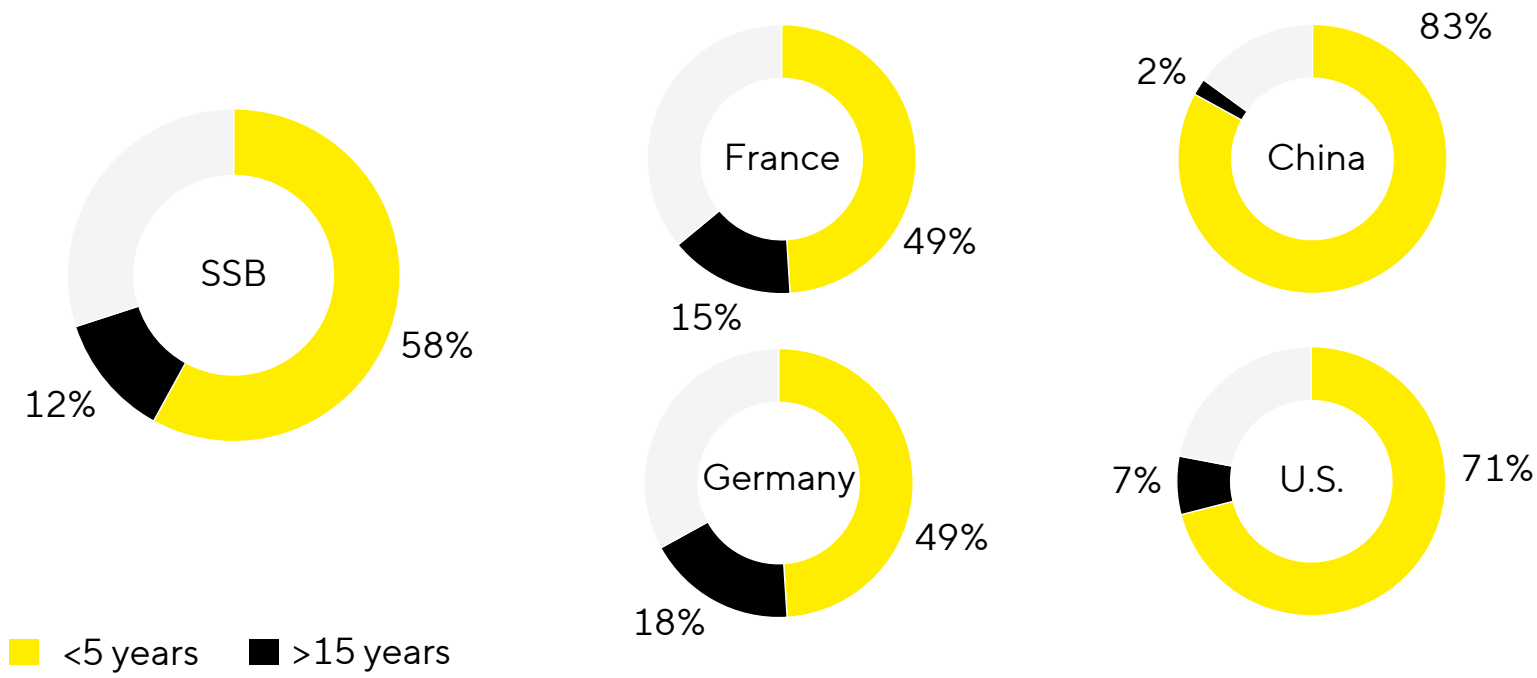


Tenure | Mixture of Experience and Fresh Ideas

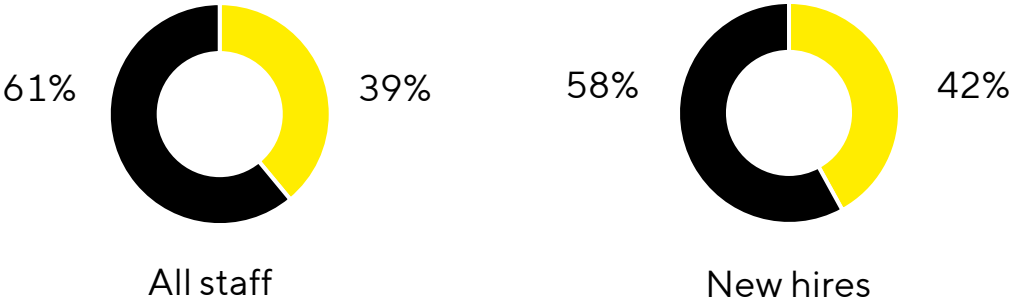
Average tenure in years



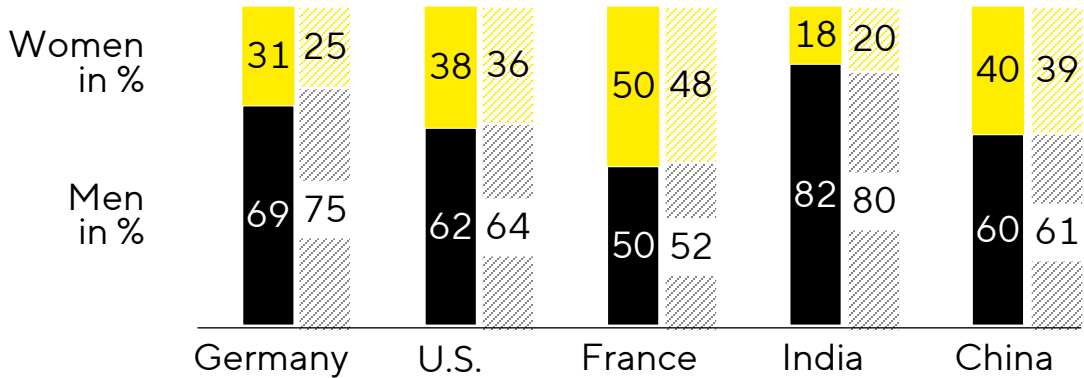
Share of staff with tenure higher than 15 years and lower than 5 years



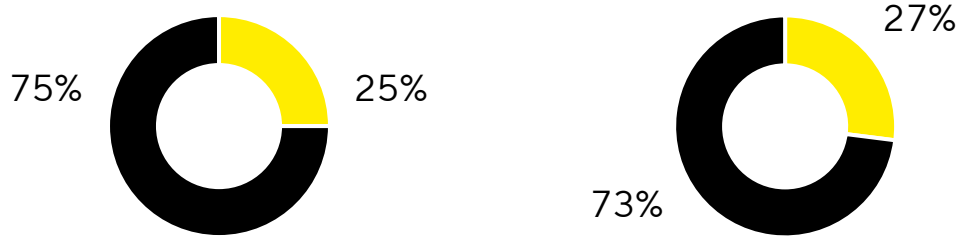
Gender | Structure Headed in the Right Direction



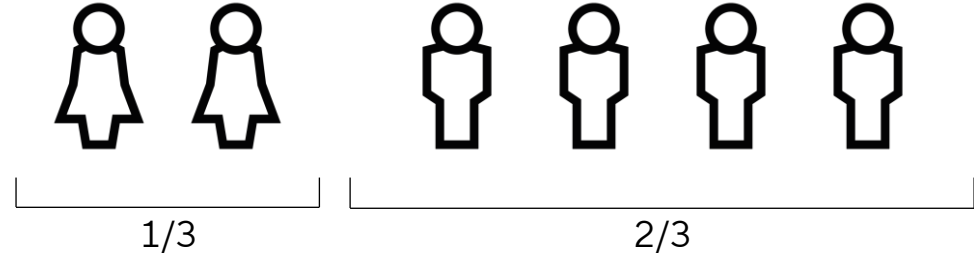
Women Men



Employees Management positions Women Men



Target 2025: One-third women in both management levels below the Executive Board



¹ As of April 1, 2024

Fostering Exchange and Dialogue for Women

Sartorius Business Women Association

The network serves to connect women across all regions and aims to support gender parity in leadership positions at Sartorius.

Partnership with global non-profit

Since 2021, Sartorius has been a corporate partner of the Healthcare Businesswomen Association, providing interested employees with access to the organizations network.



Promoting Equality and Raising Awareness of Diversity and Inclusion

North America Diversity, Equity, Inclusion and Belonging (DEIB)

The committee promotes DEIB in the U.S. through community outreach, partnerships, events, and training. So far, 800+ employees have completed Emtram’s Unconscious Bias Awareness training.

LGBT+Friends Alliance

The group strives to enhance the visibility of LGBTQIA+ individuals at Sartorius through virtual events and inclusive "Rainbow Lunches" – creating a safe space for sharing experiences.



Partnering in the Name of Diversity and Inclusion



Initiative zur Berufs- und Studienwahl