

Simplifying Progress

Diversity Fact Sheet

Sartorius Stedim Biotech 2024

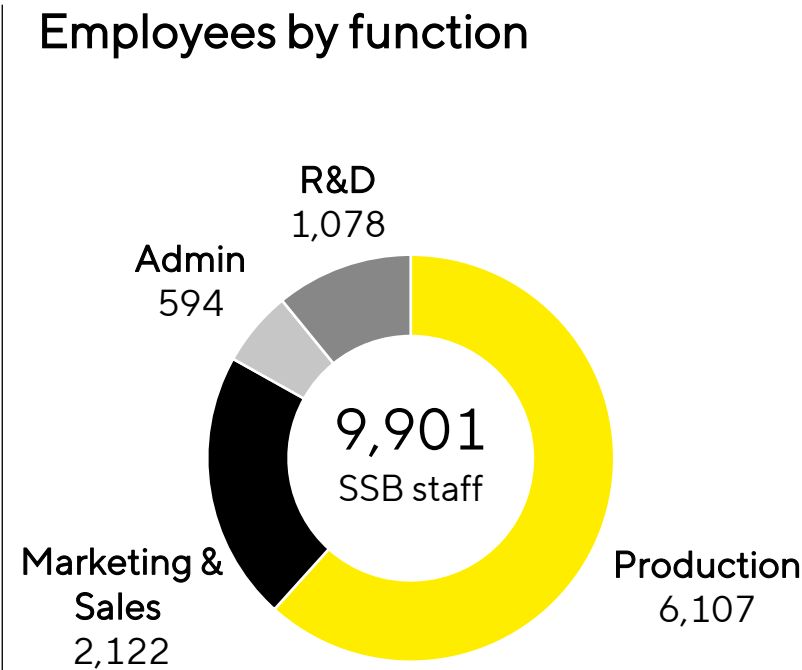
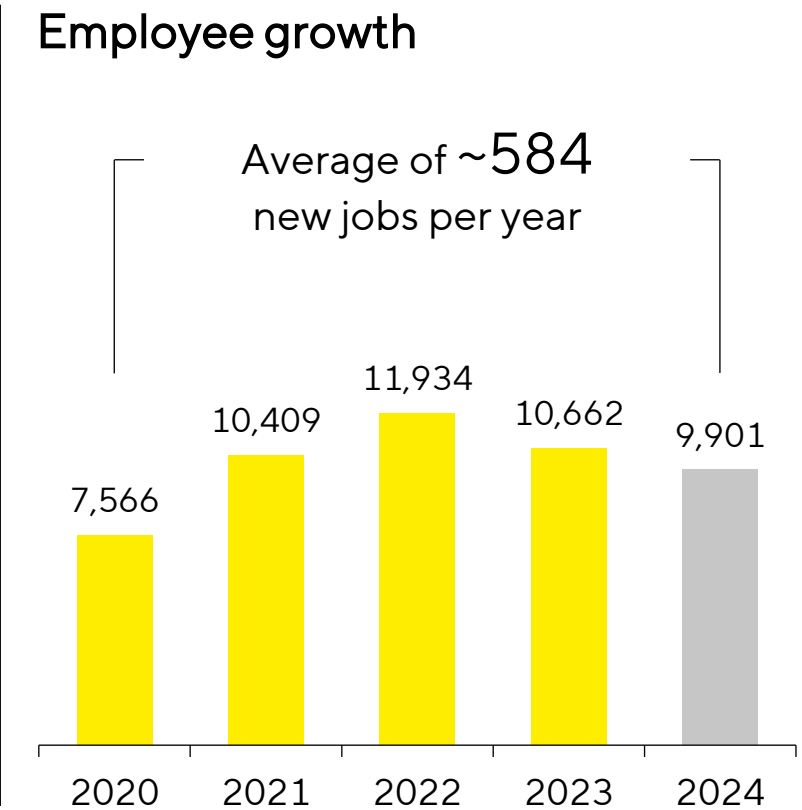
SARTORIUS

Welcome | 542 New People Joined Sartorius Stedim Biotech in 2024

542 | New hires in 2024

1,837 | Avg. hires annually 2020-2024

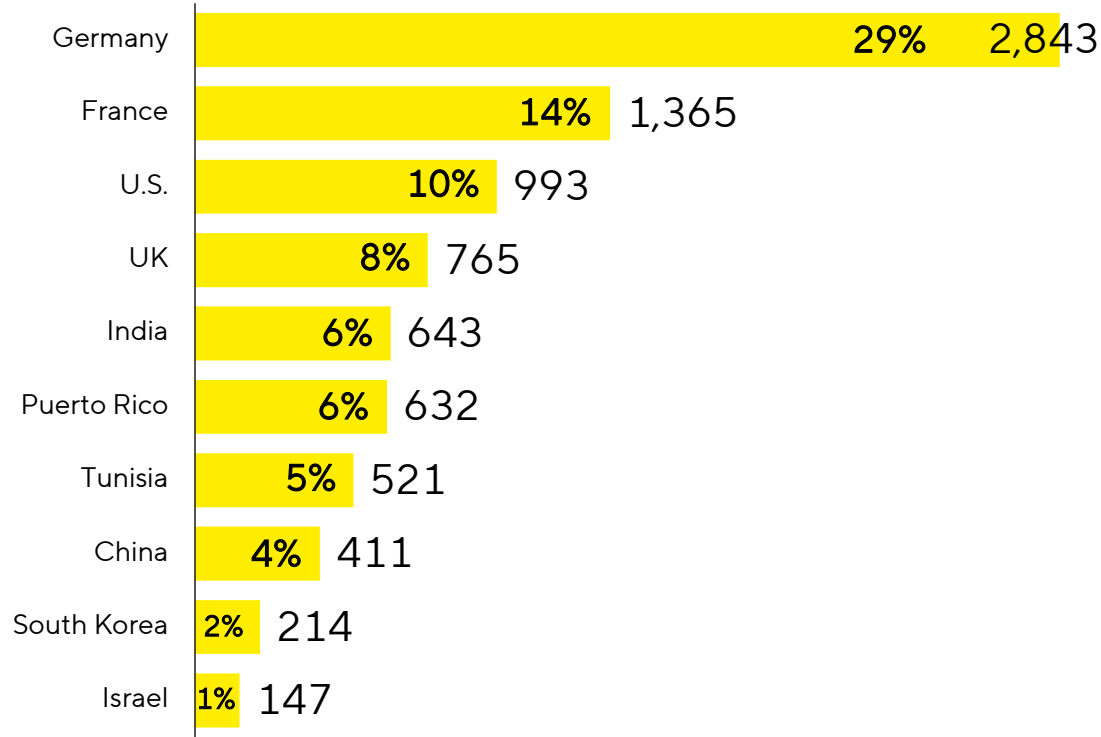
41% | Employees with higher education



Nationality | SSB Unites People from 99 Nations

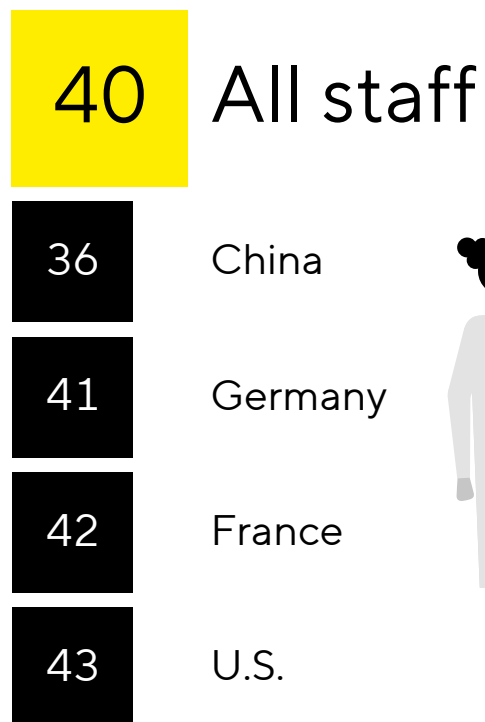


Top 10 nationalities – Share and number of staff

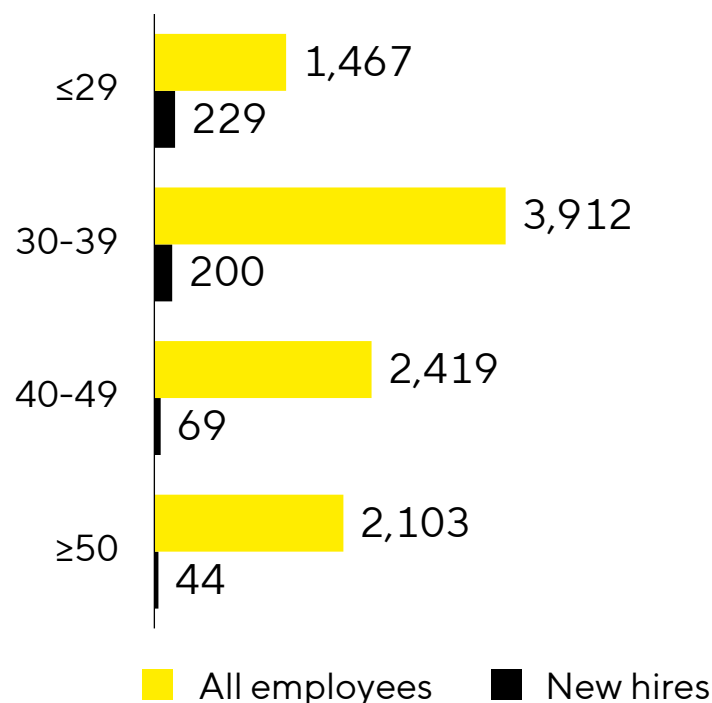


Age | Young Talent and Experienced Professionals Shape Our Workforce

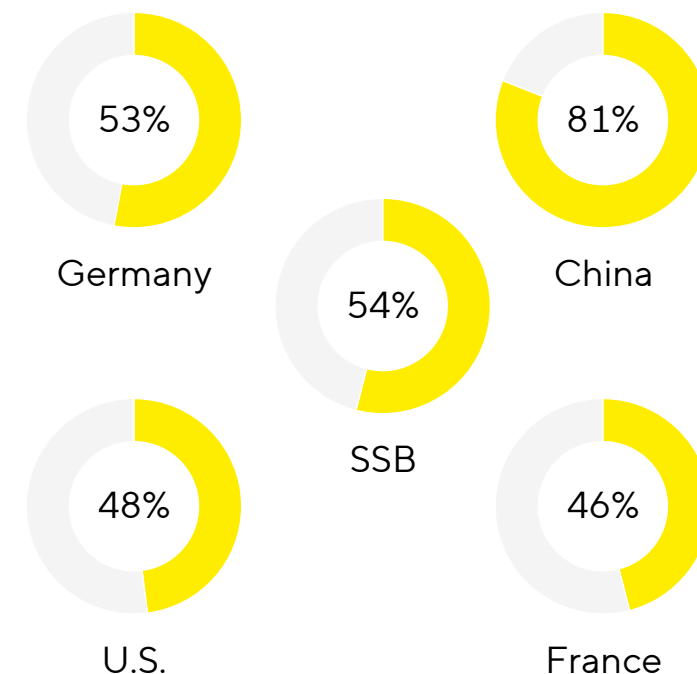
Average age



Number of staff by age

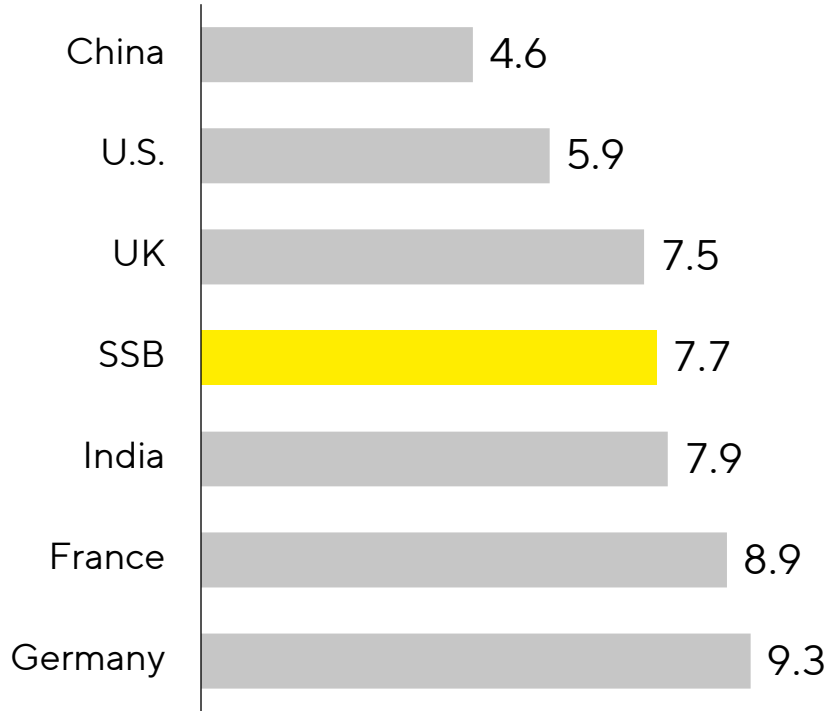


Share of staff younger than 40y

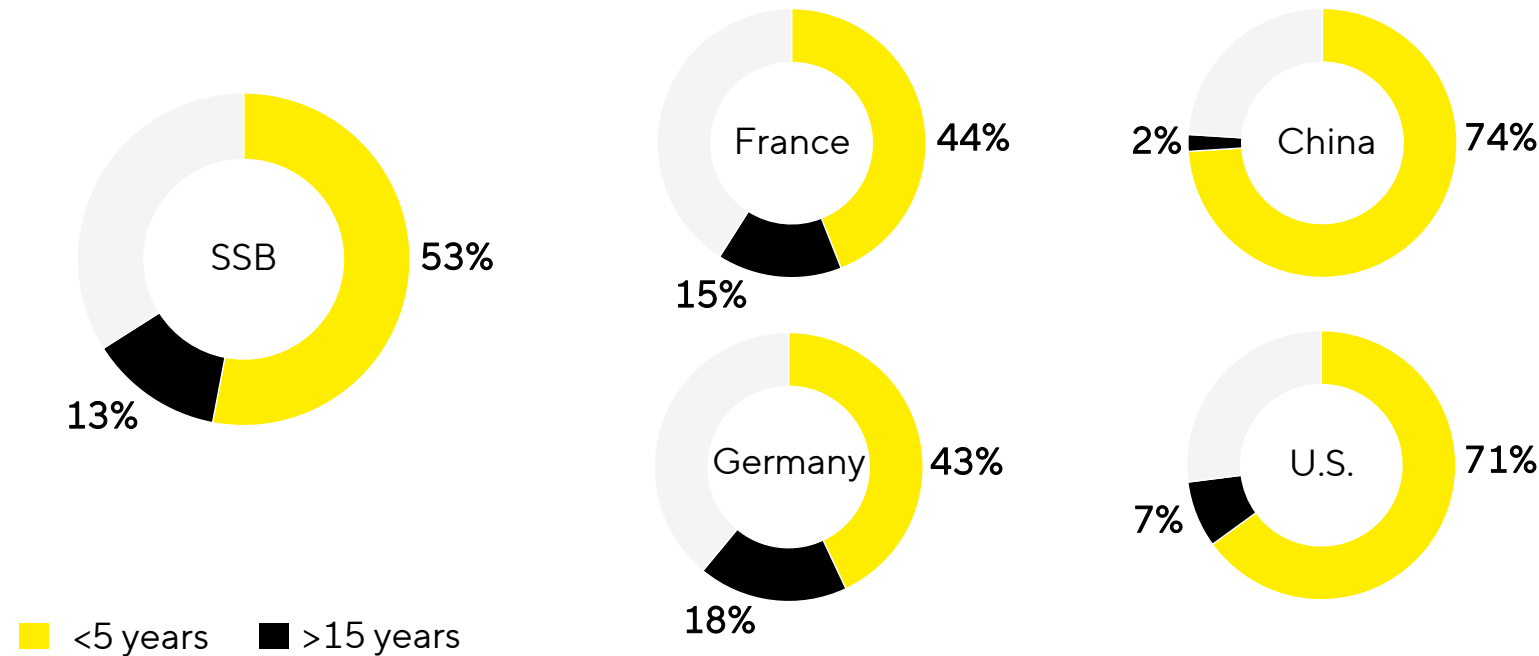


Tenure | Mixture of Experience and Fresh Ideas

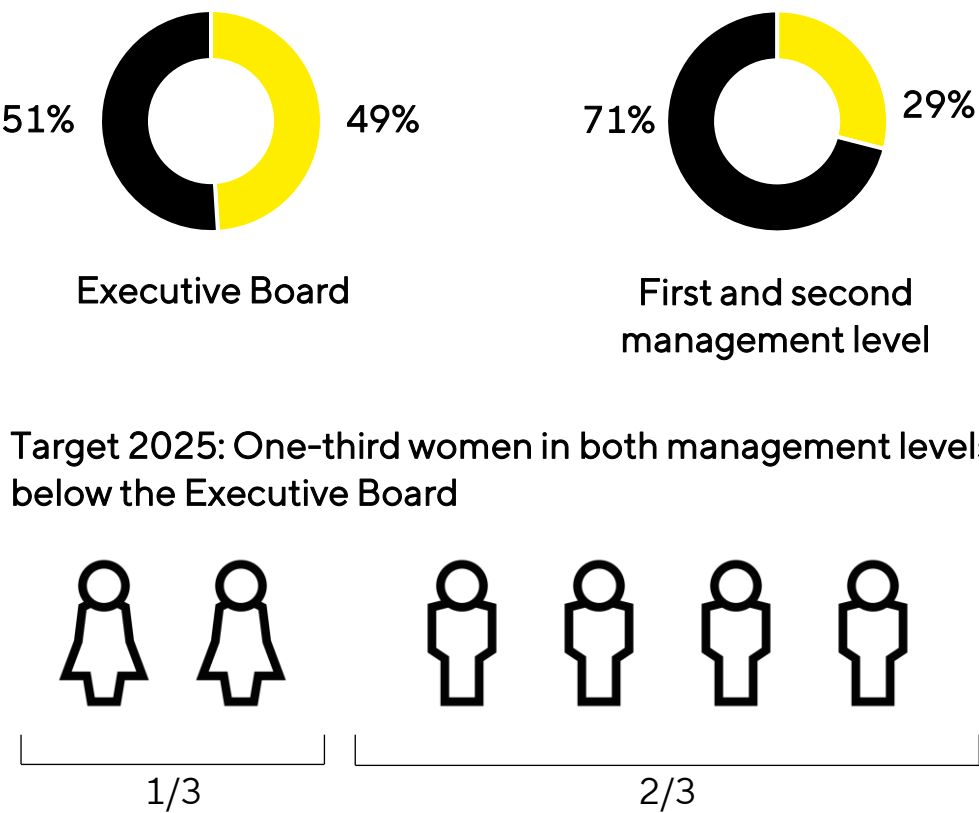
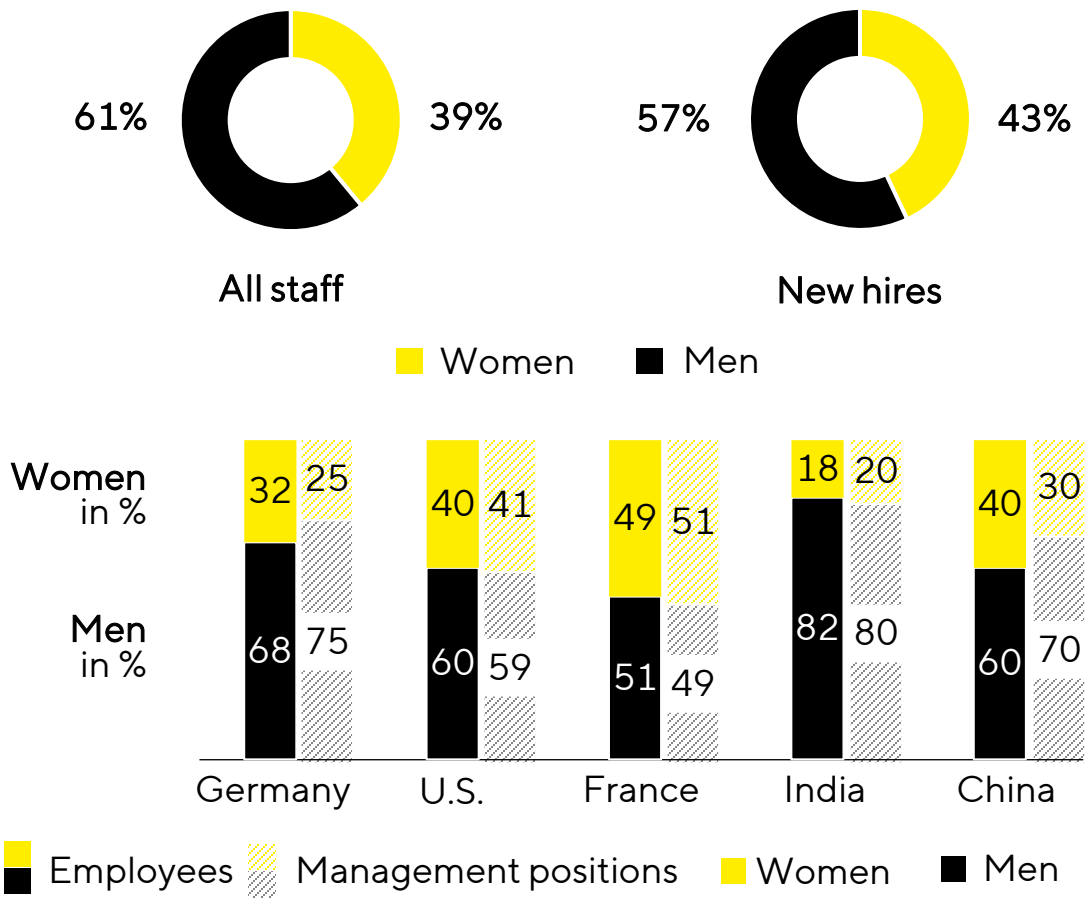
Average tenure in years



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



Gender | Structure Headed in the Right Direction



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

