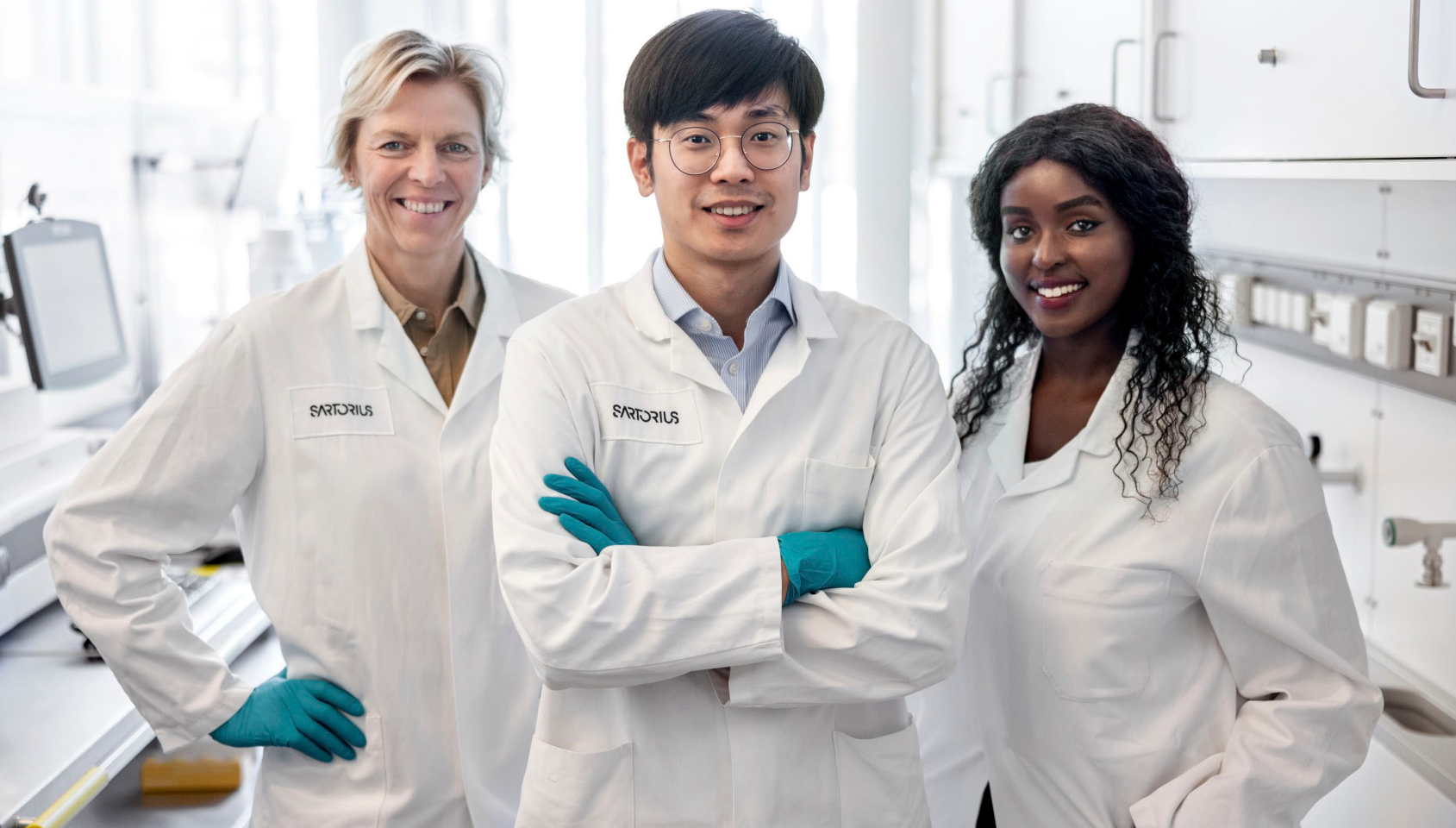




Simplifying Progress

Diversity Fact Sheet

Sartorius 2025

SARTORIUS

"Diversity is a factor of innovation. It brings together different perspectives, ways of thinking, and ideas, making real innovation possible. That's why Sartorius promotes diversity – not as an end in itself, but as a foundation for creative solutions that benefit our customers and patients worldwide."

Michael Grosse, CEO

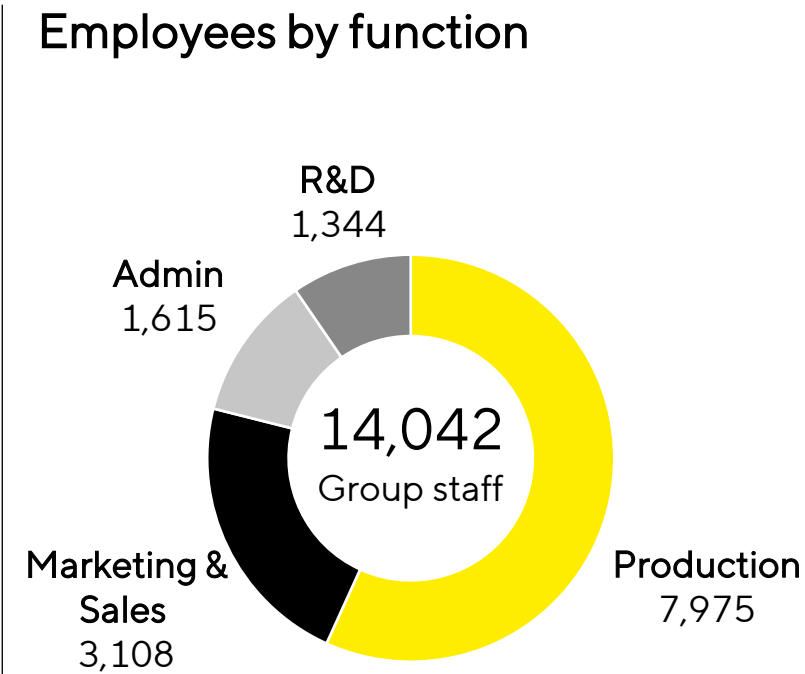
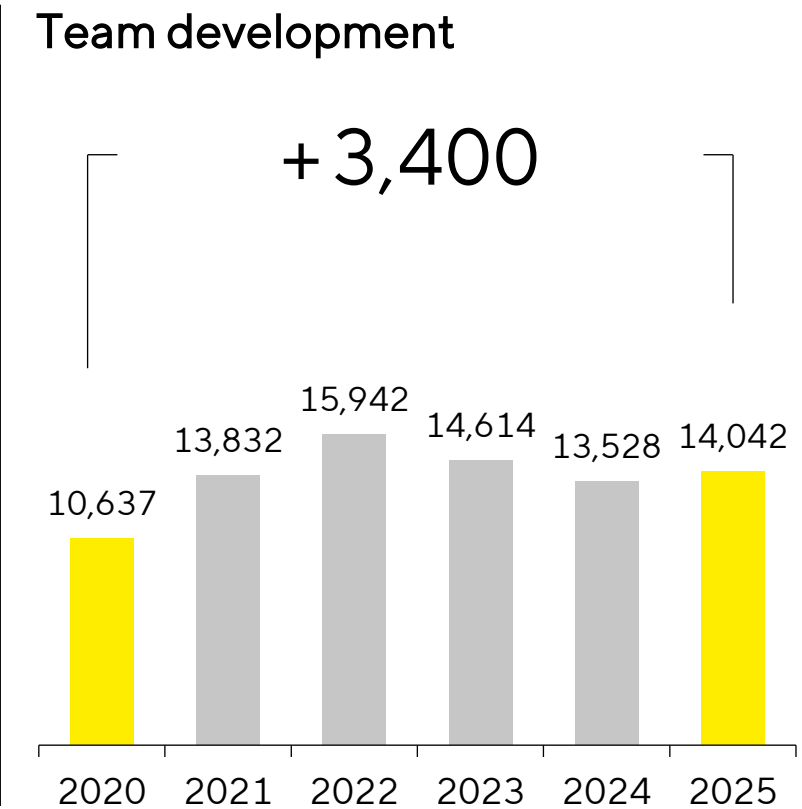


Welcome | 1,639 New People Joined Sartorius in 2025

1,639 | New hires in 2025

906 | Avg. hires annually 2020-2025

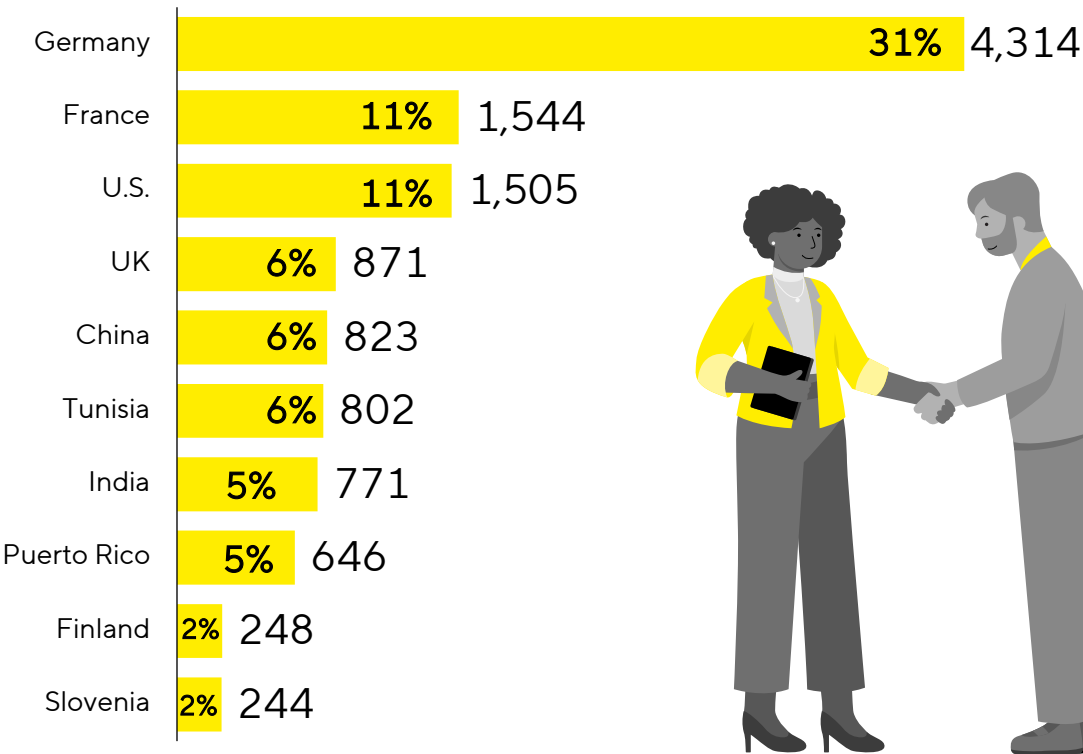
42% | Employees with higher education



Nationality | Sartorius Unites People from 108 Nations

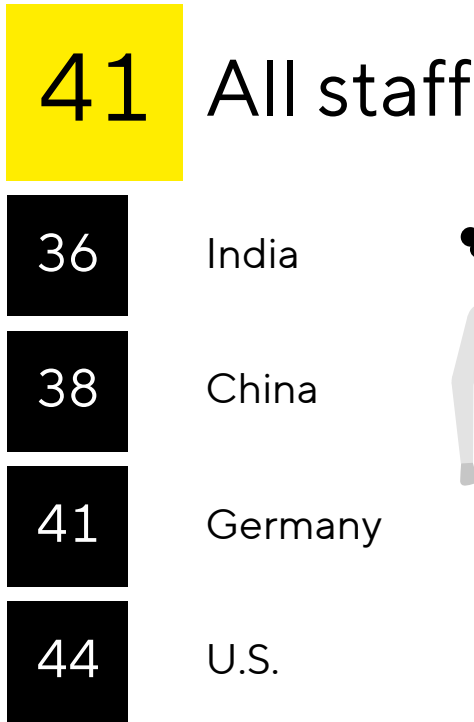


Top 10 nationalities – Share and number of staff

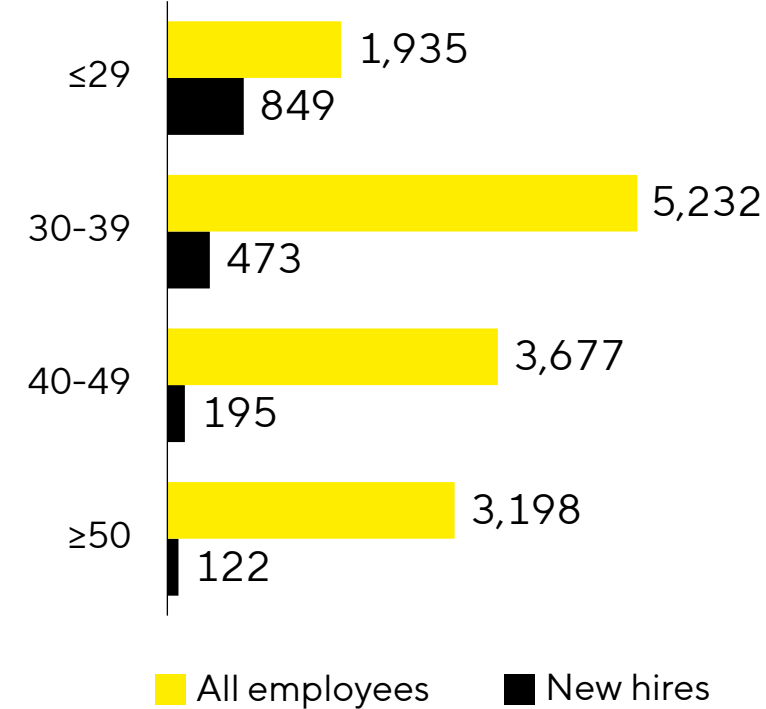


Age | Young Talent and Experienced Professionals Shape Our Workforce

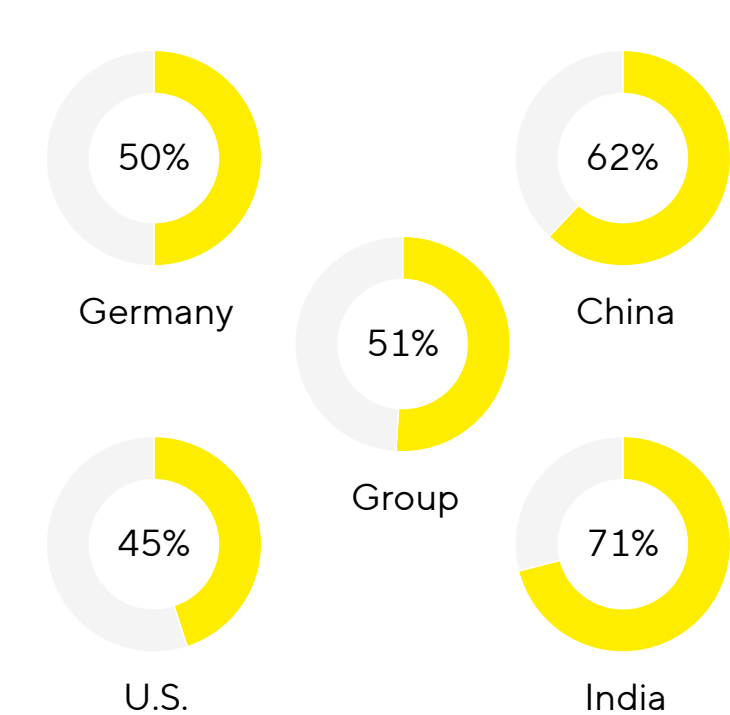
Average age



Number of staff by age

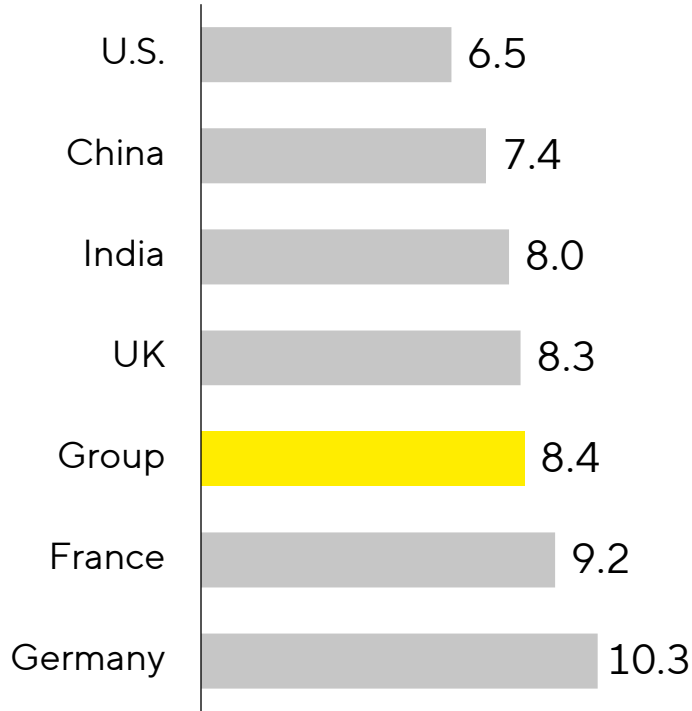


Share of staff younger than 40y

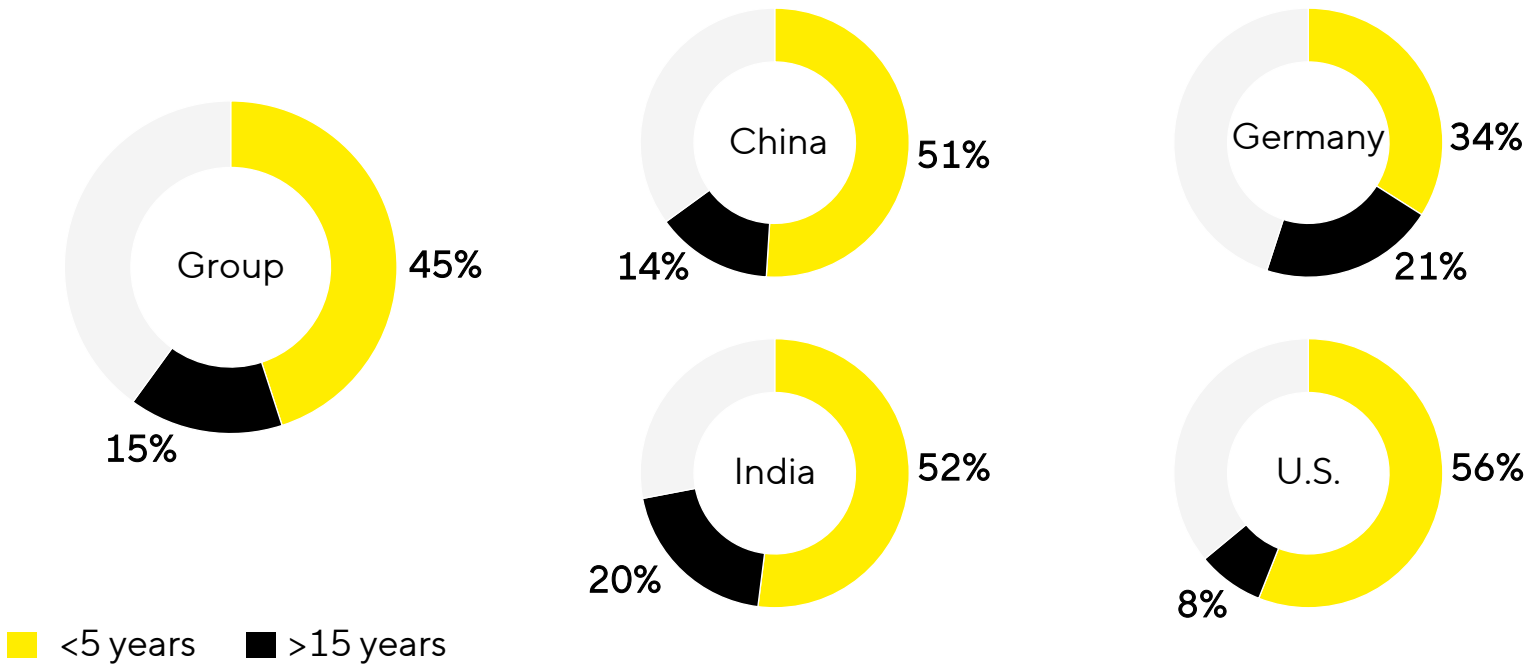


Tenure | Mix of Long-Term Experience and Fresh Ideas

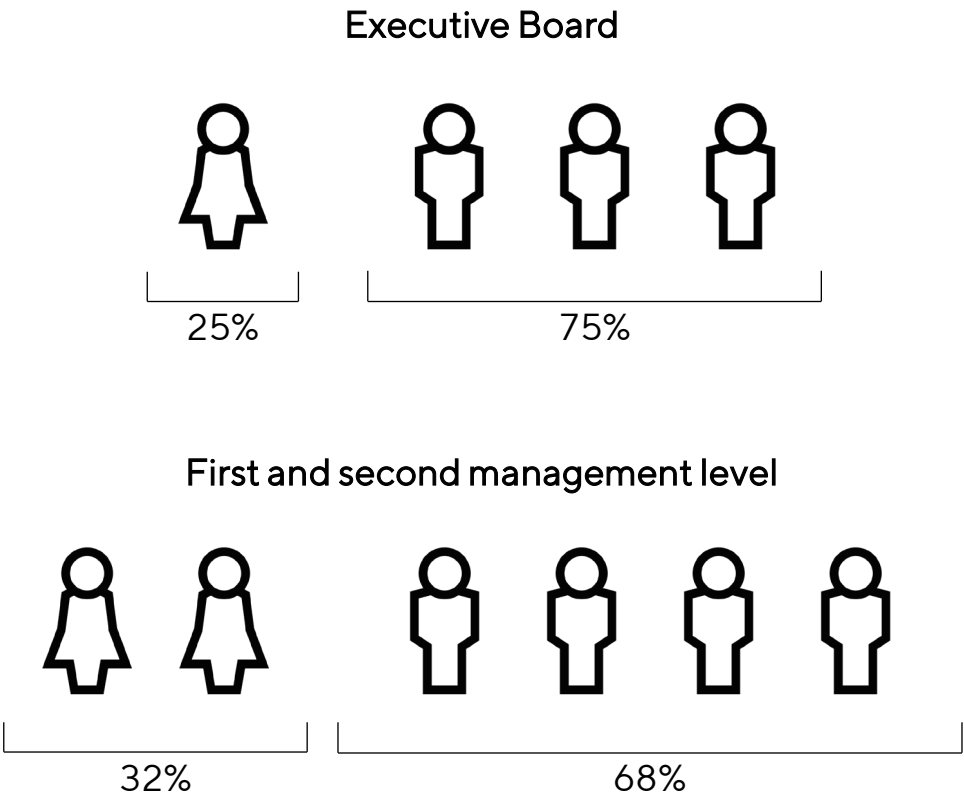
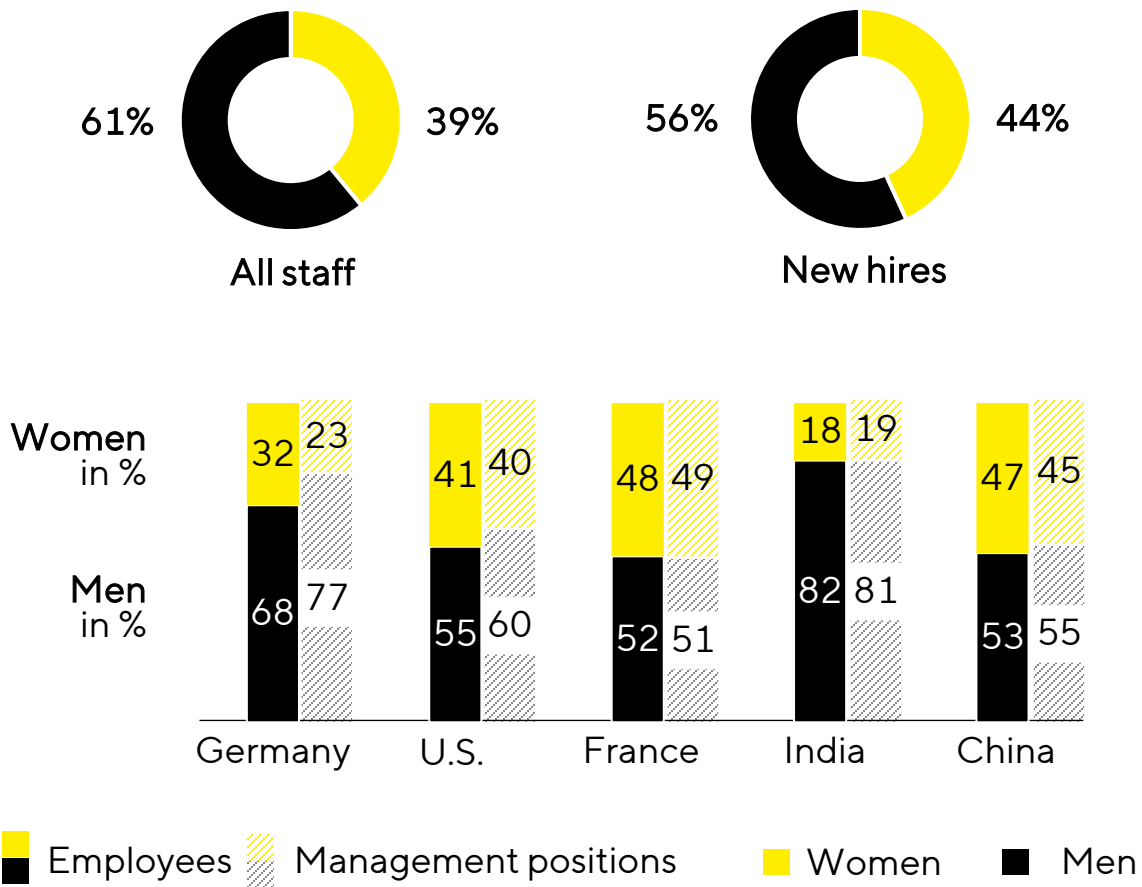
Average tenure in years



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



Gender | Men and Women Represented Across the Organization



Fostering Exchange and Dialogue for Women Internally and Externally

Sartorius Business Women Association

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

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.

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.



Sartorius' employee-led initiatives support in shaping a diverse und inclusive culture. All employees are invited to join.

Partnering in the Name of Diversity and Inclusion

