

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**

---



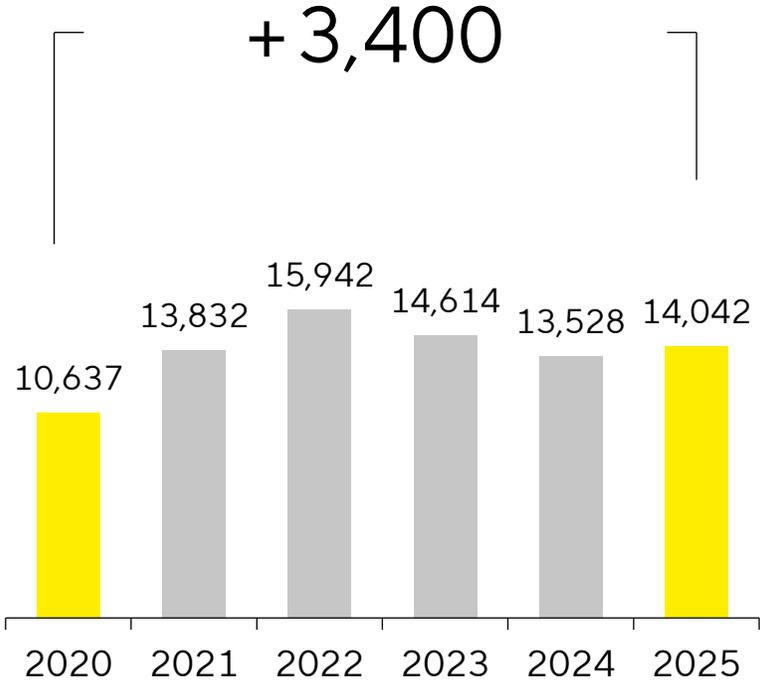
# Welcome | 1,639 New People Joined Sartorius in 2025

1,639 | New hires in 2025

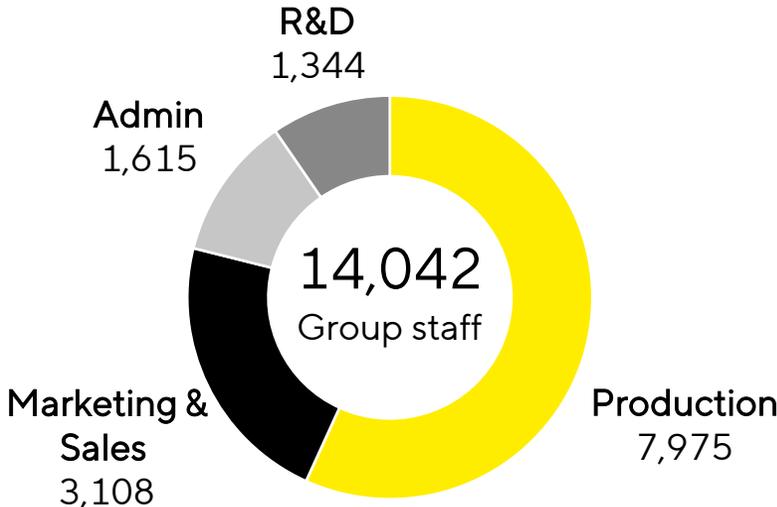
906 | Avg. hires annually 2020-2025

42% | Employees with higher education

Team development



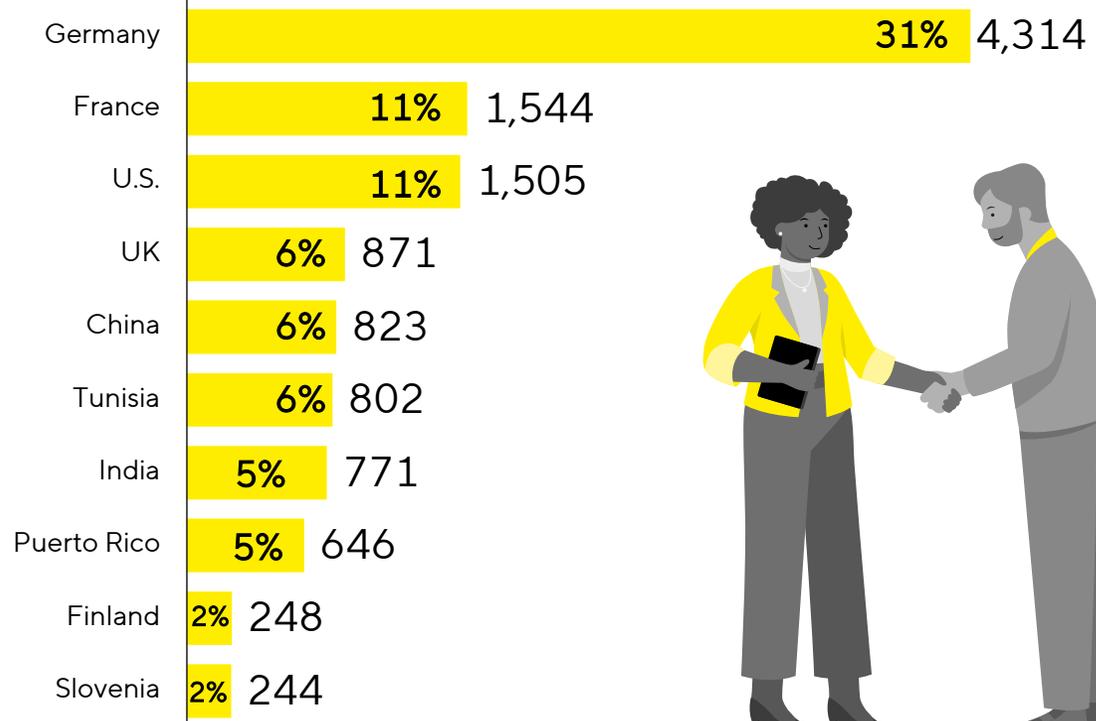
Employees by function



# Nationality | Sartorius Unites People from 108 Nations

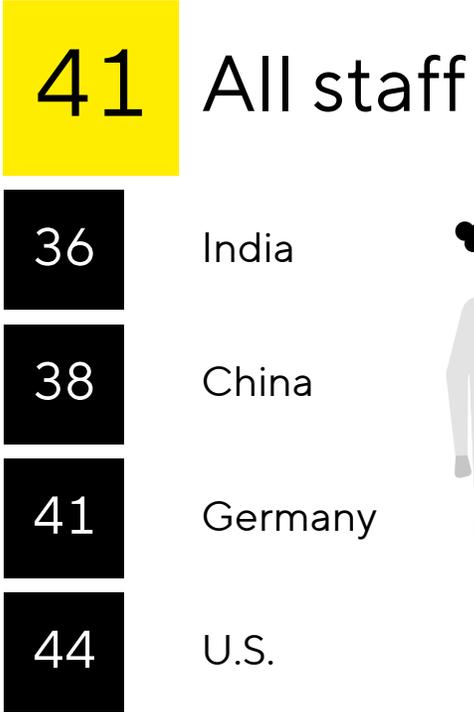


## Top 10 nationalities – Share and number of staff

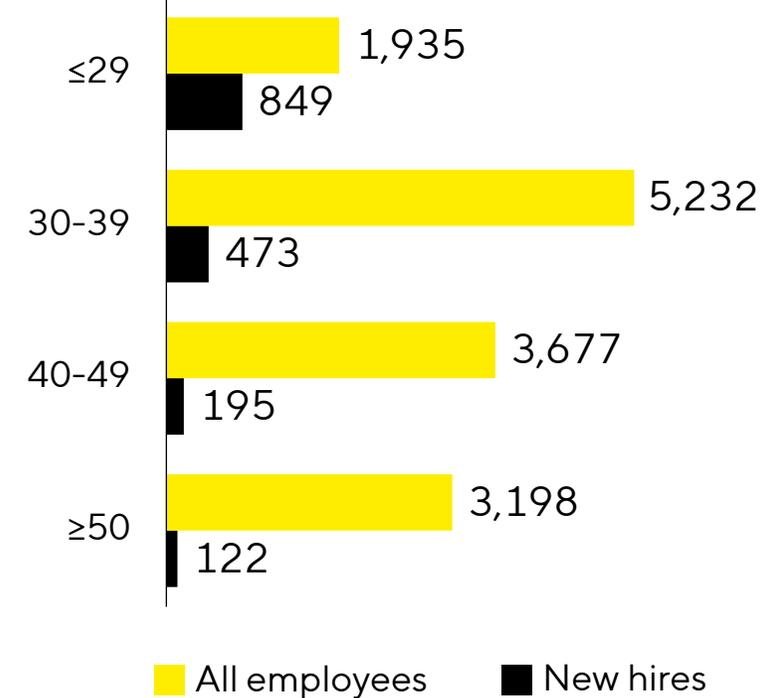


# Age | Young Talent and Experienced Professionals Shape Our Workforce

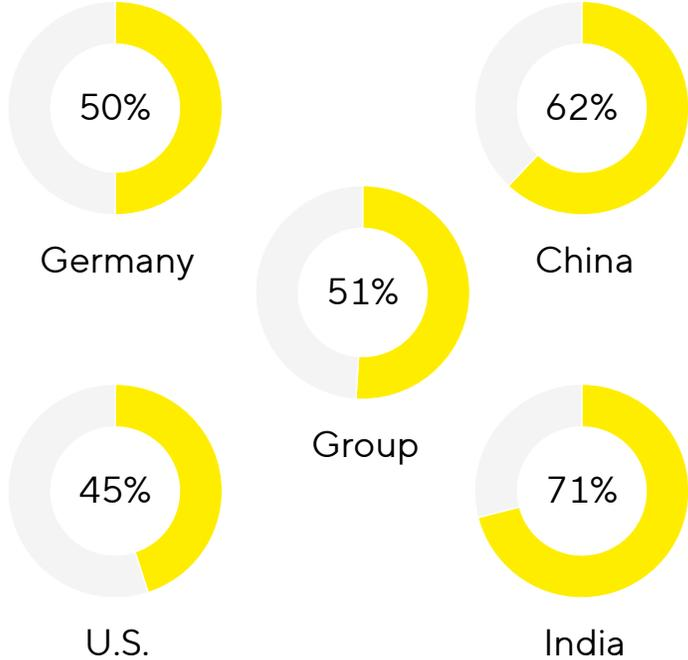
## Average age



## Number of staff by age

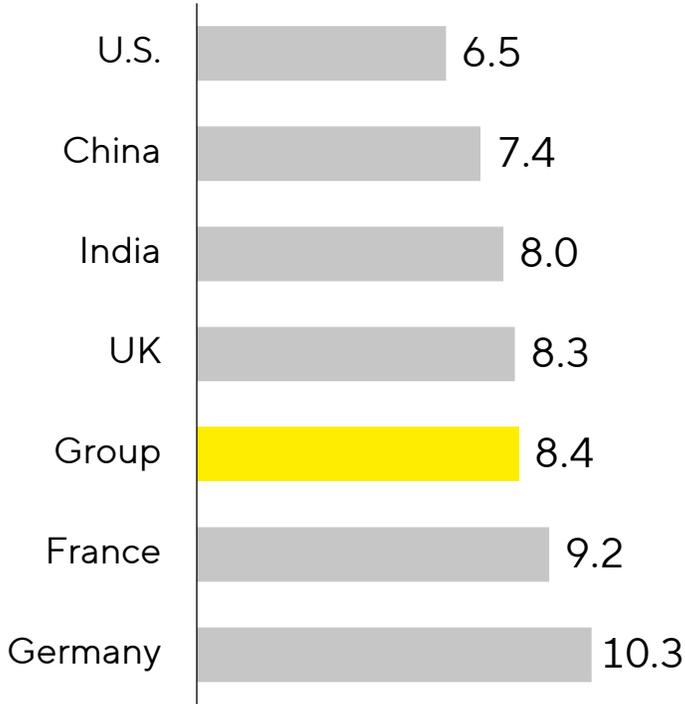


## Share of staff younger than 40y

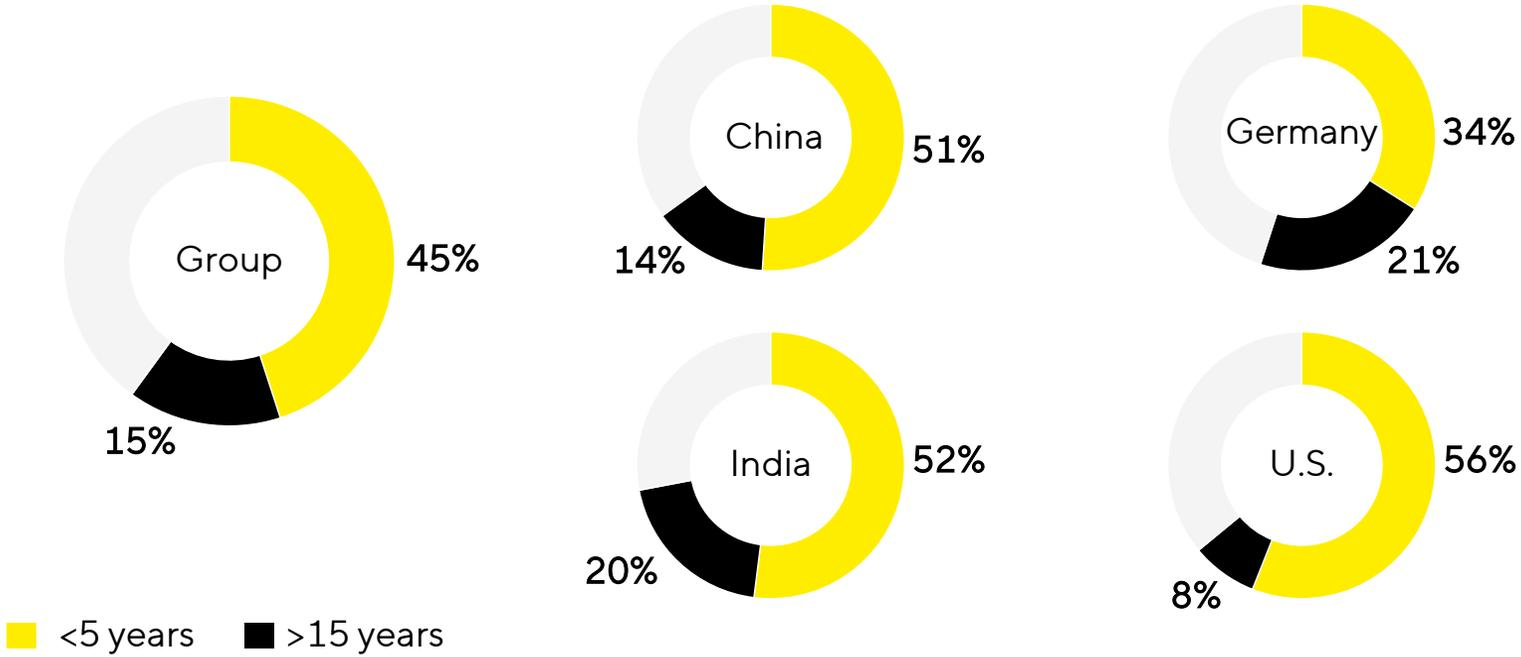


# Tenure | Mix of Long-Term Experience and Fresh Ideas

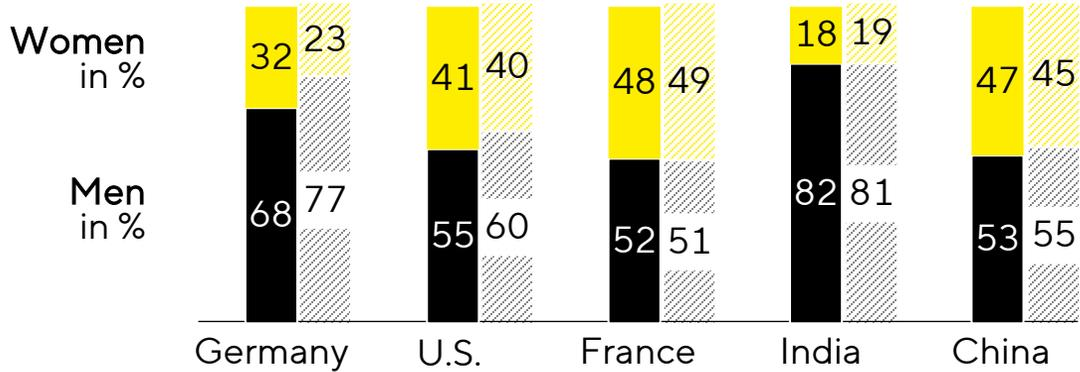
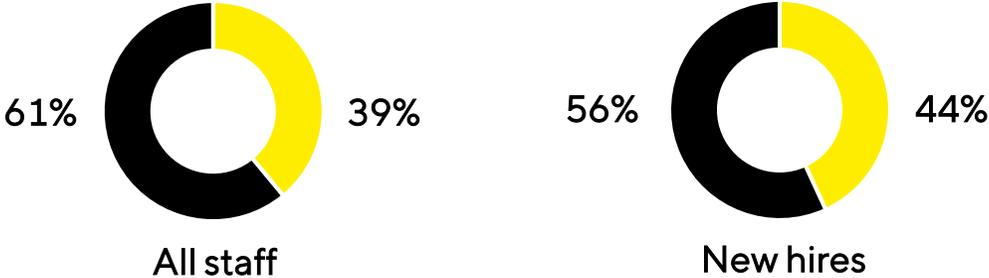
Average tenure in years



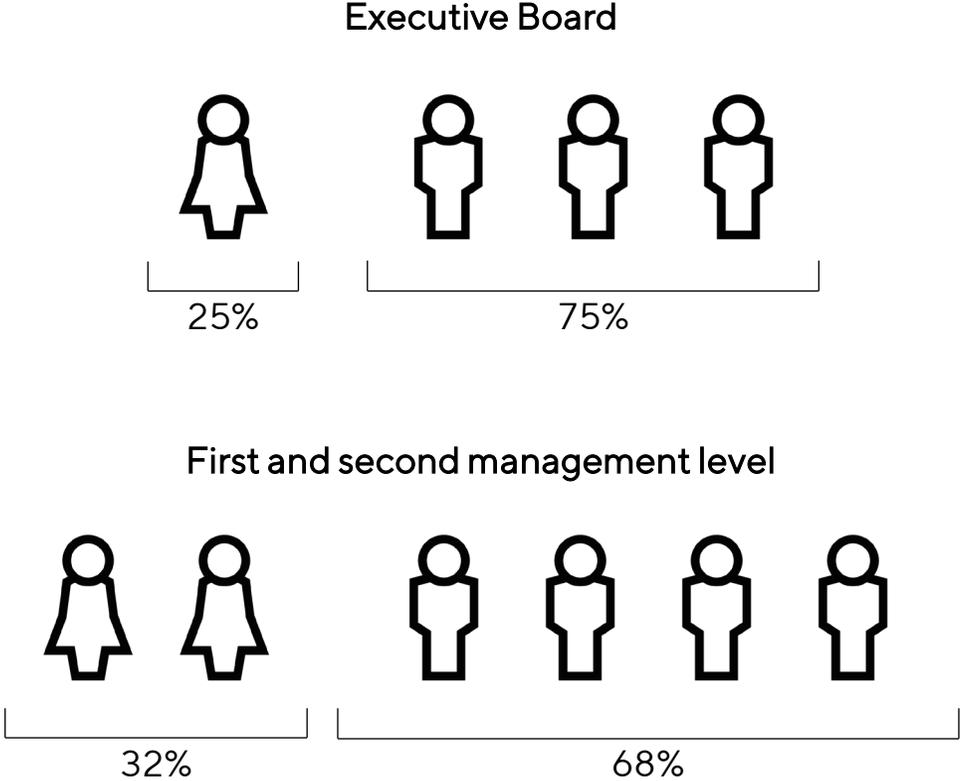
Share of staff with tenure higher than 15 years and lower than 5 years



# Gender | Men and Women Represented Across the Organization



Employees
  Management positions
  Women
  Men



# 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

