

SARTORIUS

Simplifying Progress

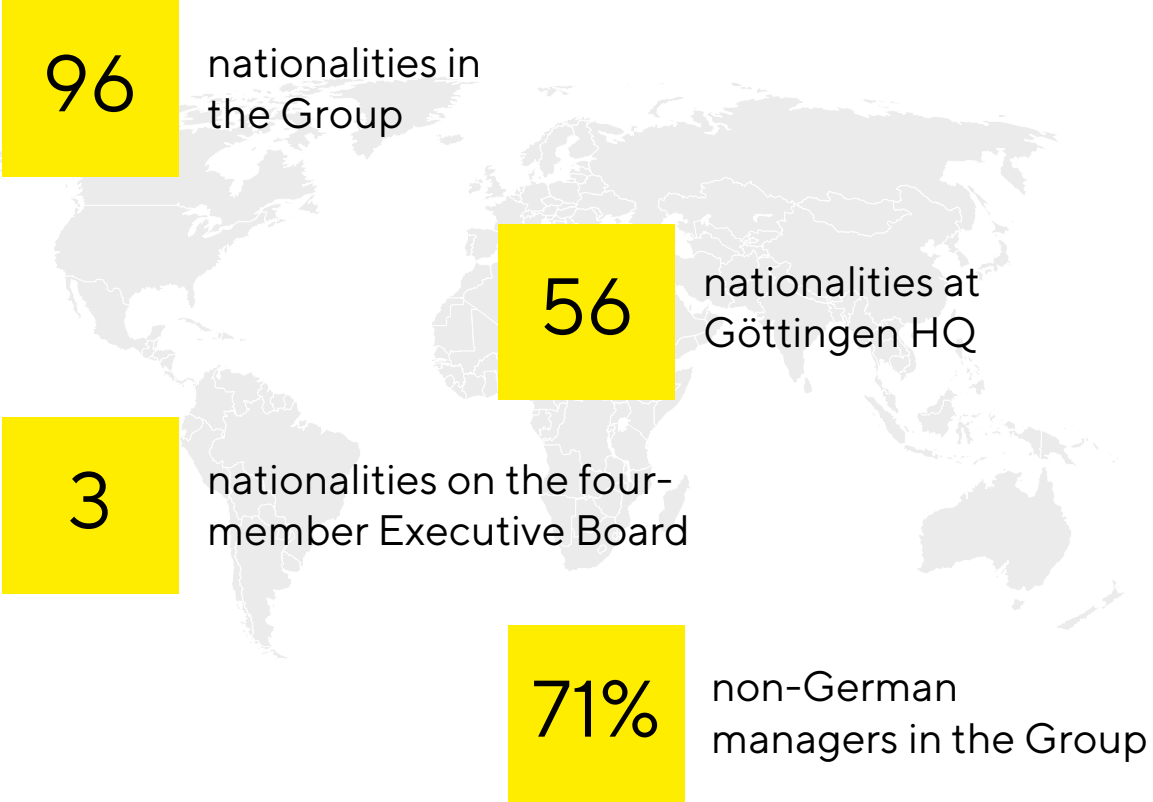
Diversity
Fact Sheet
2020

“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



Nationality Sartorius unites people from 96 nations

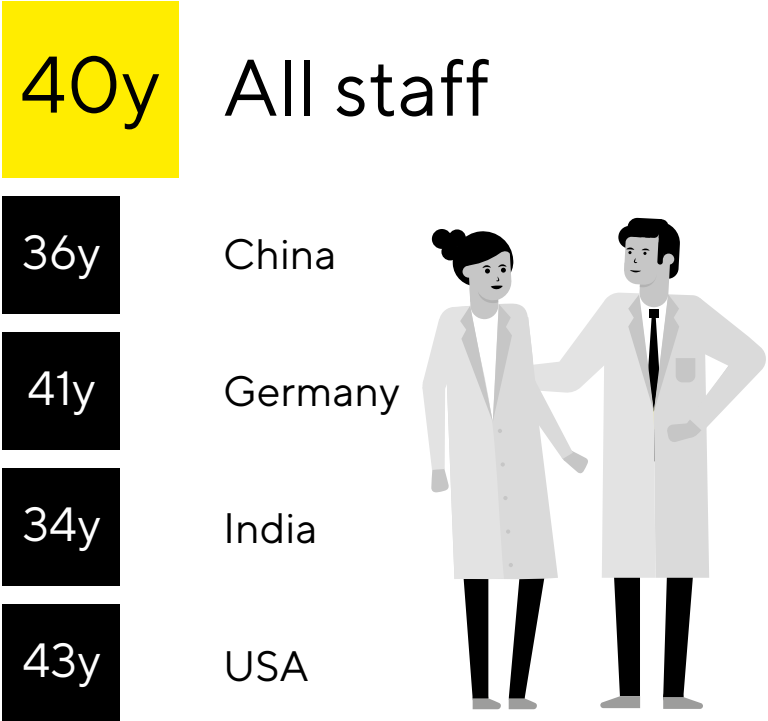


Top 10 nationalities - Share and number of staff

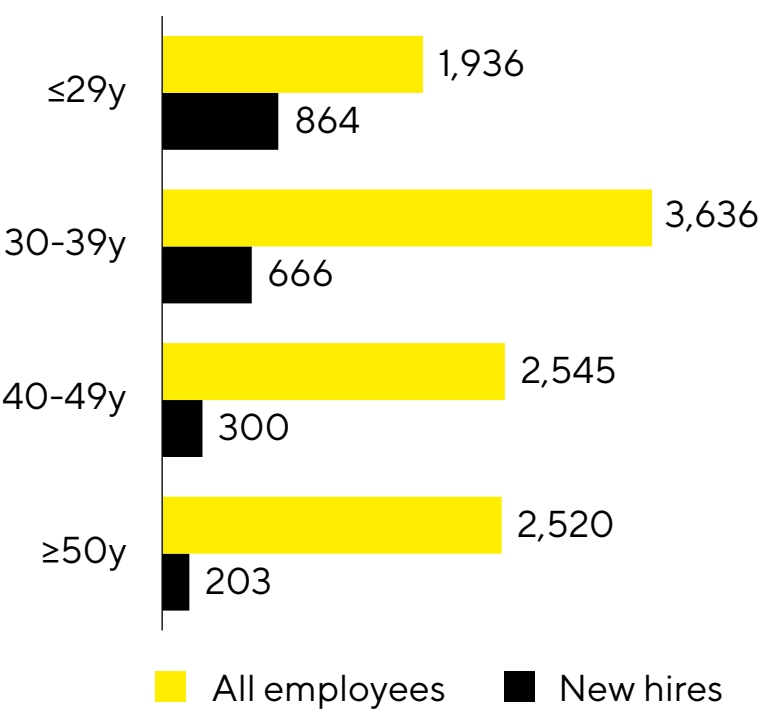
German	32%	3,445
American	16%	1,671
French	10%	1,066
British	7%	722
Chinese	6%	641
Indian	5%	577
Tunisian	4%	392
Finnish		253
Israeli		217
Korean		162

Age China and India are emerging talent markets

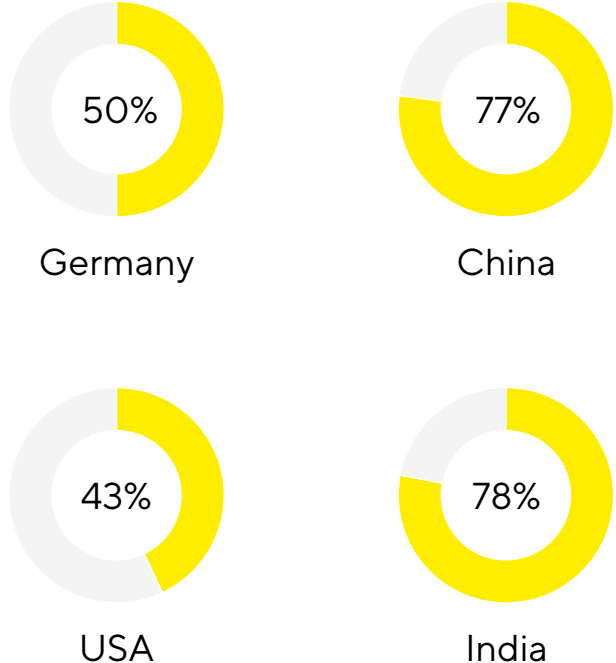
Average age



Number of staff by age

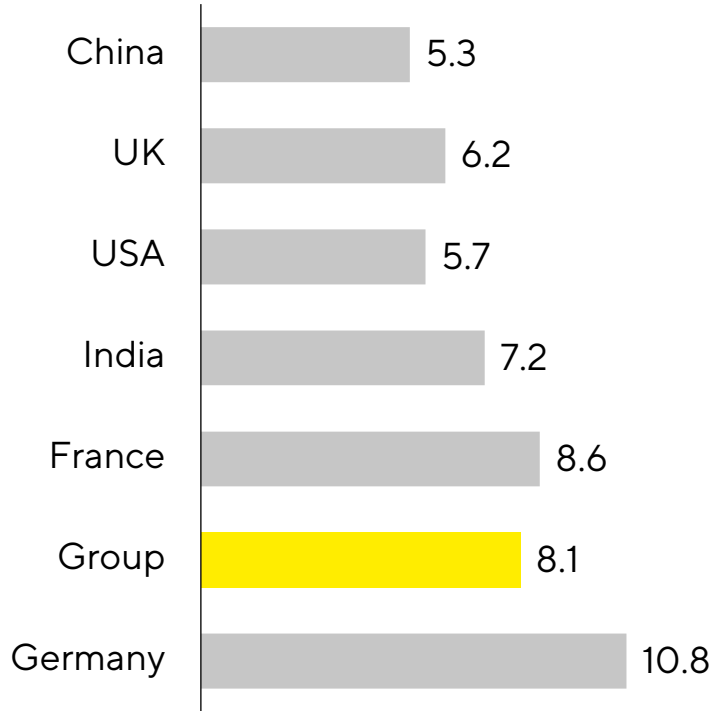


Share of staff younger than 40 yrs

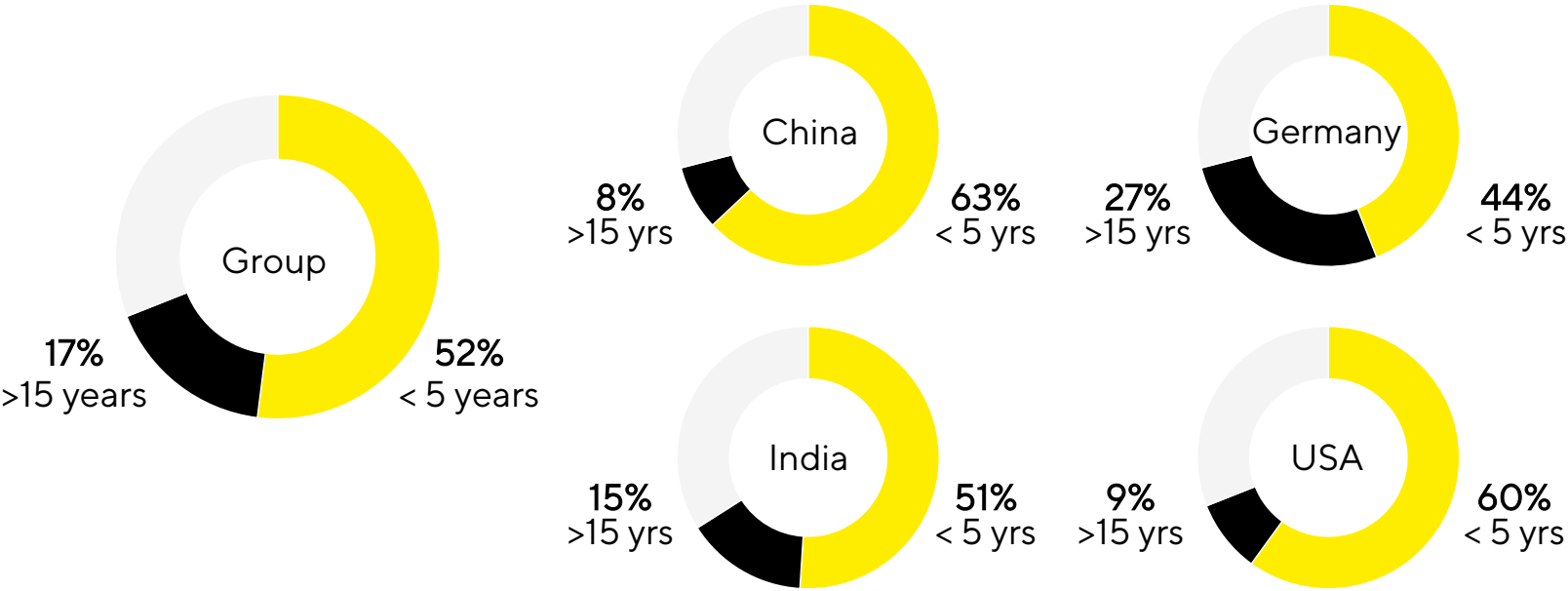


Tenure Healthy mixture of experience and fresh ideas

Average tenure in years

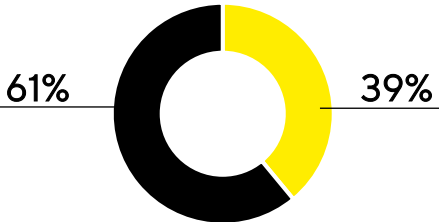


Share of staff with tenure higher than 15 yrs and lower than 5 yrs

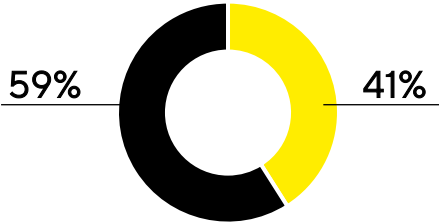


Gender 41% of new hires are women, 24% women in senior management

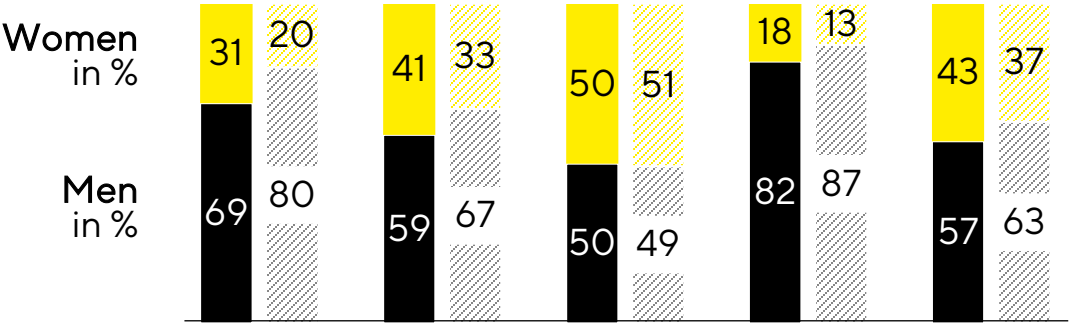
All staff



New hires

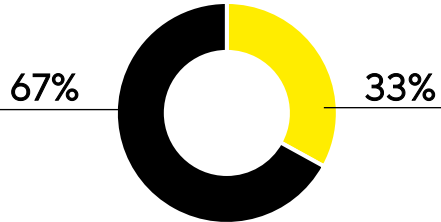


Women Men

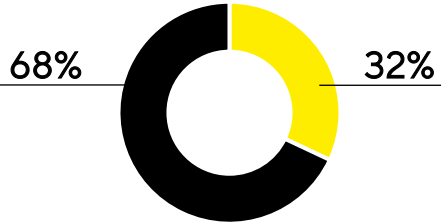


Employees Management positions

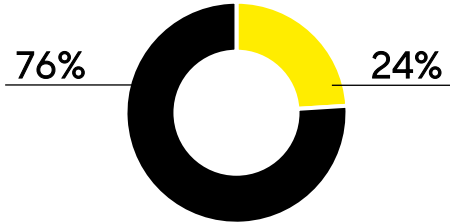
Promotions



Management



Senior Management

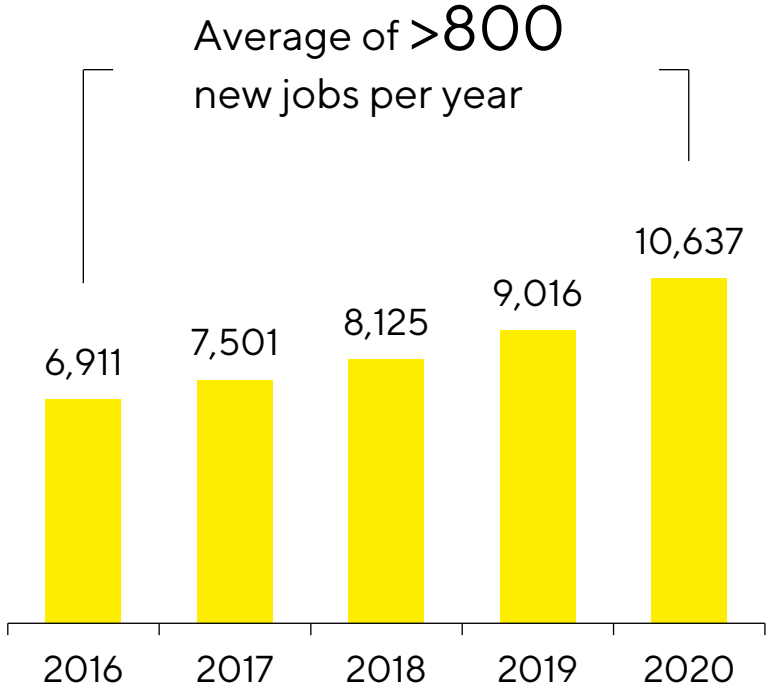


Women Men Target 2022: 30%

Welcome 2,515 new people joined Sartorius in 2020

- 2,033 | New hires in 2020
- 482 | New employees from M&A in 2020
- 1,458 | Avg. hires annually 2016-20
- 39% | Employees with higher education

Employee growth



Employees by function

