

SARTORIUS

Simplifying Progress

Diversity Fact Sheet

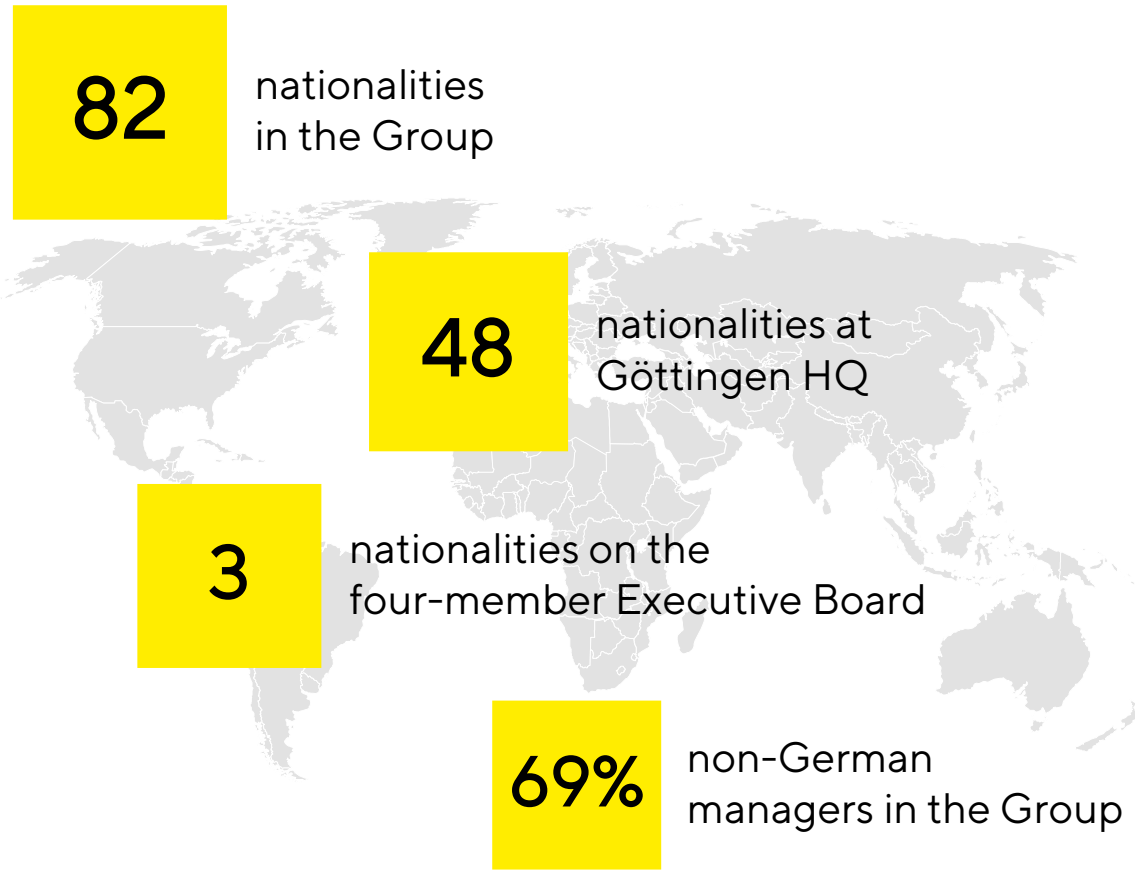
Sartorius 2019

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.

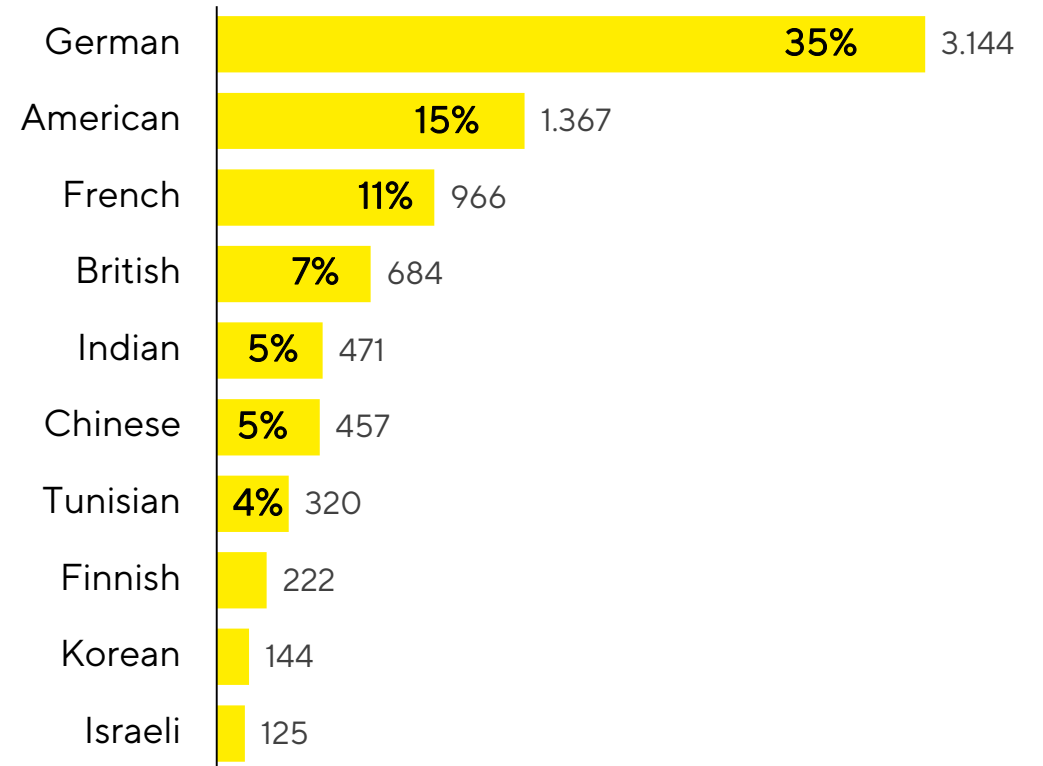
Joachim Kreuzburg, CEO



Nationality Sartorius unites people from 82 nations



Top 10 nationalities - Share and number of staff

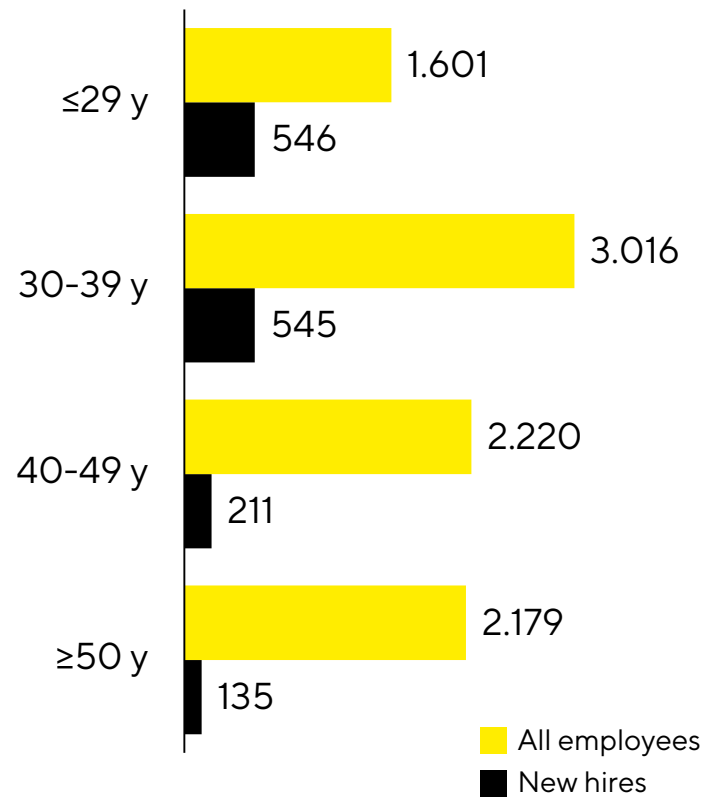


Age China and India are emerging talent markets

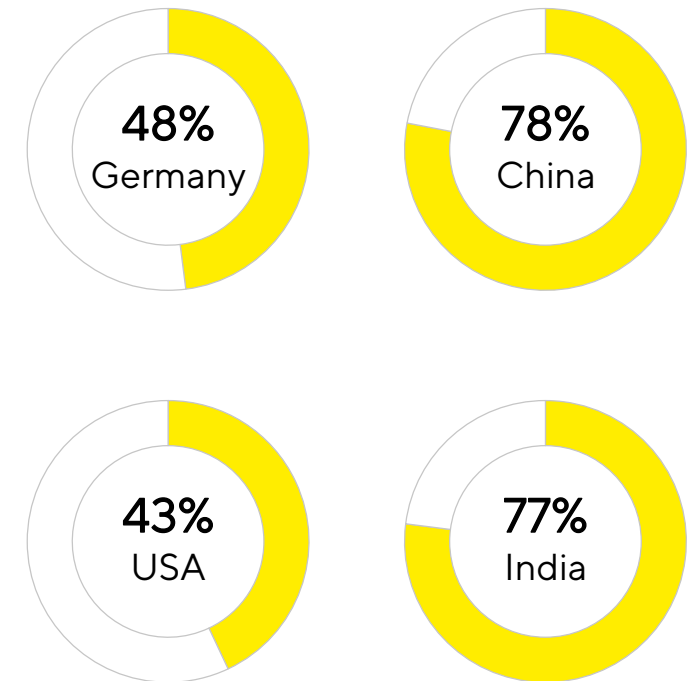
Average age



Number of staff by age

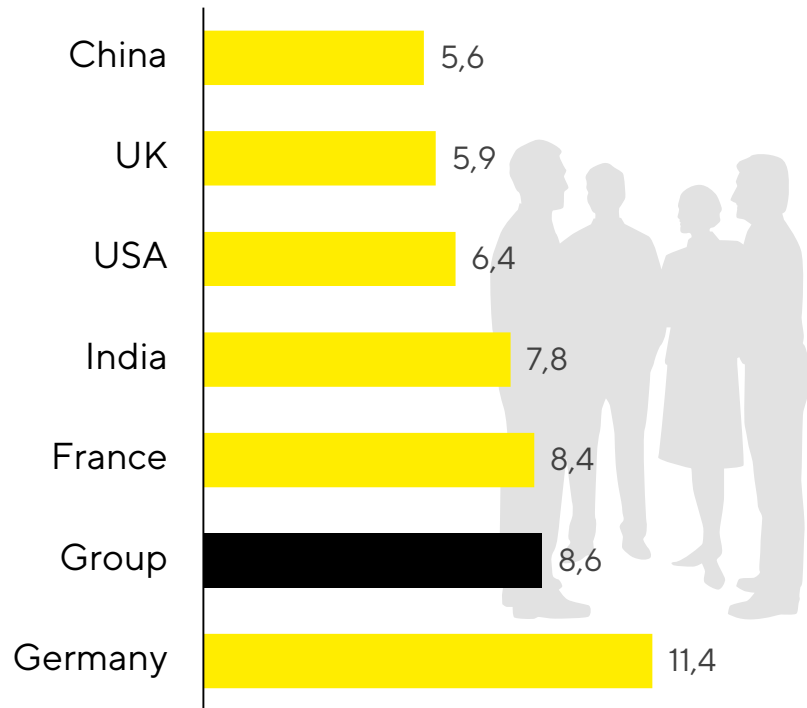


Share of staff younger than 40 yrs

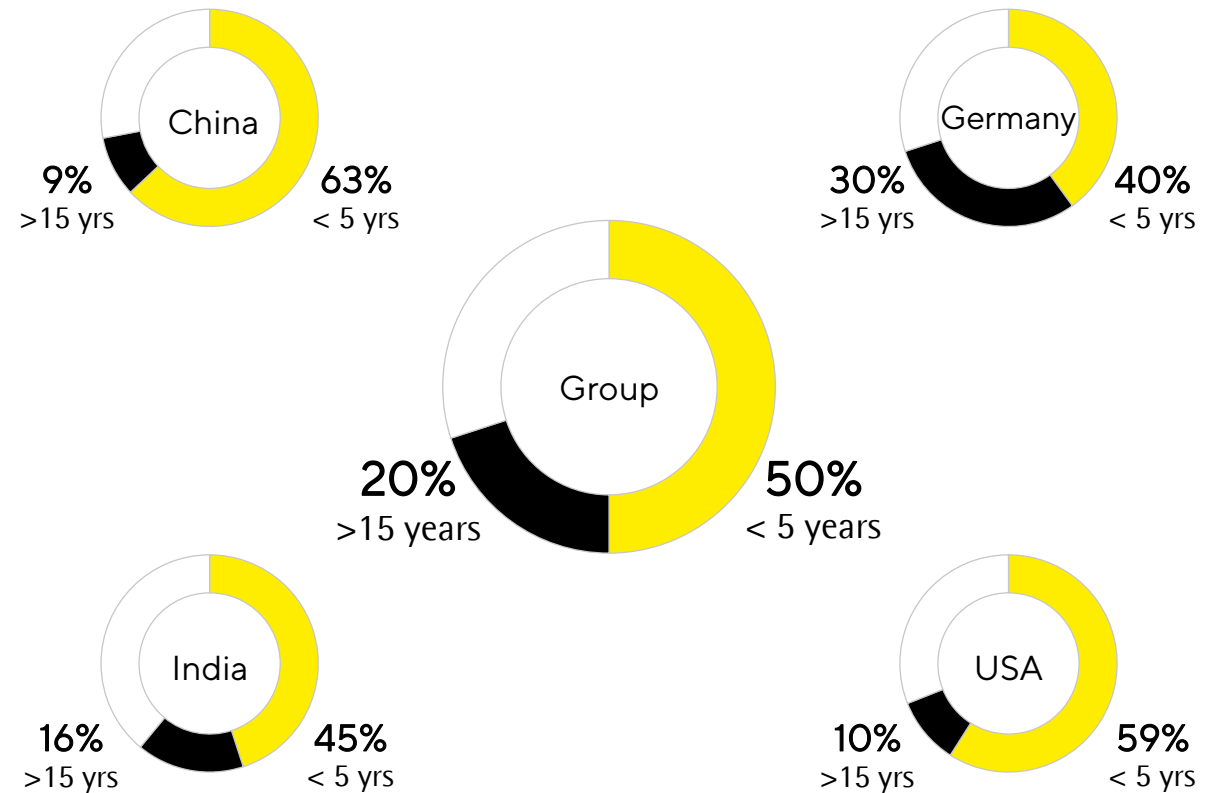


Tenure Healthy mixture of experience and fresh ideas

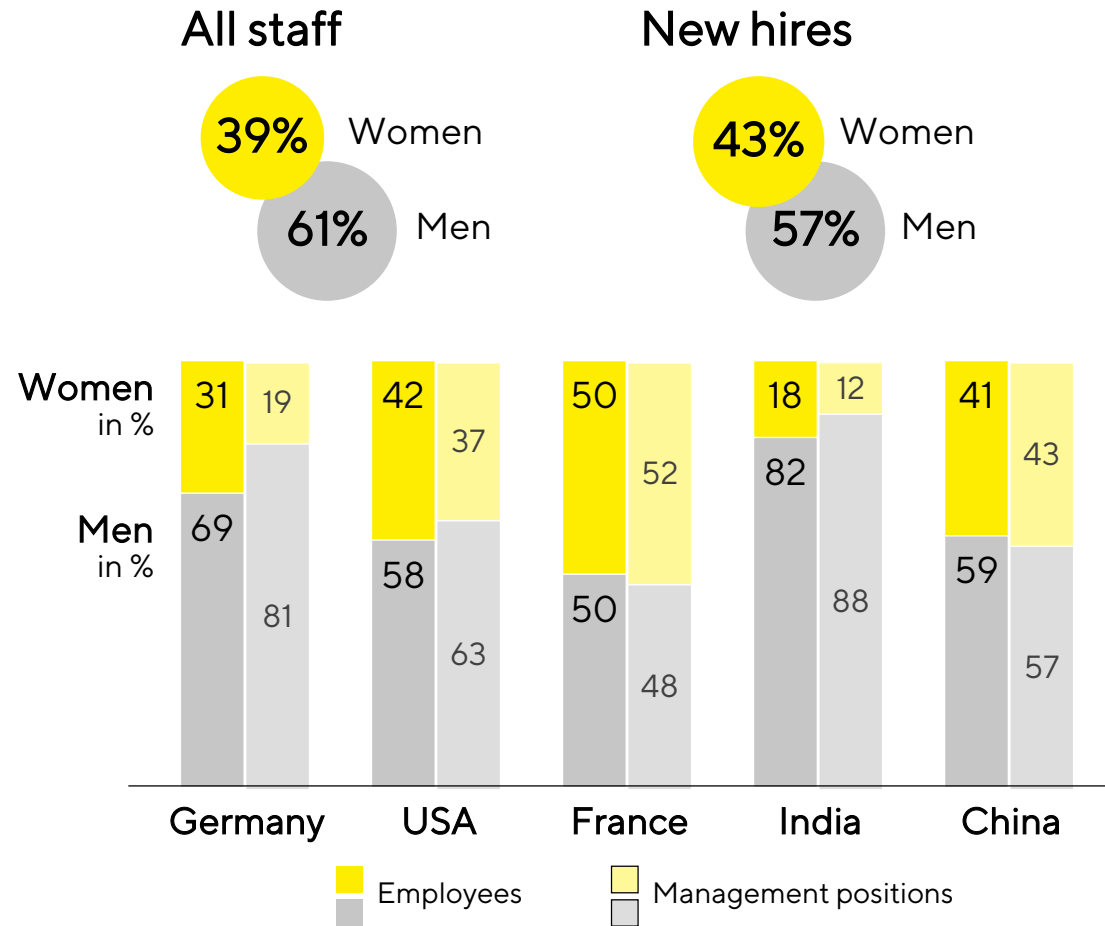
Average tenure in years



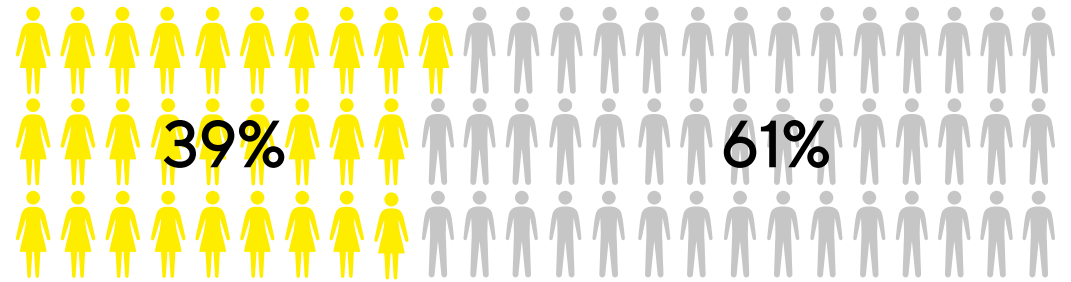
Share of staff with tenure higher than 15 yrs and lower than 5 yrs



Gender 43% of new hires are women, 28% women in senior management



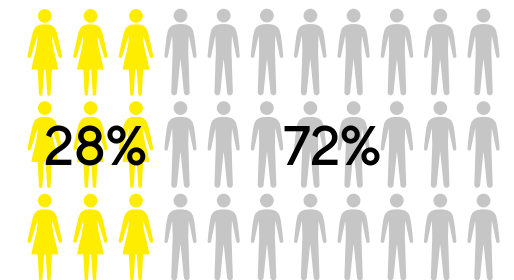
Promotions



Management

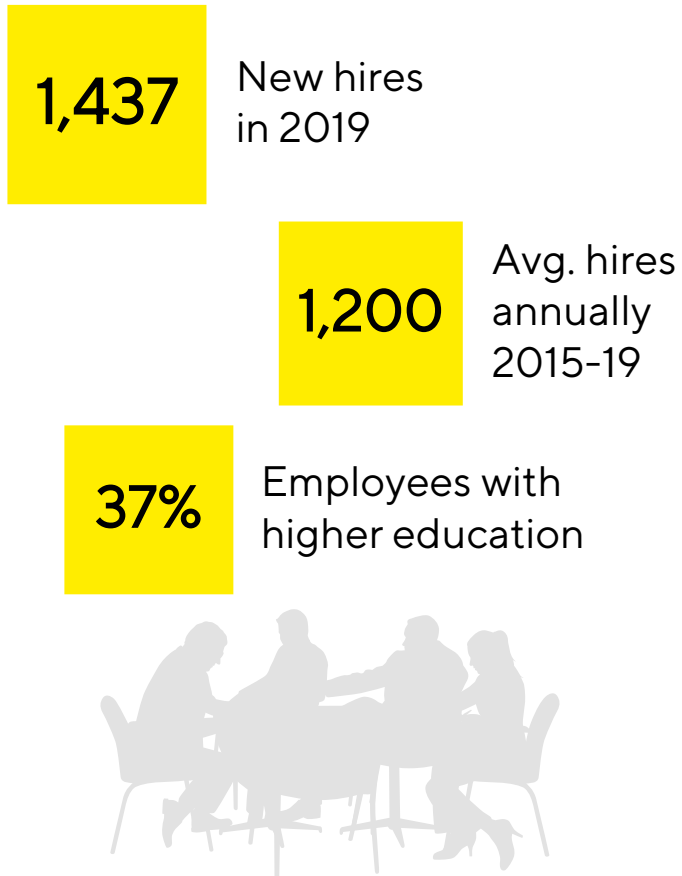


Senior Management

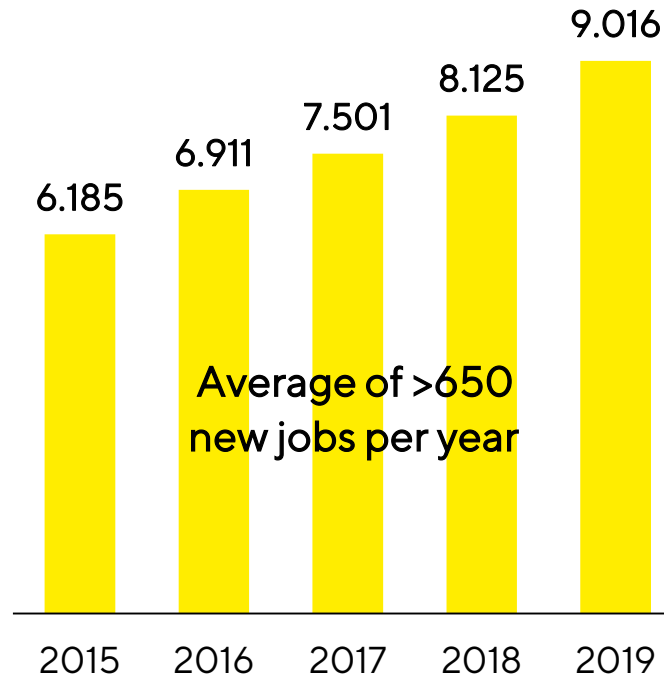


Target 2022: 30%

Welcome 1,437 new people joined Sartorius in 2019



Employee growth



Employees by function

