

# SARTORIUS

Simplifying Progress

## Diversity Fact Sheet

Sartorius 2019

---

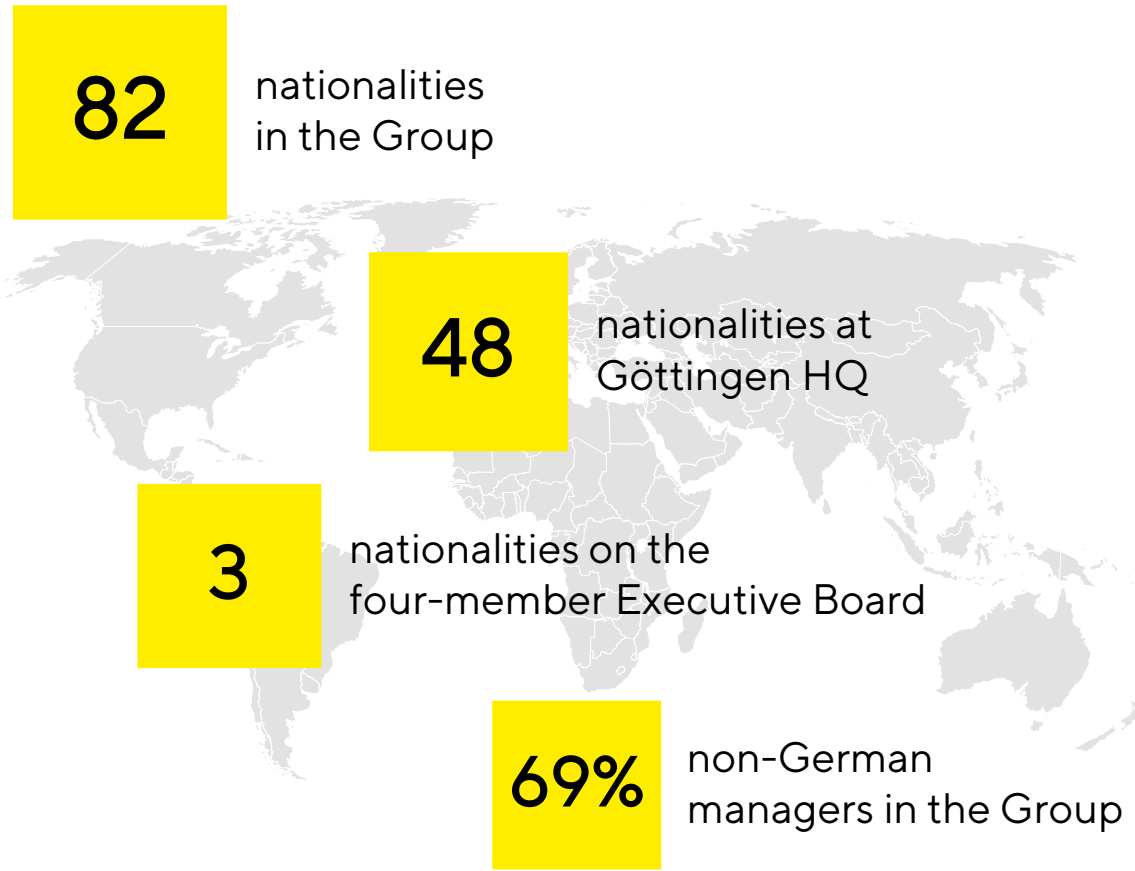
At Sartorius, diversity is not about statistics.  
It's about culture. We foster diversity  
because it ultimately teaches us to be  
more open. And we consider openness  
as a key prerequisite for success.

---

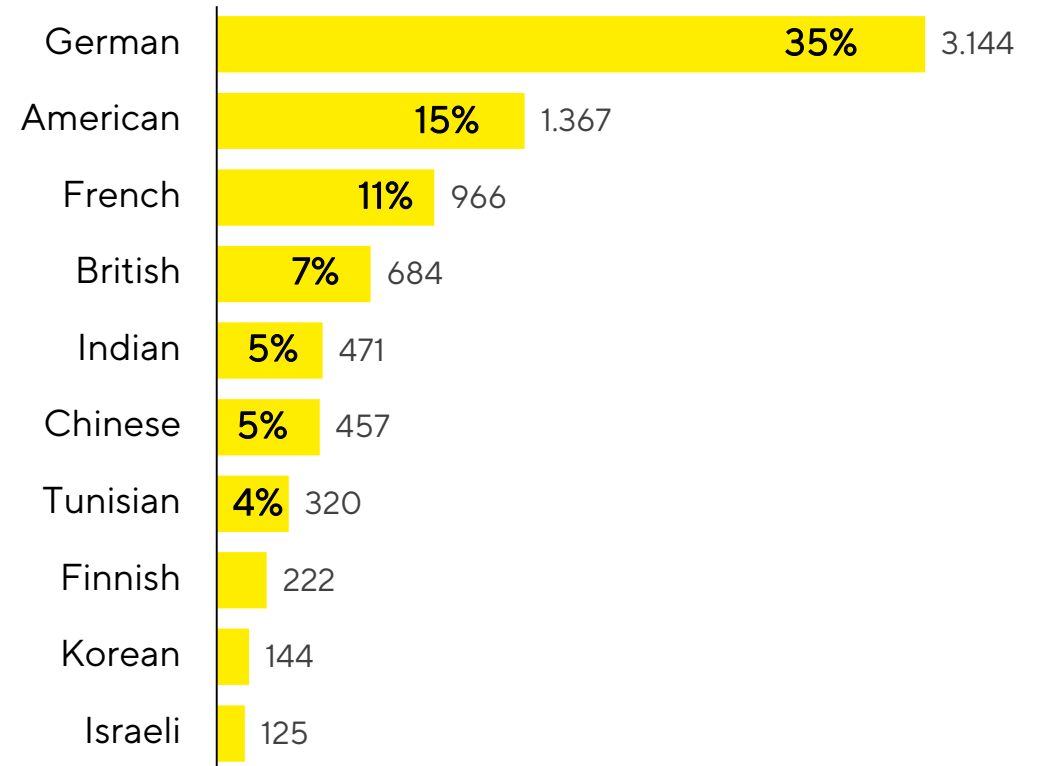
Joachim Kreuzburg, CEO



# Nationality Sartorius unites people from 82 nations



## Top 10 nationalities - Share and number of staff

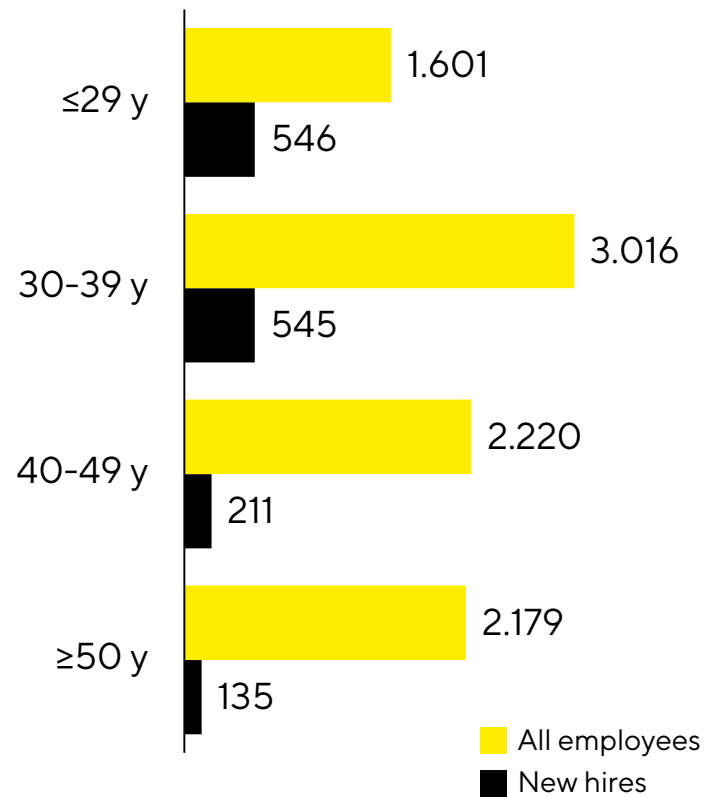


# Age China and India are emerging talent markets

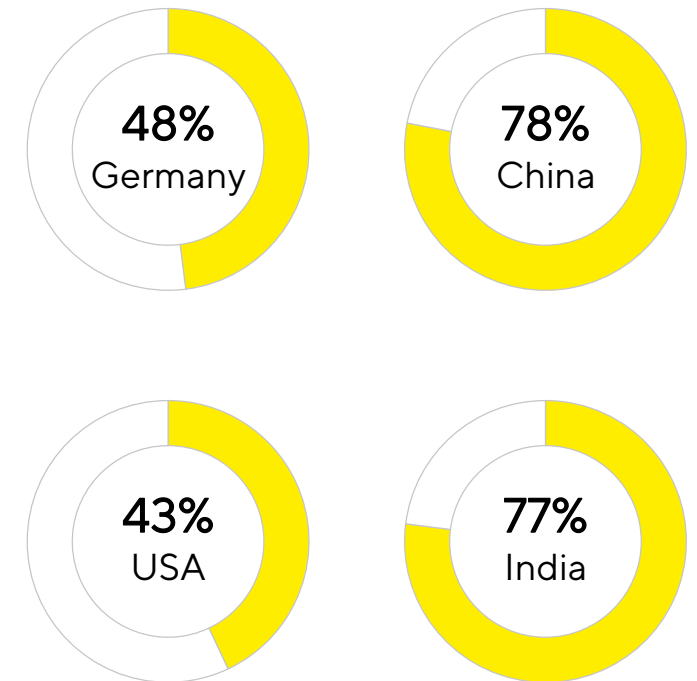
Average age



Number of staff by age

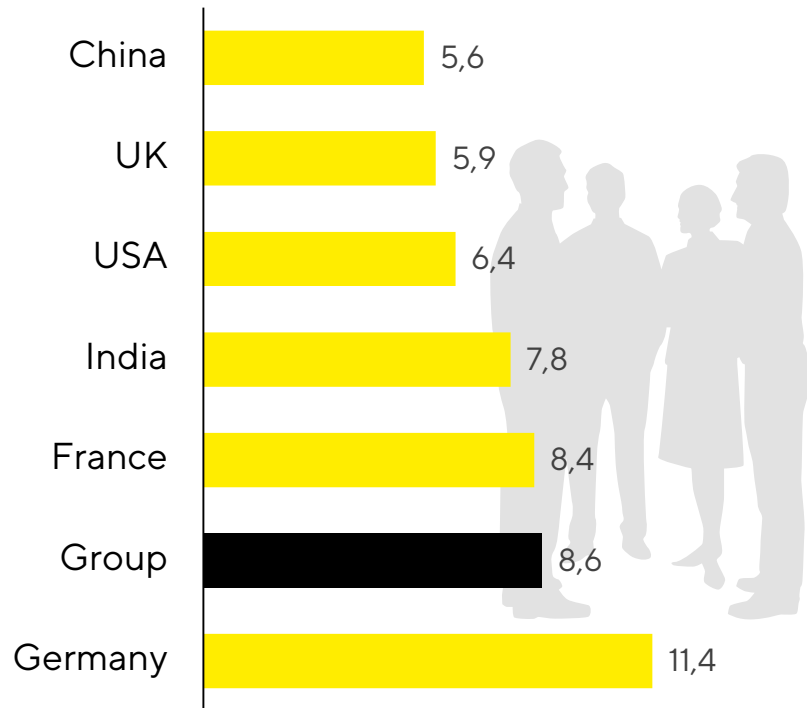


Share of staff younger than 40 yrs

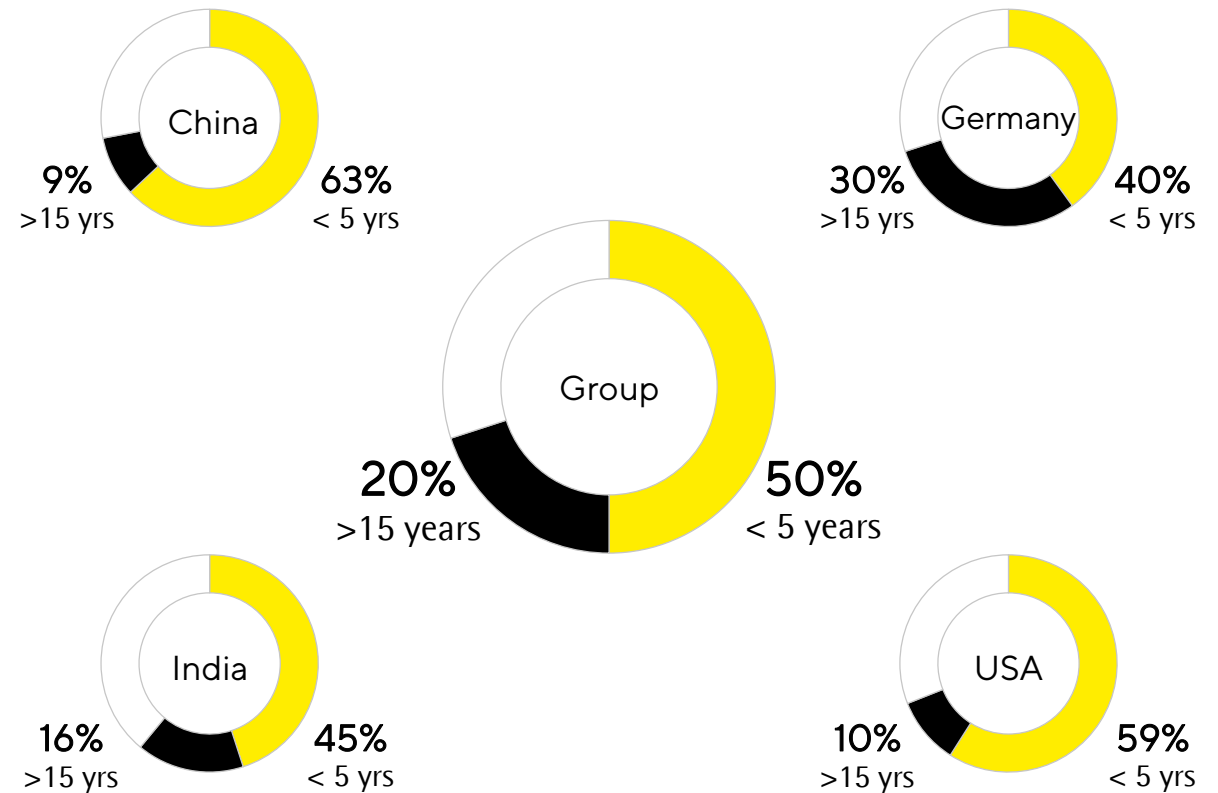


# Tenure Healthy mixture of experience and fresh ideas

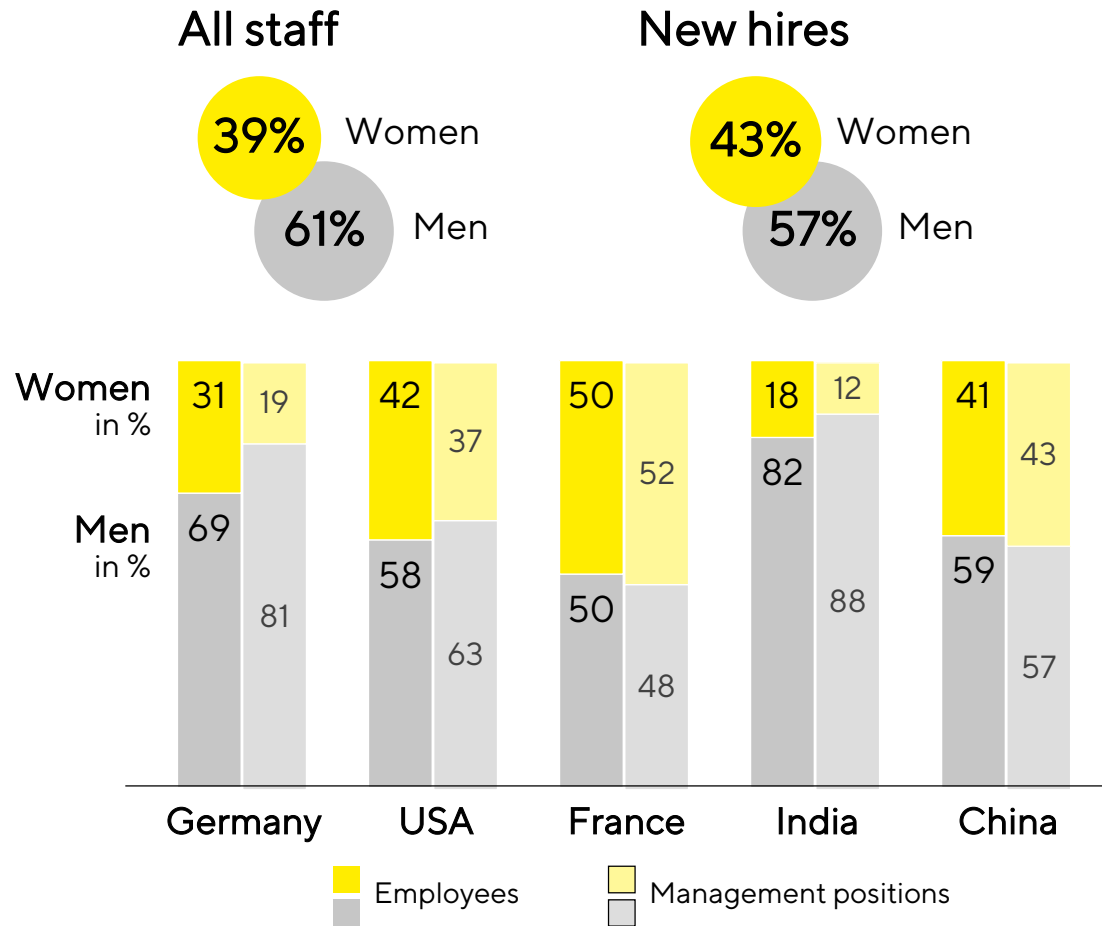
## Average tenure in years



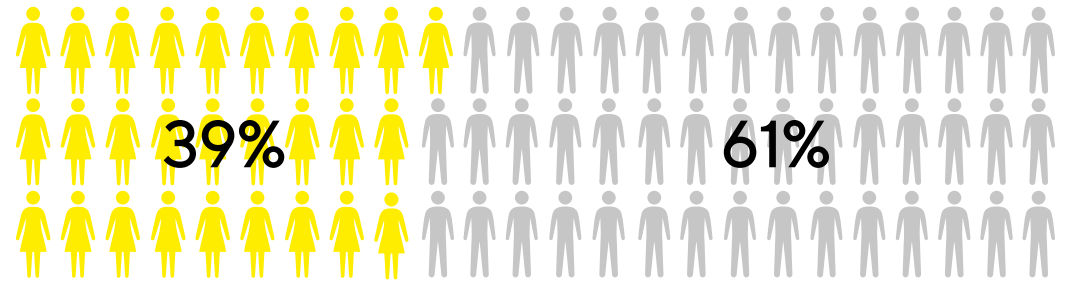
## Share of staff with tenure higher than 15 yrs and lower than 5 yrs



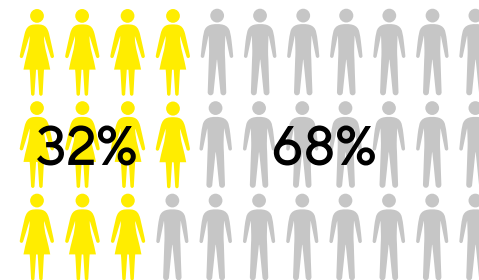
# Gender 43% of new hires are women, 28% women in senior management



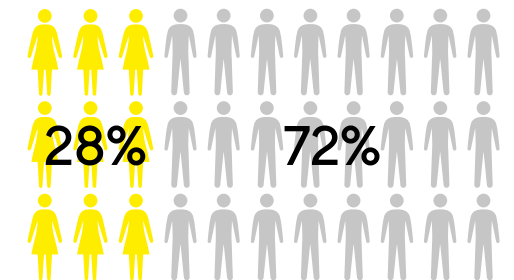
## Promotions



## Management

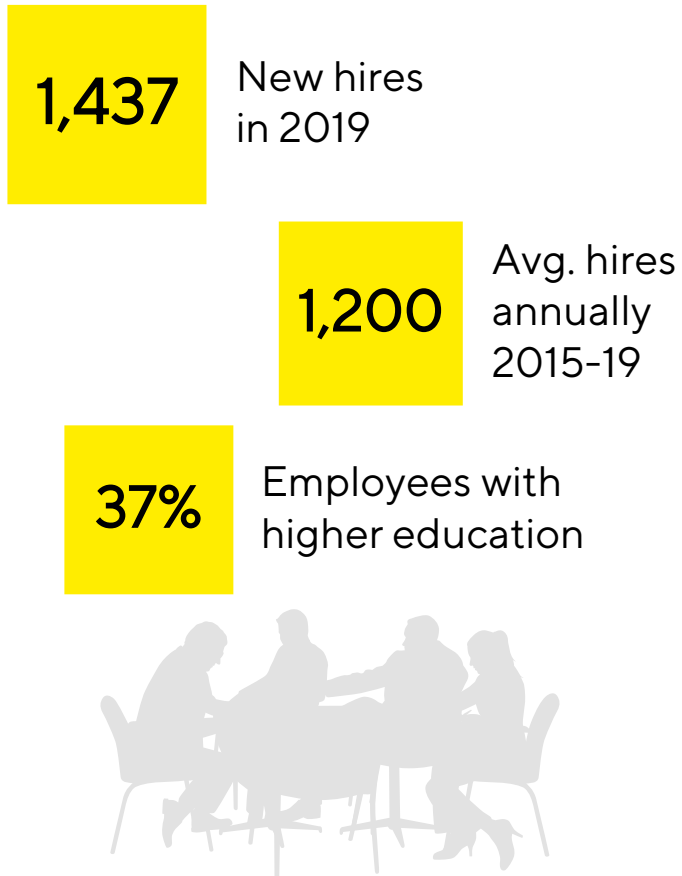


## Senior Management

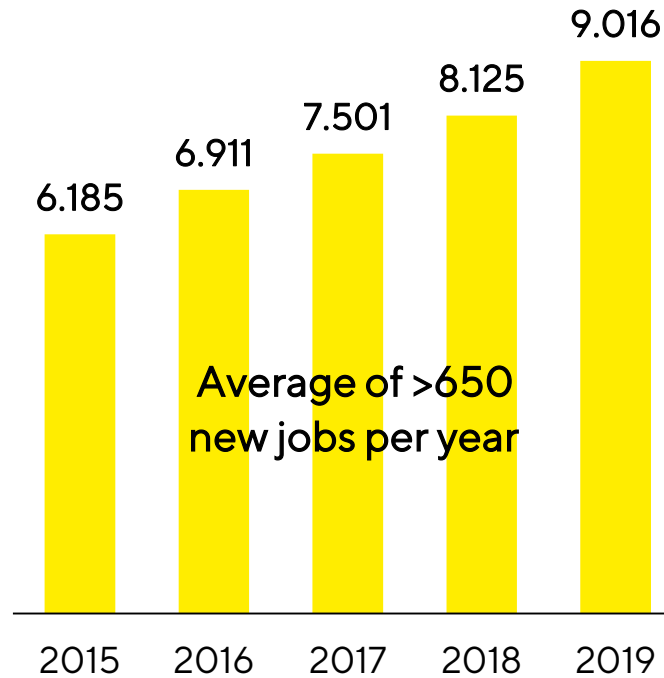


Target 2022: 30%

# Welcome 1,437 new people joined Sartorius in 2019



## Employee growth



## Employees by function

