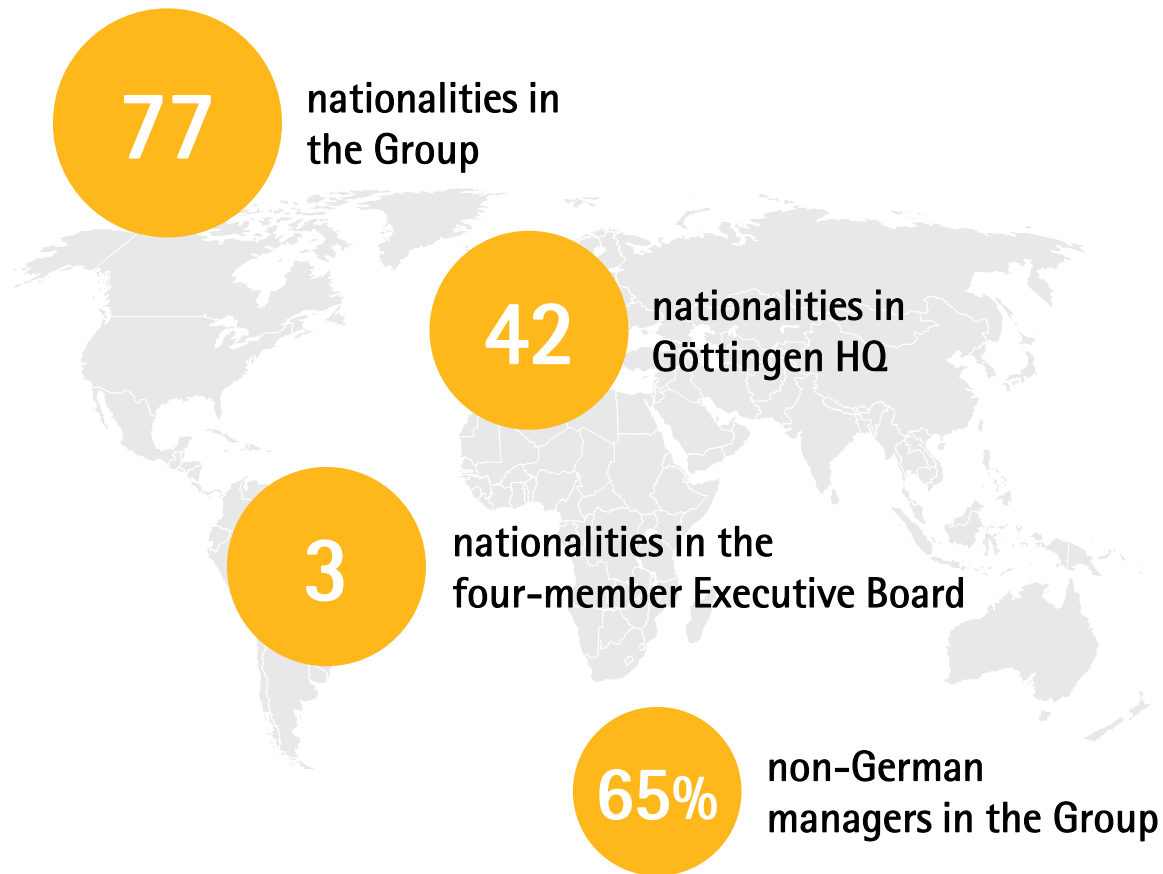




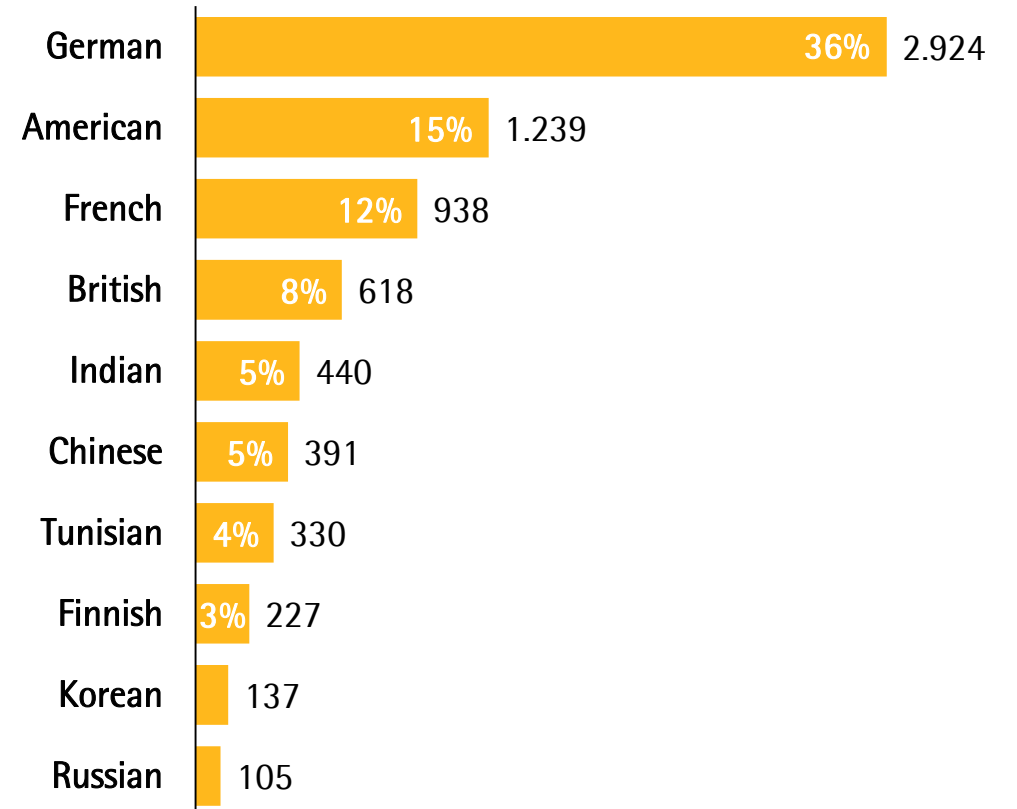
# Diversity Fact Sheet

Sartorius 2018

# Nationality Sartorius unites people from 77 nations



## Top 10 nationalities - Share and number of staff

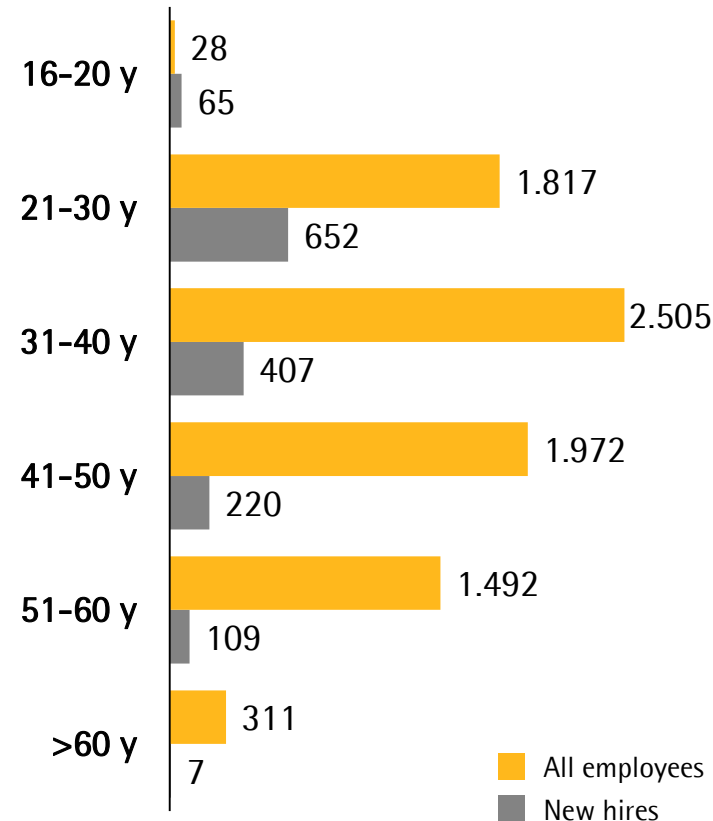


# Age China and India are emerging talent markets

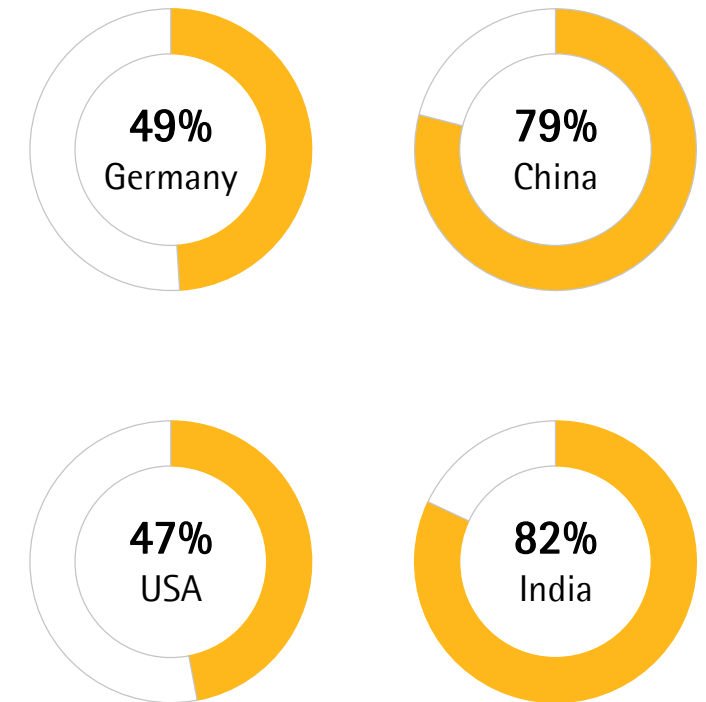
Average age



Number of staff by age

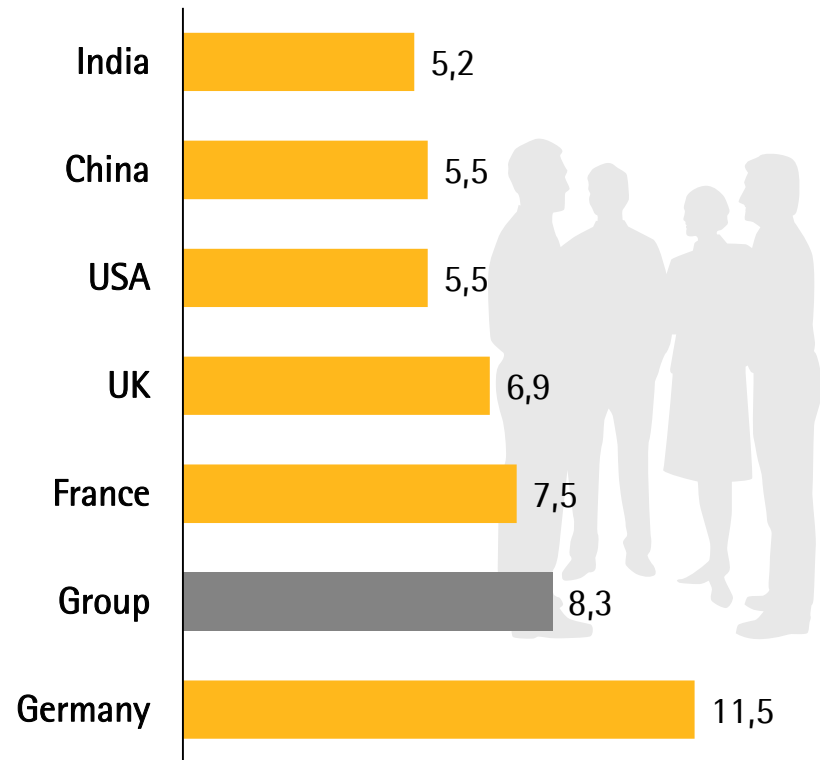


Share of staff younger than 40 yrs

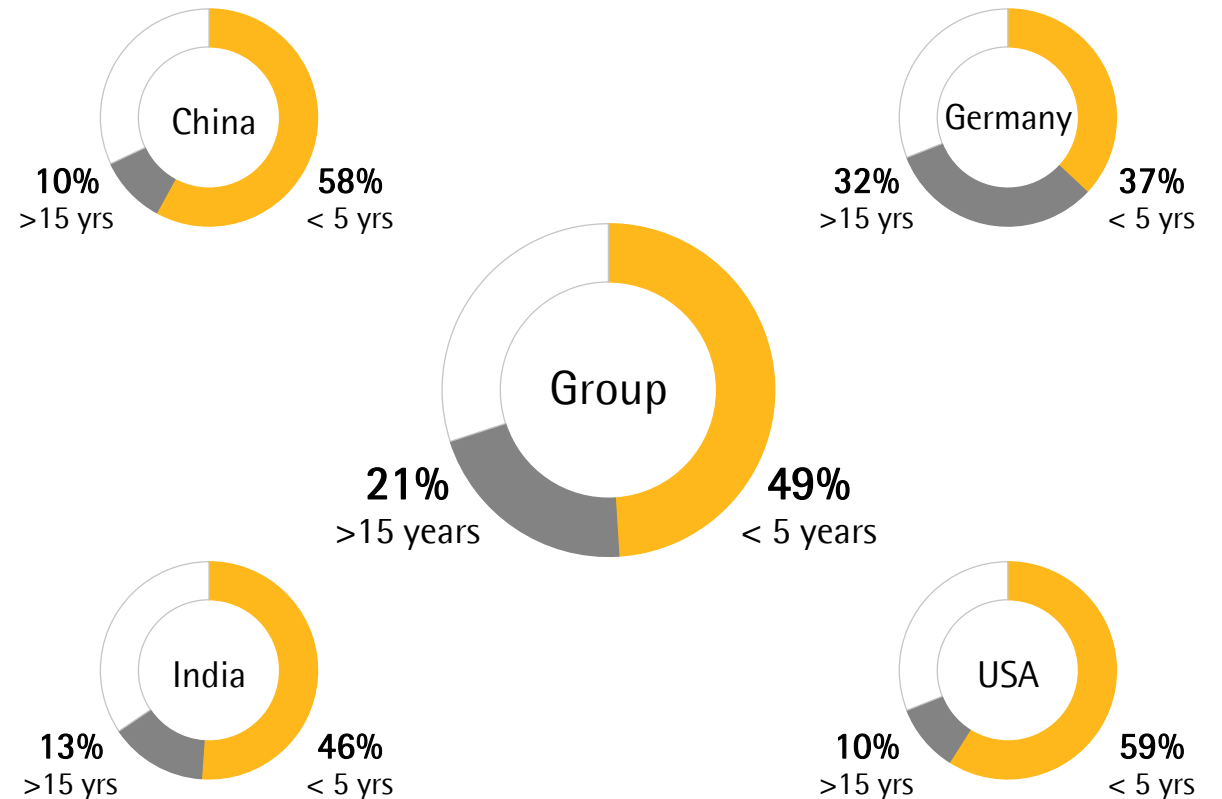


# Tenure Healthy mixture of experience and fresh ideas

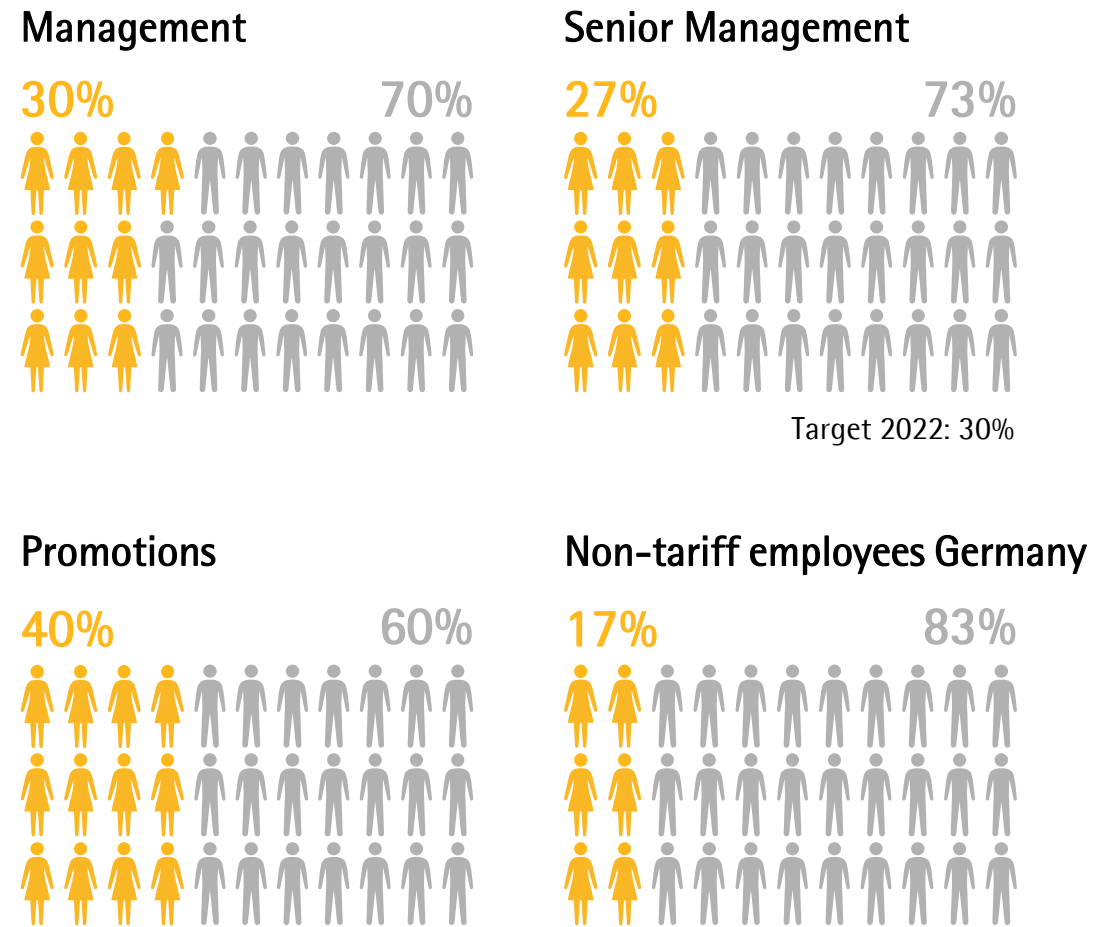
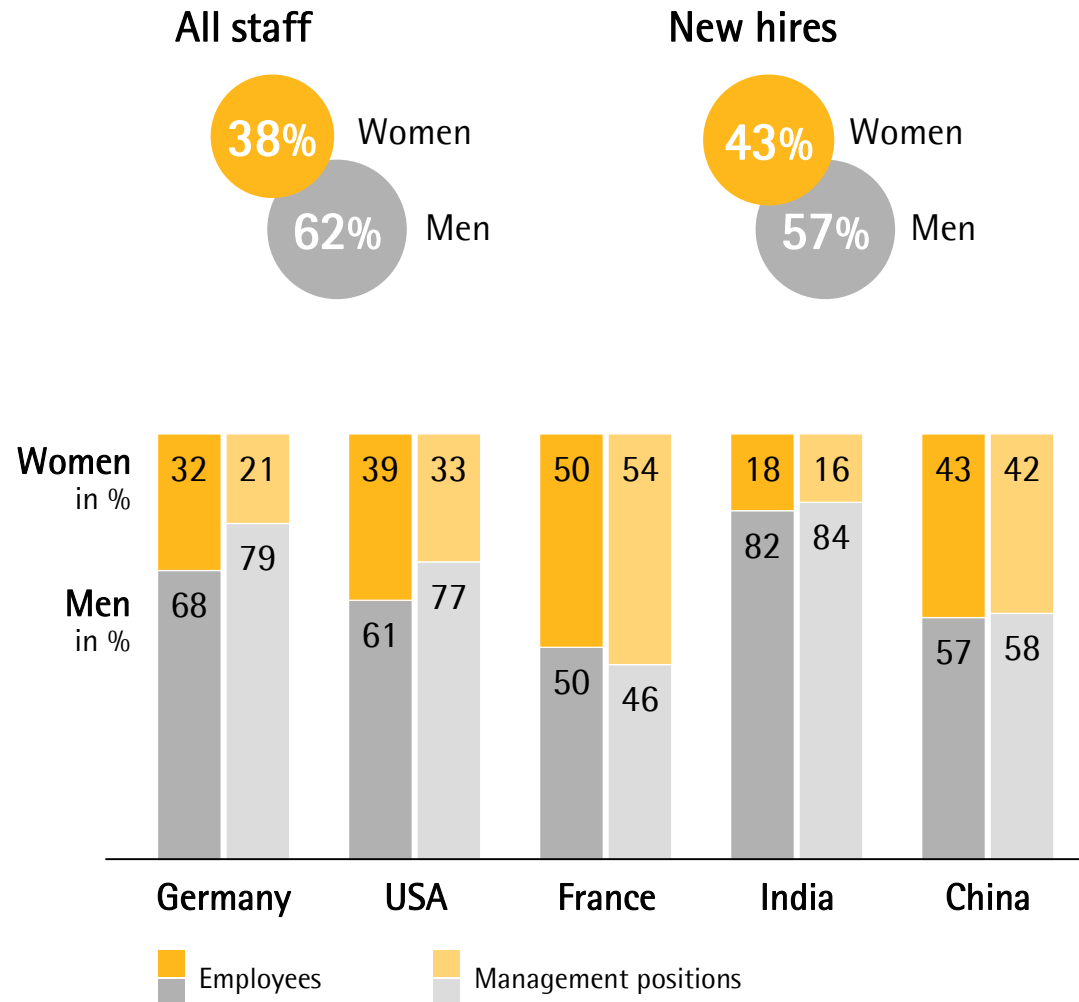
Average tenure in years



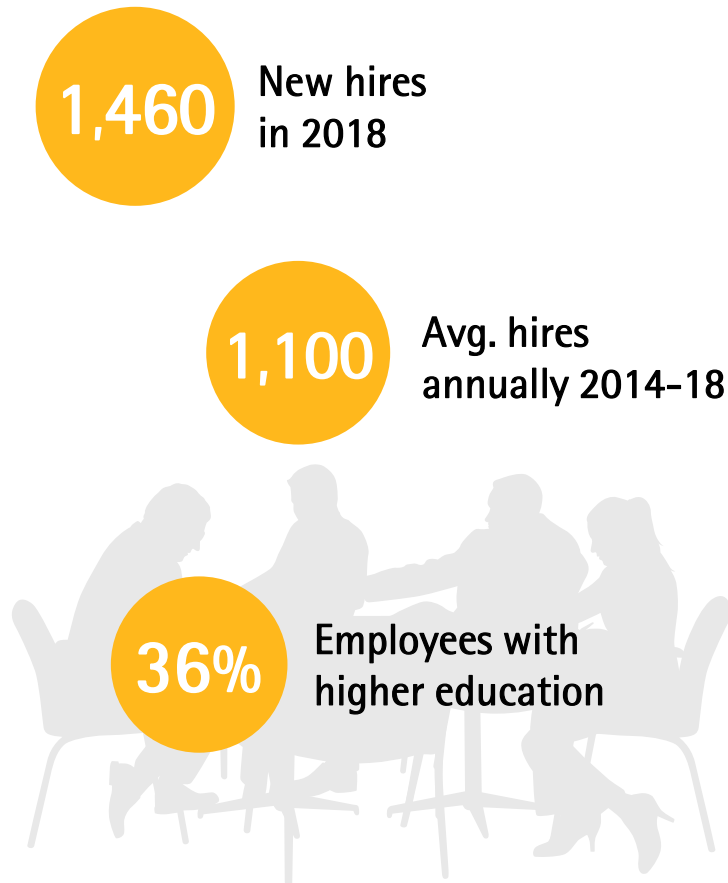
Share of staff with tenure higher than 15 yrs and lower than 5 yrs



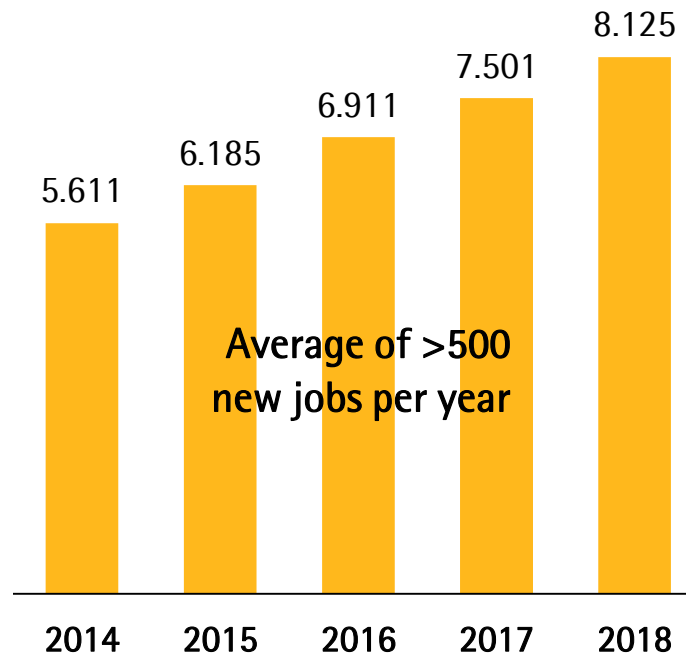
# Gender 43% of new hires are women, 27% women in senior management



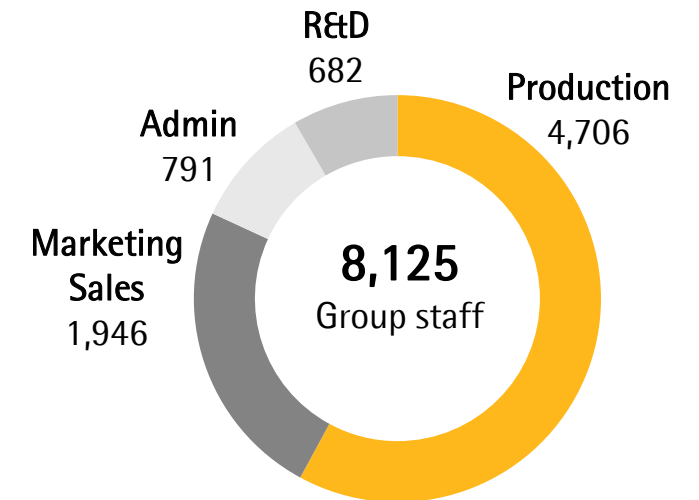
# Welcome 1,460 new people joined Sartorius in 2018



Employee growth



Employees by function



” At Sartorius, diversity is not a question of statistics. It’s a question of culture. We welcome those who live and love diversity because they see it as a chance to learn from each other, to argue constructively and to grow personally and as a team. This can be challenging. But it teaches us to see things in a new perspective in order to find the best solution on behalf of our clients.

