

Simplifying Progress

Diversity Fact Sheet

Sartorius 2023



"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



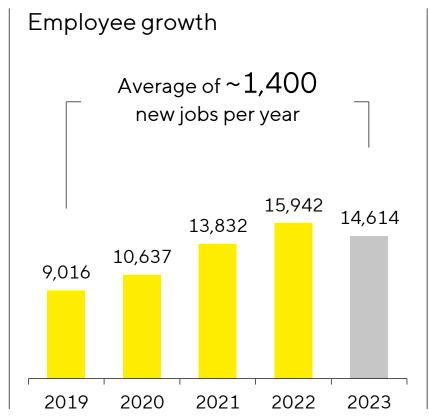
Welcome | 858 New People Joined Sartorius in 2023

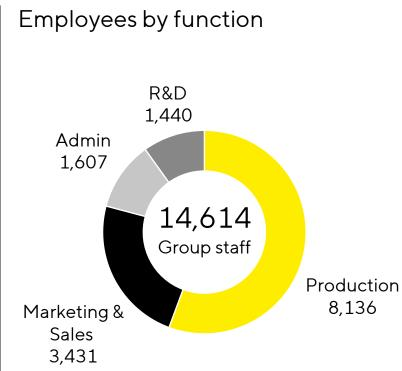
858 | New hires in 2023

New employees from M&A in 2023

2,472 Avg. hires annually 2019-2023

42% Employees with higher education

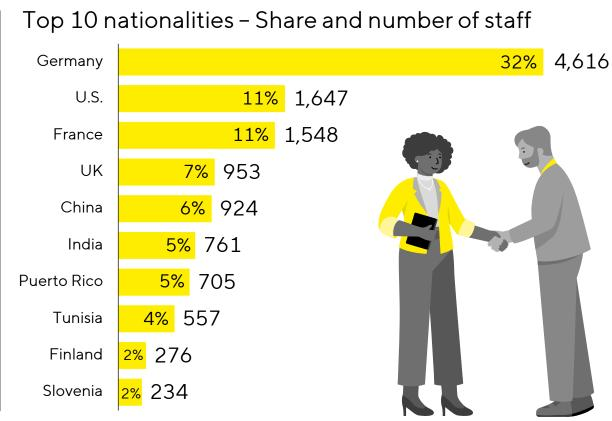






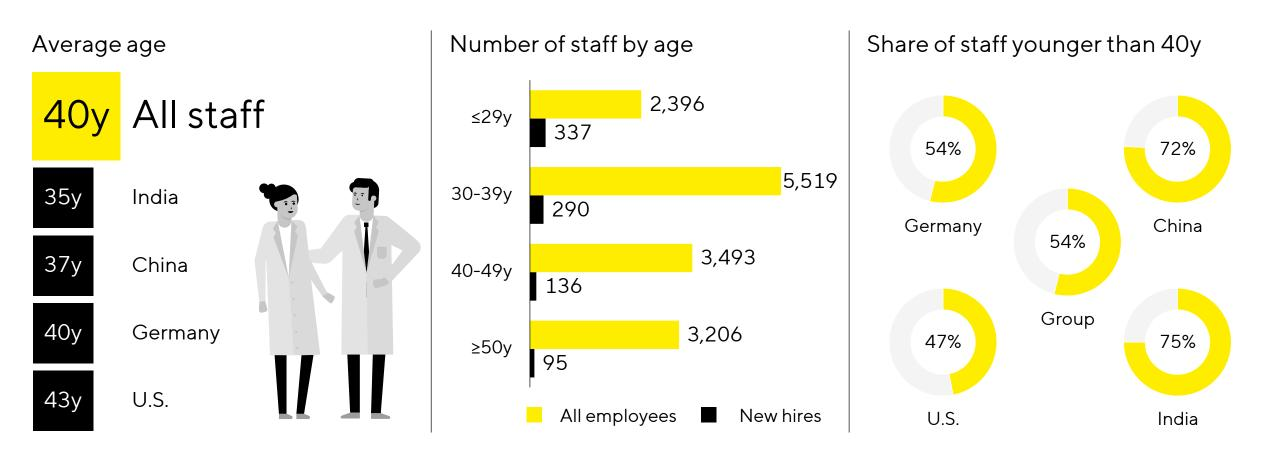
Nationality | Sartorius Unites People from 113 Nations





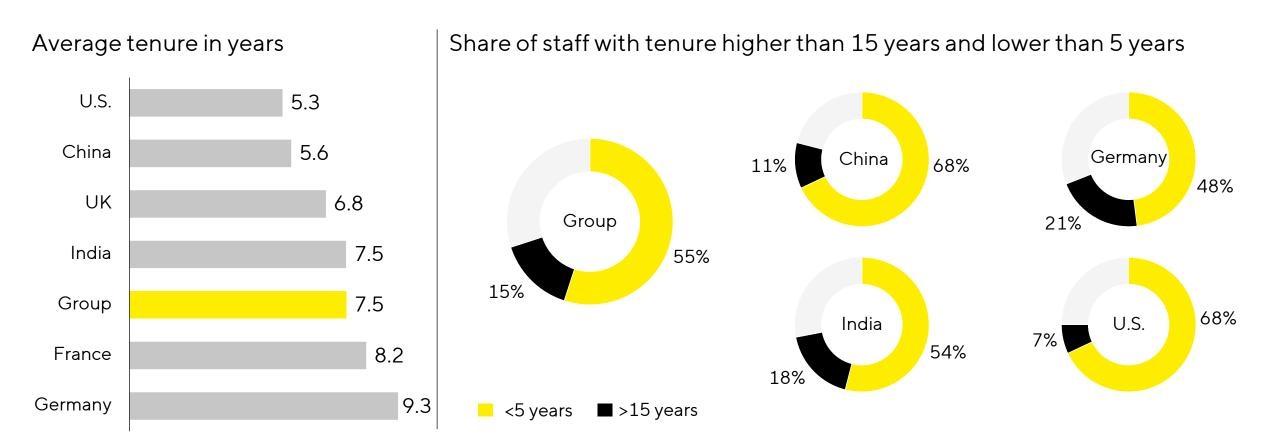


Age | Young Talent and Experienced Professionals Shape Our Workforce



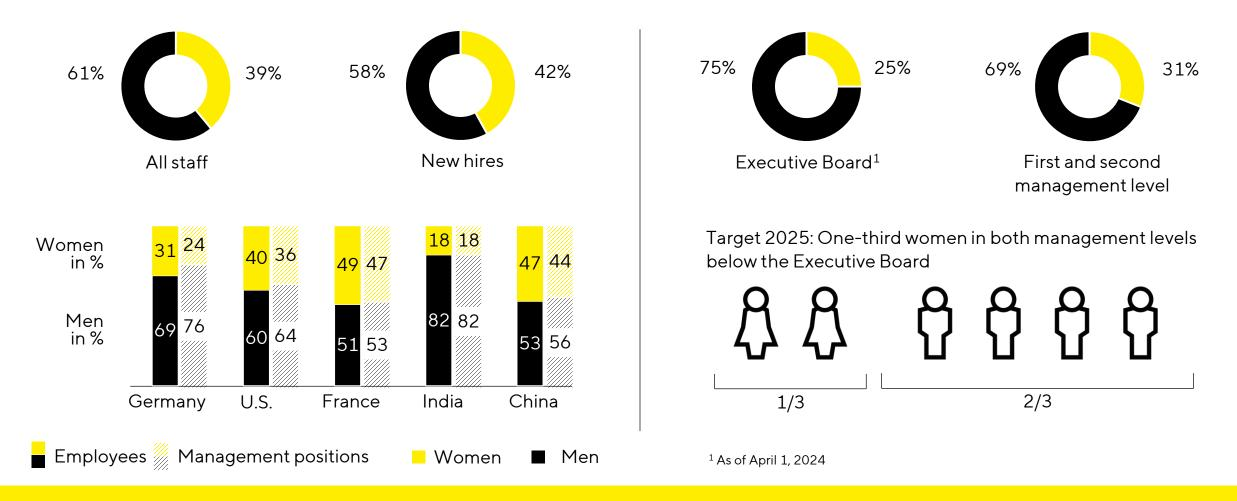


Tenure | Mix of Long-Term Experience and Fresh Ideas





Gender | Structure Headed in the Right Direction





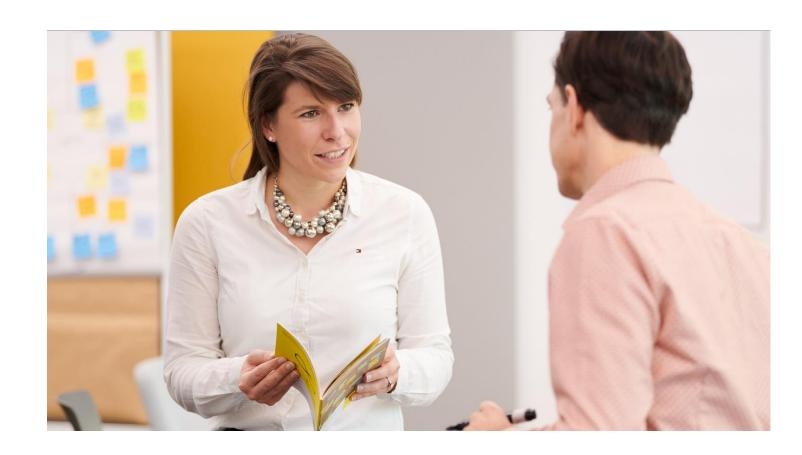
Fostering Exchange and Dialogue for Women

Sartorius Business Women Association

The network serves to connect women across all regions and aims to support gender parity in leadership positions at Sartorius.

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. So far, 800+ employees have completed Emtram's Unconscious Bias Awareness 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.



Partnering in the Name of Diversity and Inclusion













Initiative zur Berufsund Studienwahl

