

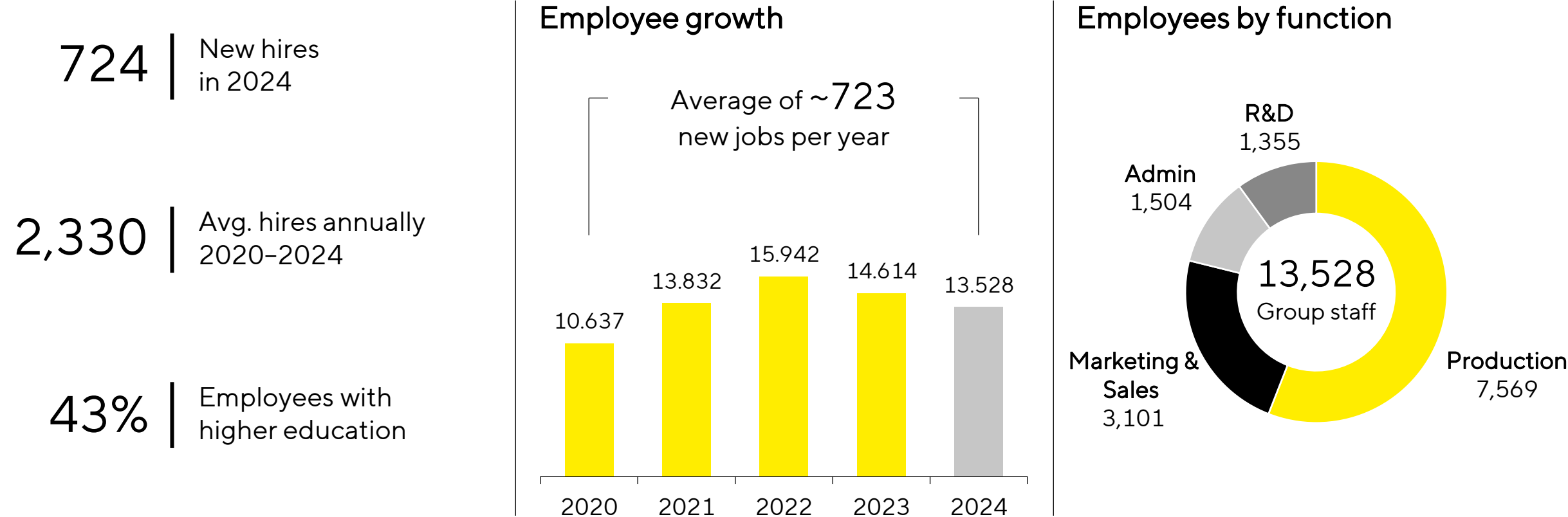
Simplifying Progress

# Diversity Fact Sheet

Sartorius 2024

**SARTORIUS**

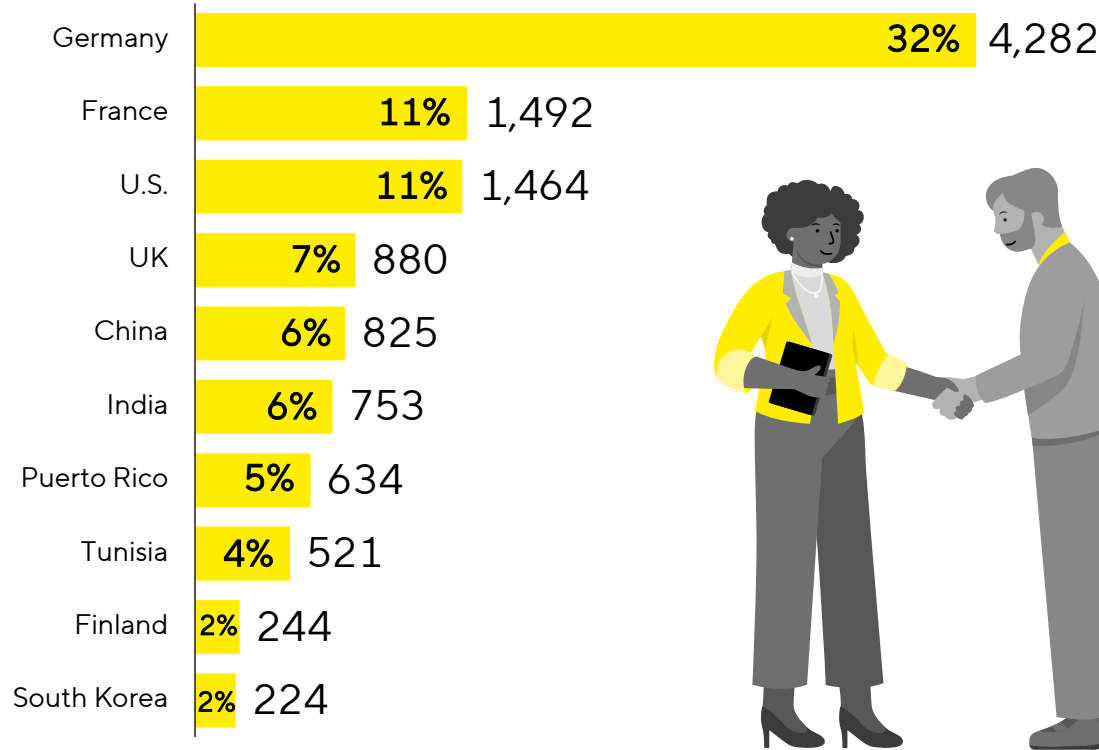
# Welcome | 724 New People Joined Sartorius in 2024



# Nationality | Sartorius Unites People from 112 Nations

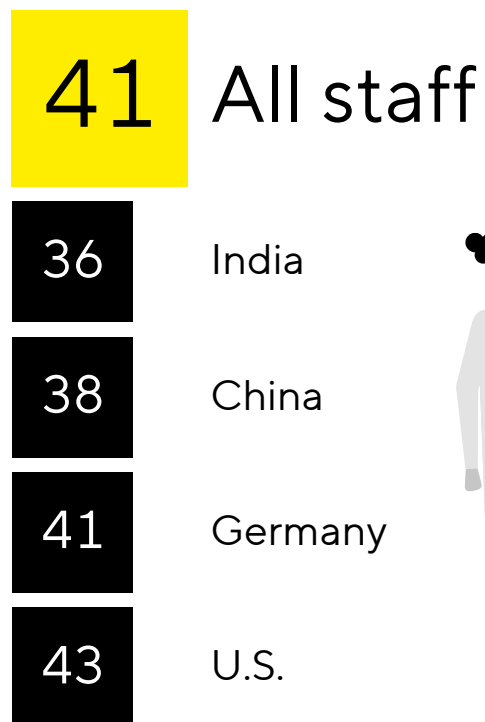


## Top 10 nationalities – Share and number of staff

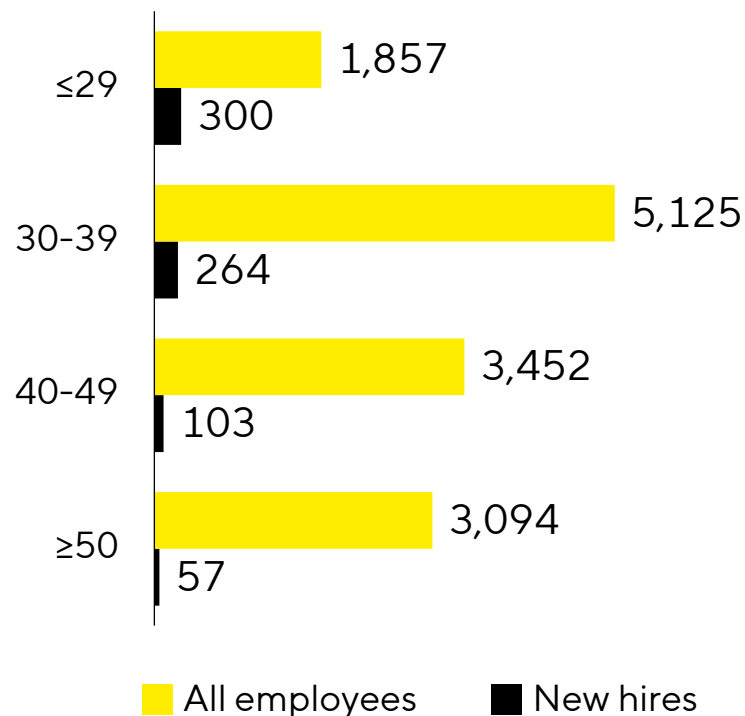


# Age | Young Talent and Experienced Professionals Shape Our Workforce

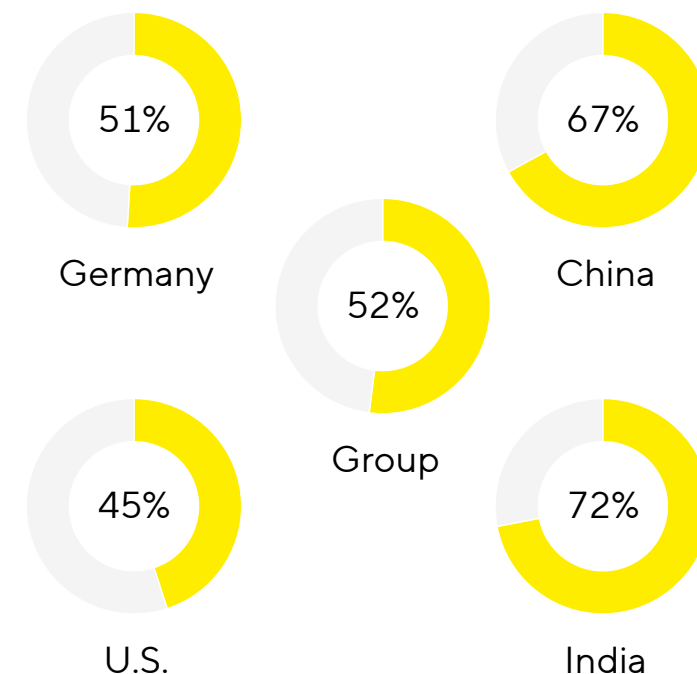
## Average age



## Number of staff by age

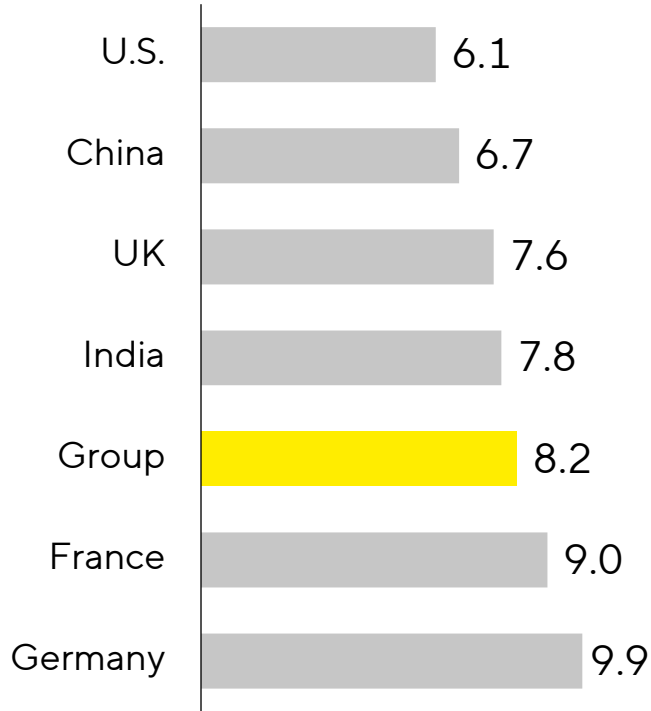


## Share of staff younger than 40y

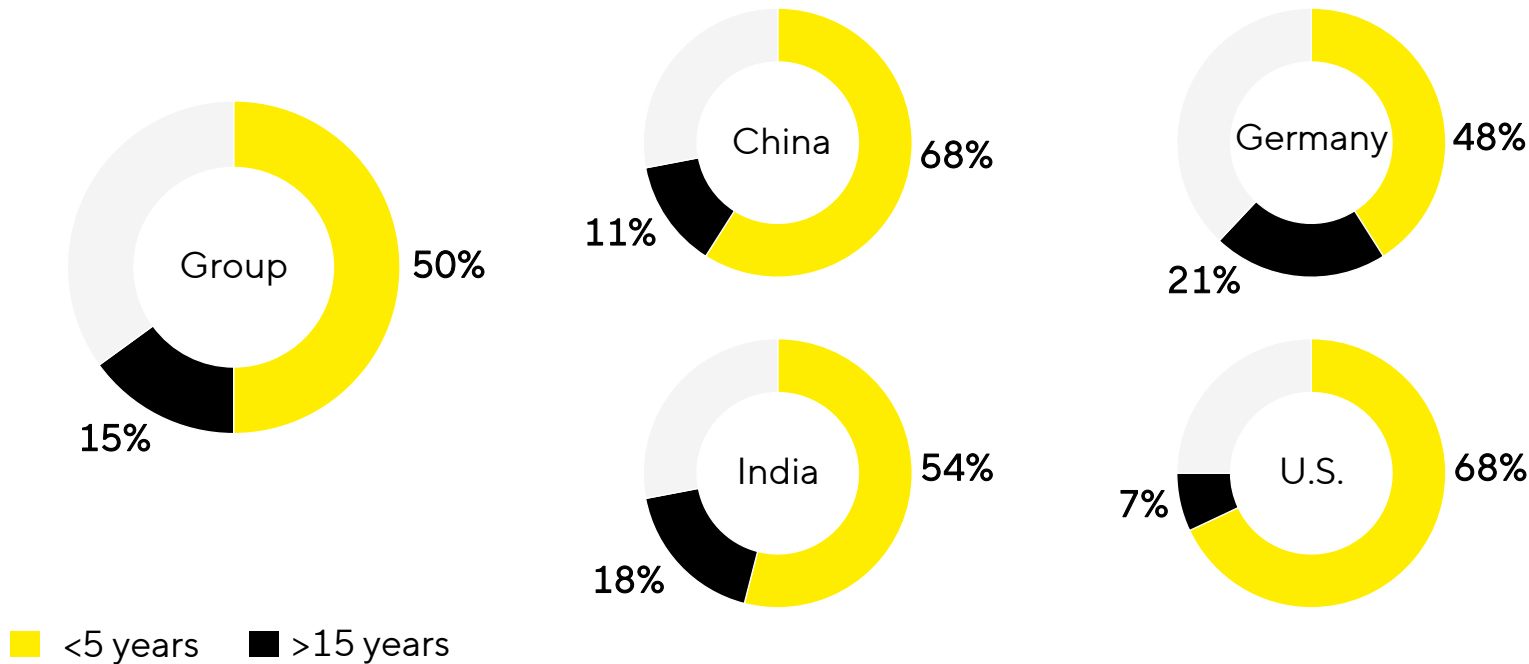


# Tenure | Mix of Long-Term Experience and Fresh Ideas

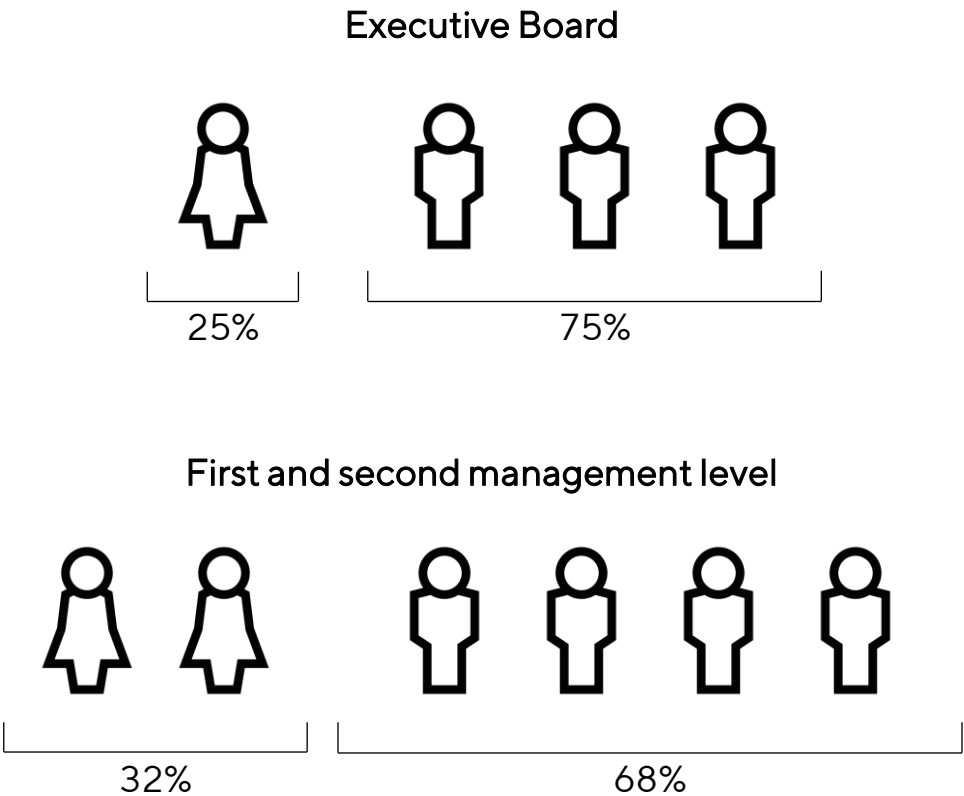
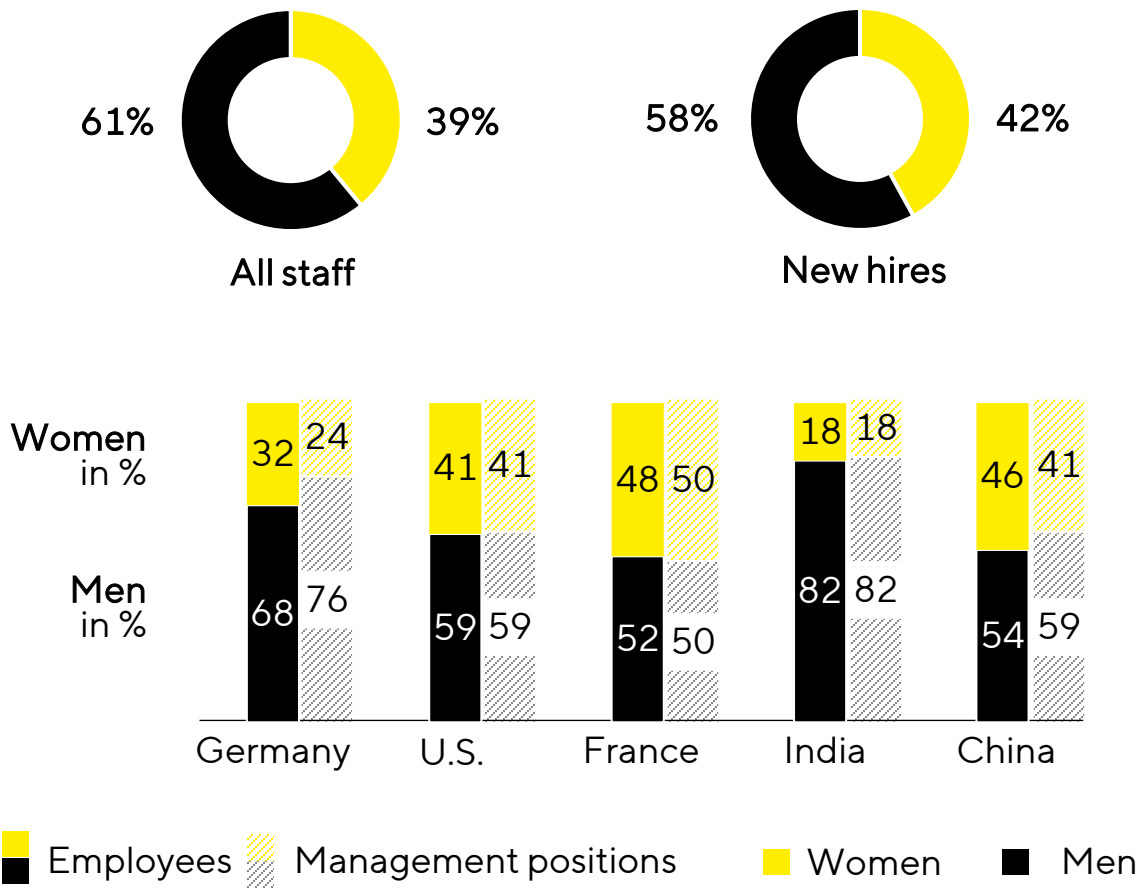
Average tenure in years



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



# 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

