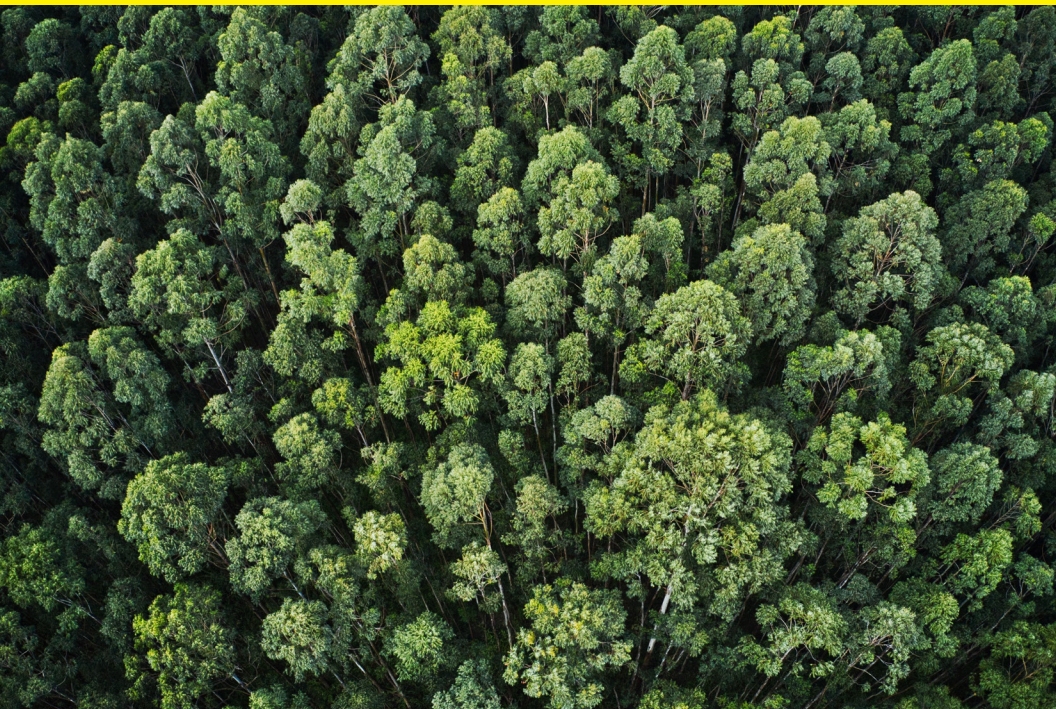




Liquid Handling

Did you know Sartorius uses 100 % renewable electricity at the production site for liquid handling? Did you know our tips, trays and boxes are 100 % recyclable? And our mechanical pipettes are up to 90 % recyclable!



Lab Filtration

Did you know that many of our filtration products (incl. packaging) are highly recyclable?



Together, we create a sustainable future

At Sartorius we aim to provide information, programs and products that support our partners in achieving sustainability goals together

CO₂ Intensity Reduction

Sartorius has set a target to reduce GHG Emission intensity by about 10 % on average per annum by 2030.



Scan the QR code to get more information

Partnership

Let's work together to solve our customer's challenge. Scan QR code and find out who your local Sartorius Account Manager is today!

Together,
We Succeed



Scan the QR code to find your local Sartorius Salesperson