

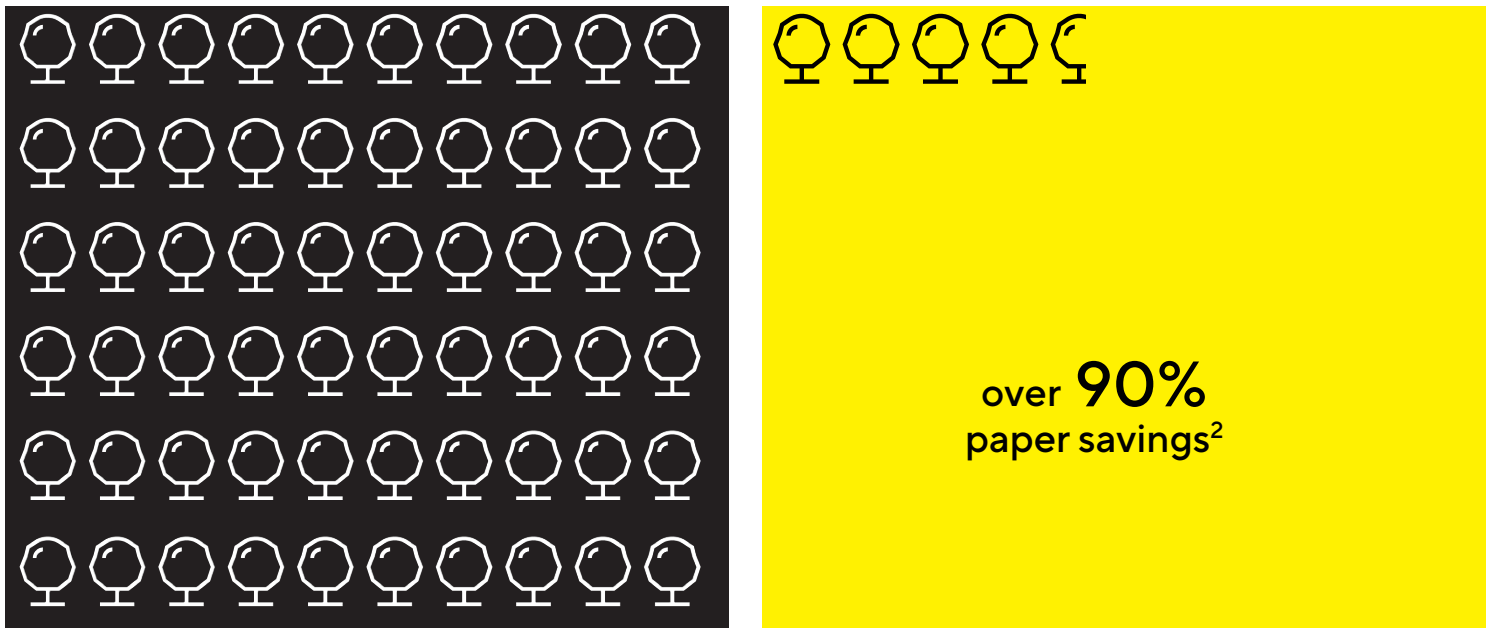
Simplified Instructions

The key information you need to use our lab ultrafiltration and chromatography products safely¹, including:

- Easy-to-follow, visual instructions
- Concise operating requirements
- Reduced environmental impact

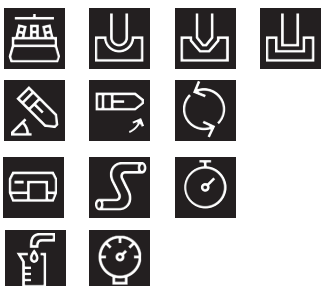
Less Paper Consumption

Number of trees required to print instruction manuals (left) and quick start guides (right):



Intuitive Layout and Symbols

Product Overview	Warnings	Visual Instructions	Specifications



Centrifuge specifications: Rotors, cavity profile (round, conical or flat base)

Centrifuge operating limits: Fixed angle rotor, swing out rotor, maximum RCF

Other specifications: Pumps, tubing, typical processing time

Other operating limits: Maximum sample volume, maximum pressure

¹ Before using our products, please download and read the full instruction manual from www.sartorius.com.

² Based on implementation of quick start guides for all lab ultrafiltration and chromatography products manufactured in the UK. Calculated in May 2023 using internal data (excluding growth) and external paper consumption calculators.