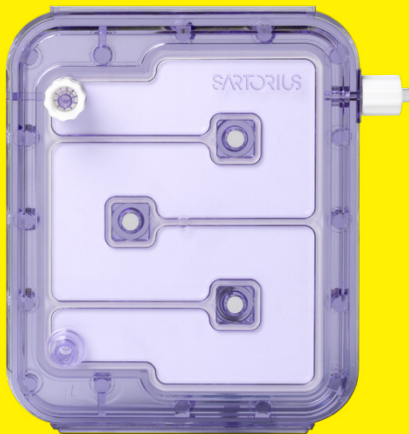


Spot the Difference\*



\* This Vivaflow® SU TFF cassette is even more sustainable.

## Vivaflow® SU Sustainability

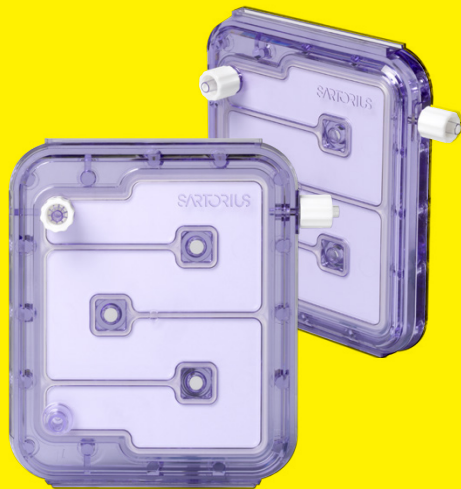
### We're Just Getting Flowing

In our mission to make TFF for the lab more sustainable, we have given Vivaflow® SU an invisible makeover.

The polycarbonate used in manufacturing Vivaflow® SU now comes certified under the International Sustainability and Carbon Certification (ISCC) PLUS program. This ensures that we have replaced a portion of the raw material, previously produced using crude oil, with the same material derived from renewable sources. Although Vivaflow® SU looks and performs the same, this change marks another step Sartorius has taken to enhance circularity and reduce the environmental impact of your research

Learn more about how Vivaflow® SU is promoting sustainable research practices:

[www.sartorius.com/en/company/sustainability/product-sustainability](https://www.sartorius.com/en/company/sustainability/product-sustainability)



# SARTORIUS