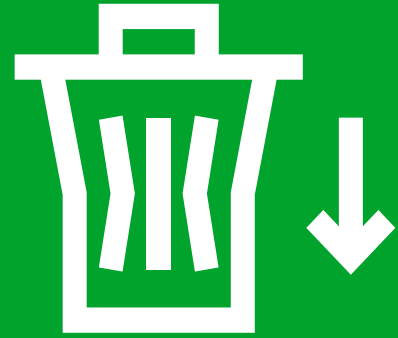


## Smarter Packaging. Smaller Footprint

Reducing waste and boosting sustainability – one product at a time.



Sartorius is committed to minimizing environmental impact in the life science and pharmaceutical industries. Beyond helping labs save energy, water, and cleanroom space, we continuously reduce our own footprint by cutting emissions, waste, and packaging materials through circular economy initiatives

### Packaging Progress Over Time



#### Cubis® II Draft Shield for high-performance balances Cubis® II

- Corrugated cardboard insert for draft shield: recyclable, plastic-free \*except pouches
- All packaging pouches, with the exception of ESD pouches with a minimum of 80 % recyclate



#### Picus® 2

- Packaging made with 100% FSC certified cardboard
- With smart packaging design packaging size reduced by 60%
- Reduced shipping emissions with compact packaging



#### Standard Lab Balances Quintix® Pro

- Plastic pouches from minimum 80% recyclate: less virgin plastic waste
- Cardboard accessory insert: Conventional PS foam parts (compared to BCA and BCE) have been replaced with cardboard inserts that are 25% lighter



#### Lab Water Systems Arium® (without Mini series)

- EPS has been replaced by multi-shock-resistant EPE cushions: more recycling streams available
- Packaging suitable for single-parcel shipping: Pallets are no longer needed.
- Improved protection: fewer damages, fewer returns



### WZB OEM Weigh Cells

- Minimum 10% less total packaging weight overall: lighter shipments, lower emissions
- Approx. 50% lighter PE foam cushion thanks to a new cushion design
- Switch to PE recyclate: less reliance on raw resources
- Bubble wrap sheets and PU foam slipcase replaced by cushioning paper material.

By rethinking packaging, Sartorius helps labs cut waste and shrink their footprint, without compromising product safety. Together, we advance science – sustainably.



### Circular Packaging Benefits

- Less plastic waste in labs
- More recyclable and renewable materials
- Reduced shipping emissions from lighter loads
- Stronger protection = fewer damaged products

## Sartorius Sustainability Initiatives

For further information, visit

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



### Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Göttingen  
Phone +49 551 308 0

### USA

Sartorius Corporation  
3874 Research Park Drive  
Ann Arbor, MI 48108  
Phone +1 734 769 1600

 For further information, visit  
[www.sartorius.com](http://www.sartorius.com)