

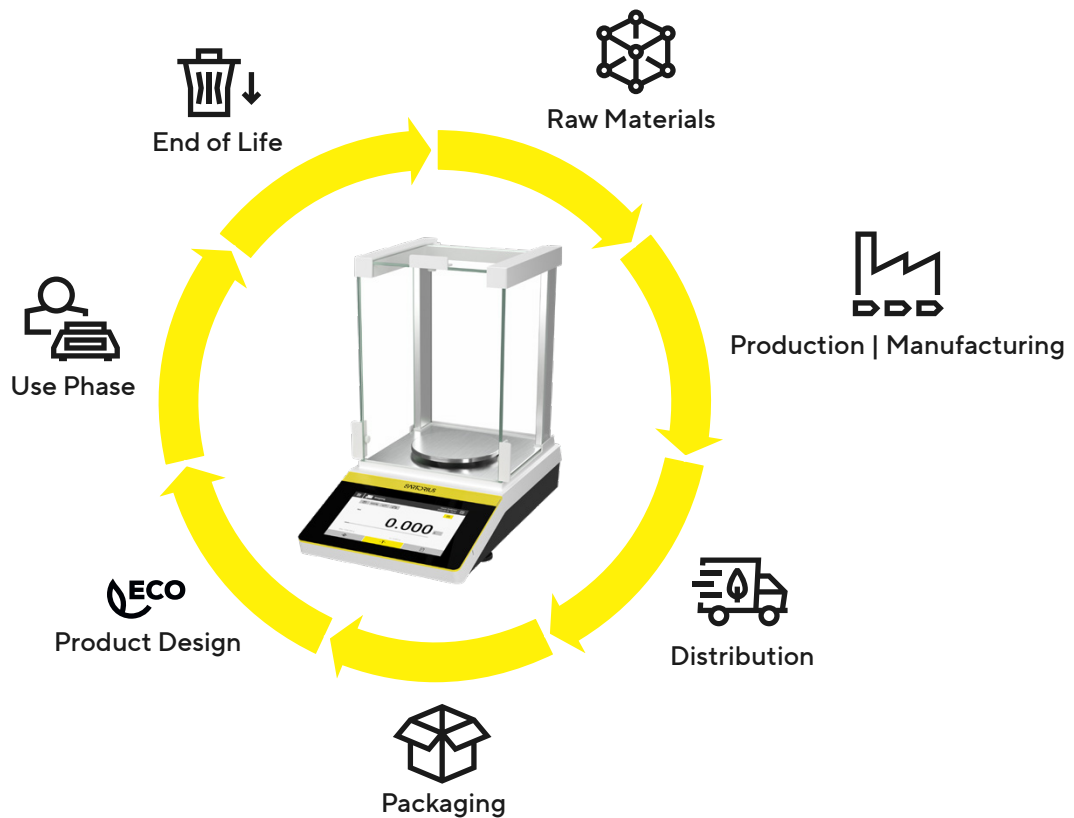
Quintix® Pro – Life Cycle Thinking

Quintix® Pro lab balances embody a significant innovation, with life cycle thinking at the core of their design and development. The early implementation of a screening Life Cycle Assessment (LCA) has been instrumental in creating a product that stands out for its technical excellence and commitment to environmental sustainability. This proactive strategy has enabled us to address potential environmental impacts from the start, positioning Quintix® Pro as a great choice in eco-friendly innovation within our industry.

Our dedication to sustainability is integral to the Quintix® Pro. We've chosen materials like aluminum with recycled content

to ensure high technical performance while embracing renewable energy in our production. We've also phased out non-recyclable packaging, reinforcing our commitment to waste reduction and the circular economy.

The design of Quintix® Pro reflects our holistic approach to product life. Energy efficiency, serviceability, and end-of-life recyclability are deliberate design choices that prolong the product's lifespan and minimize its environmental footprint. By considering the full life cycle of Quintix® Pro, sustainability remains a central focus from inception to disposal.



1. Raw Materials

The Quintix® Pro line showcases our commitment to sustainability, utilizing aluminum with a substantial recycled content sourced within Europe, reducing our carbon

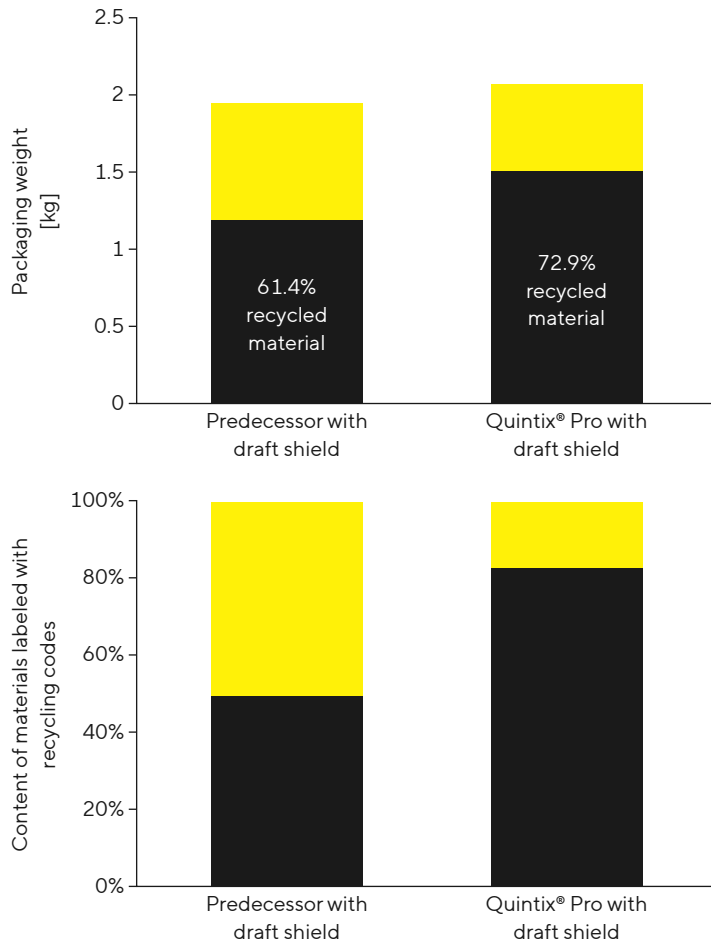
footprint. Our sourcing strategy prioritizes recycled materials to diminish environmental impact. With 20% recycled content in high-volume parts like the carrier plate, we conserve resources and support the circular economy.

2. Production | Manufacturing

At the heart of Quintix® Pro production is our Göttingen, Germany facility, where we manufacture a significant portion of the Quintix® Pro family using 100% renewable electricity. We boast a 91% recycling rate and are on track for zero-waste production by 2030. Every piece of aluminum scrap is recycled, ensuring full material utilization. Our operations, certified by the ISO 14001 Environmental Management System, are a testament to our commitment to environmental excellence.

3. Distribution

Quintix® Pro balances reach their destinations through optimized distribution channels, favoring sea transport over air to enhance environmental efficiency. Strategic distribution hubs and consolidated shipments further reduce our carbon footprint. This logistical approach ensures timely delivery and demonstrates our commitment to sustainable transportation practices.

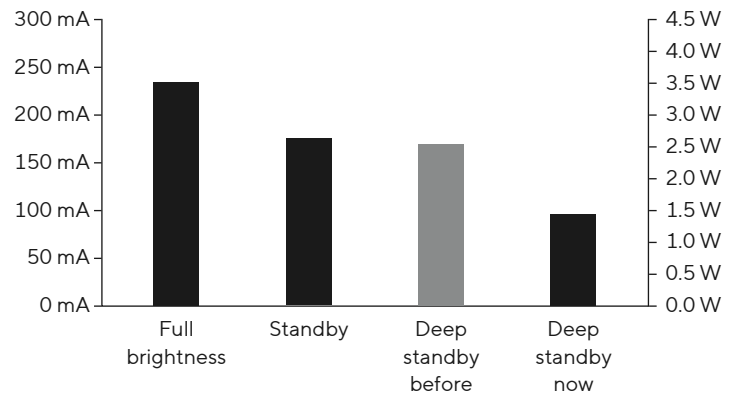


4. Packaging

We have significantly reduced our reliance on new plastic by using protective bags made from 80% post-consumer recycled material. Our cardboard boxes contain 60% to 100% recycled material, and for the Quintix® Pro with draft shield, we've increased the recycled content in packaging to over 70%. Most packaging materials are labeled with recycling codes to streamline the recycling process.

5. Product Design

Energy efficiency is a key design principle of the Quintix® Pro, featuring low power electronics and intelligent backlighting that dims to save energy and extend product life. The deep-standby mode cuts energy use by more than half for an average 8-hour, 5-day-a-week operation. We've also digitized and integrated the manual and operation instructions into the Quintix® Pro user interface, reducing the printed version by 30% to lessen our environmental impact.



6. Use Phase

The balance is designed with replaceable and repairable parts, promoting longevity and reducing the need for frequent replacements. Our service model includes a network of local service hubs close to our customers, reducing the need for extensive logistics and further cutting our carbon footprint.

7. End of Life

As a Quintix® Pro balance nears the end of its usable life, its design facilitates easy disassembly for recycling. We avoid using restrictive technologies, like adhesives, to simplify the recycling process. This design philosophy is consistent with our goal of creating products with their end-of-life disposal in mind.