

Light Pack

The Responsible Tip Box



Eco-conscious



Reduce your lab's
environmental footprint



Forward – thinking

We've created the responsible pipette tip box: a sturdy cardboard design packed with environmental benefits – without losing performance.

- Compared to the single tray plastic rack, the plastic content of the box has been reduced by 74%
- The cardboard of the Light Pack is FSC certified
- The tray is made from recyclable plastic
- The window is made from acetate cellulose
- Awarded with the My Green Lab ACT Ecolabel

All this means that you have a pipette tip box that delivers measurable impact. When you're done, the box folds flat for efficient stacking and easy integration into established recycling workflows.



my green lab.
ACT Ecolabel



reddot winner 2026

>>

Ordering Information

Volume	Purity Level	Tips/Units	Order Code
200µl	▪ Free of DNase, RNase, Human DNA and Endotoxins	▪ Optifit ▪ 10 x 96	LH-G790200
200µl	▪ Free of DNase, RNase, Human DNA and Endotoxins ▪ Pre-sterilized	▪ Optifit ▪ 10 x 96	LH-G790201
200µl	▪ Free of DNase, RNase, Human DNA and Endotoxins ▪ Pre-sterilized	▪ Safetyspace ▪ 10 x 96	LH-G790201F
120µl	▪ Free of DNase, RNase, Human DNA and Endotoxins ▪ Pre-sterilized	▪ Safetyspace ▪ 10 x 96	LH-G790101F

Optilid
Hinged lid which can be opened and closed with one hand

Foldable
At end of use, reducing waste volume by up to 50%

Cardboard box with Plastic Tray
Plastic reduction by 74%

Cellulose Acetate Window
to see the number of tips in the box

Safety Corner and Seamless Bottom
Cleanliness and reduction in contamination risk

Robust and Rigid Design
For stability during tip picking

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