



Save Up to 10 Weeks

SARTORIUS



Takeone[®] - The Time-Saving

Start immediately

Bypass parts washing, equipment preparation and assembly with Takeone[®] and break bottlenecks to your process.

Reduce costs

Take the right collection vessel for your individual samples, reduce product loss from over-sampling and minimize your costs.

Hybrid Single | Multi-Use Systems

3 min



6 min



4 min



3 min

Unpack | Connect

Parts Washing

Equipment Prep

Aseptic Sampling System

Sample safely

Every sampling line of every Takeone® device is leak tested after assembly.

Takeone®

3 min



× 1,500 Devices/Year

Unpack | Connect

Man Weeks Used



2 weeks



3 min



= 19 min



× 1,500 Devices/Year

Assembly

Disassembly

Total Time Needed



12 weeks

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For further contacts, visit
www.sartorius.com

Specifications subject to change without notice.

© 2021 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France

Publication No.: SP-1509-e | Order No.: 85037-551-69 | Status: 06 | 02 | 2021