

Quickstart Guide

Software Update BioPAT® Trace | Multi Trace



2908289-000-00



SARTORIUS

How to Update Software on BioPAT® Trace and Multi Trace

1. Unpack the zip file.
2. Copy the folder “updater” to an USB flash drive.
3. Switch off the BioPAT® Trace.
4. Insert the USB flash drive into one of the USB ports on the back site of the BioPAT® Trace.
5. Switch on the BioPAT® Trace.
6. During booting (right LED is white) the software will be updated.
7. Wait until the right LED turns red (no PC connection), which takes about 30-40 seconds.
8. Switch off the BioPAT® Trace.
9. Remove the USB flash drive.
10. Switch on the BioPAT® Trace again.
11. Connect the BioPAT® Trace to the PC according to the manual.
12. Download and install the trace_mon Software on your PC from the internal homepage of the BioPAT® Trace (connection via Internet browser und IP address 192.168.120.231 [default setting]).
13. After successful connection you see the connection information on the left top corner.
14. Open the folder information (last item on the right site).
15. Firmware version (Netus) and trace_mon version will show the actual versions and must match.

Note

If the software is **not** updated: Repeat the process starting from point 3!
Sometimes it depends on the data transfer speed of the USB flash drive!

© 2022

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

www.sartorius.com

Version: 02 | 2022
MB | DIR: 2908289-000-00