

MET ONE 3400+ Software Package Release Notes

Version: 1.0.38

Production Release Date: October 2023

Includes the following new features:

1. Added date and time to login screen.
2. Backup and restore of SOPs.
3. Added IP address to login screen.

Includes updates and fixes for the following situations:

1. Disconnecting from a network would suddenly cause the instrument to lock up.
2. Event log entries did not display the correct value when translated.
3. Negative time could appear on zero count test during calibration.
4. Audit Trail Excel Export did not reflect time zone changes.
5. Error could occur on sample export.
6. Network drive password and username inputs did not allow special characters.
7. Samples appeared in reverse order on routine monitoring report.
8. Error occurred with event log filtering when set to local language.
9. SOPs did not display consistently across different users.
10. Event log setting changes were not recorded in the audit trail.
11. Sound setting changes were not recorded in the audit trail.
12. Adding or removing a network printer was not recorded in the audit trail.
13. Some LDAP entries were not recorded in the audit trail.
14. Changes to wired network settings were not recorded in the audit trail.
15. Battery warnings were not displayed in the virtual ticker tape.
16. User could not turn on the pump during calibration.
17. User could not update the NTP server address.
18. User could not save SOPs or Buildings.
19. Updated translations for *Audit Export to PDF*.
20. Updated PDF generation.
21. Updated the calibration certificate according to A2LA standards to include correct logo and address for each site.