

# Revision History

<b>Filename</b>	Revision History IoT Client
<b>Author</b>	Syed Ghouse
<b>Storage place of valid version</b>	DHF Number 04207_00794 V1.0
<b>Corresponds to IoT Client Version</b>	IoT Client 2.8.4.276
<b>Date</b>	2024-11-28

# Contents

---

IoT Client 2.8 SP1 .....	4
Bug fixes.....	4
New Features.....	4
IoT Client 2.8.....	4
Bug fixes.....	4
New Features.....	4
IoT Client 2.7.....	5
Bug fixes.....	5
New Features.....	5
IoT Client 2.6.....	6
Bug fixes.....	6
New Features.....	6
IoT Client 2.5 SP1 .....	6
Bug fixes.....	6
New Features.....	6
IoT Client 2.5.....	7
Bug fixes.....	7
New Features.....	7
IoT Client 2.4.....	7
Bug fixes.....	7
New Features.....	7
IoT Client 2.3.....	7
Bug fixes.....	7
New Features.....	8
IoT Client 2.2.....	8
Bug fixes.....	8
New Features.....	8
IoT Client 2.1 SP1 .....	9
Bug fixes.....	9
New Features.....	9
IoT Client 2.1.....	9
Bug fixes.....	9
New Features.....	9

IoT Client 2.0 SP2 .....	10
Bug fixes.....	10
New Features.....	10
IoT Client 2.0 SP1 .....	10
Bug Fixes .....	10
New Features.....	11
IoT Client 2.0.....	11
Bug Fixes .....	11
New Features.....	11

## IoT Client 2.8 SP1

---

- Version **2.8.4.276**

### Bug fixes

ID	Title
<b>42100</b>	Impossible to register the instrument after installing IoT Client 2.8 (Installer improvement: add check for prerequisites)

### New Features

ID	Title
<b>42391</b>	Improve installer error feedback.

## IoT Client 2.8

---

- Version **2.8.2.272**

### Bug fixes

ID	Title
<b>41741</b>	IoT Client 2.8 - Upgrade installer removing Misc folder resulting in erroneous messages on restart of the IoT Client after upgrade.

### New Features

ID	Title
<b>34138</b>	IoT Client 2.8 - Remove Pairing functionalities
<b>41292</b>	IoT Client 2.8 - Cloud Registration (over new server)
<b>41302</b>	IoT Client 2.8 - Thingworx connection removal from solution.
<b>41304</b>	IoT Client 2.8 - Update Target framework.
<b>41305</b>	IoT Client 2.8 - Update IoT Client Installer.
<b>41490</b>	IoT Client 2.8 - Remove location from Registration Form and Config Tool
<b>41573</b>	IoT Client 2.8 - Remove Webcam URL
<b>41603</b>	IoT Client 2.8 - Provide more information if connection fails
<b>41623</b>	IoT Client 2.8 - Remove "Remote Update"
<b>41719</b>	IoT Client 2.8 - Remove link to Introspect manual

## IoT Client 2.7

- Version **2.7.5.254**

### Bug fixes

ID	Title
<b>33672</b>	No charts are created, but system was running all the time.
<b>34536</b>	IoT Client to accept and parse 0 and negative Introspect key values
<b>35105</b>	[CyberSecurity] Switch to Newton.Json 13.0.1
<b>35181</b>	IoT Client fails to parse log files when any custom SSL certificate is used to register instruments
<b>35348</b>	No live data visible in MissionControl with IoT Client 2.6
<b>35560</b>	UpdateModule is not deleting Updater_ReadyForUpdate.txt file & files of Updater folder after successfully package update
<b>35562</b>	Package is not getting downloaded when Software Update is triggered by Configuration Management
<b>37013</b>	CompressionRatioThreshold to be changed for IoT 2.7 for log file parsing
<b>37014</b>	When invalid SSL certificate is used, IoT Client 2.7 gets frozen while instrument registration and in configuration tool
<b>37016</b>	IoT Client Freezes when registering a new instrument via configuration tool by changing the size of an instrument
<b>37031</b>	IoT Client stops parsing when Software entry in registry is missing information
<b>37119</b>	Parsing performance dropped substantially to very slow parsing
<b>37130</b>	Serial number check missing in IoT Client for the hardware file
<b>37187</b>	IoT-Client is sending property update Type as "0" to IoT Hub instead it should send Type as "Alias"
<b>37125</b>	Release Notes: Open release notes when user clicks on hyperlink in About tab Configuration Tool
<b>37307</b>	Parse Interval change after Remote update of IoT Client fails
<b>37321</b>	Instrument not created in Configuration management when registered using IoT 2.7 client
<b>36897</b>	No additional data uploaded, only possible to see logfiles older than 1 year
<b>37359</b>	Include run name in live message
<b>37364</b>	Restarting of IoT Client in Configuration tool freezes for Universal platform

### New Features

ID	Title
<b>9691</b>	[Fluent] Track environmental data from T-GIO2
<b>29090</b>	IoT Client parses logfiles for Introspect and encounters Parse errors
<b>33504</b>	Reduce log entries for LiveData
<b>33587</b>	[Service Cockpit] Create interface to Control Application for Alias handling
<b>33592</b>	[Service Cockpit] Collect operating system information
<b>34441</b>	Upgrade .Net version, build system and used packages
<b>34924</b>	[Raptor] Parser: Tips are already Mounted
<b>35390</b>	Upgrade setup failed with IoT Client 2.6
<b>35576</b>	Sonar-cube Issues
<b>35958</b>	Update ThingWorx SDK to fix known vulnerabilities
<b>36502</b>	IoT Client does not collect hardware info file.

## IoT Client 2.6

---

- Version **2.6.2.167**

### Bug fixes

ID	Title
31083	IoTClient connection timeout
29528	Universal IoT Client Field validation on SetupConfig File
32547	Encryption of IConnectionSettings doesn't work
31685	Fix SonarQube issues
31817	Instrument state contains "idle" and "Idle" (capital I)
34426	IoT Hub message sending throws null reference exception
34885	Installation of IoTClient fails

### New Features

ID	Title
32391	Upgrade Tecan.MAP.Services.Logging to support .NETCoreApp3
32556	Change registration dialog to reflect the Freemium signup
30023	Preparation to Support new hardware Device on Fluent
30960	Switch IoTClient to .NET Core / .NET Standard
30961	Transition to TBS 6.2
30962	Transition to GIT
32827	Change Link to Manual to Tecan URL
33592	[Service Cockpit] Collect operating system information
33594	[Service Cockpit] Send move counter to IoT Hub
33595	[Service Cockpit] Send QC data to IoT Hub
33596	[Service Cockpit] Send HW data to IoT Hub
33597	[Service Cockpit] Send SW data to IoT Hub
33612	[Service Cockpit] Send initialization cycles to IoT Hub
33635	[Service Cockpit] Upload config files to IoT Hub
31508	Harmonize versions of referenced packages

## IoT Client 2.5 SP1

---

- Version **2.5.1.32637**

### Bug fixes

ID	Title
30866	Scan of too long serial number not working

### New Features

None

## IoT Client 2.5

---

- Version **2.5.0.31826**

### Bug fixes

ID	Title
<b>27894</b>	Fixed an issue where the Timezone was only set once during registration.
<b>27764</b>	Fixed an issue where the Pick&Place errors were showing up as MCA errors in the dashboard
<b>28002</b>	Fixed an issue where the RGA errors were not displayed in Introspect

### New Features

ID	Title
<b>27670</b>	The IoT Client now supports parsing of environmental data
<b>27671</b>	The IoT Client now parses deck check errors
<b>27191</b>	the IoT Client delivers a generic parser and interface that allows to parse available properties for Introspect without any platform specific implementation.
<b>28984</b>	Sample Status Rule

## IoT Client 2.4

---

- Version **2.4.1.30960**

### Bug fixes

ID	Title
<b>None</b>	None

### New Features

ID	Title
<b>25420</b>	IoT client now supports the Mission Control

## IoT Client 2.3

---

- Version **2.3.1.30399**

### Bug fixes

ID	Title
<b>24439</b>	Fixed an issue where the “ <i>FluentControl Preinstaller</i> ” was shown as only “ <i>FluentControl</i> ” in the Configuration Management.
<b>25017</b>	Fixed an issue where too much information of a barcode error was parsed by the IoT Client.
<b>25762</b>	Fixed an issue where log files that were being archived by the AuditArch were not parsed.
<b>25778</b>	Fixed an issue where the IoT Client would endlessly restart when parsing log files with FRIDA measurements.

<b>25807</b>	Fixed an issue where the IoT Client applied incorrect parsing rules for FluentControl version 3.0 and above.
<b>25955</b>	Fixed an issue where a manually replaced SSL certificate was removed during the upgrade process.
<b>26275</b>	Fixed an issue where the CPU usage of the IoT Client was too high on systems with registered EVO instruments and no running control application.
<b>26287</b>	Fixed an issue where custom SLL certificates could not be set in the configuration tool.
<b>26346</b>	Fixed an issue where the IoT Client finalizes the registration of an instrument while an invalid custom SSL certificate was selected.

## New Features

ID	Title
<b>25419</b>	The IoT Client now supports the registration and historical data parsing from D300e devices.

## IoT Client 2.2

- Version **2.2.2.29404**

## Bug fixes

ID	Title
<b>21748</b>	Fixed an issue where only the first run after the startup of FluentControl or EVOware has the Software Name set.
<b>23186</b>	Fixed an issue where obfuscation was shown as deactivated in the Configuration Tool, although obfuscation was enabled.
<b>23828</b>	Fixed an issue where the Run Count value was too high, if the run contained a pause.
<b>23884</b>	Fixed an issue where tip types for Fluent Control only contained the size not the exact type (e.g. "FCA, 1000ul" shown instead of "FCA, 1000ul SBS").
<b>24118</b>	Fixed an issue where the IoT Client was not able to start, due to an issue with setting access rights for the current user.
<b>25082</b>	Fixed an issue that caused the upgrade installation to fail when the current user %AppData%/Temp directory contained inaccessible folders.

## New Features

ID	Title
<b>22108</b>	Number of measurements for the FRIDA reader are parsed with the next release of Introspect
<b>23184</b>	Extended logging for easier debugging
<b>23480</b>	If the proxy in the local IT environment decrypts SSL traffic, the connection cannot be validated with the certificate provided by the IoT Client. It's now possible to select an export of the certificate authority to use for validation during registration or via the Configuration UI.
<b>23601</b> <b>23791</b>	Number of moves for several axis is parsed with the next release of introspect.
<b>24491</b>	Registering a Spark device is now possible to send live data to the Connect App.
<b>24931</b>	Hardware information for the registered instrument can be sent for FluentControl version 2.7 and higher.



## IoT Client 2.1 SP1

---

- Version **2.1.3.28502**

### Bug fixes

ID	Title
<b>22526</b>	The IoT Client installer now allows upgrade installations. This resolves several issues with upgrade installations of or to IoT Client versions delivered with FluentControl.
<b>23385</b>	
<b>23347</b>	Fixed an issue, where the IoT Client was not able to detect the serial number during the registration process if only a single log file exists on the system.
<b>23898</b>	Fixed an issue, that copies of parsed log files are kept on the system in the user's temp directory. Existing copies, that were created by previous versions of the IoT Client will be deleted automatically by the IoT Client for the currently logged in user.

### New Features

ID	Title
<b>23590</b>	The installation of the IoT Client now delivers the CNS Driver if an EVOware installation is detected on the system.

## IoT Client 2.1

---

- Version **2.1.2.27667**

### Bug fixes

ID	Title
<b>19352</b>	Instrument registration and sending of usage data, now use the same proxy settings. If the proxy settings are not valid, the registration will already fail.
<b>20053</b>	Proxy authentication credentials are now used during registration and sending of usage data.
<b>20601</b>	Fixed an issue, where registration was possible, but sending of usage data was not, if the environment required a proxy server and only the proxy server was allowed to send DNS queries.
<b>21409</b>	Fixed an issue, that in some cases the count of used tips was too low.
<b>22071</b>	Fixed an issue, where the IoT Client was not able to start on a system with a localized Windows user system (e.g. on a German or French Windows version).
<b>22917</b>	Fixed an issue, where the IoT Client did not send configuration data anymore, after an online update of the IoT Client from 2.0.4 to a higher version.

### New Features

ID	Title
<b>18243</b>	The IoT Client parses paused/non value added run status in addition to running or idle status.
<b>20501</b>	The value of the variable "INTROSPECT_SampleCount" is now stored to the Sample Count property in Introspect.
<b>20888</b>	QC Data is added to the configuration data and sent to the cloud.
<b>21445</b>	Enlarged the field for Webcam URL in the Configuration UI to make editing of the URL easier.
<b>21450</b>	Added additional errors to be parsed by the IoT Client.
<b>21580</b>	Extended configuration information with public IP address also for registered Freedom EVO devices.
<b>21848</b>	Uninstallation of the IoT Client no longer deletes log files.
<b>21912</b>	The IoT Client handles the sending of Live Data asynchronously.

## IoT Client 2.0 SP2

---

- Version **2.0.5.25673**

### Bug fixes

ID	Title
18696	"INTROSPECT_"-variables are now correctly tracked when repeating the value.
19191	IoT Client does not send an empty Instrument State.
19682	Scanning for the instruments serial number during registration does not crash the registration if a log file of FluentControl or EVOware is read-only. If parsing a log file causes an error, this log file is skipped, and parsing does not stop.
19919	Already used serial number cannot be applied in settings anymore
19940	Registration now logs into the same directories as other components of the IoT Client.
19944	Registration now accepts email addresses with subdomains or two-part TLDs (e.g. co.uk).
20282	The user is informed if the instrument is already registered, instead of showing the error "The service encountered an error. Please try again later."
21224	Fixed problems that failed the uninstallation of the IoT Client in some cases.
21410	Using labware in FluentControl with a label, that contains only digits, does not stop the parsing anymore.

### New Features

ID	Title
19737	Use updated file structure of Fluent 2.6 for gathering of hardware configuration.
20038	Extended parsers to cover more errors.
20743	Log additional information of IoT Client status.
20791	"EVO" prefix from instrument size dropdown values during registration removed
21245	Extended configuration information with public IP address.

## IoT Client 2.0 SP1

---

- Version **2.0.4.23795**

### Bug Fixes

ID	Title
18673	Error handling and message during setup indicates wrong entries correctly.
18828	Configuration Management – Hardware information was not sent and missing in cloud
19039	IoT Client does not send historical data for connect user, only for Introspect user.
19411	Use validated versions for test tools on build PC.
19448	Time of day chart was incorrect if run is parsed live
19494	Single FCA tip usage was not counted
19522	"Run was aborted!" was not parsed and showed no ending runtime
19630	Parsing longer runs for EVOware, while those are running, results in incorrect usage data

## New Features

ID	Title
18829	IoT Client installation and registration process are separated to allow installation without need of login to Introspect account.

## IoT Client 2.0

- Version **2.0.2.22689**

## Bug Fixes

ID	Title
9635	Connection to cloud fails with “Unable to load DLL 'twApi.dll': Attempt to access invalid address” error
9636	EVOware MCA 96 and MCA 384 tip pick-ups are always counted with maximum tips for used MCA. (Fixed for EVOware 2.8 SP2 and higher)
9637	Uninstalling the Introspect Client does not remove all installation files
9638	Entering wrong credentials, when registering an EVO instrument, forces the user to reenter all data and leads to incorrect instrument type in serial number
13552	Obfuscation is incorrectly set when run or method name are empty
15222	Tips and labware usage cannot be parsed for FluentControl 2.5
18457	Variables cannot be parsed for FluentControl 2.5
18459	IoT-Client should not run with elevated privileges if not required
18680	Operator is reset to empty after the first parsed run

## New Features

ID	Title
9652	Improved the visibility of the instrument registration window during installation
9651	Entry for Method Name is parsed for EVO log files
9667	Performance improvement: No need to parse all log files again if parsing was interrupted
9669	Reworked configuration tray tool with about box and update functionality
9676	New IoT Client that supports parsing of log files and sending live data via the same application
9677	Introspect parsers in the cloud, that allow online updates
9678	Installed version of FluentControl & EVOware is provided for Connect app
9680	IoT Client supports online updates
9683	Serial number is parsed from log files and is not editable
9688	Support for variables with user defined variable names
10562	Obfuscation is not activated by default
18456	Add link to trial account registration to instrument registration window during installation
-	Tecan Connect mobile app (CNS successor) support
-	Introduced new functionality to pair an instrument to Tecan Connect app