

Revision History

Filename	Revision History IoT Client
Author	Artur Bauer
Corresponds to IoT Client Version	IoT Client 3.1.0.326
Date	2025-06-04

Contents

IoT Client 3.1.....	3
Bug fixes.....	3
Changes	3
IoT Client 3.0.....	3

IoT Client 3.1

- Version **3.1.0.326**

Bug fixes

ID	Title
40438	[IoT Client] IoT Client doesn't start automatically when anything updated on Configuration tool
40457	[IoT Client] Remote update notification should be available all the time until it is installed in IoT Client
40603	[IoT Client] Web Cam Url not sent to IoT hub as a property update
40995	[IoT Client] Changes in configuration tab shall be logged
41912	[IoT Client] IoT Client crashes when Alias changed in OM and IoT config tool is being used when the change is in progress
43233	[IoT Client] Upload the system config and qc data, instrument info file
44038	[IoT Client] Update only "Proxy Host" and "Port" without Alias

Changes

ID	Title
40954	[IoT Client][MAP] Implement full MAPlinx compatibility
41790	[IoT Client] Update to .NET 8.0
43454	[IoT Client] Configuration tool shows errors for the ping of the URLs, but communication works fine
43690	[IoT Client] Handle case: .NET not preinstalled
44015	[IoT Client] Rename label of registration button

IoT Client 3.0

- Version **3.0.10.251**

Initial release

The IoT Client 3.0 is a newly developed component, that will be completely verified and will not replace previous versions of IoT client for legacy applications. **Legacy application support will be available with 3.1 and higher.**

Note: Support of legacy applications with IoT Client 2.8. IoT Client 2.8 is based on the old architecture. See also "04207_00794_Revision History IoT Client V2.8 SP1".