

Revision History

Filename	Revision History IoT Client
Author	Artur Bauer
Storage place of valid version	DHF Number 04207_00718 V1.0
Corresponds to IoT Client Version	IoT Client 3.0.10.251
Date	2024-07-17

Contents

IoT Client 3.0.....	3
---------------------	---

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.**