

ReadMe FluentControl 3.7 P2 Patch

2026-09-10

Included changes

PWI	Description
131074	Robustness improvement for MCA96 Tips pickup – alignment move with ejector plate
130661	Robustness improvement for FCA Tips pickup – optimized retract move
130190	Continue cLLD processing after FW error 92
132751	Reduce hold-power after doorlocks engage to reduce residual magnetism

Pre-requisites

You must have local Windows admin rights.

FluentControl 3.7.12 must be installed on the system and optionally the firmware patch 3.7 P1

Installation instructions

- (1) Shutdown FluentControl
- (2) Unzip file 3.7 P2.zip
- (3) run install.bat as windows administrator
- (4) Start FluentControl
- (5) Import updated labware definitions from FcaMca96Tips.zeia