

QPatch Compact Manifold Recovery Guide

In case of unforeseen mechanical events, the manifold may be jammed into a position where it will not release itself during normal operation. Follow this guide to recover the manifold via the built-in recovery function.



Start up and attempt homing

- 1. Read through guide** Please read the entire guide before starting the recovery process. Some steps need to be monitored manually.
- 2. Close latches**

Open the instrument lid and ensure that both latches are fully engaged.

If the manifold has been dislodged from the interface module so latches cannot fully close, make sure to push the manifold and the interface module together as much as possible without closing the latches.

Caution! Never use excessive force to attempt closing the latches.
- 3. Start up the QPC** Turn on the instrument (on the left-hand side). Log in via the touchscreen using your normal login credentials for the QPC database.
- 4. Wait for homing**

If the manifold releases itself during homing, remove the QPlate after homing is done and proceed to checking manifold latches and verifying functionality.

If the manifold is stuck, the instrument will make a low-pitched humming during homing and the "Preflight Check" status will stay in "Initializing QPlate slot". In that event proceed to the next step.



Run recovery function

- 1. Enter Maintenance menu** Click the Maintenance icon on the left-hand side to enter the maintenance menu.
- 2. Start recovery function** Press the "Recover from jam" button in the "Manifold" menu to start the recovery process. A countdown timer next to the button will show the remaining time for the process.

The recovery function will attempt moving the manifold back and forth between the intra- and extracellular positions via a forced move. The instrument may make a low-pitched humming during the process. This is normal.

The humming should stop during the process or at the latest a few seconds after the countdown has ended.

Caution! *If the humming does not stop shortly after the countdown, open the instrument lid to stop the motor.*
- 3. Monitor manifold move** Monitor the manifold through the pipetting access cut-out during the process. When the manifold starts lifting, it is recommended to very gently pushing down on the protruding QPlate handle a couple of times to aid the release process.

Caution! *Do NOT attempt to remove the QPlate during the recovery process.*

Caution! *Do NOT touch the manifold during the recovery process.*
- 4. Wait for countdown** When the countdown has stopped, remove the QPlate if possible.
- 5. Close latches** Open the instrument lid to check and close the manifold latches. See the user manual and installation manual on how to check the manifold latches.

