

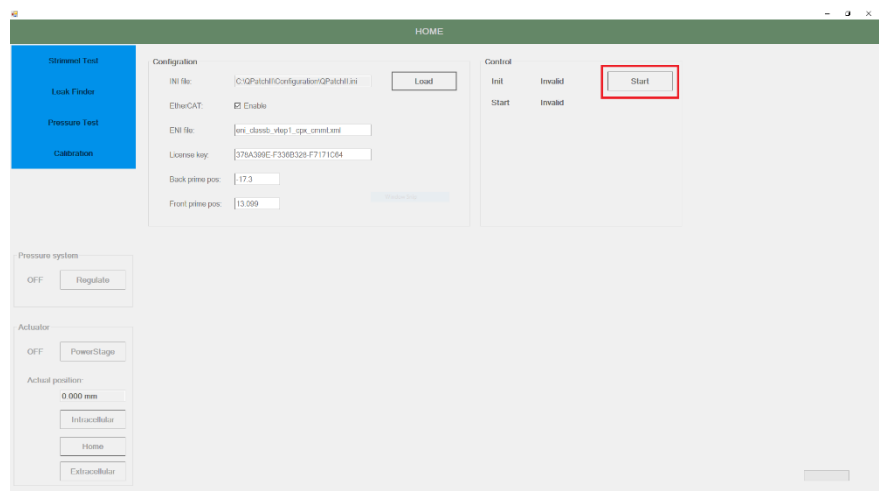
QPatch Compact pressure test

In order to make sure the QPatch Compact instrument is working as expected, e.g. after shipping or service, it is possible to execute a pressure test to ensure that pressure connections to the QPlate are leak tight within specifications.

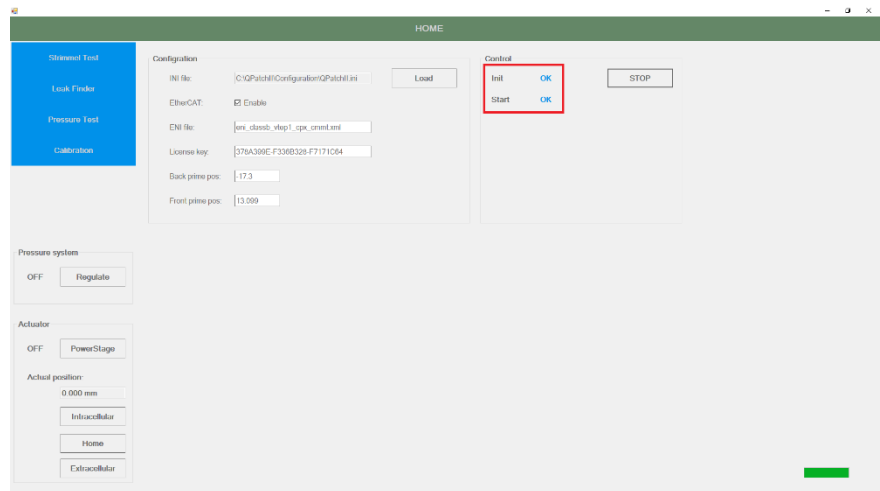


Conducting a pressure test

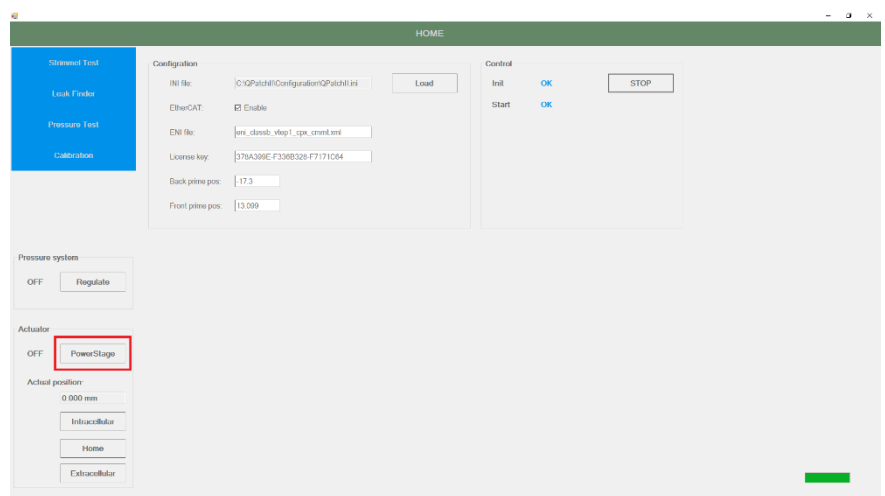
1. **Locate the pressure ARQ plate** The QPatch Compact is delivered with a special Artificial Reference QPlate (ARQ) pressure plate. It has a red frame and a label stating "Pressure ARQ".
2. **Start up the QPC** Turn on the instrument (on the left-hand side).
3. **Close main application** At the login screen of the main application, close it by pressing the "X" in the corner. It may take a while before the button is clickable due to initializing of the instrument.
4. **Open test application** Locate and start the application "VtepTest.exe" found in the C:\QPatchII\tools folder.
5. **Initialize the system** Click "Start".



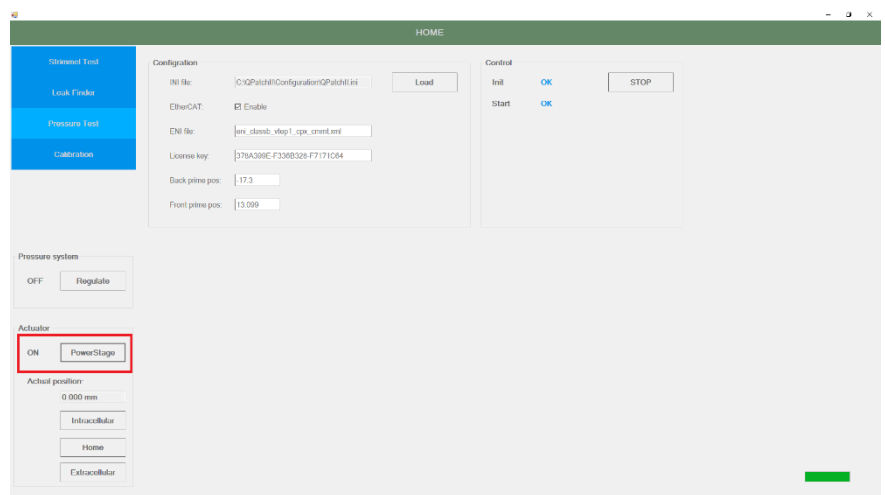
Wait for it is finished with "OK".



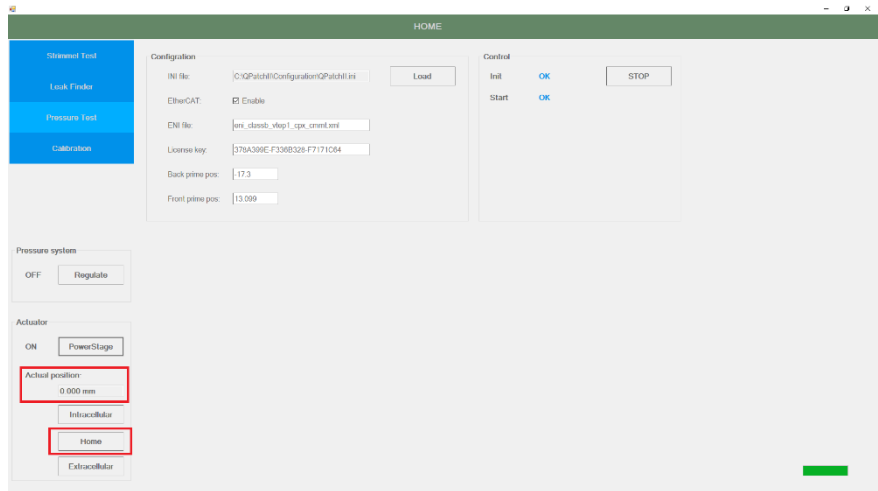
Click "PowerStage" in the "Actuator" section in bottom left side.



"Off" state will change to "On".



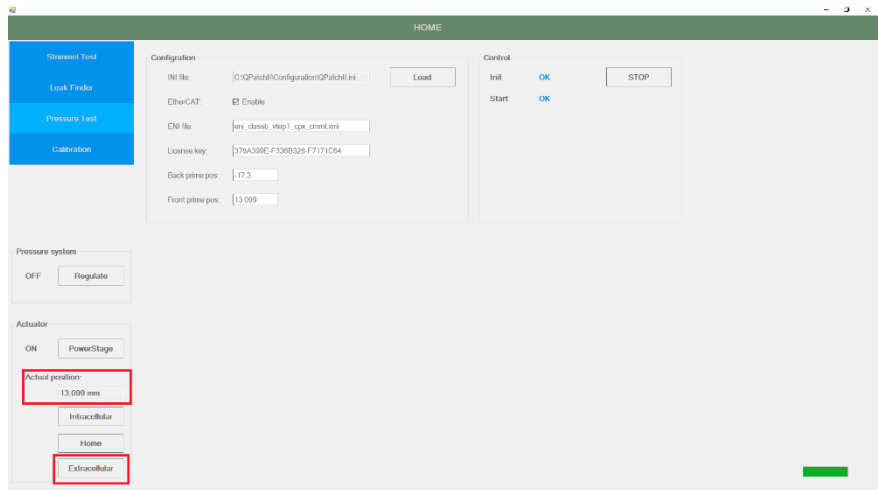
Click "Home" and wait for the actuator finishing its movement and the "Actual position" is at 0.



Prepare the test

Insert a pressure ARQ QPlate into the system.

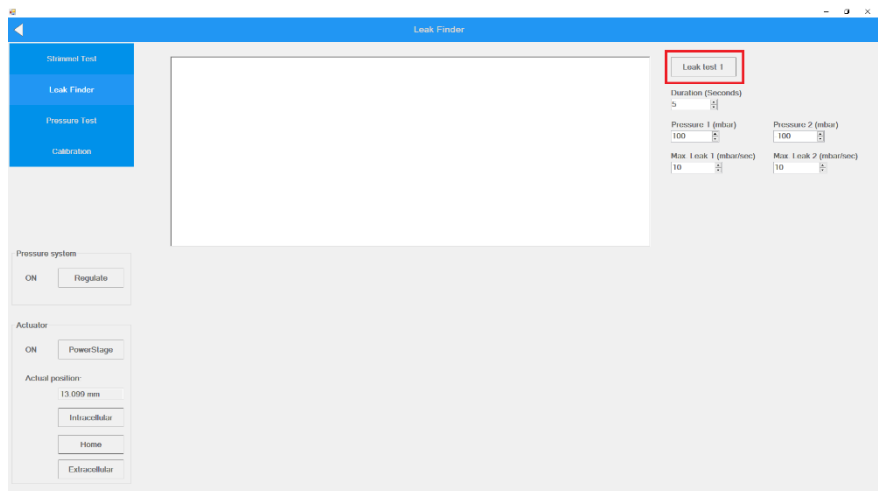
Click "Extracellular" and wait for the actuator finish its movement when the "Actual position" is idle.



Perform the test

Click on "Leak Finder" in the top left section.

Click "Leak test 1" using default parameters.



Wait for the test to perform.

Read the test outcome and expect a message saying "No leaks detected".

