


NanoDrop Ultra Spectrophotometers and Fluorometers

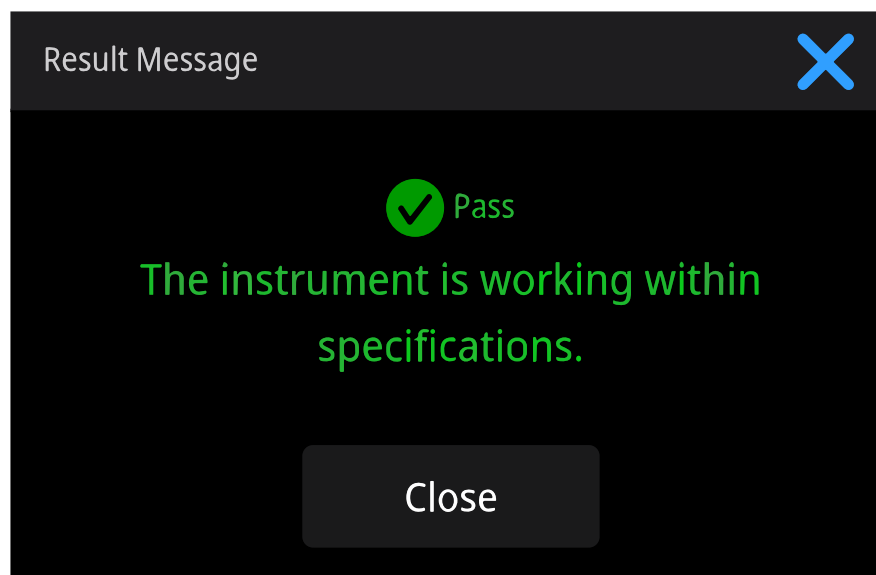
Pedestal image check


Materials needed

- Lint-free laboratory wipes

Pedestal image check procedure

1. From the home screen, select the **Diagnostics** icon , followed by **Pedestal Image Check**.
2. Clean the upper and lower pedestals using a clean, dry laboratory wipe. Visually confirm that there is no cuvette in the cuvette well.
3. Select **Measure** to begin the measurement.
4. After measurement is complete the overall result of the Pedestal Image Check will be displayed on screen. Select **OK** or **Close** to see a more detailed summary of the results.
5. Select **End Experiment** when done.



6. The experiment name can be changed at this time and up to five unique identifiers/tags can be added, once complete, select **Next**.
7. Results can be exported and printed at this time by selecting **Continue** or at a later time from the History.
 - After exporting or printing, select **OK** to go back to the Diagnostics screen.
8. If results do not need to be exported or printed, select **Finish** to return to the Diagnostics screen.
9. To review results from a previous Pedestal Image Check, select the **History** icon  from the home screen and locate the Pedestal Image Check results from the list of experiments.

Interpreting the results

1. The overall result of the test will then be displayed at the top of the screen as a green “Pass” or red “Fail”.
2. The screen will also display the parameters of the Pedestal Image Check including Empty column detection, Sharpness, Background exposure, Background quality, and Edge detection.
3. In the case of a failed Pedestal Image Check, immediately repeat the test ensuring that the image sensor is clean and not blocked. Rotating the instrument 90° or moving it away from a nearby window can also help. If the test fails again, contact NanoDrop Technical Support (nanodrop@thermofisher.com) or your local distributor.

