


Performing an Intensity Check




Materials Needed:

- Lint Free Laboratory Wipes

Pedestal Intensity Check Procedure:

1. From the Home Screen, tap the **Diagnostics** icon , followed by **Intensity Check**.
2. Clean the upper and lower pedestals using a clean, dry laboratory wipe.
3. If **Auto-Measure** feature is off, tap **Measure** to begin the measurement. If **Auto-Measure** feature is ON, the measurement will begin automatically after the arm is lowered.

Cuvette Intensity Check Procedure:

1. From the Home Screen, tap the **Diagnostics** icon , followed by **Intensity Check**.
2. Tap the **Menu** icon , followed by **Settings**.
3. From the **General** tab, tap the box next to **Use cuvette** and confirm that the box is checked.
4. Tap **Done**  when finished.
5. Visually confirm that there is no cuvette in the cuvette well.
6. Tap **Measure** to begin the measurement.

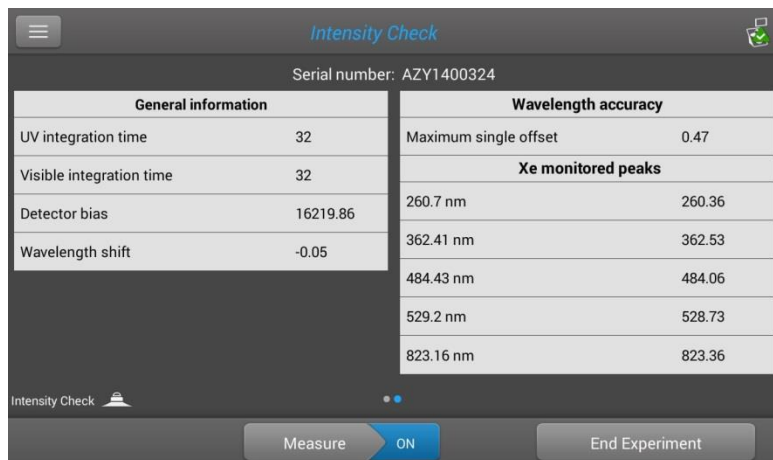
Interpreting results:

1. The results will display as green and orange spectra.
2. A green check mark will be seen adjacent to the “UV”, “Visible”, and “Bias” notations if they are within specification. If they are not in specification, a yellow check mark will display.



3. Slide your finger to the left across the screen to display a second screen showing more information including: Integration times, detector bias, and the Xe monitored peaks. See image below.

Performing an Intensity Check



4. Tap **End Experiment** when done
 - A. Results can be exported and printed at this time or at a later time from the Data Viewer
 - B. Experiment name can be changed at this time and Identifiers can be added
5. Tap **End Experiment** when done
6. To review results from a previous intensity check, select Data Viewer from the Main Menu and locate the intensity check results from the list of experiments