


## Performing a Pedestal Image Check

### Materials Needed:

- Lint Free Laboratory Wipes

### Pedestal Image Check Procedure:

1. From the Home Screen, tap 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. 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.

### Interpreting results:

1. The results will show an image of the upper and lower pedestals. To the right of the image, the instrument will indicate whether the diagnostic passed or failed.
2. A green triangle with a check indicates that the instrument passed the Pedestal Image Check. A yellow triangle with an exclamation point indicates that the instrument failed the Pedestal Image Check. See Figure 1 below.
3. Slide your finger to the left across the screen to display a second screen showing the parameters of the Pedestal Image Check including: Sharpness, Background exposure, Background quality, and Edge detection. See Figure 2 below.

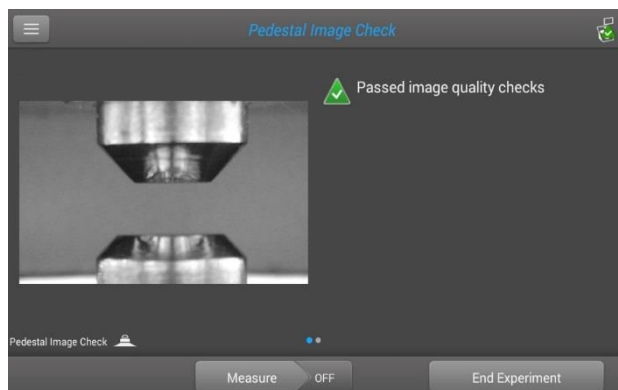


Figure 1



Figure 2

S083 Rev 18 January 2018