

Mettler Toledo Calibration Instructions (For imPULSE Systems) Quick Start Guide



This guide provides a brief overview of calibration for the Mettler Toledo™ IND331 and IND560 models used on the Thermo Scientific™ imPULSE™ Single-Use Mixers (S.U.M.s). Please take the time to review the imPULSE S.U.M. User's Guide (DOC0043) for important information on safety and operation.

Mettler Toledo IND331 Calibration steps

IND331 control descriptions and navigations:

Figure 1 includes a description of the scale function keys. In the Setup Menu, the scale function keys are used for navigation as described in Figure 2.

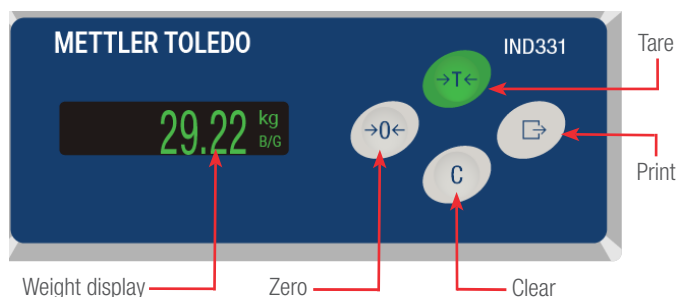


Figure 1. Functions of the scale function keys in the Setup Menu.

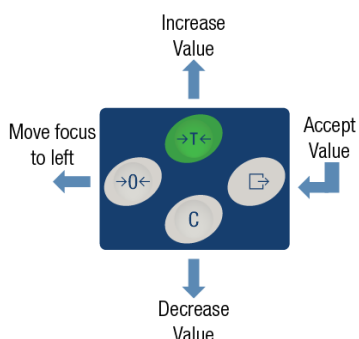


Figure 2. Description of navigation using the scale function keys.

Process steps

1. Press and hold the **Print** key (see Figure 1) for three seconds. The display will change from showing a normal weight to the information recall icon: .
2. Press the **Clear** key (Figure 1) until the Setup icon appears , then press the **Print** key to enter the Setup Mode. **Note:** This gives access to all setup parameters for the terminal.
3. Use the scale function keys to navigate (see Figure 2) to Set Zero in the setup menu. See Figure 3 for the path in the menu tree.

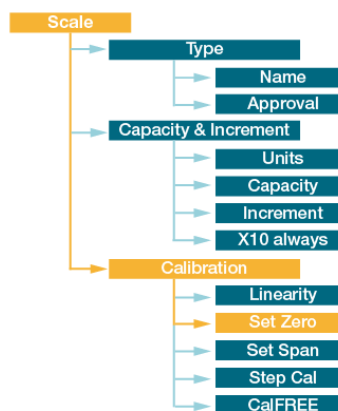


Figure 3. Location of Set Zero in the Setup Menu.

- Follow the prompts and remove all weight. Press **Print**. The display will indicate “Capture Zero Ok”.
- After Set Zero is completed, press the **Print** key. Use the scale function keys to navigate through the calibration menu again to select **CalFREE** (Figure 4) and press **Print**.

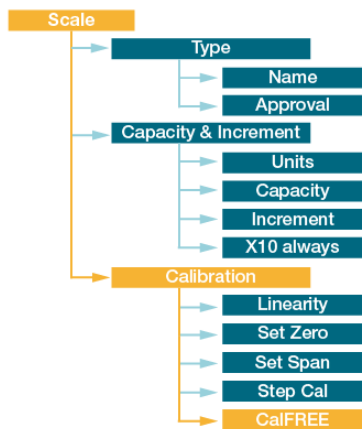


Figure 4. Location of CalFREE in the Setup Menu.

- After selecting **CalFREE**, follow the prompts and either input the full scale weight (using the navigation keys), or accept the current value and press the **Print** key.
- Press the **Print** key again to accept the rated cell output. The display will indicate “Perform CalFREE?”
- Press **Print**, and after cycling through the percentage completed, the display will read: “CALIBRATION OK.”
- Press **Print** to return to the scale menu.
- Press the **Zero** key twice to exit the Setup Menu and then press the **Tare** key until you return the the weight indication display.

Mettler Toledo IND560 Calibration steps

IND560 control descriptions and navigations:

Figure 5 includes a description of the scale function keys. Use the **Up** and **Down** navigation (nav.) keys to navigate the menu tree. Use the **Right** nav. key to expand a branch, and the **Left** nav. key to collapse a branch.

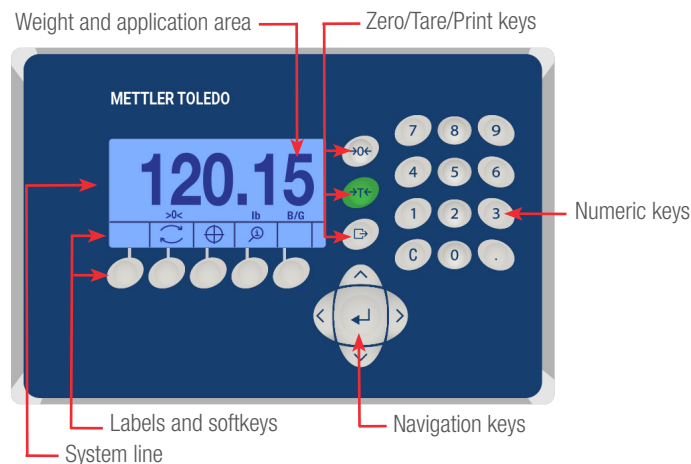


Figure 5. IND560 control descriptions.

Process steps

- Press the **Down** nav. key (Figure 5) to view the Setup Menu softkey label \diamond .
- Press the **Setup Menu** \diamond softkey (Figure 6).

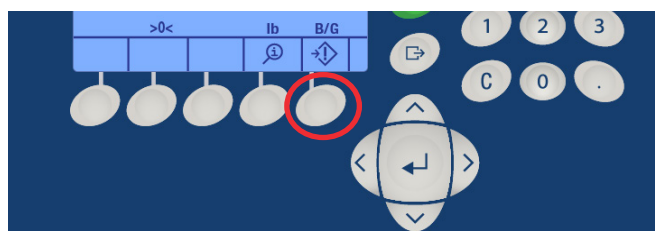


Figure 6. Softkey that corresponds with Setup Menu icon.

- Navigate to **Calibration** (Figure 7).
- Once the desired menu option is highlighted, press the **Enter** key (center nav. key).

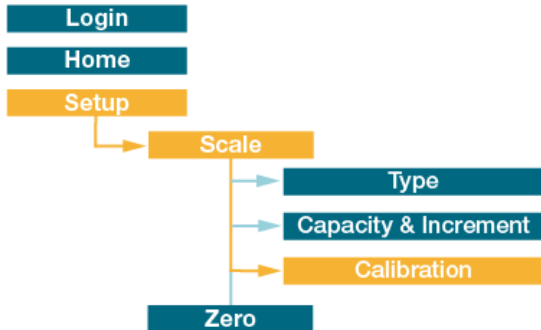


Figure 7. Location of Calibration in the Setup Menu.

- The calibration menu has multiple pages similar to those shown in Figures 8 and 9. Press the **Down** nav. key to navigate through the calibration menu pages until the Softkey Calibration Labels are visible.

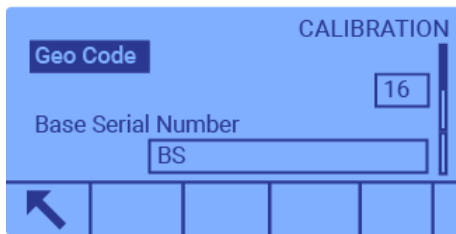


Figure 8. Softkey Calibration Labels not visible on page 1.

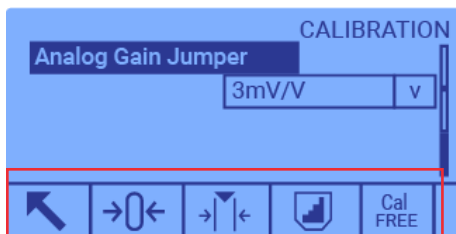


Figure 9. Softkey Calibration Labels visible on page 3.

- Press the **Capture Zero** $\rightarrow 0 \leftarrow$ softkey.

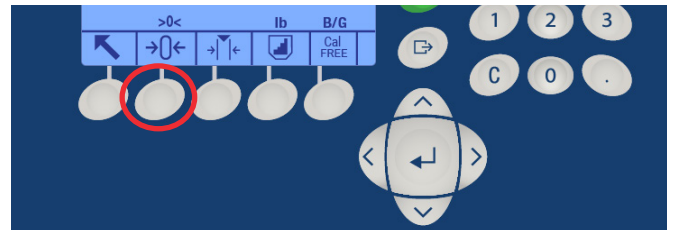


Figure 10. Softkey that corresponds with the Capture Zero icon.

- Remove any weights and follow the prompts. Press the **Start** \diamond softkey (Figure 11). The screen will read "Capture Zero Ok."

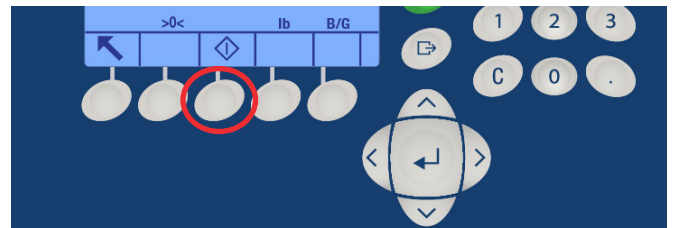


Figure 11. Softkey that corresponds with the Start icon.

- After Set Zero is complete, press the **Exit** \swarrow softkey (Figure 12) to exit the current menu.

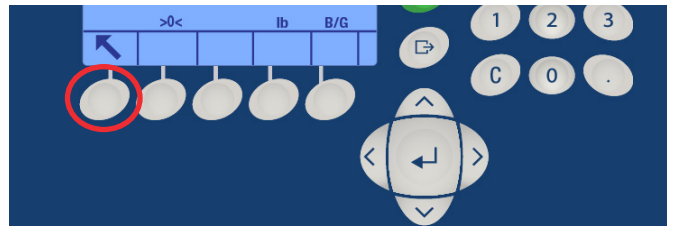


Figure 12. Softkey that corresponds with the Exit icon.

- Press the softkey that corresponds with the **CalFREE** softkey label (Figure 13).

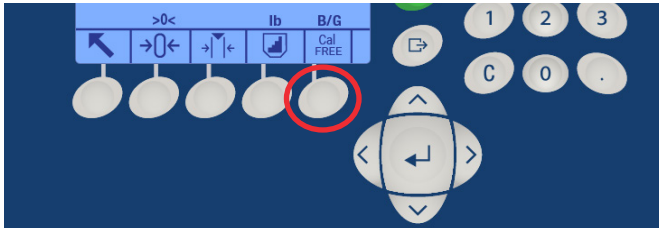


Figure 13. Softkey that corresponds with the CalFREE icon.

- Navigate to Home in the Setup Menu tree and press **Enter** (center nav. key) to return home.

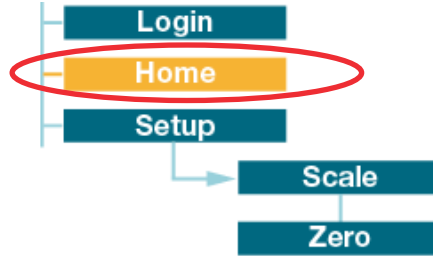


Figure 16. Location of Home in the Setup menu.

- Press the **Down** nav. key to navigate through the CalFREE menu pages until the **OK** softkey label is visible (Figures 14 and 15).

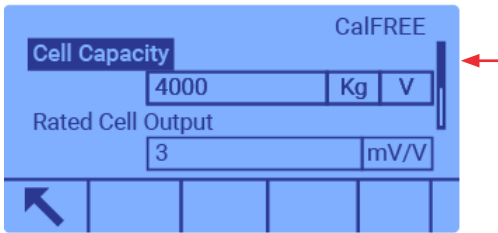


Figure 14. OK softkey label not visible on page 1 of the CalFREE menu.

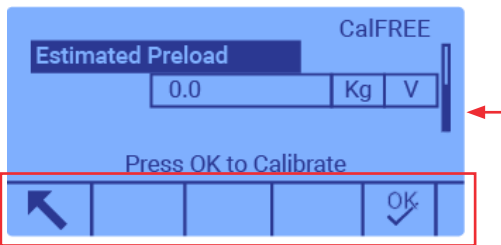


Figure 15. OK softkey label visible on page 1 of the CalFREE menu.

- Follow the prompts and press the **OK** softkey. The screen should read “Calibration OK.”
- Press the softkey that corresponds with the **Exit** softkey twice to return to the navigation menu.

Find out more at thermofisher.com/sut