

TC10™ Verification Slide

Catalog #145-0014

Product Description

Printed glass slide

For research purposes only. Store at room temperature.

Product Content

The TC10 verification slide is a glass slide with a defined number of circles simulating live and dead cells printed on sides A and B of the slide. The verification slide is used to validate functionality of the TC10 automated cell counter. The TC10 automated cell counter is calibrated at the time of manufacture and does not require further adjustment by the customer.

Use

Materials needed:

TC10 verification slide

TC10 automated cell counter

Procedure

To perform the TC10 system test:

1. Power-on the TC10 automated cell counter.
2. From the Home screen, use the down arrow key to select **Options**, then press **Enter**.
3. From the Options screen, use the down arrow key to select **System test**, then press **Enter**.
4. “Insert verification slide, side A first, then press Enter” appears on the screen. Insert side A of the verification slide into the slide port of the instrument. Press **Enter** to continue.
5. “Pull out verification slide and insert side B into the instrument, then press Enter” displays on the screen. Remove the slide from the instrument and insert side B of the verification slide into the slide port. Press **Enter** to continue.

6. Pull out the verification slide and press **Enter**. (This is a test step without the slide to check for debris on the optics.)

7. The results of the system test are displayed on the System test screen. For each sample, a Pass or Fail result is reported, along with the number of objects found, the number of objects expected, and the ratio of live objects (example: Pass 218 of 221, 0.50). The number expected to be found can change each time the test is repeated due to slight position variations of the slide in the TC10 cell counter.

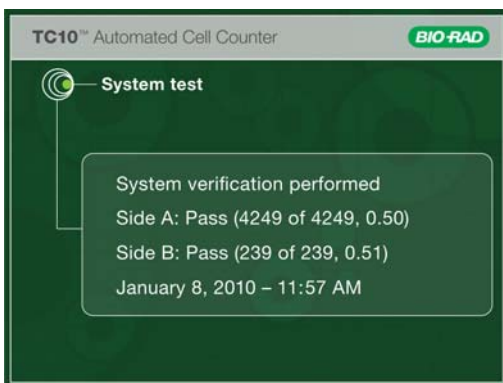


Fig. 1 System verification results

8. Results of the TC10 system test can be exported via a USB device using the Export previous counts function.

9. Save the TC10 verification slide for future use.

Important: If the TC10 cell counter fails the system test procedure, a reason for the failure will be displayed on the System test screen.

The TC10 counter will indicate that you should either clean the verification slide and repeat the test or that you should contact Bio-Rad Technical Support.

Do not insert verification slide if it is damaged. It could damage the instrument.