

# Agilent 708-DS Dissolution Apparatus: Pre-Test Preparation in Under a Minute

## Pre-Test Preparation

The following checklist can be completed in less than a minute. It will help to ensure that the 708-DS is properly configured and setup to run your dissolution test. Making this quick and easy procedure part of your routine can help to eliminate testing errors and prevent potential investigations.

## Pre-Test Diagnostics

1. Once the method is loaded – either from the 708-DS, 850-DS, or software – examine your environment. Make sure that the area is free of any objects that could contribute to interference when running the dissolution test (e.g., vibration).
2. Complete the following visual checks:
  - A. Are the paddles or baskets clean, in good condition, and set at the proper height?
  - B. Are the vessels clean and in good condition (e.g., free from cracks)?
  - C. Have the vessels been filled with the proper media and volume?
  - D. If applicable, have the correct filters been properly installed on the sample cannulas?
  - E. If performing media replacement, has an appropriate length of tubing been connected to the proper 850-DS port and properly submerged in the media container? If using vessel position 7 or 8, ensure the tubing is free from any shaft interference when rotating.
  - F. If applicable, has the 850-DS sample tray been loaded with an appropriate number of tubes/vials for sample collection?

3. From the Main screen of the 708-DS, press **Menu**.
4. From the Menu screen, press **Diagnostics**. The following screen displays:
5. In the Spindle box, press Run. Confirm that all shafts are secure and operational. Press Stop.
6. Verify that the bath temperature is set to the appropriate value.
7. In the DDM control box, press All to start all DDMs. Ensure that each DDM opens and closes freely.
8. Press Finish to return to the Main Menu.
9. Press Main to return to the Main screen.
10. In the Manifold box, press the down arrow to lower the manifold. Ensure the manifold and all tubing lines move freely.
11. If applicable, verify that the vessel temperatures are visible, uniform, and within tolerance before beginning your test.
12. Press the up arrow to raise the manifold.

## Key Reminders

- If manual sampling, it is a good idea to prepare the area around the apparatus with all necessary items – syringes, cannulas, filters, etc. A sample timer is available within the 708-DS or an external timer could also be used to ensure that sampling is not missed.
- If an 850-DS is configured, ensure any solutions required – for media replacement or cleaning – are sufficiently filled and connected.
- For any troubleshooting assistance, you may contact [dissolution.hotline@agilent.com](mailto:dissolution.hotline@agilent.com)

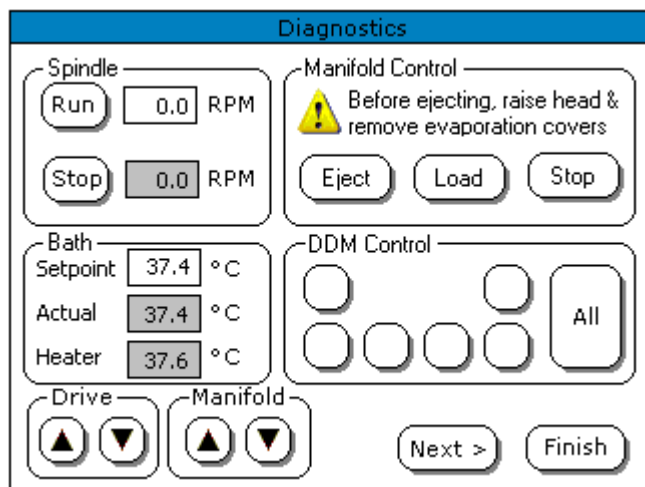


Figure 1. Diagnostics box on 708-DS.

[www.agilent.com/chem/dissolution](http://www.agilent.com/chem/dissolution)

DE44474.4604861111

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021  
Published in the USA, October 01, 2021  
5994-4248EN