

xCELLigence RTCA HT—High Throughput

System maintenance

This quick reference guide describes common procedures for maintenance of the Agilent xCELLigence RTCA HT Station (part number 5919878001). It is suggested that maintenance be performed at least annually or as directed below.

General cleaning (excluding the contact pins)

Using a soft cloth or fiber-free tissue paper moistened slightly with 80% alcohol, gently wipe off any dust or contaminants from the surface of the HT Station. A small air blower can also be used to help remove any dust.

Operational check

In the System menu, click "System -> QC Test" and select "Operational Check". Run the test with a resistor plate 384 (part number 5867878001). Upon completion of the test, the test results will show either pass or fail. If any of the wells failed, the contact pins should be cleaned or replaced as shown below.



Regular cleaning of the RTCA contact pins 384

If a well is out of specification, cleaning of the Agilent xCELLigence RTCA contact pins is recommended. It's recommended to clean the RTCA contact pins 384 after every experiment or when contaminants are present. Power off and unplug the HT Analyzer. Carefully remove the RTCA HT Heating Plate from the HT Station. Clean the entire surface of the RTCA HT Station using the dust blower supplied in the RTCA HT cleaning kit. Brush the RTCA contact pin 384 with the nylon brush provided with the system. Brush gently across the tips of the contact pins approximately 10 times.

Extended cleaning of the RTCA contact pins 384

If a well is consistently out of specification even after a regular cleaning procedure, an extended cleaning procedure is recommended. It's recommended to perform extended cleaning of RTCA contact pins 384 once a month or after 2,000 cycles. See the RTCA HT Operator's Guide under the "Maintenance and Care" chapter for detailed instructions.

Cleaning the RTCA HT Heating Plate

If the contact pin cleaning procedures didn't solve the contact issue, cleaning the RTCA HT Heating Plate is recommended. Refer to the xCELLigence RTCA HT Operator's Guide "Maintenance and Care" chapter for detailed instructions.

Exchange the RTCA contact pins 384

If the test result for any well from the Operational Check shows "F" (fail) even after the contact pin and Heating Plate cleaning procedure, the corresponding contact pin should be replaced (part number 380601270). See the RTCA HT Operator's Guide "Maintenance and Care" chapter for detailed instructions.

Inspection of the RTCA cables

The two cables connecting the RTCA HT Station, Analyzer, and control unit (computer) should be inspected annually. If any sign of breakage or corrosion is observed, the cables should be replaced.

Decontamination of the RTCA HT System and RTCA control unit

The following compositions have been tested for compatibility with the surfaces of the RTCA HT Station, Analyzer, control unit, and cables:

- 80% Ethanol
- A mixture containing propan-1-ol (450 mg/g), propan-2-ol (250 mg/g), and ethanol (47 mg/g). This mixture is available from Bode Chemie under the tradename Bacillol AF. It is fungicidal, tuberculocidal, mycobactericidal, and virucidal against enveloped viruses (including HBV, HIV, HCV), adenovirus, FCV, papovaviruses, and rotavirus when used as described in the supplier's instructions.

Ordering information

Product	Description	Part Number
Consumables		
xCELLigence RTCA E-Plate 384	2 x 5 plates	5867681001
xCELLigence RTCA E-Plate 384	8 x 5 plates	5867673001
IMT assay (anti-CD40) tethering kit	Tethering reagent (250 µL), buffer and cytolysis reagent	8100005
IMT assay (anti-CD29) tethering kit	Tethering reagent (250 µL), buffer and cytolysis reagent	8100008
IMT assay (ant-CD19) tethering kit	Tethering reagent (250 µL), buffer and cytolysis reagent	8100011
IMT assay (anti-CD71) tethering kit	Tethering reagent (250 µL), buffer and cytolysis reagent	8100017
Maintenance items		
xCELLigence RTCA resistor plate 384		5867878001
xCELLigence RTCA contact pins 384, rounded (20 units)		380601270
EPS cable RTCA HT		6204821001



www.agilent.com

RA45089.5580671296

This information is subject to change without notice.



© Agilent Technologies, Inc. 2023, 2024
Printed in the USA, March 14, 2024
5994-6620EN

 **Agilent**
Trusted Answers