



# Seahorse XFp Extracellular Flux Analyzer

## User Manual



# Notices

© Agilent Technologies, Inc. 2020

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

## Manual Part Number

S7802-90001

## Edition

Second Edition, November 2020

Printed in USA

Agilent Technologies, Inc.  
121 Hartwell Ave.  
Lexington, MA 02466



Manufactured by:  
Agilent Technologies, Inc.  
300 Griffith Rd,  
Chicopee, MA 01022

## Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data -Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A CAUTION notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a CAUTION notice until the indicated conditions are fully understood and met.

### WARNING

A WARNING notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a WARNING notice until the indicated conditions are fully understood and met.

# Contents

## 1 Introduction

- Safety Considerations **5**
- Overview **6**
- Technical Specifications **7**

## 2 Installation

- Unpacking and Component Identification **9**
- Suitable Locations for the XFp Analyzer **10**

## 3 Basic Operations

- Power and Warm Up **11**
  - Power up **11**
  - Welcome screen on first time power up **11**
- XFp Analyzer Communications and Network Connection Setup **12**
- Turning the XFp Analyzer Off **12**

## 4 Navigating the XFp Analyzer

- Setting Up an XFp Assay **14**
- Running an XFp Assay **19**
- Modifying XFp Analyzer Settings **24**
- XFp Analyzer System Files **25**
  - System files **25**
  - Diagnostics tests **26**
  - System settings **29**
  - Assay results **36**
  - Template management **37**
  - Help **38**

|   |  |    |
|---|--|----|
|   | Using Safe Lock (Trend Micro)                        | 40 |
|   | Cleaning and Routine Maintenance                     | 41 |
|   | Air filter replacement                               | 41 |
|   | Analyzing XFp Analyzer Data Using Seahorse Analytics | 42 |
| 5 | <b>XFp Analyzer Network Setup</b>                    |    |
|   | Introduction   | 48 |
|   | Why should I network my XFp Analyzer?                | 48 |
|   | Network access features on the XFp Analyzer          | 50 |
|   | XFp Analyzer System Information                      | 51 |
|   | Wired Network Setup                                  | 53 |
|   | Wired connection                                     | 53 |
|   | Wireless Network Setup                               | 55 |
|   | Wireless connection                                  | 55 |
|   | Joining a wireless network                           | 55 |
|   | Shared folder setup                                  | 57 |
|   | Email configuration                                  | 58 |
|   | Network Checklist                                    | 62 |
| 6 | <b>Support</b>                                       |    |
|   | Troubleshooting Guide                                | 63 |
|   | Technical Support and Ordering Information           | 64 |
|   | Worldwide technical support                          | 64 |
|   | Ordering   | 64 |
|   | Online help and support                              | 65 |
|   | Additional Resources                                 | 67 |

# 1

## Introduction

Safety Considerations 5

Overview 6

Technical Specifications 7

## Safety Considerations

### **WARNING**

**The protection provided by this instrument may be compromised if it is used in a manner not specified by Agilent.**

Safe operation of the Agilent Seahorse XFp Analyzer requires all covers to be securely attached and the plate tray door to be closed. This also prevents heat loss and system cooling, which can affect data quality.

The door opens automatically when the tray is extended, allowing the operator to insert or remove the assay consumables. Exercise caution when loading the miniplate/cartridge to avoid the possibility of a pinch hazard. After the miniplate/cartridge is placed securely on the tray, the remove your hand from the tray area before continuing the assay. After the command is given to continue the assay through the user interface, the tray will move slowly back into the instrument, and the door will close. An optical sensor is used to determine the status of the door.

The XFp Analyzer has heaters around the miniplate that maintain a stable tray temperature. The tray temperature is maintained at 37 °C, and monitored by temperature sensors and controllers embedded above the tray. A thermal fuse disables the heater when it reaches an abnormally high tray temperature.

Do not replace the power cord provided with any other power cord that is rated at less than what is specified in **“Technical Specifications”** on page 7. This depends on the power mains of the country where the instrument will be used.

## Overview

The XFp Analyzer measures the rate of change in dissolved oxygen and pH in the media immediately surrounding living cells cultured in a miniplate. Changes in the extracellular media are caused by cell consumption or production of analytes. A sensitive measurement of the media flux can be used to determine rates of cellular metabolism with precision and in a noninvasive, label-free manner.

A unique feature of XF technology is its ability to make accurate and repeatable measurements in as little as five minutes. This is accomplished by isolating an extremely small volume (approximately 2  $\mu$ L) of media above the cell monolayer. Cellular metabolism causes rapid, easily measured changes to the microenvironment in this small volume.

A measurement cycle is performed in six minutes. The media is gently mixed and the analyte levels are then measured until the oxygen concentration drops approximately 20 to 30%, and media pH declines approximately 0.1 to 0.2 pH units. The measurement is performed using optical fluorescent biosensors embedded in a disposable cartridge. The cartridge is placed into the XFp Analyzer cell culture miniplate.

Baseline metabolic rates are typically measured three to four times, and are reported in pmol/min for oxygen consumption rate (OCR) and in mpH/min for extracellular acidification rate (ECAR). The compound is then added to the media and mixed, and then the post-treatment OCR and ECAR measurements are made and repeated. As cells shift metabolic pathways, the relationship between OCR and ECAR changes.

The XFp Analyzer system includes:

- Benchtop analyzer
- Disposable sensor/compound
- Delivery cartridges
- Tissue-culture treated miniplates for seeding cells for analysis
- Calibration solution
- Seahorse Analytics for XF data analysis

Consumables are sold separately and include FluxPaks (sensor cartridges, cell plates, and calibrant) as well as various assay kits, reagents, and media. Seahorse XFp sensor cartridges are available exclusively from Agilent.

# Technical Specifications

## **REF** S7802A

|   |  |
|---|--|
| <b>Model</b>                              | Seahorse XFp Analyzer (model number S7802A)  |
| <b>Dimensions</b>                         | Width × height × depth<br>12" × 17" × 23"<br>30 cm × 43 cm × 58 cm   |
| <b>Weight</b>                             | 33 lb /15 kg   |
| <b>Power requirements</b>                 | 100-240 VAC 50/60Hz<br>6A/3A   |
| <b>Power cord rating</b>                  | 3- wire (grounded) AC power cord rated 10A or greater  |
| <b>Power fuse ratings</b>                 | 250V/5A time lag (2 fuses) 5 mm × 20 mm  |
| <b>Environmental operating range</b>      | +39 °F - 86 °F / +4 °C - 30 °C<br>No direct sunlight<br>Humidity 20 to 70% RH, noncondensing   |
| <b>Sample temperature and environment</b> | Controlled to user-selected temperature between 16 and 40 °C, but at least 10 °C above ambient temperature<br>No gas or humidity control |
| <b>Data interface</b>                     | TCP/IP (external)<br>USB type A (one in front, two in back)  |
| <b>Software OS</b>                        | Windows 10 embedded  |

**1 Introduction**  
Technical Specifications

## 2

# Installation

Unpacking and Component Identification 9



Suitable Locations for the XFp Analyzer 10

This chapter provides unpacking and installation instructions for the Agilent Seahorse XFp Analyzer.

## Unpacking and Component Identification

After receiving the package, immediately check each box for damage. Report any shipping damage to the transportation company, and Agilent using **“Technical Support and Ordering Information”** on page 64. **Table 1** shows the Seahorse XFp Analyzer system components.

**Table 1** Seahorse XFp Analyzer system components

| Instrument/Controller                           | Quantity | Image  |
|---|----------|--|
| XFp Analyzer                                    | 1        |  |
| Power cord<br>(region-specific)                 | 1        |  |
| Alternate wireless adapter<br>(p/n S7802-80000) | 1        |   |

## 2 Installation

### Suitable Locations for the XFp Analyzer

# Suitable Locations for the XFp Analyzer

The XFp Analyzer is designed for laboratory use. The internal environment for the XFp Analyzer cell culture miniplate is controlled to a preset tray temperature. The laboratory room temperature must therefore be maintained within the range listed in **“Technical Specifications”** on page 7. The miniplate tray temperature can be monitored using the status display in the upper right portion of the user interface.

The XFp Analyzer uses optical detection technology to measure extremely low levels of fluorescent emission from analyte sensors. Although the instrument has been designed to shield room light, avoid excessive light (such as direct sunlight).

## 3

# Basic Operations

Power and Warm Up 11

XFp Analyzer Communications and Network Connection Setup 12

Turning the XFp Analyzer Off 12

This chapter provides basic operating procedures for the Agilent Seahorse XFp Analyzer.

## Power and Warm Up

### Power up

To power on the XFp Analyzer:

- 1 Press the power switch on the back of the instrument. (**Figure 1**)



Figure 1. XFp Analyzer power switch.

- 2 Allow at least 20 minutes for the instrument to fully warm and equilibrate to the set temperature.
- 3 Check the status icon in the upper right corner of the screen. When ready the tray temperature and a green check mark is displayed.

### Welcome screen on first time power up

During the first power up, a welcome screen displays initial diagnostic tests on the instrument. Follow the on-screen directions to run them.

Run these tests at any time using the Diagnostics menu selection on the instrument user interface. For more details, see **“Navigating the XFp Analyzer”** on page 13.

## XFp Analyzer Communications and Network Connection Setup

The USB connectors on the instrument transfer template and result files between the XFp Analyzer and a computer running Seahorse Analytics. Agilent recommends connecting the XFp Analyzer to a local network directory to ease file movement and aid Cell Analysis Technical Support in servicing the instrument. Please contact Cell Analysis Technical Support, **“Worldwide technical support”** on page 64, with any questions.

For instructions on setting up a wired network connection, see **“Wired Network Setup”** on page 53. For instructions on setting up a wireless network connection using the included wireless adapter, see **“Wireless Network Setup”** on page 55.

#### NOTE

**Only the included wireless adapter has been qualified to work with the XFp Analyzer.**

## Turning the XFp Analyzer Off

To turn off the instrument, press the power button in the lower-left corner of the Home screen to first shut down the XFp Analyzer screen. After the screen turns black, use the power switch on the back of the instrument to completely shut down power.

#### NOTE

**Shut down the software before powering off using the rear power switch.**

## 4

# Navigating the XFp Analyzer

Setting Up an XFp Assay 14

Running an XFp Assay 19

Modifying XFp Analyzer Settings 24

XFp Analyzer System Files 25

Using Safe Lock (Trend Micro) 40

Cleaning and Routine Maintenance 41

Analyzing XFp Analyzer Data Using Seahorse Analytics 42

This chapter provides navigation information for the Agilent Seahorse XFp Analyzer.

## Setting Up an XFp Assay

### Select a template from local, USB, or network drive

Agilent provides default templates for most Seahorse XF assay kits. Templates can also be created in Wave desktop on a desktop or laptop computer and then transferred to the XFp Analyzer through a network connection or USB flash drive.

- 1 Click **Start** from the Home screen. See **Figure 2**.

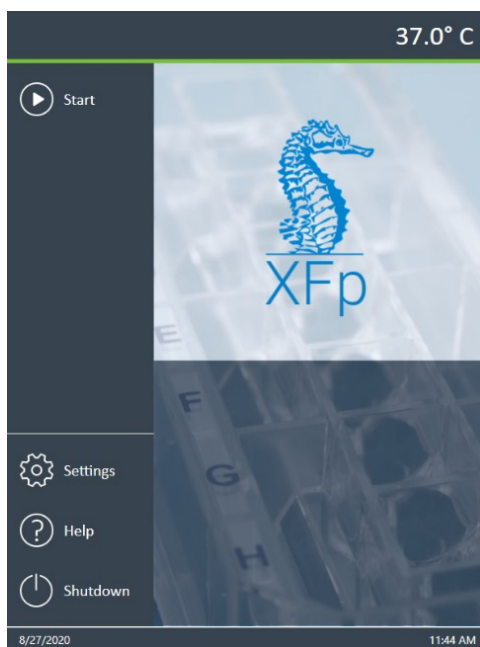


Figure 2. Home screen.

## 4 Navigating the XFp Analyzer

### Setting Up an XFp Assay

- 2 Select a template provided by Agilent to perform a specific assay. These templates can be found in the **Local** tab. See **Figure 3**.

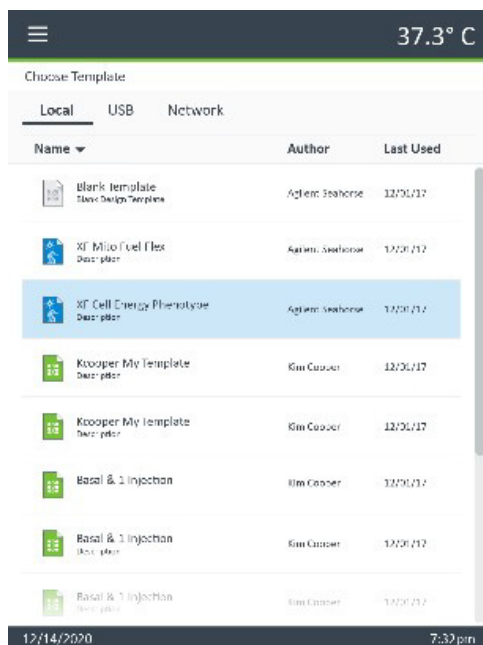


Figure 3. Local tab.

- 3 Templates designed in Wave can be transferred to the XFp Analyzer through a USB or network location.
  - If the instrument is networked, click the **Network** tab, select and run the template from the networked location.
  - If a USB drive containing a valid XFp Analyzer template is inserted into the instrument, click the **USB** tab, select and run the template directly from the USB drive.

#### NOTE

**Only XFp Analyzer assay template files (\*.asyt) are recognized by the XFp Analyzer. Templates must reside on the root of the USB flash drive and not in a subfolder.**

#### Verify groups and plate map

After selecting the template file, make any necessary modifications to the plate map and groups to be analyzed.

## 4 Navigating the XFp Analyzer

### Setting Up an XFp Assay

- 1 To see the conditions defined for a group, click the group name and look at the header information, as shown in **Figure 4**.

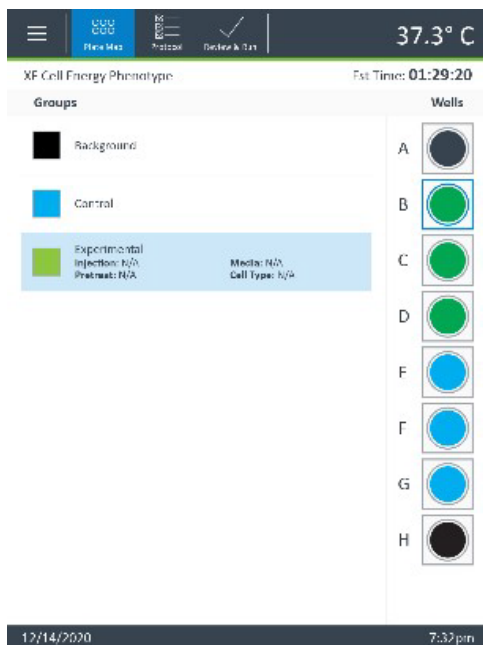


Figure 4. Groups.

- 2 To change the wells assigned to each group, click the group name followed by the well or wells to be included.
- 3 For all other changes to **Groups**, use Seahorse Analytics.

### Review/edit instrument protocol

- 1 Check that all desired steps are selected, check boxes checked (**Figure 5** on page 17). Agilent strongly recommends performing equilibration for all cell-based XF assays.
- 2 If needed, increase or decrease the number of measurements that will be performed during the assay on the **Instrument Protocol** page.
  - a Click the number in the cycle column that corresponds with the step to modify.

## 4 Navigating the XFp Analyzer

### Setting Up an XFp Assay

b Use the arrows to adjust the number of cycles up or down. See **Figure 5**.

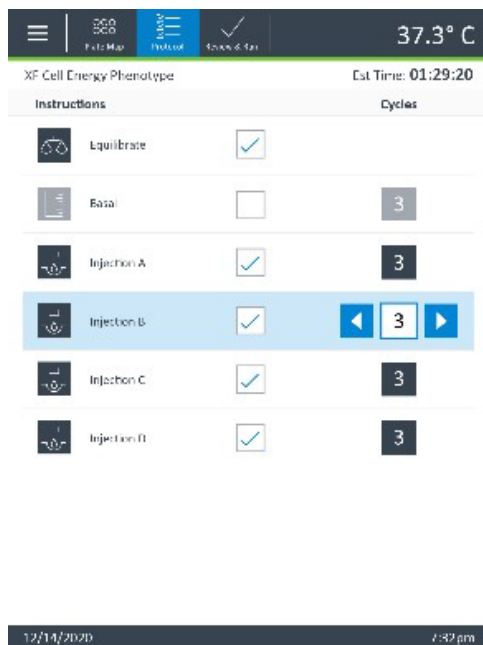


Figure 5. Instrument protocol page.

## 4 Navigating the XFp Analyzer

### Setting Up an XFp Assay

#### Review summary and start assay

- 1 Review the “**Safety Considerations**” on page 5 to verify the desired settings.
- 2 Before beginning an assay, the following optional steps may be taken:
  - a Click the **Assay Result File Name** text box to customize the name of the assay result file. See **Figure 6**.

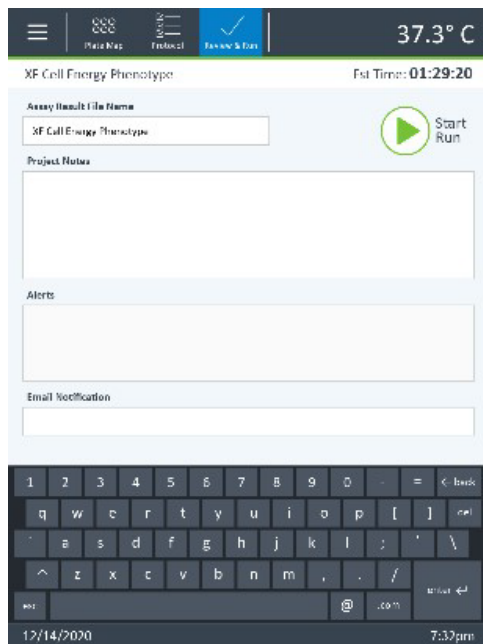


Figure 6. Assay name.

- b Click **Project Notes** to add any notes related to the assay or protocol being performed.
  - c Click **Email Notification** to add email addresses for recipients to be notified when the assay is complete (requires an active network connection).
- 3 Click **Start Run** then follow the onscreen instructions.

#### NOTE

The XFp Analyzer will save assay results to a USB flash drive by default (if available). Agilent recommends inserting a USB flash drive into the instrument at the beginning of assay setup so that results will automatically be saved. If the XFp Analyzer does not detect a USB flash drive after completing an assay, the system will prompt the user to insert a USB to save their assay result file.

## Running an XFp Assay

- 1 Load the cartridge (hydrated and loaded with compounds) and utility plate onto the tray when prompted.
- 2 Ensure the cartridge fits properly on the utility plate, the lid is removed from the cartridge, and the direction of the cartridge matches the image on the screen. See **Figure 7**.



Figure 7. Load cartridge screen.

- 3 The XFp Analyzer will perform calibration of the sensor cartridge. This takes approximately 20 minutes.

### NOTE

**Although the sensor calibration may appear complete after 15 minutes, the instrument will be busy for a few more minutes. Wait for the user prompt to appear.**

- 4 The XFp Analyzer tray will open and present the utility plate. Remove the utility plate and load the cell plate. Ensure the lid is removed from the cell plate before loading onto the XFp Analyzer tray.

## 4 Navigating the XFp Analyzer

### Running an XFp Assay

- 5 Click **Continue** to begin the equilibration step. The progress bar in the upper-left area of the screen displays the progress of the equilibration step.
- 6 Click either the **Overview** or **OCR vs. ECAR** tab to toggle between the views during the assay. See **Figure 8**. Overview and OCR versus ECAR are the two run-time views available as data is acquired by the XFp Analyzer.

### Overview

The **Overview** tab displays both OCR and ECAR as a function of time. Red vertical lines indicate the injections and are labeled by injection port letter. See **Figure 8**.

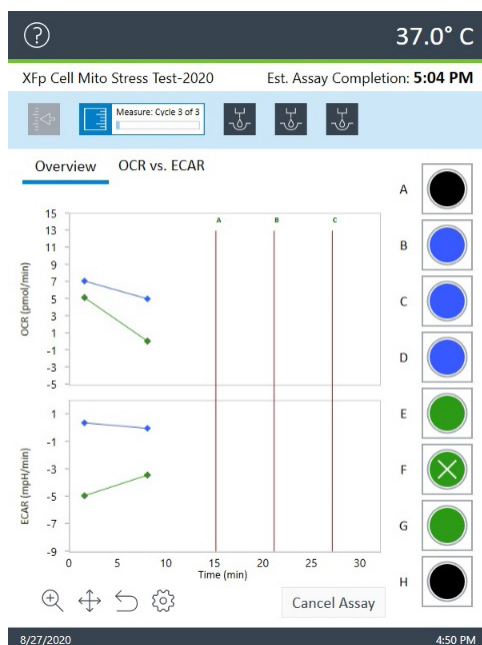


Figure 8. Overview tab.

In this view the charts can be zoomed and scaled by choosing one of the tools below the charts:

- Zoom
- Move
- Reset

## 4 Navigating the XFp Analyzer

### Running an XFp Assay

- Options. By default, rate data is displayed in **Group** mode. Click the Options icon to change display modes and turn on error bars. See **Figure 9**.



Figure 9. Options icon.

- **Show StdDev** turns on display of error bars for each **Group**.
- **Level Data** shows the concentration data from which the rate data is derived.
- **Well Mode** shows the data from the individual wells instead of the average of the wells in each group.

### OCR versus ECAR

The **OCR vs. ECAR** tab displays OCR on the Y-axis and ECAR on the X-axis.

Choose a single time point at which to examine OCR versus ECAR using the **Measurement** selector below the chart in **Figure 10**. Data is displayed in Group mode with standard deviations.

The scale of this chart is fixed to allow easy comparison of values across measurements. Quadrants are labeled **Aerobic**, **Energetic**, **Glycolytic**, or **Quiescent** to show the energy preference of the cell during the assay at a given measurement. See **Figure 10**.

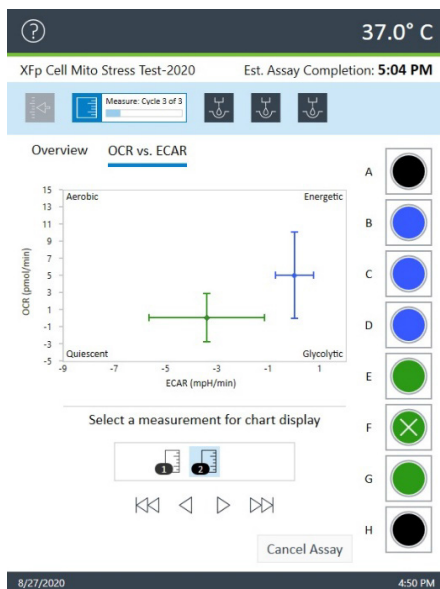


Figure 10. OCR vs. ECAR tab.

## 4 Navigating the XFp Analyzer

### Running an XFp Assay

#### NOTE

In both the Overview and OCR vs. ECAR views, the display of individual wells can be turned off by touching the well graphic on the right. Background correction is always performed while the experiment is running.

- 7 Once the assay has finished, the XFp Analyzer prompts the user to eject the cartridge and remove the cartridge and cell plate. See [Figure 11](#).



Figure 11. Remove cartridge and cell plate prompt.

- 8 After the assay is complete, the data is automatically saved to a USB flash drive (if inserted), the local drive, or the networked drive (if configured). A message appears indicating the location of the save assay result File (\*.asyr). See [Figure 12](#).

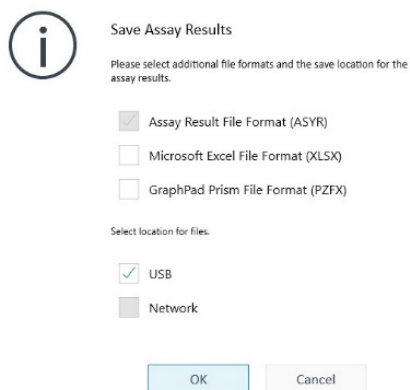


Figure 12. Save results.

## 4 Navigating the XFp Analyzer

### Running an XFp Assay

#### NOTE

Canceling during an assay may cause the XFp Analyzer to stall. If this occurs, shut down the instrument by pressing the power switch on the back. Use the same power switch to turn the instrument ON.

- 9 After the assay is complete, data is automatically saved on the assay result file format to the location specified at the start of your assay (USB flash drive, network location, or locally). It is mandatory to save result data in the assay result file format after completing the assay and will always be checked ON. You can also save result data as an Excel file (.xlsx) and GraphPad Prism file (.pzfx) directly from the XFp Analyzer as well. Check the box next to the desired file formats and click **OK** (Figure 12 on page 22).

Assay result files saved locally on the XFp Analyzer will be stored for 60 days after the assay. Saving assay result files to a USB or network drive is encouraged.

## Modifying XFp Analyzer Settings

From the Start page, click **Settings** to access the XFp Analyzer settings, see **Figure 13**. The XFp Analyzer settings page contains:

- System Files
- Diagnostic Tests
- System Settings
- Assay Results
- Template Management

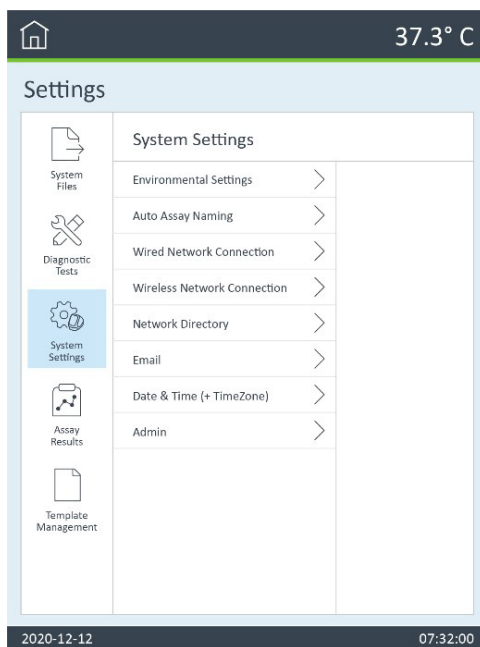


Figure 13. XFp Analyzer settings page.

# XFp Analyzer System Files

## System files

To access the **System Files** page, go to the Home page and select **Settings > System Files**. (Figure 14) Cell Analysis Technical Support routinely requests that System Files be sent if an XFp Analyzer encounters an error.

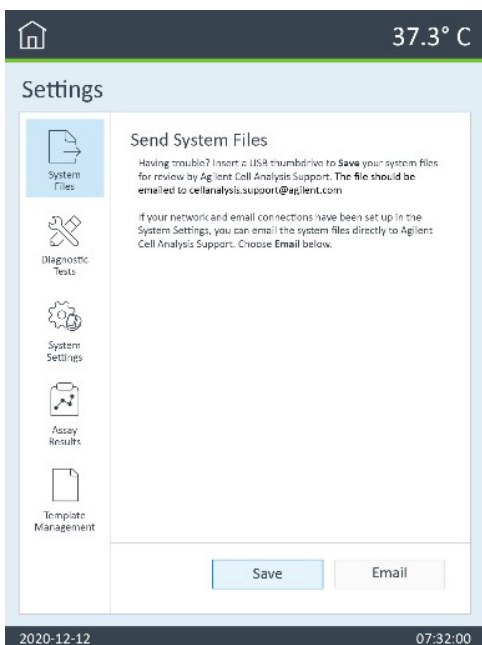


Figure 14. Send system files.

### Option 1 - Send system files to a USB flash drive

Save System Files to a USB flash drive as a compressed (zipped) folder. This can then be attached to an email to Cell Analysis Technical Support. **Save** will become active after a USB is connected to one of the three USB ports. (Figure 14)

#### NOTE

The file will be named: **Diagnostic Report\_Date\_Time** (Example: **Diagnostic Report\_2020\_05\_19\_15\_34\_14.zip**).

## 4 Navigating the XFp Analyzer

### Diagnostics tests

#### Option 2 - Send system files directly to Cell Analysis Technical Support

The diagnostic report compressed folder may also be emailed directly to Cell Analysis Technical Support from the XFp Analyzer. The XFp Analyzer must have an active wired or wireless network connection to directly send system files to Cell Analysis Technical Support.

## Diagnostics tests

The XFp Analyzer **Diagnostic Tests** page contains various functions that assist Cell Analysis Technical Support in debugging any issues or errors encountered during normal operation. Perform the **System Check** diagnostic function at the initial start of the XFp Analyzer and when Cell Analysis Technical Support request.

To access the **Diagnostic Tests** page, go to the Home page and select **Settings > Diagnostic Tests**. (Figure 15)

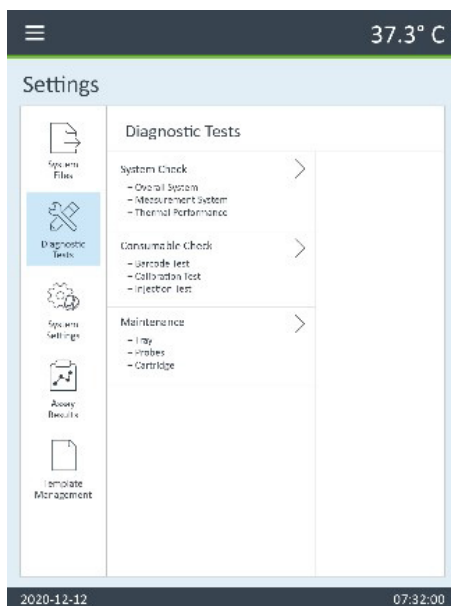


Figure 15. Diagnostic tests.

The Diagnostic Tests page has the following sections:

- System Check
- Consumable Check
- Maintenance

## 4 Navigating the XFp Analyzer

### Diagnostics tests

#### System check

##### Overall system

- **Communication test** - Verifies all system modules are communicating appropriately.
- **Voltage test** - Ensures voltage levels are within specified ranges for operation.
- **Motor test** - Verifies the movements and positions of the tray and probes.
- **Holding pressure test** - Ensures the injection system does not have any leaks.
- **Injection zero test** - Ensures function of the zero pressure sensor.
- **Injection solenoid test** - Verifies proper open/close operation of internal valves.

##### Measurement system

- **LED reference test** - Verifies that the system records LED intensity signal when ON.
- **Autozero test** - Analyze and record any offset values found within the channels on each board.
- **Noise test** - Analyze and record the amplitude of the noise level on each channel.
- **High-Gain rise time test** - Conduct an amplifier rise time test.

##### Thermal performance

- **Thermal test** - Verifies that the temperature remains within tolerance after a "Wait" period.

##### Consumable check

- **Barcode test** - Verifies XFp Analyzer can read barcodes (requires user interaction to complete).
- **Calibration test** - Performs calibration procedure and will verify that all wells are calibrated correctly.
- **Injection test** - Verifies proper performance of injection step.

## 4 Navigating the XFp Analyzer

### Diagnostics tests

#### Maintenance

The XFp Analyzer is designed to require minimal cleaning and maintenance. All consumables are disposable, and none of the instrument components are exposed to the cell plate at any time, preventing cross-contamination of biological or chemical materials. However, it may be necessary to perform some tasks. (Figure 16)

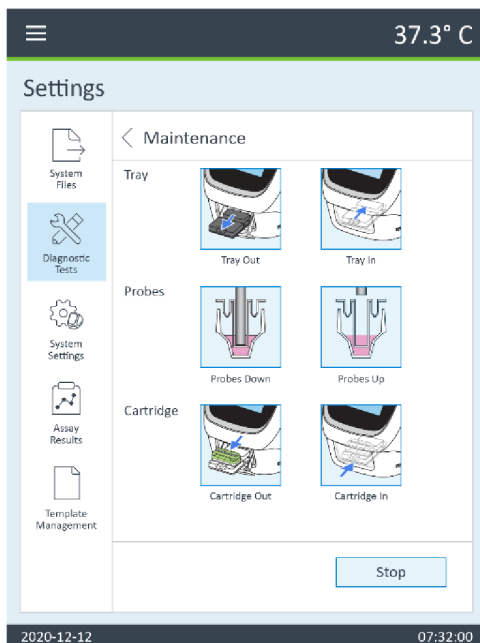


Figure 16. Maintenance window.

- **Tray** - Press **Tray Out** to open the door and eject the tray. Press **Tray In** to return the tray inside the XFp Analyzer.
- **Probes** - Probe movement will adjust the probes within the XFp Analyzer up or down.
- **Cartridge** - Use this function to take a **Cartridge Out** or put a **Cartridge In** the XFp Analyzer.

#### NOTE

If a cartridge is suspected to be inside the XFp Analyzer, use this function to remove the cartridge before beginning another assay.

## 4 Navigating the XFp Analyzer

### System settings

# System settings

The XFp Analyzer System Settings page contains various functions that permit assay customizations and IT integration options.

To access the **System Settings** page, go to the Home page and select **Settings > System Settings**. (Figure 17)

The **System Settings** page has the following sections:

- Environmental Settings
- Auto Assay Naming
- Wired Network Connection
- Wireless Network Connection
- Network Directory
- Email
- Date and Time (plus TimeZone)
- Admin

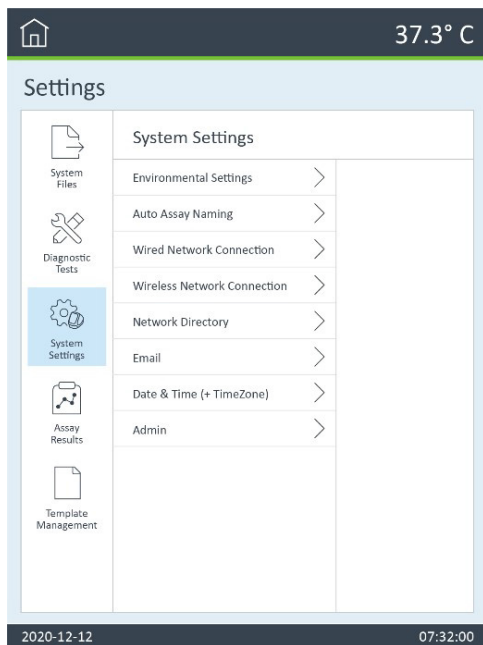


Figure 17. System Setting page.

## 4 Navigating the XFp Analyzer

### System settings

#### Environmental settings

Temperature target: Adjust the tray temperature on the XFp Analyzer. See **Figure 18**.

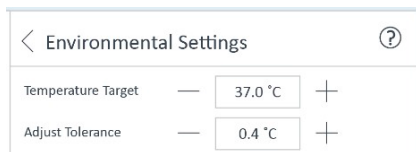


Figure 18. Temperature target setting.

Temperature Control on XFp Analyzer: Users may adjust the tray temperature of the XFp Analyzer for a wider array of assays and applications. The validated tray temperature range of the XFp Analyzer is 16 to 40 °C. See **Figure 19** for complete details of the environmental settings required to accurately achieve the desired tray temperature setting.

Temperature specifications: Users can change the sample (tray) temperature to 16 to 40 °C; there is no internal cooling function.

- Lower tray temperatures require the instrument to be placed in a cold room or refrigerated unit that is 8 to 20 °C cooler than the desired sample tray temperature.
- Users can set the tray temperature to any value (in increments of 0.1 °C) within the allowed range. **Figure 19** shows the range of supported ambient and sample tray temperatures.

| XFp 1.1                    |    | sample (tray) temperature |    |    |    |    |    |    |    |    |    |    |    |    |  |  |  |  |  |  |
|----------------------------|----|---------------------------|----|----|----|----|----|----|----|----|----|----|----|----|--|--|--|--|--|--|
| ambient (room) temperature | 4  | 16                        | 18 | 20 | 22 | 24 |    |    |    |    |    |    |    |    |  |  |  |  |  |  |
|                            | 6  | 16                        | 18 | 20 | 22 | 24 | 26 |    |    |    |    |    |    |    |  |  |  |  |  |  |
|                            | 8  | 16                        | 18 | 20 | 22 | 24 | 26 | 28 |    |    |    |    |    |    |  |  |  |  |  |  |
|                            | 10 |                           | 18 | 20 | 22 | 24 | 26 | 28 | 30 |    |    |    |    |    |  |  |  |  |  |  |
|                            | 12 |                           |    | 20 | 22 | 24 | 26 | 28 | 30 | 32 |    |    |    |    |  |  |  |  |  |  |
|                            | 14 |                           |    |    | 22 | 24 | 26 | 28 | 30 | 32 | 34 |    |    |    |  |  |  |  |  |  |
|                            | 16 |                           |    |    |    | 24 | 26 | 28 | 30 | 32 | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 18 |                           |    |    |    |    | 26 | 28 | 30 | 32 | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 20 |                           |    |    |    |    |    | 28 | 30 | 32 | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 22 |                           |    |    |    |    |    |    | 30 | 32 | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 24 |                           |    |    |    |    |    |    |    | 32 | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 26 |                           |    |    |    |    |    |    |    |    | 34 | 36 | 37 | 40 |  |  |  |  |  |  |
|                            | 28 |                           |    |    |    |    |    |    |    |    |    | 36 | 37 | 40 |  |  |  |  |  |  |
| 30                         |    |                           |    |    |    |    |    |    |    |    |    | 37 | 40 |    |  |  |  |  |  |  |

Figure 19. Ambient room temperature (yellow) required to achieve the desired XFp Analyzer tray temperature setting (green).

## 4 Navigating the XFp Analyzer

### System settings

#### Cartridge hydration and equilibration steps for temperatures of 28 °C and below

##### Day before Assay:

Hydrate the XFp Cartridge in a non-CO<sub>2</sub> 37 °C incubator overnight.

##### Day of Assay:

The utility plate + cartridge require 40 minutes of equilibration time inside the XFp when running at temperatures below 20 °C. Approximately 40 minutes before starting the assay:

- 1 Place the cell plate in an incubator at the desired XFp operational temperature.
- 2 Bring the hydrated cartridge loaded with compounds and paired with the utility plate (with calibrant solution) to the XFp Analyzer.
- 3 Click **Diagnostics**, then click **Maintenance**.
- 4 Press **Tray Out**.
- 5 Once the tray fully extends, remove the lid from the cartridge and place the utility plate and cartridge on the tray.
- 6 Press **Tray In**, then click the lower-left arrow twice to return to the Home screen.
- 7 Wait 40 minutes.
- 8 Click **Start** on the home screen, then select the **Assay Template** to run.
- 9 Click the lower-right arrow twice to navigate to the **Summary** screen, then click **Start Assay**.
- 10 The **Load Cartridge** message appears and the tray extends with the utility plate and cartridge. *Do not* remove the utility plate and cartridge (ensure the lid is removed from the cartridge).
- 11 Click **Continue** to start calibration.
- 12 After completing calibration, the XFp will prompt the user for the cell plate. Load the cell plate and begin assay.

#### Returning the XFp Analyzer to room temperature after cold room use

To use the XFp Analyzer at room temperature after use in the cold room, complete the following steps:

## 4 Navigating the XFp Analyzer

### System settings

- 1 Unplug the analyzer and move it to the desired location.
  - 2 Leave the instrument unplugged in the destination overnight.
  - 3 Resume normal use.
    - a Adjust tolerance:** The temperature tolerance can be set from 0.2 to 9.9 °C. The tolerance defines the acceptable range in which the tray temperature may deviate from the target tray temperature.
    - b Temperature alarm:** When enabled, the XFp Analyzer will automatically notify users (**Figure 20**) when the tray temperature of the XFp Analyzer is out of the acceptable range, as specified by the temperature tolerance above (only while running an assay on the XFp Analyzer).
- Example: At a target tray temperature of 37 °C and a tolerance of 0.3 °C, the acceptable tray temperature range is 36.7 to 37.3 °C.
- c Atmospheric pressure:** Customize the atmospheric pressure set on the XFp Analyzer to match that of the lab where the instrument is located. This value must be entered manually, the local atmospheric pressure is not automatically detected by the XFp Analyzer.

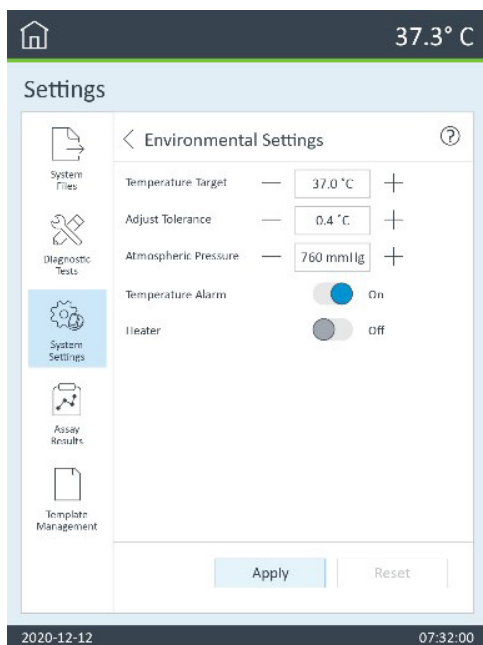


Figure 20. Temperature alarm window.

## 4 Navigating the XFp Analyzer

### System settings

#### Auto assay naming

Auto assay naming allows users to specify the default name for each Assay Result File (\*.asyr) that is created on the XFp Analyzer. See **Figure 21**.

- 1 Adjust the order of each variable by dragging and dropping the selection in the order desired. **Figure 21** shows the configuration (template name) - (Timestamp).
- 2 The **Custom** field can be used to add a keyword, instrument name, or other items to all assay results files.
- 3 Click **Apply** to complete the process of setting up a custom name.

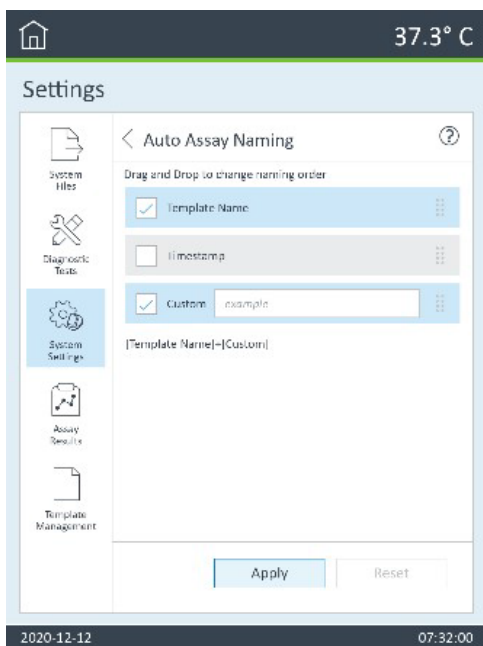


Figure 21. Auto assay naming window.

## 4 Navigating the XFP Analyzer

### System settings

#### Wired network

An ethernet cable is required to set up wired network access. The ethernet cable port is located on the lower backside of the XFP Analyzer.

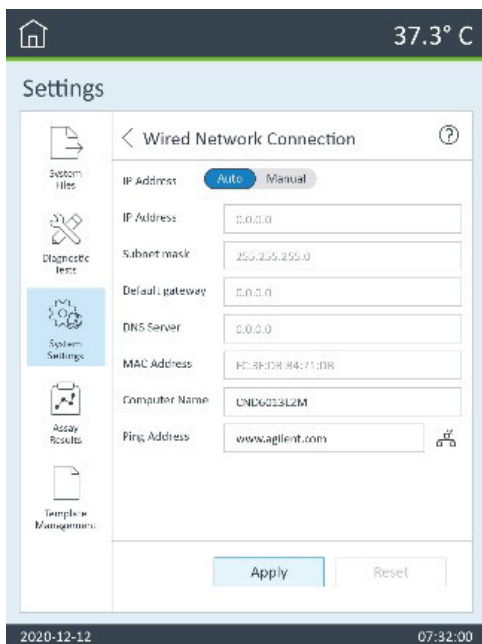


Figure 22. Wired network.

#### NOTE

Agilent recommends that the device settings be configured by the institution IT department using the “[XFP Analyzer Network Setup](#)” on page 47 as a reference.

#### Wireless network

A wireless USB adapter can be plugged into one of the available USB ports (recommended on lower back side of XFP Analyzer) to gain wireless access. Once finished, select an available **SSID** from the list and click **Connect**. If there are no visible networks available, click **Refresh**.

## 4 Navigating the XFP Analyzer

### System settings

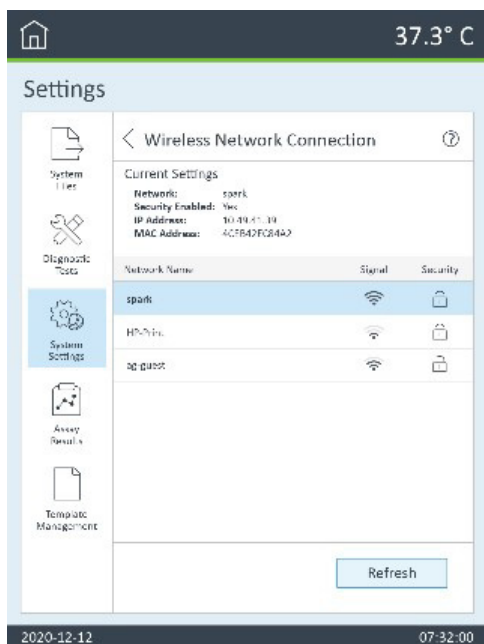


Figure 23. Wireless network.

### NOTE

The XFP Analyzer must be restarted after the wireless USB adapter is plugged in.

Saving assay template or result files to a shared network directory requires an active wired or wireless network connection.

### Network directory

Network configuration is covered in “[XFP Analyzer Network Setup](#)” on page 47.

### Email configuration

Configure an email account to automatically send 'Assay Complete' email notification as well as the Assay Result file to specified email addresses entered before starting an assay (see “[Review summary and start assay](#)” on page 18 for more information).

### Date and Time (plus Time Zone)

Adjust the date or time. This can be done automatically by connecting to a time server (requires an active network connection) or set manually.

## 4 Navigating the XFp Analyzer

### Assay results

Set the time zone for the XFp Analyzer. Once set, the correct time will be visible on the screen as well as when using the time stamp function for auto assay naming.

### Admin

Exit and re-enter Kiosk mode to perform Windows Updates or install security software such as Trend Micro's Safe Lock application. The Using Safe Lock section provides the steps to reach the Windows desktop and return to XFp application.

## Assay results

The XFp Analyzer Assay Results page contains a copy of each assay performed within the previous 60 days on the local storage within the XFp Analyzer. After 60 days, Assay Result files are deleted from the XFp Analyzer.

To access the **System Settings** page, go to the Home page, then select **Settings > Assay Results**.

- 1 **Assay results:** Individual assay results can be viewed by clicking the small check box next to the result file to view and click **View**. See [Figure 24](#).

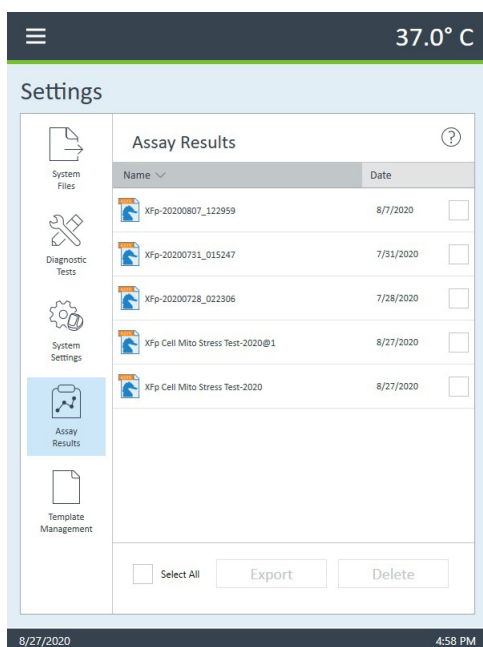


Figure 24. Assay results view.

## 4 Navigating the XFp Analyzer

### Template management

- 2 Export options:** Assay result files may be exported to Microsoft Excel or GraphPad Prism file formats (requires a USB flash drive or shared network directory). See **Figure 25**. To export these files (either individually or collectively) to a network location or USB, click the individual check box or click **Select All**, and then click **Export**.

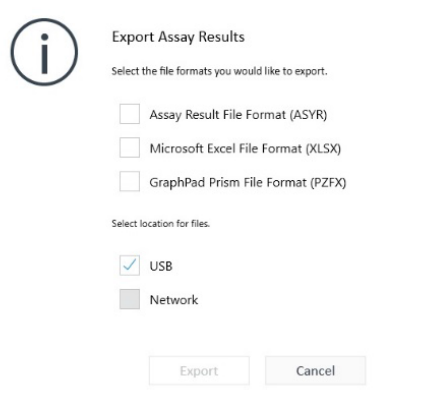


Figure 25. Export options dialog box.

- 3 Delete:** To remove assay result files from the XFp Analyzer, click the check boxes next to all result files to remove, and click **Delete**.

## Template management

Template management provides a simple way to import or export assay template files to/from the XFp Analyzer.

### Import assay template file (or files):

- 1** First, transfer assay template file to a USB flash drive or shared network directory location.
- 2** On the XFp Analyzer, select the location where the assay template file will be imported from (insert the USB flash drive, then select the tab when it appears).
- 3** Check the box next to an individual assay template, or click **Select All** to select all the assay template files.
- 4** To import assay template files to the **Local** tab, click **Import**.
- 5** A successful assay template import results in an **Import Complete** message, click **OK**.

## 4 Navigating the XFp Analyzer

### Help

#### Export assay template files:

- 1 On the **Local** tab, click the check box next to the assay template file to be exported.
- 2 Select the desired location to export the assay template file (USB flash drive or shared network directory).
- 3 If the assay template has exported successfully, an **Export Complete** message will appear. Click **OK**.

## Help

From the Home page Click **Help**. The Help page has the following sections:

- Support (see “**Support**” on page 65)
- Software version (see “**Support**” on page 65)
- Getting started
- License agreement (see “**Support**” on page 65)
- Remote assistance

#### Getting started

Provides useful tips and information regarding the setup and operation of the XFp.

## 4 Navigating the XFp Analyzer Help

### Remote assistance

Cell Analysis Technical Support has the ability to remotely access each XFp Analyzer that has an active network connection.

To access the Remote Assistance page, go to the Home page, then select **Help > Remote Assistance**. See **Figure 26**.

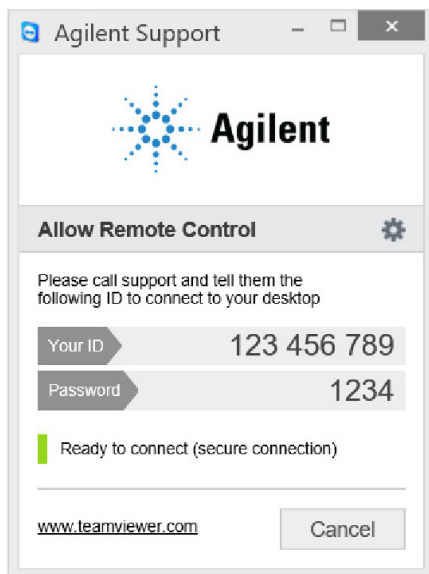


Figure 26. Remote assistance window.

- 1 To begin a remote session with Cell Analysis Technical Support, call the appropriate geographically located support line then click **Start**.
- 2 An ID and password will be presented, Cell Analysis Technical Support will request this information to connect to the XFp Analyzer.

## 4 Navigating the XFp Analyzer

### Using Safe Lock (Trend Micro)

# Using Safe Lock (Trend Micro)

The computer security application Safe Lock by Trend Micro has been validated on the XFp Analyzer. Safe Lock provides several levels of security while allowing the XFp Analyzer to perform its functions. More information can be found at [www.trendmicro.com](http://www.trendmicro.com). To reach the Windows 10 desktop to install application, follow the steps below:

- 1 From the Home page tap **Settings > System Settings > Admin**.
- 2 Click **Exit Kiosk Mode**.
- 3 The Windows 10 desktop is now available
- 4 Follow the installation instructions for the Safe Lock application
- 5 Once all the steps and configurations are complete, double-click the **Exit Maintenance Mode** icon on the XFp Desktop
- 6 The system will reboot and start the XFp controller application

## Cleaning and Routine Maintenance

The XFp Analyzer is designed to require minimal cleaning and maintenance. All consumables are disposable and none of the instrument components are exposed to the cell plate at any time, preventing cross-contamination of biological or chemical materials.

### Air filter replacement

At approximately one-year intervals, the air filter at the upper rear of the analyzer should be replaced. The filter retainer is held in place by magnets and can be pulled away from the housing by hand. (Replacement filters can be ordered using part number: 102799-000.)

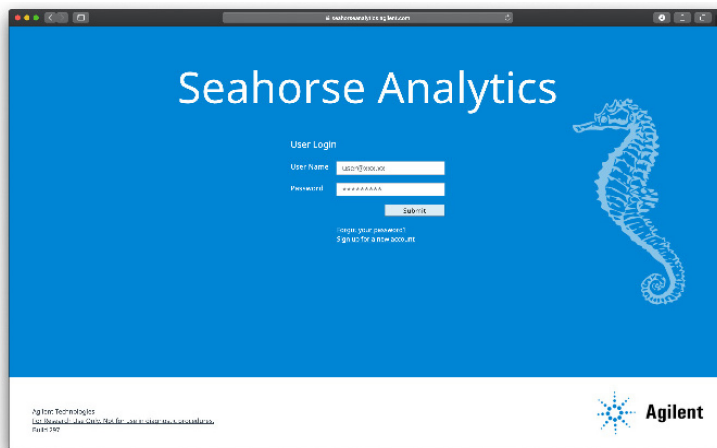
## 4 Navigating the XFp Analyzer

### Analyzing XFp Analyzer Data Using Seahorse Analytics

# Analyzing XFp Analyzer Data Using Seahorse Analytics

Agilent Seahorse Analytics is a web-based analysis and file management software application for XFp assay data. Use the integrated assay kit companion analysis views to quickly report and share result data, or create a custom analysis view using the 30+ graphing options available. Create a Seahorse Analytics account: <https://seahorseanalytics.agilent.com> Basic data analysis:

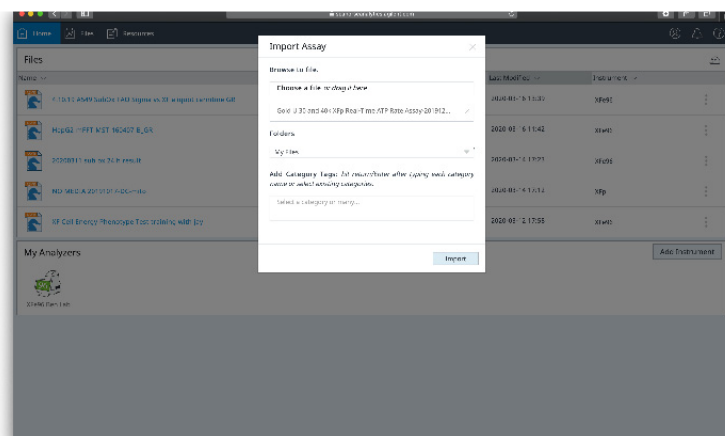
- 1 After the assay is completed, transfer your assay result file to your personal computer using a USB drive or network drive.
- 2 Go to <https://seahorseanalytics.agilent.com> to register or log in to your Seahorse Analytics account.



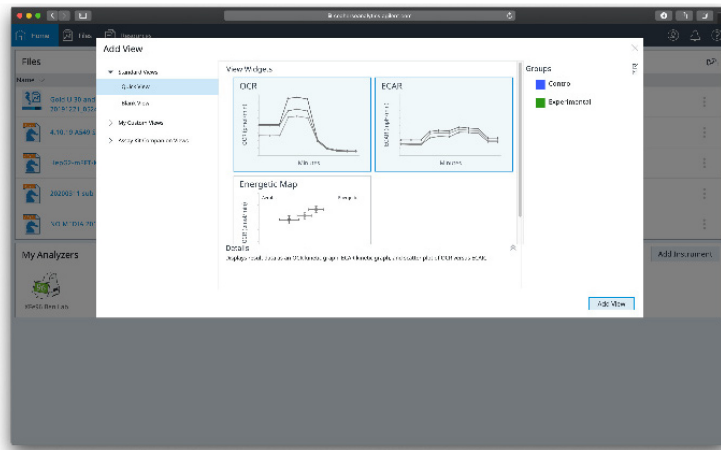
## 4 Navigating the XFp Analyzer

### Analyzing XFp Analyzer Data Using Seahorse Analytics

- 3 From the home view, import your XFp assay result file to your account.



- 4 Click the file to open it. You will be presented with various analysis view options that you can apply to your data file, including kinetic graphs or bar charts as well as analysis specific for an assay kit.

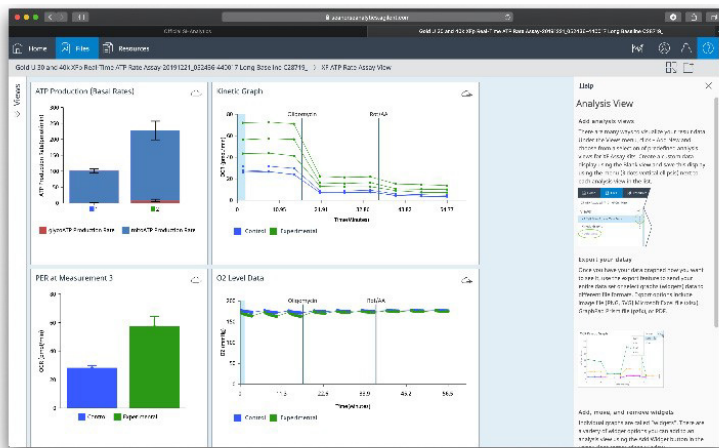




## 4 Navigating the XFp Analyzer

### Analyzing XFp Analyzer Data Using Seahorse Analytics

- 6 Use the embedded help seen on each software screen to learn about other features available.



## 4 Navigating the XFp Analyzer

Analyzing XFp Analyzer Data Using Seahorse Analytics

# 5

## XFp Analyzer Network Setup

|                                 |    |
|---------------------------------|----|
| Introduction                    | 48 |
| XFp Analyzer System Information | 51 |
| Wired Network Setup             | 53 |
| Wireless Network Setup          | 55 |
| Network Checklist               | 62 |

This chapter provides network setup information for the Agilent Seahorse XFp Analyzer.

## Introduction

### Why should I network my XFp Analyzer?

Connecting your XFp Analyzer to a network provides unique workflow advantages to users, such as:

- Immediate acquisition of Assay Results.
- Simple File Transfer for Assay Templates and Results, no need for USB drives.
- Remote access for Cell Analysis Technical Support.

The XFp Analyzer supports two types of network connections:

- **Wide Area Network (WAN):** A computer network that covers a broad area, such as the internet.
- **Local Area Network (LAN):** Typically found in offices or schools, a computer network that connects multiple devices (computers, printers, lab devices, etc.) to enable file sharing between those networked devices in addition to internet access. LANs also have some level of security to access the shared file directories.

Both a WAN and LAN can be configured through a wired (Ethernet cable) or wireless (USB WiFi adapter) connection on the XFp Analyzer. Depending on your company's IT infrastructure, configuring WAN or LAN access may require additional support from your local IT department to complete set up.

Give the **"Network Checklist"** on page 62 to your IT department to use during the network setup of the XFp Analyzer.

5 XFp Analyzer Network Setup  
Why should I network my XFp Analyzer?



Figure 27. Networked XFp Analyzer.

## Network access features on the XFp Analyzer

### Immediate data delivery



Before starting an assay, users enter in one or multiple email addresses to receive a copy of the Assay Result file (\*.asyr) for data analysis. The email service also informs users when they can begin another assay.

### File transfer made simple



Use a shared network directory to simplify data transfer to and from the XFp Analyzer. A shared network directory allows users to easily transfer Assay Template files to the XFp Analyzer to perform an assay, then retrieve the Assay Results for analysis using **Agilent Seahorse Analytics**.

### Rapid remote assistance



An active network connection on the XFp Analyzer allows a Cell Analysis Technical Support representative to diagnose and troubleshoot potential issues quickly by allowing remote access to view and control the XFp Analyzer.

### Send 'System Files' directly to cell analysis technical support



When an issue is encountered, Cell Analysis Technical Support routinely requests System Files from the XFp Analyzer. System Files assist Cell Analysis Technical Support in identifying the root cause of an issue. With network access, users are able to send System Files directly to Cell Analysis Technical Support.

# XFp Analyzer System Information

## Required materials

- Approved wireless USB adapter or wired connection (Ethernet cable)
- IT department network connection settings

### NOTE

**XFp Analyzer instruments use Windows Defender with default settings. These include Quick Scan, real-time protection and cloud-protection turned on.**

- The XFp Analyzer can be connected to any Microsoft Windows compatible network and the Local Area connections can be configured as required by the network.
- Complete the **“Network Checklist”** on page 62 to ensure you have everything needed to successfully connect to a new network.
- The XFp Analyzer ships with an integrated 100Mbps Ethernet network adapter.
- The XFp Analyzer ships with a Netgear Wireless AC Adapter AC600 Dual Band (only the provided adapter is qualified for use). See **Figure 28**.

### NOTE

**The software required to use the USB WiFi Adapter is installed on the XFp Analyzer by default.**

- There is no internal WiFi adapter in the XFp Analyzer.
- An Ethernet cable is using a wired network setup. The Ethernet (RJ-45) jack is located at the base of the instrument in the back. See **Figure 29** on page 52.



**Part number S7802-80000**

Figure 28. USB WiFi adapter for the XFp Analyzer.

## 5 XFp Analyzer Network Setup

### XFp Analyzer System Information



Figure 29. Ethernet (RJ-45) jack location outlined in red.

## Wired Network Setup

### Wired connection

- 1 Plug the wired network connection (Ethernet cable) into the Ethernet port (**Figure 30**), located on the back of the XFp Analyzer. Ensure that the connection is firmly seated.



Figure 30. Ethernet port with wired network connection cable.

- 2 Power XFp Analyzer **OFF**.
- 3 Power XFp Analyzer **ON**.
- 4 Wait for the temperature to display in the upper right-hand corner of the home screen before proceeding with the Wired Setup instructions, see **"Option 2: Manual IP address assignment"** on page 54.
- 5 From the **Home** screen, click **Settings**.
- 6 Click **System Settings** to access the Wired Network Connection Settings screen.

There are two options depending on the network settings required by your facility's IT department:

#### **Option 1: Automatic IP address assignment (DHCP)**

Dynamic Host Configuration Protocol (DHCP) is enabled by default on the XFp Analyzer and automatically uses a DHCP server on the network to retrieve IP address values. The XFp Analyzer is set to obtain the IP address and subsequent information (Subnet mask, Default gateway, etc.). If these fields do not automatically populate, follow the steps in **"Option 2: Manual IP address assignment"** on page 54 or contact your local IT administrator.

## 5 XFp Analyzer Network Setup

### Wired connection

#### Option 2: Manual IP address assignment

- 1 In the Wired Network Connection screen, switch the toggle from **Auto** to **Manual**.
- 2 Manually enter the information provided by the local IT department in each field.
- 3 After setting the IP options, click **Save**.
- 4 Power XFp Analyzer **OFF**.
- 5 Power XFp Analyzer **ON**.
- 6 Once the temperature is displayed in the upper right-hand corner, click **Settings**, and then click **System Settings**.
- 7 To ensure you have a working connection, click **Ping**. A **Pass** message appears if the setting are working and the connection is active.



- 8 If the connection fails, reconfirm the IP address setting with the IT department, and confirm the information in the hardware setup steps.
- 9 Use the left and right arrows to continue editing instrument settings or exit the settings options.

## Wireless Network Setup

### Wireless connection

- 1 Plug the USB WiFi Adapter into one of the available USB ports on the back of the XFp Analyzer. See **Figure 31**.
- 2 Power XFp Analyzer **OFF**.
- 3 Power XFp Analyzer **ON**.
- 4 Wait for the temperature to display in the upper right-hand corner of the home screen before proceeding with the Wireless Setup instructions.



Figure 31. USB WiFi adapter inserted into an available USB port on the back of the XFp Analyzer.

### Joining a wireless network

- 1 From the Home screen click **Settings**.
- 2 Click **System Settings**.
- 3 Click the right arrow (bottom right) once to access the Wireless Network Connection Settings screen shown in **Figure 32** on page 56.
- 4 Select the network you want to connect to, and click **Connect**.

#### NOTE

If you do not see your wireless network on the list of available networks, click **Refresh**.

## 5 XFp Analyzer Network Setup

### Joining a wireless network

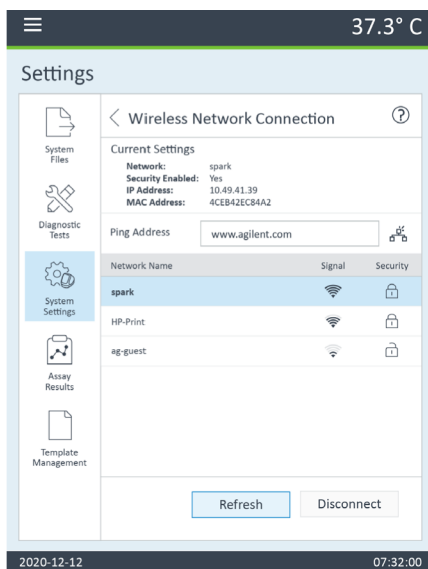


Figure 32. Example wireless network selection table.

- 5 Enter the password when prompted by the system.
- 6 If no errors are received, select the desired network again and verify that the connect button now displays disconnect.
- 7 Power XFp Analyzer **OFF**.
- 8 Power XFp Analyzer **ON**.
- 9 After the XFp Analyzer starts up and the temperature value in the upper right-hand corner is displayed, click **System Settings**.
- 10 Click **Ping** in the Wired Network Connection Settings screen.



- 11 If you received a **Pass** message and icon, your wireless setup is complete. If not, confirm the information in the wireless networking setup steps and reattempt connecting to a wireless network.
- 12 Use the left and right arrows to continue editing instrument settings or exit the settings options.

#### NOTE

**An active Internet connection is required for this feature.**

## Shared folder setup

This procedure requires an active wired or wireless network connection.

- 1 From the Home screen, click **Settings**, then **System Settings**.
- 2 Select the **Network Directory** tab.
- 3 Enter the information for the shared network drive into the dialogue boxes. Refer to the **“Shared directory settings”** on page 62 in the **“Network Checklist”** on page 62 for the correct information. (IT should provide this information.)
  - **Shared Directory:** The desired location on the LAN where all Assay Template and Assay Result files will be saved for access.
  - **Domain:** Name of the LAN.
  - **User Name:** User name of the Windows account that has read/write permission to access the shared directory location.
  - **Password:** Password for the Windows account.

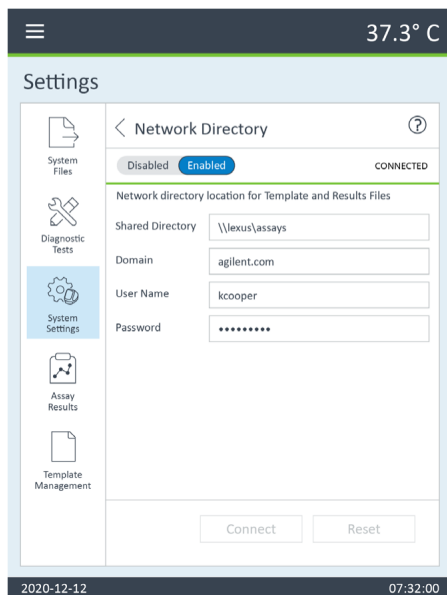


Figure 33. Example of a complete Network Directory configured on the XFp Analyzer.

### NOTE

**Accuracy is essential, double check your entry before checking access.**

## 5 XFp Analyzer Network Setup

### Email configuration

- Once all information has been entered, click **Enable**. Successfully configuring a LAN results in a **CONNECTED** message on the XFp Analyzer. (Figure 34)

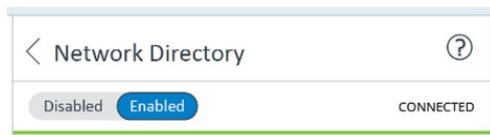


Figure 34. Successful network connection status.

- Use the left and right arrows to continue editing instrument settings or exit the settings options.

## Email configuration

- From the Home screen, click **Settings**, then **System Settings**.
- Select the **Email Configuration** tab. (Figure 35).

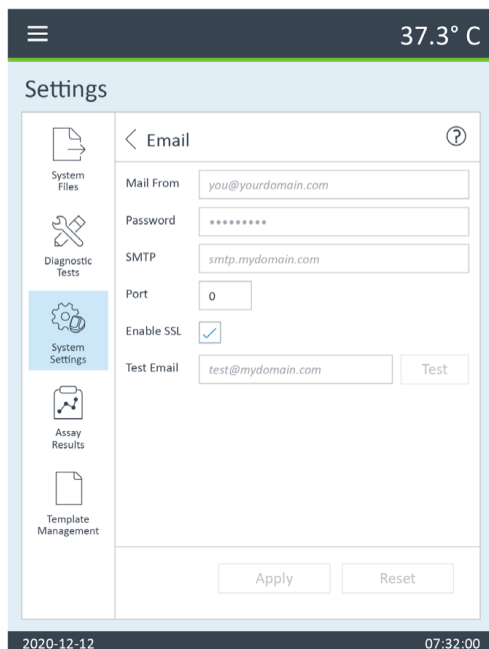


Figure 35. Example of a completed Email Configuration setup.

## 5 XFp Analyzer Network Setup

### Email configuration

- 3 Enter the information below:
  - **Mail From:** The email address that will send Assay Result files to email recipients.
  - **Password:** Password for the email address account in the **Mail From** field.
  - **Port:** Contact your local IT administrator for the correct port.
  - **Enable SSL:** Typically required by email providers, SSL protects data transmission between devices.
- 4 Verify the email configuration is completed by entering an email address to send a test email message from the XFp Analyzer. Click **Test** once an email address has been entered into the field. If an email is not received, ensure the information provided is correct.
- 5 Use the left and right arrows to continue editing instrument settings or exit the settings options.

### Time zone setup

Time can be set manually, or if you are connected to a network, it can be auto set. First, you must choose a Time Zone.

- 1 From the Home screen click **Settings**, then click **Go to Setup**.
- 2 Select the **Date & Time** tab. (**Figure 36** on page 60)
- 3 Select the **Time Zone** from the list provided.

Optional: Check the box 24 Hour Clock to activate the 24 hour clock setting.

## 5 XFp Analyzer Network Setup

### Email configuration

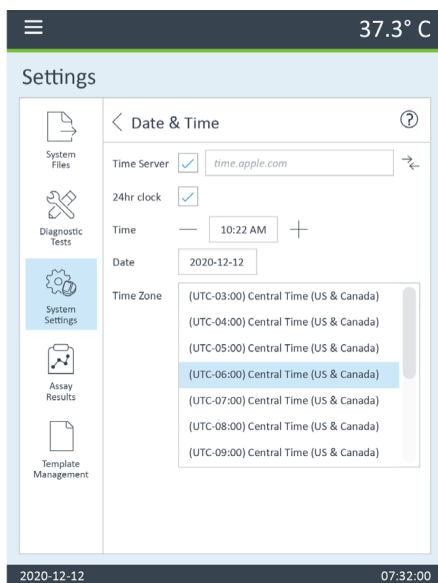


Figure 36. Select the local time zone for your XFp Analyzer.

### Setting the time

- 1 From the Home screen click **Settings > System Settings > Data & Time (+ Time Zone)**.
- 2 Select the **Date & Time** settings tab. (Figure 37 on page 61)
- 3 Manually enter the appropriate date and time.

## 5 XFp Analyzer Network Setup

### Email configuration

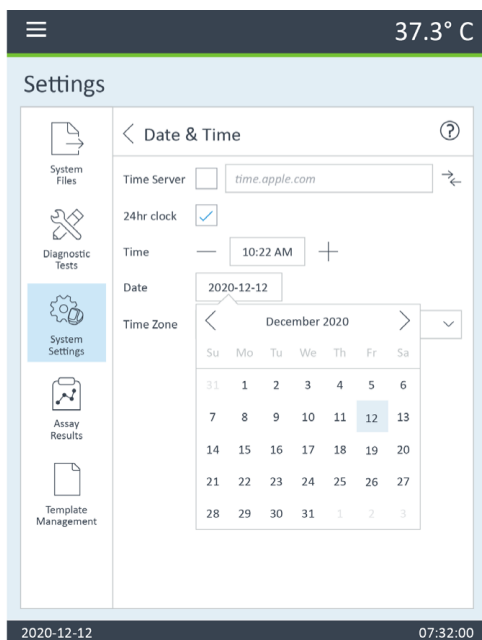


Figure 37. Configure date and time for the XFp Analyzer.

### Time server (optional)

For XFp Analyzers with an active network connection, the XFp Analyzer can sync to a network's time server, if provided. (Figure 38)

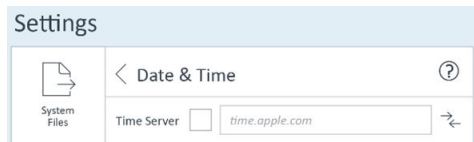


Figure 38. Configuring a time server on the XFp Analyzer.

**Optional:** A national time server (example: time.nist.gov) can be used to automatically set the time and account for daylight savings adjustments, or any time server provided by your local IT department. Once entered, click **Sync**, then click **Apply** for the changes to take place. Use the top left menu button to exit the settings options and return to Home screen.

# Network Checklist

Only complete the section that pertains to the desired network connection type.

## **Wired network settings**

- IP address
- Subnet mask
- DNS server
- Computer name

## **Wireless network settings**

- Wireless network name
- Wireless network password

## **Email settings**

- Email address
- Password
- SMTP address
- Email port
- SSL required

## **Shared directory settings**

- Shared directory address
- Shared directory domain
- Shared directory user name
- Shared directory password
- Time server address

# 6

# Support

Troubleshooting Guide 63

Technical Support and Ordering Information 64

Additional Resources 67

This chapter provides support and troubleshooting information for the Agilent Seahorse XFp Analyzer.

## Troubleshooting Guide

| Problem   | Possible Cause   | Solution  |
|---|--|---|
| Instrument gives a “load error message” and it is not clear whether a cartridge has already been loaded     |  | <ul style="list-style-type: none"><li>• Click <b>Settings</b> from Home Screen.</li><li>• Click <b>Diagnostic tests</b>.</li><li>• Click <b>Maintenance</b>.</li><li>• Click <b>Cartridge out</b>.</li><li>• The XFp Analyzer will eject a cartridge if one is left inside the instrument</li></ul> |
| Intermittent barcode read errors  | USB Wi-Fi adapter plugged into back USB port not actively being used                         | <ul style="list-style-type: none"><li>• Unplug the USB Wi-Fi adapter from the XFp Analyzer and store in a safe location</li></ul>   |
| Low rates   | Not enough cells   | <ul style="list-style-type: none"><li>• Consult the <b>Cell Reference Database</b> to identify a working range or perform a dilution series of your cell type. The wells of the XFp miniplate are the same dimensions as those of the XFe96 standard plate.</li></ul>                               |
| “An error has occurred” window pops up and displays the message: “A Task’s Exception(s) were not observed.” | Instrument was not shut down properly. See <b>“Turning the XFp Analyzer Off”</b> on page 12. | <ul style="list-style-type: none"><li>• Contact technical support. See <b>“Technical Support and Ordering Information”</b> on page 64.</li></ul>  |

# Technical Support and Ordering Information

## Worldwide technical support

For questions about XF technology, the XFp Analyzer, XF experimental design, data analysis, troubleshooting, and other information, contact Agilent Cell Analysis Technical Support:

Email: [cellanalysis.support@agilent.com](mailto:cellanalysis.support@agilent.com)

Phone:

|                           |                 |
|---------------------------|-----------------|
| Global/United States      | +1 719 528 7500 |
| United States (toll free) | +1 800 227 9770 |
| United Kingdom            | 0800 096 7632   |
| Germany                   | 0800 180 6678   |
| Europe                    | +45 3136 98 78  |
| China (toll free)         | 800 820 3278    |

## Ordering

Link to online store: <https://www.chem.agilent.com/store/>

US direct ordering:

- Email: [css\\_afo\\_fax@agilent.com](mailto:css_afo_fax@agilent.com)
- Phone: 1.800.227.9770 option #1 #1
- Fax purchase orders to: 302.633.8901

Europe:

Contact your local Customer Care Center  
<https://www.agilent.com/en-us/contact-us/page>

## Online help and support

To access the Help page, go to the Home screen then click **Help**.

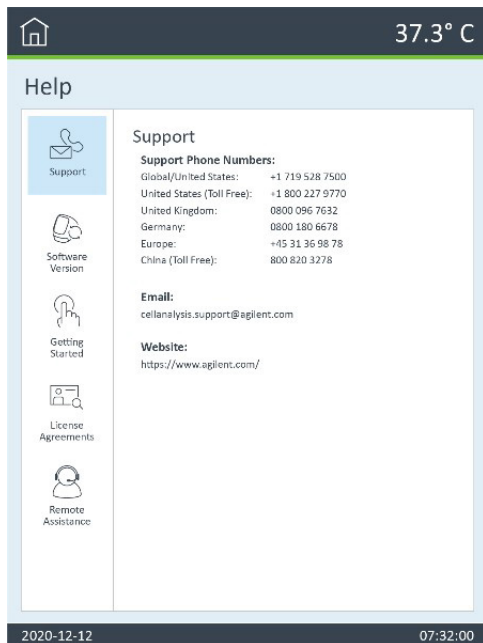


Figure 39. Help window.

### Support

Worldwide Agilent Support contact information.

## Version of software/hardware

Currently installed software and firmware versions on XFp Analyzer are displayed on this screen. If the XFp Analyzer is networked, software updates (Product Update Available) or firmware updates (Firmware Update Available) will be selectable. See **Figure 40**.

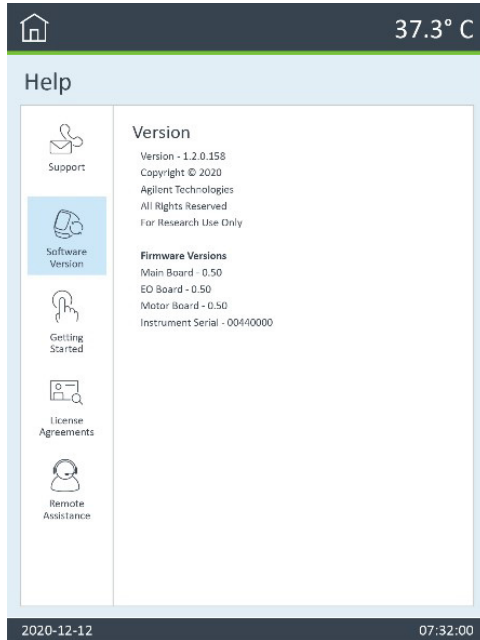


Figure 40. Software version.

## “Help” text info

Most screens on the XFp Analyzer have a small **Help** button that will display help text related to the current screen. See **Figure 41**.



Figure 41. Help button.

## Additional Resources

---

|   |   |
|---|---|
| XFp data analysis software registration | <a href="https://www.agilent.com/en/products/cell-analysis/cell-analysis-software/data-analysis/seahorse-analytics">https://www.agilent.com/en/products/cell-analysis/cell-analysis-software/data-analysis/seahorse-analytics</a> |
| XFp Consumables webpage                 | <a href="https://www.agilent.com/en/products/cell-analysis/seahorse-xfp-consumables">https://www.agilent.com/en/products/cell-analysis/seahorse-xfp-consumables</a>   |
| Cell Reference Database                 | <a href="https://www.agilent.com/cell-reference-database/">https://www.agilent.com/cell-reference-database/</a>   |
| Links to other useful information       | <a href="https://www.agilent.com/en/promotions/cell-analysis-technology">https://www.agilent.com/en/promotions/cell-analysis-technology</a>   |

---

**6 Support**  
Additional Resources



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

For Research Use Only. Not for use in diagnostic procedures.

DE.4990046296

© Agilent Technologies, Inc. 2020

Second Edition, November 2020



S7802-90001\_REV B

