

Smart-Vue Pro LoRaWAN Enabled Receiver

User Guide

331676H02 • Revision A • 05/14/2020

IMPORTANT Read this user guide. Failure to follow the instructions in this user guide can result in damage to the unit, injury to operating personnel, and poor equipment performance.

CAUTION All internal adjustments and maintenance must be performed by qualified service personnel.

Material in this manual is for informational purposes only. The contents and the product it describes are subject to change without notice. Thermo Fisher Scientific makes no representations or warranties with respect to this manual. In no event shall Thermo be held liable for any damages, direct or incidental, arising from or related to the use of this manual.

© 2020 Thermo Scientific. All rights reserved. Thermo Scientific, the Thermo Scientific logo, Smart-View Pro Duo and Smart-View Pro Quattro are the exclusive property of Thermo Scientific. All other brands mentioned are the property of their respective owners. This is a non-contractual document. Features and specifications mentioned here are subject to change.

Contents

Regulatory Information	1
Legal Information	1
FCC Certification	1
FCC Interference Notice	1
Industry Canada Class B Notice	1
Safety Notices	2
Lithium Battery	2
User Responsibility	2
Power Supply Caution	2
Device Maintenance	2
Introduction	3
Product Overview	3
Main Features	3
Architectures and Technologies	4
Setting Up your Receiver	5
Getting Started	5
LoRaWAN Configuration	7
Setting Up an Ethernet Connection	8
Setting Up a Wi-Fi Connection to your Network	10
Configuring “Failover” Order for Network Access	12
Troubleshooting	14
Firmware Upgrade from Version 1.4.17 to 5.1.2	18
Set Up Procedure	19
Smart-View Pro LoRaWAN Set Up Procedure with Smart-View Pro Duo/Quatro and Web app	19

Appendix - Product Specifications	28
Smart-Vue Pro LoRaWAN Enabled	
Receiver Features	28
Related Products	28
Installation Details	29
WEEE Compliance	30
Contact Information	31

Regulatory Information

Legal Information

47 CFR Part 15 Regulation Class B Devices

This equipment has been tested and found to comply with the limits for a Class B digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation. This equipment generates, uses and can radiate radio frequency energy if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

- Reorient or relocate the receiving antenna.
- Increase the separation between the equipment and receiver.
- Connect the equipment into an outlet on a circuit different from that to which the receiver is connected.
- Consult the dealer or an experienced radio/TV technician for help.



CAUTION: Changes or modifications to this unit not expressly approved by the party responsible for compliance could void the user's authority to operate the equipment.

FCC Certification

The antenna(s) used for this transmitter must be installed to provide a separation distance of at least 20 cm from all users and must not be co-located or operated in conjunction with any antenna or transmitter not described under this FCC id, except in accordance with FCC multi-transmitter product procedures.

FCC Interference Notice

This device complies with part 15 of the FCC Rules. Operation is subject to the following two conditions:
(1) This device may not cause harmful interference
(2) This device must accept any interference received, including interference that may cause undesired operation.

Industry Canada Class B Notice

This Class B digital apparatus meets all requirements of the Canadian Interference-Causing Equipment Regulations.

This device complies with Industry Canada license-exempt RSS standard(s). The operation is permitted for the following two conditions:

1. The device may not cause interference.
2. This device must accept any interference, including interference that may cause undesired operation.

Safety Notices

Lithium Battery

- A lithium battery (3 V, coin cell, CR1632) located within the product provides backup power for the timekeeping. This battery has an estimated life expectancy of ten years.
- When this battery starts to weaken, the date and time may be incorrect.
- Battery is not user replaceable. If the battery fails, the device must be sent back to **MultiTech** Systems for battery replacement.



CAUTION: Risk of explosion, if this battery is replaced by an incorrect type. Dispose off batteries according to instructions.

User Responsibility

Respect all local regulations for operating your wireless device. Use the security features to block unauthorized use and theft.

Power Supply Caution



CAUTION: Do not replace the power supply with one designed for another product; it can damage the modem and void your warranty. Adapter is installed near the equipment and is easily accessible.

Device Maintenance

When maintaining the device:



CAUTION: Do not attempt to disassemble the device. There are no user serviceable parts inside.



CAUTION: Do not misuse the device. Follow instructions on proper operation and only use as intended. Misuse could make the device inoperable, damage the device and/or other equipment or harm users.



CAUTION: Do not apply excessive pressure or place unnecessary weight on the device. This could result in damage to the device or harm to users.



CAUTION: Do not use this device in explosive or hazardous environments unless the model is specifically approved for such use. The device may cause sparks. Sparks in explosive areas could cause explosion or fire and may result in property damage, severe injury and/or death.



CAUTION: Do not expose the device to any extreme environment where the temperature or humidity is high. Such exposure could result in damage to the device or fire.



CAUTION: Do not expose the device to water, rain or spilled beverages. It is not waterproof. Exposure to liquids could result in damage to the device.



CAUTION: Do not place the device alongside computer discs, credit or travel cards or other magnetic media. The information contained on discs or cards may be affected by the device.



CAUTION: Using accessories, such as antennas, that **MultiTech** has not authorized or that are not compliant with **MultiTech**'s accessory specifications may invalidate the warranty.

Introduction

This document describes the set up and use of the Smart-Vue Pro LoRaWAN™ enabled receiver, a key component for collecting data in your monitoring solution.

Product Overview

It is designed for the use with Thermo Scientific wireless modules featuring Smart-Vue Pro LoRaWAN connectivity.

The Smart-Vue Pro LoRaWAN receiver is a wireless receiver that leverages new-generation Smart-Vue Pro LoRaWAN Internet of Things technology to offer exceptionally long-range wireless communication.

This receiver can be connected to your local network via a wired Ethernet or a Wi-Fi connection.



Figure 1. Smart-Vue Pro LoRaWAN Enabled Receiver

Main Features

Technical Highlights

- Up to 10 miles (16 km) line-of-sight wireless range.
- 2-ways wireless communications.
- Available Smart-Vue Pro LoRaWAN protocol frequencies:
 - 868 MHz
 - 915 MHz
- Wired Ethernet or Wi-Fi network connection.

In the Box

- Smart-Vue Pro LoRaWAN enabled receiver
- AC adapter (100-240 V)
- Antenna
- Ethernet cable

Architectures and Technologies

Installed locally at your site, the Smart-Vue Pro LoRaWAN enabled receiver collects data from compatible Thermo Scientific modules within wireless range.

The receiver is connected permanently via internet to the server where data logger data is pushed by the receiver and accessed using the Smart-Vue Pro web application. The following diagram shows the overall solution organization with the receiver collecting data from data loggers and forwarding it to the cloud platform.

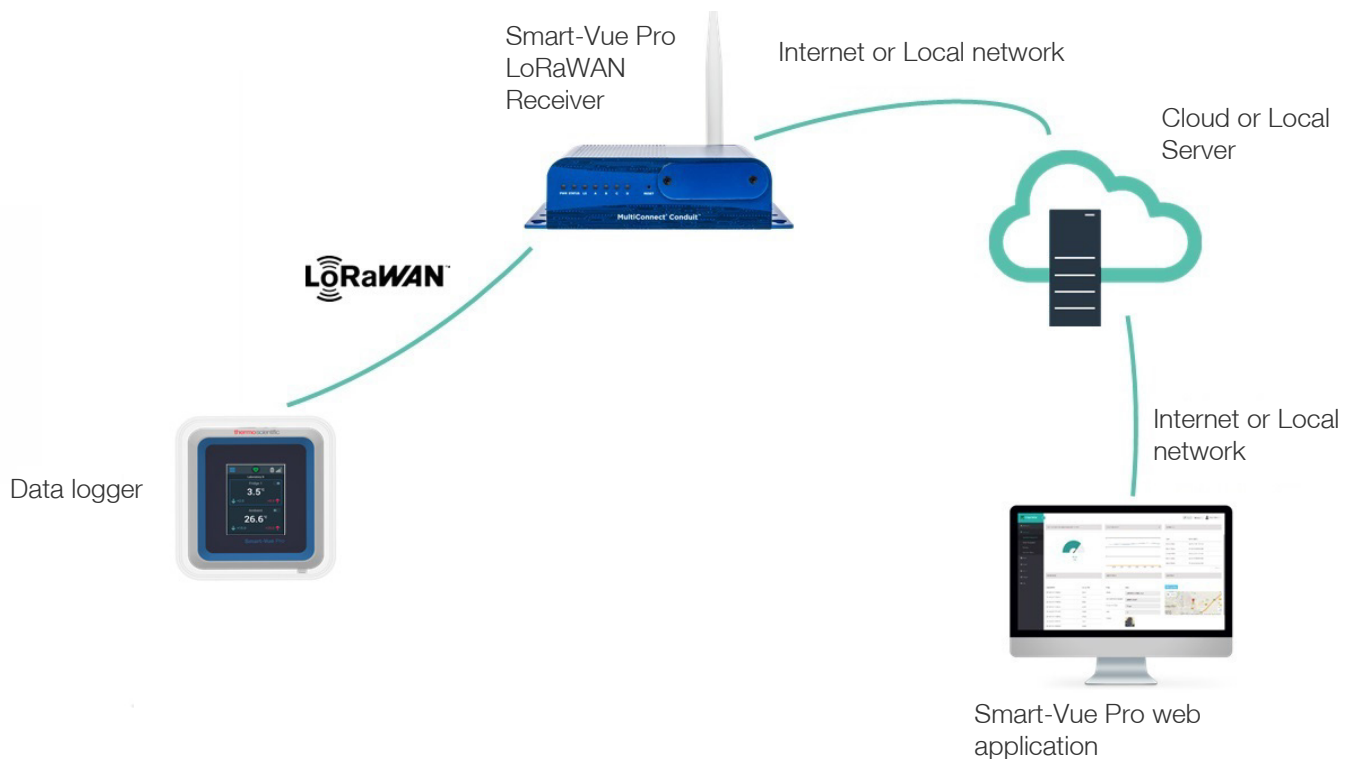


Figure 2. Receiver connecting Smart-Vue Pro Quatro/Duo data logger to the Cloud

Setting Up your Receiver

The Smart-Vue Pro LoRaWAN enabled receiver is configured using an integrated web interface. You must use that interface to determine how your receiver connects to the internet, that is, via an Ethernet or Wi-Fi. This chapter describes how to connect to your receiver and enter the appropriate setting for your solution.

Getting Started

Plug in the Smart-Vue Pro LoRaWAN Antenna

1. Plug the white Smart-Vue Pro LoRaWAN radio antenna into the “RF” connector (the location may vary according to your unit’s options).
2. Hand-tighten the antenna by rotating the ring clockwise.



Figure 3. Attach Smart-Vue Pro LoRaWAN Antenna to connector labeled “RF”

Plug in the Power Cable

1. Plug the power cable into the stainless-steel connector behind the receiver. The cable fits correctly onto one of the connectors.
2. Tighten the cable by rotating the steel ring clockwise to attach it firmly to the unit.



Figure 4. Plug in the power cable to the stainless steel connector

3. Plug the AC adapter into a power socket to boot the receiver. The startup process may take 2-3 minutes to complete. When the unit is ready for use, the left-hand LED remains lit as shown here (the status LED continues to blink):

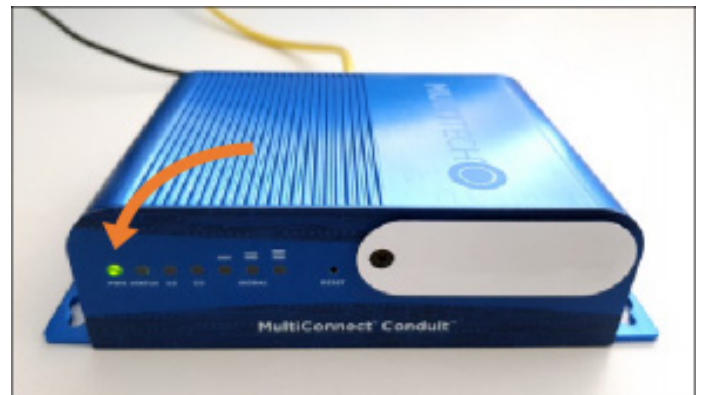


Figure 5. Power indicator remains on when receiver is ready to use

Connect to the Configuration Interface

1. Plug the ethernet network cable into the ethernet port behind the receiver.



Figure 6. Receiver with ethernet cable

2. Plug the other end of the cable into either your:

- a. Local Area Network

If your Local Area Network uses the following IP addresses: 192.168.2.x, you may skip to step 3. Otherwise or if you are unsure, see step (b).

- b. Computer

The receiver does not have a DHCP server which means you must manually set an IP address on your computer to communicate with the receiver.

In your computer's network settings, assign the following IP information temporarily (ensure you change it back when done configuring your receiver):

IP: 192.168.2.199

Mask: 255.255.255.0

Do not change the Default Gateway or DNS

3. Use your web browser to connect to the receiver's default IP address: <http://192.168.2.1>



IMPORTANT NOTE: We recommend to use Google Chrome as your web browser. Internet Explorer is not supported.

Upon first boot, the gateway is in "commission mode" and you will be prompted to create a new user-name and password. There is no default user-name or password.

4. Assign user-name in commissioning mode and press **OK**.



Figure 7. Assign user-name in commissioning mode

5. Assign a password for that user-name and press **OK**:



Figure 8. Assign a password for the user-name

6. You may then login using the information you just entered.



Figure 9. Enter Login Credentials

Note: User-name and Password

While performing the initial configuration of the device for the network, an administrative user-name and password is set.

Ensure that this information complies with your organization's user account provisioning policies and follows best security practices.

Rather than setting a password for the administrative account, consider setting a pass phrase. A combination of 4 or more unrelated words with spaces is statistically more secure than assigning a password, regardless of complexity.

Store this user-name and password information in the designated repository defined by your organization's policies. A digital password vault that has the ability for multiple users is one option, while storing it in a fireproof safe is another. It is recommended that the user-name and password should not be stored in a location easily accessible by those not requiring access because of the access privileges allowed to the device with these credentials.

First-Time Setup

Setting Up Your Device Using Setup Wizard (After Choosing Reset and Factory Default Settings)

When the power is turned on the device, set the device to factory default settings to see the first-time setup. This wizard helps you configure the main features of your device.

Here are the steps for first-time setup:

1. When the power is turned on for the first time or after you set factory default settings, the device goes into commissioning mode. The system requires you to set up an admin user. Enter your user-name and click **OK**.
2. Enter a password for the admin user and click **OK**. This password must be of sufficient length and strength (with a mix of character classes such as letters, numbers and symbols). Enter the password again to confirm. Click **OK**.
3. On the first page, you see the below screen. Click **Next**.



Figure 10. First Time setup wizard

4. Configure **Call Home**.

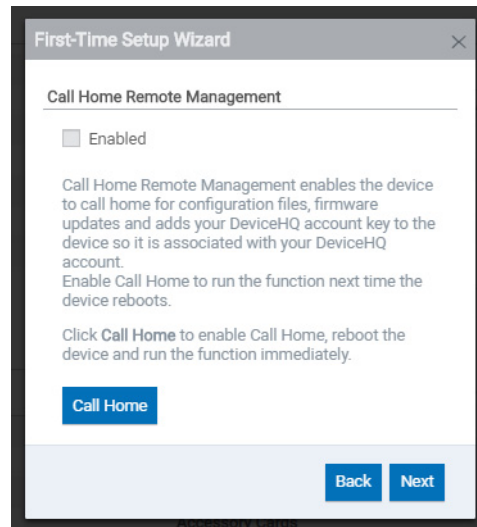


Figure 11. Call home configuration

- a. This feature is not supported for the Smart-Vue Pro solution so click **Next** to go to the next screen without changing the settings.
5. Set the date, time and time zone.

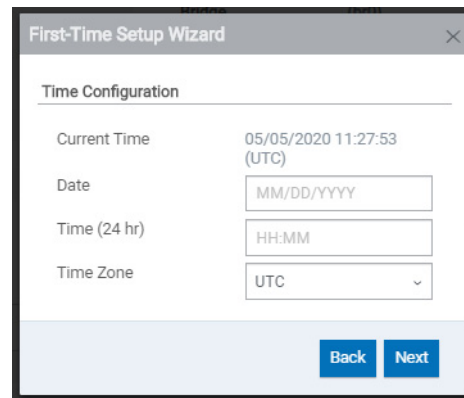


Figure 12. Time configuration

- a. Enter the **Date**.
- b. Enter the **Time**.
- c. Select the **Time Zone** in which the device operates.
- d. Click **Next**.
- e. By default the date and time are in UTC format. If you do not wish to change the current settings of date and time simply click **Next** to go the next screen.
6. Configure LAN network interfaces Eth0 and Br0.

First-Time Setup Wizard

Network Interface Configuration - eth0

Bridge: --

IPv4 Address: 192.168.2.1

Mask: 255.255.255.0

Network Interface Configuration - br0

IPv4 Address: 192.168.99.1

Mask: 255.255.255.0

Back **Next**

Figure 13. Network interface configuration

- a. Leave the settings to default and click **Next** to go to the next screen. The detailed procedure to configure the Ethernet is mentioned in “**Setting Up an Ethernet Connection**”.
7. Set up **Remote Management**.

First-Time Setup Wizard

Remote Management

☐ Enabled ☒ SSL Enabled

Server Name: ds.devicehq.com

Server Port: 5798

App Store URL: https://www.devicehq.com

Account Key:

Back **Next**

- a. This feature is not supported for the Smart-Vue Pro solution so simply click **Next** to go to the next screen without changing the settings.
8. Configure **HTTP/HTTPS Access**.

First-Time Setup Wizard

Access Configuration

HTTP Redirect to HTTPS

☒ Enabled

Port: 80

☒ Via LAN

☒ Via WAN

HTTPS

☒ Via WAN

Port: 443

Back **Next**

Figure 14. Access configuration

- a. This feature is not supported for the Smart-Vue Pro solution so click **Next** to go to the next screen without changing the settings.
9. Set up **Bootloader Protection** by setting a u-boot password.

First-Time Setup Wizard

Bootloader Protection

Authentication Status: Disabled

Set Password

Password:

Enable

Back **Finish**

Figure 15. Bootloader Protection

- a. This feature is not supported for the Smart-Vue Pro solution so click **Next** to go to the next screen without changing the settings.
10. Click **Finish**.
11. To save your changes, click **Save** and **Restart**.

LoRaWAN Configuration



IMPORTANT NOTE: When you connect to the receiver for the first time, a First-Time Setup Wizard runs automatically. We recommend that you close the wizard and set up the receiver manually as described here.

Start your receiver configuration by verifying LoRaWAN networking settings.

1. Click **LoRaWAN (Network Settings)**.

The screenshot shows the 'LoRaWAN NETWORKING' configuration page. The left sidebar contains navigation links: Home, Save And Restart, LoRaWAN (selected), Network Settings, Setup, Wireless, Firewall, Tunnels, Administration, Status & Logs, Commands, Apps, and Help. The main content area is titled 'LoRaWAN NETWORKING' and includes a 'LoRa Mode' section with fields for Mode (PACKET FORWARDER), Packet Forwarder (4.0.1-r26.0), Network Server (2.3.0), and Lens Server (2.3.0). Below this are status indicators for 'Restart LoRa Services', 'RUNNING', 'DISABLED', and 'DISABLED'. The 'LoRa Card Information' section shows Frequency Band (868) and FPGA Version (31). The 'LoRa Packet Forwarder Configuration' section includes a 'Channel Plan' dropdown (EU868), an 'Intervals' section with 'Keep Alive Interval (s)' set to 10, and a 'Server' section with 'Server Address' (smartvueconnect.com) and 'Port' (1700). At the bottom are 'Submit', 'Reset To Default', and 'Test LoRa Server' buttons.

Figure 16. LoRaWAN configuration screen

2. Ensure the following default values are entered.
 - a. For Smart-Vue Pro Quatro/Duo data loggers:
 - Server Address: **smartvueconnect.com**
 - Port (up & down): **1700**
 - Keep Alive Interval: 10 seconds
3. Click **Test LoRa Server** in the lower right-hand corner to verify the connection.



CAUTION: If you enter a domain name rather than an IP address, the DHCP client in the receiver must be able to resolve the name to connect to the internet (described in the next section).



CAUTION: Test Lora Server will only work once you configure your network properly (see **Setting Up an Ethernet Connection to Setting Up a Wi-Fi Connection to your Network**).

4. When the test run has succeeded, click **Submit** to apply your changes.

5. You must restart the gateway services for the settings to be saved. Click **Save And Restart** on the left side to save and restart the gateway.
6. Once the services are restarted, LoRaWAN networking status is updated.

Setting Up an Ethernet Connection

This section describes how to configure your receiver with an Ethernet wired connection. The receiver does not require to be placed in its final location for this operation, but it does need to be accessible over your network.

1. Click **Setup → Network Interfaces → Options** (Edit) for “eth0” **(1)**.

mPower™ Edge Intelligence Conduit - Application Enablement Platform
MTC01-247A Firmware 5.1.2-OCEA-r50

admin as administrator

Home
Save And Restart
LoRaWAN ®
Setup

Network Interfaces
WAN Configuration
Global DNS
DDNS Configuration
DHCP Configuration
GPS Configuration
SMTP Configuration
SNMP Configuration
Time Configuration

Wireless
Firewall
Tunnels
Administration
Status & Logs
Commands
Apps
Help

NETWORK INTERFACES CONFIGURATION ⓘ

Reset To Default

Name	Direction	Type	IP Mode	IP Address	Bridge	Options
eth0	WAN IPv4	ETHER	Static	192.168.0.250/24		
wlan0	WAN IPv4	WIFI_AS_WAN	DHCP Client	192.168.225.45/16		
wlan1	LAN IPv4	WIFI_AP	Static		--	
br0	LAN IPv4	BRIDGE	--	--	br0	

Copyright © 1995 - 2020 by Multi-Tech Systems, Inc. - All rights reserved.

mPower™ Edge Intelligence Conduit - Application Enablement Platform
MTC01-247A Firmware 5.1.2-OCEA-r50

admin as administrator

Home
Save And Restart
LoRaWAN ®
Setup

Network Interfaces
WAN Configuration
Global DNS
DDNS Configuration
DHCP Configuration
GPS Configuration
SMTP Configuration
SNMP Configuration
Time Configuration

Wireless
Firewall
Tunnels
Administration
Status & Logs
Commands
Apps
Help

NETWORK INTERFACE CONFIGURATION - ETH0 ⓘ

Direction
WAN

IPv4 Settings

Mode
Static

IP Address
192.168.0.250

Mask
255.255.255.0

Gateway
192.168.0.1

Primary DNS Server
8.8.8.8

Secondary DNS Server

Submit Cancel

Copyright © 1995 - 2020 by Multi-Tech Systems, Inc. - All rights reserved.

Figure 17. Ethernet adapter settings for your network

Smart-Vue Pro LoRaWAN Enabled Receiver

Setting Up your Receiver | 11

2. Adjust the various settings in the **Network Interface Configuration** window to match your network:

Table 1. Network Interface Configuration

Option	Settings
Direction	WAN
Mode	Static (in which case you must set the IP Address in the field below) or DHCP Client (the receiver gets its IP address from your network's DHCP server).
IP Address, Mask, Gateway, Primary DNS Server, Secondary DNS Server	Enter the IP addresses according to your network configuration.

3. Click **Submit** when you are done.

If you connected your computer directly to the receiver with an ethernet cable remember to plug the receiver into your Local Area Network.



CAUTION: It is essential to leave the DHCP server option disabled on your receiver (**Setup → DHCP**). Enabling that option could create a conflict with your network's DHCP server. This is not to be confused with the DHCP client option described above. Its not required to configure other options in the Network Interfaces (described in **Setting Up a Wi-Fi Connection to your Network**).

Note: Ethernet Network Address Settings

If you wish to specify what devices can communicate with the system by way of Access Control Lists (ACLs), the network address should be set to a static IP. Your IT team will assist in ensuring the routes required are open for each device to connect through any firewalls.

Setting Up a Wi-Fi Connection to your Network

This section describes how to configure your receiver with a Wi-Fi wireless connection using DHCP automatic IP address assignment.



CAUTION: If you need to use your LoRaWAN router in Wi-Fi mode with a static IP address, See **Workaround for Wi-Fi Connection with Static IP Address**.

To set up Wi-Fi wireless access for your receiver:

1. Plug the provided Wi-Fi antenna into the Wi-Fi connector and rotate the ring clockwise to fasten the antenna securely:



Figure 18. Attaching antenna for Wi-Fi network

2. Click **Wireless → Wi-Fi as WAN**, as shown below:

Figure 19. Adding a Wi-Fi network

3. Click **Add Network (1)** and fill in the information as required:

Figure 20. Wi-Fi Network Settings

Table 2. Network Information

Option	Setting
Network Name	Assign a name for the network.
SSID	Enter the exact network name.
Security Mode	Choose the security and key encryption methods used by your Wi-Fi network.

4. Click **Finish → Submit (2)** to save your settings, then click **Save and Restart** to complete configuration.

Note: Wi-Fi Network Connection Settings

All configuration to connect to your local WiFi network is set internal to your organization. You may have a captive portal or pre-shared key required to obtain access to your wireless network. If this is the case, follow the internal process to receive an exception for the LoRaWAN receiver for approval and to maintain this connection.

Workaround for Wi-Fi Connection with Static IP Address

As of this writing, the Smart-Vue Pro LoRaWAN receiver does not support the use of a static IP addresses for Wi-Fi communications.

To use your Smart-Vue Pro LoRaWAN receiver in Wi-Fi mode with a static IP address, you must implement a workaround solution using an intermediary Wi-Fi repeater or router whose setup allows a fixed IP address, as shown below, to connect to your network.



CAUTION: The Wi-Fi router or repeater must also support being configured as a Wi-Fi client in order to connect to your organization's Wi-Fi access point.

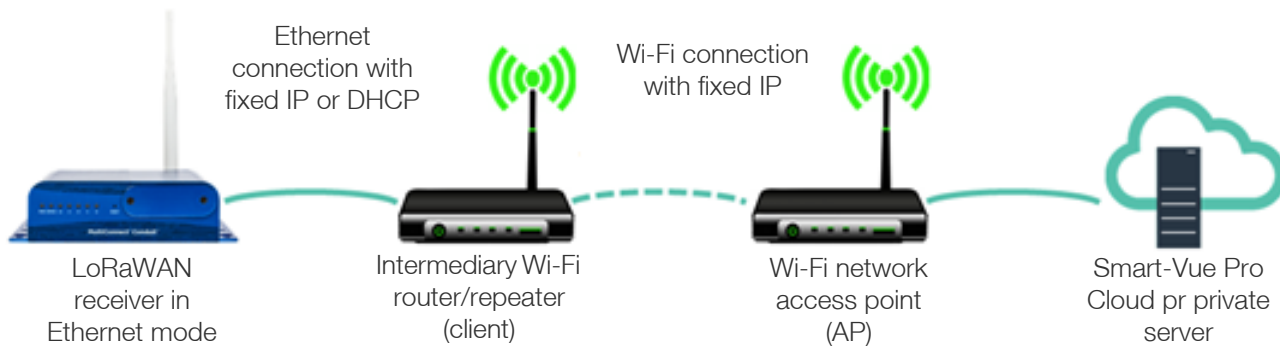


Figure 21. Using an intermediary Wi-Fi router or repeater to provide a static IP address

To implement this workaround:

1. Set up an intermediary Wi-Fi router/repeater whose configuration options allow you to set a static IP address for connecting to your network. That device must also be configurable as a Wi-Fi client (and not just as an access point (AP)).
2. Connect the router to your network.



Note: We recommend that you test the router using a laptop computer plugged into an Ethernet port (as the Smart-Vue Pro LoRaWAN receiver will be).

3. Set up the Smart-Vue Pro LoRaWAN receiver to use an Ethernet connection (see **Setting Up an Ethernet Connection**) with either a static IP address or using DHCP according to your requirements.
4. Connect the Smart-Vue Pro LoRaWAN receiver to the Wi-Fi router using an Ethernet cable.

Configuring “Failover” Order for Network Access

If you have more than one internet connectivity option installed in your receiver, the receiver can switch from one to another in case of failure. For example, if you choose to run ethernet as your primary connection and the connection fails at some point, you could backup the communication channel by also configuring the Wi-Fi data connection. To do this:

1. Click **Setup → WAN Configuration**.

The list below shows the priority order in which connectivity options are used. In the case below, Ethernet is used in priority. The receiver would switch automatically to Wi-Fi data if communication via Ethernet fails.

2. Click the up/down arrows **(1)** to change network priority order.

WAN CONFIGURATION ⓘ

General Configuration

Mode

FAILOVER

WANS

Priority	State	Name	Type	Options
1	Enabled	eth0	ETHERNET	^ v ✎
2	Enabled	wlan0	WIFI	^ v ✎

Reset To Default

Figure 22. Determine network “failover” order

Note: Network Configuration

Upon completion of the network configuration of the LoRaWan receiver and to support ongoing availability of the device and its functionality, Thermo Fisher recommends configuring the device for a “Failover” network configuration. This helps to ensure that variables in network performance (wired or wireless) are not impacting the data flows, thus supporting the availability of the device.

Troubleshooting

How do I access the Smart-Vue Pro LoRaWAN configuration page?

For the first time when you plug in your Smart-Vue Pro LoRaWAN, use your web browser to connect to the receiver's default IP address: `http://192.168.2.1`.

Once Wi-Fi or ethernet is successfully configured and saved then use the IP address assigned to the Smart-Vue Pro LoRaWAN to connect.

The assigned IP address for your Smart-Vue Pro LoRaWAN can be seen on the home page of Smart-Vue Pro LoRaWAN. Note the IP address for later use.

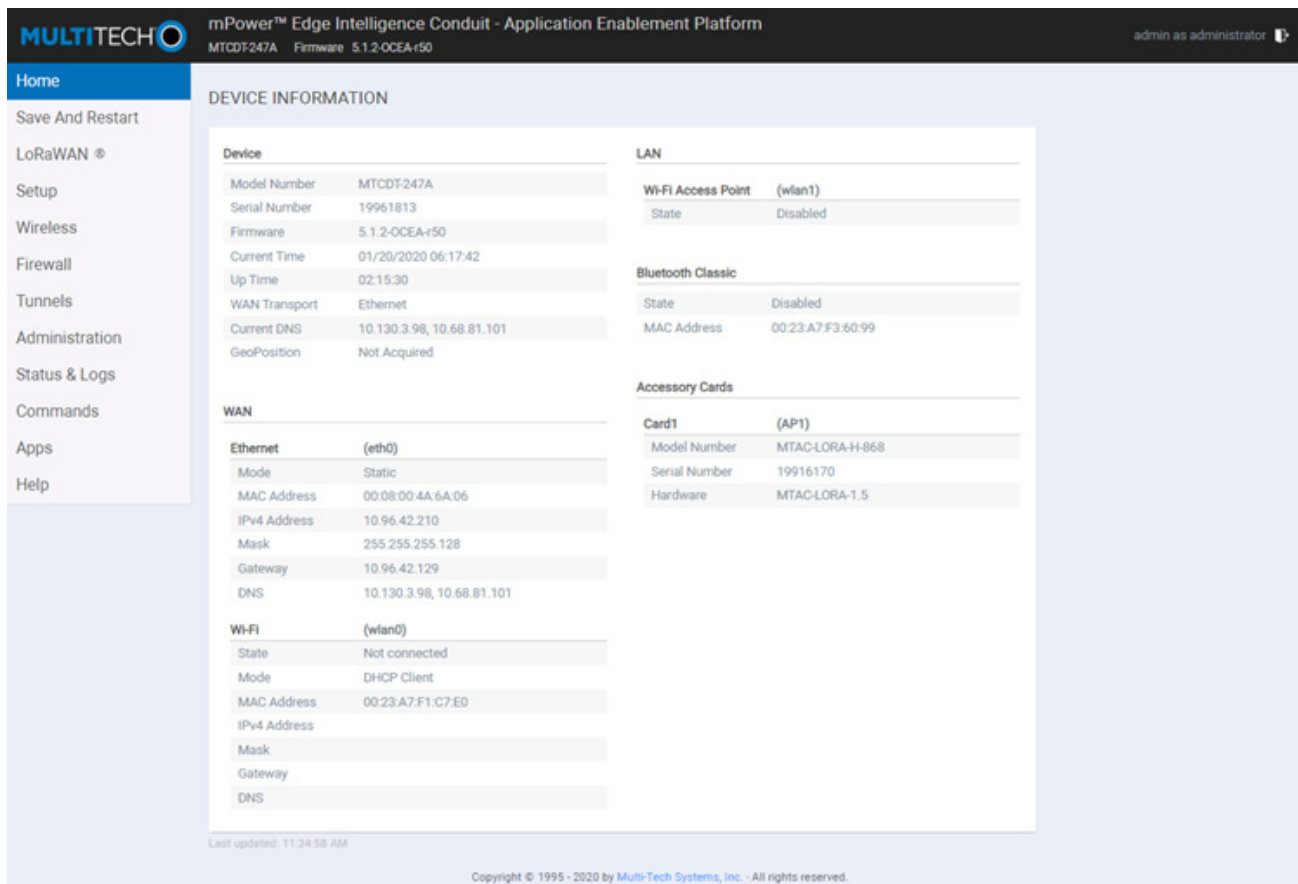


Figure 23. Smart-Vue Pro LoRaWAN configuration page

Can I test my Smart-Vue Pro LoRaWAN network connection?

In the **Setup → LoRa** window (described in **Smart-Vue Pro LoRaWAN Configuration**), is a **Test LoRa Server** button. When you Click that button, you receive an acknowledgment that the connection is up and running.

Why sensors connected to Smart-Vue Pro Quatro does not get detected in the web application?

Sensors attached on Smart-Vue Pro Quatro sometimes may not be detected due to the latency in the network which is configured in Smart-Vue Pro LoRaWAN. If the latency is greater than 200 mSec, then this issue occurs. Use the 'Test LoRa Server' option to measure the latency.

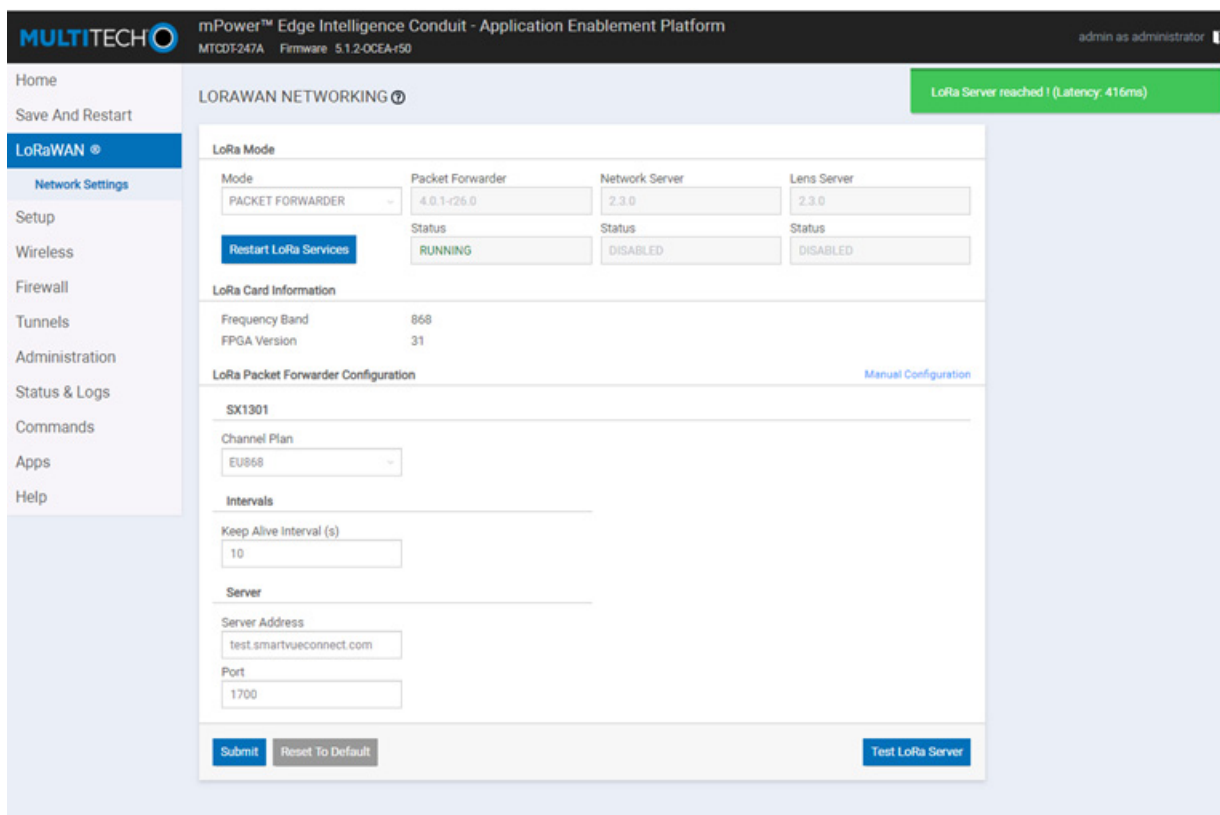


Figure 24. Smart-Vue Pro LoRaWAN network connection

How can I be confirmed that communication between the Smart-Vue Pro LoRaWAN and the webapp is established?

This can be confirmed by the latency time. The latency (time taken for the data logger to communicate with the web application via the Smart-Vue Pro LoRaWAN and receive a response) should be less than 200 ms. Latency details are visible when the 'Test LoRa server' option is clicked on the Smart-Vue Pro LoRaWAN web interface page. (See **Figure 24**).

I would like to see how well the LoRaWAN connection works at my site. Is there an easy way to test this?

You may test the wireless connection between your receiver and a Smart-Vue Pro LoRaWAN data logger as described in the data logger user guide. For more information, or to conduct a more complete site survey, contact your Thermo Scientific representative.

How can I check the coverage status of the Smart-Vue Pro LoRaWAN signal between Smart-Vue Pro Quatro and the Smart-Vue Pro LoRaWAN?

On Smart-Vue Pro Quatro, the 'Range Test' feature helps you know if there is Smart-Vue Pro LoRaWAN signal coverage between Smart-Vue Pro Quatro and the Smart-Vue Pro LoRaWAN.

A good coverage (100%) will be shown on the device.

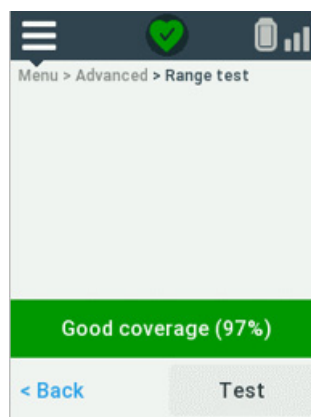


Figure 25. Smart-Vue Pro LoRaWAN signal

Why is the telemetry data from the sensors (connected to Smart-Vue Pro Quatro) not showing on the web app?

Verify the below:

- Check if the option “**Smart-Vue Pro LoRaWAN ON/OFF**” on Smart-Vue Pro Quatro is enabled.
- Ensure the sensor is disconnected from earlier data logging sessions completed.
- Smart-Vue Pro LoRaWAN is switched **ON** and configured network latency is less than 200 mSec.

- Switch off the Smart-Vue Pro LoRaWAN Quatro by disconnecting the AC Adapter (If plugged in) and remove the batteries and then replace batteries and connect back AC Adapter (if required). Press show sensor on device.
- Disconnect sensor and connect the sensor back. Press “Refresh sensors” and ensure the sensor value is seen in “Show Sensor” screen of Smart-Vue Pro LoRaWAN Quatro.

What is the default time to wait for complete functionality to start on Smart-Vue Pro LoRaWAN after restart / power on restart?

After a power on restart/ restart waits for at least 2 minutes for the complete setup to be up.

Why is data logging not starting/stopping on web app?

- This would be due to bad network. If the latency is more than 200 msec, then the selected command will not reach the Smart-Vue Pro LoRaWAN and Smart-Vue Pro LoRaWAN Quatro.
- Ensure that the network used for communication from Smart-Vue Pro LoRaWAN has good internet speed.
- Stop data logging takes certain time to stop based on the data which need to be synced from Smart-Vue Pro LoRaWAN Quatro to cloud. Audit trails in web app shall show the time of the event happened.

What to do when it shows “Unable to reach Smart-Vue Pro LoRaWAN server”?

Ensure port 1700 is open and inform IT team that it should communicate via UDP.

I'd like to test LoRaWAN wireless coverage at my site. The receiver is in place, but I don't want to use it on the internet and just use LoRaWAN to check coverage. Is it possible?

There is an auto-acknowledgment wireless testing feature in SVP Quatro data logger firmware. If you want to test on-site coverage without an internet connection (that is, just the LoRaWAN part), you may configure the SVP Quatro data logger as follows:

Select **Menu → Advanced → LoRaWAN → Network → Custom**.

Select the region and enter this 8-digit number via the keypad: “00000000”. On gateway side, click **LoRaWAN (Network Settings)**, change the “Server Address” to 127.0.0.1. Click to use the **Range test** feature on the SVP Quatro data logger to test LoRaWAN connectivity with your receiver.

How to reset the login password if password is blocked or forgotten?

The LoRaWAN receiver need to be reset, to set a new password.

How do you reset the LoRaWAN receiver?

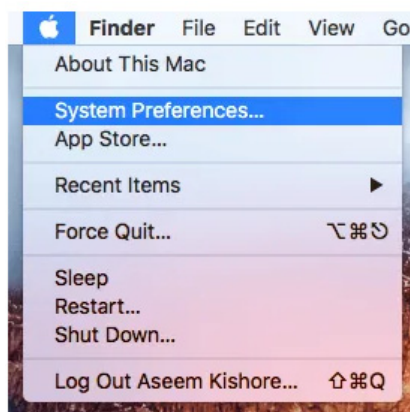
Switch on the power on the Lora receiver and then press and hold the reset button on front panel for more than 30secs and then release the button. The status LED on front panel will be steady for more than 2mins and then blink. Once the device is ready, as explained in section “**Connect to the Configuration Interface**” configure the device.

How to assign a Static IP address in Windows 7, 8, 10, XP or Vista?

In the “Network Connections” window, right-click the adapter for which you want to set a static IP address, and then select the “Properties” command.

Step by step illustration:

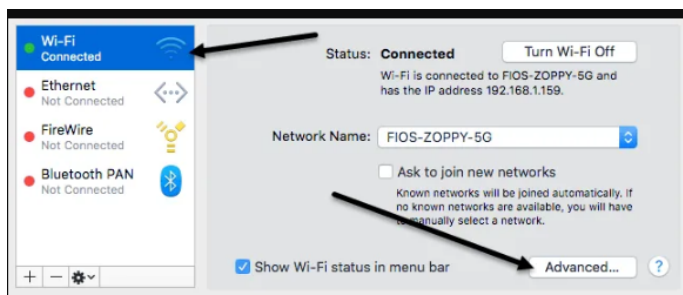
1. Go to **System Preferences**.



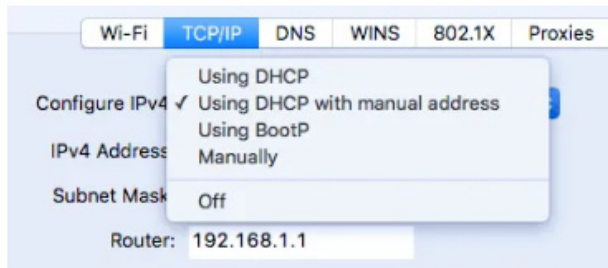
2. Click the **Network icon**.



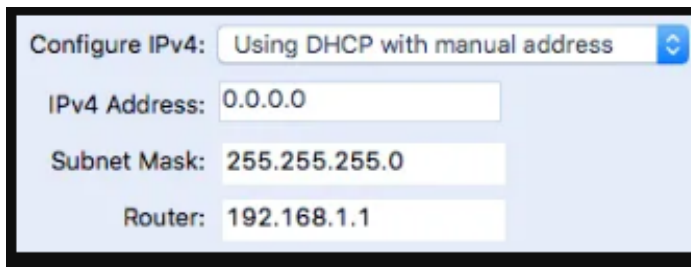
3. You will see a list of network connections on the left. If the connection is green, that means it is active. Click the connection and then click the **Advanced** button at the bottom right.



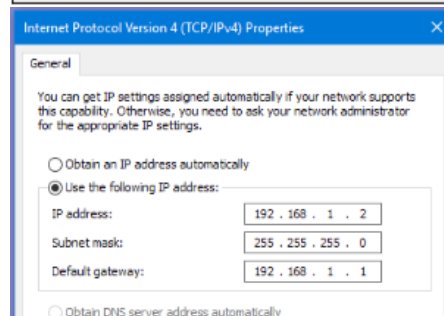
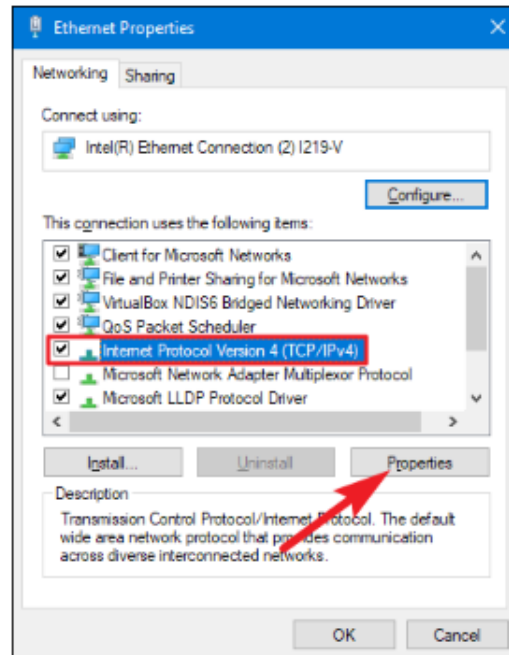
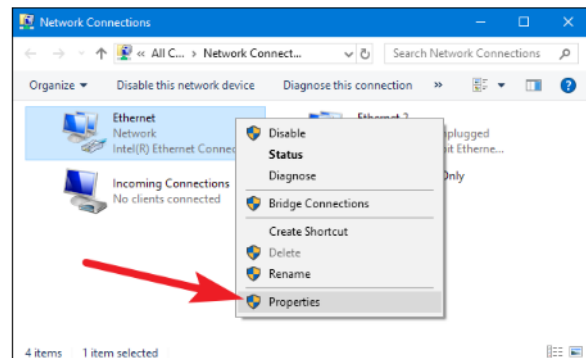
4. This will bring up all the advanced settings for the network connection. Click the TCP/IP tab and you'll see a drop-down next to Configure IPv4.



5. This will bring up all the advanced settings for the network connection. Click the TCP/IP tab and you'll see a drop down next to Configure IPv4.



For Windows OS:



Firmware Upgrade from Version 1.4.17 to 5.1.2

This section provides specific information when upgrading your gateway's firmware from version 1.4.17 to version 5.1.2.

To upgrade the firmware in your gateway, typically follow notification from thermo fisher scientific or contact with our technical support team, follow these steps:

1. Connect to your gateway with your login name and password.
2. Select **Administration** → **Firmware upgrade**.
3. Select **Choose Firmware Upgrade** File and browse your computer to locate the firmware file provided to you by Smart-Vue Pro.



CAUTION: Use official firmware files provided by thermo scientific.

5. When upgrading from 1.x.x firmware, this process resets your previous LoRaWAN configuration. You must then click **LoRaWAN** in the main menu, then click the **Normal Configuration** link (1) on the right-hand side of the screen to open **Manual Configuration**.
6. You may then proceed with LoRaWAN configuration as described earlier. This specific behavior is caused by the migration from firmware 1.x.x to 5.1.2. After this upgrade, you should not have to perform this procedure again in the future.

4. Select Start upgrade and continue the process as directed on the screen.

The screenshot displays the MULTITECH mPower Edge Intelligence Conduit - Application Enablement Platform interface. The top header shows the MULTITECH logo and the text 'mPower™ Edge Intelligence Conduit - Application Enablement Platform'. Below the header, the page is titled 'LoRaWAN NETWORKING'. The interface includes a sidebar with navigation links: Home, Save And Restart, LoRaWAN (selected), Network Settings, Setup, Wireless, Firewall, Tunnels, Administration, Status & Logs, Commands, Apps, and Help. The main content area shows the 'LoRa Mode' section with fields for Mode (PACKET FORWARDER), Packet Forwarder (4.0.1-r26.0), Network Server (2.3.0), and Lens Server (2.3.0). Below these are status indicators: Status (RUNNING), Status (DISABLED), and Status (DISABLED). A 'Restart LoRa Services' button is also present. The 'LoRa Card Information' section shows Frequency Band (868) and FPGA Version (31). The 'LoRa Packet Forwarder Configuration' section includes a 'Config (examples)' text area with a code snippet. A blue link labeled 'Normal Configuration' is highlighted with an arrow and the number 1. At the bottom, there are 'Submit', 'Reset To Default', and 'Test LoRa Server' buttons.

Figure 26. Normal Configuration Link

Set Up Procedure

Smart-Vue Pro LoRaWAN Set Up Procedure with Smart-Vue Pro Duo/Quatro and Web app

1. Ensure Smart-Vue Pro Quatro is connected with sensors and sensor value is shown in show sensor screen.

Example: Smart sensor value displayed in show sensor screen.

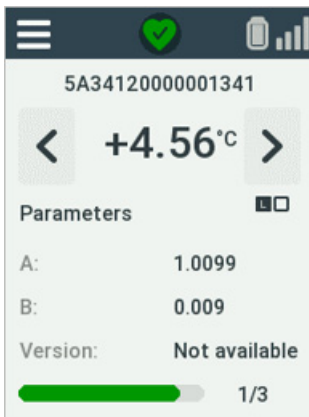


Figure 27. Smart Sensor value display

2. Ensure LoRaWAN is ON in Smart-Vue Pro Quatro.

In Smart-Vue Pro Quatro device,
Click **Menu** → **Advanced** → give the code which is set in webapp and added by user during configuring the user setting i.e., pin code set in webapp by the user.

Then in Advanced, once a valid code is entered, select the **LoRaWAN** → select **ON** option from **ON/OFF** to enable Smart-Vue Pro LoRaWAN connectivity. After selecting the LoRaWAN as ON, go back to LoRaWAN → Network → ThermoScientific → select your region and Save.

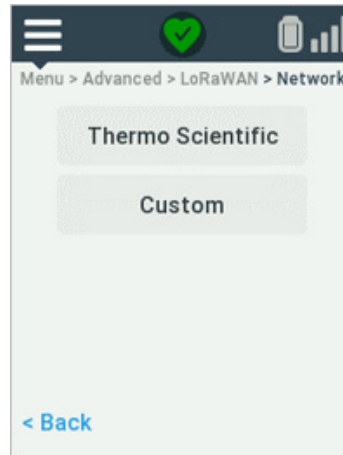


Figure 28. In Smart-Vue Pro Quatro device

3. Press “Test” in Smart-Vue Pro Quatro device to ensure the Smart-Vue Pro LoRaWAN coverage is good.

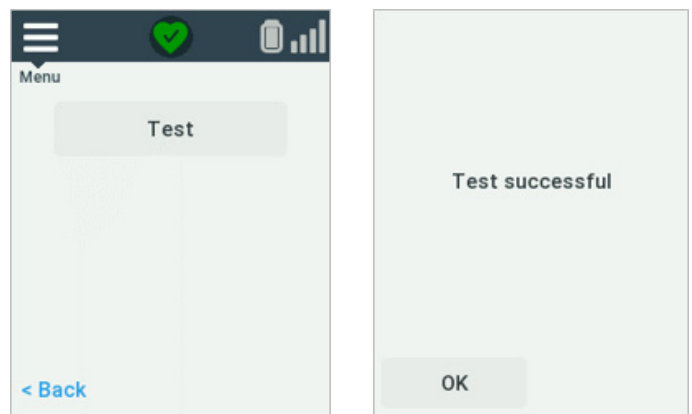


Figure 29. Smart-Vue Pro LoRaWAN coverage

4. Open Smart-Vue Pro LoRaWAN web browser page, Configure the Network in Smart-Vue Pro LoRaWAN settings to communicate to cloud.
Refer **Setting Up an Ethernet Connection** and **Setting Up a Wi-Fi Connection to your Network** which explains the setting up of different networks like Wi-Fi and Ethernet.
 - a. Ethernet Configuration:

Click **Setup** → **Network Interfaces** → **Options** (Edit) for “eth0” (1).

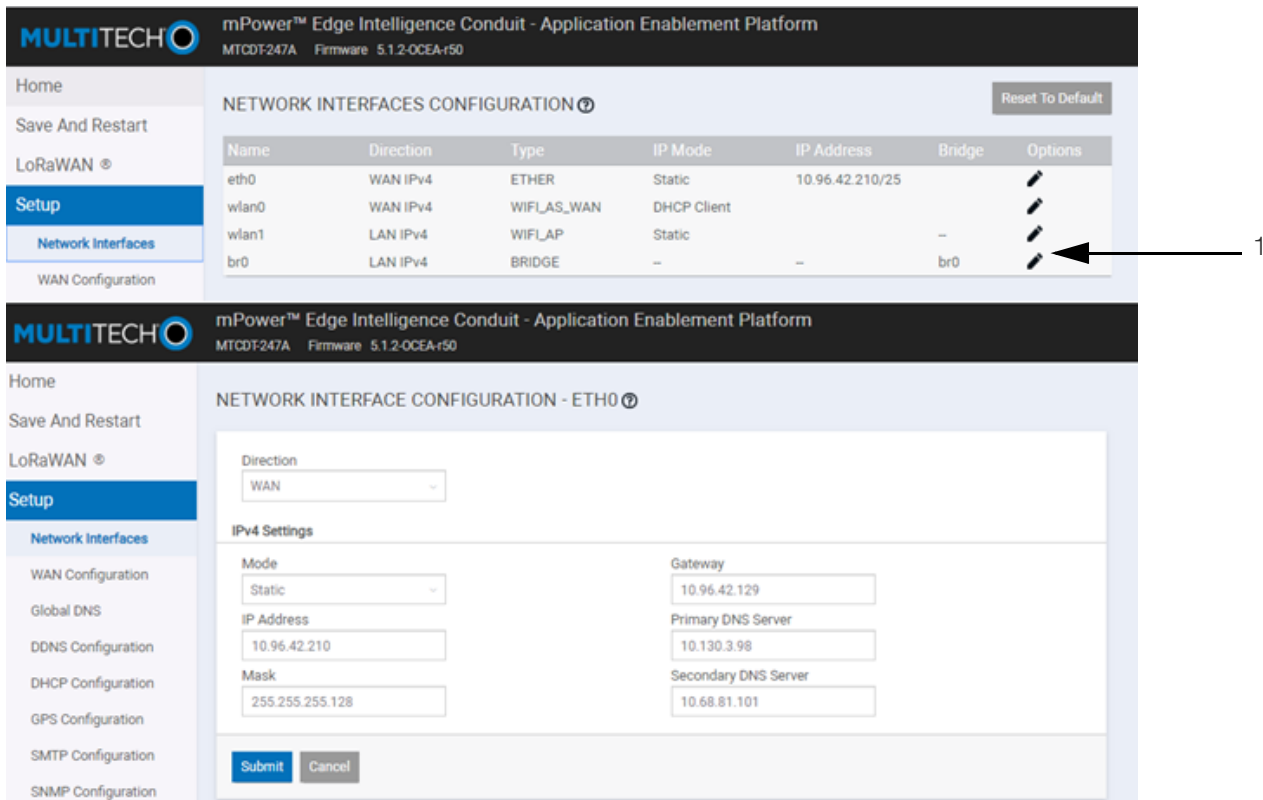


Figure 30. Ethernet Configuration

- b. Wi-Fi onfiguration:
 - To save the settings, save and restart to be selected.
 - **Wireless → Wi-Fi as WAN →** add the network and select the enabled option and submit the settings to save.

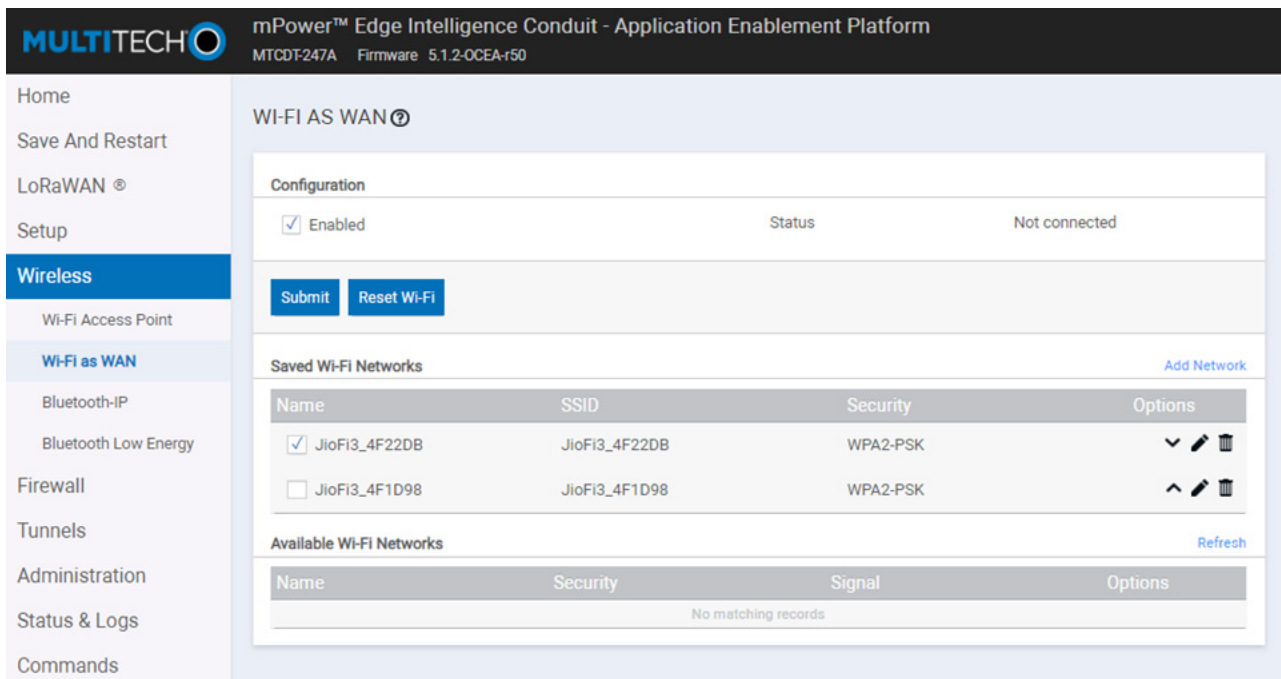


Figure 31. Wi-Fi as WAN

- c. Press “Test Lora Server” in Smart-Vue Pro LoRaWAN page after infrastructure is added in webapp (This only works once you configure your network properly (as described in the **Setting Up an Ethernet Connection** and **Setting Up a Wi-Fi Connection to your Network**).

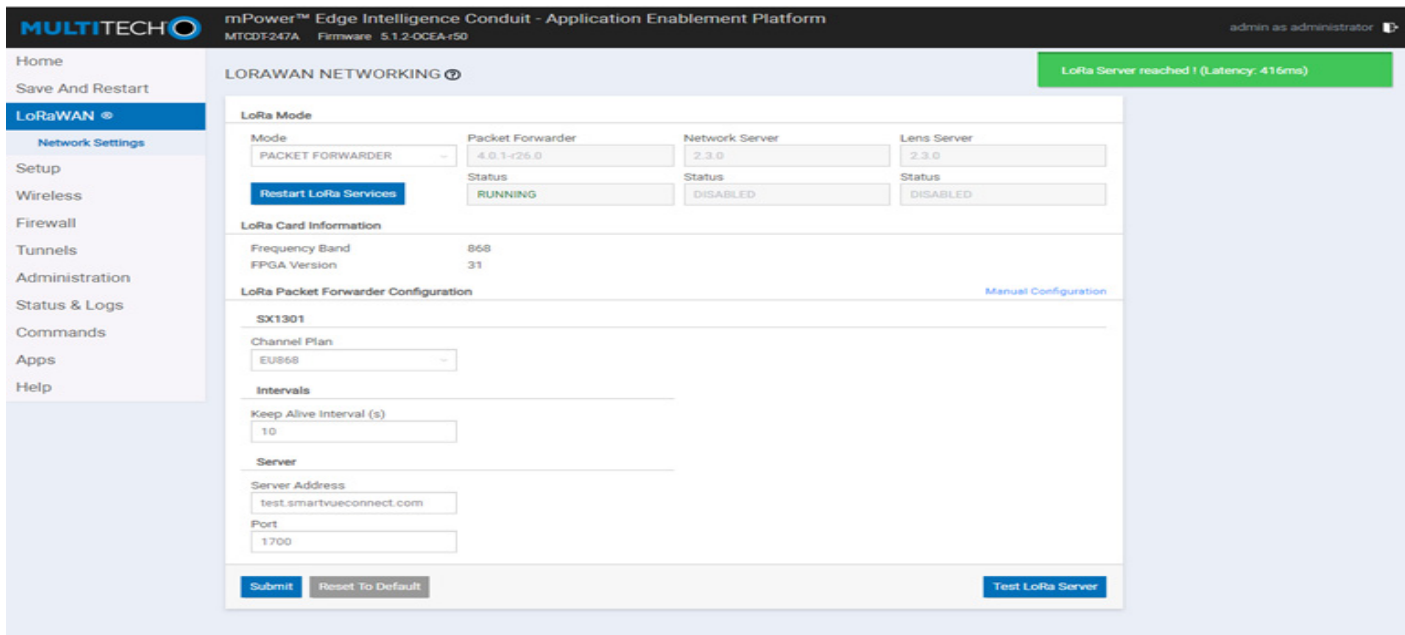


Figure 32. Smart-Vue Pro LoRaWAN Networking

5. Configuration in webapp.
 - a. In devices page, add the device name (Smart-Vue Pro Quatro) which is under test.

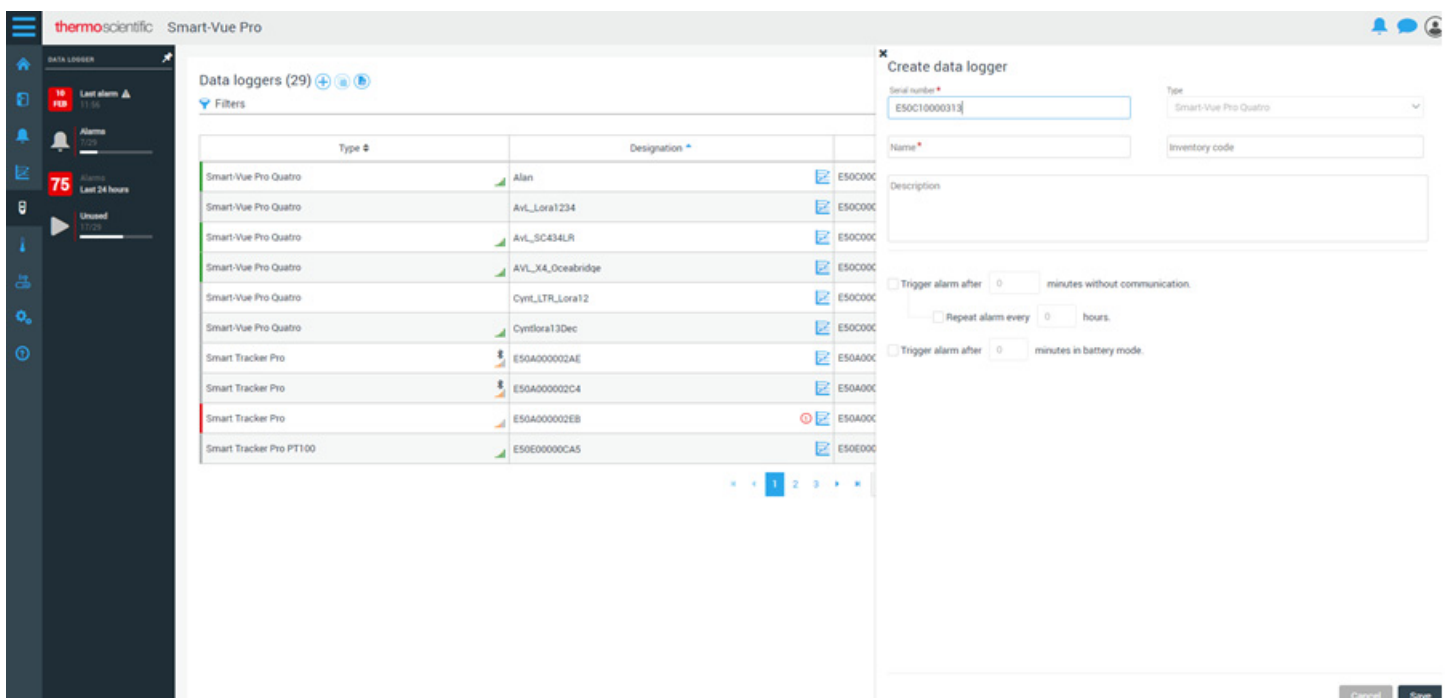


Figure 33. Add the device name

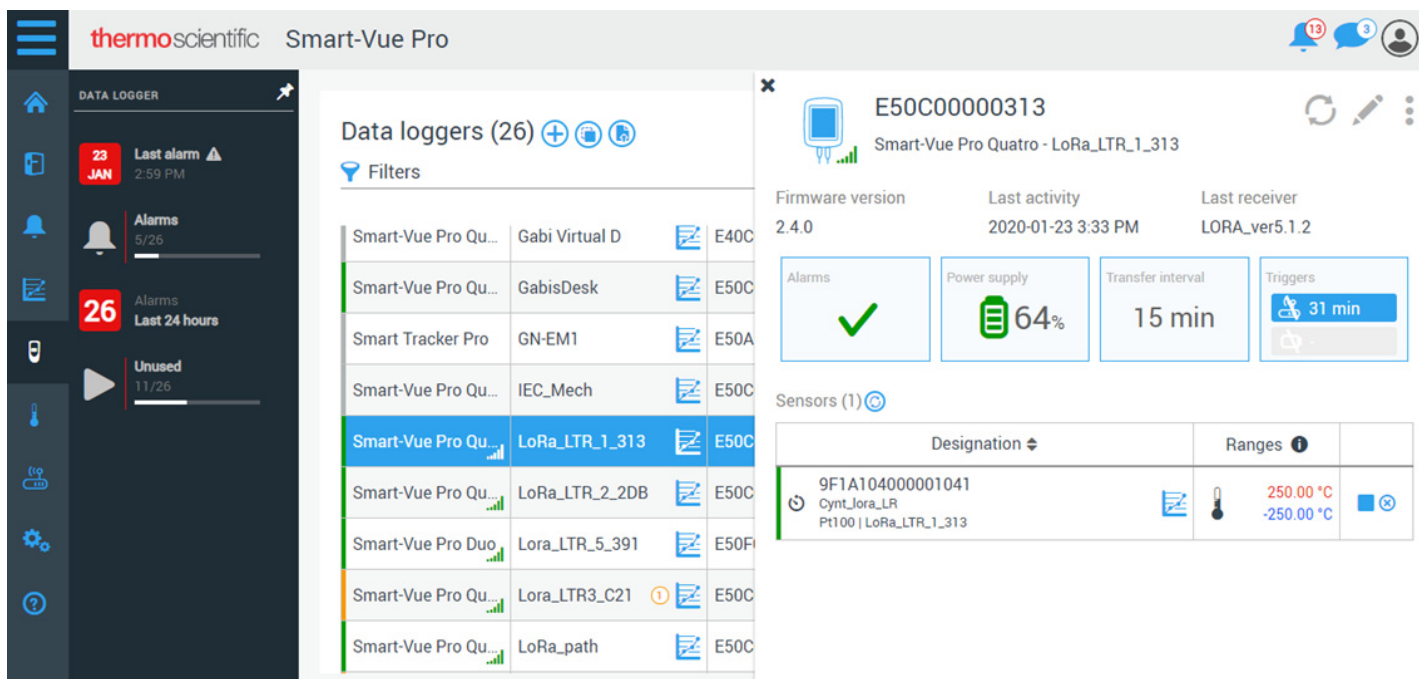


Figure 34. After adding the device

- In sensor page, verify the associated sensor serial number shown.
- In infrastructure, add the Smart-Vue Pro LoRaWAN details.

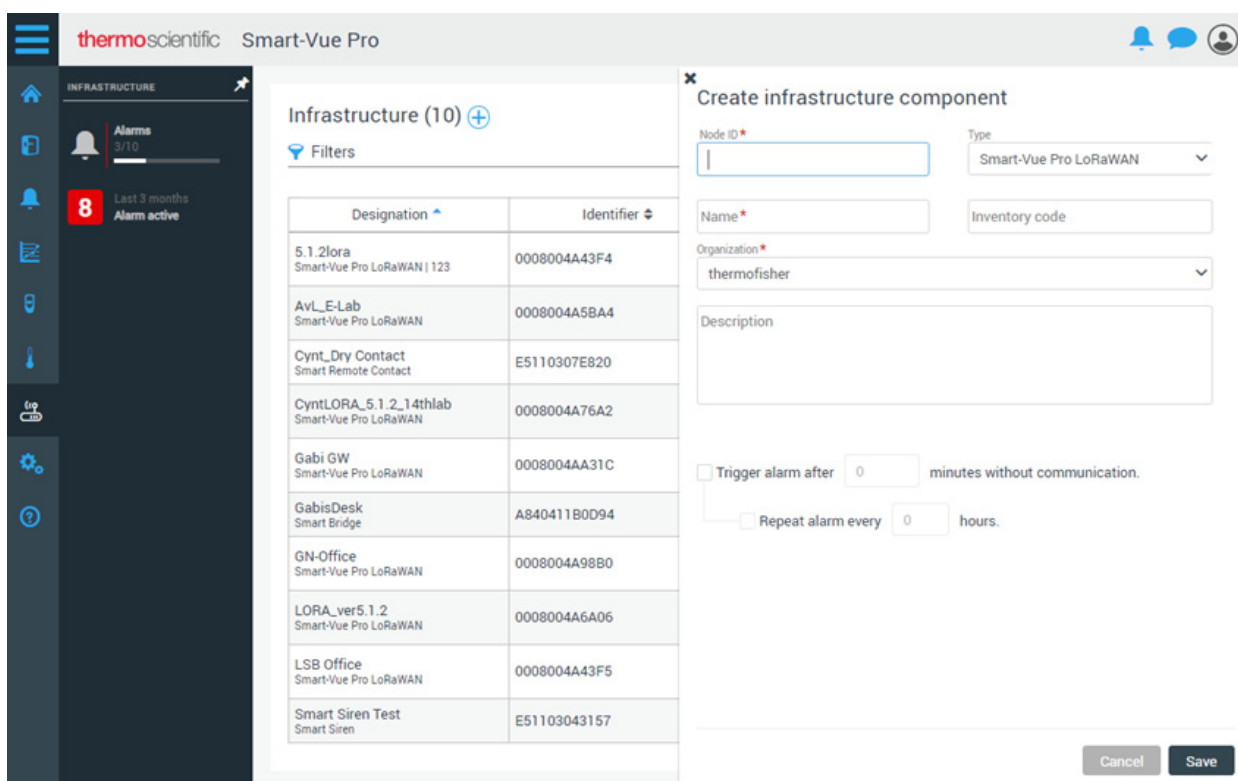


Figure 35. Add Smart-Vue Pro LoRaWAN details in Infrastructure

- d. To add a Smart-Vue Pro LoRaWAN, in serial number field, add the node id mentioned on the Smart-Vue Pro LoRaWAN label.

Infrastructure (9) +

Filters

Designation	Identifier
5.1.2lora Smart-Vue Pro LoRaWAN 123	0008004A43F4
AvL_E-Lab Smart-Vue Pro LoRaWAN	0008004A5BA4
Cynt_Dry Contact Smart Remote Contact	E5110307E820
CyntLORA_5.1.2_14thlab Smart-Vue Pro LoRaWAN	0008004A76A2
GabisDesk Smart Bridge	A840411B0D94
CN-Office Smart-Vue Pro LoRaWAN	0008004A98B0
LORA_ver5.1.2 Smart-Vue Pro LoRaWAN	0008004A6A06
LSB Office Smart-Vue Pro LoRaWAN	0008004A43F5

Edit infrastructure component

Node ID * 0008004A6A06 Type Smart-Vue Pro LoRaWAN

Name * LORA_ver5.1.2 Inventory code

Organization * AlertsTesting1

Description

Firmware version: - Last activity: 2020-01-23 3:38 PM

☒ Trigger alarm after 15 minutes without communication.

☐ Repeat alarm every 0 hours.

Cancel Save

Figure 36. Add the node ID in Smart-Vue Pro LoRaWAN label

- e. Create an equipment and then add the sensor to equipment which is recently added to device.

Equipment (28) +

Filters

Designation	Type
#116_equip_lora	box
Alan Office _E50C000c	
Alan Office _E50C000c	
AVL_ELAB_451	fridge
AvL_Lora	E-Lab Bench
CX2 50C000004C5	
CX2 50C000005B0	
Cynt_lora	normal testing 3
Cynt_Lora_331	
Gabi Digital	

Create equipment

Name *

Inventory code

Icon Type

Organization * thermofisher

Critical Not critical

Mobile Static

Cancel Save

Figure 37. Adding Sensor to equipment (1 of 2)

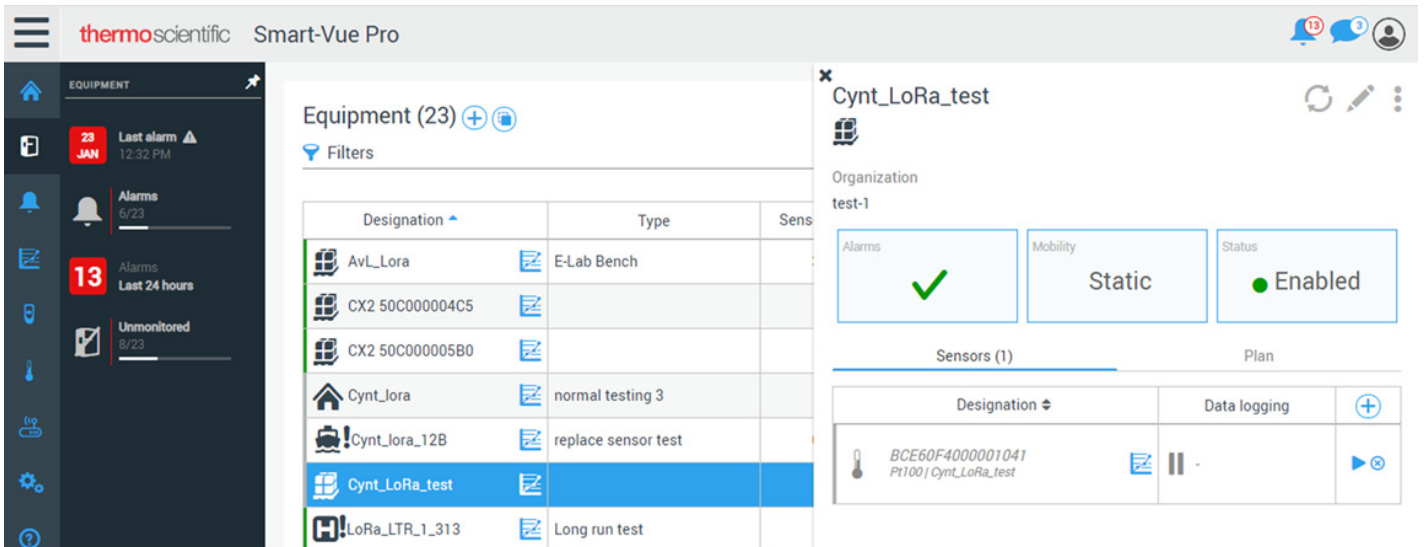


Figure 38. Adding Sensor to equipment (2 of 2)

- f. For adding sensor to equipment, open the particular equipment created and click + in equipment details page under sensor.

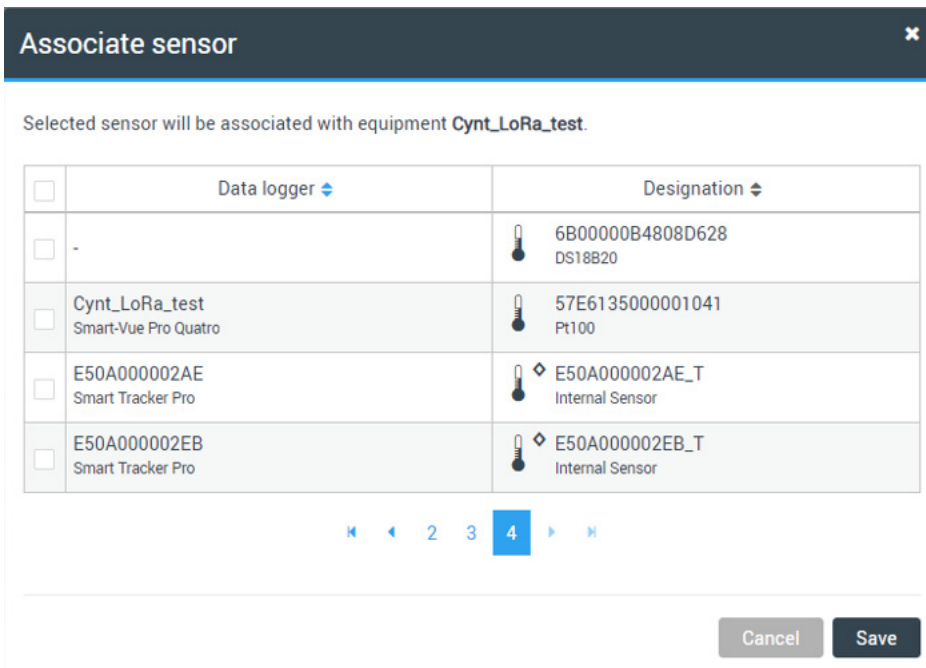


Figure 39. Associate sensor

- g. After successfully adding the equipment, device and testing the Smart-Vue Pro LoRaWAN coverage on Smart-Vue Pro Quatro, start the data logging from either Equipment page or Device page by clicking on the play icon.

Equipment (23)

Designation	Type	Sens
AvL_Lora	E-Lab Bench	
CX2 50C000004C5		
CX2 50C000005B0		
Cynt_lora	normal testing 3	
Cynt_lora_12B	replace sensor test	
Cynt_LoRa_test		
LoRa_LTR_1_313	Long run test	

Cynt_LoRa_test

Organization: test-1

Alarms: ☒ Mobility: Static Status: ● Enabled

Sensors (1)

Designation	Data logging
BCE60F4000001041 P1100 Cynt_LoRa_test	II

Figure 40. Equipment Page

Data loggers (26)

Type	Designation	Serial n
Smart-Vue Pro Quatro	AvL_Lora1234	E50C000004
Smart-Vue Pro Quatro	AvL_SC434LR	E50C000004
Smart-Vue Pro Quatro	AVL_X4_Oceabridge	E50C000004
Smart-Vue Pro Quatro	Cynt_LoRa_test	E50C00000B
Smart-Vue Pro Quatro	Cynt_LTR_Lora12	E50C000003
Smart-Vue Pro Quatro	Cyntlora13Dec	E50C000003
Smart Tracker Pro	E50A000002AE	E50A000002

E50C00000BCB
Smart-Vue Pro Quatro - Cynt_LoRa_test

Firmware version: 2.4.0 Last activity: 2020-01-23 4:05 PM Last receiver: LORA_ver5.1.2

Alarms: ☒ Power supply: 96% Transfer interval: 20 min Triggers: ☐

Sensors (2)

Designation	Ranges
57E6135000001041 P1100 Cynt_LoRa_test	250.0 °C -250.0 °C
BCE60F4000001041 Sc-Demo-BcB P1100 Cynt_LoRa_test	250.0 °C -250.0 °C

Figure 41. Data loggers page

- h. We can see the equipment thumbnail in home screen when data logging is in progress.



Note: For more detail refer the user guides of the data logger.

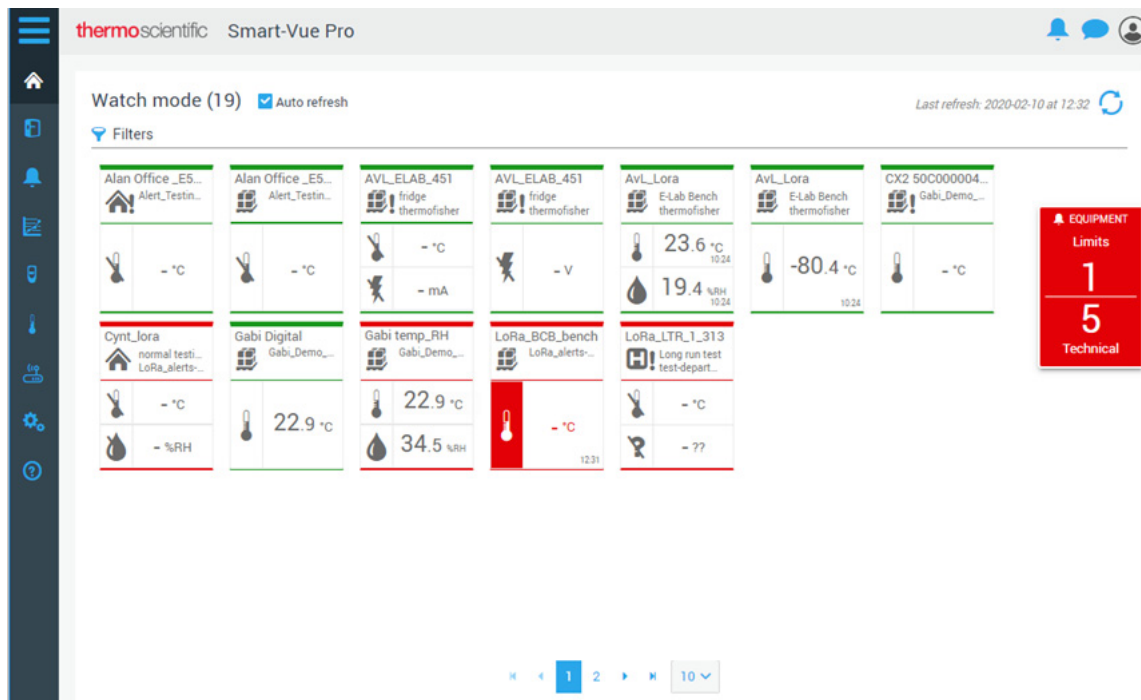


Figure 42. Watch Mode

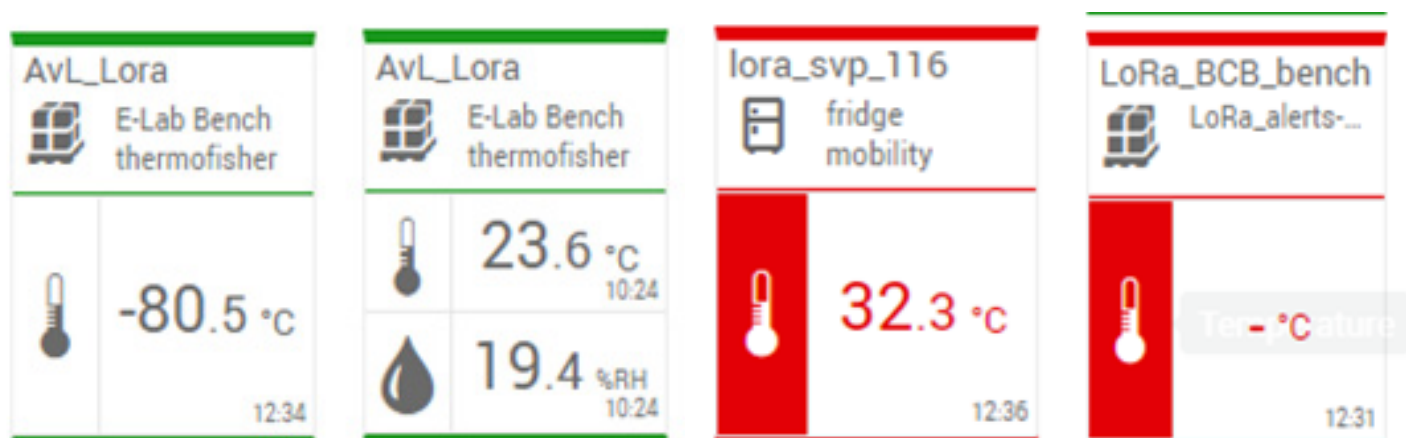


Figure 43. Tiles on watch mode

Appendix - Product Specifications

Smart-Vue Pro LoRaWAN Enabled Receiver Features

- Internet of Things connectivity featuring Smart-Vue Pro LoRaWAN wireless connectivity
- Long asset management range – up to 10 miles/16 km line-of-sight; 1-3 miles/2 km within buildings.

Note: This represents ideal network configuration and equipment set up. Results vary depending on various technical considerations, as well as RF interference and obstruction type (e.g., metal, cement, etc.)


- Ethernet: One RJ-45 Ethernet 10/100 port
- Wi-Fi (optional) via integrated data logger
- Input voltage: 9 V - 32 V

- Dimensions (L x W x H): 6.35" x 4.23" x 1.69" (161.3 mm x 107.4 mm x 42.8 mm)
- Weight: About 16.2 oz. (459 g)
- Chassis Type: Metal
- Operating Temperature: -30°C to +70°C
- Storage Temperature: -40°C to +85°C
- Relative Humidity: 20% to 90%, non-condensing

Related Products

- Smart-Vue Pro sensor monitoring solution (web application).
- Compatible Smart-Vue Pro LoRaWAN enabled wireless sensor modules.

Table 3. Back Panel Connectors

Label	Description	Remarks
CELL, AUX	Cellular antenna inputs CELL: Primary AUX: Rx Diversity/MIMO	Not used.
AP1, AP2	Slots for MultiTech accessory cards.	AP1 slot used for LoRa radio antenna. AP2 slot not used.
USB DEVICE	User-defined, high-speed 480 Mbps, standard USB 2.0 Micro B Connector.	Not used.
E-NET	RJ-45 receptable for standard Ethernet 10/100 Base-T.  CAUTION: Ethernet ports and command ports are not designed to be connected to a public telecommunication network.	Used for connecting to the ethernet cable.
USB HOST	High-speed, standard USB 2.0 Type A connector. 500 mA maximum current draw.	Not used.
POWER	9.32 V DC power receptable for provided power cord.	Used for connecting to the power supply.

Note: Power Source

When setting up the Smart-Vue Pro LoRaWAN Enabled Receiver (antenna), connect to a power source that is grounded with surge protection as well as on a battery backup, if at all possible. This ensures the availability of the functionality of the device.

Table 4. Front Panel Connectors

Label	Name	Description
PWR	Power	Solid (constant green) if unit is on, indication that power is present.
STATUS	Power Status	Default Condition: LED blinks when mLinux is fully loaded.
LS	Link Status	Varies with radio model.
CD	Carrier Detect	This LED is on when a cellular data connection is made. Present on the Conduit Application model only.
Signal	Signal Strength	These 3 LEDs display the strength of the cellular signal. Present on the Conduit application model only.

Installation Details

The LoRaWAN module is a critical device in your wireless monitoring system. Therefore, we recommend that you keep it plugged it into an interruptible power supply or safety plug.

Smart-Vue Pro LoRaWAN enabled receiver may be placed on a flat surface or mounted using the mounting holes located at bottom of the casing.



Figure 44. Back/bottom view of LoRaWAN gatewaydata logger

WEEE Compliance

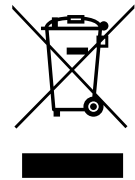
WEEE Compliance. This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at www.thermofisher.com under Services & Support.

WEEE Konformität. Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliedstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter www.thermofisher.com unter Services & Support.

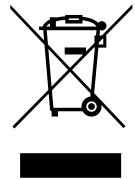
Conformità WEEE. Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2012/19/EU in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito www.thermofisher.com in Servizi e Supporto.

Conformité WEEE. Ce produit doit être conforme à la directive européenne (2012/19/EU) des Déchets d'Équipements Électriques et Électroniques (DEEE). Il est marqué par le symbole suivant. Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur www.thermofisher.com sous Services et Assistance.

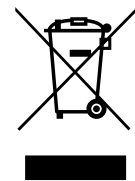
Great Britain



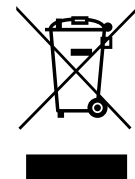
Deutschland



Italia



France





IF YOU NEED ASSISTANCE:

Thermo Fisher Scientific products are backed by a global technical support team ready to support your applications. We offer cold storage accessories, including remote alarms, temperature recorders, and validation services.

Visit [thermofisher.com](https://www.thermofisher.com) or call:

Countries	Sales	Services
North America	+1 866 984 3766	(800) 438-4851
India	1800 22 8374, +91 22 6716 2200	+91 22 6716 2200
China	+800 810 5118, +400 650 5118	+8621 68654588
Japan	+81 3 5826 1616	+81 3 3816 3355
Australia	+61 39757 4300	1 300 735 292
Austria	+43 1 801 40 0	+43 1 801 40 0
Belgium	+32 53 73 42 41	+32 2 482 30 30
France	+33 2 2803 2180	+33 2 2803 2180
Germany	0800 1 536 376, +49 6184 90 6000	0800 1 536 376
Italy	+32 02 95059 552	+39 02 95059 552, 432 254 375
Netherlands	+31 76 579 55 55	+31 76 571 4440
Nordic/Baltic/CIS	+358 9 329 10200	+358 9 329 100
Russia	+7 812 703 4215	+7 812 703 4215
Spain/Portugal	+34 93 223 09 18	+34 93 223 09 18
Switzerland	+41 44 454 12 22	+41 44 454 12 12
UK/Ireland	+44 870 609 9203	+44 870 609 9203
New Zealand	+64 9 980 6700	+64 9 980 6700
Other Asian Countries	+852 2885 4613	+852 2885 4613
Countries not listed	+49 6184 90 6000	+49 6184 90 6000

Thermo Fisher Scientific Inc.
275 Aiken Road
Asheville, NC 28804
United States

Find out more at [thermofisher.com/cold](https://www.thermofisher.com/cold)

ThermoFisher
S C I E N T I F I C