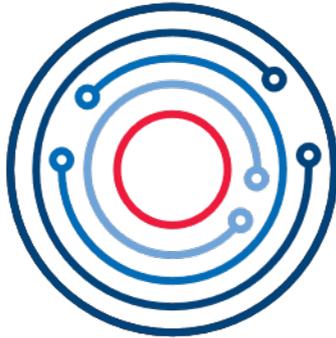




January 2025

QIAsphere[®] Connectivity Package Quick-Start Guide



QIAsphere

Table of Contents

Introduction	3
Safety Information	5
External Feature of QIAsphere Base	6
Site Requirements	7
Installation Procedure	8
QIAsphere Setup	8
Requirements on site	8
Connect QIAsphere Base to Local Network	9
Wireless network connection	9
Wired Ethernet connection (with DHCP) – recommended	9
Connect instrument to local network and QIAsphere Base	10
Setup QIAsphere Workgroup	11
Document Revision History	12

Introduction

QIASphere connects users with QIAGEN's comprehensive digital ecosystem to deliver a unique user experience and improve laboratory efficiency and safety through cloud-based connectivity. The QIASphere Base (IoT* gateway device) of the QIASphere Connectivity Package connects QIASphere-ready instruments to the QIASphere App running on mobile devices or desktop PCs, enabling the user for an increasing number of applications such as remote instrument monitoring.

Using QIASphere, you can immediately enhance your lab experience:

- **Stay connected** – Monitor your instrument(s) while being network independent. The QIASphere App can access the QIASphere Cloud from anywhere.
- **Monitor instrument status with mobile devices or from desktop PC** – either via the mobile app for iOS and Android devices or the web app version for your desktop.
- **Full walk-away convenience** – Keep track of instrument status, such as run completion, through push notifications.
- **Faster service response** – Creating and sending service packages via e-mail is no longer necessary for instruments connected to QIASphere (for devices that support it).
- **QIASphere Insights** – Explore cloud-based analytic capabilities by means of epidemiology monitoring reports that deliver statistics on pathogens' detection and system usage reports that include kit consumption and instrument usage of your instruments connected to QIASphere.
- **Connect to multiple QIASphere Bases** – Monitor instruments in your QIASphere App that are connected to different local networks and to QIASphere Bases in the same networks.

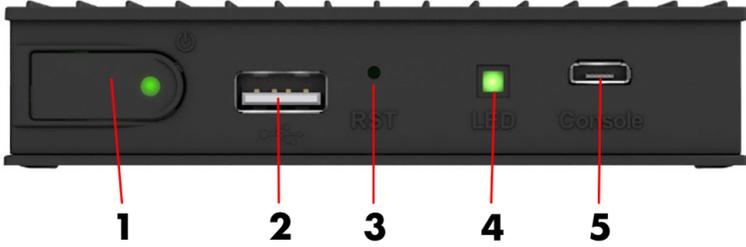
*Internet of Things

Note: QIAGEN® does not collect any personal identification or sensitive data such as patient health data when using QIASphere cloud-connected mode. QIASphere Cloud is used only to provide you with fully functional instrument monitoring and other cloud services. For more information on cyber security, see our cyber security technical information to be downloaded from www.qiagen.com/qiasphere.

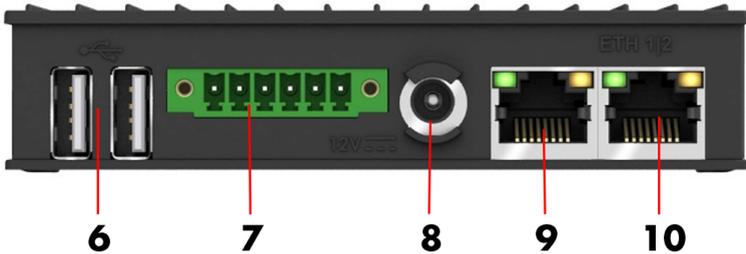
Safety Information

Before installing or using the QIAsphere Base, it is essential that you carefully read the user manual provided by the hardware supplier, CompuLab, and pay attention to their safety information. The instructions and safety information in this user manual must be followed to ensure safe operation of the instrument and to maintain the instrument in a safe condition. The safety information for QIAsphere Base (CompuLab – IOT-GATE-iMX8) can be found at <https://www.complab.com/products/iot-gateways/iot-gate-imx8-industrial-arm-iot-gateway>.

External Feature of QIAsphere Base



Front view of QIAsphere Base.



Rear view of QIAsphere Base.

- | | | | |
|---|--|----|--|
| 1 | Power button with LED power indicator and factory reset button | 6 | Two USB ports |
| 2 | USB port | 7 | RS485/RS232 serial interface |
| 3 | Reset button | 8 | Power cord socket |
| 4 | Status three-color LED | 9 | ETH1 Ethernet port to connect personal computer to QIAsphere Base – denoted as a service port |
| 5 | Debug console | 10 | ETH2 Ethernet port to connect QIAsphere Base to local network – denoted as a connectivity port |

Site Requirements

QIAsphere Base must be connected to the same network to which QIAsphere-ready instruments are to be connected. QIAsphere Base does not need to be located in close proximity to ensure a proper connection.

Installation Procedure

This section provides you important information on how to install the QIAsphere Base and to set up QIAsphere. Depending on your preference, you can connect QIAsphere Base wireless or via Ethernet cable to your local network. The described installation procedure in this Quick Start Guide comprises the configuration of QIAsphere Base using a personal computer.

If you need detailed installation information, or if you want to use QIAsphere App to configure QIAsphere Base, refer to the QIAsphere user manual. Please check and update your QIAsphere Base to the current software using update packages and the Update Guide provided at www.qiagen.com/qiasphere.

Depending on your local network settings, you may need to consult your IT department. If required, please share the *QIAsphere User Manual* and our technical information about cybersecurity with your IT specialist for detailed information on network architecture and IT prerequisites, see www.qiagen.com/qiasphere.

QIAsphere Setup

Requirements on site

- My QIAGEN account: Sign up on www.qiagen.com/myqiagen
- QIAsphere License: On **My QIAGEN** account, go to **QIAsphere Administration** widget.

Note: If a license already exists, the main screen is displayed. If no license is available, you can request for it.

- Download QIAsphere App.

Note: Mobile app can be downloaded from the App Store or Google Play Store.

Connect QIAsphere Base to Local Network

Wireless network connection

1. Plug in Wi-Fi USB adapter into QIAsphere Base. Power on the QIAsphere Base by plugging in the QIAsphere Base power cord and pressing the power button.
2. Wait for approximately 2 minutes until User 1 LED is on. Then, connect your PC to the QIAsphere Base **ETH1** Port via Ethernet cable.
3. On your PC web browser (e.g., Chrome™ or Mozilla® Firefox) open QBSP* by entering this URL: **https://172.16.0.1**
4. Log in with default user name admin and default password **Qiagen@123**.
5. Change password as prompted.
6. Go to **Network configuration**. Select the desired Wi-Fi network and enter the Wi-Fi password. Then, click **Connect**.
7. QIAsphere Base is successful connected. Note down QIAsphere Base **IP address** from the QBSP main screen, for example, 192.168.1.10.
8. Click **Log out** on the side menu and close the browser.
9. Remove the cable between PC and **ETH1** port.

Wired Ethernet connection (with DHCP) – recommended

Note: Most Ethernet networks support automated IP address assignment using DHCP. In some special cases, networks do not support DHCP. In this case, static IP settings can be configured. See the *QIAsphere User Manual* for details.

*QIAsphere Base Setup Portal. QIAsphere Base uses self-assigned certificate. Therefore, the browser may issue a security warning. Ignore the warning or add an exception to access QSBP.

1. Connect QIAsphere Base ETH2 port to a wired network using the Ethernet cable. Power on the QIAsphere Base by plugging in the QIAsphere Base power cord. Power on the QIAsphere Base by plugging in the QIAsphere Base power cord and pressing the power button.
2. Wait for approximately 2 minutes until **User 2** LED is on. Then, connect your PC to the QIAsphere Base ETH1 Port via Ethernet cable.
3. On your PC web browser (e.g., Chrome or Mozilla Firefox) open QBSP* by entering this URL: <https://172.16.0.1>
4. Log in with default user name **admin** and default password **Qiagen@123**.
5. Change password as prompted.
6. Follow instructions from the installation wizard to connect QIAsphere Base to cloud.
7. After finishing the set-up, note down QIAsphere Base IP address from the QBSP main screen, for example, 192.168.1.10.
8. Click **Log out** on the side menu and close the browser.
9. Remove cable between PC and **ETH0** port.

Connect instrument to local network and QIAsphere Base

- QIAsphere-ready instrument must be connected to the same local network as the QIAsphere Base.
- See instrument user manual for instructions on how to connect to local network and to the QIAsphere Base (QIAsphere Base IP address is required).

*QIAsphere Base Setup Portal. QIAsphere Base uses self-assigned certificate. Therefore, the browser may issue a security warning. Ignore the warning or add an exception to access QBSP.

Setup QIASphere Workgroup

1. Log in to **My QIAGEN**, and select the **QIASphere Administration** widget.
2. Go to **Workgroups** menu, and click **Add Workgroup** button.
3. Enter the workgroup name as prompted, and click **Add**.
4. Go to **Instruments** menu, and select a relevant instrument or QIASphere Base.
5. Click on the pencil icon to edit instrument details.
6. Ensure that the QIASphere connection shows **Enabled**. Click the pencil icon in the Workgroup name field, then select a relevant workgroup and click **Save**.
7. Go to Users menu and click **Add User** button.
8. Enter required information about the new user as prompted. Then, select a workgroup and click Invite.
9. New user receives invitation email. By clicking Activate your **QIASphere Account** button in the email, the user is added to QIASphere. The user can launch now the QIASphere App and see instruments that are assigned to the user's workgroup (My QIAGEN account required).

Document Revision History

Revision	Description
January 2025	Updated the document to new hardware of QIASphere Base. Cleaned up the Local Mode.
August 2022	Updated the document based on the new wizard process. Deleted the Local Mode section.

Trademarks: QIAGEN®, Sample to Insight®, QIASphere® (QIAGEN Group); Chrome™ (Google LLC); Mozilla® (The Mozilla Foundation). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

HB-2842-004 01/2025 © 2025 QIAGEN, all rights reserved.

