

## Connecting Imaging Instruments with Image Lab Touch to a Local Computer

Introduction.....	1
Required Components.....	1
Specifying a Link-Local IP Connection on the Instrument.....	1
Specifying a Link-Local IP Address for the Computer.....	2
Creating a Shared Folder on the Computer.....	3
Assigning Specific Users to the Shared Folder.....	4
Exporting Files to the Shared Folder.....	4
Legal Notices.....	5
Contacting Technical Support.....	5

### Introduction

This document explains how to connect your imaging instrument, with Image Lab Touch, **v3.0 and later**, installed, to a shared folder on a computer that does not have a network connection. Once your local connections are established, you can export images from the instrument to the shared folder.

### Required Components

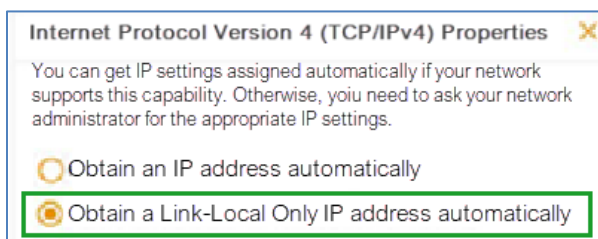
- ChemiDoc MP or GelDoc Go instrument
- Windows 10 computer
- Ethernet cable

### Specifying a Link-Local IP Connection on the Instrument

Before you can connect a computer to the instrument, you must specify a Link-Local IP address as part of the instrument's network configuration.

#### To specify a Link-Local IP address

1. Turn on the instrument and log in as an administrative user.
2. Tap the gear icon to open the Settings menu.
3. Tap Network Configuration and then tap Obtain a Link-Local Only IP address automatically.



4. Close the settings screens and turn off the instrument.

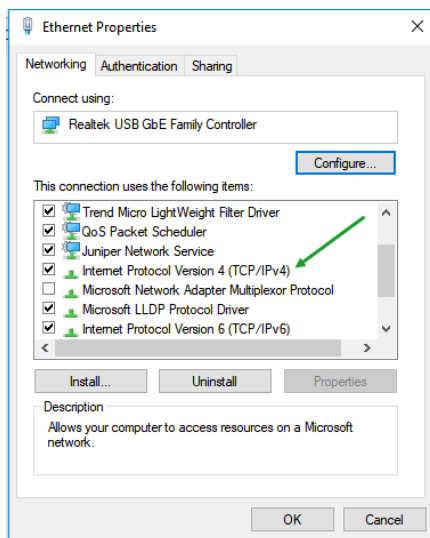
After you complete the remaining setup and turn on the instrument again, a Link-Local IP address will appear in the About This Imager screen.

## Specifying a Link-Local IP Address for the Computer

After setting up the instrument for a Link-Local IP address, complete the steps below to create an IP address for the computer.

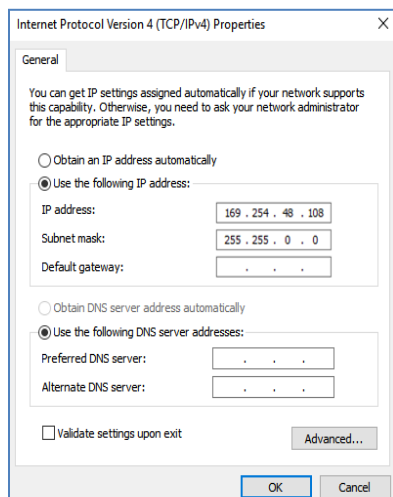
### To create a Link-Local IP address on the Windows computer

1. Turn on the computer and log into Windows using an administrator account.
2. Click the Windows icon, and then click the gear icon to open the Settings.
3. Click Network and Internet to display the Status window (default).
4. Click Change adapter options to open the Network Connections window, and then double-click Ethernet.
5. In the Ethernet Properties dialog box, scroll to and select Internet Protocol Version 4 (TCP/IPv4), and then click Properties.



6. In the Internet Protocol Version 4 (TCP/IPv4) Properties dialog box, select **Use the following IP address** and then enter an IP address beginning with 169.254.

**Tip:** You can use any numbers for the remaining parts of the address, since the connection is not part of your institution's network. The following example uses 169.254.48.108.



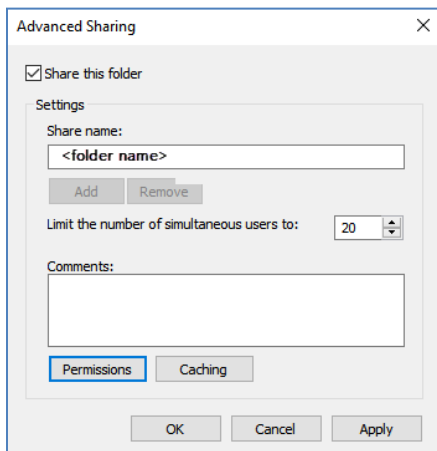
7. For the Subnet mask, enter 255.255.0.0.
8. Click OK and then click OK again to close the dialog boxes.
9. Close the Network Connections window.

## Creating a Shared Folder on the Computer

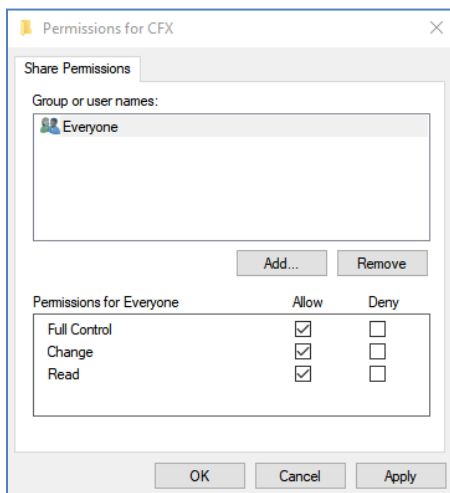
You must also create a shared folder on the computer, where the image files exported from the instrument will be stored.

### To create the folder to contain exported images

1. On the computer, create a folder in the location where you want to store your images.
2. Select the folder and right-click, and then select Properties to display the folder Properties dialog box.
3. Select the Sharing tab and click Advanced Sharing.



4. Select the Share this folder checkbox.
5. Enter a folder name in the Share name field and click Permissions.
6. Select the Full Control checkbox.



7. Do one of the following:

- If everyone will require access to the folder, leave the Everyone default and click OK, and then click OK again twice to close the remaining dialog boxes. Go to [Exporting Files to the Shared Folder](#).
- If only certain users require access, click Add and continue to [Assigning Specific Users to the Shared Folder](#).

## Assigning Specific Users to the Shared Folder

**Important:** The users requiring access must be set up as Windows accounts before you can add them as users with folder permissions. For information, contact a system administrator.

### To add Windows users

1. For each user, enter the name in the Object Names field, and then click Check Name.  
If the Windows account has been set up and the user name is valid, it appears as verified in the field.
2. When you are finished, click OK, and then click OK twice to close the remaining dialog boxes.

## Exporting Files to the Shared Folder

### To export files to the shared folder

1. Connect the computer to the instrument using an Ethernet cable.
2. Turn on the instrument and log in as an administrator.
3. Tap the gear icon.
4. Tap About This Imager and verify that a 169.254.x.x IP address appears for IPv4.  
**Note:** This IP address is the Link-Local address for the instrument, and is different than the address you configured earlier on the computer.
5. Close the settings and tap Gallery.
6. Select the images to export and tap Send/Save.
7. Click Save to Network to open the corresponding dialog box.
8. Enter the computer IP address in the field and tap Connect.
9. In the resulting dialog box, enter your Windows user name and password.  
**Note:** Enter your user name in the following format:  
<computer name\user name>  
The shared folder appears in the Save to Network dialog box.
10. Select the folder and tap Save to Folder.  
For future exports, the folder appears for selection in the Recents list.

## Legal Notices

No part of this publication may be reproduced or transmitted in any form or by any means, electronic or mechanical, including photocopy, recording, or any information storage or retrieval system, without permission in writing from Bio-Rad Laboratories, Inc.

Bio-Rad reserves the right to modify its products and services at any time. This guide is subject to change without notice. Although prepared to ensure accuracy, Bio-Rad assumes no liability for errors or omissions, or for any damage resulting from the application or use of this information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

All trademarks used herein are the property of their respective owner.

Copyright © 2020 Bio-Rad Laboratories, Inc. All rights reserved.

## Contacting Technical Support

Bio-Rad Technical Support in the United States is open Monday–Friday, 5:00 AM to 5:00 PM, Pacific Time.

**Phone:** 1-800-424-6723, option 2

**E-mail:** Support@bio-rad.com (U.S./Canada Only)

For technical assistance outside the U.S. and Canada, contact your local technical support office or click the Contact us link at [www.bio-rad.com](http://www.bio-rad.com).