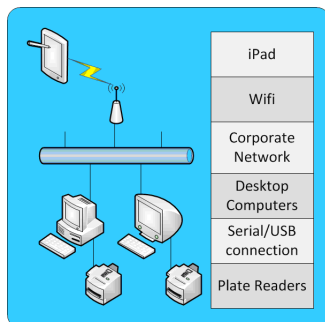


Setting Up Access to SoftMax Pro Software from an iPad Mobile Digital Device

Introduction

You can use an iPad® Mobile Digital Device and wireless network access to remotely control SoftMax® Pro Microplate Data Acquisition and Analysis Software that is installed on a computer on the network. You can create protocols, run experiments, and analyze acquired data from anywhere within range of your wireless access point.



Requirements

To remotely control SoftMax Pro Software from an iPad Device, you need the following items:

- iPad Mobile Digital Device
- iTunes® account (credit card required)
- Jump Desktop App (purchased separately) installed on the iPad Device
- Wireless network access
- Host computer with network access
- SoftMax Pro Software installed on the host computer

The iPad Device accesses the network through a wireless access point set up by your system administrator. Your networked host computer needs to be set up to allow remote access or screen sharing.

- For information on network access and permissions, contact your system administrator.
- For information on using the iPad Device, contact Apple support at <http://www.apple.com/support/ipad>.
- For information on using the Jump Desktop App, contact Jump Desktop support at <http://jumpdesktop.com>.

Setting Up the Host Computer

The Jump Desktop App on an iPad Device can remotely access an Apple or Windows computer. Follow the setup instructions for the computer that has the SoftMax Pro Software installed on it.

Setting Up Access to an Apple Computer

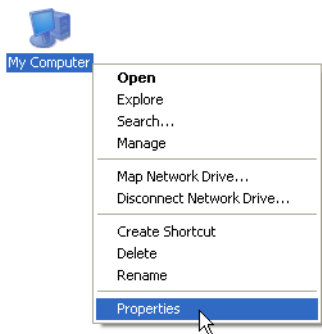
1. Open **System Preferences**.
2. Click the **Sharing** preference.
3. Select the **Screen Sharing** service.
4. Close the Sharing preference screen.

To get the IP address of your Apple Computer

1. Open **System Preferences**.
2. Click the **Network** preference.
3. Make a note of the **IP Address**.
4. Close the Network preference screen.

Setting Up Access to a Windows Computer

1. Right-click **My Computer**.
2. Click **Properties**.
3. Click the **Remote** tab.
4. Select **Allow user to connect remotely to this computer**.
5. Click **OK**.



To get the name of your Windows Computer

1. Right-click **My Computer**.
2. Click **Properties**.
3. Click the **Computer Name** tab.
4. Make a note of the **Full computer name**.
5. Click **OK**.

Setting Up the iPad Device

Connecting the iPad Device to the Network

1. Start the **Settings** App.
2. From the menu on the left, tap the **Wi-Fi** option.
3. Tap the wireless network and type the access keys provided by your system administrator.
4. Return to the iPad Desktop.

Accessing or Creating an iTunes Account

1. Start the **Settings** App.
2. From the menu on the left, tap the **Store** option.
3. Tap **Sign In**.
4. If you already have an iTunes account, tap **Use Existing Apple ID**. Otherwise, tap **Create New Apple ID**.
5. Follow the on-screen instructions.

Getting the Jump Desktop App

1. Start the **App Store** App.
2. In the upper-right corner of the AppStore screen, tap the **Search** box and search for the **Jump Desktop** App.
3. Purchase and install the Jump Desktop App.

Accessing the Host Computer

1. Start the **Jump Desktop** App.
2. In the upper-right corner of the Jump Desktop screen, click the **+** button.
 - ♦ To access an Apple computer, tap the **VNC** option.
 - ♦ To access a Windows computer, tap the **RDP** option.
3. Type the IP address or name of the computer. See [Setting Up the Host Computer on page 2](#).
4. Tap **Save**.
5. Tap **Done**.
6. Tap **Jump**.

Controlling the Software from the iPad Device

Connecting to the Host Computer

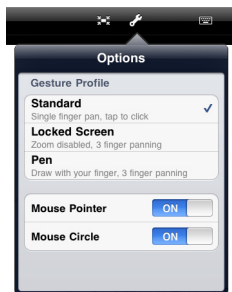
1. From the iPad Device, start the **Jump Desktop** App.
2. Tap the host computer IP address or computer name.
3. Logon to the host computer, if required.

Running the SoftMax Pro Software

1. Start the SoftMax Pro Software.
2. Use pinch and other iPad gestures to zoom and navigate the SoftMax Pro Software interface.

Most of the time, you can use the **Standard** Gesture Profile.

To make it easier to select a group of wells in a plate template, you can temporarily switch to the **Pen** Gesture Profile.



Disconnecting from the Host Computer

1. From the top of the Jump Desktop screen, tap the white **X**.
2. Tap **Disconnect**.



For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, Inc. or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, Inc.

iPad and iTunes are a registered trademarks of Apple Inc.

© 2011 Molecular Devices, Inc.

All rights reserved
