



Luminometers

Fluorometers

Multimode
ReadersDNA
PurificationApplication
MethodsProduct
Selector ToolProduct
VideosFind A
DistributorOEM
OpportunitiesCustomer
Care Center

Modulus Microplate Multimode Reader

This Technical Note provides instruction for using an iPod as a mass storage device with the Modulus™ Microplate Multimode Reader

iPod™ configuration

1. Download and install the latest version of iTunes™ from <http://www.apple.com/itunes/>
2. Connect the iPod to your computer using the supplied USB cable.
3. Open iTunes if it does not automatically open.
4. On the left hand menu bar, under devices, click on the iPod icon.
5. Choose the device summary tab.
6. **CHECK** the box next to **ENABLE DISC USE**.
7. Click on **APPLY** in the bottom right corner.
8. **EJECT** the iPod from your computer by selecting the eject arrow to the right of the iPod name on the menu bar or by selecting the eject icon on the bottom right of the screen.

Transferring data from the Modulus™ Microplate Multimode Reader to the iPod

1. Using the supplied USB cable, attach the iPod to the USB drive on the left of the touch screen of the Modulus™ Microplate Multimode Reader.
2. Wait for the USB icon to appear on the instrument status bar (depending on the storage device, this may take several minutes).
3. Touch the **RESULTS** button on the **INSTRUMENT CONTROL** screen.
4. Touch the **FILES** button and highlight the file(s) to be exported.
5. Touch the **USB** button under the **TOOLS** button.
6. Touch the **OK** button to export only the selected file(s).
7. Touch the **COPY ALL FILES** button to export all of the results files.
8. Disconnect the iPod by removing the cable from the USB drive.

Transferring data from the iPod to your computer

1. Move the iPod to your local PC or Mac computer.
2. Find the folder by opening the **MY COMPUTER** system folder.
3. Click on the iPod name under the Devices with Removable Storage heading.
4. The .csv formatted file(s) will be saved onto the iPod in the folder named Modulus™ Microplate.
 1. For Windows, double-click to open the file in Excel.
 2. For Mac, click to open the file and import the data into Excel.

TRADEMARKS

Modulus™ is a trademark of Turner BioSystems, Inc.
iPod™ and iTunes™ are registered trademarks of Apple, Inc.

TURNER BIOSYSTEMS CONTACT INFORMATION

Phone: (408) 636-2400 or
Toll Free: (888) 636-2401 (US & Canada)
Fax: (408) 737-7919
Web Site: www.turnerbiosystems.com

[About Promega](#) • [Support](#) • [Site Map](#) • [Distribution](#) • [Grants Programs](#) • [Sunnyvale Office Career Opportunities](#)

Promega Corporation
Sunnyvale Office • Detection Instruments
645 N Mary Avenue • Sunnyvale CA 94085