



MetaXpress[®] 6 FAQ

Changing the Fluidics Tip Adapter and Type

Date Revised 07/07/15 Version B



Frequently Asked Questions Overview

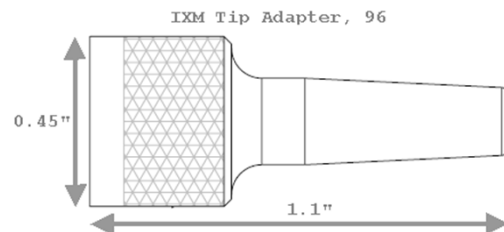
The purpose of this FAQ guide is to help the user change between 96-format and 384-format pipette tips for the ImageXpress Micro with Fluidics (pipetting) chapter.



Fluidic Tip Types and Adapters

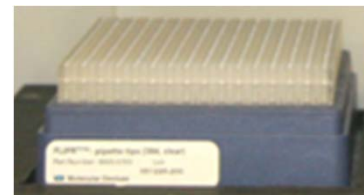
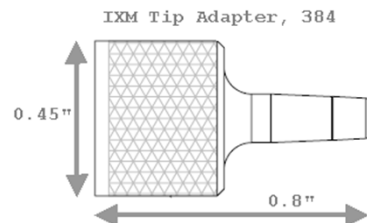
96-format tips

- 5-200 ml
- Tip Part **9000-0761** FLIPR Tetra Pipette tips, clear, non-sterile, 96-well 50 racks/case
- Replacement Adapter Part **2300-2460** ImageXpress Micro Fluidics Tip Adapter, 96-well



384-format tips

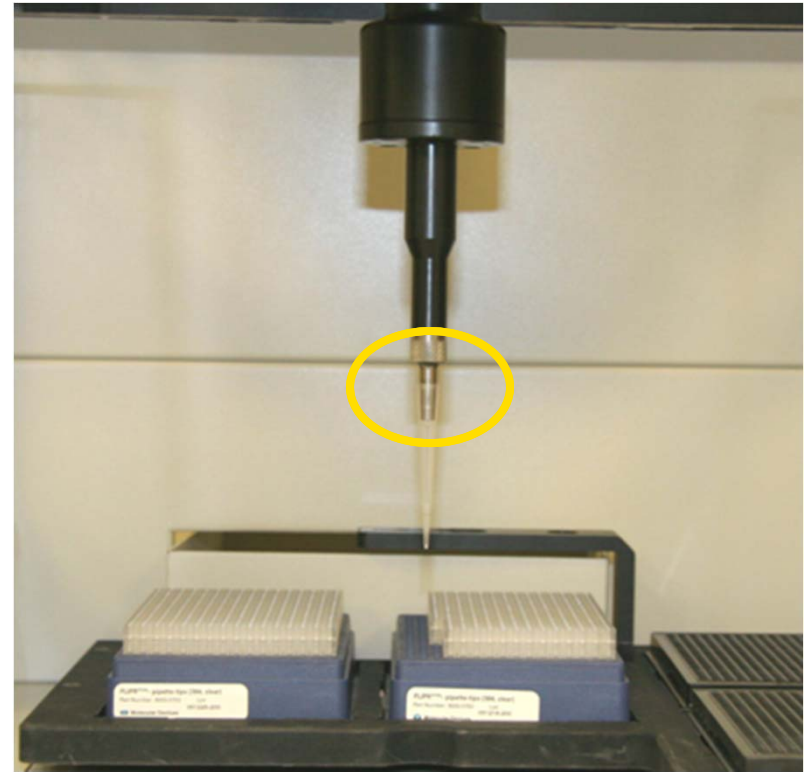
- 3-25 ml
- Tip Part **9000-0763** FLIPR Tetra Pipette tips, clear, non-sterile, 384-well 50 racks/case
- Replacement Adapter Part **2300-2280** ImageXpress Micro Fluidics Tip Adapter, 384-well



Changing Fluidics Tip Types: Summary

There are 5 basic steps for changing tip types in the ImageXpress Micro Fluidics option:

1. Hardware: change the tip adapter
2. Hardware: load the appropriate tips racks on the system
3. Software: select the correct tip adapter
4. Software: select the correct tip type
5. Software: reset the tips

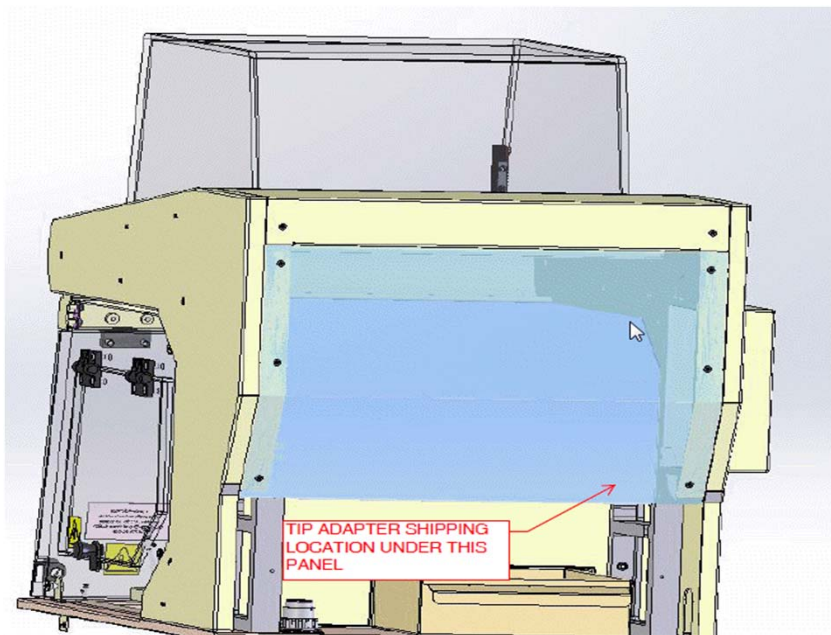


Change Tip Adapter

1. Change the tip adapter

- Locate tip adapter on the back of the instrument or in the IXM accessory kit
- Open access door and remove previous tip adapter (untwist with fingers)
- Screw the new tip adapter on until finger-tight
- Store the previous tip adapter

NOTE Do not change while an experiment is running

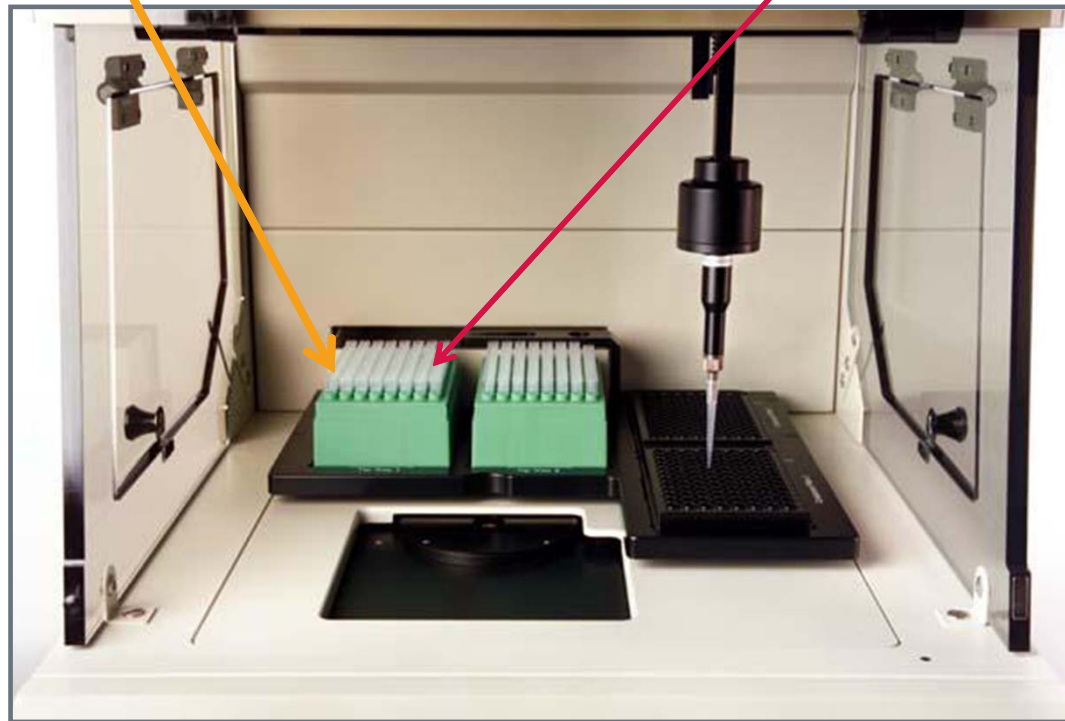


Load Appropriate Tips

2. Load instrument deck with tip racks
 - Remove any existing tip racks
 - Insert new tip type rack [insert **right side** first, then push down on **left side**]
 - Remove tip rack lid

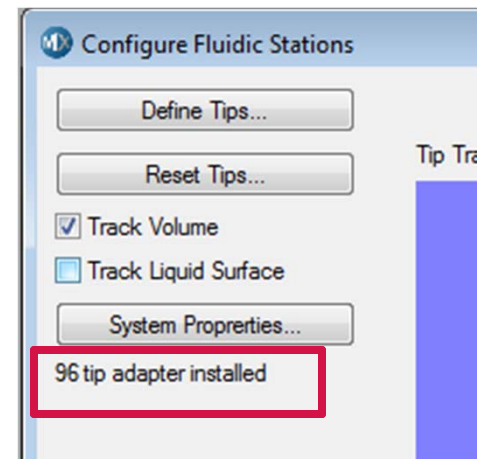
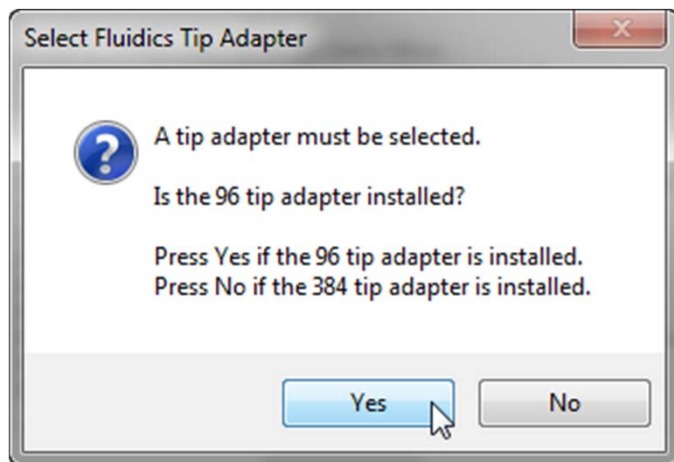
2. Push down on left side

1. Insert right side



Select Tip Adapter in Software – Ask Mode

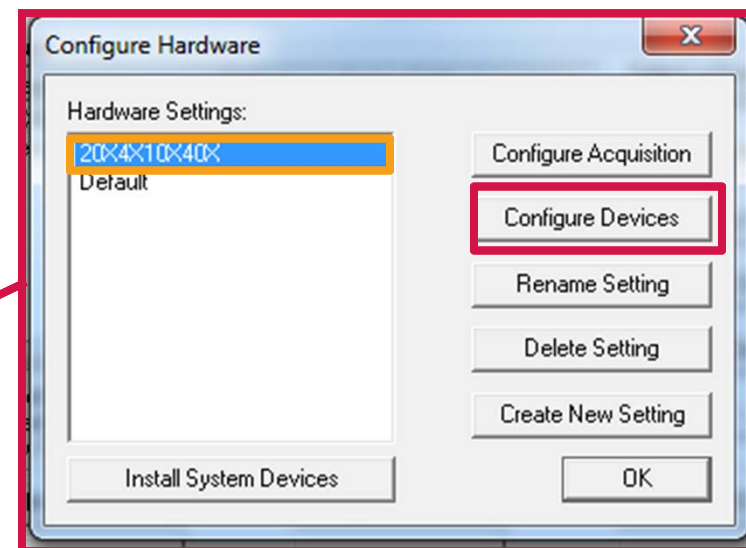
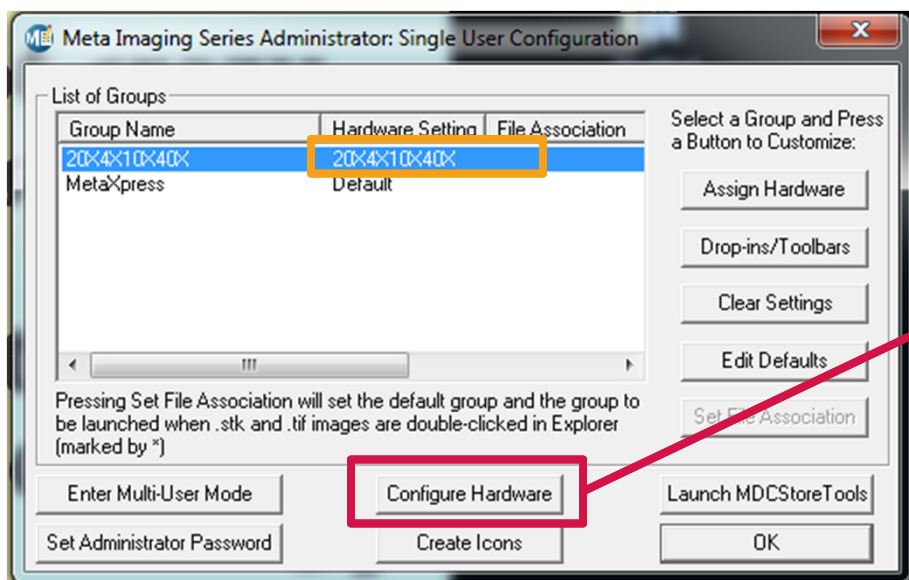
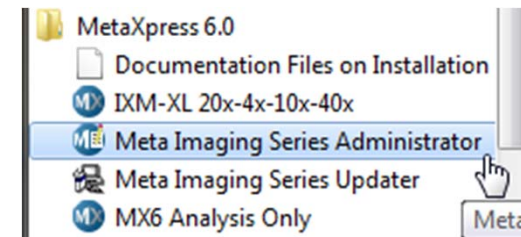
3. If Ask Mode is activated in **MetaXpress**, you will be asked to verify which tip adapter is installed when starting the software
 - Select the appropriate option for the installed tip adapter
 - Verify the current tip adapter selection in the **Plate Acquisition Setup** dialog (**Configure > Fluidics > Configure Stations**)



Select Tip Adapter in Software – Not In Ask Mode

3. If Ask Mode is deactivated in MetaXpress, you will need to exit the software and change the setting in **Meta Imaging Series Administrator**

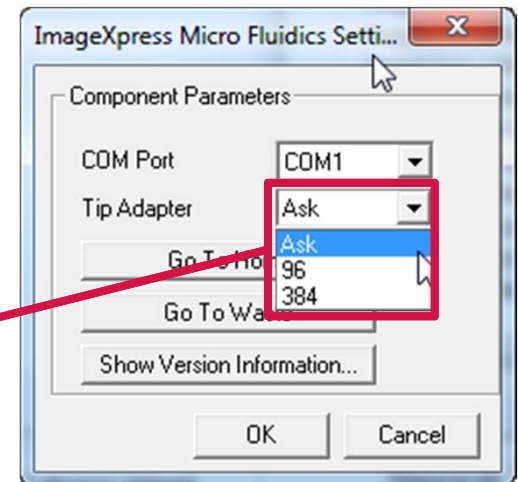
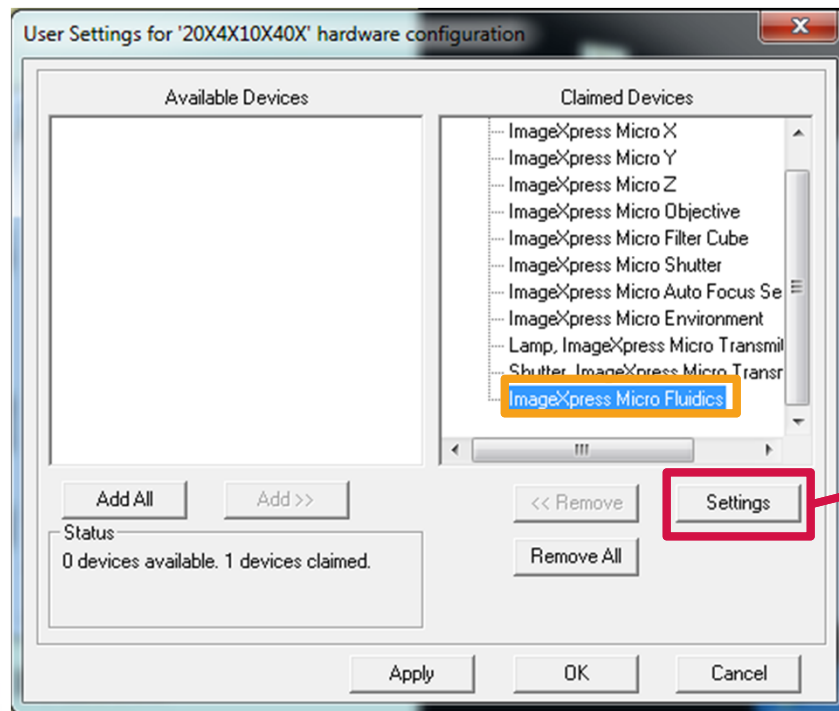
- Open **Meta Imaging Series Administrator**
- Identify the correct hardware setting and click on **Configure Hardware**
- Highlight the correct hardware setting in the **Configure Hardware** dialog, then click **Configure Devices**



Select Tip Adapter in Software – Not In Ask Mode

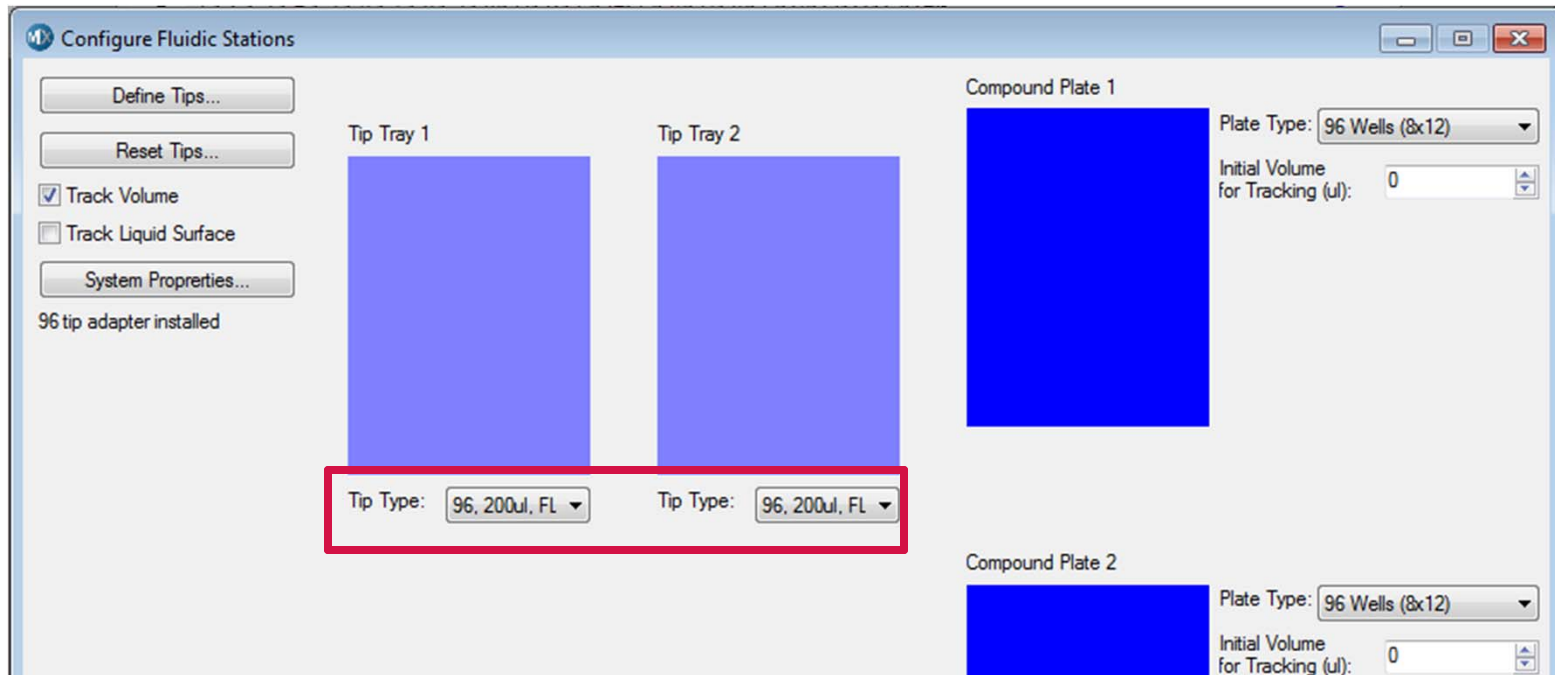
3. In **Meta Imaging Series Administrator** (cont.)
 - Highlight **ImageXpress Micro Fluidics** under the **Claimed Devices** section and click **Settings**
 - Select the appropriate option from the **Tip Adapter** drop-down menu
 - Click **OK** until you have exited **Meta Imaging Series Administrator**

NOTE By selecting **Ask** in the ImageXpress Micro Fluidics Settings dialog, you will activate Ask Mode for verifying tip adaptor selection every time MetaXpress is started.



Select Correct Tip Type in Software

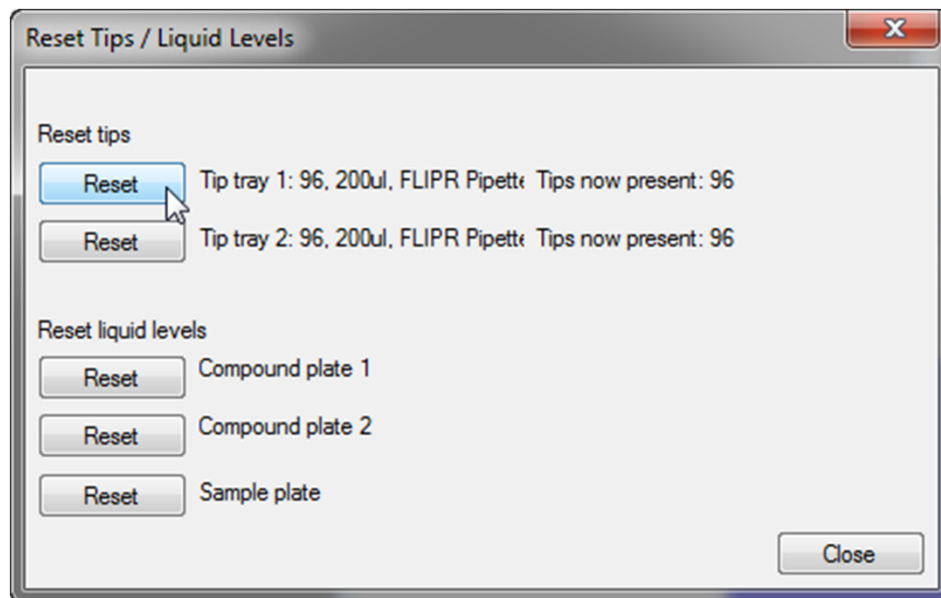
4. Select the correct tip type in **Plate Acquisition Setup**
 - Sign back into **MetaXpress** if necessary
 - Open the **Plate Acquisition Setup** dialog
 - **Configure > Fluidics > Configure Stations**
 - Change the **Tip Type** from the drop-down menus



Reset Tips in Software

5. Reset the tips

- Verify that tip rack(s) on system deck are full
- In the **Plate Acquisition Setup** dialog, select the **Fluidics** tab, click on the **Reset Tips** button
- Reset the appropriate tip track locations and click **Close**



Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- User Forum: <http://metamorph.moleculardevices.com/forum/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments





ADVANCING PROTEIN AND CELL BIOLOGY