

# ClonePix 2 Training Guide

## Picking Head Maintenance



Date Revised 2/27/2017 Version A

# Index

- [Index](#)
- [Chapter Purpose](#)
- [Removing the Picking Head](#)
- [Removing Picking Pins from Picking Head](#)
- [Picking Pin Cleaning Procedure](#)
- [Picking Head Cleaning Procedure](#)
- [Replacing Picking Pins in Picking Head](#)
- [Replacing the Picking Head](#)
- [Pin Fire Test](#)
- [Select Source Plate Type](#)
- [Load Source Plate Type](#)
- [Adjusting Well Alignment](#)
- [Adjusting Primary Picking Pin Z Datum Position](#)
- [Confirm Plate Type](#)
- [Adjusting Primary Picking Pin XY Alignment](#)
- [Adjusting Picking Pins #2-8 XY Alignment](#)
- [Filling the Ethanol Wash Bath](#)
- [Sanitizing Picking Pins](#)
- [UV Sanitizing ClonePix 2 Deck](#)
- [Check Wash & Waste Bottles](#)
- [Rinsing Picking Pins](#)
- [Conclusion of Process](#)
- [Support Resources](#)

# Chapter Purpose

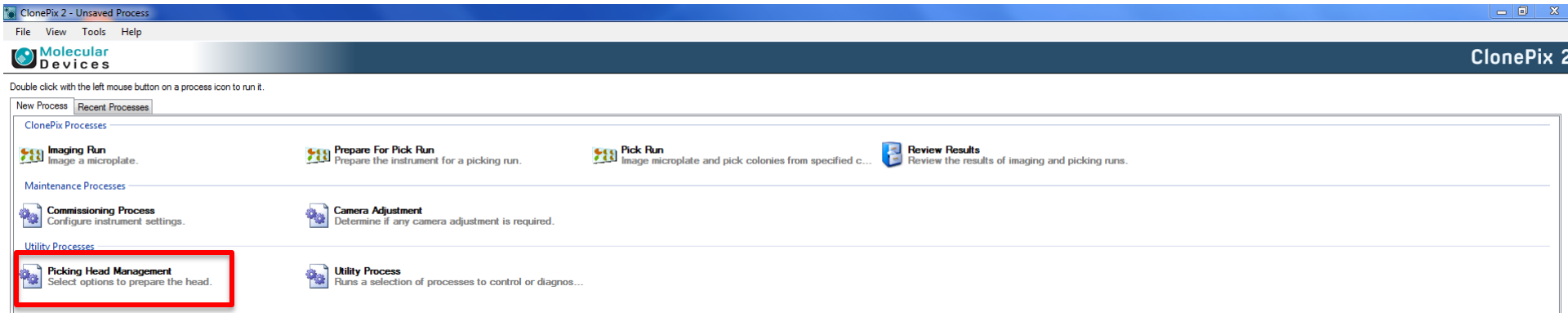
The purpose of this chapter is to guide the user through the process for **maintaining the picking head** on the **ClonePix 2**.

The Prepare for Pick Run and Pick Run setup processes are not covered in this document. Please consult the dedicated training modules for detailed information on these topics.



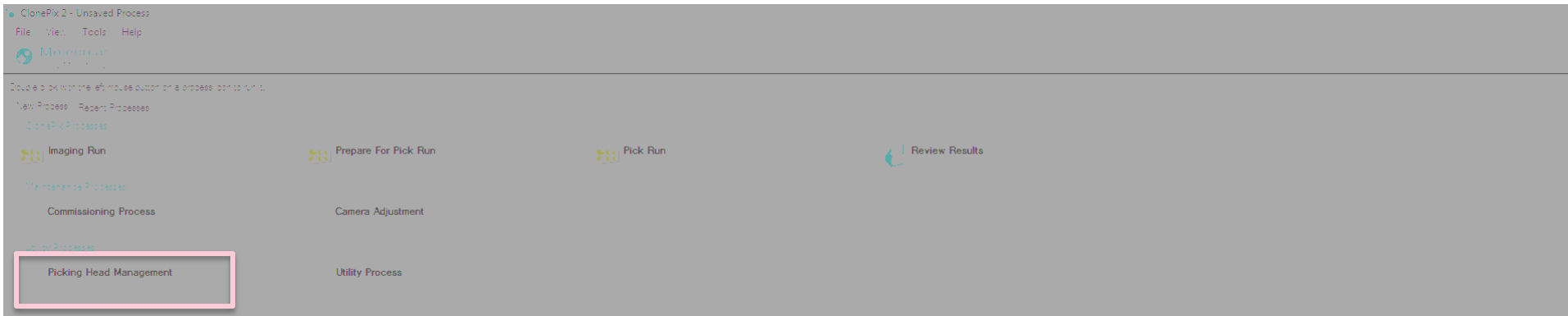
# Removing the Picking Head

1. From the **Main Menu**, under **Utility Processes** select **Picking Head Management**.

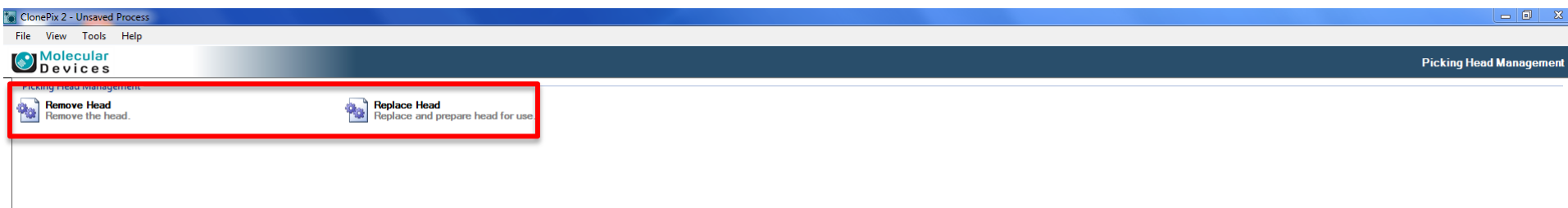


# Removing the Picking Head

1. From the **Main Menu**, under **Utility Processes** select **Picking Head Management**.

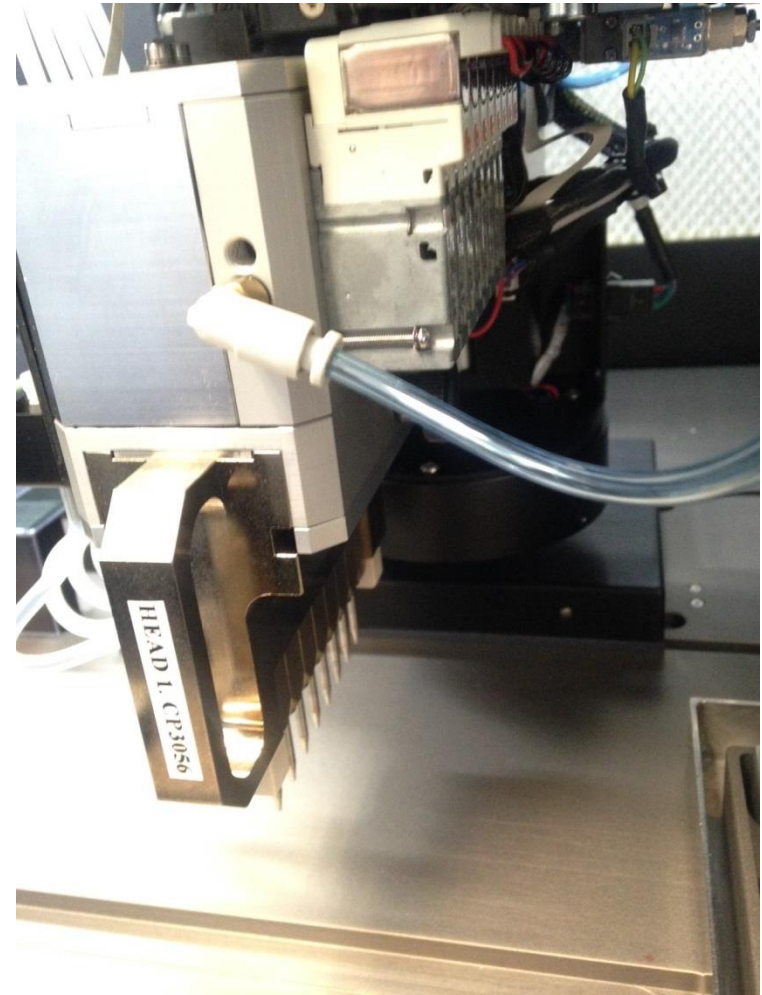


2. From the **Picking Head Management** menu:
  - a. If you **are not** planning to replace with an alternate picking head immediately: select **Remove Head**.
  - b. If you **are** planning to replace the head with an alternate picking head: select **Replace Head**.



# Removing the Picking Head

3. The ClonePix 2 will now home the **picking head** to a **safe position** for removal – it will be moved to the **front right** of the **instrument deck**.

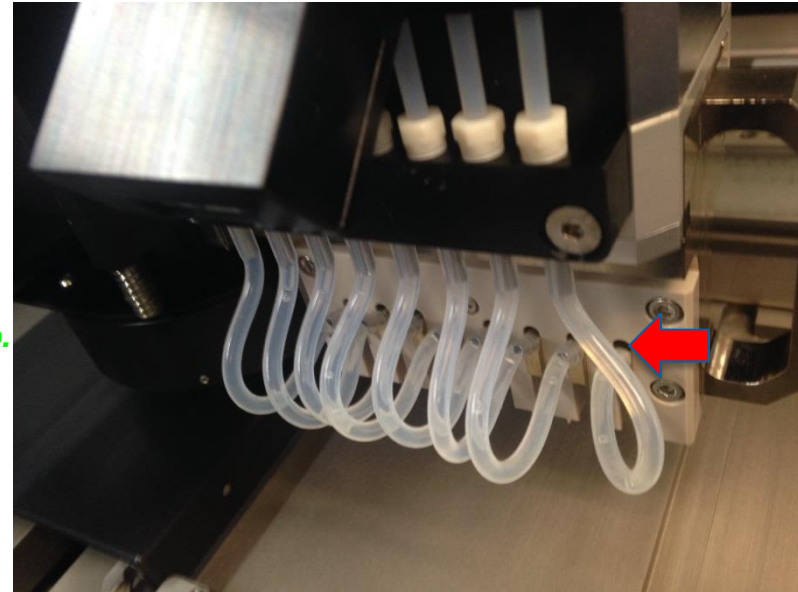


# Removing the Picking Head

3. The ClonePix 2 will now home the **picking head** to a **safe position** for removal – it will be moved to the **front right** of the **instrument deck**.
  - a. Using gloved hands, **detach** the **tubing** from **each pin** of the **picking head**.

**NOTE:** *If you wish to clean the tubing, you can also detach from the head carrier.*

**TIP:** *Place a sterile plate lid beneath the head to catch any liquid that may be released by the head or tubing at this step.*



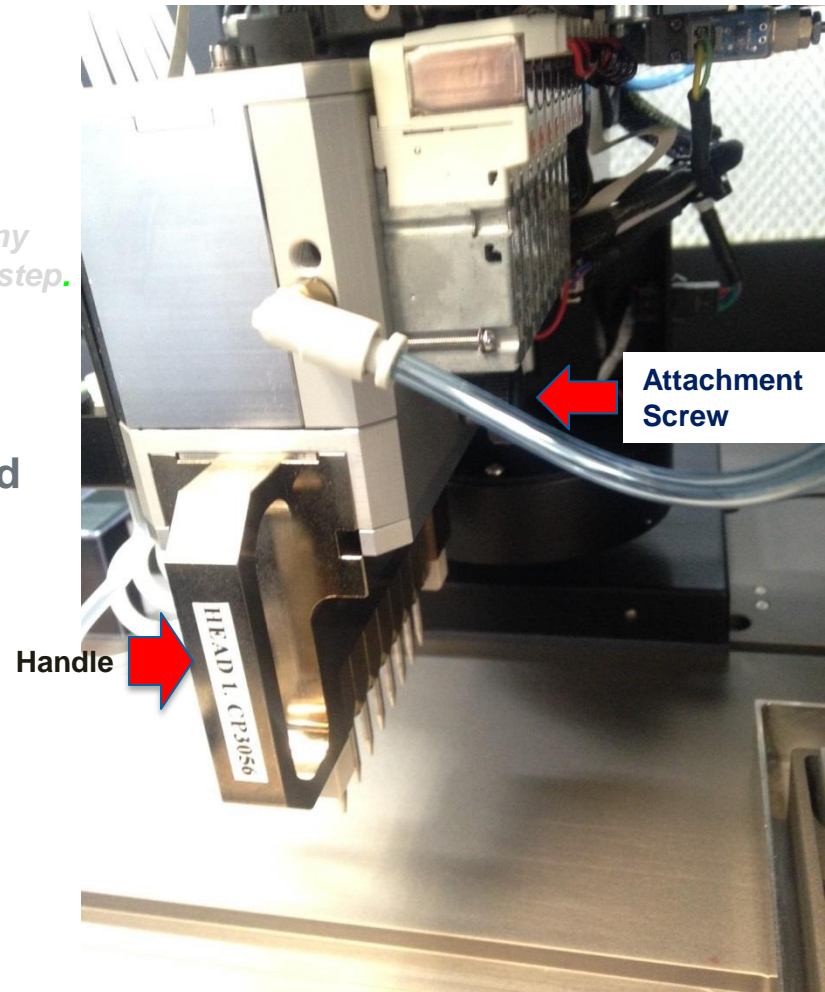
# Removing the Picking Head

- The ClonePix 2 will now home the **picking head** to a **safe position** for removal – it will be moved to the **front right** of the **instrument deck**.
  - Using gloved hands, **detach** the **tubing** from **each pin** of the **picking head**.

*NOTE: If you wish to clean the tubing, you can also detach from the head carrier.*

*TIP: Place a sterile plate lid beneath the head to catch any liquid that may be released by the head or tubing at this step.*

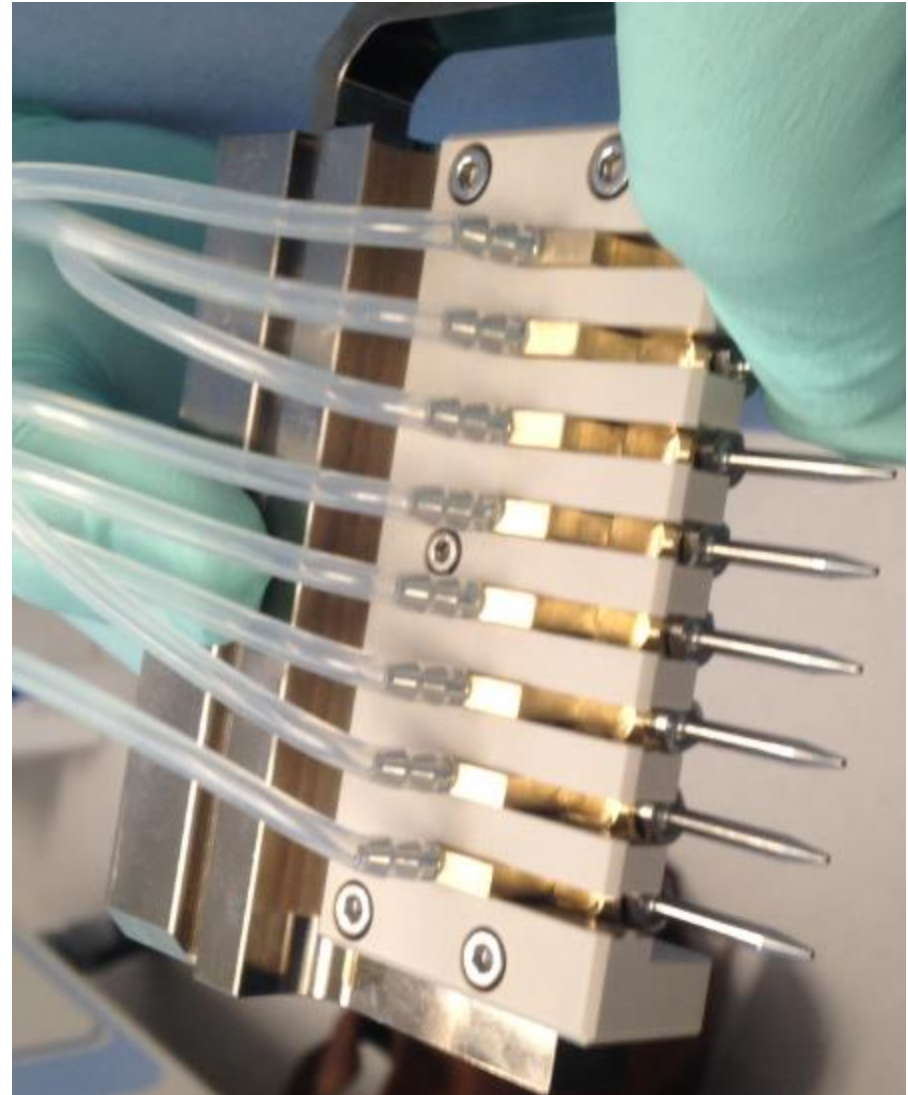
- While holding the **picking head** firmly by the **handle** at the front, locate and **loosen** the **black attachment screw**. Slide the **picking head** out of the **carrier**, taking care to support the head (it is heavy!).





# Removing Picking Pins from Picking Head

1. Once the picking head is removed, you **MUST** remove all picking pins from the head for cleaning and sterilization as follows:



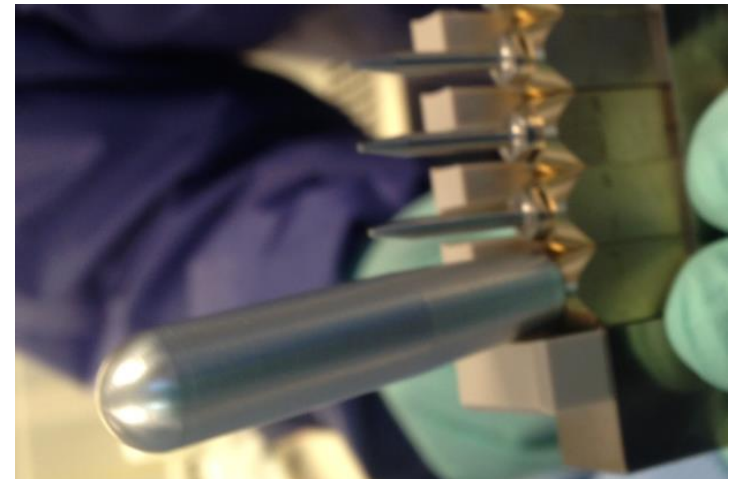
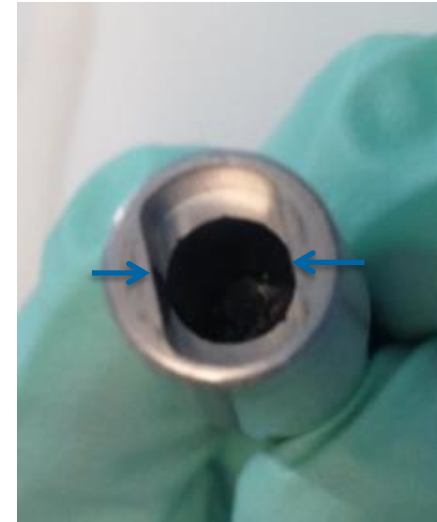
# Removing Picking Pins from Picking Head

1. Once the **picking head** is removed, you must **remove all picking pins** from the head for **cleaning and sterilization** as follows:
  - a. Locate the **Pin Removal Key** provided with your instrument.



# Removing Picking Pins from Picking Head

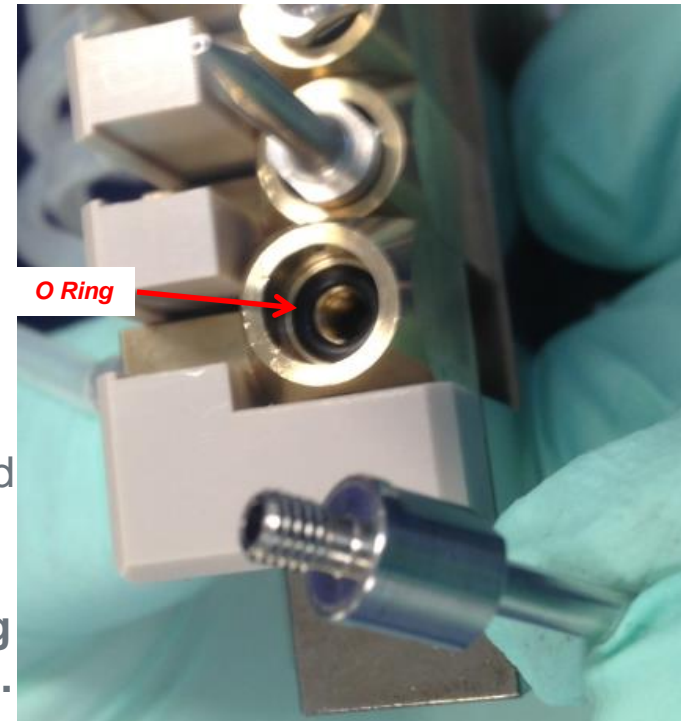
1. Once the **picking head** is removed, you must **remove all picking pins** from the head for **cleaning and sterilization** as follows:
  - a. Locate the **Pin Removal Key** provided with your instrument.
  - b. To remove **picking pins** from the **head**, slide the key over each pin, taking care to **align the flat edges** of the **pin removal key opening** with the corresponding edges at the **base** of the **pin**.



# Removing Picking Pins from Picking Head

1. Once the picking head is removed, you must remove all picking pins from the head for cleaning and sterilization as follows:
  - a. Locate the **Pin Removal Key** provided with your instrument.
  - b. To remove **picking pins** from the **head**, slide the key over each pin, taking care to **align the flat edges** of the pin removal key opening with the corresponding edges at the **base** of the **pin**.
  - c. While firmly holding the **pin removal key**, **twist counter-clockwise** to unscrew the pin from the head. Once the pin is loose, you can **remove by hand**.

**NOTE:** You will see a small black rubber “O” ring either still inside the head OR attached to the pin. **DO NOT autoclave the O ring** – this will compromise the integrity of the seal.



# Picking Pin Cleaning Procedure

1. **Picking Pins** should be cleaned by **sonication** in **aQuClean (K2505)** and **autoclaved** prior to use.
  - a. Prepare **50 mL** of a **2% solution** of **aQuClean** in **di-ionized water**.

A **sample** of **aQuClean** is provided with your instrument. To **order more**, please follow the **link** below and click on **Request Quote** or **contact your local sales representative** :  
<https://www.moleculardevices.com/reagents-supplies/cell-biology-reagents/cleaning-and-sterilizing-solutions>
  - b. Immerse **picking pins** in **aQu Clean** and **sonicate** for **10 minutes**.
  - c. **Rinse pins** thoroughly in **de-ionized water**.
  - d. Place **pins** in **de-ionized water** and **sonicate** for **10 minutes**.
  - e. **Autoclave pins** and **pin removal key**.



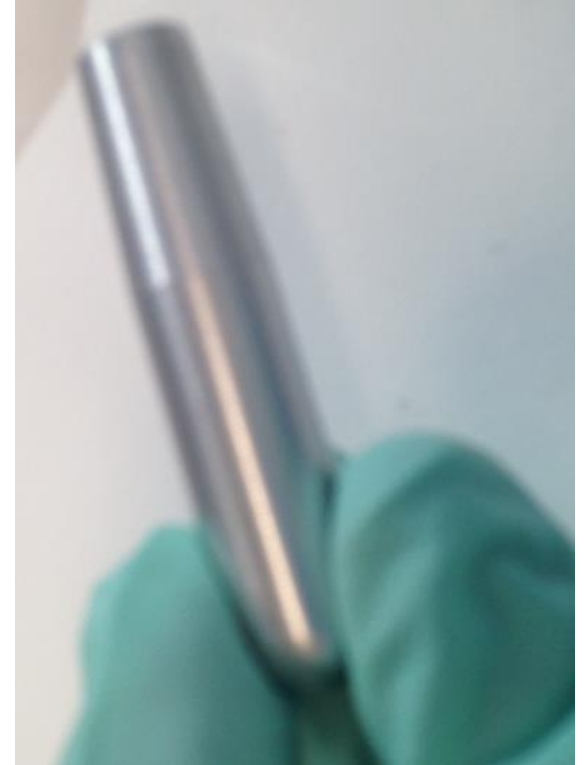
# Picking Head Cleaning Procedure

1. The Picking Head & O rings can be cleaned using the following procedure:
  - a. Soak the **picking head/O rings** in **100% Ethanol**.
  - b. If you suspect that the **head is clogged**, you may **flush** the head using a **sterile 5 mL syringe** filled with **100% Ethanol**.
  - c. Allow **head to completely dry** in a **sterile tissue culture hood** (several hours to overnight).
  - d. **DO NOT autoclave the picking head or O rings** – this will cause **damage** to the head and O rings and affect picking capability.



# Replacing Picking Pins in Picking Head

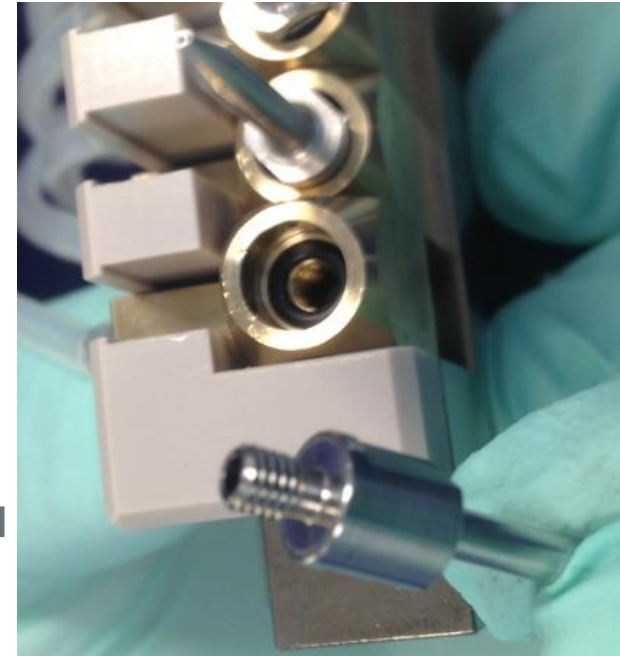
1. Once the picking head is cleaned and the pins autoclaved, you can replace the pins in the picking head as follows:
  - a. Autoclave the **Pin Removal Key** provided with your instrument.



# Replacing Picking Pins in Picking Head

1. Once the picking head is cleaned and the pins autoclaved, you can replace the pins in the picking head as follows:
  - a. Autoclave the **Pin Removal Key** provided with your instrument.
  - b. To replace **picking pins** in the **head**, in a sterile tissue culture hood with gloved hands, insert each pin into the head by holding the base of the pin and twisting clockwise.

NOTE: At this point you should also inspect the **rubber O ring** in each position and ensure that it is **present and intact before** inserting the pin. If the **O ring** is **damaged** or **missing**, you can **replace** with a **new ring** (contact your regional sales rep to order more if needed – part # X1036).





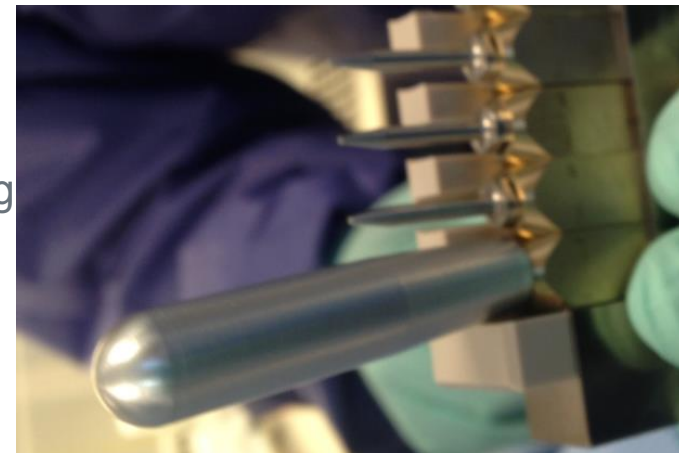
# Replacing Picking Pins in Picking Head

1. Once the picking head is cleaned and the pins autoclaved, you can replace the pins in the picking head as follows:

- a. Autoclave the **Pin Removal Key** provided with your instrument.
- b. To replace **picking pins** in the **head**, in a sterile tissue culture hood with gloved hands, insert each pin into the head by holding the base of the pin and twisting clockwise.

NOTE: At this point you should also inspect the **rubber O ring** in each position and ensure that it is **present and intact before** inserting the pin. If the **O ring** is **damaged or missing**, you can **replace** with a **new ring** (contact your regional sales rep to order more if needed – part X1036).

- c. Next, place the **pin removal key** over each **pin**, taking care to **align the flat edges of the key opening** with the corresponding edges at the **base** of the **pin**. **Twist the key** firmly in the **clockwise direction** until the **pin** is seated **finger-tight**.



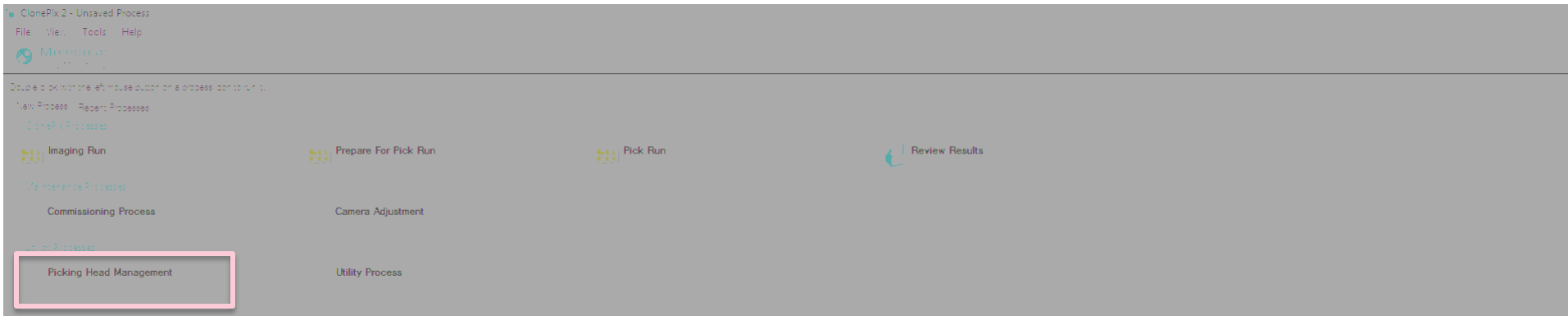
# Replacing the Picking Head

1. From the **Main Menu**, under **Utility Processes** select **Picking Head Management**.

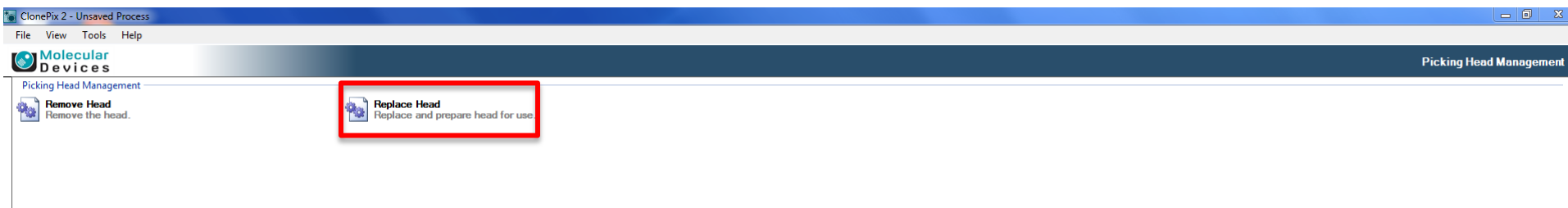
The screenshot shows the ClonePix 2 software interface. The window title is "ClonePix 2 - Unsaved Process". The menu bar includes "File", "View", "Tools", and "Help". The Molecular Devices logo is in the top left, and "ClonePix 2" is in the top right. Below the menu bar, there is a instruction: "Double click with the left mouse button on a process icon to run it." There are two tabs: "New Process" and "Recent Processes". The main area is divided into sections: "ClonePix Processes", "Maintenance Processes", and "Utility Processes". Under "ClonePix Processes", there are four icons: "Imaging Run" (Image a microplate.), "Prepare For Pick Run" (Prepare the instrument for a picking run.), "Pick Run" (Image microplate and pick colonies from specified c...), and "Review Results" (Review the results of imaging and picking runs.). Under "Maintenance Processes", there are two icons: "Commissioning Process" (Configure instrument settings.) and "Camera Adjustment" (Determine if any camera adjustment is required.). Under "Utility Processes", there are two icons: "Picking Head Management" (Select options to prepare the head.) and "Utility Process" (Runs a selection of processes to control or diagnos...). The "Picking Head Management" icon is highlighted with a red rectangular box.

# Replacing the Picking Head

1. From the **Main Menu**, under **Utility Processes** select **Picking Head Management**.

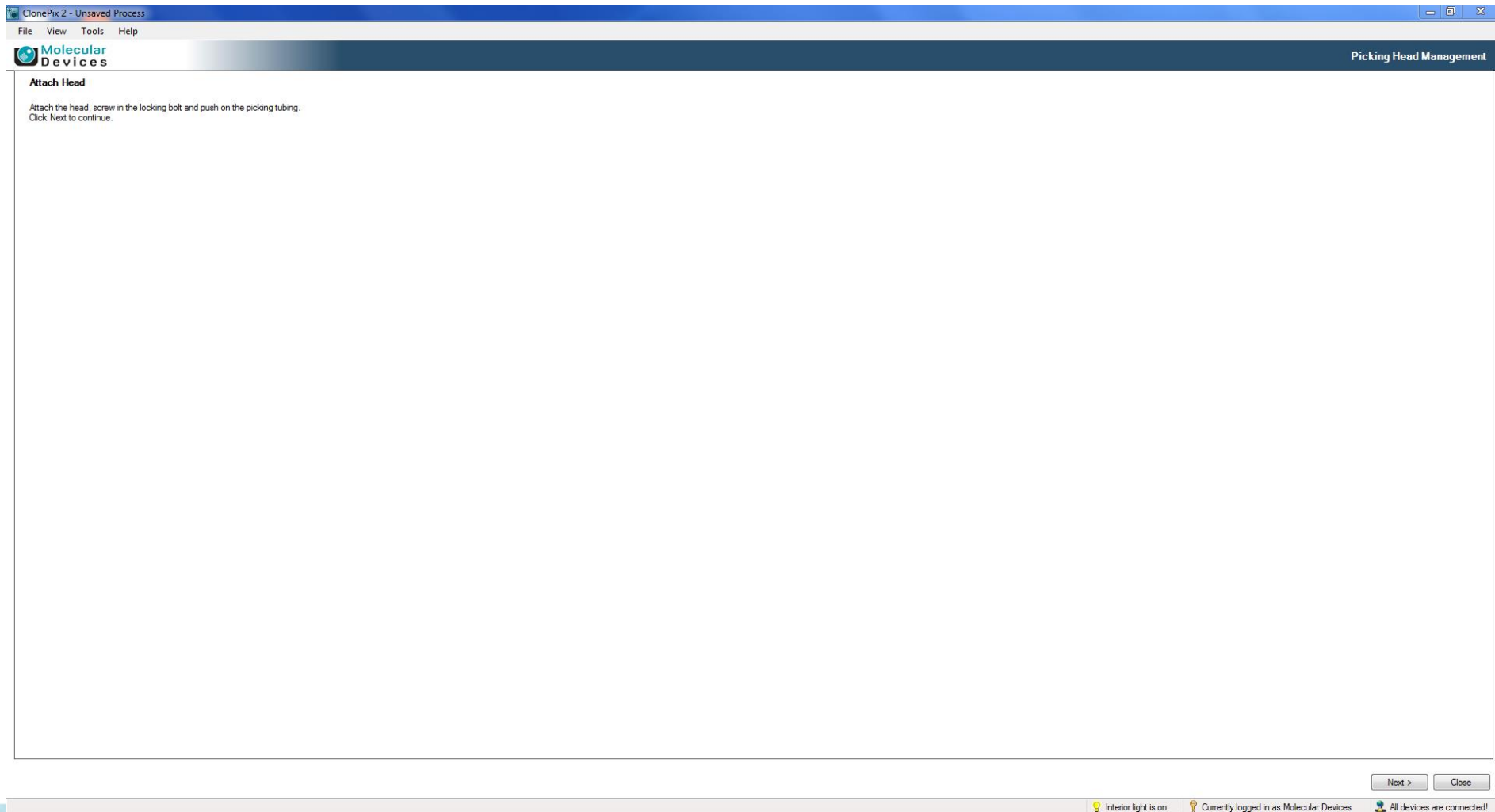


2. From the **Picking Head Management** menu, select **Replace Head**.



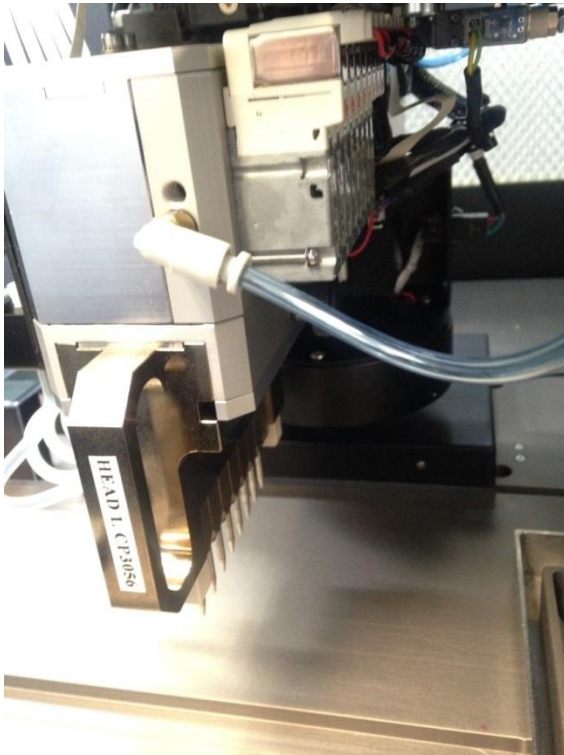
# Replacing the Picking Head

3. The **Remove Head** dialog appears – click **Next** to continue. Then the **Attach Head** dialog will appear.



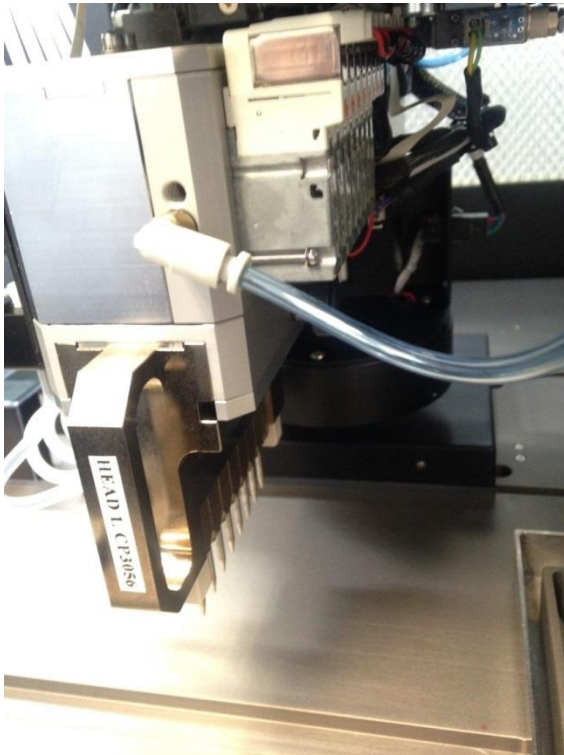
# Replacing the Picking Head

4. The picking head carrier (actuator) will now be in a position where you can easily insert the head.
  - a) Using two gloved hands, carefully slide the picking head into the actuator.



# Replacing the Picking Head

4. The picking head carrier (actuator) will now be in a position where you can easily insert the head.
  - a) Using two gloved hands, carefully slide the picking head into the actuator.
  - b) Reattach the feed tubing to each picking pin, starting from back to front as you are facing the head. Ensure that tubing is fully attached to each pin before moving on to the next one.



# Replacing the Picking Head

## 5. Click Next.

The screenshot shows a software window titled "ClonePix 2 - Unsaved Process" with a menu bar (File, View, Tools, Help) and the Molecular Devices logo. The main window is titled "Picking Head Management" and contains the following text:

**Attach Head**

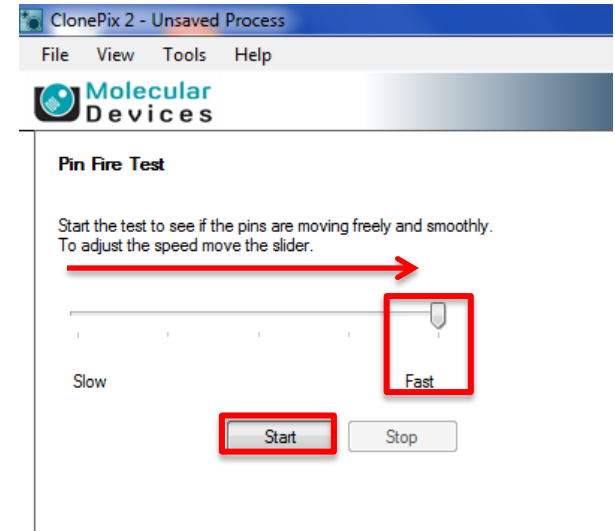
Attach the head, screw in the locking bolt and push on the picking tubing.  
Click Next to continue.

At the bottom right of the window, there are two buttons: "Next >" (highlighted with a red box) and "Close".

At the bottom of the screen, there is a status bar with the following text: "Interior light is on.", "Currently logged in as Molecular Devices", and "All devices are connected!".

# Pin Fire Test

6. The **Pin Fire Test** dialog now appears.
  - a) **Click and drag the speed slider** to select the **Fast** setting, then click **Start**.





# Pin Fire Test

6. The **Pin Fire Test** dialog now appears.
  - a) **Click and drag the speed slider** to select the **Fast** setting, then click **Start**.
  - b) You will now see that each **picking pin** in the **head** is fired (**extended**), then **retracted sequentially**.

Watch carefully for **3-4 full cycles** and ensure that **each of the 8 picking pins** is moving **smoothly**. If you note any **sticking** or **significant hesitation** of the **pin** movement, contact **Molecular Devices Technical support** for further assistance (**800-635-5577** or [support@moldev.com](mailto:support@moldev.com))

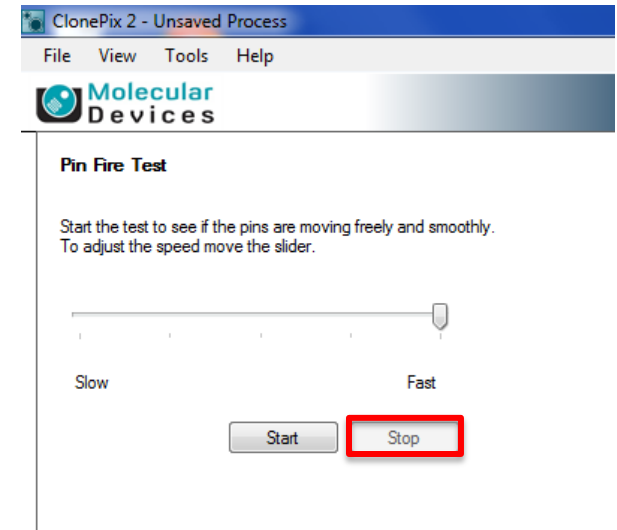


# Pin Fire Test

6. The **Pin Fire Test** dialog now appears.
  - a) Click and drag the **speed slider** to select the **Fast** setting, then click **Start**.
  - b) You will now see that each **picking pin** in the **head** is fired (**extended**), then **retracted sequentially**.

Watch carefully for **3-4 full cycles** and ensure that **each of the 8 picking pins** is moving smoothly. If you note any **sticking or significant hesitation of the pin movement**, contact **Molecular Devices Technical support** for further assistance (800-635-5577 or [support@moldev.com](mailto:support@moldev.com))

- c) Click the **Stop** button to conclude the test, then click **Next** to proceed.



# Select Source Plate Type

7. You will now see a screen where you are prompted to **select your source microplate type**. **Select your source microplate type** from the **dropdown list**, then click **Next**.

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Picking Head Management

Select the microplate type that you wish to use from the list:

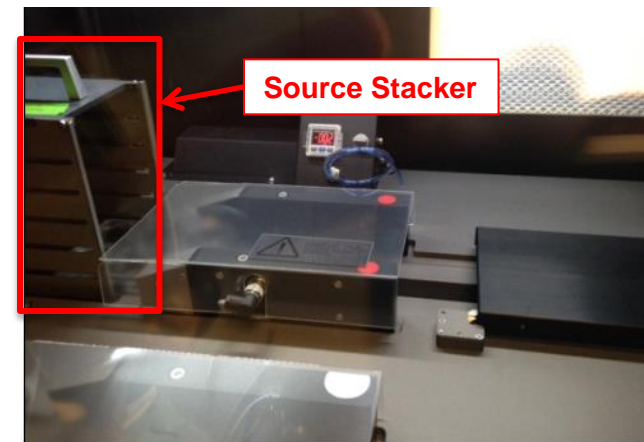
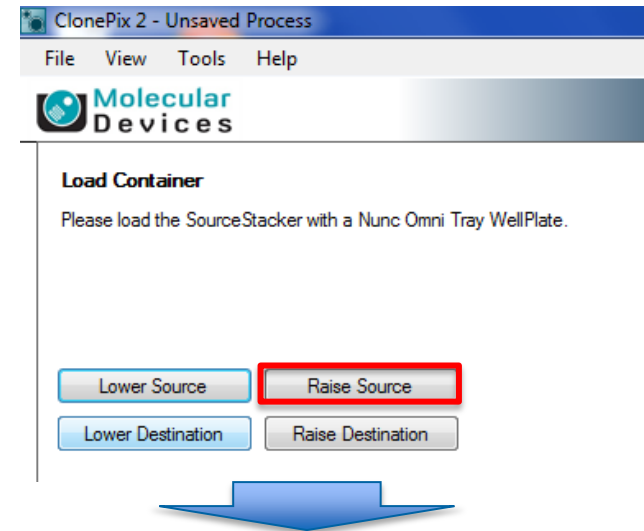
Nunc Omni Tray

Next > Close

Interior light is on. Currently logged in as Molecular Devices All devices are connected!

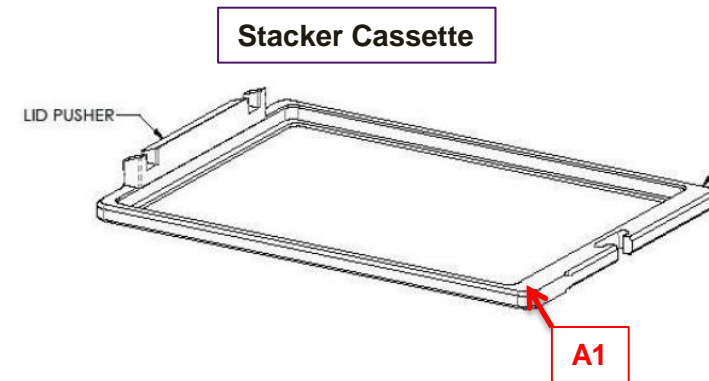
# Load Source Plate Type

8. The Load Container dialog appears, where you are prompted to **load the source stacker** with your selected **source plate type**.
  - a) Select **Raise Source** - the **Source Stacker** will now be **raised** within the **ClonePix 2** instrument.



# Load Source Plate Type

8. The **Load Container** dialog appears, where you are prompted to **load the source stacker** with your selected **source plate type**.
  - a) Select **Raise Source** - the **Source Stacker** will now be **raised** within the **ClonePix 2** instrument.
  - b) Place an **empty** (no media, no cells) **source plate** in the **appropriate stacker cassette** (see chart for details), with well **A1** facing the **lower right hand corner** (away from the color-coded lid pusher).



Source Plate	Lid Pusher Color	Lid Lifter Assembly
Greiner 6 well	Green	Red
Nunc Omni -	Blue	Silver
Nunc - 6 Well	Green	White

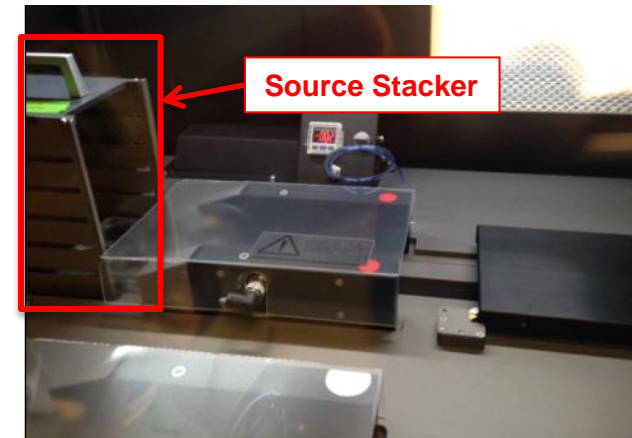


# Load Source Plate Type

Source Plate + Stacker Cassette



Lid Pusher



Source Stacker

8. The **Load Container** dialog appears, where you are prompted to **load the source stacker** with your selected **source plate type**.
  - a) Select **Raise Source** - the **Source Stacker** will now be **raised** within the **ClonePix 2** instrument.
  - b) Place an **empty** (no media, no cells) **source plate** in the **appropriate stacker cassette** (see chart for details), with well A1 facing the lower right hand corner (away from the color-coded lid pusher).
  - c) Insert your **source plate + stacker cassette** into **any position** within the **source stacker** with the **lid pusher** facing the **back** of the **stacker** (until you hear the **magnet engage**), then click **Next**.

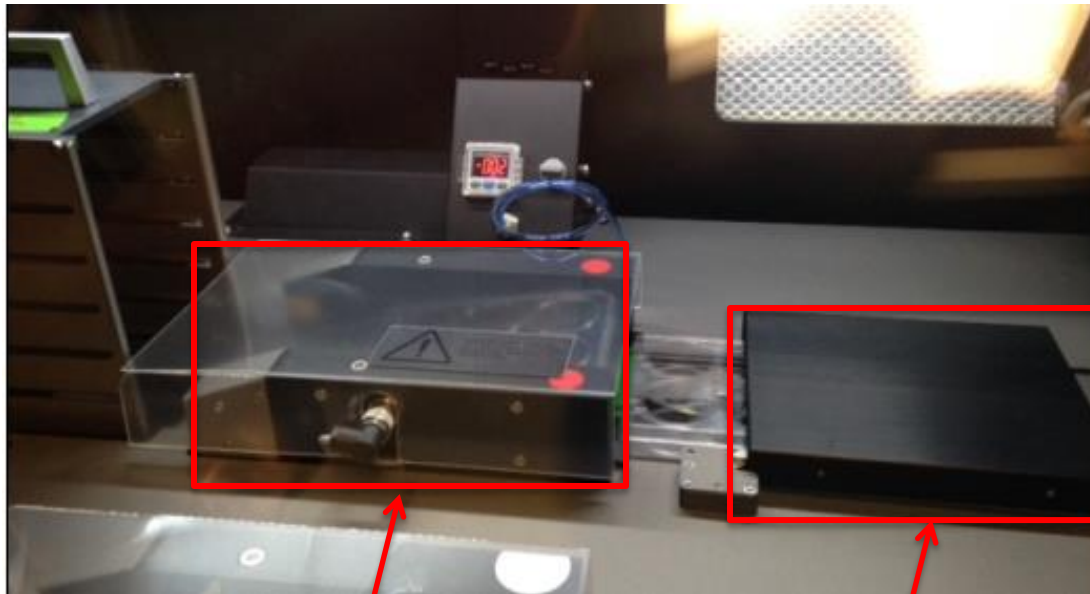
# Load Source Plate Type

8. The **Load Container** dialog appears, where you are prompted to **load the source stacker** with your selected **source plate type**.
  - a) Select **Raise Source** - the **Source Stacker** will now be **raised** within the **ClonePix 2** instrument.
  - b) Place an **empty** (no media, no cells) **source plate** in the **appropriate stacker cassette** (see chart for details), with well **A1** facing the **lower right hand corner** (away from the color-coded lid pusher).
  - c) Insert your **source plate + stacker cassette** into **any position** within the **source stacker** with the **lid pusher** facing the **back** of the **stacker** (until you hear the **magnet** engage), then click **Next**.
  - d) The **stacker** will now **automatically** move through the stack and **detect the position of your loaded plate** (this could take up to a minute). You will see the **Loading holders** indicator on your computer screen during this process.



# Load Source Plate Type

8. The **Load Container** dialog appears, where you are prompted to **load the source stacker** with your selected **source plate type**.
  - e) The **loaded plate** will next be pulled from the **stacker** into the **de-lidding station**, then to the **imaging station** for the next step. You will see the **Loading microplate** indicator on your computer screen during this process.



**De-Lidding  
Station**

**Imaging  
Station**

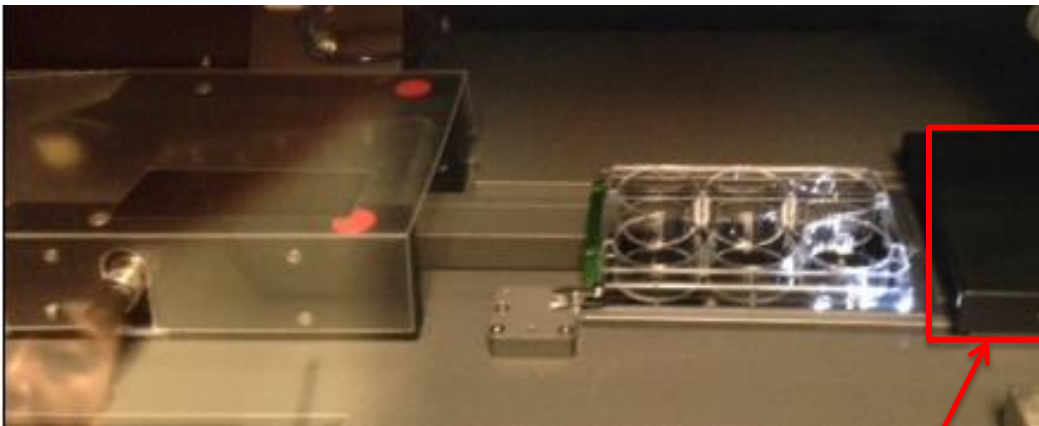




# Adjusting Well Alignment

9. The **ClonePix 2 system** will now capture **white light images** of a **well** of your **source plate** and the software will enable any necessary **adjustment to alignment of the well** as follows:

- a) The **illumination cover** will automatically **retract** and the system will capture **4 images** of your **well edges** (north, south, east, and west sides). You will see the **Acquiring Alignment Images** indicator appear on your computer screen during this process.



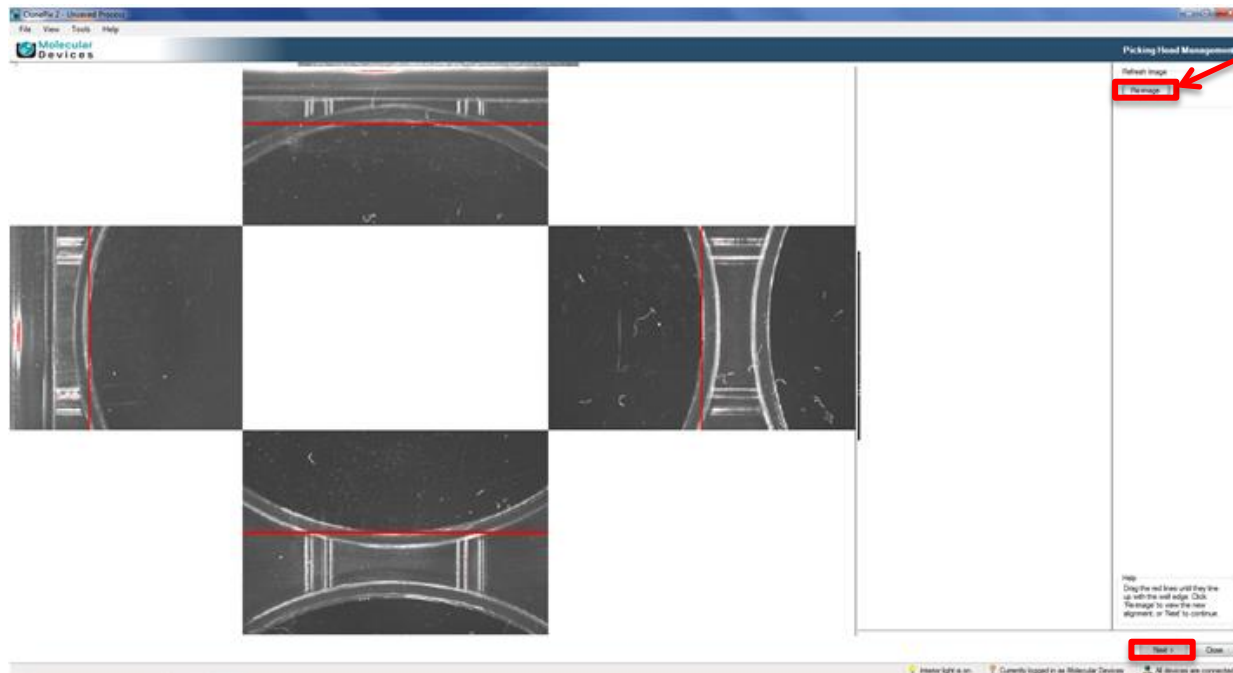
**Illumination  
Cover**



# Adjusting Well Alignment

9. The **ClonePix 2** system will now capture **white light images** of a well of your **source plate** and the software will enable any necessary **adjustment to alignment of the well** as follows:

- The **illumination cover** will automatically **retract** and the system will capture **4 images** of your **well edges** (north, south, east, and west sides). You will see the **Acquiring Alignment Images** indicator appear on your computer screen during this process.
- You will now see the **4 well edge images** appear on your screen. The **red lines** superimposed on the image set should **align** closely with the **well edges**. **Adjust** the line positioning by **clicking and dragging** – note that the top/bottom (north/south) lines will move together, same for left/right (east/west). If you make an adjustment, click the **Re-image** button to confirm alignment. Click **Next** to proceed.

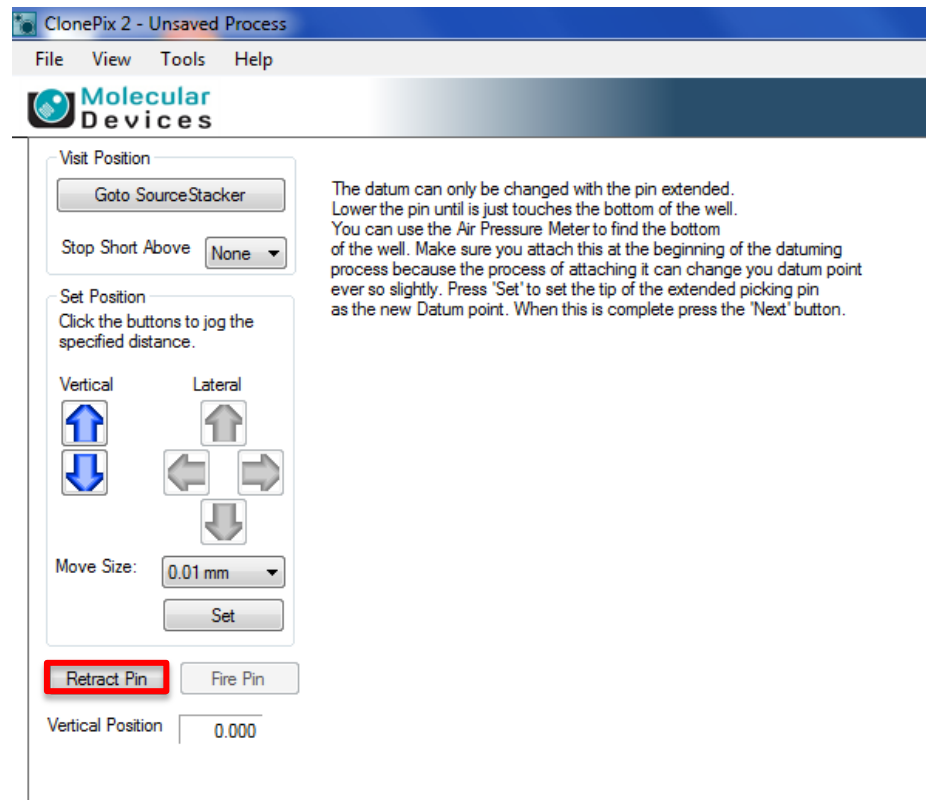


Re-image  
Button

# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

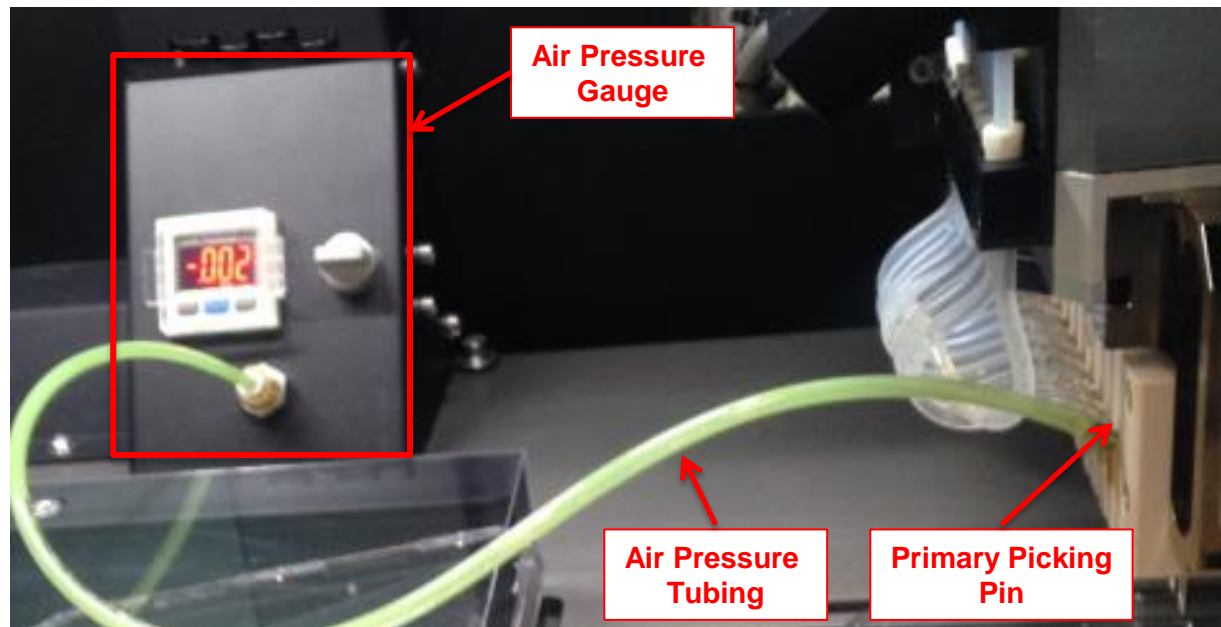
- The **illumination cover** will remain **retracted** and the **primary picking pin** (#1, closest to you as you face the instrument) will be **extended** to the center of a well in your loaded source plate. You will see the **adjustment dialog** appear (screenshot below) – click on **Retract Pin** before proceeding to the next step.



# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

- a) The **illumination cover** will remain **retracted** and the **primary picking pin** (#1, closest to you as you face the instrument) will be **extended** to the center of a well in your loaded source plate. You will see the **adjustment dialog** appear (screenshot below) – click on **Retract Pin** before proceeding to the next step.
- b) **Disconnect the picking tubing** from the **primary picking pin**, then connect the **free end** of the **air pressure tubing** to the **primary picking pin** as shown below.



# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

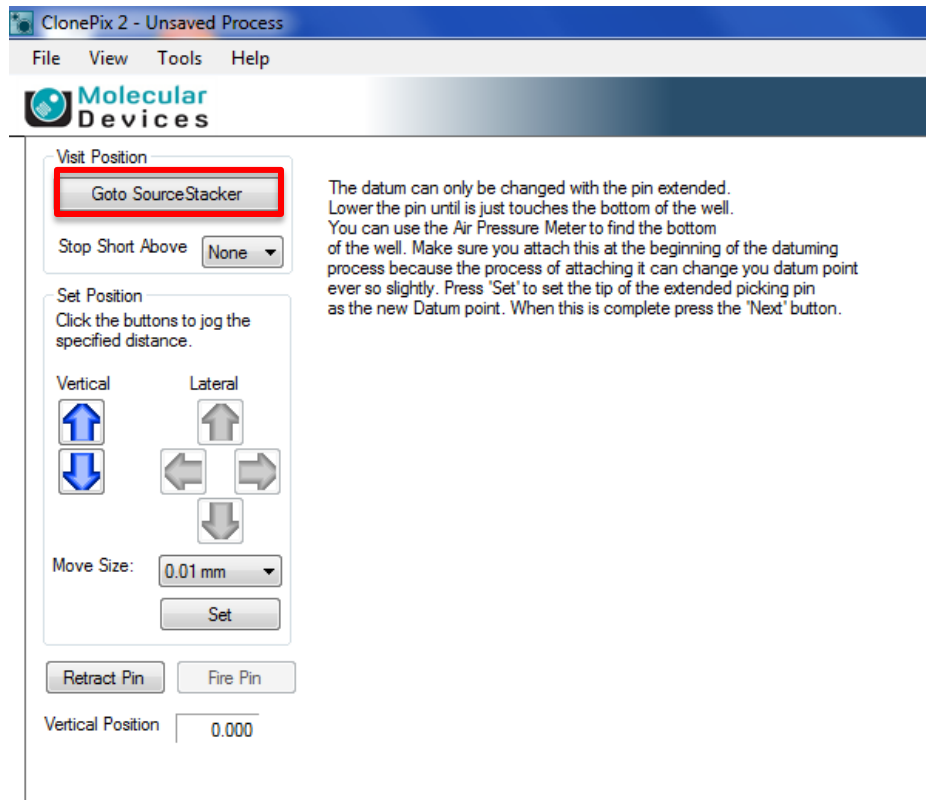
c) **Switch on the air pressure gauge** by twisting the **grey knob counterclockwise** so that the **handle is vertical** as pictured.



# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

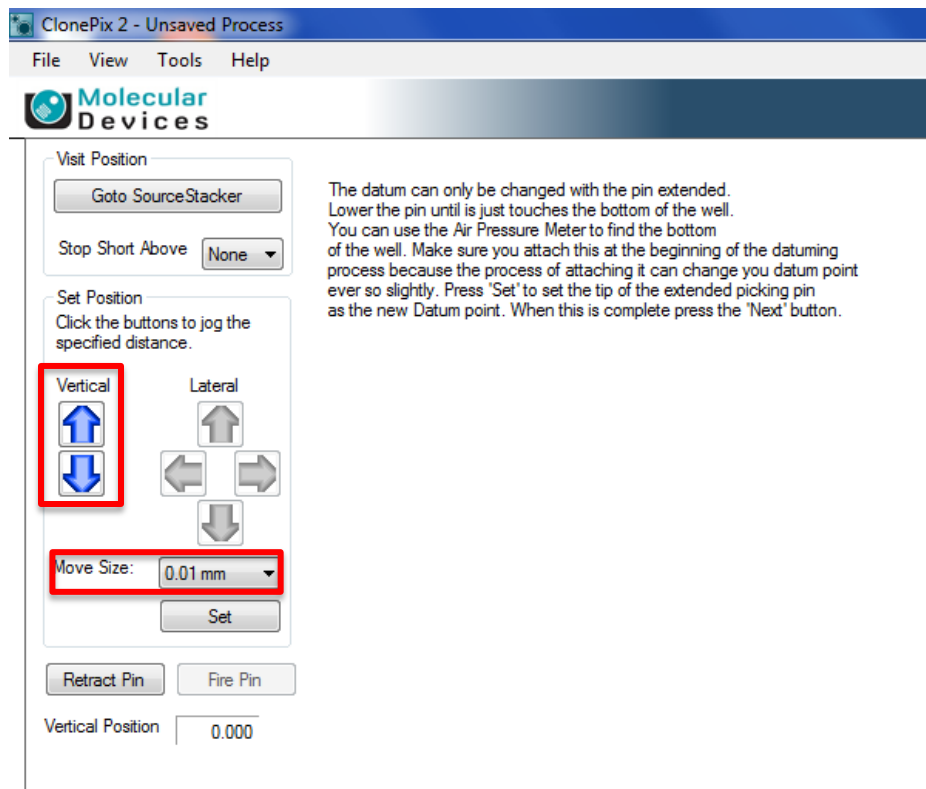
d) Click the **Goto SourceStacker** button - you will now see the **primary picking pin extend** to the bottom of the well and the **air pressure gauge readout value** will increase ( $>0.25$ ).



# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a reference point for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

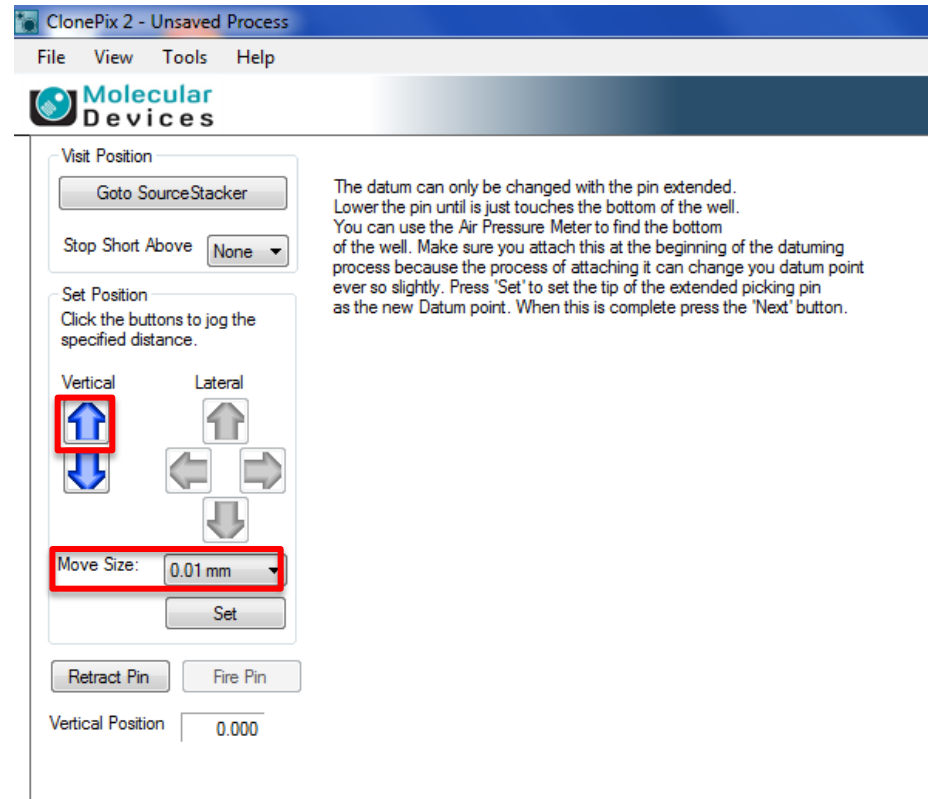
e) Use the **vertical arrow keys** to adjust the **Z position** of the **primary picking pin** until the **air pressure gauge readout color** changes to **green**. Select **move size increments** from the **dropdown** to refine the positioning.



# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

f) Once you have refined the position such that the indicator just turns **green** at the *smallest move increment (0.01 mm in the dropdown)*, **move the pin up** by selecting the **1 mm move size setting** and clicking the **up key once** (the **air pressure gauge readout** will now change back to **red**).

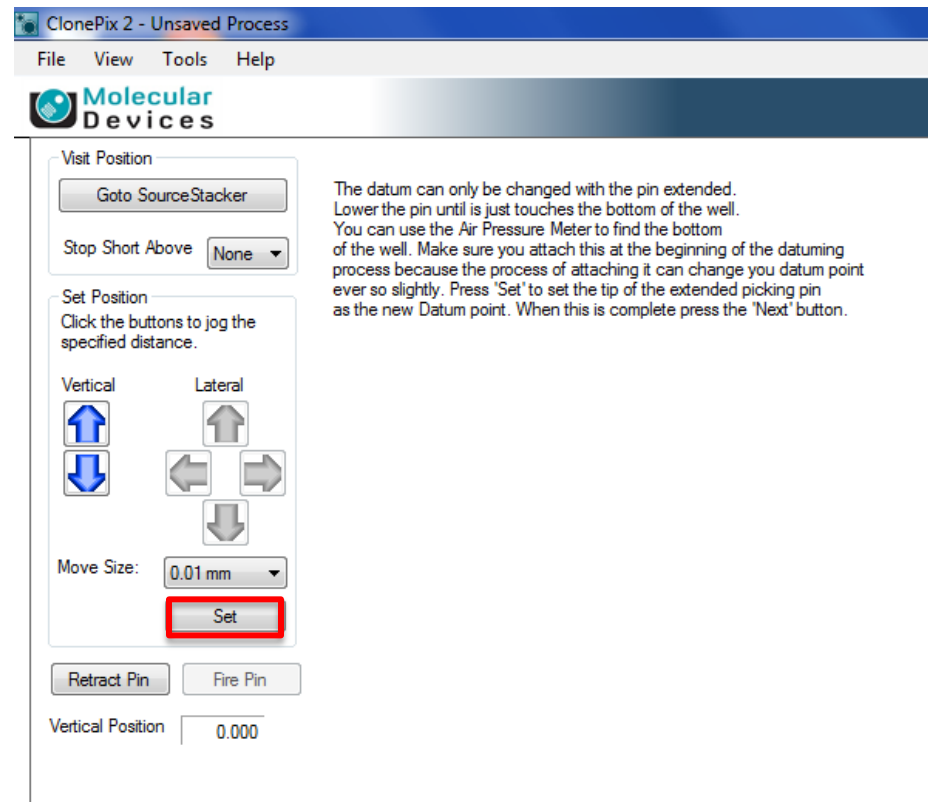




# Adjusting Primary Picking Pin Z Datum Position

10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

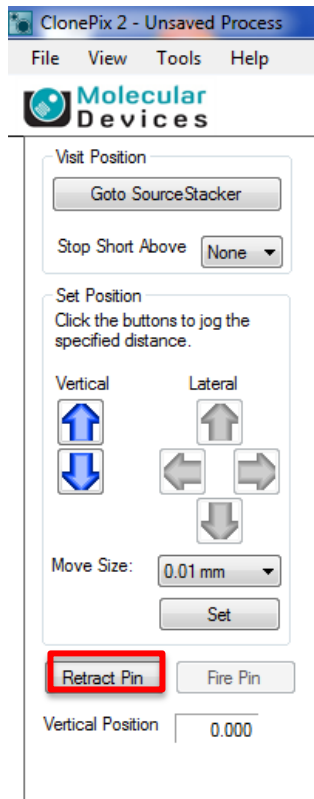
**g)** Click the **Set** button to save the current **Z datum position**. You will be prompted to confirm your choice to save the position – click **Yes**.



# Adjusting Primary Picking Pin Z Datum Position

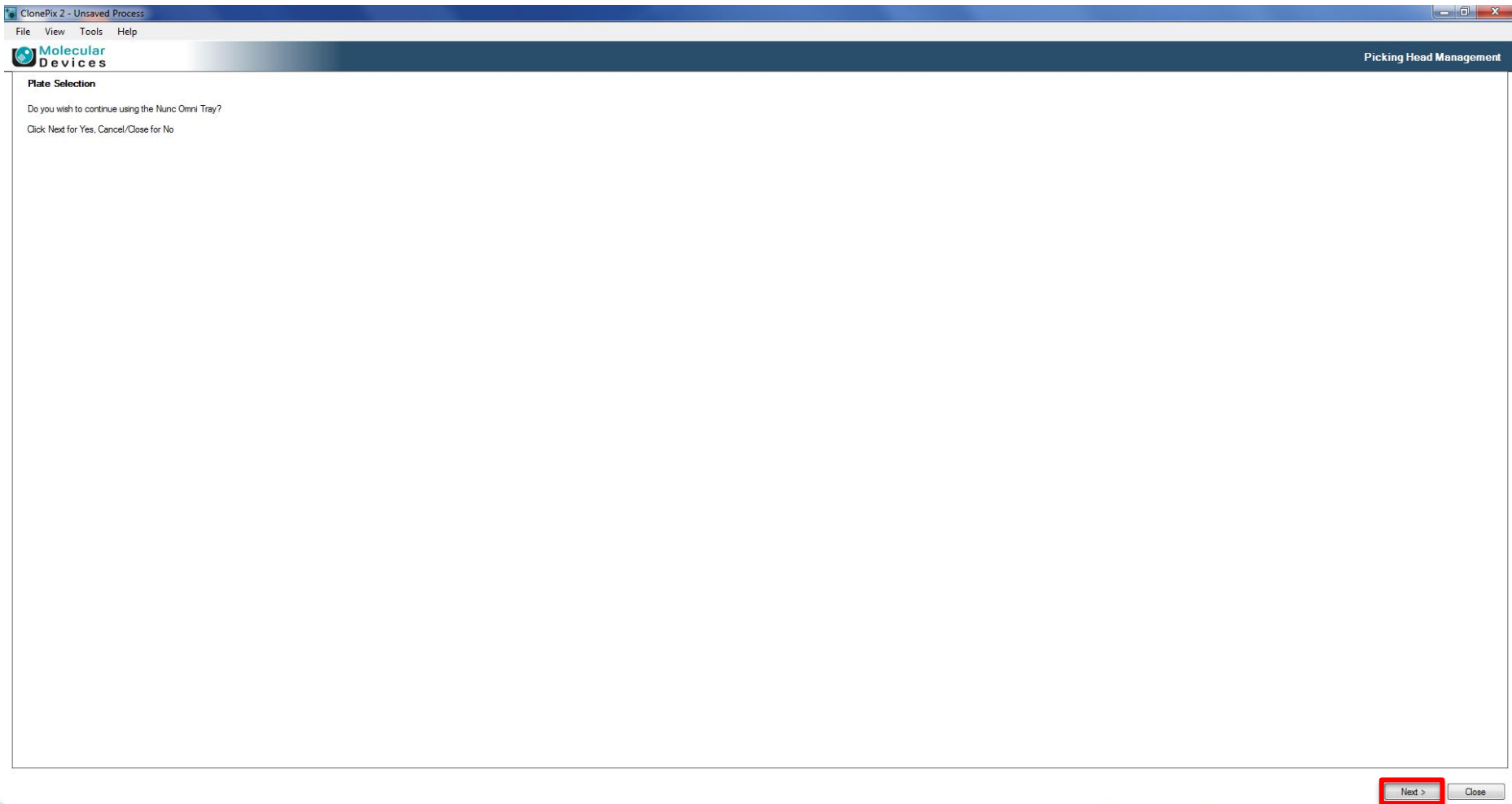
10. The **primary picking pin Z datum position** provides a **reference point** for the instrument to accurately pick your colonies while preventing collision with the plate bottom. This adjustment step ensures proper configuration of this position:

**h) Click on Retract Pin. Disconnect the air pressure tubing from the primary picking pin, then reconnect the picking tubing to the pin. Click Next to proceed.**



# Confirm Plate Type

11. The **Plate Selection** dialog now appears. Click the **Next** button to **confirm** that you want to **continue** to the next step using the **same source plate type** that you have been working with up to this point in the process.

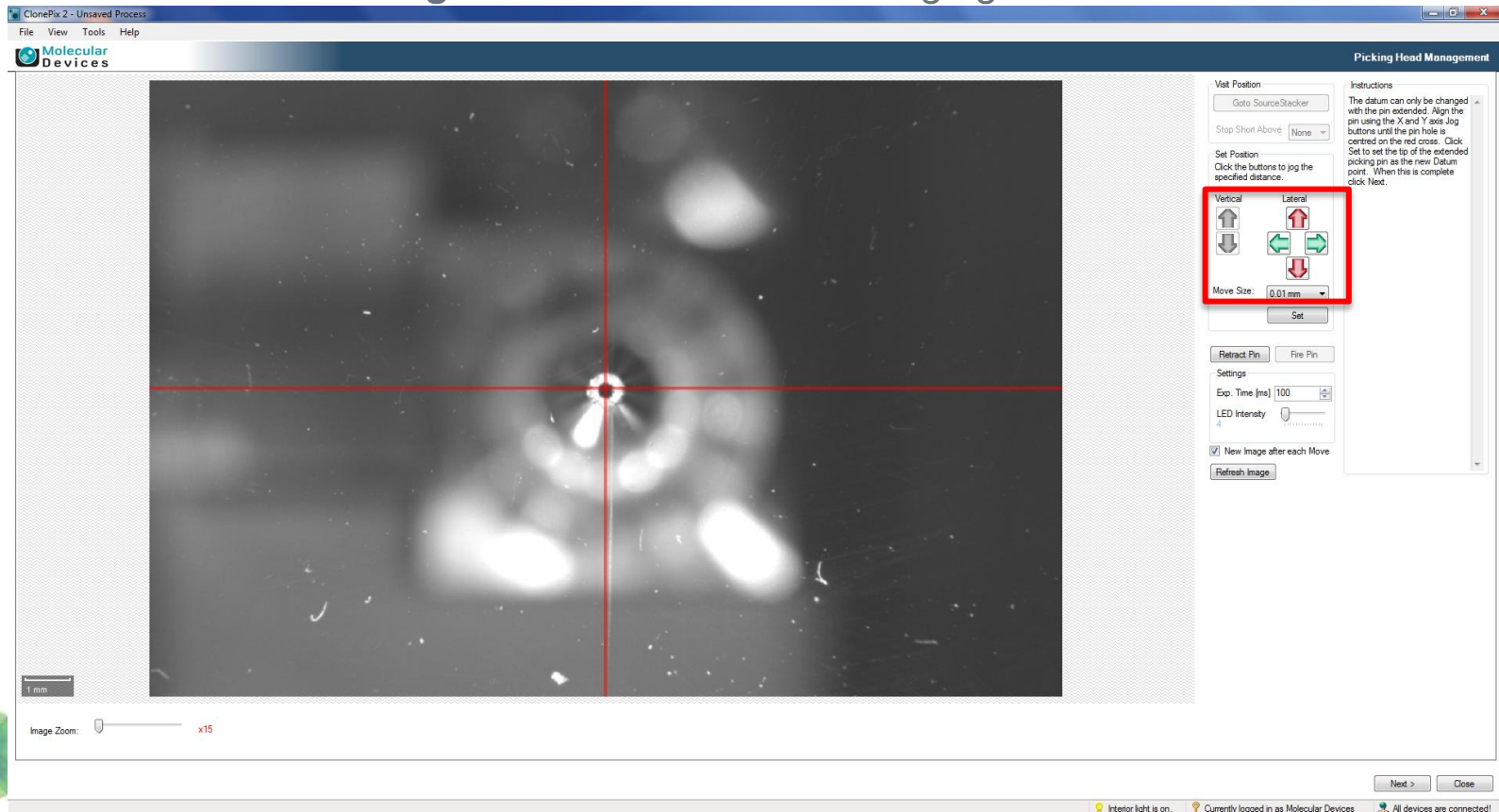


# Adjusting Primary Picking Pin XY Alignment

12. This next step allows for **adjustment** of the **primary picking pin XY alignment**. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

- a) Click on the **lateral arrow keys** to adjust the position of the **primary picking pin** so that the **center** of the **pin** aligns as closely as possible with the **intersection** of the **red crosshairs**. Change the **Move Size** via the **dropdown** as needed to refine positioning.

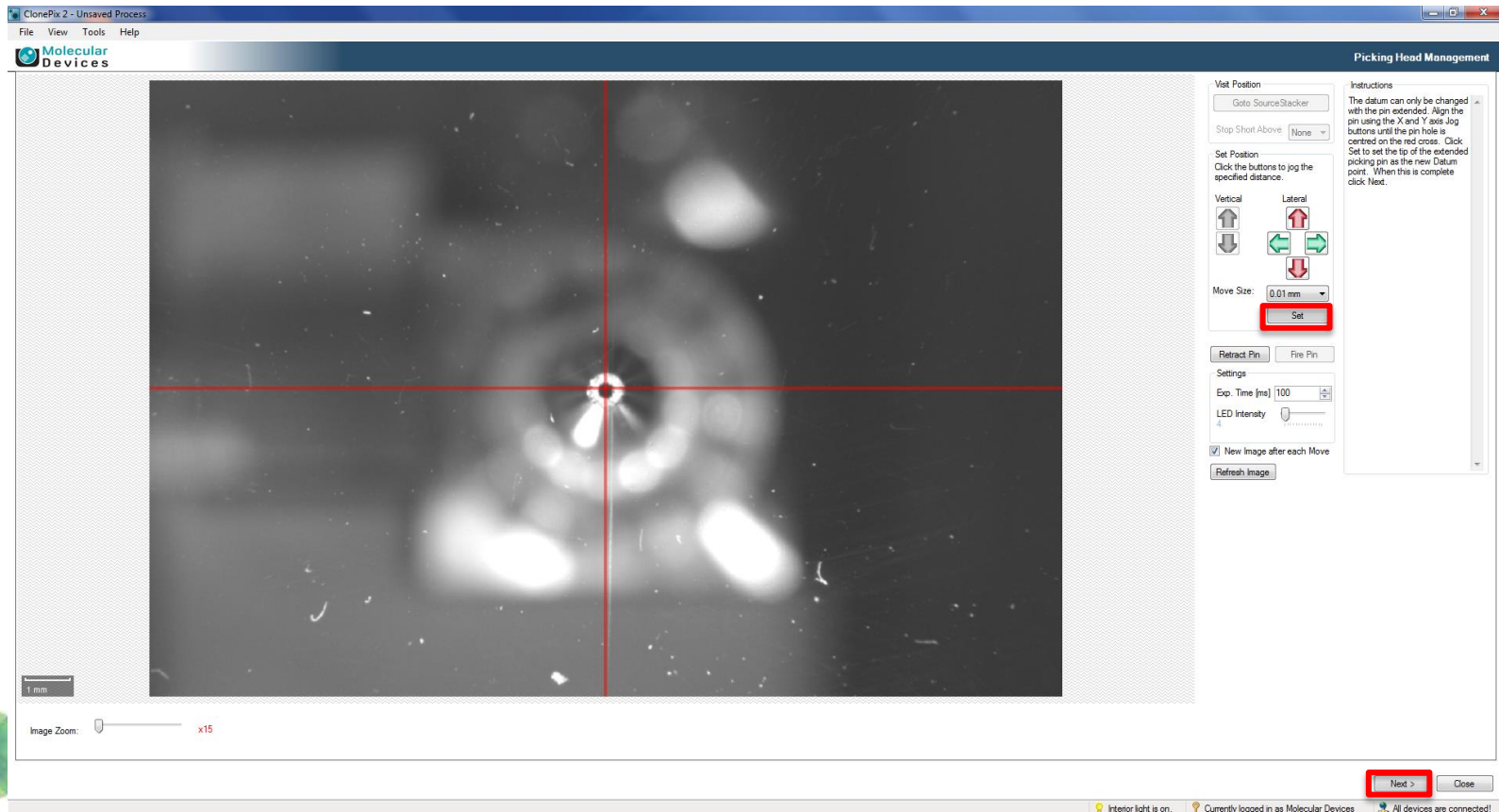
**NOTE:** The **arrow keys** move the **pin** itself (NOT the crosshairs). Please click **once**, then **wait** for the **image** to **refresh** before clicking again.



# Adjusting Primary Picking Pin XY Alignment

12. This next step allows for adjustment of the primary picking pin XY alignment. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

- b) Once the positioning is correct, click the **Set** button. You will then see a **dialog** that will ask you to **confirm save** of the **datum point** – click **Yes**.
- c) Click the **Next** button to proceed to the next step.



# Adjusting Picking Pins #2-8 XY Alignment

13. This next step allows for **adjustment** of the **#2-8 picking pins XY alignment**. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

- Click on the **lateral arrow keys** to adjust the position of the **current picking pin** so that the **center** of the **pin** aligns as closely as possible with the **intersection** of the **red crosshairs**. Change the **Move Size** via the **dropdown** as needed to refine positioning.

**NOTE:** The **arrow keys** move the **pin** itself (**NOT** the crosshairs). Please click **once**, then **wait** for the **image** to **refresh** before clicking again.

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Align the Primary Pin  
Ensure the pin is extended. Jog the Primary Pin around until the red cross is aligned over the centre. When this is complete you can proceed to Datum the other pins. You can not set the datum for the primary pin here, this is performed elsewhere. Here we are setting the datum points of the other pins relative to the primary pin. Use the Previous/Next Pin buttons to move through the picking pins.

Retract Pin  
Extend Pin  
Previous Pin Pin 1 Next Pin

Picking Head Management

Vial Position  
Gate Indexed/NozzleHead  
Stop Shot Above None

Set Position  
Click the buttons to jog the specified distance.

Vertical Lateral  
Move Size: 0.01 mm  
Set

New Image after each Move  
Refresh Image

Settings  
Exp. Time [ms] 100  
LED Intensity

Pin	Datum Adj. (X, Y)
1	Primary Pin
2	(-0.03413427, -0.1906661)
3	(-0.01951265, -0.05644594)
4	(0.0110141, -0.2349299)
5	(-0.08171952, -0.1827612)
6	(-0.002403617, -0.2120091)
7	(0.07930684, -0.2756613)
8	(-0.01829326, -0.1512002)

Image Zoom: x12

Next > Close

Interior light is on. Currently logged in as Molecular Devices All devices are connected!

ECULAR VICES

# Adjusting Picking Pins #2-8 XY Alignment

13. This next step allows for adjustment of the #2-8 picking pins XY alignment. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

- b) Once the positioning of the **current pin** is correct, click the **Set** button. You will then see a **dialog** that will ask you to **confirm save** of the **datum point** – click **Yes**.

**NOTE:** Changing the position of the **primary pin (Pin 1)** in this dialog will **not** affect its **XY datum**.

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Align the Primary Pin  
Ensure the pin is extended. Jog the Primary Pin around until the red cross is aligned over the centre. When this is complete you can proceed to Datum the other pins. You can not set the datum for the primary pin here, this is performed elsewhere. Here we are setting the datum points of the other pins relative to the primary pin. Use the Previous/Next Pin buttons to move through the picking pins.

Retract Pin  
Extend Pin  
Previous Pin Pin 1 Next Pin

Vial Position  
Goto Indexed/NozzleHead  
Stop Shot Above: None

Set Position  
Click the buttons to jog the specified distance.

Vertical Lateral  
Move Size: 0.01 mm  
Set

New Image after each Move  
Refresh Image

Settings  
Exp. Time [ms]: 100  
LED Intensity

Pin	Datum Adj. (X, Y)
1	Primary Pin
2	(-0.03413427, -0.1906661)
3	(-0.01951265, -0.0564694)
4	(0.0110141, -0.2349299)
5	(-0.08171952, -0.1827612)
6	(-0.002403617, -0.2120091)
7	(0.07930684, -0.2756613)
8	(-0.01829326, -0.1512002)

Image Zoom: x12

Next > Close

Interior light is on. Currently logged in as Molecular Devices. All devices are connected!

ECULAR VICES

# Adjusting Picking Pins #2-8 XY Alignment

13. This next step allows for adjustment of the #2-8 picking pins XY alignment. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

- b) Once the positioning of the **current pin** is correct, click the **Set** button. You will then see a **dialog** that will ask you to **confirm save** of the **datum point** – click **Yes**.

*NOTE: Changing the position of the primary pin (Pin 1) in this dialog will **not** affect its XY datum.*

- c) Click **Next Pin** to adjust the positioning of the remaining pins.

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Align the Primary Pin  
Ensure the pin is extended. Jog the Primary Pin around until the red cross is aligned over the centre. When this is complete you can proceed to Datum the other pins. You can not set the datum for the primary pin here, this is performed elsewhere. Here we are setting the datum points of the other pins relative to the primary pin. Use the Previous/Next Pin buttons to move through the picking pins.

Retract Pin  
Extend Pin  
Previous Pin Pin 1 **Next Pin**

Vial Position  
Goto Indexed/NozzleHead  
Stop Shot Above: None

Set Position  
Click the buttons to jog the specified distance.

Vertical Lateral  
Move Size: 0.01 mm  
Set

New Image after each Move  
Refresh Image

Settings  
Exp. Time [ms] 100  
LED Intensity

Pin	Datum Adj. (X, Y)
1	Primary Pin
2	(-0.03413427, -0.1906661)
3	(-0.01951265, -0.0564694)
4	(0.0110141, -0.2349299)
5	(-0.08171952, -0.1827612)
6	(-0.002403617, -0.2120091)
7	(0.07930684, -0.2756613)
8	(-0.01829326, -0.1512002)

Image Zoom: x12

Next > Close

Interior light is on. Currently logged in as Molecular Devices. All devices are connected!

MOLLECULAR DEVICES



# Adjusting Picking Pins #2-8 XY Alignment

13. This next step allows for adjustment of the #2-8 picking pins XY alignment. This step is important to ensure accurate colony picking. Follow these steps to adjust the alignment:

d) You will see that any updates to the picking pin XY datum points will be highlighted in red in the table below the Set Position dialog. Click the Next button to proceed to the system sanitizing steps.

ClonePix 2 - Unsaved Process

File View Tools Help

Molecular Devices

Align the Primary Pin

Ensure the pin is extended. Jog the Primary Pin around until the red cross is aligned over the centre. When this is complete you can proceed to Datum the other pins. You can not set the datum for the primary pin here, this is performed elsewhere. Here we are setting the datum points of the other pins relative to the primary pin. Use the Previous/Next Pin buttons to move through the picking pins.

Retract Pin

Extend Pin

Previous Pin Pin 1 Next Pin

Vial Position

Goto Indexed/NozzleHead

Stop Shot Above None

Set Position

Click the buttons to jog the specified distance.

Vertical Lateral

Move Size: 0.01 mm

Set

New Image after each Move

Refresh Image

Settings

Exp. Time [ms] 100

LED Intensity

Pin	Datum Adj. (X, Y)
1	Primary Pin
2	(-0.03413427, -0.1906661)
3	(-0.01951265, -0.0564594)
4	(0.0110141, -0.2349299)
5	(-0.08171952, -0.1827612)
6	(-0.002403617, -0.2120098)
7	(0.07930684, -0.2756613)
8	(-0.01829326, -0.1512002)

Image Zoom: x12

Next > Close

Molecular Devices

# Filling the Ethanol Wash Bath

14. The **Wash Bath** dialog now appears.

- Ensure that your **Wash Bottle** is at least  $\frac{3}{4}$  full with **70% Ethanol** (second bottle from **left front** side of instrument).
- Ensure that the **Auto Replenish Wash Bath** checkbox is checked, and keep the **default settings** (30 seconds for pump duration, 30 minutes for intervals between durations)

## Wash Bath

Click the 'Start' button to start filling the Wash Bath.

Click the 'Stop' button to stop filling the Wash Bath.

Wash Bath Pump

Auto Replenish Wash Bath

Duration the pump will be on for:  seconds

Interval between durations:  minutes

# Filling the Ethanol Wash Bath

14. The **Wash Bath** dialog now appears.

- a) Ensure that your Wash Bottle is at least  $\frac{3}{4}$  full with 70% Ethanol (second bottle from left front side of instrument).
- b) Ensure that the **Auto Replenish Wash Bath** checkbox is checked, and keep the default settings (30 seconds for pump duration, 30 minutes for intervals between durations) for the **Wash Bath Pump**.
- c) Click the **Start** button to initiate filling of the **Ethanol Wash Bath**.

## Wash Bath

Click the 'Start' button to start filling the Wash Bath.

Click the 'Stop' button to stop filling the Wash Bath.

Wash Bath Pump

Auto Replenish Wash Bath

Duration the pump will be on for:  seconds

Interval between durations:  minutes

# Filling the Ethanol Wash Bath

14. The **Wash Bath** dialog now appears.

- Ensure that your Wash Bottle is at least  $\frac{3}{4}$  full with 70% Ethanol (second bottle from left front side of instrument).
- Ensure that the **Auto Replenish Wash Bath** checkbox is checked, and **keep the default settings (30 seconds for pump duration, 30 minutes for intervals between durations)** for the **Wash Bath Pump**.
- Click the **Start** button to **initiate filling** of the **Ethanol Wash Bath**.
- Observe the **filling** of the **ethanol wash bath** in the **ClonePix 2** – allow the fill to **continue** until the **fluid reaches the exit tube** at the **back** of the bath (it is okay if you still see bubbles in the exit tube). Click the **Stop** button, then click **Next** to proceed.



Ethanol Wash Bath

## Wash Bath

Click the 'Start' button to start filling the Wash Bath.

Click the 'Stop' button to stop filling the Wash Bath.

Wash Bath Pump

Auto Replenish Wash Bath

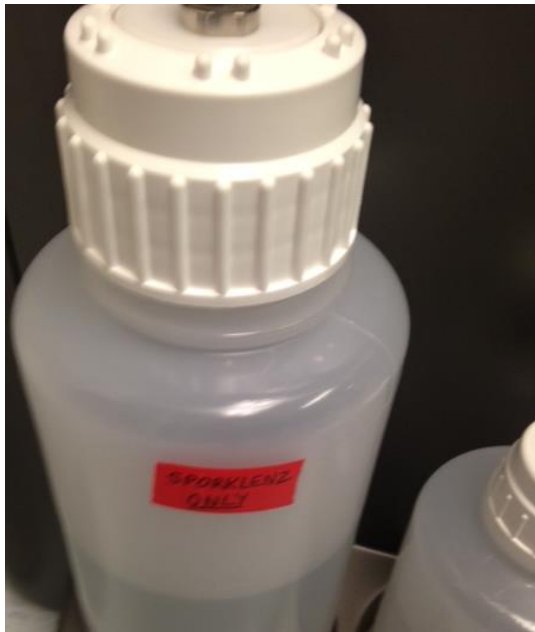
Duration the pump will be on for:  seconds

Interval between durations:  minutes

# Sanitizing Picking Pins

15. This next step ensures that the **picking pins & fluidics** are **flushed with sanitizing solution** (important for maintaining sterility during picking):

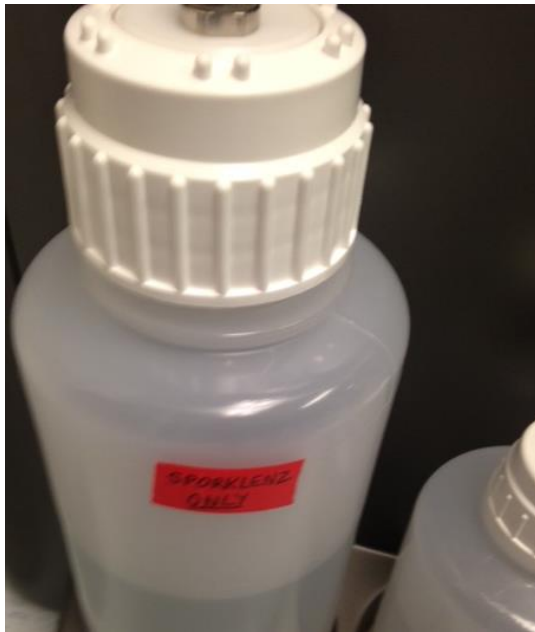
- a) Ensure that the **5L wash bottle** containing approximately **4L of SporKlenz solution** is attached to the **leftmost fluidics connection** on the ClonePix 2 instrument.



# Sanitizing Picking Pins

15. This next step ensures that the **picking pins & fluidics** are flushed with **sanitizing solution** (important for maintaining sterility during picking):

- a) Ensure that the **5L wash bottle** containing approximately **4L of SporKlenz solution** is attached to the **leftmost fluidics connection** on the ClonePix 2 instrument.
- b) Leave the **default** settings in the **Sanitise Options** dialog (**30/5/10** as pictured).



**Sanitise Pins**

Purge the pump system with sanitising agent

Sanitise Options

Number of cycles for the head to purge	30
Number of cycles the pins will be scrubbed in the Wash Bath	5
Number of seconds the Dryer will be on for	10

Start

# Sanitizing Picking Pins

15. This next step ensures that the **picking pins & fluidics** are flushed with **sanitizing solution** (important for maintaining sterility during picking):

- Ensure that the **5L wash bottle** containing approximately **4L of SporKlenz solution** is attached to the **leftmost fluidics connection** on the ClonePix 2 instrument.
- Leave the **default** settings in the **Sanitise Options** dialog (30/5/10 as pictured).
- Click the **Start** button in the **Sanitise Pins** dialog to begin **flushing** the ClonePix 2 with **SporKlenz**. You will see the **picking head** move to the **wash station** and the **pins** inserted into the **8 ports** within. You will see the **Purge Progress** dialog appear as well. Once this step is complete, click **Next**.

## Sanitise Pins

Purge the pump system with sanitising agent

### Sanitise Options

Number of cycles for the head to purge

30

Number of cycles the pins will be scrubbed in the Wash Bath

5

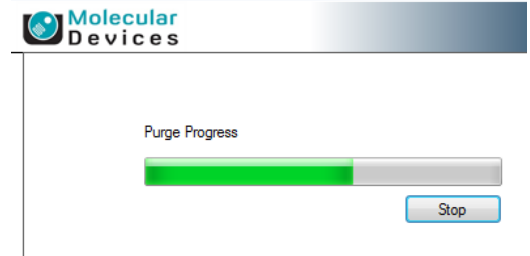
Number of seconds the Dryer will be on for

10

Start



Wash Station



# UV Sanitizing ClonePix 2 Deck

16. This next step ensures that the **ClonePix 2 instrument deck** is **sterilized** with **UV light** (important for maintaining sterility during picking):

- a) The **Ultra Violet Sanitise** dialog appears. Leave the **Duration of Ultra Violet exposure** setting at the **default** (600 seconds).

## Ultra Violet Sanitise

1. Check the bed of the instrument is clear of plates.
2. Perform any manual cleaning tasks.
3. Set the number of seconds for the UV light to be on.
4. Click the 'Begin' button to turn the light on. You can stop at any time by clicking the 'Stop' button.

While the UV sanitise is in progress you can check the fluid level in the waste bottles and make sure the feed bottles are full.

Ultra Violet Sanitise

Duration of Ultra Violet exposure:  seconds



# UV Sanitizing ClonePix 2 Deck

16. This next step ensures that the **ClonePix 2 instrument deck** is **sterilized with UV light** (important for maintaining sterility during picking):

- a) The **Ultra Violet Sanitise** dialog appears. Leave the **Duration of Ultra Violet exposure** setting at the **default (600 seconds)**.
- b) Click **Begin** to start **UV sterilization** – this step takes **10 minutes**. You will see the **UV light illuminate the deck**, as well as the **UV light indicator** on the instrument front panel illuminate.

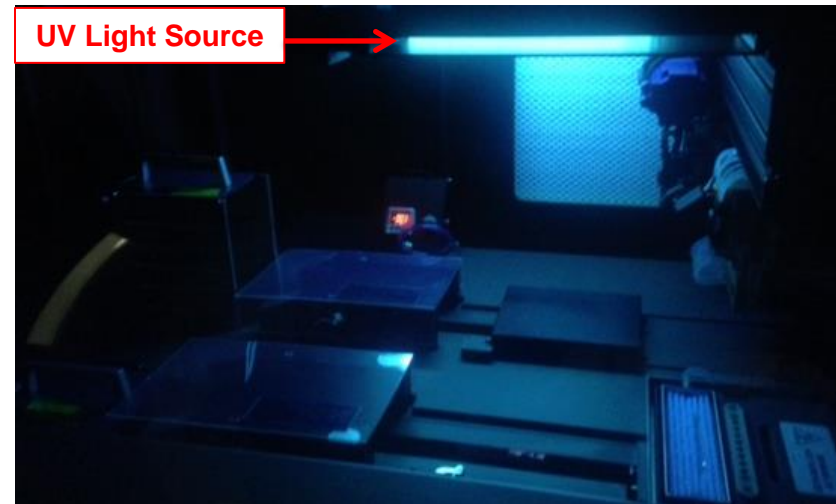
## Ultra Violet Sanitise

1. Check the bed of the instrument is clear of plates.
2. Perform any manual cleaning tasks.
3. Set the number of seconds for the UV light to be on.
4. Click the 'Begin' button to turn the light on. You can stop at any time by clicking the 'Stop' button.

While the UV sanitise is in progress you can check the fluid level in the waste bottles and make sure the feed bottles are full.

Ultra Violet Sanitise

Duration of Ultra Violet exposure:  seconds



# Check Wash & Waste Bottles

17. This next step is a reminder to check that the **wash bottles** are **full** and that the **waste bottles** are **empty**.

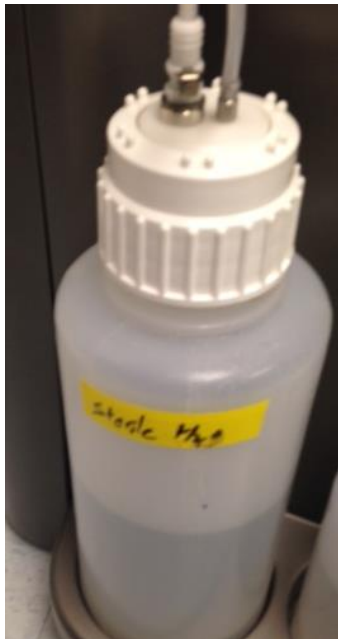
- a) Ensure that the **wash bottles (SporKlenz 5L, Sterile Water 5L, 70% Ethanol 2L)** are at least **75% full** and that the **waste bottles (wash & picking waste)** are **empty**.
- b) Click the **Next** button to proceed.

The screenshot shows a software window titled "ClonePix 2 - Unsaved Process" with a menu bar (File, View, Tools, Help) and a toolbar with a "Prepare For Pick Run" button. The main interface has a sidebar with a list of steps: Pin Fire Test, Alignment, Check Wash Bath, Wash Bath Utility, Sanitise Pins In, Sanitising Agent, UV Sanitise, **Check Bottles** (highlighted), and Sanitise Pins. The main content area displays the "Check Bottles" step with the instruction: "Check the fluid level in the wash bottle and make sure the waste bottle is empty." At the bottom of the window, there is a status bar with a "Start Time: 9:14:28 AM" on the left and a row of buttons: "< Back", "Next >" (highlighted with a red box), and "Cancel". To the right of the buttons are three status indicators: "Interior light is on.", "Currently logged in as Molecular Device", and "All devices are connected!". The "CULAR ICES" logo is visible in the bottom right corner.

# Rinsing Picking Pins

18. This next step ensures that the **picking pins & fluidics** are **flushed** with **sterile water** (important for removing sanitizing solution & maintaining sterility during picking):

- a) **Disconnect the 5L wash bottle containing SporKlenz** and then **connect the 5L wash bottle containing approximately 4L of Sterile Water** is **attached to the leftmost fluidics connection** on the ClonePix 2 instrument.
- b) Leave the **default** settings in the **Sanitise Options** dialog (30/5/10 as pictured).



**Sanitise Pins**

Purge the pump system to clear any air bubbles or sanitising agent

Sanitise Options

Number of cycles for the head to purge	30
Number of cycles the pins will be scrubbed in the Wash Bath	5
Number of seconds the Dryer will be on for	10

Start

# Rinsing Picking Pins

18. This next step ensures that the **picking pins & fluidics** are flushed with **sterile water** (important for removing sanitizing solution & maintaining sterility during picking):

- Disconnect the 5L wash bottle containing SporKlenz and then connect the 5L wash bottle containing approximately 4L of Sterile Water is connected to the leftmost fluidics connection on the ClonePix 2 instrument.
- Leave the default settings in the Sanitise Options dialog (30/5/10 as pictured).
- Click the **Start** button in the **Sanitise Pins** dialog to begin flushing the ClonePix 2 with **Sterile Water**. You will see the **picking head** move to the **wash station** and the **pins** inserted into the **8 ports** within. You will see the **Purge Progress** dialog appear as well.

## Sanitise Pins

Purge the pump system to clear any air bubbles or sanitising agent

### Sanitise Options

Number of cycles for the head to purge

30

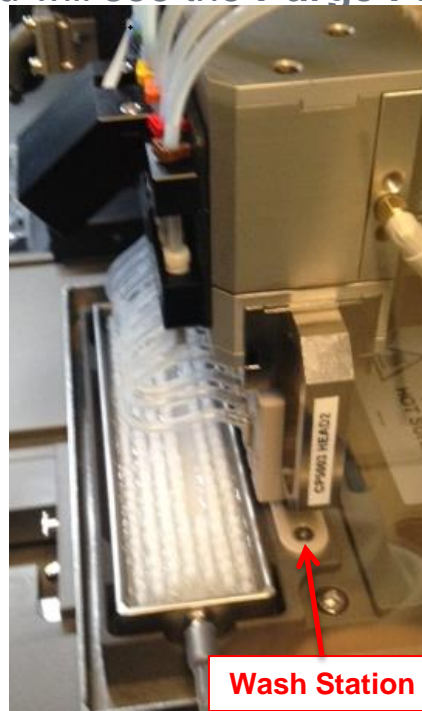
Number of cycles the pins will be scrubbed in the Wash Bath

5

Number of seconds the Dryer will be on for

10

Start



Molecular  
Devices

Purge Progress



Stop

# Rinsing Picking Pins

18. This next step ensures that the **picking pins & fluidics** are flushed with **sterile water** (important for removing sanitizing solution & maintaining sterility during picking):

d) The **picking head** will now **automatically** move to the **ethanol wash bath** and **scrub** the **picking pins**, followed by **drying** of the pins in the **halogen drying station**. You will see the **Washing Pins** indicator followed by the **Drying Pins** indicator on your **computer screen** during these steps. At the conclusion of the drying step, click **Next** to continue.



Ethanol Wash Bath



Halogen Drying Station



# Conclusion of Process

19. This final step will result in **completion** of the **Replace Head** process:

- a) You will see the **Process Completed** screen appear. Click the **Finish** button to return to the **main menu**. Click the **Finish** button.

## Process Completed

Click the Finish button to return to the process properties.



- b) Within the instrument, the following will occur **automatically**:

- **Re-homing** of the **picking head**
- **Re-lidding and return** of the **source plate** to the **source stacker**
- **Closing** of the **illumination cover** (at the imaging station).

- c) Be sure to **remove** your **empty source plate** from the **source stacker** before proceeding with setting up a **Pick Run**.

# Support Resources

- Go to the HELP menu within ClonePix 2 Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- Request Support: <http://mdc.custhelp.com/app/ask> or via email [support@moldev.com](mailto:support@moldev.com)
- Technical Support can also be reached by telephone:
  - 1 (800) 635-5577
  - Select options for Tech Support → Biotherapeutics Products → ClonePix Instruments





# **MOLECULAR** DEVICES

ADVANCING PROTEIN AND CELL BIOLOGY