

# QPix QUICK GUIDE – Pin Calibration

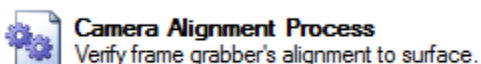
For more complete information, refer to page 244 of the User Manual

## Required Materials

- QTray with Black Foam
- Light colored lab tape
- 96 pin E. coli picking head (recommended)

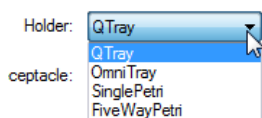
Objective: Align the picking head with the camera to allow for accurate colony picking.

1. Under “QPix Utility Processes,” double click “Camera alignment process”

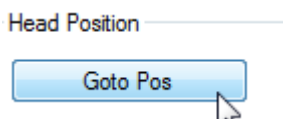


2. Select “QTray” from the holder list.

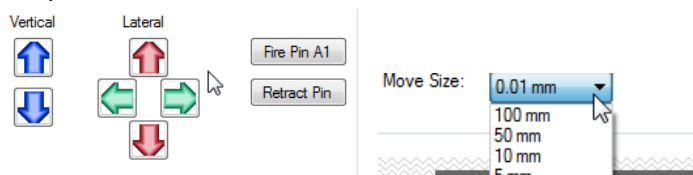
*NOTE: if you are not using the QTray provided by Molecular Devices, make sure you select the appropriate container from the menu.*



3. Make sure you have installed a picking head. (Molecular Devices recommends using the 96 pin E. coli picking head. Click “Go To Pos” to send head to the picking position.



4. Click “Fire Pin A1” and use the blue arrows to lower the pin until it pokes a hole in the lab tape. You can adjust how far the pin moves down at every step with the “Move Size” drop down menu.



5. Click “Retract Pin” and then “Go to Camera”



6. Use the red and green arrows labeled “Lateral” to center the hole left in the lab tape in the red crosshairs in the image (see picture above). You can adjust how far the camera moves each click with the “Move Size” drop down menu.
7. Once you are satisfied, hit “Set” and “Next” to finish the process.

