

# Rearranging

## Application Guide



# Contents

<b>Table of Figures</b> .....	<b>4</b>
<b>Rearranging</b> .....	<b>6</b>
What is Rearranging? .....	6
Preparation .....	7
Fitting the Head .....	7
Maintaining a Picking Head .....	7
Replacing Pins .....	8
Loading Source Plates .....	8
Loading Destination Plates .....	9
Deep Well Plates .....	9
Alternative Locations for Source and Destination Plates .....	10
<b>Defining Well Data</b> .....	<b>12</b>
Note .....	12
Define Source Data Only .....	12
Importing a Source Text File.....	12
Selecting Wells on a Plate Representation .....	16
Recording Barcodes .....	17
<b>Rearranging Software</b> .....	<b>18</b>
Overview .....	18
Sign On .....	19
Rearranging Setup .....	20
Description .....	20
Head .....	20
Source.....	21
Destination .....	23
Sterilize .....	24
Barcodes .....	25
Start .....	27
Running a ReArraying Routine.....	28
<b>Biology Guide</b> .....	<b>32</b>
Preparing Media .....	32
Luria-Bertani medium (LB) – per litre.....	32
LB + 8% Glycerol – per litre .....	32
LB Agar – per litre .....	32
Antibiotic Preparation.....	32
<b>Glossary of Terms</b> .....	<b>33</b>



<b>Appendix A</b> .....	<b>35</b>
Description of Toolbar and Menu Items .....	35
<b>Appendix B</b> .....	<b>38</b>
GTX Messaging Server .....	38
Set the Delay Time .....	42
<b>Index</b> .....	<b>43</b>
<b>Contact Details</b> .....	<b>44</b>



# Table of Figures

Figure 1: The Principle of Rearranging.....	6
Figure 2: QPix Bed Layout.....	8
Figure 3: Bioassay Tray Locator.....	9
Figure 4: An Example Import File.....	13
Figure 5: Rearranging Data Selection Dialog.....	14
Figure 6: Import Data Message.....	14
Figure 7: Rearranging Data Selection Dialog.....	15
Figure 8: Import Data Message.....	15
Figure 9: Import Rearranging Data.....	16
Figure 10: Rearranging Data Plate Representation.....	17
Figure 11: Routine Prompt.....	18
Figure 12: Sign on dialog.....	19
Figure 13: Welcome.....	19
Figure 14: Head.....	20
Figure 15: Source Settings.....	21
Figure 16: Control Data.....	22
Figure 17: Destination.....	23
Figure 18: Sterilize.....	24
Figure 19: Barcodes.....	25
Figure 20: Invalid Barcode Dialog.....	26
Figure 21: Start.....	27
Figure 22: Prepare Script.....	28
Figure 23: Wells Found.....	28
Figure 24: View Script.....	28
Figure 25: Example Script Code.....	29
Figure 26: Plates Required.....	29
Figure 27: Change Head.....	29
Figure 28: Destination Plate Locations.....	30
Figure 29: Source Plate Locations.....	30
Figure 30: Set Up and Ready.....	30
Figure 31: Script Progress.....	30
Figure 32: Script Complete.....	31
Figure 33: Carry Over Pin and Well Settings.....	31
Figure 34: Outlook Express Tools Menu.....	39
Figure 35: Outlook Express Security Settings.....	39
Figure 36: Messaging Server Configuration.....	40
Figure 37: New Contact.....	40



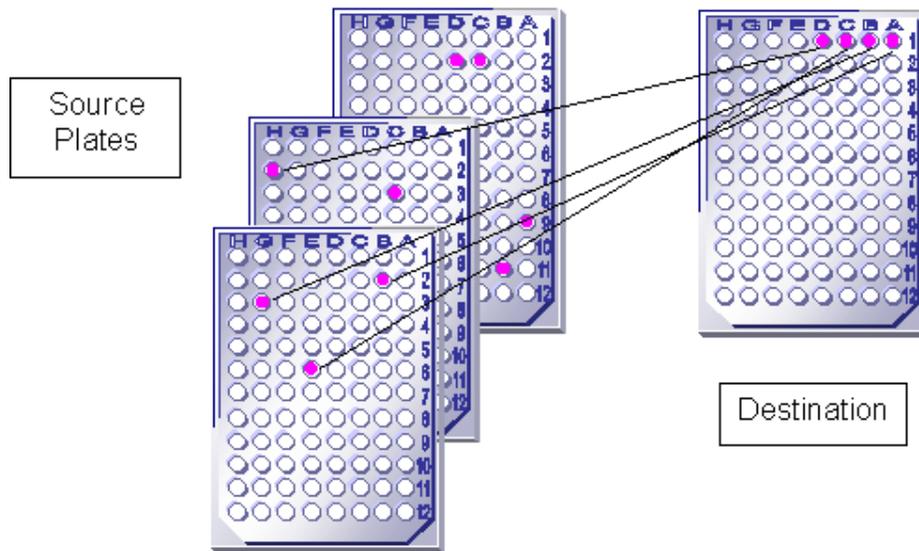
**Figure 38: Configure Email** ..... 41  
**Figure 39: Windows System Properties** ..... 41  
**Figure 40: Contact Details** ..... 42



# Rearranging

## What is Rearranging?

The objective of rearranging is to **REDISTRIBUTE** collected samples into **NEW** well plates. That is, to take user-selected samples from wells on source WellPlates and deposit them into wells on a destination well plate as shown below.



**Figure 1: The Principle of Rearranging**

Of course the robot can accommodate, simultaneously, many more source well plates and destination well plates in one rearranging sequence than that shown in the diagram.

**QPix** can accommodate up to 21 well plates. Usually 6 source well plates and 15 destination well plates.

If more destination or source plates are required to complete a rearranging sequence, you will be prompted at the relevant time to replace the well plates.

## Preparation

Rearranging uses the following robot accessories:

- Replicating/rearranging table (**QPix**)
- Picking head

Before preparing the robot for Rearranging, check that the workbed area of the base plate is clean and free from debris. The working volume of the machine should also be UV sterilised before setting up the bed for any routine, (see **General Maintenance** section in main robot manual).

## Fitting the Head

If the robot has been used previously for any other application, the head may need to be changed. The head must be removed and replaced with the correct head (see **General Maintenance** section in main robot manual).

## Maintaining a Picking Head

Cleaning the head can be a long process, but it is vital for good results. The head should be cleaned every time a picking routine is completed. Handle all parts with care when cleaning to avoid bending any pins or losing springs.

- Use a flat bladed screwdriver to unscrew the 10 screws. A support is needed that allows for stability of the head but keeps the pins suspended (e.g. a pipette tip box top, or the Robot wash bath).
- Very carefully remove the top plate to expose the springs and the tops of the pins. **Be very careful not to lose the springs!**
- Remove the pins from the main body and place in a container suitable for sonic cleansing. Sonicate the pins, body and springs for 10 minutes in a 1% aQu Clean.
- Remove the pins, plate and springs from the sonicator and rinse thoroughly in distilled water.
- Blow through the plate with an airline and dry thoroughly, along with the pins and springs.
- Insert the pins into the holes of the body, followed by the springs (all pins should fall down under their own weight).
- Place the top plate over the pins (aligned with the main body) and screw into place with the 10 screws. **Do not over tighten.**



## Replacing Pins

Occasionally colonies may be missed. This is most commonly due to bent or sticky pins.

Bent pins can be easily identified by carefully checking the head before each use. Hold the head so that the tips of the pins are at eye level and look along each row of pins from each side of the head, the problem pin should be easily spotted.

Remove and replace the damaged pin as follows:

- Use a flat bladed screwdriver to unscrew the screws. A support is needed that allows for stability of the head but keeps the pins suspended (e.g. a pipette tip box top).
- Very carefully remove the top of the head to expose the springs and the tops of the pins.
- Using the end of an Allen key, push the damaged pin up from the bottom of the head until. Carefully remove the pin.
- Place a new pin into the hole that has been vacated, ensuring that the pin does not stick (as above). Place the top of the head over the pins and tighten the screws. **Do not over tighten.**

## Loading Source Plates

The source plates may contain the colony libraries that are to be rearrayed, they may also contain just DNA, for example PCR products.

### QPix

Traditionally the 2 right-most containers are used for the source plates and the 3 left-most containers are used for the destination plates. For possible variations see the section "**Alternative locations for Source and Destination plates**" (page 10).

On **QPix**, the source plates are loaded on to the source plate holders. The source plate holders are fixed to a removable carrier which fits onto the robot bed.

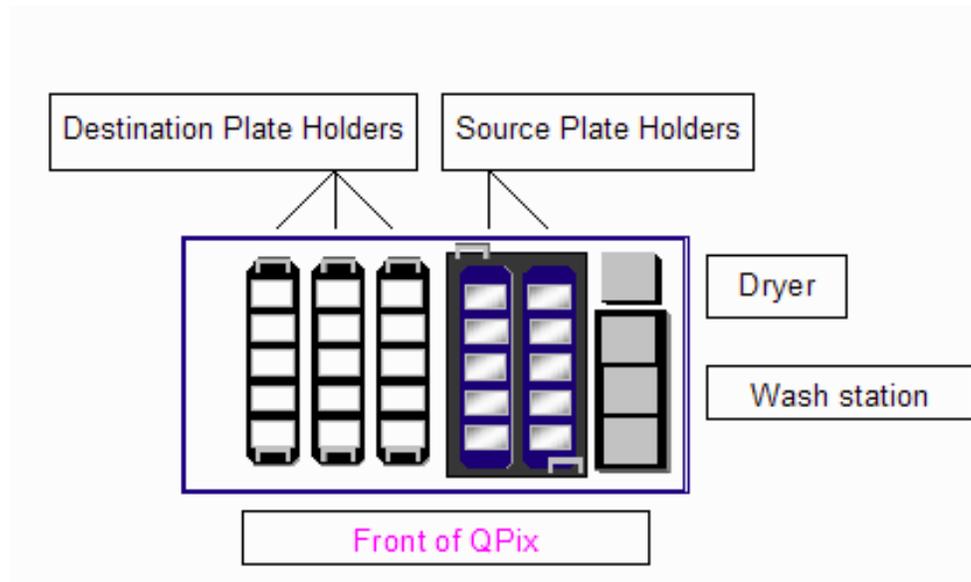


Figure 2: QPix Bed Layout

To fit the source plate carrier, take out the black thumbscrews that hold the gridding table. Remove the bioassay tray locator from the back left corner of the light table, which is held by the white thumbscrew.



**Figure 3: Bioassay Tray Locator**

Position the source plate carrier so that the four locating lugs are seated over the locating posts on the **QPix** bed and the bioassay tray holder/imaging area is obscured.

Replace the black thumbscrews.

Plates should be loaded **with the lids off** and well A1 should be facing the front right of the machine. Make sure that all plates are correctly inserted into each location.

## Loading Destination Plates

### QPix

Traditionally the 2 right-most containers are used for the source plates and the 3 left-most containers are used for the destination plates. For possible variations see the section "**Alternative locations for Source and Destination plates**" (page 10).

The destination plates are loaded directly on to plate holders which fit on to the robot bed.

The well plates are loaded without lids and aligned so that well A1 is in position front-right. The well plates are secured in place by the notches in the plate holders.

### Deep Well Plates

When using deep well plates you must ensure that the following plate properties are correctly set:

**InkingDepth** – must be set so that the pins reach no more than 5mm into the media. This is to ensure that media isn't drawn back into the head when the pins are retracted.

**UseLowerBedLevel** – (QPix only) must be set to TRUE (-1)

### QPix

**QPix** has a section that lifts out to reveal a recess which will accommodate deep well plates if required. To allow access to the lower bed level undo the four holding screws on the removable plate.



# Alternative Locations for Source and Destination Plates

Please Note: Although the physical objects are actually referred to as Source **objectname** and Destination **objectname**, it is possible to designate which area is to be defined as the source and destination. For example:

## On a QPix....

This is slightly more complicated because there are more permutations.

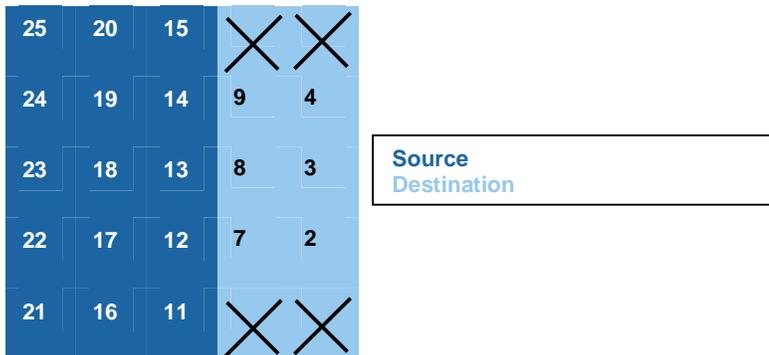
When picking source data in the ReArraying module on a **QPix**, the head cannot address locations on the back row, the front row or the leftmost well plate holder. The unavailable locations are shown in the diagrams with a **X** through them.

When depositing the destination data the head will align directly above well plates and can therefore address any location on the bed.

### Standard Well Plates

#### Option 1

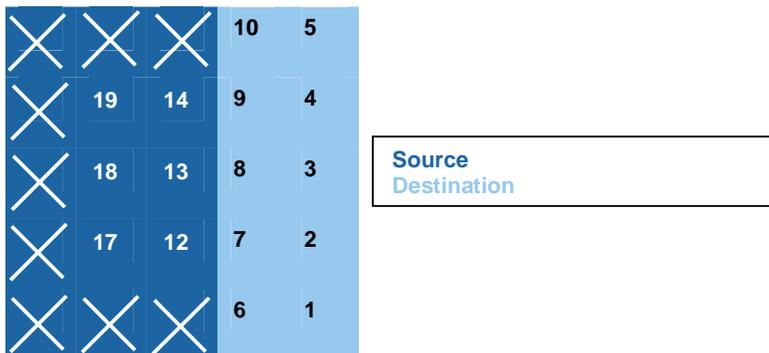
Up to 6 source well plates and up to 15 destination well plates.



**Note:** Make sure that you select **Source Plate Holder (1X3)** as the **Source** container.

#### Option 2

Up to 6 source well plates and up to 10 destination well plates.



**Note:** When using locations 1 - 10 as the destination, a **Source Plate Holder (1X5)** as the Destination container and a **Destination Plate Holder (1X3)** as the Source Container must be selected.



### Deep well Plates

**Note:** Deep well plates (indicated by white boxes) can only be used as **either** source **or** destination not as both.

#### Option 1

Up to 6 **standard** well plates as source and up to 15 **deep well** plates as destination.

25	20	15	×	×
24	19	14	9	4
23	18	13	8	3
22	17	12	7	2
21	16	11	×	×

Destination (Deep well plates)  
Source (Deep well plates)

**Note:** Make sure that you select **Source Plate Holder (1X3)** as the Source container.

#### Option 2

Up to 6 deep well plates as source and up to 10 standard well plates as destination

×	×	×	10	5
×	19	14	9	4
×	18	13	8	3
×	17	12	7	2
×	×	×	6	1

Source (Standard well plates)

**Note:** When using locations 1 - 10 as the destination, a **Source Plate Holder (1X5)** as the Destination container and a **Destination Plate Holder (1X3)** as the Source Container must be selected.



# Defining Well Data

---

## Note

Well data can be either "Source Data Only" or "Source **and** Destination Data".

## Define Source Data Only

"Source Data Only" can be defined in one of two ways:

Importing a Source text file.

or

Selecting wells on a plate representation.

## Importing a Source Text File

A data file defining the selection of source sample wells could be created *outside* the Rearranging module. This can be done using any software that produces output in text format e.g. Windows<sup>®</sup> Notepad or Microsoft<sup>®</sup> Excel<sup>®</sup>.

**Note:** If the data file is created using an application other than QSoft, you must ensure that the file is saved as a **comma delimited text** file. It may have the extension TXT or IMP

The data file would then be imported into the rearranging module at run time. You may need to change the "Files of Type" list box option to "All Files (\*.\*)" in order to locate your file.



### Rearranging Import File Format (Source only)

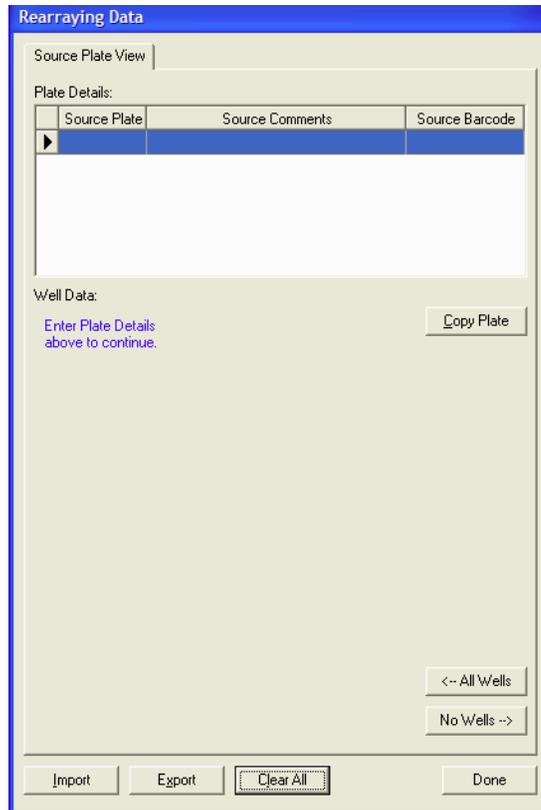
The Rearranging Import File is made up of a list of keywords, well references and plate numbers. All keywords should be uppercase and suffixed with a colon and a space (i.e. ": "). The default import file extension is IMP, however, any ASCII file can be read.

For each plate to be rearranged, there must be a **PLATE** keyword followed by a list of well references. The import or export file specification is as follows:

PLATE	Required – Source plate number to be rearranged. The plate numbers must be contiguous.
BARCODE	Optional - The barcode on the plate to be checked when the plate is loaded.
COMMENT	Optional - any comments for/about this plate.
<p><b>Figure 4: An Example Import File</b></p>	<pre> PLATE: 1 BARCODE: 12345 COMMENT: Some Comments A3 B3 B9 C1 C5 D5 E7 F5 F11 G9 PLATE: 2 B3 B7 B9 PLATE: 3 COMMENT: Some notes about this plate B2 C3 D4 E5                     </pre>

To import the file - when the Rearranging software is running - select the Source tab and click the Source Data button. The following screen is displayed:

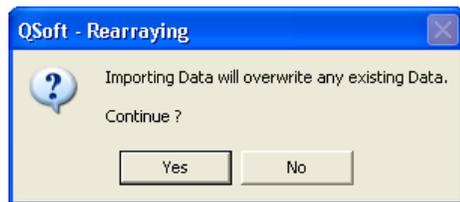




**Figure 5: Rearranging Data Selection Dialog**

Click the Import button.

If there is rearranging data already defined for the current routine, the following screen will appear.



**Figure 6: Import Data Message**

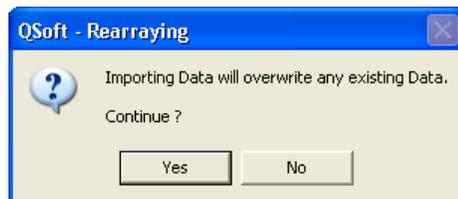
Click **Yes** to display the **Import ReArraying Data** dialog (Figure 9).



**Figure 7: Rearranging Data Selection Dialog**

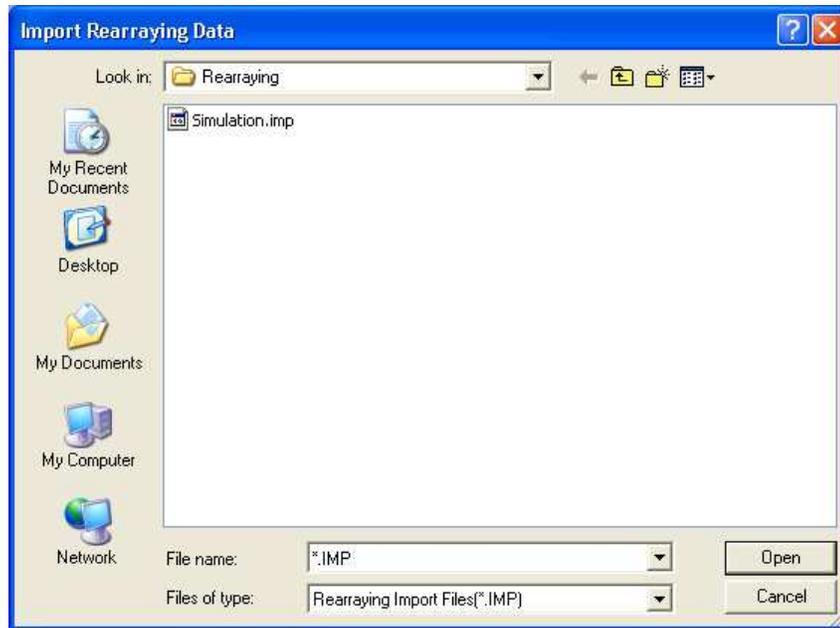
Click the Import button.

If there is rearranging data already defined for the current routine, the following screen will appear.



**Figure 8: Import Data Message**

Click **Yes** to display the **Import ReArraying Data** dialog (Figure 9).



**Figure 9: Import Rearranging Data**

This is a standard Windows **File Open** dialog and is used to locate your prepared source data file. Highlight the filename and Click on Open.

This will import all the information into the Rearranging Data Input screen. It is then possible to check the data for each plate if necessary - by clicking the plate number in the Plate Details list, every imported plate will be shown and the chosen wells are highlighted in red.

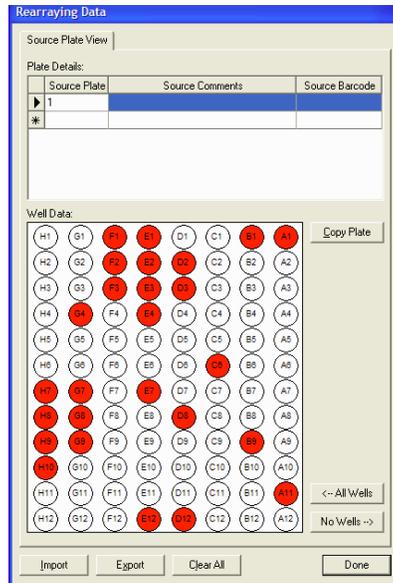
## Selecting Wells on a Plate Representation

When the Rearranging software is running, select the Source tab and click the Source Data button.

The screen shown in Figure 7 on page 15 is displayed.

**Note:** If 'Source and Destination Data' was previously imported, the existing data will need to be cleared using the 'Clear All' button in order to see the Plate Representation.

Enter the plate number (and optional plate barcode and any comments) for the plate to be rearranged from. The data entry screen for this plate will then appear on screen.



**Figure 10: Rearranging Data Plate Representation**

To view the wells better, you can expand the "Rearranging Data" window by dragging one of its corners. Individual wells can then be select/deselect by clicking on them. Selected wells are displayed with a red background.

Having chosen the required wells for the first plate, click the next blank row of the **Plate Details** section and the second plate can then be set up. Continue until all required plates have been entered.

## Recording Barcodes

Barcodes can be recorded to the log file for both source and destination. To enable barcodes to be recorded, appropriate selections must be made in the barcode options screen. The barcodes are recorded during the program run.

Barcodes can be input either directly from the keyboard or using the in-line barcode reader.

For detailed information about recording barcodes see the section headed Barcodes later in this manual.



# Rearranging Software

## Overview

Double click on the **QSoft** Rearranging Icon on the desktop, the application splash screen will appear.

A prompt will appear asking for a new routine to be created or to load a previously saved routine, any existing routines will be listed in the lower window.

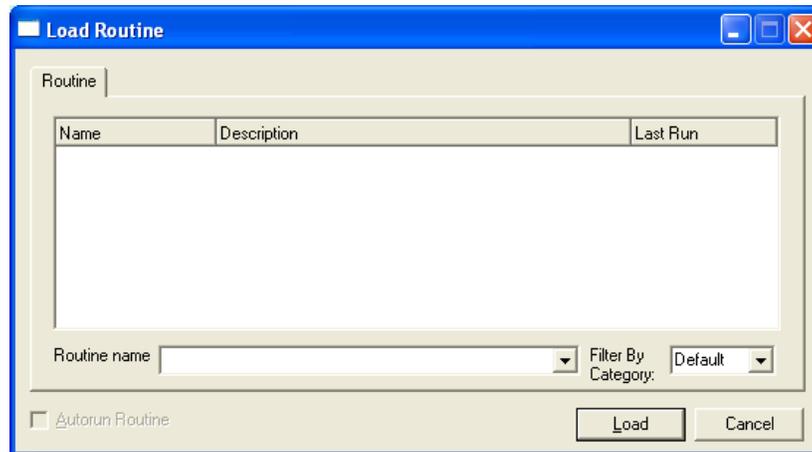


Figure 11: Routine Prompt

### Create a New Routine

Select this option then click OK. The default routine settings will be loaded, this routine can be edited then saved if required.

### Load an Existing Routine

Previously saved routines are listed here. Select this option then highlight the required routine name. Click OK to start the application with the routine settings loaded.

### Autorun Routine

If this box is checked and OK is clicked the selected routine is started automatically.

### Filter by category

Previously saved routines can be filtered by the category added when the routine was created.

## Sign On

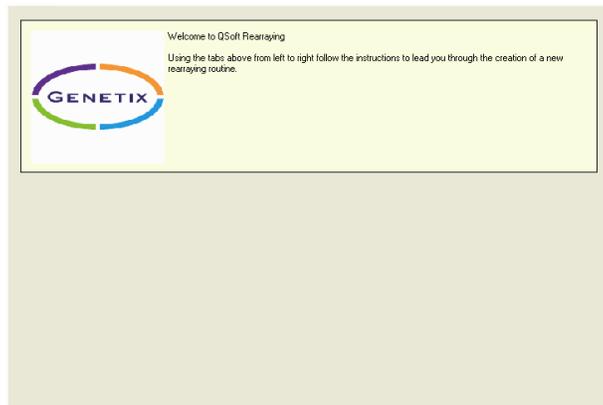
If the **Show at Start-up** box has been checked, the **Sign On** screen will be displayed.



The Sign on dialog box is titled "Sign on" and contains a "Run Details" tab. It features several input fields: "Run Number" (with a spinner box showing "1" and a "(Last Run No. shown)" label), "Operator", "Library", "Set Number", "Replica Number", and "Description". At the bottom, there are two checkboxes: "Show at Startup" (checked) and "Show for Each Run" (unchecked). "Cancel" and "OK" buttons are also present.

**Figure 12: Sign on dialog**

Complete the Sign On screen and then click OK. The Rearranging Welcome screen will be displayed.



**Figure 13: Welcome**

## The Menu and Toolbar Options

There is a reference to all of the menu items in Appendix A of this manual. More detailed information about the use of some of these can be found in the robot manual.



## Rearranging Setup

The Rearranging setup screen is split into tabbed dialogs. Each tab contains instructions to guide you through setting up a routine.

Select each tab in turn from left to right and fill in appropriate fields or select required options.

### Description

Use the text boxes here to enter information about the routine you are creating.

### Head



**Figure 14: Head**

**Head** - Gives the option to choose between the different heads available. Currently there is only 1 Head option for rearranging: 96-Pin Picking Head.

**Pin Order** - This option allows the order in which the pins in the Picking Head will fire to be chosen. The options are:

A1-H1, H2-A2... (snaking)

A1-A12, B12-B1... (snaking)

A1-H1, A2-H2, A3-H3...

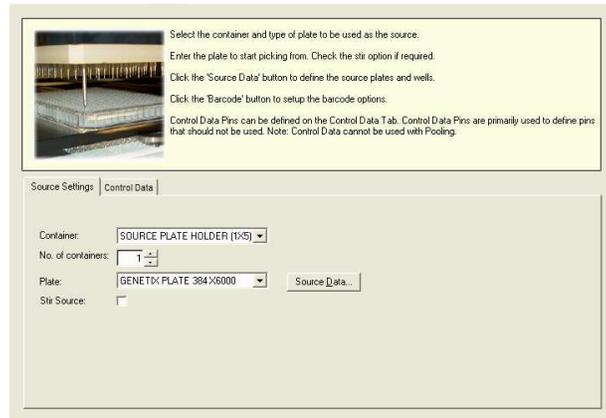
A1-A12, B1-B12, C1-C12

**First Picking Pin** - Allows the starting pin to be specified. If the number of wells to be rearranged from 1 well plate is less 96 then **First Picking Pin** shows which will be the first pin used to rearrange from the next well plate.



## Source

### Source Settings



**Figure 15: Source Settings**

**Container** - This refers to the Container that is to hold the source plates – i.e. Well plates that are to be rearrayed from. There are several options available, these vary according to which robot being used. For example...

#### QPix

Source Plate holder (1 x 3)

Source Plate holder (1 x 5)

Destination Plate holder (1 x 3)

Destination Plate holder (1 x 5)

**No. of Containers** – Refers to the number of addressable locations depending on the type of Plate selected (below).

**Plate** - This field allows the type of source plate that is to be rearrayed from to be defined. These plates will be located in the container specified above.

Note: It is essential that the correct type of microplate be selected. Severe damage to the Picking pins can result if incorrectly set.

**Stir Source** – This allows the source to be stirred prior to inking. The stir option moves back and forth then side to side within a known region as defined by the following plate properties.

- WellDiameter – Well diameter (microns)
- StirAreaPC – percentage area of the well to stir in (ensure WellDiameter is correct)
- NoStirs – how many times to stir



**Source Data** - Click on this to open the Rearranging Data dialog (see Figure 10 above).

**Copy** - Will copy the selected wells to another plate - a prompt to enter a new plate number to copy to will appear.

**<--All Wells** - Will select all wells on current plate.

**No Wells -->** - Clears all selected wells on current plate.

**Import** - Imports predefined source data from a text file.

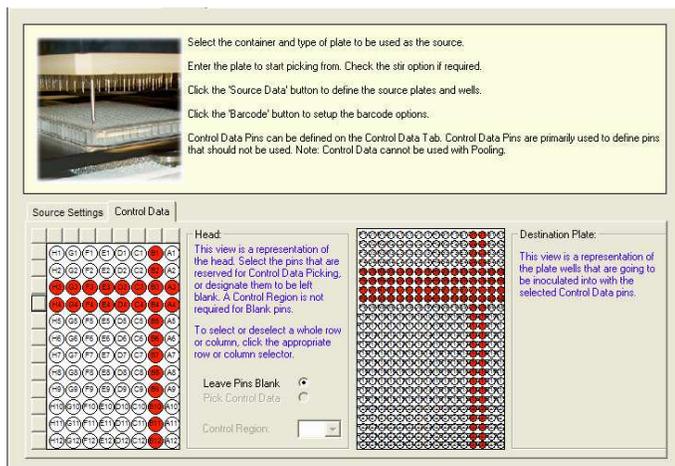
**Export** - Allows the plate configuration data in a text file to be saved. The file is automatically given an extension .exp. The file format conforms to the same standards as import data files and can be viewed using text editing software such as Windows Notepad.

**Clear All** - Clears all information related to all plates.

**Done** - Click this button when satisfied with the data.

Prior to carrying out a Rearranging run, information relating to the wells that are to be rearranged must be entered. The methods for doing this are described in detail in the section headed "Defining Well Data" on page 12.

## Control Data



**Figure 16: Control Data**

Please Note: Control data cannot be used when destination well data has been defined (for example in a Source and Destination data import file).

### Head

This diagram is a representation of the head with pin locations labelled. Click to select/deselect pins to be left blank.

Note: A complete row or column can be selected by clicking the relevant selector button.

### Destination Plate

This diagram represents the selected destination plate, as pins to leave blank are selected, the corresponding plate wells are highlighted in red.

### Leave Pins Blank

This is the only available option in Rearranging.



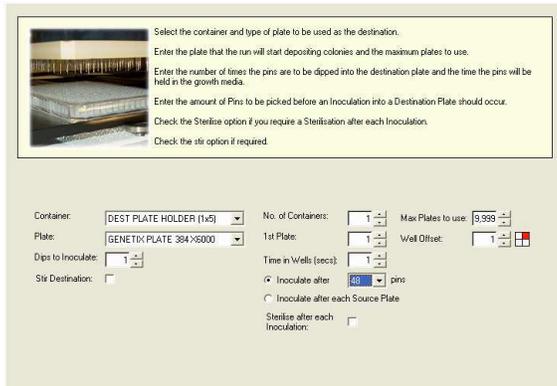
### Pick Control Data

This is currently not available in Rearranging.

### Control Region

Works in conjunction with the **Pick Control Data** option.

### Destination



**Figure 17: Destination.**

**Container** - This refers to the Container that is to hold the destination plates – i.e. Well plates that are to be picked into. There are several options available. For example...

**QPix**

- Destination Plate holder (1 x 3)
- Destination Plate holder (1 x 5)
- Source Plate holder (1 x 3)
- Source Plate holder (1 x 5)

**No. of Containers** - Refers to the number of addressable locations depending on the type of Plate selected (below).

**Plate** - This field allows the type of destination plate that is to be picked into to be selected. These plates will be located in the container specified above. There are several options available, these will vary depending on which robot you are using.

Note: It is essential that the correct type of microplate be selected. Severe damage to the Picking pins can result if incorrectly set.

**1st Plate** - This determines the well plate into which the picking head will start depositing colonies. It is normally set at 1 but can be set at any value up to the maximum number of plates.

**Max Plates to use** - This specifies the maximum number of well plates available. If this value is set to a higher number of plates than the total Destination Holder capacity a **pop-up** screen will appear and prompt for "Change WellPlates" before continuing.

**Dips to Inoculate** - In order to control the inoculation of the Destination plates, you can choose the number of dips that the Picking Head will make into the well plates can be chosen. This setting would normally be in the range of 1 to 3.

**Time in Wells (secs)** - Varies the time that the Picking pins are held in the Growth media in the well plates in seconds. This parameter is used in conjunction with the number of Dips to Inoculate.

**Well Offset** – The well offset corresponds to one of the 4 inoculation offsets (wells A1, A2, B1



and B2) when using 384 well plates. This is not applicable when using 96 well plates. Default value - 1.

**Stir** – This allows the source to be stirred prior to inking. The stir option moves back and forth then side to side within a known region as defined by the following plate properties.

- WellDiameter – Well diameter (microns)
- StirAreaPC – percentage area of the well to stir in (ensure WellDiameter is correct)
- NoStirs – how many times to stir

**Inoculate After nn pins** – Will pick colonies sequentially from all available plates and will only deposit after the set number of pins have been used.

**Inoculate After each Source Plate** – Will pick colonies sequentially from each plate and will deposit *either* when all pins have been used *or* when all colonies from the current plate have been picked.

**Sterilise After Each Inoculation** – Check this box if you want to sterilise the head after every deposit.

## Sterilize



Figure 18: Sterilize

**Bath Cycles** – The number of cycles in the wash bath. Usually set to 3 or 4.

**Dry Time** – The length of time the pins are dried in the dryer. If the number of bath cycles is set to zero this box is greyed out.

**Wait After (drying)** – A time delay (in milliseconds) can be introduced to allow the head to cool after drying.

If the robot uses a halogen dryer, QSoft automatically adds to this wait time in order to allow enough time for the pins to cool properly.

This additional wait time is based on the following calculation:

$$3 \text{ Seconds} + (1.5 \times \text{Dry Time}).$$

Thus even if Wait Time was set to 0 and Dry Time was set to 5000 ms, the head would remain in the dryer for 10.5 seconds after drying.



**Wash Solutions** - A single wash bath containing 80% ethanol will ensure sterility when handling *E. coli*. However, some organisms (e.g. yeast) form 'sticky' colonies that can build up on the pins and others are particularly robust (e.g. spore-forming organisms). To ensure sterility in these cases it is recommended to use all three wash baths with the following solutions:

Bath 1 – 1% sodium hypochlorite

Bath 2 – deionised water

Bath 3 – 80% ethanol

The washes should be performed in the order 1% sodium hypochlorite, water then 80% ethanol followed by drying and the wait after drying.

**Note:** Use of sodium hypochlorite at concentrations greater than 1% may cause damage to the instrument.

## Barcodes

This area allows you to set the Barcoding options for the Run.  
A Separate License may be required.  
Select either Manual or Automatic Barcoding.

780747521256

Source | Destination

Enable Logging:

Input Method:  
 Manual  
 Automatic

Behaviour Options:  
Barcode not Found:  
 Manual  
 Automatic

Validate:   
Barcode not Validated:  
 Manual  
 Automatic

Figure 19: Barcodes

**Important Note:**

The thickness of some barcode labels can affect the fit of the lid so that it becomes too tight to remove the lid. (Only applies to robots that have Stacker Units and/or Lid Lift mechanisms).

This problem does not arise so much with Genetix plates as they have been designed with special ribs on the lid which do not obstruct the barcode label.

To enable barcodes to be recorded, the appropriate options must be set in the Barcodes setup dialog. Select the Enable Logging option on the Source and/or Destination tabs of the Barcodes dialog.

Barcodes can be input either directly from the keyboard or using the in-line barcode reader.



## Enable Logging

Check this box to enable barcode reading.

QSoft will automatically generate a unique key in the barcode field for any plate that is used during a run and the logs will display these keys in the barcode field. The 'Enable Logging' barcode reading options will overwrite this key. The keys are generated based on the current date and time and in the format UID-YYMMDDHHNNSSss-X where ss is milliseconds and YY is a two character ascii representation of the year.

**Input Method** – Choose Manual if scanning barcodes with a hand held barcode reader is required or to input barcodes at the keyboard. Automatic barcode reading is possible if the robot has a barcode reader installed. If automatic barcode reading is required contact Genetix Ltd for a licence.

**The following options are enabled when Automatic barcode reading is selected.**

## Behavior Options

This section determines how QSoft will behave in the specified circumstances.

### Barcode Not Found

Set required behaviour if a barcode is not present.

- Manual will produce a prompt and will wait for a barcode to be input (either via the keyboard or using a hand held barcode scanner).
- Automatic causes QSoft to generate a unique barcode based on the system date and time.

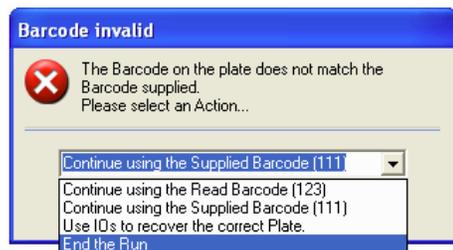
### Validate

Check this box to enable barcode validation.

### Barcode Not Validated

- Manual will pause allowing the order of the plates being processed to be verified before continuing the run.
- Automatic, in the event of an invalid barcode the run will continue automatically replacing the expected barcode with the read barcode. These actions will be recorded in the log.

**Note:** If barcode validation fails a set of options as shown in Figure 26 below will be displayed.



**Figure 20: Invalid Barcode Dialog**

**Add** – Creates a new row for entry of the next barcode

**Remove** – Deletes the highlighted entry

**Import** – Prompts for the name of the text file that contains the validation barcodes

**Export** – Creates a text file of the current list

**Clear All** – Removes all entries from the list

### Disable Barcoding for source or destination

If Datatracking is selected, Barcoding is automatically enabled. Barcoding for a given container can be disabled – for example if barcoding the source bioassay trays is not required – by

changing the **BarcodeReaderType** property this can be achieved. Do this as follows:

- Highlight the container in the **Hardware Configuration** tab
- Click the **Properties** button
- Select **BarcodeReaderType** and click the **Edit** button
- Enter **None-AutoGenerate** and click **OK**

## Start

Having set all of the variables, save the routine by clicking the Save or Save As button on the toolbar.



**Figure 21: Start**

### Enable Data Tracking

Selecting this option will generate a Data Tracking XML file, which can then be imported into the Data Tracking database.

This is only applicable if barcoding options have been set for both source and destination plates.

The data tracking file is stored in the following location:

**C:\Program Files\Genetix\Logs\**

To view the xml file, double-click the filename in Windows Explorer.

For more information on Data Tracking, please refer to the **QSoft Data Tracking** manual.

**Start in Slow Motion** – The robot will run at a slower speed for diagnostic purposes.

**Run** – Click this button to start the routine.



## Running a ReArraying Routine

Once all the parameters have been set and plates have been loaded, the Rearraying Routine can begin.

Click the **Run** button at the bottom of the setup screen.

The following screen appears:



**Figure 22: Prepare Script**

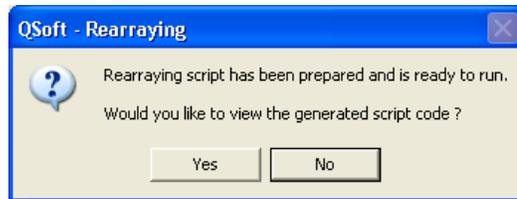
Click **Yes** to continue **No** to abort.

The following screen will appear if continuing, which displays the total number of wells to rearray.



**Figure 23: Wells Found**

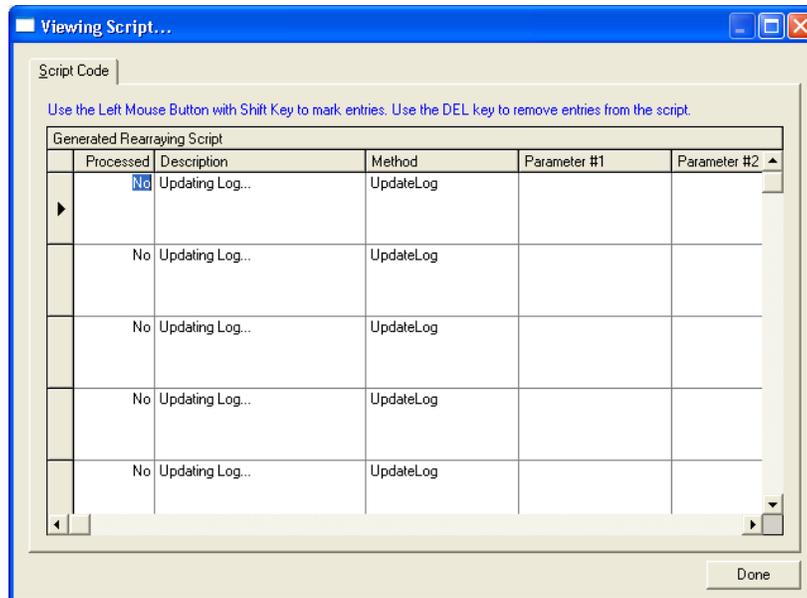
Click OK. And the following screen will appear.



**Figure 24: View Script**

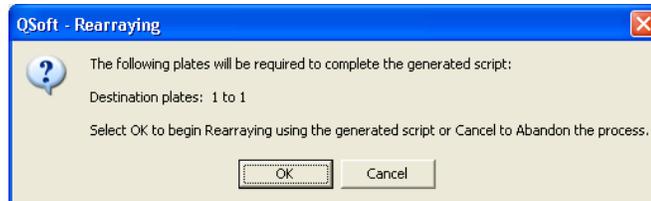
To view the script code click on **Yes** otherwise click **No**.





**Figure 25: Example Script Code**

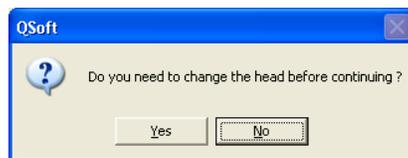
Click on Done to continue, the following screen will appear:



**Figure 26: Plates Required**

This screen will indicate how many destination plates are required for the rearraying run. Ensure that these have been loaded. Click OK to carry on.

The screen will now ask to ensure that the correct head has been fitted to the robot.

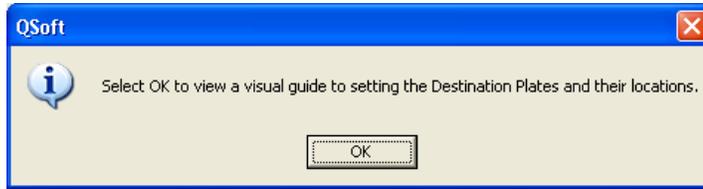


**Figure 27: Change Head**

**Remember that for a rearraying run the Picking Head must be fitted.**

Once this has been checked click on **Yes** to continue. The following screen will now appear.

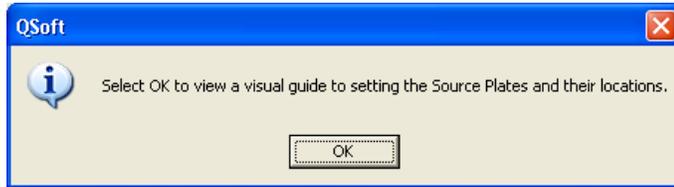




**Figure 28: Destination Plate Locations**

This prompt is for the destination plate holders to be fitted on the bed of the robot. When you click **OK** you will be presented with a visual display that indicates the positioning of the destination well plates in the holders.

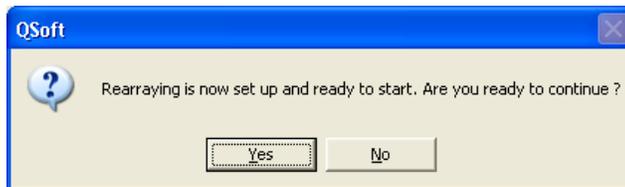
Click **Done** to continue, the following screen appears:



**Figure 29: Source Plate Locations**

Click **OK** and the screen will show the required source well plates.

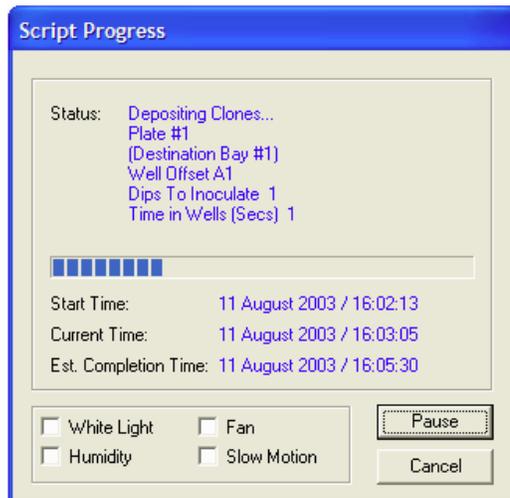
Click **Done**.



**Figure 30: Set Up and Ready**

This screen displays that the ReArranging is now ready to start. To commence the run click **Yes**. To abort click **No**.

During the run the progress screen (shown below) will be displayed. This screen displays the status of the robot, the start time of the run, current time and expected time of completion.



**Figure 31: Script Progress**

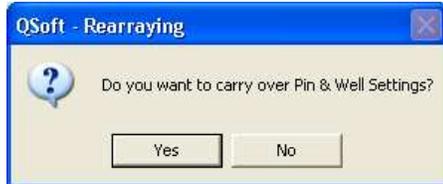


If at any time the run needs to be stopped, the **Pause** button can be clicked.  
When the run has finished, the following screen will appear.



**Figure 32: Script Complete**

Clicking **OK** will generate the following screen.



**Figure 33: Carry Over Pin and Well Settings**

The last well and pin settings are stored, so that more data can be loaded to carry out a further run, utilising any pins on the head that have not been used and any destination plates that have not been filled.

As you become more familiar with the robot some of the above screens need not be shown. Under **Additional Options** in the ReArraying setup screen, there is a command button – **Script Options**, allows which screens will appear and which will not to be chosen.

Click the **Exit** button when ReArraying is finished.



# Biology Guide

## Preparing Media

### Luria-Bertani medium (LB) – per litre

To 1 litre of de-ionised H<sub>2</sub>O add 25g of pre-prepared LB (Sigma, Gibco, LAB3). LB can also be made by adding (per litre) 10g tryptone, 5-10g yeast extract, 5g NaCl (pH 7.2) and stir on a magnetic stirrer until the powder has dissolved. Sterilise by autoclaving at 121°C for 15 minutes.

### LB + 8% Glycerol – per litre

As above but replace 80 ml of water with glycerol (80 ml glycerol + 920 ml deionised H<sub>2</sub>O). Sterilise by autoclaving at 121°C for 15 minutes.

### LB Agar – per litre

To 1 litre of de-ionised H<sub>2</sub>O add appropriate amount of pre-prepared LB agar (Sigma, Gibco) and stir until the powder has dissolved. If making your own LB add 16g of agar per litre of LB. Sterilise by autoclaving at 121°C for 15 minutes.

## Antibiotic Preparation

### Ampicillin

**Stock solution:** Dissolve 1g of ampicillin in 20ml of sterile distilled water. Filter sterilise using a 0.2 µm syringe filter and dispense 1ml aliquots into 1.5ml Eppendorf tubes. Store at -20°C. (50mg/ml stock solution).

**Working solution:** Add 1ml of stock solution per litre of medium.

### Chloramphenicol

**Stock solution:** Dissolve 1.25g of chloramphenicol in 100ml of ethanol. Store in 1.5ml Eppendorf tubes at -20°C. (12.5 mg/ml stock solution).

**Working solution:** Add 1ml of stock solution per litre of medium.

### Kanamycin

**Stock solution:** Dissolve 1g of kanamycin in 20ml of sterile distilled water. Sterile filter and store in 1.5ml Eppendorf tubes at -20°C. (50 mg/ml stock solution)

**Working solution:** Add 1ml of stock solution per litre of medium.



# Glossary of Terms

---

## Arrayed

Distribution of colonies or samples into known positions from 96 or 384 WellPlates

## Base Class

Blueprint for the properties of an object

## Bioassay Tray (QTray)

22x22 cm clear plastic tray from which colonies/phage are picked

## Bioassay Tray Holder

Perspex holder that drops into the robot bed for holding two Bioassay trays in place whilst carrying out a Picking routine

## Compressing

Converting 4 x 96 well plates into 1 x 384 well plate etc.

## Datum Point

A series of X, Y, Z co-ordinates that define a set position on the Robot bed

## Destination Plate Holders

Holders for microplates located on the bed of the robot. There are 3 destination plate holders on QPix. They are used for replicating, rearraying and picking.

## DMF

Dimethyl formamide

## Expanding

Converting 1 x 384 well plate into 4 x 96 well plates etc.

## I/O

Inputs / Outputs

## IPTG

Isopropyl-thio- $\beta$ -D-galactoside

## LB

Luria-Bertani Medium



**Phage**

Bacteriophage

**QSoft.DLL**

ActiveX software component housing all the functionality of robot software

**Rearranging**

Redistribution of selected colonies into new plates performed with picking head

**Replicating**

To copy, compress or expand 96 or 384 well plates

**Script**

Listing of all moves needed to complete a routine

**SDS**

Sodium Dodecyl Sulphate

**SSC**

Sodium Chloride/Sodium Citrate buffer

**X Drive**

Axis running from back to front of the QBot bed or right to left on QPix

**Y Drive**

Axis running from left to right across the QBot bed or back to front on QPix

**Z Drive**

Axis running vertically on the Robot bed



# Appendix A

## Description of Toolbar and Menu Items

	Add Sample to Rearranging Checkout	Add the highlighted sample(s) from the currently selected plate to the Rearranging checkout.	Data Tracking
	Align	Align camera to pin	Picking Excision Excellerate Rearranging
	Analysis path	Allows you to select the Gel Analysis program to be used	Excision
	Calibrate	Calibrate the camera	Picking Excision Excellerate Rearranging
	Change head	Moves the actuator into an accessible position to allow you to change the head	All
	Chiller	Turns the chiller on or off	Robot Specific
	Clear Rearranging Checkout	Will clear all samples from the rearranging checkout.	Data Tracking
	Configuration	Displays the robot configuration dialog which gives you access to all the hardware settings, datum point settings and database facilities	All
	Configure Messaging Server	Messaging Server provides a means for you to remotely monitor your robot. For example: if a robot run is interrupted for any reason, one or more contacts can be notified thus eliminating the need for constant supervision of your robot.  Click this button to set messaging server preferences. For detailed information about setting up Messaging Server see Appendix A of this manual.	All
	Create Rearranging file	Write the contents of the Rearranging checkout to a text file.	Data Tracking
	Diagnostics	Displays an animated representation of the robot bed layout for diagnostics purposes or for use when running the software in simulation mode. Click <b>Stay on Top</b> to keep the dialog visible while the application is running.	All
	Exit	Exits the application	All
	Fan	Turns the fan on or off	Robot Specific
	Humidity	Turns the humidity system on or off	Robot Specific
	Import Custom Properties	Import properties to use in searches to identify samples of interest	Data Tracking
	Import Process file	Import a Data Tracking process file	Data Tracking



	IO's	Displays the robot Inputs and Outputs to allow you to manually control hardware components	All
	Logs	Displays the QSoft Logs dialog which gives you access to the Text, XML and DataTracking log files. Highlight the log file name and click Open to view it.	All
	Park head	Parks the head in the wash bath	All
	Pin fire test	Displays a dialog whilst continuously firing all pins in the head. The purpose of the Pin Firing Test is to check, by sound, that each pin is firing correctly. <b>Continuous Test</b> – When checked will continuously fire pins until it is either unchecked or the Cancel button is selected. <b>Control Valve</b> – When checked, slows down the pin retraction. This is useful for first time use or for demonstration purposes to allow you to observe the mechanical actions of the pins. <b>Cancel</b> – Ends the Pin Firing test.	Picking Excision Excellerate Rearranging
	Purge	Purges an Excision head to clean the pins	Excision Excellerate
	Remove Sample from Rearranging Checkout	This will remove selected samples from the rearranging checkout.	Data Tracking
	Reset toolbars	You can rearrange the buttons on the toolbar if necessary, this option will set them back to the default order.	All
	Routine close	Closes the current routine	All
	Routine delete	Deletes the specified routine	All
	Routine Export	Permits you to save routines as XML files in a user-defined location. This enables the transfer of routines between robots.	All
	Routine Import	Allows externally created routines to be used.	All
	Routine new	Create a new routine	All
	Routine open	Opens a previously saved routine	All
	Save	Saves the current routine	All
	Save as ...	Saves the current routine with a new name	All
	Script options	Allows you to select which script options to display during the application run	All
	Select Database	Displays the login dialog to allow you to connect to a different database or login as a different user.	Data Tracking
	Set pin height	Provides the facility for setting the picking height	Robot Specific



	Sign on	Displays the Sign on dialog to allow you to enter specific information about the current run	All
	Switch User	<p>Displays the Login prompt. There are currently 3 levels of user, permissions are as follows:</p> <p><b>Operator</b>                      Load routines                      Run routines</p> <p><b>Creator</b>                      Create routines                      Load routines                      Run routines                      Save routines                      Save other user's routines with a new name</p> <p><b>Admin</b>                      No restrictions on use</p>	All
	System info	Displays information about your computer system	All
	Test Image	Displays the test image window so that you can set criteria for picking colonies	Robot Specific
	UV light	Turns the UV light on or off	Robot Specific
	Vacuum	Turns the vacuum on or off	Robot Specific
	White light	Turns the white light on or off	All



# Appendix B

---

## GTX Messaging Server

Messaging Server provides a means for you to remotely monitor your robot. So that if a robot run is interrupted for any reason, one or more contacts can be notified thus eliminating the need for constant supervision of your robot.

In order to use Messaging Server you need to allocate one or more contacts and ensure that at least one of those contacts is enabled before the run starts.

Once QSoft has started to process the script, Messaging Server will notify the contact(s) whenever the software requires a response. For example any of the standard message boxes that appear during the run and also "Timeout" messages that are displayed when a problem occurs.

The messaging server currently supports dial-up (without voice) and email notification.

### Dial-Up Settings

The dial up facility is implemented through the use of a Hayes compatible modem installed in the robot PC. A phone line is also required. Upon receiving a message, the modem will dial the appropriate number, and ring it for a predefined period of time. If the receiving contact's number is a mobile phone, a phone book entry can be made for the robot. Then when the robot rings the number, the caller-id facility will display to the operator which number is called.

### Email Settings

The email facility requires that a MAPI compliant mail application (such as Microsoft Outlook<sup>®</sup> Express) has been configured on the robot PC.

The steps for configuring Outlook Express are as follows:

- Set up the email account that is required to send mail with. If unsure how to do this then look in the Outlook Express help or contact the system administrator.
- Change the security settings to allow the messaging server to automatically send an email unattended:
  - Open the Tools menu, then select Options



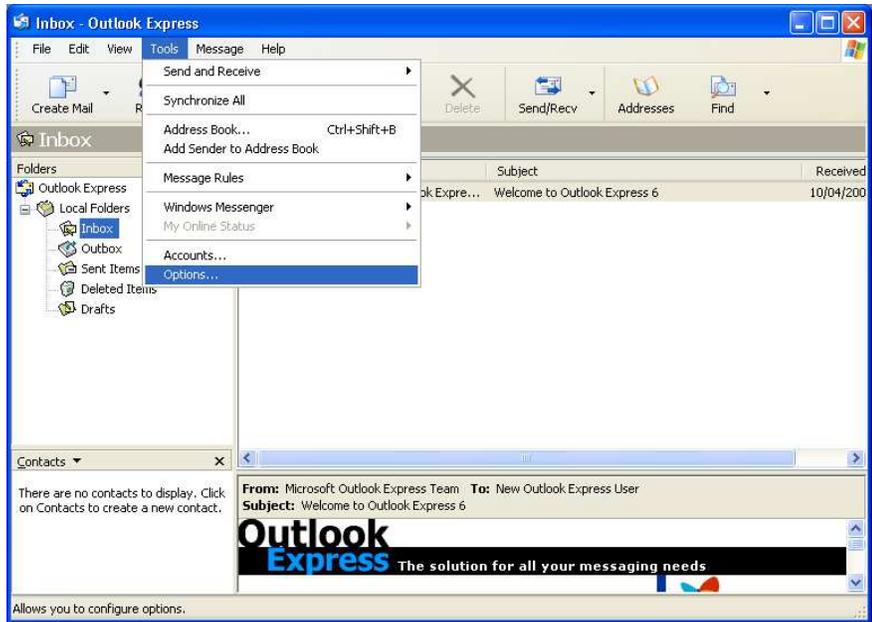


Figure 34: Outlook Express Tools Menu

Uncheck the option that says:

**'Warn me when other applications try to send mail as me'**



Figure 35: Outlook Express Security Settings

Click **OK**



## Configure Messaging Server

Click the **Configure Messaging Server** button on the toolbar.

The following screen will be displayed

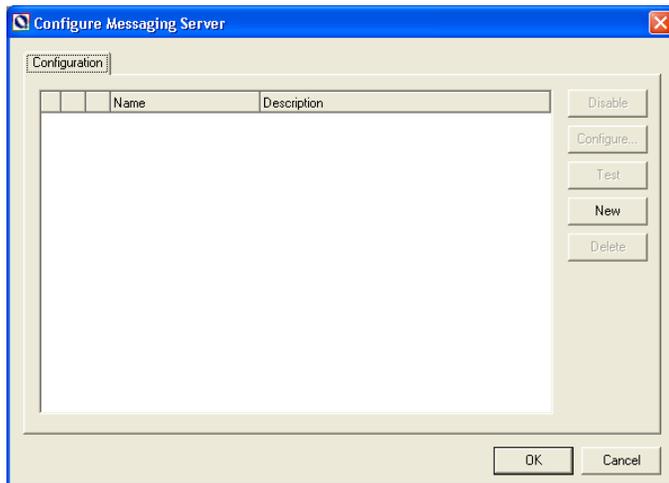


Figure 36: Messaging Server Configuration

Click the **New** button to add a new contact.

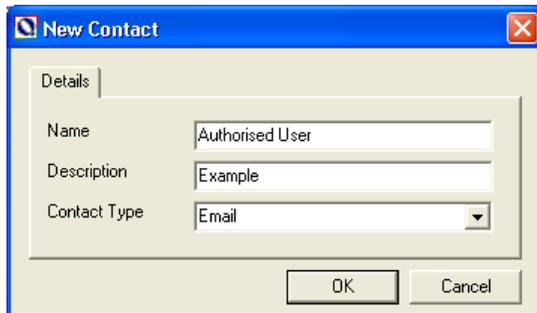


Figure 37: New Contact

All the parameters here are required.

### Name

Type in a name for the contact

### Description

Type in any descriptive text.

## Contact Type

Select the required option from the drop-down list

Click OK when you have entered the required details. The following dialog is displayed.

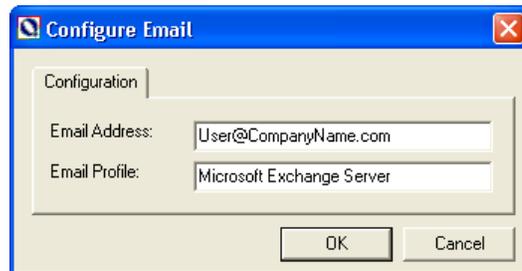


Figure 38: Configure Email

## Email Address

Enter the contact's valid email address.

## Email Profile

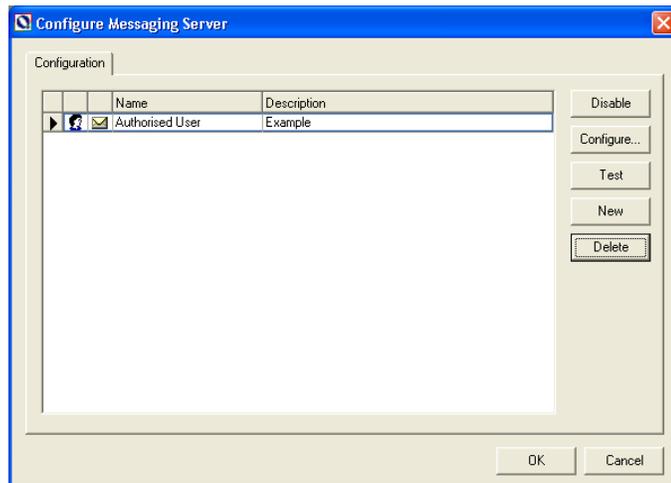
If there is no specific profile set up (such as Microsoft Exchange Server) enter the name of the registered user of the copy of Windows that is installed on the robot PC. To identify the registered user, locate System Properties in Windows Control Panel (see below).



Figure 39: Windows System Properties

Click **OK** to store these settings.

The Configuration screen will show the details of the contact you have just added.



**Figure 40: Contact Details**

### Disable

This allows the selected (highlighted) contact to be disabled. The button's state will change to say **Enable** when a contact with **Disabled** status is highlighted. Operators can be enabled or disabled on a run-by-run basis.

### Configure

Allows the configuration settings of the selected contact to be changed.

### Test

Use this to check that the email and/or dial-up settings work as expected.

### New

Allows you add a new contact and configure their settings.

### Delete

Will delete the highlighted contact.

## Set the Delay Time

By default the delay time is set in QSoft to 10 minutes. This is the time that will elapse before either a mail message is sent or the dial-up will be activated. If necessary, you can change the delay time as follows:

Click the **Configuration** button on the toolbar.

In the **Hardware** tab select the robot name.

Select the property named **MessagingSendingDelayMinutes**.

Click **Edit** to change the value.

# Index

<b>A</b>		MicroArraying .....	27
Actuator .....	35	Microns .....	21, 24
Addressable locations.....	10, 21, 23	Monitor .....	35
Airline.....	7	<b>O</b>	
Antibiotic Preparation.....	32	Options .....	26
<b>B</b>		<b>P</b>	
Barcode .....	13, 16, 17, 25, 26	Picking.....	7, 20, 29
Barcode reader .....	17, 25, 26	Picking head.....	7, 20, 23, 29
Barcodes .....	17, 25, 26	Pin .....	8, 20, 31, 36
Base plate.....	7	Pin(s).....	8, 20, 36
Bath .....	7, 8, 24	Pipette .....	7, 8
Bed .....	7, 10, 30	Plate holder .....	21, 23
<b>C</b>		Preparing Media .....	32
Cleaning .....	7	Progress screen .....	30
Colonies.....	8	Properties .....	21, 24
Container .....	7, 21, 23	<b>Q</b>	
<b>D</b>		QArray .....	27
Data tracking .....	27	QPix .....	6, 7, 8, 9, 10, 21, 23
Description.....	13	QSoft .....	18, 26, 27
Destination.....	25	<b>R</b>	
Destination plate holder.....	30	ReArray .....	20, 28
Distilled water .....	7	Rearranging	
DNA .....	8	Import File Format.....	13
<b>E</b>		Source Data .....	22
Export .....	26	What is Rearranging?.....	6
<b>F</b>		Remove .....	7, 8, 26
File extension .....	13	Replacing Pins .....	8
Fitting the head.....	7	Replicating	
<b>G</b>		Preparation .....	7
Gel Analysis.....	35	Run .....	27
Genetix .....	27	Run.....	12, 17, 22, 26, 27, 28, 29, 30, 31, 35
<b>H</b>		<b>S</b>	
Head .....	7, 8, 10, 20, 24, 29, 31	Scanner .....	26
Hotel .....	21, 23	Screen .....	18, 19, 25, 28
<b>I</b>		Setup.....	18, 25
Import .....	26	Setup screen .....	25
<b>L</b>		Sign on .....	19
Licence .....	26	Sonic cleansing .....	7
Log barcodes .....	25	Sonicate .....	7
<b>M</b>		Sonicator .....	7
Maintenance .....	7	Source plates .....	6, 8, 9, 21
Media .....	23	Stir source .....	21, 24
Menu.....	19	<b>U</b>	
Menu Options .....	19	UV sterilise .....	7
		<b>W</b>	
		Wash bath .....	7, 8, 24
		Wellplate.....	20
		Workbed.....	7



# Contact Details

---

## Corporate Headquarters

### Genetix Ltd

Queensway, New Milton  
Hampshire BH25 5NN, UK

Tel: +44 (0) 1425 624 600

Fax: +44 (0) 1425 624 700

Visit [www.genetix.com](http://www.genetix.com) for contact details of your nearest Genetix sales and support.

### Trademarks

ClonePix™, CloneSelect™, CellReporter™, HalfBD®, 'NRich®, SlidePath™, Data Arena™ and Image Arena™ are trademarks of Genetix Ltd., UK

CytoVision®, Ariol®, CytoNet® and CytoTalk® are registered trademarks of Genetix Corp

All third party trademarks are the property of their respective owners.

### Copyright © 2010 by Genetix Ltd

All rights reserved. No part of this publication may be reproduced, stored in a retrieval system, or transmitted, in any form by any means, electronic, mechanical, by photocopying, recording, or otherwise, without the prior written permission of Genetix Ltd.

Information furnished by Genetix Ltd is believed to be accurate and reliable; however, no responsibility is assumed by Genetix Ltd, for its use; nor for any infringements of patents or other rights of third parties which may result from its use. No license is granted by implication or otherwise under any patent rights of Genetix Ltd.

Revised January 2010

