

Screen Acquisition Option is not present

Abstract

Reasons and solutions for the absence of the Screen Acquisition option from the Apps menu.

Instructions

Symptoms

The Screen Acquisition option does not appear under the Apps menu.

Solution

There are three reasons why Screen Acquisition would not be listed in the Apps menu:

1. A stage MetaDevice is not open.
2. The dropin **htacquir** is not being loaded.
3. No camera / acquisition device is being used.

Diagnosing and solving the problem

The first possibility may be diagnosed and solved within the Discovery-1 application.

1. Choose *Install and Configure Devices* from the Devices menu. The *Install and Configure Devices* dialog will appear.
2. The third button down should say "Close Stage MetaDevice" and a metadvice will be listed. If the button says "Open Stage MetaDevice", then push the button to open a stage metadvice. If more than one stage metadvice is available you will be asked to choose one. Consult your sales rep. or dealer for more information on which one to choose.

The second possibility requires the use of the Discovery-1 software administrator (called the Meta Imaging Series Administrator) application.

1. Start the Meta Imaging Series Administrator program.
2. Determine whether you are in Single User Mode or Multiple User Mode by looking at the application window title bar. If you are in Multiple User Mode skip to step 6.
3. Select the application you are using from the list of Group names and choose the *Configure Drop-ins* button. The *Configure Drop-ins* dialog will appear.
4. The **htacquir** drop-in should be listed in the Installed Drop-ins list (right column). If **htacquir** is not in the installed list make sure that the Category: Special is checked, select the drop-in from the Available Drop-ins list (left column) and hit the *Add* button.
5. Choose *Close* to return to the main Meta Imaging Series Administrator window. Choose *OK* to exit the Meta Imaging Series Administrator application. (Do not follow steps 6-9 as they apply to the multi-user mode).

Document ID
D20002

Product
Discovery-1

Platform
Windows 95/98
Windows NT
Windows 2000

Created
02-Jan-2002

Last Reviewed
14-Jul-2003

Keywords: hcs discovery-1 screen acquisition dialog missing

Issue Type: geninfo admn

Discovery-1 Online Support

6. Select the application or the user within the application from the list of Group names and Users. Choose the *Edit Group* or *Edit User* button. The Edit group or Edit User dialog will appear.
7. Choose the *Configure Drop-ins* button. The Configure Drop-ins dialog will appear.
8. The **htacquir** drop-in should be listed in the Installed Drop-ins list (right column). If htacquir is not in the installed list make sure that the Category: Special is checked, select the drop-in from the Available Drop-ins list (left column) and hit the *Add* button.
9. Choose *Close* to return to the Edit group or Edit User dialog. If you were editing a group you will be asked whether the drop-ins for all users should be updated - choose Yes. Choose OK to return to the main Meta Imaging Series Administrator window. Choose OK to exit the Meta Imaging Series Administrator application.

You can check this by clicking the camera icon in the standard software toolbar. If an image appears, the camera is loaded and working properly. If an image does not appear, or the camera icon is unavailable, check that the camera is on and that the cables are connected, then restart your software. If this does not help, use of the Meta Imaging Series Administrator application.

1. Start the Meta Imaging Series Administrator program.
2. Find the Group you are using from the list of Group names and note the Hardware Setting assigned to the group.
3. Choose the *Configure Hardware* button. The *Configure Hardware* dialog will appear.
4. Select the *Hardware Setting* you are using and choose the *Configure Acquisition* button. The *Configure Acquisition* dialog will appear.
5. Confirm that the appropriate driver for your camera is installed. If it is not, contact Universal Imaging or your local representative for assistance in adding and configuring the appropriate driver.
6. Choose *OK* to return to the *Configure Hardware* dialog.

Choose *OK* to return to the main Meta Imaging Series Administrator window. Choose *OK* to exit the Meta Imaging Series Administrator application.