

Release Notes
Molecular Analyst / PC Molecular Imager
PC version 1.5

INSTALLATION INSTRUCTIONS FOR THE MOLECULAR ANALYST VERSION 1.5

New Item for Windows 95 Users:

After installation, the Molecular Analyst program group contains an additional icon, "Install HASP Hardware Key". Window 95 users must double click on the "Install HASP Hardware Key" icon to install necessary system drivers to properly access the hardware key.

In Windows 3.1 or Windows 95:

1. Turn on your computer and start Microsoft Windows.
2. Insert the "Molecular Analyst Setup" diskette in the A: drive.
3. In Windows 3.1, select "Run" from the "File" menu in Windows' Program Manager.
In Windows 95, select "Run" from the start menu.
4. Type a:setup and press the OK button. The Molecular Analyst Setup program will display its window.

The setup program is self explanatory. It will examine your video system software driver to be sure that it is set to 256 color mode and will then determine whether there is enough free space (about 4 Mbytes) on the hard disk. The "Preload SCSI Drivers" checkbox should be checked during setup by most users. This will insure that SCSI drivers are loaded into conventional memory on the PC. If you later get the error message "Not enough conventional memory to scan" you must rerun setup and set this option.

The setup program will, by default, install the software in the C:\MA directory. Any other directory may be specified instead. If it does not exist, it will be created.

If a previous version of the Molecular Analyst software is already installed in the same "C:\MA" directory, it will be overwritten. User created data will not be disturbed. Sample images included with the software will be updated.

When the installation is complete:

In Windows 3.1, activate the Windows Program Manager application.

In Windows 95, the Molecular Analyst program in the start menu will be open on the desktop.

You will see a new program group called the Molecular Analyst. Within this group are three icons: Molecular Analyst , Read Me Imager and Read Me Copy Protection.

Double click on the "Read Me Imager" icon to review the document you are reading. Double click on the "Read Me Copy Protection" icon for information on the Copy Protection key installation.

ASPI Installation Instructions

For both Windows 3.1 and Windows 95, your CONFIG.SYS file must load the ASPI device manager in order for the Molecular Analyst software to communicate with the Molecular Imager. This is true in Windows 95 even though Windows may be loading its own protected-mode driver for the SCSI adapter.

This ASPI driver is usually called ASPI4DOS.SYS.

Important note: The Molecular Analyst does not support the Adaptec 2940 SCSI PCI card.

The recommended SCSI adapters for the Molecular Analyst are the Adaptec 1540B/42B, 1540C/42C or 1540CF/42CF.

If you are running Windows 3.1 and you are already using a CD-ROM or optical disk drive on the same SCSI bus to be used for the Molecular Imager, the CONFIG.SYS most likely already has the proper entry.

In Windows 3.1, if the Imager will be the only device on this SCSI card, or in all cases for Windows 95, the ASPI driver must be installed by using the Adaptec EZSCSI diskette included with your SCSI card.

GENERAL INFORMATION

RECOMMENDED 'autoexec.bat' FILE ENTRIES for Windows 3.1

```
REM DOS Version
VER
REM Set DOS prompt to <DRIVE:DIR>
PROMPT $P$G
REM Load disk cache for drive C with a size of 4096k bytes.
C:\WINDOWS\smartdrv.exe C+ 4096 4096
REM Set path variable to DOS and Windows directories.
PATH=C:\WINDOWS;C:\DOS
REM Set TEMP variable to the correct directory for temporary file.
REM Make sure directory exists.
SET TEMP=C:\TEMP
REM Check the integrity of your hard disk
CHKDSK /F
REM Launch Windows
WIN
```

RECOMMENDED "config.sys" FILE ENTRIES for Windows 3.1

```
REM Load High memory manager device driver
DEVICE=C:\WINDOWS\himem.sys
REM Load DOS in high memory.
DOS=HIGH
REM Set maximum files to 30
FILES=30
REM Set maximum file buffers to 20
BUFFERS=20
REM Set system stacks to 9,256
STACKS=9,256
REM Install SCSI Manager
DEVICE=aspi4dos.sys /D
```

RECOMMENDED 'autoexec.bat' FILE ENTRIES for Windows 95

```
REM Set TEMP variable to the correct directory for temporary file.  
REM Make sure directory exists.  
SET TEMP=C:\TEMP
```

RECOMMENDED "config.sys" FILE ENTRIES for Windows 95

```
REM Install SCSI Manager  
DEVICE=aspi4dos.sys /D
```

IMPORTANT REQUIREMENTS

The environment variable 'TEMP' should be set as is recommended. The application relies on it to create temporary files.

When printing in Windows 3.1 with the HP LaserJet 4 family, use raster mode, not HP-GL/2 mode for graphics.

For Systems with Multiple SCSI Cards

Previous versions of Molecular Analyst could communicate successfully with a Molecular Imager *only if* the SCSI card it is connected to had a host adapter number equal to 0. It is now possible to specify the host adapter number for the SCSI card. To specify the host adapter number, add the following lines to your MA.INI file:

```
[Imager]  
HostAdapterNum=x (depends on which SCSI card the Imager is attached to)
```

To determine which SCSI bus the Imager is attached to, add the /D switch to the ASPI4DOS.SYS command line. Reboot the PC with the scanner on, and note on which host adapter the scanner is attached. We recommend the following configuration:

```
DEVICE=ASPI4DOS.SYS /p<port address> /d
```

The card at IO port xxx will be configured as host adapter number 0. The card at IO port yyy will be host adapter number 1. Make sure to configure the IRQ and DMA settings of any additional cards to prevent conflicts. This can be done through the program SCSISelect, activated by hitting Ctrl-A during SCSI driver initialization. This is only available however when the card's BIOS is enabled.

Molecular Analyst / Densitometry Version 1.5

New Item for Windows 95 Users:

After installation, the Molecular Analyst program group contains an additional icon, "Install HASP Hardware Key". Window 95 users must double click on the "Install HASP Hardware Key" icon to install necessary system drivers to properly access the hardware key.

Addition of Freehand Draw Volume Analysis Tool

Molecular Analyst now has a freehand draw volume analysis tool. This tool provides users a method to define a number of irregularly shaped objects for volume analysis.

TIFF Export

In some cases, the previous version of Molecular Analyst incorrectly exports the calibration table during TIFF export. When the error occurs the calibration table values are off by a factor of 10. This would cause an O.D. value of 1.0 to be incorrectly exported as 10.0. This has been fixed in Molecular Analyst 1.5.

Fixed Windows 95 Image Printing Problem:

Printing images from Molecular Analyst, running under Windows 95, to Hewlett-Packard Laser Jet printers now works correctly.

VERSION HISTORY

Version 1.4.1 Revisions

Support for Windows 95 Operating System

Molecular Analyst now runs under Windows 95.

Version 1.4 Revisions

Scanner Support

Now supports 525 Molecular Imager

Saving TIFF Images

Images can now be saved in 3 different TIFF formats.

1. Raw - saves raw data only
2. View - saves 8-bit display data with a table of palette indices
3. Calibrate - saves raw data with calibration table

Reading TIFF Images

TIFF images can now be read using the IMPORT function in the FILE Menu. Support is available for reading 8-bit and 16-bit raw data TIFF files, 8-bit display data with a table of palette colors, and 8-bit and 12-bit data with a calibration table. RGB data import is *not* supported.

This TIFF import/export is compatible with most popular third-party packages.

Version 1.3 Revisions

.IMG and .SCN Association

.IMG and .SCN files are automatically associated with Molecular Analyst/PC during the setup process. When files with these extensions are double clicked, Molecular Analyst/PC is executed and opens the associated file.

Table Report Heading

The printed Table Report now has a heading that displays the file name and page number.

Object Table Save and Export

The table save and user interface has changed. The table is now saved with the image. The table may be saved in the old manner to a text file via the *File/Export* menu item.