

NOTICE: Varian, Inc. was acquired by Agilent Technologies in May 2010. This document is provided as a courtesy but is no longer kept current and thus will contain historical references to Varian. For more information, go to [www.agilent.com/chem](http://www.agilent.com/chem).



Varian, Inc.  
2700 Mitchell Drive  
Walnut Creek, CA 94598-1675/USA

---

# Star Chromatography Workstation

## Version 6.41

# Upgrade Notes

If you have Internet access and a web browser installed on your computer, visit the Star Chromatography Workstation Product Support Web Site at [www.chromatography.varian.com](http://www.chromatography.varian.com). Here you will find the latest software and documentation updates for the Star Chromatography Workstation suite of products, along with additional notes, tips, and answers to frequently asked questions. You may wish to visit this site periodically to see if new information is available that may be pertinent to you.

---

## What is New in Version 6.41

Version 6.41 is a maintenance release of Version 6.30, containing the following fixes and improvements:

- Integrates the drivers for the ProStar 325 UV-Vis detector (in the LC Control) and the ProStar 335 UV-Vis PDA detector (in PolyView 2000).
- Added capability to quickly update a Recalc list when the data files it refers to has been moved to another location.
- Added search capability to the message log viewer.
- Added ability to save size and settings to some of the file selection dialogs.
- Improvements to the Combi PAL driver: eliminated occasional communication errors, added the ability to agitate without heating, other minor improvements.
- And many other minor corrections and improvements...

---

## What was New in Version 6.30

Version 6.30 is a maintenance release of Version 6.20, containing the following fixes and improvements:

- Corrections to the Access Control and Audit Trail option,
- A fix for a problem limited to file access rights under Windows XP when using an account without local Administrator Privileges,
- The ability to add the Combi PAL AutoSampler to instruments containing a 3900 or 39XL GC and no other AutoSampler without resorting to STAR32.INI file switches.
- Usability improvements in System Control: AutoControl module is highlighted in Instrument Status Window, double-clicking on a module in the Instrument Status Window restores the module window to its most recent size, instead of showing it maximized.

---

## What was New in Version 6.20

Version 6.20 is a maintenance release of Version 6.0, containing corrections to the Access Control and Audit Trail option, support for Windows XP Professional, improvements to several drivers (Combi PAL, ProStar 500, ProStar 701, ProStar/Dynamax System) and the ability to acquire data from 2 separate CIMs in the ProStar/Dynamax system.

---

## What was New in Version 6

**‘21CFR Part 11’:** Star Workstation Version 6 supports a new add-on ‘Access Control and Audit Trail option’, which provides native tools to facilitate your organization’s compliance with the requirements of the Code of Federal Regulations, Title 21, Part 11. Refer to the [Access Control and Audit Trail documentation](#) for details.

### **New and Updated Drivers**

- ProStar 310 / 9050 UV-Vis detector driver has new digital filtering capabilities
- ProStar 364 4-channel Fluorescence Detector
- ProStar 500 Column Valve Module
- ProStar 701 Fraction Collector driver now supports method download
- 39XL GC
- CP-4900 Micro-GC driver
- Combi PAL AutoSampler

**Permanent Message Logs:** Message Logs documenting the automation processes are now saved for each automation process under unique names, and a new application lets you review and print past message logs.

**StarSD and StarDHA** are no longer on the Star Workstation CD, separate media will be provided with the products.

---

## Read Before Installing Star Workstation Version 6

Refer to the following documents, also accessible from the Star Workstation CD ROM, for additional information important to customers upgrading to Star Workstation Version 6.

### **Star Chromatography Workstation Release Notes**

These notes contain instructions for customers running Star Workstation under Windows NT, as well as important notes and safeguards to ensure the most reliable system operation.

### **LC Control Release Notes**

Information for Star Workstation customers running LC Control software.

### **GC Control Release Notes**

Information for Star Workstation customers running GC Control software.

## Upgrading to Star WS Ver. 6.41 from Ver. 6.x

This is an automated process. Use the 'Update Version 6.x to Version 6.41' button displayed in the CD Browser after you select 'Install → Install Upgrades and Updates'. Existing Star WS Ver. 6.x serial numbers will be used to install the Star WS Ver. 6.41 software. Please have your Star WS Ver. 6.x serial numbers available in case you need to manually enter the serial numbers.

## Upgrading to Star WS 6 from Star 5.x

### ***Prior to upgrading to Star WS Ver. 6 from Ver. 5.x, you may need to save your Star32.ini***

Upgrading from Star WS Ver. 5.x will uninstall Star WS Ver. 5.x. Uninstalling Star WS Version 5.x will also remove the star32.ini file located in the Windows directory. This file contains all of the preference settings used by Star Workstation. Installing Star Workstation Version 6 installs a new default version of star32.ini. In some cases, you may have edited star32.ini manually and you may want to use these same settings in the new star32.ini. If so, you should make a copy of the original star32.ini so that you can refer to it as you make changes to the new one. **Do not, however, simply replace the new one with your old star32.ini.**

In particular, you may want to verify the following settings:

- The "directory" field in the [Varian Star Workstation] section.
- The "GPIBAddress" field in the [System Control] section.
- The "ADC\_Board\_Base\_Address" in the [System Control] section.

### ***Performing the upgrade to Star WS 6 from Star 5.x***

This is an automated process. Use the 'Upgrade from Ver.5.x to Version 6' button displayed in the CD Browser after you select 'Install → Install Upgrades and Updates'. Existing 5.x serial numbers will be converted to Version 6 serial numbers, and a one page document will be composed showing your new serial numbers. Please print and save this document with the license cards for the software used on this workstation. These serial numbers will be needed should if you have to reinstall the Star Workstation Version 6 software. The installation will proceed, re-installing the existing options and automatically providing the new serial numbers; you may add new options (such as 'Access Control and Audit Trail') if you purchased them at the same time as your upgrade.

**NOTE: On some Windows 2000 and Windows NT installations, the Star Workstation Version 6 installation engine may not be able to automatically enter the Star WS Version 6 serial numbers for your existing installed Star 5.x applications. In this case, you will need to manually enter the Star WS Version 6 serial numbers that you printed after Star WS 5.x was removed.**

## Removing a Version Prior to Star 5

Prior to installing Version 6 on a computer already containing a version of the Star Workstation prior to Star 5, it is strongly recommended that you remove the previous version. You may use the Data Files, Methods, SampleLists, SampleLogs, and Sequences created with the previous version of the Star Workstation with Version 6 (see upgrade notes below). To remove the previous version, perform the following steps. In this procedure, it is assumed that the previous version of the Star Workstation is installed in the directory C:\STAR.

1. Close all Star Workstation applications that may be running.
2. Copy the Data Files, Methods, SampleLists/SampleLogs, and Sequences that you wish to upgrade to a backup directory (for example, C:\STAR45).
3. Delete the STAR directory using the Windows Explorer.
4. Run the System Configuration Editor by selecting Run... from the Windows Start button and typing sysedit.exe. The System Configuration Editor allows you to edit autoexec.bat and win.ini.
5. Edit your autoexec.bat file and remove C:\STAR from your path.
6. Edit the file win.ini. Find the line starting with "run=" in the file. Remove start.exe from the line. For example, if the line was 'run=start.exe', replace it with 'run='.
7. Delete the file star.ini from your Windows directory.
8. Reboot your computer.

You should also remove the *Varian Star 4.5* item from the Windows Start Menu. To do so, use the following procedure:

1. Right-click on the Windows taskbar (where the Start button is located) and select *Properties* from the popup menu.
2. Select the Start Menu Programs tab in the Taskbar Properties dialog box.
3. Click on the *Remove* button and select the *Varian Star 4.5* item from the Programs folder.
4. Click on the *Remove* button to remove the item.

If you have made associations between files that end with .RUN and the Interactive Graphics/Data Handling application in versions of the Star Workstation prior to 5.0, you will need to remove this association before installing Version 6. Similarly, associations between files that end with .MTH and the Method Editor application will also need to be removed. *This is not needed if you upgrade from 5.x.*

To remove an association between an application and a file type, use the following procedure:

1. In My Computer or Windows Explorer, click the View menu, and then click Options.
2. Click the File Types tab.
3. In the list of file types, click the one you want to remove (for example RUN file or MTH file).
4. The settings for that file type are shown in the File Type Details box.
5. Click Remove.

If you do not remove these associations before installing Star Workstation Version 6, you may see inappropriate selections available on Data File or Method Quick Link buttons. If you select one of these inappropriate entries, you may get error messages stating that the file you are attempting to access is corrupt.

## Upgrading Data Files, Methods, SampleLists and Sequences

Data Files, Methods, SampleLists, and Sequences are automatically upgraded from versions prior to Version 6 as soon as they are accessed by the Star Workstation Version 6. There have been no changes to the data file format from Star 5.x to Star 6, so files created with Version 5.x will not be upgraded. A message box will prompt you to upgrade the file. Note that if you cancel the upgrade operation, you may be prompted several times with this message box—once for each section of the file requiring the upgrade. To preserve the contents of the file in the previous version's format, it is important that each message box is canceled.

If you wish to automatically upgrade all files you access without being prompted, use the following procedure:

1. Edit the STAR32.INI file in your Windows directory using a text editor such as Notepad.
2. Find the line starting with "AutoUpgrade=" in the file. Change the line so that it reads **AutoUpgrade=1**.
3. Save the file.