

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 This document is provided as a courtesy but is no longer kept current and thus will contain historical references to Varian.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).



**Agilent Technologies**

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

---

## Star Chromatography Workstation Version 6

# Star Custom Report Writer Upgrade and Release Notes

The following information can be found in this document:

### **Upgrading from Previous Versions of Star Custom Report Writer**

For customers who are installing Version 6 on a computer where a version prior to version 6.0 is installed. Also includes operational differences between Version 6 and previous versions not covered in other documents.

### **Miscellaneous Notes**

Additional information for all Star Report Writer customers not covered in other documents.

The following information can be found in other related documents:



### **Star Chromatography Workstation Upgrade Notes**

For customers who are installing Version 6 on a computer where a previous version is installed. Also includes operational differences between Version 6 and previous versions not covered in other documents.



### **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 from Previous Versions of the Star Custom Report Writer

## Report Template upgrades

Report templates generated using versions of the Star Report Writer older than 5.0 are automatically upgraded when they are accessed by Version 6. You will be prompted to save the upgraded template even if you made no other changes to it. Once it is saved, it can no longer be used by the previous version. Templates generated by version 5.x are directly compatible with version 6 and do not need to be upgraded.

## Upgrading Old Report Templates containing bitmaps

You may get an "Unexpected File Format" error when old (Pre 5.0) report templates containing bitmaps are opened in Version 6. To avoid this situation, it is recommended that you remove bitmaps embedded in 'Pre 5.0' report templates before opening the templates in Version 6. After the templates have been upgraded, you can then embed the bitmaps back into the templates.

---

## Miscellaneous Notes

### Embedding OLE objects into Report Templates

If you are embedding an OLE object into a Star Report Writer Template (by using the *OLE Object* icon in the *General* tab of the Report Sections window), do not have the Palette dialog box displayed in Star Report Writer. If you do, an application error will result and Star Report Writer will close. Once an OLE object has been embedded into a Report Template, you may open and use the Palette dialog box *as long as the OLE object is not activated*. Be sure to close the Palette Dialog box *before* activating your OLE object.

### Overlaying Report Objects

If report objects such as the chromatogram object are overlaid on top of each other and *Overflow Printing* is checked in the File menu, the overlaid objects may be offset vertically in the report printout. The amount of the offset varies depending upon the number of objects appearing above the overlaid objects in the report template. If the overlaid objects are at the top edge of the report, no vertical offset will occur. To avoid the offset, you can disable overflow printing, but make sure that all report objects in the template are sized appropriately for the data that they will display.

## Displaying Report Templates without a data source specified

If you open a Report Template with no data source specified, chromatogram objects that specify a fixed data source will not be drawn until you activate them (by double-clicking on the object) or select a data source.

## Automation Files Object fields

The *RecalcList* field in the Automation Files Object will display the old SampleLog name and won't be updated to display the RecalcList (if any) used in the most recent recalculation of the Data File from System Control. When the Data File is recalculated from Interactive Graphics, both of these fields correctly display "Interactive Graphics".

## Localization of Objects

With the exception of the chromatogram object, all of the report objects are designed to use the decimal point, date and time conventions specified in the Windows Control Panel Regional Settings. Numbers displayed in the chromatogram object always use a "." for the decimal point.

## Missing Status Codes

Some Star Custom Report Writer objects have a key to the status codes used in the data table and their meanings printed at the end of the object. This status code key resides in the first column of the table and will NOT be printed if the first column is not printed. Whenever status codes are used in an object and they are not printed at the end of the object, check to make sure that the first column of the table is enabled. If it is not selected, enabling its printing will restore the printing of the status code key for that object.

## Solvent Gradient Display

If you are using the ProStar / Dynamax HPLC System and are running a solvent gradient, plots of the gradient profiles will be added to the chromatogram display. When viewed on the screen or printed on a color printer, each of the profiles will be plotted in a different color. A Solvent %s scale will also be displayed at the right of the chromatogram.

To learn more about this feature, or to turn it off, please consult the [Star Custom Report Writer Operation Manual](#)

## Access Control and Audit Trail support

Version 6.0 supports the Access Control and Audit Trail option if it is installed. Access to the application requires the operator to log-in and to have certain rights, and the Revision Log object will display the Audit Trail for the data file.