

# Overview of SpectrAA-300/400 Software Capability

## Application Note

Atomic Absorption

### Author

Jonathan Moffett

### Introduction

It is probable that operators of the SpectrAA-300/400 are more instrument orientated than computer orientated. Indeed it is possible to operate the SpectrAA-300/400 purely as an instrument. However operators wishing to make advantage of the controlling computer can refer to this summary.

There has been an increasing trend for scientific instruments to be designed as "black boxes" controlled by an external microcomputer. Except for basic instruments and special requirements, there are a number of reasons for this.

Microcomputers have become increasingly smaller, more powerful and lower priced. The average personal computer (PC) can potentially provide much greater automation of instruments than previously possible except for expensive dedicated instruments. The PC also provides powerful data manipulation and so can produce final answers rather than raw data from the instrument. Additionally it is possible to transmit data between computers and thus a variety of instruments can send their results to a central data collection point to be integrated into a complete report.

The SpectrAA series was the first commercial atomic absorption spectrometer to use a central computer/ controller to operate both the base instrument and its accessories. SpectrAA was originally designed around an Agilent Data Station (DS-15). However SpectrAA can also be controlled by an IBM PC. This configuration is designated 300/400.

DS-15 is a dedicated instrument controller and has the facility to transfer data to a computer for data manipulation. However the use of a PC as an instrument controller has advantages. Hardware improvements include fixed magnetic data storage devices that allow rapid storage and retrieval of large amounts of data and monitors that provide high resolution color displays. There is a wide range of software available that allow data manipulation within the PC environment itself.



**Agilent Technologies**

## Hardware Configuration

While the IBM PCs have many clones and compatibles, not all of these behave in a manner identical to the IBM. The computer systems known to operate as per the SpectrAA 300/400 Operation Manual are listed in Table 1 although performance is guaranteed on IBM computers only.

Table 1. Recommended Computer Systems for SpectrAA

Computer	Recommended memory (RAM)	Comments
EPSON AX-2	640 KB	80286 CPU
IBM PS/2 Model 30	640 KB	8086 CPU; superseded
IBM PS/2 Model 30/286	1024 KB	80286 CPU

NOTES: A fixed (hard) disk (20 MB or greater) is required. Display must be equivalent to 128 KB EGA or better (built-in IBM Model 30/286 display card with the IBM 8513 monitor is suitable). SpectrAA IDB card must be installed.

## Brief Overview of DOS

Each computer requires an operating system (OS) to be loaded into its memory. IBM PCs and compatibles use a Disk Operating System (DOS). All DOS information is kept in files which must have unique names. DOS nomenclature allows up to eight alphanumeric characters for the name and up to three characters for an extension (for example, FILENAME.EXT). Files are stored on (magnetic) media called disks. Disks can either be removable (floppy) or fixed (hard). DOS assigns letters to the disk drives. The computers recommended for SpectrAA have one floppy drive (A:) and one hard drive (C:). The names of all the files present on a disk are stored in a directory. Floppy disks with their limited storage are generally assigned a single (root) directory called A:\. Hard disks have a much larger capacity and for convenience and good house-keeping are divided into multiple directories and subdirectories (which are all chained or PATHed back to the root directory, C:\). Directory names also follow DOS nomenclature rules. Hence the SpectrAA software is installed on the C: drive under the AGILENT directory in a subdirectory called SPECTRAA. This path is designated as C:\AGILENT\SPECTRAA\, which is displayed if the DOS command PROMPT \$PSG is used.

## SpectrAA Software

There are two forms of SpectrAA software - flame/furnace (SPECTRAA) and Zeeman (ZEEMAN). Where appropriate in examples below both names are shown. In examples that only show SPECTRAA, ZEEMAN may be used interchangeably. Each form is distributed on one 3.5" 720 KB floppy disk or one 5.25" 1.2 MB high density floppy disk depending on the floppy drive of the controlling computer.

The distribution disk contains four files:

ARCE.COM  
INSTALL.BAT  
INSTALL.DOC  
SPECTRAA.ARC (or ZEEMAN.ARC)

To install, type A: [ENTER] to change to A drive. Type INSTALL [ENTER] to run the INSTALL.BAT file. The SpectrAA files are all compressed into one archive file (.ARC) and the program ARCE.COM "unpacks" the files which are then copied into the appropriate directories. The final file INSTALL.DOC contains information for use at installation. This may be read by the command

```
TYPE INSTALL.DOC [ENTER]
```

which displays it on the screen or

```
TYPE INSTALL.DOC > PRN: [ENTER]
```

which directs the text to a printer.

After installation, the distribution disk should be stored away. It is recommended a copy be made of it for additional security.

To run the SpectrAA instrument, type SPECTRAA (or ZEEMAN) [ENTER]. Within a few seconds a welcome screen (with a version number) is displayed and then the instrument software PROGRAM MODES page. If this does not happen, turn the computer OFF and follow the troubleshooting directions below.

## Troubleshooting

- Program does not load but returns immediately to DOS command line.  
There is insufficient RAM memory available. SpectrAA requires about 400 KB. Check that memory is not being occupied by a program which terminates but stays resident (TSR) for example, Borland Sidekick or by a RAM disk. Also check the CONFIG.SYS file for excess open files or buffers.
- Program loads but only displays the welcome page.  
The IDB card is not installed. This card must be present for the software to function.

The SpectrAA software does not address any memory above the 640 KB DOS limit. Operating SpectrAA should therefore not affect any customization of additional RAM. Once the software has been loaded, the computer acts as a controller for the SpectrAA instrument.

## Review of DOS Data Output

The data generated by a software package such as SpectrAA should be recorded in a permanent method as well as being displayed. The simplest method is to direct output to a printer which is what SpectrAA as supplied does. In addition to a printer, DOS is able to address a number of devices and the three most commonly used devices are outlined in Table 2.

Table 2. DOS Devices

### Device–DOS Name–Example

File on disk drive:\path\filename.ext A:\TEST\DATA.TXT

Parallel port PRINTER:, PRN:, LPT1: Printer (Centronics)

Serial port COM1:, COM2:, AUX: Printer (Serial) or another computer.

#### NOTES:

1. The drive need only be specified if it is not the current drive (usually C:).
2. The colon (:) is essential as it differentiates between a physical device and a file.
3. SpectrAA is supplied configured for an EPSON compatible parallel printer.

It is possible to redirect SpectrAA's output if required. When sending data to an external device such as a printer or another computer, care is needed to place it in a known condition. For serial devices, the transmission rate must be specified. For parallel devices, special ASCII escape code sequences must be transmitted along with the data.

An operator can select most of SpectrAA's communication options with the SYSTEM CONFIGURATION page (page 20) shown in Figure 1. Under this page is the EXTERNAL CONTROL subpage shown in Figure 2. OUTPUT CONTROL (Figure 3) is used to select the device and PRINTER CONTROL (Figure 4) to specify the escape codes. Table 4 summarizes the escape codes most often used with SpectrAA. Figure 4 shows the default printer conditions supplied with the SpectrAA software.

PROGRAM 52	Cu EPA Test QC ON4	FURNACE
BLANK	REPLICATE 1	
ABSORBANCE	CONCENTRATION	SAMPLER AUTOMIX
PEAK HEIGHT	1.0 (sec)	BC OFF
PRINTING		
SYSTEM CONFIGURATION		
COLOR SELECTION	FOR COLOR SELECTION	
EXTERNAL CONTROL	FOR SETTING COMMUNICATION PARAMETERS	
		F COLOR F EXTERNAL 5 SELECTION 6 CONTROL

Figure 1. Main SYSTEM CONFIGURATION page screen.

PROGRAM 52	Cu EPA Test QC ON4	FURNACE
BLANK	REPLICATE 1	
ABSORBANCE	CONCENTRATION	SAMPLER AUTOMIX
PEAK HEIGHT	1.0 (sec)	BC OFF
PRINTING		
EXTERNAL CONTROL		
INPUT CONTROL	TO DEFINE THE INPUT CHANNEL CHARACTERISTICS	
OUTPUT CONTROL	TO DEFINE THE OUTPUT CHANNEL CHARACTERISTICS	
PRINTER CONTROL	TO DEFINE THE CONTROL AND INITIALIZATION STRINGS	
SYNCHRONIZATION	TO DEFINE SYNCHRONIZATION AND HANDSHAKING CHARACTERISTICS FOR EXTERNAL CONTROL	
LEARN CONTROL	TO CREATE A FILE OF RECORDED KEYSTROKES TO BE REPLAYED LATER	
		F INPUT F OUTPUT F PRINTER F SYNCH F LEARN 2 CONTROL 3 CONTROL 4 CONTROL 5 CONTROL 6 CONTROL

Figure 2. EXTERNAL CONTROL option selection page.

PROGRAM 52	Cu EPA Test QC ON4	FURNACE
BLANK	REPLICATE 1	
ABSORBANCE	CONCENTRATION	SAMPLER AUTOMIX
PEAK HEIGHT	1.0 (sec)	BC OFF
PRINTING		
OUTPUT CONTROL		
OUTPUT CHANNEL NAME	_MOFF\SSC.PRN	
		F OUTPUT F INPUT F PRINTER F SYNCH F LEARN 1 SETUP 2 CONTROL 4 CONTROL 5 CONTROL 6 CONTROL

Figure 3. Example of SpectrAA's output being directed to a file SSC.PRN in the author's directory on drive C:.

PROGRAM 52	Cu EPA Test QC ON4	FURNACE
BLANK	REPLICATE 1	
ABSORBANCE	CONCENTRATION	SAMPLER AUTOMIX
PEAK HEIGHT	1.0 (sec)	BC OFF
PRINTING		
PRINTER CONTROL		
PRINTER INITIALIZATION STRING (13) (27) (10) (27) (10) (10) (13)		
REPORT START STRING (14)		
REPORT TERMINATOR STRING (26)		
OUTPUT RECORD TERMINATOR (10) (13)		
		F INPUT F OUTPUT F SYNCH F LEARN 2 CONTROL 3 CONTROL 5 CONTROL 6 CONTROL

Figure 4. Default PRINTER CONTROL screen.

## Output Data to Printer

The default condition as supplied on the distribution disk is to output the report to an Epson printer which is Centronics compatible. The necessary escape codes are shown in Table 3. Table 4 shows how to configure SpectrAA for the default PRINTER: output.

Table 3. Selected SpectrAA ASCII Codes  
(To be entered on PRINTER CONTROL page)

Field string	Printer (default)	File
PRINTER		(leave blank)
INITIALIZATION	{13}{27}@{27}O{27}H{18}{13}	or optional
REPORT START	{14}	message)
REPORT TERMINATOR	{26}	{13}{10}
OUTPUT TERMINATOR	{10}{13}	{13}{10}
NOTE:	{10} - Line feed (LF)	
	{13} - Carriage Return (CR)	
	{14} - Double width (Epson)	
	{18} - Cancel condensed mode (Epson)	
	{26} - End of file (EOF)	
	{27} - Escape (ESC)	
	{27}@ - Initialize printer (Epson)	
	{27}O - Cancel bottom margin (Epson)	
	{27}H - Cancel double strike (Epson)	

Table 4. Configuring SpectrAA for PRINTER:

Keypresses	Comments
[F10]	Go to INDEX page
20 [ENTER]	Go to SYSTEM CONFIGURATION page ([F6] or [F10] also accepts the page number)
[F6]	Go to EXTERNAL CONTROL page
[F3]	Go to OUTPUT CONTROL page
[DELETE]	Clear device name.
[F1]	Setup OUTPUT CHANNEL NAME The default PRINTER: appears
[F4]	Go to PRINTER CONTROL page
Enter escape codes	Refer to Table 4 (Figure 4)
[F10]	Return to INDEX page.

## Output Data to File

SpectrAA can be configured to output its report to a file on disk (Table 5) for later use. Figure 3 shows an example of file output. Such an option has two major uses, spooling and data manipulation.

Spooling is simply storing the printer output until it can be printed out at a later time. The printer codes can be left as the default ones supplied. Graphics data such as calibration graphs and screen prints can be included. Printing out this file is discussed in the review.

The most common use of a data file is incorporating the output in a spreadsheet or text editor program. Including graphics data is not recommended. This is done by changing the CALIBRATION GRAPH option to NO on REPORT FORMAT page (SpectrAA page 13 as shown in Figure 6). Do not press [PRINT SCREEN] key. The PRINTER option LX-800 (monochrome printer) should be used to suppress color codes.

If the file name used in Table 5 is not changed then successive reports are chained together. It is recommended that the ASCII character 26 (EOF) not be used to terminate a report. If the file is to be imported into another software package the EOF character may cause only the first report to be recognized. Custom messages or strings can be inserted in the first two printer control fields if desired.

Table 5. Configuring SpectrAA for File Output

Keypresses	Comments
[F10]	Go to INDEX page
20 [ENTER]	Go to SYSTEM CONFIGURATION page
[F6]	Go to EXTERNAL CONTROL page
[F3]	Go to OUTPUT CONTROL page
[DELETE]	Clear device name
filename.ext	Follow DOS rules, 16 characters allowed
[F1]	Setup OUTPUT CHANNEL NAME
[F4]	Go to PRINTER CONTROL page
Enter escape codes	Refer to Table 3
[F10]	Return to INDEX page

## Output Data to Serial Port

SpectrAA can be configured to output the report through the serial communication port (Table 6). In this option, data is sent out as if going to a printer (live). It is suggested that the file escape codes of Table 4 be used and that no graphics data are sent as discussed in Output Data File. It is also possible to transmit a report file by this method.

The recommended computer systems come with one serial port, designated COM1:. If it is to be used, this port must first be initialized at the DOS command level. The parameters used must match those of the receiving device and the appropriate operating manuals must be consulted.

To use an example from another application, the Varian Report Manager program used for trapping data from a DS-15 via the RS-232A channel uses 9600 baud, even parity, 8 databits and 1 stopbit through serial port COM1:. The DOS initialization is then

```
MODE COM1:96,E,8,1
```

Initialization of the communications port must be done before running SpectrAA. It is recommended that the MODE command be incorporated in the AUTOEXEC.BAT file so that the port is initialized on power up. Note however that SpectrAA's output must still be set to COM1: as shown in Table 6.

The most common application is a one way data communication to another computer. The receiving computer must have appropriate software to scan the serial port continuously for incoming data. There are many commercial and public domain programs available for this purpose.

For the serial link between the computers a three wire "null modem" connection with appropriate plug jumpers must be used. A configuration known to work is detailed in Figure 5.

For two way communication, the EXTERNAL CONTROL option must be used.

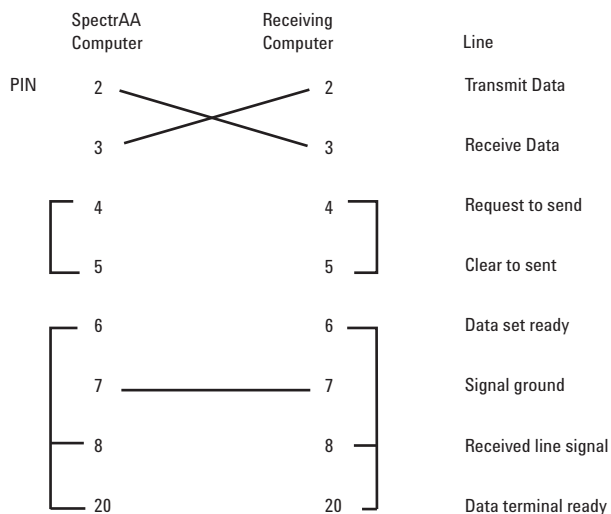


Figure 5. Serial computer linking (PC to PC) using a standard D25 connector.

## Save Current Results on Drive:\Path\

When [F11] (START) is pressed to start an automatic run all readings and instrument settings are stored in a current results file in the data file directory shown on the REPORT FORMAT page (Figure 6). When the run is completed as many reports as desired may be generated from the current results file. Note that when [F11] is pressed this file is automatically cleared for the new automatic run results.

Table 6 Configuring SpectrAA for COM1:

Keypresses	Comments
Initialize serial port	Use MODE command at DOS level
[F10]	Go to INDEX page
20 [ENTER]	Go to SYSTEM CONFIGURATION page
[F6]	Go to EXTERNAL CONTROL page
[F3]	Go to OUTPUT CONTROL page
[DELETE]	Clear device name.
COM1:	Do not omit the colon (:), otherwise a file named COM1 will be created.
[F1]	Setup OUTPUT CHANNEL NAME
[F4]	Go to PRINTER CONTROL page
Enter escape codes	Refer to Table 3
[F10]	Return to INDEX page.

Different operators using the same spectrometer may wish to keep their current results file in separate data file directories (as the author did in Figure 6). Changing the data file directory is outlined in Table 7. If the directory name field is blank, then the default directory (C:\AGILENT\SPECTRAA\ or C:\AGILENT\ZEEMAN\) is used.

PROGRAM 52		Cu EPA Test QC 0N4		FURNACE
BLANK	REPLICATE	1		
ABSORBANCE	CONCENTRATION		SAMPLER AUTOMIX	
PEAK HEIGHT	1.0	(sec)	BC OFF	
<b>PRINTING</b>				
<b>REPORT FORMAT</b>				
OPERATOR	-			
DATE	-			
BATCH NAME	-			
FORMAT	SEQUENTIAL			
DATA PRINTED	EACH READING			
PRINT	DURING RUN			
INSTRUMENT STATUS	YES			
NOTES	YES			
CALIBRATION RESULTS	YES			
CALIBRATION GRAPH	YES			
SAMPLE LABELS	YES			
LINES PER REPORT PAGE	70			
PRINTER TYPE	LX800			
DATA FILE DIRECTORY	C:\MOFF\			
F PRINT	Z STATUS	F SAMPLE	S LABELS	

Figure 6. Example where the current results file is in the author's directory on drive C:.

Table 7. Changing Location of Current Results File (Data File Directory)

Keypresses	Comments
[F10]	Go to INDEX page
13 [ENTER]	Go to REPORT FORMAT page
[Up arrow]	Move cursor to DATA FILE DIRECTORY Either select default directory
[DELETE] [ENTER]	(C:\AGILENT\SPECTRAA\ by having a blank line. Or select another directory following
Drive:\path\ [F10]	DOS rules for naming the directory Return to INDEX page (up to 30 characters).

NOTE: This can be incorporated into setting up an automatic run as SpectrAA goes through the REPORT FORMAT page.

### Archive Results on Drive:\Path\

To store a current results file permanently, the file must be archived. This is done in the Utilities section of SpectrAA in the ARCHIVE SET DIRECTORY page shown in Figure 7. The current results directory (DATA FILE DIRECTORY) and the archive directory (ARCHIVE FILE DIRECTORY) may be specified. If blank, the default directory (C:\AGILENT\SPECTRAA\ or C:\AGILENT\ZEEMAN\) is used.

PROGRAM 52	Cu EPA Test QC OM4	FURNACE
BLANK	REPLICATE 1	
ABSORBANCE	CONCENTRATION	SAMPLER AUTOMIX
PEAK HEIGHT	1.0 (sec)	BC OFF
PRINTING		
ARCHIVE SET DIRECTORY		
1.	11.	
2.	12.	
3.	13.	
4.	14.	
5.	15.	
6.	16.	
7.	17.	
8.	18.	
9.	19.	
10.	20.	
DATA FILE DIRECTORY C:\MOFFN		
ARCHIVE FILE DIRECTORY A:\		
REQUIRED SET NO. 1		
F 1 PURGE	F 4 CREATE R/OR READ	F 5 ARCHIVE F 6 UTILITIES INDEX

Figure 7. Example in which the file from Figure 6 is being archived onto a floppy disk.

Table 8. Archiving Current Results

Keypresses	Comments
[F10]	Go to INDEX page
21 [ENTER]	Go to UTILITIES INDEX page
2 [ENTER]	Go to ARCHIVE SET DIRECTORY page
	Optional
	Change data file: Move cursor to DATA FILE DIRECTORY line
drive:\path\ [F4]	Give valid DOS drive and path (up to 30 characters allowed) Read the file.
	Optional
	Change archive file: Move cursor to ARCHIVE FILE DIRECTORY line.
drive:\path\ [F4]	Give valid DOS drive and path (up to 30 characters allowed) Read the archive set or prepare for one
	Save current results file: Move cursor (if necessary) to REQUIRED SET NO. line
Enter valid number [F5]	Number from 1 to 20 Archive current results data file Optional Purge data set
Enter valid number [F1]	Enter number from 1 to 20 of existing set Remove the existing archived data file.

NOTE: Sample weight and/or volume correction values will be archived provided they are entered using page 5 of the UTILITIES INDEX before archiving results.

### Record LEARN Files

No matter how flexible a software package is, using the same sequence of keystrokes frequently can become tedious. SpectrAA has the facility to record keypresses and to replay these memorized keypresses. For example if an operator wishes to print out results during an automatic run and then to place the report in a file after the run, the printer codes that need to be entered can all be done using a LEARN file. Naming the file and starting the recording process is shown in Figure 8. Note that when recording/replaying a LEARN file, the INDEX page is the first page to be used.

It must be noted that the LEARN mode does not record settings, only keypresses. To avoid unexpected outcomes input fields that are being changed must be initialized. For numeric field entries press [DELETE] and the required digit keys. For [HOME] selectable fields first press [CONTROL] [HOME] to reset the field and then press [HOME] until the desired option is seen.

In LEARN mode most keys behave as they do in the normal mode. However some keys do have special functions and these are summarized in Table 9.

Table 9 Special LEARN Mode Commands

Key	Function
[INSERT]	Replay a LEARN file
[END]	Terminate LEARN mode
[CONTROL] [HOME]	Reset menu choice to default condition
[CONTROL] [PAGE UP]	Rewind/Terminate LEARN mode
[CONTROL] [PAGE DOWN]	Pause three seconds
[CONTROL] [DELETE]	Clear NOTES page (from any page)

## Replay LEARN File

A LEARN file whose name appears on the LEARN CONTROL subpage (Figure 8) can be replayed any time the [INSERT] key is pressed. If another file is to be used, its name must first be entered as shown in Table 10.

The screenshot shows a terminal window with the following text:

```

PROGRAM 52      Cu EPA Test QC 0M4      FURNACE
BLANK          REPLICATE 1
ABSORBANCE    CONCENTRATION 1      SAMPLER AUTOMIX
PEAK HEIGHT   1.0 (sec)      BC OFF
PRINTING
LEARN CONTROL

LEARN FILE NAME      LEARN.KEY
    
```

At the bottom, there is a status bar with the following text:

```

F START  P INPUT  F OUTPUT  F PRINTER  F SYNCH
1 LEARN  2 CONTROL 3 CONTROL 4 CONTROL 5 CONTROL
    
```

Figure 8. The LEARN CONTROL screen showing the file either where new keystrokes will be recorded or from which stored keystrokes will be replayed. The file would be in the C:\VARIAN\SPECTRAA\subdirectory.

Table 10. Recording a LEARN file

Keypresses	Comments
[F10]	Go to INDEX page
20 [ENTER]	Go to SYSTEM configuration page
[F6]	Go to EXTERNAL CONTROL page
[F6]	Go to LEARN CONTROL page
drive:\path\filename.ext	Use a valid DOS file name (up to 16 characters). It is suggested that the extension .KEY be used
[F1]	Start LEARN mode - INDEX page is shown automatically
Enter program	Configure SpectrAA as desired. Note that all keys behave as normal except for special ones listed in Table 9
[END]	Terminate LEARN mode

### NOTES –

- For numeric entries first press [DELETE] and enter the number.
- For [HOME] selected entries press [CONTROL] [HOME] and then [HOME] the required number of times for the desired option.

## External Control

SpectrAA has the capability of being operated from a remote computer. This is a specialized option and it is recommended only computer specialists implement this. Refer to the SpectrAA Operation Manual (Section 15 - External Control). For further information, contact your Agilent sales office.

## At DOS Command Line

### Review

Some DOS commands have already been outlined,

PROMPT \$P\$G

and

MODE COM1: [baud rate],[parity],[databits],[stopbits].

As mentioned above (Output Data to File) the printer output can be directed to a file for later printing. To send a file to a printer, use the DOS command

COPY filename PRN:

(note that DOS on occasion may have trouble handling graphics codes and output may not be what is expected).

There are some other SpectrAA options that are possible only at the DOS command level. As the commands may be used frequently and to avoid tedious typing, it is suggested that batch files for each option be created. A text editor or the DOS editor, EDLIN, may be used to create an ASCII text file containing the necessary DOS commands. This file is then given the extension .BAT and when run performs the commands. Some names for batch files are suggested below but of course other names could be used.

## Copy Method Files

SpectrAA can store up to 90 flame methods and up to 90 furnace methods. SpectrAA Zeeman can store up to 90 furnace methods. It is recommended that back-up copies of methods should be made in case the hard disk files are damaged.

CAUTION: This is not recommended as a means of transferring methods between spectrometers unless it is known they have the same software version and same hardware options. Otherwise it is suggested a LEARN file be used.

Each set of methods (flame, furnace and Zeeman) are stored in two files per set. The two files for each set must be copied as a pair otherwise the methods are not usable. Table 11 suggests batch file formats to ensure the files are copied as a pair to and from floppy disk.

Table 11. Copy SpectrAA Method Files

Copy flame and furnace methods to floppy disk (SAVEALL.BAT)	
CD\AGILENT\SPECTRAA	(Change directory)
COPY IOSPGRM.* A:	(Copy flame and furnace methods programs)
COPY IOSDIR.* A:	(Copy flame and furnace methods directory)
CD\	(Change to root directory)
Copy flame methods to floppy disk (SAVEFLAM.BAT)	
CD\AGILENT\SPECTRAA	(Change directory)
COPY IOSPGRM.I A:	(Copy flame method programs)
COPY IOSDIR.I A:	(Copy flame method directory)
CD\	(Change to root directory)
Copy furnace methods to floppy disk (SAVEFURN.BAT)	
CD\AGILENT\SPECTRAA	(Change directory)
COPY IOSPGRM.K A:	(Copy furnace method programs)
COPY IOSDIR.K A:	(Copy furnace method directory)
CD\	(Change to root directory)
Copy Zeeman methods to floppy disk (SAVEZEEM.BAT)	
CD\AGILENT\ZEEMAN	(Change directory)
COPY IOSPGRM.L A:	(Copy Zeeman method programs)
COPY IOSDIR.L A:	(Copy Zeeman method directory)
CD\	(Change to root directory)
Restore flame/furnace methods from floppy disk (GETMETH.BAT)	
CD\AGILENT\SPECTRAA	(Change to SpectrAA directory)
COPY A:\IOSPGRM.*	(Copy methods programs from disk)
COPY A:\IOSDIR.*	(Copy method directory from disk)
CD\	(Change to root directory)
Restore Zeeman methods from floppy disk (GETZMETH.BAT)	
CD\AGILENT\ZEEMAN	(Change to SpectrAA directory)
COPY A:\IOSPGRM.L	(Copy methods programs from disk)
COPY A:\IOSDIR.L	(Copy method directory from disk)
CD\	(Change to root directory)

For example, an operator of a flame system would create the SAVEFLAM and GETMETH batch files. To store the methods the operator places a formatted floppy disk in drive A: and at the C> prompt types SAVEFLAM. To reinstall these methods, the operator inserts the same floppy disk in drive A: and at the C> prompt types GETMETH.

## Copy Archive Files

It is suggested that the SpectrAA utilities be used to store the files onto a separate floppy disk (see Archive Results Drive:\Path). However an operator may wish to utilize the speed of a hard disk when archiving and editing data and then transfer files when the archive directory is full. In this case, it is easier to create a special directory at the DOS level using the make directory command MKDIR or MD. For example, assuming AGILENT\SPECTRAA exists, a subdirectory USER1 can be created as shown below:

```
MKDIR \AGILENT\SPECTRAA\USER1
```

Use this directory only for archiving data (see Archive Results Drive:\Path). Then at a suitable time such as when the archive data directory has all 20 sets filled, use the following batch file:

```
CD \AGILENT\SPECTRAA\USER1 (Change to example archive directory)
```

```
COPY *.* A: (Copy all files to drive A:)
```

```
DEL *.* (Delete existing files to clear directory)
```

```
CD\ (Change back to root directory)
```

## Move SPECTRAA.BAT/ZEEMAN.BAT Files

The SPECTRAA.BAT and ZEEMAN.BAT files are read-only protected so that they cannot be deleted inadvertently. These files are installed in the root directory. Operators may prefer to keep the root directory for essential files only and so create a separate directory for batch files (for example, C:\BAT).

Before a read-only file can be erased, the DOS command ATTRIB must be used to remove the read-only attribute. An example procedure is given below. As this should only be done once, creating a batch file would not be necessary.

```
CD\ (Change to root directory)
```

```
ATTRIB -R SPECTRAA.BAT (Remove read-only attribute)
```

```
COPY SPECTRAA.BAT C:\BAT (Copy to another directory)
```

```
ATTRIB +R C:\BAT\SPECTRAA.BAT (Set read-only attribute of copied file)
```

```
DEL SPECTRAA.BAT (Remove original file)
```

## Interaction with Third Party Software

### Export ASCII Files

Many programs generate files that can be read only by that program. Hence the current results file of SpectrAA could not be read by any other program. The most common way to pass information between different programs is to use an ASCII text file. To export SpectrAA current results in an ASCII report file refer to Output Data to File. This ASCII file can then be imported into a variety of programs.

The two most useful types of programs using the report file tend to be a text editor and a spreadsheet.

### Import ASCII Files Into Text Editor

A text editor would be used to create a report with headings, comments, conclusions which would need to incorporate all or part of a SpectrAA report file. Importing a file is usually very simple and the program documentation would explain how this is done. The report file can then be modified as required.

### Import ASCII Files Into Spreadsheet

A spreadsheet program is more complex to use than a text editor because numbers within the SpectrAA report would be required for numerical operations such as plotting results, trends or comparing values and so on. In an ASCII text file, there are no numerical values only strings of text which cannot be used immediately for numerical manipulations. The numbers imbedded in the text must therefore be converted to numerical values. To explain the process, Lotus 1-2-3 is used as an example but other programs such as Borland Quattro and Microsoft Excel operate in a similar manner.

Lotus 1-2-3 reads in ("imports") an ASCII file in two ways. One option is to import the file as a set of values (or numbers) using the File Import Numbers command (/FIN). The other option is to import the file as a set of labels (or text) using the File Import Text command (/FIT).

Using the /FIN command causes all non-numerical text to be stripped out and the remaining numbers loaded into individual cells. These numbers can be used immediately for calculations and graphing.

If it is necessary to retain, say, instrument parameters or sample labels then the whole file must be imported as text using the /FIT command. Each line of the ASCII file is then loaded

as a label in the first column of the spreadsheet. All numbers are still part of a text string and cannot be used immediately. To use the concentrations and absorbance values as numbers, the relevant lines must be converted using the data parse command (/DP). This breaks up or parses the text line into numbers and labels. To do this a template line must first be created. This line indicates how all the lines below are to be parsed. For SpectrAA report files printed with the each reading option, the following format line is suggested.

```
:L>>>>>>>>>>V>>>>>*****V>>***V>>>>>***V>>>>>>>>>>
***V>>>>>***V>>>>>***V>>>>>***
```

Using this template line means that parsing the SpectrAA report gives eight columns. Table 12 shows the format and the appropriate SpectrAA value of each column.

Table 12. Parsed SpectrAA Data Values

Column	Type	Width	SpectrAA value
1	Label	12	Sample labels
2	Value	6	Concentration/emission
3	Value	3	%RSD
4	Value	6	Mean absorbance/emission
5	Value	6	Readings
6	Value	6	Readings
7	Value	6	Readings
8	Value	6	Readings

## Conclusion

Operating the SpectrAA-300/400 atomic absorption spectrometer can be done with only a minimal knowledge of computers. However the SpectrAA software allows easy interaction with the PC environment if the operator wishes it. If desired, the standard SpectrAA report can be enhanced by text and numerical manipulations without compromising the original automatic run results which may be permanently archived and recalled if required.

## For More Information

For more information on our products and services, visit our Web site at [www.agilent.com/chem](http://www.agilent.com/chem)

[www.agilent.com/chem](http://www.agilent.com/chem)

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 1990  
Printed in the USA  
November 1, 2010  
AA094



**Agilent Technologies**