



Agilent Technologies

Agilent Chemstation Sequence Summary Report

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material.
The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.



Title Agilent ChemStation summary reports

After each sequence of runs a sequence summary report can be printed. Typically this is done to obtain statistical results and determine system suitability. In addition to the entries in the sequence table and before the report can be calculated and printed, several data inputs for sequence parameter and sequence output are required, see figure1.

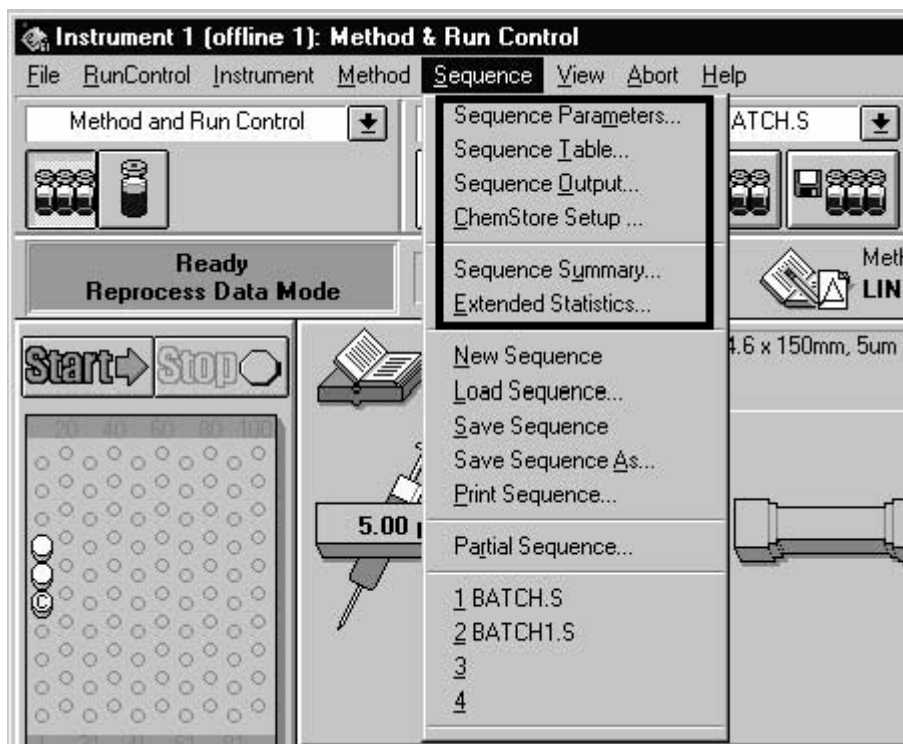


Figure 1
Entries need to be made in these sections to obtain automatically a sequence Summary report at the end of a sequence

Sequence Parameters: Instrument 1

Operator Name:

Data File

☒ Auto ☐ Prefix/Counter

Prefix: Counter:

Subdirectory:

Path:

Bar Code Reader

☐ Use In Sequence

On a bar code mismatch

☐ Inject anyway

☒ Don't inject

Part of methods to run

(after loading a new method)

Shutdown

☐ Post-Sequence Cmd / Macro

nRdy Timeout: min

Sequence Comment:

OK Cancel Help

Figure 2
Sequence parameters screen

In the *Sequence Parameters* screen (figure 2) the item *Parts of Method to Run* must be set to *According to Runtime Checklist*. This entry determines which part of a method is executed during a sequence and *According to Runtime Checklist* refers to the run-time checklist configuration that was previously edited as part of the method in order to obtain integration and quantitative results.

If data acquisition is completed and the user wants to reanalyze a sequence of samples without data acquisition, the option *Reprocessing Only* allows to recalculate the sequence summary report easily.

In the *Sequence Output* screen the report destination and the content of a sequence summary report are defined by selecting the appropriate check boxes, see figure 3.

The content of the sequence summary report is defined by the items on the right side of the screen shown in figure 3. Selecting *Setup* in the *Sequence Output* dialog box accesses this configuration screen. The sequence summary report allows a variety of information to be printed in one continuously enumerated report.

In addition to a wide selection of statistical results from sample and/or calibration runs, other items can be selected such as sample summary reports that list all

acquired samples, complete printouts of all parameters in the methods that were used, and printouts of sequence logbooks and so on.

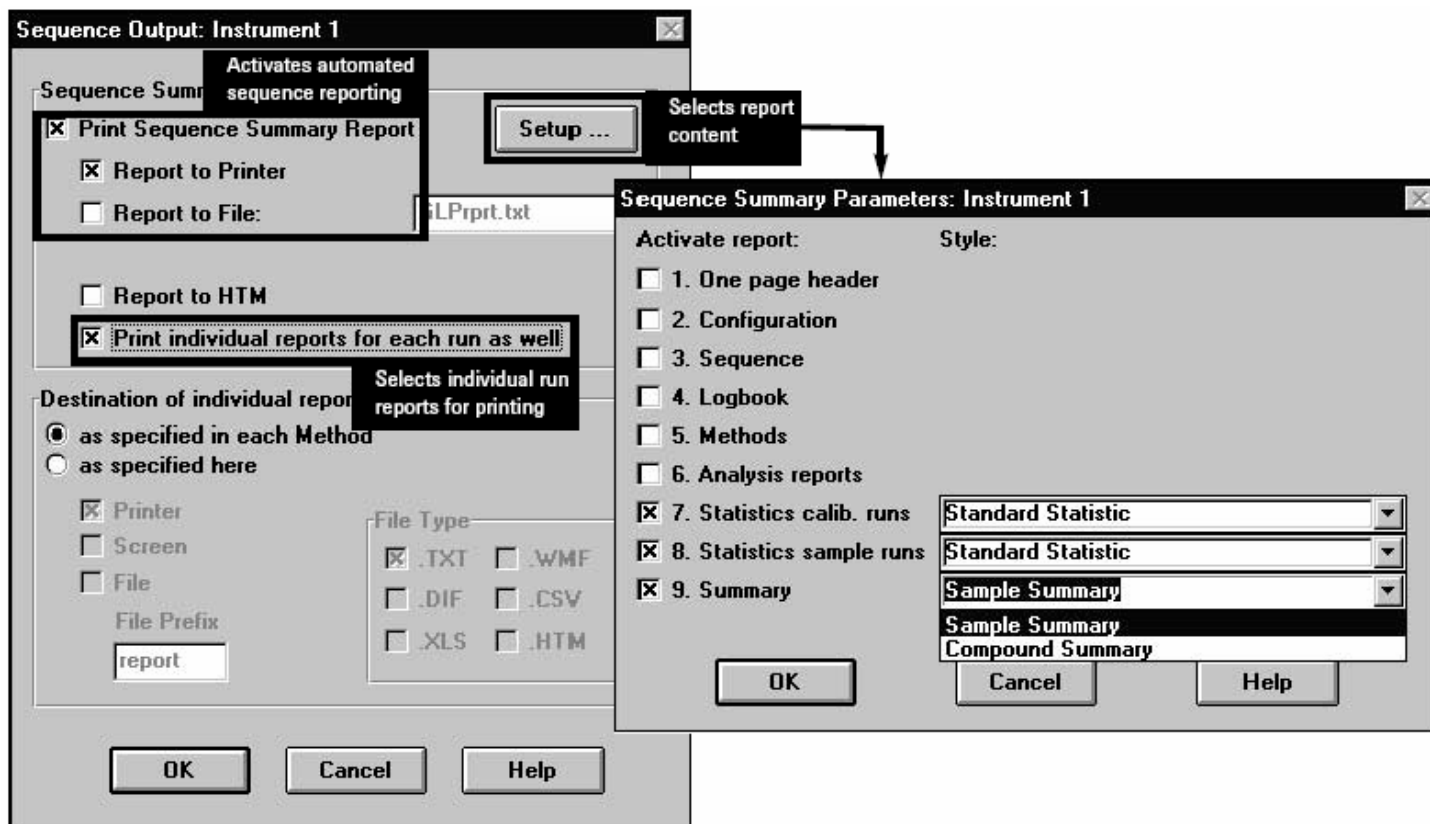


Figure 4
Selection of report destination and content of a sequence summary report

It is also possible to include the individual result reports for each run as part of the summary report instead of individual printouts after the end of each run.

The statistical evaluation of sequence runs is defined in the *Extended Statistic Parameter* screen, see figure 4. Statistical results can be obtained for all parameter shown in this dialog box. Either standard deviation or relative standard deviation or 95% confidence interval can be applied and upper/lower limits for each parameter can be specified.

A calibrated method is necessary to be able obtain statistical results.

Figure 5 shows the *Sequence Table* screen, in which it is important to ensure that the sample type is correctly set to *Sample*, *Calibration* or *Control Sample*, because statistical calculations can be selected based on sample type.

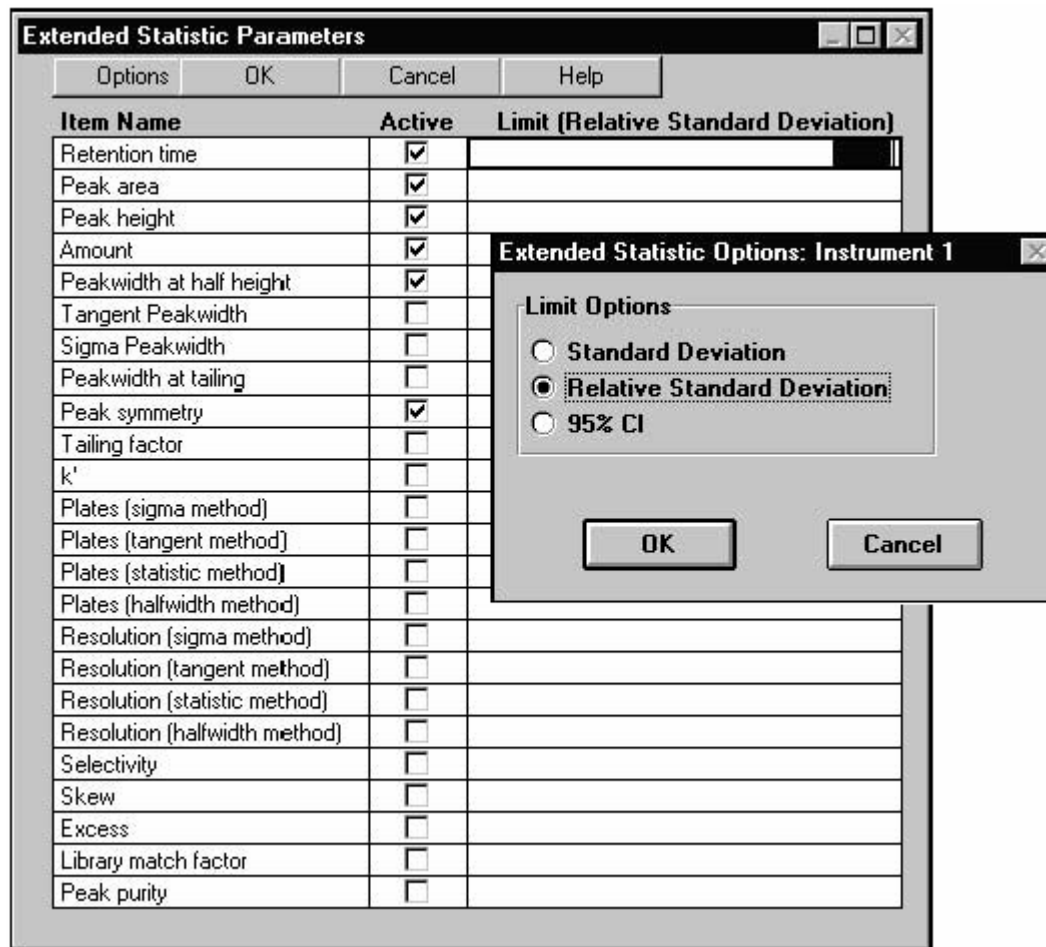


Figure 5
Setup of statistical calculations for sequence runs

Figure 6 shows an example of a sequence summary report. It contains information about the analyzed samples such as location, sample name, filename, and so on. The header includes information such as operator name, the used chromatographic method, and date of acquisition.

Sequence Table: Instrument 1

Currently Running
 Line: Method: Location: Inj:

Sample Info for P1-D-12:
 Lib105

Plate 1 ID:
 Plate 2 ID:

Line	Location	Sample Name	Method Name	Inj/Location	Sample Type	Cal Level	Update RF	Update RT	Interval	Sam
1	P1-D-12	Glucocorticoide+	96WELL	2	Sample					
2	P1-A-01	Theophyllin	96WELL	1	Sample					
3	P1-A-02	Theobromine	96WELL	1	Sample					
4	P1-A-03	Caffeine	96WELL	1	Sample					
5	P1-A-08	Phenobarbital	96WELL	1	Sample					
6	P1-A-09	Diazepam	96WELL	1	Sample					
7	P1-A-10	Paracetamol	96WELL	1	Sample					
8	P1-A-11	Phenacetin	96WELL	1	Sample					
9	P1-B-03	Antipyrin	96WELL	1	Sample					
10	P1-C-01	Minocycline	96WELL	1	Sample					
11	P1-C-02	Tetracycline	96WELL	1	Sample					
12	P1-C-03	Doxycycline	96WELL	1	Sample					
13	P1-D-01	Amoxicillin	96WELL	1	Sample					
14	P1-D-02	Ampicillin	96WELL	1	Sample					
15	P1-D-03	PenicillineG	96WELL	1	Sample					
16	P1-D-04	PenicillineV	96WELL	1	Sample					
17	P1-E-01	Triphenelamine	96WELL	1	Sample					
18	P1-E-02	Chlorpheniramine	96WELL	1	Sample					
19	P1-E-03	Promethazine	96WELL	1	Sample					
20	P1-F-01	Dextromethorphan	96WELL	1	Sample					
21	P1-F-02	Verapamil	96WELL	1	Sample					

Sample Type must be filled in appropriately as Sample, Calibration or Control

Figure 6
The Sequence Table screen

S a m p l e S u m m a r y

Sequence table: C:\HPCHEM\1\SEQUENCE\96WELL.S
 Data directory path: C:\HPCHEM\1\DATA\96WELL
 Logbook: C:\HPCHEM\1\DATA\96WELL\96WELL.LOG
 Sequence start: 10/18/2002 5:25:06 PM
 Operator: agratz

Method file name: C:\HPCHEM\1\METHODS\96WELL.M

Run #	Location #	Inj	Sample Name	Sample Amt	Multip.* [ng]	File name	Cal #	Cmp
1	P1-D-12	2	Glucocorticoide+	-	1.0000	lib10002.D	3	
2	P1-A-08	1	Phenobarbital	-	1.0000	lib10006.D	1	
3	P1-A-09	1	Diazepam	-	1.0000	lib10007.D	1	
4	P1-A-11	1	Phenacetin	-	1.0000	lib10009.D	1	
5	P1-B-03	1	Antipyrin	-	1.0000	lib10010.D	1	
6	P1-C-03	1	Doxycycline	-	1.0000	lib10013.D	1	

Figure 7
 Example of a sequence sample summary report

