
Software Update Notification



Filename: QSpec Software Update Notification for version 2.18 Doct. Version No: 2
software Issue 2.doc
Date: 14th June 1999 Author: R . M. Jones

Mastersizer (S, X, Micro, Microplus) version 2.18 Software
for use with (English) Microsoft Windows version 3.1,
Windows for Workgroups version 3.11, Windows95,
Windows98 and WindowsNT v4.0

Mastersizer S

(Part Number PSW0002/2.18)

Mastersizer X

(Part Number PSW0003/2.18)

Mastersizer Micro

(Part Number PSW0004/2.18)

Mastersizer Microplus

(Part Number PSW0005/2.18)

Contents

Mastersizer S	1
(Part Number PSW0002/2.18)	1
Mastersizer X.....	1
(Part Number PSW0003/2.18)	1
Mastersizer Micro.....	1
(Part Number PSW0004/2.18)	1
Mastersizer Microplus.....	1
(Part Number PSW0005/2.18)	1
Contents.....	2
1. Introduction.....	2
2. Instrument compatibility	2
3. Functional Features	3
3.1 Additions.....	3
3.2 Deletions.....	3
3.3 Changes.....	3
3.4 Table 1 - Bug Fixes	3
3.5 Table 2 - Known Bugs.....	3
3.6 Distribution Media	5
4. Installing Software.....	5
5. Directory Structure.....	5
7. Appendix A.....	6
8. Appendix B.....	11
9. Appendix C.....	16
10. Appendix D.....	22

1. Introduction

- 1.1** This update notification describes the version 2.18 release of Mastersizer software for Microsoft Windows and describes the software bugs addressed by this release. When ordering this version of software you will be supplied with a copy of this 'QSpec Update Notification' document and the software on floppy disks. No user manuals are supplied (as the originals remain valid and unchanged) when updating from any prior version (2.xx) of software. Issue 2 states support for Windows98 & WindowsNT v4.0

2. Instrument compatibility

- 2.1** This release requires that that the IHSD card in the Mastersizer (S, X, Micro or Microplus) instrument is fitted with the kernal EPROM version 1.1a or above and should have version 2.04 firmware or above. (Select "About" from the Help menu and click on the "Instrument connected" check-box in your current software to check this information)

The application software can only be run from a computer with a hard disc drive and will require about 7 Mbytes of free space.

The installation program will check your disk drive for the space required. If you do not have sufficient disk space, the installation programme will advise you.

The software is distributed on 5 high density 3.5" disks and comprises the following:

Installation Program: INSTALL.EXE

The Mastersizer Application: Sizer.EXE

A Presentation Generator SMG.EXE

Help File: Sizer.HLP - and various ancillary files. (See paragraph 3.5 for a detailed list)

3. Functional Features

3.1 Additions

See Changes (3.3)

3.2 Deletions

There are no deletions.

3.3 Changes

This is a maintenance issue of the Mastersizer software for Windows and, as such, the changes are limited to fixing the following bugs:

3.4 Table 1 - Bug Fixes

The following single bug fix has been carried out on this release of the software.

Malvern Software Change Request Number (SCR)	Version in which Bug was spotted	Description
160	2.15, 2.17	Small differences in the result distribution were noted and reported by two users when a sample record was re-analysed.

3.5 Table 2 - Known Bugs

The following is a list of known bugs (to date) in this and previous releases of software. The bug is generic to all four versions of product (Mastersizer S, X, Micro and Microplus) unless the entry states otherwise.

SCR Number	Description
D046-2	If kill cursors are used to reset the kill channels to zero then the checkbox in the Setup-Result Modification dialog remains checked but disabled. The table shows no channels killed but does include a "Killed Results Channels" message on the modification line.
D046-3	If Kill Result is set in the Setup-Result Modification dialog, but the checkbox is not set to enable kill values, the red cursors still appear on the result graph (the result channels are, correctly, not zeroed.)
D046-4	On the MAM5000 or MAM5002 (Mastersizer X) 45mm range lens the lowest kill cursor does not go down to zero (the bottom size band at the extreme left hand edge of the graph.)
D046-5	When changing the shape file values from within the Setup-Result Modification dialog the user must remember to save the edited values before leaving the Edit Shape Factors dialog and returning to the Setup-Modification dialog.
D046-6	On the MAF5000 (Mastersizer Micro) the set graph to auto-scaling, accumulated mode. Load DEFAULT, 1 then (separately) load DEFAULT, 3. The area where the Open File and Load Record dialog was has a different scale to the rest of the graph, (not instrument dependent).
D046-7	With 100 as the start size the 100.0 label is missing from the graph axes.
D046-9	If all spaces are filled on a notes line and the documentation is saved to the current result and the result is saved. The information display from the Open File and Load Record dialog does not show the correct separation into lines. If one character is removed from the lines then the notes are correct.

D046-10	At maximum scale (20000), the left hand graph labels (View Data and View Fit) are slightly clipped.
D046-11	Using generated sizes (auto scale, 20 bands), the histogram graph & table are updated. If you try to load LOG100.SIZ using Edit-Input Sizes, you will have to load it twice.
D046-12	Save Configuration - if no path is given then the file is saved in the Windows directory. No default extension is added.
D046-15	DTLastElectBackground\$ is the date and time of the last <u>default</u> electrical background, and is a value stored on the IHSD - not as in documentation.
D046-16	If the file is on a floppy disk and there is no disk in the drive and the user persists in cancelling the request for the floppy to be inserted then, eventually, the open file dialog will appear with all the listboxes empty. The only way to close the dialog is to type in a path such as "C:*.SAM"
D048	Graph title incorrectly re-loaded - last character missing.
D055	Auto sample dispersion unit has wrong label in Setup Hardware dialog box.
D057	Error trap for DDE Connect needed.
D058	No saving of blended results.
D059	If drive B: exists, software will not load from A: unless there is a floppy disk in B:.
D060	Possible to perform an align on an empty cell?
D061	Print settings not saved in configuration.
D062	When using File-Save As, if the file already exists, the message "File *.* already exists is displayed.
D065	Information dialog should shown all result modification settings.
D066	Distribution stats. of averaged results do not agree with averaged stats. of the components.
D067	Alert BASIC function returns 2 for both "no" and "cancel".
D068	Measure BASIC command always measures 100 experiments.
D069	Concentration value does not change when result channels killed.
D070	BASIC does not allow "\" as the last character in a string.
SCR 37-1	Erratic behaviour of error message when number of sweeps or experiments is out of range.
SCR 37-2	Sieve tables show % below size, not % above as in documentation.
SCR 37-3	DIF2002 (MSX17) or the DIF2012 (Jacuzzi) sample handling unit dialog box sometimes closes after dispersant or sample valves are opened or closed.
SCR 37-4	Bottom half of second line of text in "Printing" message box is cut off.
SCR 37-5	Manual align not permitted in measurement sequence.
SCR 37-6	"STOP" bitmap ("STOP.BMP") appears inverted.
SCR 37-7	Only 4 levels of program nesting are possible.
SCR 37-8	Setup Page BASIC command always sets up draft table printing.
SCR 37-9	With security disabled, SysSecurityLevel returns -1, not 255 as in documentation.

3.6 Distribution Media

The full complement of files is listed in the following Appendices:

For Mastersizer S (Part Number PSW0002/2.18)	Appendix A
For Mastersizer X (Part Number PSW0003/2.18)	Appendix B
For Mastersizer Micro (Part Number PSW0004/2.18)	Appendix C
For Mastersizer Microplus (Part Number PSW0005/2.18)	Appendix D

4. Installing Software

For installation of the software, refer to the appropriate user manual.

5. Directory Structure

Once this software is installed, you should have one of the following directory structures:-

For Mastersizer S

\SIZERS	Malvern application code and data files
\SIZERS\DATA	User data
\SIZERS\PAGES	Malvern and User report and table definitions
\SIZERS\PRESENT	Malvern Presentation Matrices
\SIZERS\PROGRAMS	Malvern and User measurement and analysis procedures

For Mastersizer X

\SIZERX	Malvern application code and data files
\SIZERX\DATA	User data
\SIZERX\PAGES	Malvern and User report and table definitions
\SIZERX\PRESENT	Malvern Presentation Matrices
\SIZERX\PROGRAMS	Malvern and User measurement and analysis procedures

For Mastersizer Micro

\SIZERMU	Malvern application code and data files
\SIZERMU\DATA	User data
\SIZERMU\PAGES	Malvern and User report and table definitions
\SIZERMU\PRESENT	Malvern Presentation Matrices
\SIZERMU\PROGRAMS	Malvern and User measurement and analysis

For Mastersizer Microplus

\SIZERMP	Malvern application code and data files
\SIZERMP\DATA	User data
\SIZERMP\PAGES	Malvern and User report and table definitions
\SIZERMP\PRESENT	Malvern Presentation Matrices
\SIZERMP\PROGRAMS	Malvern and User measurement and analysis

7. Appendix A

Complete listing of disks and files for version 2.18 Software
(Part Number PSW0002/2.18) for the Mastersizer S

Disk 1

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,113	02-19-98	11:14a	SIZER.INF
SIZER	<DIR>		02-19-98	4:02p	sizer
		2 file(s)			30,801 bytes

Directory of A:\sizer

.	<DIR>		02-19-98	4:02p	.
..	<DIR>		02-19-98	4:02p	..
SIZER	EXE	651,728	01-26-98	12:06p	sizer.exe
SMG	EXE	109,008	11-06-95	5:12a	smg.exe
PRESENT	DLL	48,000	11-06-95	4:52a	present.dll
MSYAN	DLL	53,920	01-26-98	11:44a	msyan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	8,358	02-19-98	11:16a	SIZER.INI
PAINT	EXE	86,528	03-11-92	8:50a	PAINT.EXE
USR_RI	INI	532	04-26-95	10:50a	USR_RI.INI
STD_RI	INI	438	09-22-95	5:20p	STD_RI.INI
DIMSZS	DD	1,040	07-14-93	12:37p	DIMSZS.DD
R0300F	SD	206	03-10-93	8:31a	R0300F.SD
R0300RF	SD	266	03-10-93	8:31a	R0300RF.SD
R1000F	SD	186	03-10-93	8:33a	R1000F.SD
LOG100	SIZ	1,420	11-14-94	8:58a	LOG100.SIZ
STND300R	SHA	2,206	11-03-95	10:25a	STND300R.SHA
STND300	SHA	1,695	11-03-95	10:25a	STND300.SHA
STND1000	SHA	1,526	11-03-95	10:25a	STND1000.SHA
CTL3D	DLL	14,448	02-08-93	4:04p	CTL3D.DLL
DEFAULT	SAM	1,536	11-11-97	9:52a	DEFAULT.SAM
CEMENTM	BSC	3,012	10-28-92	2:09p	CEMENTM.BSC
DDELINK	BSC	268	10-28-92	2:19p	DDELINK.BSC
LOGFILEH	BSC	947	10-28-92	2:25p	LOGFILEH.BSC
LOGFILEL	BSC	772	10-28-92	2:30p	LOGFILEL.BSC
LOGREMH	BSC	651	10-28-92	2:32p	LOGREMH.BSC
LOGREML	BSC	652	10-28-92	2:35p	LOGREML.BSC
PCDEMOC	BSC	1,717	10-28-92	2:44p	PCDEMOC.BSC
SETUPALL	BSC	35	12-09-91	6:54a	SETUPALL.BSC
SPRAY	BSC	1,218	11-23-92	9:28a	SPRAY.BSC
MTROMP	BSC	2,104	11-11-94	11:13a	MTROMP.BSC
8255	BSC	1,182	10-28-92	1:54p	8255.BSC
SHAPES	BSC	19,578	02-17-97	4:16p	SHAPES.BSC
1	BMP	502	01-20-92	9:25p	1.BMP
2	BMP	502	01-20-92	9:29p	2.BMP
3	BMP	502	01-20-92	9:31p	3.BMP
4	BMP	502	01-20-92	9:35p	4.BMP
5	BMP	502	01-20-92	9:36p	5.BMP
6	BMP	502	01-20-92	9:37p	6.BMP
7	BMP	502	01-20-92	9:38p	7.BMP
8	BMP	502	01-20-92	9:38p	8.BMP
ALIGN	BMP	502	11-02-91	5:14p	ALIGN.BMP
BACKGND	BMP	502	01-07-92	11:04p	BACKGND.BMP

Appendix A to QSpec Software Update Notification for version 2.18 software

CALCULA2	BMP	502	11-19-91	11:00p	CALCULA2.BMP
CALCULAT	BMP	502	11-02-91	5:23p	CALCULAT.BMP
CHECK	BMP	502	01-17-92	10:54p	CHECK.BMP
CLEARGRF	BMP	502	11-02-91	5:01p	CLEARGRF.BMP
DISKLOAD	BMP	502	11-12-91	11:00p	DISKLOAD.BMP
DISKSAVE	BMP	502	11-12-91	11:00p	DISKSAVE.BMP
DOCUMENT	BMP	502	11-03-91	11:27a	DOCUMENT.BMP
EXIT	BMP	502	04-07-93	2:48p	EXIT.BMP
GO	BMP	502	11-03-91	11:46a	GO.BMP
GRFDOWN	BMP	502	11-02-91	4:56p	GRFDOWN.BMP
GRFLEFT	BMP	502	11-02-91	4:39p	GRFLEFT.BMP
GRFRIGHT	BMP	502	11-02-91	4:37p	GRFRIGHT.BMP
GRFUP	BMP	502	11-02-91	4:55p	GRFUP.BMP
HARDWARE	BMP	502	01-08-92	8:32a	HARDWARE.BMP
HELP	BMP	502	01-17-92	10:37p	HELP.BMP
INSPECT	BMP	502	11-02-91	5:42p	INSPECT.BMP
LOADREC	BMP	502	11-03-91	12:36p	LOADREC.BMP
MIX	BMP	502	11-03-91	12:11p	MIX.BMP
OPENFIL2	BMP	502	11-19-91	10:55p	OPENFIL2.BMP
OPENFILE	BMP	502	11-03-91	12:46p	OPENFILE.BMP
PAGE	BMP	502	01-07-92	11:00p	PAGE.BMP
PRESENT	BMP	502	01-07-92	10:54p	PRESENT.BMP
PRINT	BMP	502	11-12-91	11:09p	PRINT.BMP
PROGRAM	BMP	502	11-22-91	7:53p	PROGRAM.BMP
REPORT	BMP	502	01-07-92	10:58p	REPORT.BMP
RR	BMP	1,870	02-17-97	2:55p	RR.BMP
SAMPLE	BMP	526	07-20-93	11:56a	SAMPLE.BMP
SAMPLER	BMP	502	01-07-92	11:18p	SAMPLER.BMP
SAVEREC	BMP	502	11-03-91	12:36p	SAVEREC.BMP
SEQUENC	BMP	502	11-03-91	11:40a	SEQUENC.BMP
SETUP	BMP	502	11-26-91	9:03p	SETUP.BMP
SPACE	BMP	502	05-06-93	11:22a	SPACE.BMP
STOP	BMP	502	08-03-93	5:25p	STOP.BMP
TEMPLATE	BMP	502	11-02-91	2:14p	TEMPLATE.BMP
VDATA	BMP	502	11-17-91	11:01a	VDATA.BMP
VRESULT	BMP	502	11-17-91	11:03a	VRESULT.BMP
WASH	BMP	502	11-03-91	12:06p	WASH.BMP
ZMLEFT	BMP	502	11-02-91	4:46p	ZMLEFT.BMP
ZMRIGHT	BMP	502	11-02-91	4:47p	ZMRIGHT.BMP
MDATA	PAG	6,138	01-02-96	10:18a	MDATA.PAG
MDIFF	PAG	6,791	12-15-95	9:43a	MDIFF.PAG
MFIT	PAG	7,095	12-15-95	9:41a	MFIT.PAG
MPARAM	PAG	10,997	12-15-95	9:43a	MPARAM.PAG
MRESHIST	PAG	12,112	12-15-95	9:34a	MRESHIST.PAG
MRESMEAS	PAG	11,710	12-15-95	9:30a	MRESMEAS.PAG
MRESSIVA	PAG	12,026	12-15-95	9:39a	MRESSIVA.PAG
MRESSIVB	PAG	12,035	12-15-95	9:39a	MRESSIVB.PAG
MRESSIVI	PAG	11,951	12-15-95	9:44a	MRESSIVI.PAG
MRESSTAT	PAG	12,841	12-15-95	9:38a	MRESSTAT.PAG
MSHAPE	PAG	3,388	12-14-95	5:30p	MSHAPE.PAG
MSTATS	PAG	4,359	12-15-95	9:42a	MSTATS.PAG
MTROMP	PAG	11,039	12-15-95	9:47a	MTROMP.PAG
COLOURS	PAG	510	07-21-91	10:17p	COLOURS.PAG
MDATA	REP	6,815	01-02-96	10:19a	MDATA.REP
MDIFF	REP	8,142	08-07-95	1:28p	MDIFF.REP
MFIT	REP	8,352	08-09-95	4:20p	MFIT.REP
MPARAM	REP	21,440	11-01-95	11:37a	MPARAM.REP
MRESHIST	REP	13,362	08-07-95	1:35p	MRESHIST.REP

Appendix A to QSpec Software Update Notification for version 2.18 software

```

MRESMEAS REP          12,693  08-07-95  1:39p MRESMEAS.REP
MRESSIVA REP          13,239  08-07-95  2:10p MRESSIVA.REP
MRESSIVB REP          13,308  11-01-95 10:53a MRESSIVB.REP
MRESSIVI REP          13,249  08-07-95  2:13p MRESSIVI.REP
MRESSTAT REP          14,148  08-07-95  2:19p MRESSTAT.REP
MSHAPE  REP           5,499  12-14-95  5:22p MSHAPE.REP
MSTATS  REP           4,675  02-13-95  2:59p MSTATS.REP
MTROMP  REP           14,183  08-07-95  2:22p MTROMP.REP
COLOURS REP            510  07-21-91 10:17p COLOURS.REP
HEADER  WMF           11,126  07-05-94  1:21p HEADER.WMF
SIZER   WMF            7,766  08-04-91  7:39p SIZER.WMF
IHSD204 IH           24,078  12-07-95  4:17p IHSD204.IH
      111 file(s)          1,402,218 bytes
Total files listed:
      113 file(s)          1,433,019 bytes
      3 dir(s)             4,096 bytes free

```

Disk 2

```

Directory of A:\
SIZER          <DIR>          02-19-98  4:07p sizer
      0 file(s)             0 bytes

```

```

Directory of A:\sizer
.          <DIR>          02-19-98  4:07p .
..         <DIR>          02-19-98  4:07p ..
SIZER     HLP           1,229,802  09-04-97  4:08p sizer.hlp
DE1000    SHA            1,526  02-17-97  2:54p DE1000.SHA
DE300     SHA            1,695  02-17-97  2:54p DE300.SHA
DE300RF   SHA            2,206  02-17-97  2:54p DE300RF.SHA
ASPECT    PAG           11,748  02-17-97  2:55p ASPECT.PAG
ASPECT    REP           12,791  02-17-97  2:55p ASPECT.REP
RRT       PAG            5,758  02-17-97  2:55p RRT.PAG
RRT       REP            5,758  02-17-97  2:55p RRT.REP
ASPECT2   BSC            1,456  02-17-97  2:55p ASPECT2.BSC
RORA2     BSC            2,818  02-17-97  2:55p RORA2.BSC
      10 file(s)          1,275,558 bytes
Total files listed:
      10 file(s)          1,275,558 bytes
      3 dir(s)             179,712 bytes free

```


Disk 3

Directory of A:\

```
SIZER          <DIR>          02-19-98  4:08p  sizer
                0 file(s)              0 bytes
```

Directory of A:\sizer

```
.              <DIR>          02-19-98  4:08p  .
..            <DIR>          02-19-98  4:08p  ..
3$$1         PM           111,485  07-06-94  1:06p  3$$1.PM
3$$A         PM           111,485  07-04-94  5:12p  3$$A.PM
3$$D         PM           111,485  07-04-94  5:02p  3$$D.PM
3_METAL     PM           111,485  11-15-94  12:15p  3_METAL.PM
3JAD        PM           111,485  06-24-94  11:25p  3JAD.PM
3JHD        PM           111,485  11-14-94  3:23p  3JHD.PM
3KAF        PM           111,485  06-25-94  2:48a  3KAF.PM
3KHF        PM           111,485  06-25-94  6:07a  3KHF.PM
3LAD        PM           111,485  11-14-94  4:13p  3LAD.PM
3MAD        PM           111,485  06-25-94  9:12a  3MAD.PM
3NAD        PM           111,485  06-25-94  12:16p  3NAD.PM
3NAE        PM           111,485  06-25-94  3:28p  3NAE.PM
3NDD        PM           111,485  11-14-94  5:04p  3NDD.PM
                13 file(s)        1,449,305 bytes
Total files listed:
                13 file(s)        1,449,305 bytes
                3 dir(s)          6,144 bytes free
```

Disk 4

```
Directory of A:\
SIZER          <DIR>          02-19-98  4:11p  sizer
                0 file(s)                0 bytes
```

```
Directory of A:\sizer
.              <DIR>          02-19-98  4:11p  .
..            <DIR>          02-19-98  4:11p  ..
3NHD         PM           111,485  06-25-94  6:29p  3NHD.PM
3NHE         PM           111,485  06-25-94  9:38p  3NHE.PM
3NJE         PM           111,485  11-14-94  5:57p  3NJE.PM
3OAD         PM           111,485  06-24-94  8:20p  3OAD.PM
3ODD         PM           111,485  09-23-94  10:12a 3ODD.PM
3OFD         PM           111,485  06-26-94  12:40a 3OFD.PM
3OHD         PM           111,485  06-24-94  3:50p  3OHD.PM
3OHE         PM           111,485  11-14-94  6:50p  3OHE.PM
3OJD         PM           111,485  11-14-94  7:42p  3OJD.PM
3OJE         PM           111,485  11-15-94  9:20a  3OJE.PM
3PAD         PM           111,485  06-26-94  3:46a  3PAD.PM
3PDD         PM           111,485  11-15-94  10:10a 3PDD.PM
3PFD         PM           111,485  06-26-94  6:49a  3PFD.PM
                13 file(s)        1,449,305 bytes
Total files listed:
                13 file(s)        1,449,305 bytes
                3 dir(s)          6,144 bytes free
```

Disk 5

```
Directory of A:\
SIZER          <DIR>          02-19-98  4:13p  sizer
                0 file(s)                0 bytes
```

```
Directory of A:\sizer
.              <DIR>          02-19-98  4:13p  .
..            <DIR>          02-19-98  4:13p  ..
3PHD         PM           111,485  06-26-94  9:52a  3PHD.PM
3QAA         PM           111,485  09-30-94  11:55a 3QAA.PM
3QHD         PM           111,485  10-14-94  12:57a 3QHD.PM
3RAA         PM           111,485  06-26-94  12:26p  3RAA.PM
3RHA         PM           111,485  06-26-94  2:57p  3RHA.PM
3RHD         PM           111,485  10-14-94  7:32a  3RHD.PM
3RJA         PM           111,485  06-26-94  5:35p  3RJA.PM
3SAA         PM           111,485  06-26-94  8:13p  3SAA.PM
3SHA         PM           111,485  06-26-94  10:48p 3SHA.PM
3SJA         PM           111,485  06-27-94  1:30a  3SJA.PM
3THD         PM           111,485  06-27-94  4:55a  3THD.PM
3TJD         PM           111,485  06-27-94  8:23a  3TJD.PM
3TKD         PM           111,485  11-15-94  11:16a 3TKD.PM
                13 file(s)        1,449,305 bytes
Total files listed:
                13 file(s)        1,449,305 bytes
                3 dir(s)          6,144 bytes free
```

8. Appendix B

Complete listing of disks and files for version 2.18 Software
(Part Number PSW0003/2.18) for the Mastersizer X

Disk 1

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,575	02-19-98	11:43a	SIZER.INF
SIZER	<DIR>		02-20-98	9:18a	sizer
		2 file(s)			31,263 bytes

Directory of A:\sizer

.	<DIR>		02-20-98	9:18a	.
..	<DIR>		02-20-98	9:18a	..
SIZER	EXE	651,728	01-26-98	12:06p	sizer.exe
SMG	EXE	109,008	11-06-95	5:12a	smg.exe
PRESENT	DLL	48,000	11-06-95	4:52a	present.dll
MSXAN	DLL	39,504	11-07-97	8:41a	msxan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	8,316	02-19-98	11:44a	SIZER.INI
PAINT	EXE	86,528	03-11-92	8:50a	PAINT.EXE
USR_RI	INI	533	05-30-95	2:33p	USR_RI.INI
STD_RI	INI	438	09-22-95	5:20p	STD_RI.INI
DIMSZX	DD	626	12-12-91	9:43a	DIMSZX.DD
R0045	SD	138	11-07-91	1:19p	R0045.SD
R0300	SD	138	11-07-91	1:19p	R0300.SD
R1000	SD	138	11-07-91	1:20p	R1000.SD
R0100	SD	138	11-07-91	1:19p	R0100.SD
LOG100	SIZ	1,420	06-09-95	11:23a	LOG100.SIZ
CTL3D	DLL	14,448	02-08-93	4:04p	CTL3D.DLL
DEFAULT	SAM	1,536	01-11-93	12:29p	DEFAULT.SAM
STND45	SHA	1,116	11-03-95	10:19a	STND45.SHA
STND100	SHA	1,117	11-03-95	10:19a	STND100.SHA
STND300	SHA	1,117	11-03-95	10:20a	STND300.SHA
STND1000	SHA	1,118	11-03-95	10:20a	STND1000.SHA
CEMENTM	BSC	3,012	10-28-92	2:09p	CEMENTM.BSC
DDELINK	BSC	268	10-28-92	2:19p	DDELINK.BSC
LOGFILEH	BSC	947	10-28-92	2:25p	LOGFILEH.BSC
LOGFILEL	BSC	772	10-28-92	2:30p	LOGFILEL.BSC
LOGREMH	BSC	651	10-28-92	2:32p	LOGREMH.BSC
LOGREML	BSC	652	10-28-92	2:35p	LOGREML.BSC
PCDEMOC	BSC	1,717	10-28-92	2:44p	PCDEMOC.BSC
SETUPALL	BSC	35	12-09-91	6:54a	SETUPALL.BSC
MTROMP	BSC	2,104	11-11-94	11:13a	MTROMP.BSC
SPRAY	BSC	1,218	11-23-92	9:28a	SPRAY.BSC
8255	BSC	1,182	10-28-92	1:54p	8255.BSC
SHAPEX	BSC	19,527	02-17-97	2:54p	SHAPEX.BSC
1	BMP	502	01-20-92	9:25p	1.BMP
2	BMP	502	01-20-92	9:29p	2.BMP
3	BMP	502	01-20-92	9:31p	3.BMP
4	BMP	502	01-20-92	9:35p	4.BMP
5	BMP	502	01-20-92	9:36p	5.BMP
6	BMP	502	01-20-92	9:37p	6.BMP
7	BMP	502	01-20-92	9:38p	7.BMP
8	BMP	502	01-20-92	9:38p	8.BMP

Appendix B to QSpec Software Update Notification for version 2.18 software

ALIGN	BMP	502	11-02-91	5:14p	ALIGN.BMP
BACKGND	BMP	502	01-07-92	11:04p	BACKGND.BMP
CALCULA2	BMP	502	11-19-91	11:00p	CALCULA2.BMP
CALCULAT	BMP	502	11-02-91	5:23p	CALCULAT.BMP
CHECK	BMP	502	01-17-92	10:54p	CHECK.BMP
CLEARGRF	BMP	502	11-02-91	5:01p	CLEARGRF.BMP
DISKLOAD	BMP	502	11-12-91	11:00p	DISKLOAD.BMP
DISKSAVE	BMP	502	11-12-91	11:00p	DISKSAVE.BMP
DOCUMENT	BMP	502	11-03-91	11:27a	DOCUMENT.BMP
EXIT	BMP	502	04-07-93	2:48p	EXIT.BMP
GO	BMP	502	11-03-91	11:46a	GO.BMP
GRFDOWN	BMP	502	11-02-91	4:56p	GRFDOWN.BMP
GRFLEFT	BMP	502	11-02-91	4:39p	GRFLEFT.BMP
GRFRIGHT	BMP	502	11-02-91	4:37p	GRFRIGHT.BMP
GRFUP	BMP	502	11-02-91	4:55p	GRFUP.BMP
HARDWARE	BMP	502	01-08-92	8:32a	HARDWARE.BMP
HELP	BMP	502	01-17-92	10:37p	HELP.BMP
INSPECT	BMP	502	11-02-91	5:42p	INSPECT.BMP
LOADREC	BMP	502	11-03-91	12:36p	LOADREC.BMP
MIX	BMP	502	11-03-91	12:11p	MIX.BMP
OPENFIL2	BMP	502	11-19-91	10:55p	OPENFIL2.BMP
OPENFILE	BMP	502	11-03-91	12:46p	OPENFILE.BMP
PAGE	BMP	502	01-07-92	11:00p	PAGE.BMP
PRESENT	BMP	502	01-07-92	10:54p	PRESENT.BMP
PRINT	BMP	502	11-12-91	11:09p	PRINT.BMP
PROGRAM	BMP	502	11-22-91	7:53p	PROGRAM.BMP
REPORT	BMP	502	01-07-92	10:58p	REPORT.BMP
RR	BMP	1,870	02-17-97	2:55p	RR.BMP
SAMPLE	BMP	526	07-20-93	11:56a	SAMPLE.BMP
SAMPLER	BMP	502	01-07-92	11:18p	SAMPLER.BMP
SAVEREC	BMP	502	11-03-91	12:36p	SAVEREC.BMP
SEQUENC	BMP	502	11-03-91	11:40a	SEQUENC.BMP
SETUP	BMP	502	11-26-91	9:03p	SETUP.BMP
SPACE	BMP	502	05-06-93	11:22a	SPACE.BMP
STOP	BMP	502	08-03-93	5:25p	STOP.BMP
TEMPLATE	BMP	502	11-02-91	2:14p	TEMPLATE.BMP
VDATA	BMP	502	11-17-91	11:01a	VDATA.BMP
VRESULT	BMP	502	11-17-91	11:03a	VRESULT.BMP
WASH	BMP	502	11-03-91	12:06p	WASH.BMP
ZMLEFT	BMP	502	11-02-91	4:46p	ZMLEFT.BMP
ZMRIGHT	BMP	502	11-02-91	4:47p	ZMRIGHT.BMP
MDATA	PAG	6,138	01-02-96	10:18a	MDATA.PAG
MDIFF	PAG	6,791	12-15-95	9:43a	MDIFF.PAG
MFIT	PAG	7,095	12-15-95	9:41a	MFIT.PAG
MPARAM	PAG	10,997	12-15-95	9:43a	MPARAM.PAG
MRESHIST	PAG	12,112	12-15-95	9:34a	MRESHIST.PAG
MRESMEAS	PAG	11,710	12-15-95	9:30a	MRESMEAS.PAG
MRESSIVA	PAG	12,026	12-15-95	9:39a	MRESSIVA.PAG
MRESSIVB	PAG	12,035	12-15-95	9:39a	MRESSIVB.PAG
MRESSIVI	PAG	11,951	12-15-95	9:44a	MRESSIVI.PAG
MRESSTAT	PAG	12,841	12-15-95	9:38a	MRESSTAT.PAG
MSHAPE	PAG	3,388	12-14-95	5:30p	MSHAPE.PAG
MSTATS	PAG	4,359	12-15-95	9:42a	MSTATS.PAG
MTROMP	PAG	11,039	12-15-95	9:47a	MTROMP.PAG
COLOURS	PAG	510	07-21-91	10:17p	COLOURS.PAG
MDATA	REP	6,815	01-02-96	10:19a	MDATA.REP
MDIFF	REP	8,142	08-07-95	1:28p	MDIFF.REP
MFIT	REP	8,352	08-09-95	4:20p	MFIT.REP

Appendix B to QSpec Software Update Notification for version 2.18 software

```

MPARAM    REP          21,440  11-01-95  11:37a  MPARAM.REP
MRESHIST  REP          13,362  08-07-95  1:35p  MRESHIST.REP
MRESMEAS  REP          12,693  08-07-95  1:39p  MRESMEAS.REP
MRESSIVA  REP          13,239  08-07-95  2:10p  MRESSIVA.REP
MRESSIVB  REP          13,308  11-01-95  10:53a  MRESSIVB.REP
MRESSIVI  REP          13,249  08-07-95  2:13p  MRESSIVI.REP
MRESSTAT  REP          14,148  08-07-95  2:19p  MRESSTAT.REP
MSHAPE    REP           5,499  12-14-95  5:22p  MSHAPE.REP
MSTATS    REP           4,675  02-13-95  2:59p  MSTATS.REP
MTROMP    REP          14,183  08-07-95  2:22p  MTROMP.REP
COLOURS   REP           510    07-21-91  10:17p  COLOURS.REP
HEADER    WMF          11,126  07-05-94  1:21p  HEADER.WMF
SIZER     WMF           7,766  08-04-91  7:39p  SIZER.WMF
IHSD204   IH          24,078  12-07-95  4:17p  IHSD204.IH
          113 file(s)      1,386,231 bytes
Total files listed:
          115 file(s)      1,417,494 bytes
           3 dir(s)        18,432 bytes free

```

Disk 2

```

Directory of A:\
SIZER     <DIR>          02-20-98  9:25a  sizer
          0 file(s)      0 bytes

```

```

Directory of A:\sizer
.         <DIR>          02-20-98  9:25a  .
..        <DIR>          02-20-98  9:25a  ..
SIZER     HLP          1,229,802  09-04-97  4:08p  sizer.hlp
          1 file(s)      1,229,802 bytes
Total files listed:
          1 file(s)      1,229,802 bytes
           3 dir(s)        227,328 bytes free

```

Disk 3

```

Directory of A:\
SIZER     <DIR>          02-20-98  9:26a  sizer
          0 file(s)      0 bytes

```

```

Directory of A:\sizer
.         <DIR>          02-20-98  9:26a  .
..        <DIR>          02-20-98  9:26a  ..
2$$1     PM           50,045  03-29-93  2:00a  2$$1.PM
2$$A     PM           50,045  02-21-92  9:52a  2$$A.PM
2$$D     PM           50,045  02-20-92  5:27p  2$$D.PM
2AKD     PM           50,045  02-18-92  2:05p  2AKD.PM
2GAD     PM           50,045  01-29-92  4:44a  2GAD.PM
2GDD     PM           50,045  03-03-92  3:52a  2GDD.PM
2GFD     PM           50,045  01-28-92  5:54a  2GFD.PM
2GHD     PM           50,045  01-28-92  5:13p  2GHD.PM
2IAD     PM           50,045  03-11-92  4:12a  2IAD.PM
2IDD     PM           50,045  03-10-92  5:42p  2IDD.PM
2IFD     PM           50,045  03-10-92  7:25a  2IFD.PM
2IHD     PM           50,045  03-09-92  9:08p  2IHD.PM
2IHF     PM           50,045  09-29-93  9:01p  2IHF.PM
2JAD     PM           50,045  01-25-92  4:47p  2JAD.PM
2JDD     PM           50,045  03-02-92  5:34p  2JDD.PM

```

Appendix B to QSpec Software Update Notification for version 2.18 software

```

2JFD      PM          50,045  02-02-92  2:26p  2JFD.PM
2JHD      PM          50,045  02-01-92  4:35a  2JHD.PM
2LAD      PM          50,045  12-15-91  2:26p  2LAD.PM
2LAF      PM          50,045  12-23-93  9:59p  2LAF.PM
2LDD      PM          50,045  02-26-92  5:24a  2LDD.PM
2LDF      PM          50,045  12-24-93  1:59a  2LDF.PM
2LFD      PM          50,045  02-02-92  3:09a  2LFD.PM
2LFF      PM          50,045  12-23-93  11:59p 2LFF.PM
2LHD      PM          50,045  02-01-92  3:52p  2LHD.PM
2LHF      PM          50,045  12-23-93  7:58p  2LHF.PM
2MAD      PM          50,045  02-04-92  5:33a  2MAD.PM
2MDD      PM          50,045  02-25-92  7:06p  2MDD.PM
2MFD      PM          50,045  02-04-92  4:51p  2MFD.PM
2MHD      PM          50,045  02-05-92  4:08a  2MHD.PM
      29 file(s)          1,451,305 bytes
Total files listed:
      29 file(s)          1,451,305 bytes
      3 dir(s)            1,536 bytes free

```

Disk 4

```

Directory of A:\
SIZER      <DIR>          02-20-98  9:28a  sizer
      0 file(s)          0 bytes

```

```

Directory of A:\sizer
.           <DIR>          02-20-98  9:28a  .
..          <DIR>          02-20-98  9:28a  ..
2NAD      PM          50,045  01-22-92  5:39a  2NAD.PM
2NDD      PM          50,045  03-04-92  6:18p  2NDD.PM
2NFD      PM          50,045  01-25-92  5:15a  2NFD.PM
2NFE      PM          50,045  07-04-92  1:37p  2NFE.PM
2NGE      PM          50,045  11-08-93  9:13p  2NGE.PM
2NHD      PM          50,045  12-14-91  3:37p  2NHD.PM
2NHE      PM          50,045  10-13-93  12:37a 2NHE.PM
2NJF      PM          50,045  11-30-93  9:46p  2NJF.PM
2OAD      PM          50,045  12-17-91  1:09a  2OAD.PM
2ODD      PM          50,045  03-04-92  5:21a  2ODD.PM
2OFD      PM          50,045  01-22-92  4:59p  2OFD.PM
2OHD      PM          50,045  12-13-91  4:02a  2OHD.PM
2OHE      PM          50,045  10-12-93  9:19p  2OHE.PM
2OID      PM          50,045  07-24-93  5:33a  2OID.PM
2OKD      PM          50,045  07-24-93  9:25a  2OKD.PM
2PAD      PM          50,045  12-13-91  3:36p  2PAD.PM
2PDD      PM          50,045  03-03-92  6:59p  2PDD.PM
2PFD      PM          50,045  02-06-92  8:09p  2PFD.PM
2PFF      PM          50,045  06-30-93  12:31a 2PFF.PM
2PHD      PM          50,045  12-15-91  2:58a  2PHD.PM
2PHF      PM          50,045  06-29-93  9:06p  2PHF.PM
2PJD      PM          50,045  07-24-93  2:24a  2PJD.PM
2PKD      PM          50,045  07-04-94  8:25p  2PKD.PM
2QAA      PM          50,045  02-19-92  2:10p  2QAA.PM
2QAD      PM          50,045  02-07-92  7:55a  2QAD.PM
2QDD      PM          50,045  03-06-92  11:07a 2QDD.PM
2QFD      PM          50,045  02-08-92  2:05p  2QFD.PM

```

Appendix B to QSpec Software Update Notification for version 2.18 software

```

2QHD      PM          50,045  12-17-91 10:51p 2QHD.PM
2QJD      PM          50,045  07-23-93 11:07p 2QJD.PM
          29 file(s)      1,451,305 bytes
Total files listed:
          29 file(s)      1,451,305 bytes
          3 dir(s)        1,536 bytes free
    
```

Disk 5

```

Directory of A:\
SIZER      <DIR>          02-20-98  9:31a  sizer
          0 file(s)      0 bytes
    
```

```

Directory of A:\sizer
.          <DIR>          02-20-98  9:31a  .
..         <DIR>          02-20-98  9:31a  ..
2RAA      PM          50,045  02-24-92  5:19p 2RAA.PM
2RAD      PM          50,045  02-09-92  2:08a 2RAD.PM
2RAE      PM          50,045  12-02-93 12:33a 2RAE.PM
2RDD      PM          50,045  03-06-92 12:33a 2RDD.PM
2RFD      PM          50,045  02-16-92 11:38p 2RFD.PM
2RHA      PM          50,045  12-15-91 11:46p 2RHA.PM
2RHD      PM          50,045  02-03-92  2:14a 2RHD.PM
2RID      PM          50,045  09-02-93  8:44p 2RID.PM
2RJA      PM          50,045  11-15-93 11:42p 2RJA.PM
2SAA      PM          50,045  01-21-92 12:13a 2SAA.PM
2SFA      PM          50,045  02-08-92  2:32a 2SFA.PM
2SFD      PM          50,045  09-28-94  7:04p 2SFD.PM
2SHA      PM          50,045  12-16-91  9:23a 2SHA.PM
2SHF      PM          50,045  06-24-93 10:03p 2SHF.PM
2SKE      PM          50,045  05-19-93 12:53p 2SKE.PM
2SKF      PM          50,045  06-29-93  1:41a 2SKF.PM
2TDD      PM          50,045  07-07-93 12:47a 2TDD.PM
2TFD      PM          50,045  07-06-93  9:12p 2TFD.PM
2THA      PM          50,045  02-20-92  3:36a 2THA.PM
2THD      PM          50,045  12-14-91  4:21a 2THD.PM
2TJD      PM          50,045  05-17-93  3:18p 2TJD.PM
2TKD      PM          50,045  02-19-92  4:51a 2TKD.PM
2UHD      PM          50,045  08-16-93  9:41p 2UHD.PM
2UJA      PM          50,045  11-15-93  8:56p 2UJA.PM
2UJD      PM          50,045  08-17-93  1:51a 2UJD.PM
2UKA      PM          50,045  11-29-93  9:13p 2UKA.PM
2VHF      PM          50,045  12-07-93  8:32p 2VHF.PM
2VJF      PM          50,045  12-07-93 11:11p 2VJF.PM
2VKF      PM          50,045  12-08-93  2:02a 2VKF.PM
          29 file(s)      1,451,305 bytes
Total files listed:
          29 file(s)      1,451,305 bytes
          3 dir(s)        1,536 bytes free
    
```

9. Appendix C

Complete listing of disks and files for version 2.18 Software
(Part Number PSW0004/2.18) for the Mastersizer Micro

Disk 1

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,852	02-19-98	11:42a	SIZER.INF
SIZER	<DIR>		02-25-98	3:45p	sizer
		2 file(s)			31,540 bytes

Directory of A:\sizer

.	<DIR>		02-25-98	3:45p	.
..	<DIR>		02-25-98	3:45p	..
SIZER	EXE	651,728	01-26-98	12:06p	sizer.exe
SMG	EXE	109,008	11-06-95	5:12a	smg.exe
PRESENT	DLL	48,000	11-06-95	4:52a	present.dll
MSYAN	DLL	47,728	11-07-97	8:49a	msyan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	8,001	02-19-98	11:42a	SIZER.INI
PAINT	EXE	86,528	03-11-92	8:50a	PAINT.EXE
USR_RI	INI	532	04-26-95	10:50a	USR_RI.INI
STD_RI	INI	413	11-08-95	11:09a	STD_RI.INI
DIMSZS	DD	1,040	07-14-93	12:37p	DIMSZS.DD
R0100RF	SD	190	11-11-93	3:47p	R0100RF.SD
LOG100	SIZ	1,420	04-25-95	12:28p	LOG100.SIZ
CTL3D	DLL	14,448	02-08-93	4:04p	CTL3D.DLL
DEFAULT	SAM	4,608	11-19-93	11:10a	DEFAULT.SAM
STANDARD	SHA	1,559	11-30-93	2:23p	STANDARD.SHA
DDELINK	BSC	268	10-28-92	2:19p	DDELINK.BSC
LOGFILEH	BSC	947	10-28-92	2:25p	LOGFILEH.BSC
LOGFILEL	BSC	772	10-28-92	2:30p	LOGFILEL.BSC
LOGREMH	BSC	651	10-28-92	2:32p	LOGREMH.BSC
LOGREML	BSC	652	10-28-92	2:35p	LOGREML.BSC
SETUPALL	BSC	35	12-09-91	6:54a	SETUPALL.BSC
MTROMP	BSC	2,104	11-11-94	11:13a	MTROMP.BSC
8255	BSC	1,182	10-28-92	1:54p	8255.BSC
PROMSAM	BSC	62	02-19-97	5:27p	PROMSAM.BSC
SHAPEM	BSC	19,284	02-17-97	2:54p	SHAPEM.BSC
1	BMP	502	01-20-92	9:25p	1.BMP
2	BMP	502	01-20-92	9:29p	2.BMP
3	BMP	502	01-20-92	9:31p	3.BMP
4	BMP	502	01-20-92	9:35p	4.BMP
5	BMP	502	01-20-92	9:36p	5.BMP
6	BMP	502	01-20-92	9:37p	6.BMP
7	BMP	502	01-20-92	9:38p	7.BMP
8	BMP	502	01-20-92	9:38p	8.BMP
ALIGN	BMP	502	11-02-91	5:14p	ALIGN.BMP
BACKGND	BMP	502	01-07-92	11:04p	BACKGND.BMP
CALCULA2	BMP	502	11-19-91	11:00p	CALCULA2.BMP
CALCULAT	BMP	502	11-02-91	5:23p	CALCULAT.BMP
CHECK	BMP	502	01-17-92	10:54p	CHECK.BMP
CLEARGRF	BMP	502	11-02-91	5:01p	CLEARGRF.BMP
DISKLOAD	BMP	502	11-12-91	11:00p	DISKLOAD.BMP
DISKSAVE	BMP	502	11-12-91	11:00p	DISKSAVE.BMP

Appendix C to QSpec Software Update Notification for version 2.18 software

DOCUMENT	BMP	502	11-03-91	11:27a	DOCUMENT.BMP
EXIT	BMP	502	04-07-93	2:48p	EXIT.BMP
GO	BMP	502	11-03-91	11:46a	GO.BMP
GRFDOWN	BMP	502	11-02-91	4:56p	GRFDOWN.BMP
GRFLEFT	BMP	502	11-02-91	4:39p	GRFLEFT.BMP
GRFRIGHT	BMP	502	11-02-91	4:37p	GRFRIGHT.BMP
GRFUP	BMP	502	11-02-91	4:55p	GRFUP.BMP
HARDWARE	BMP	502	01-08-92	8:32a	HARDWARE.BMP
HELP	BMP	502	01-17-92	10:37p	HELP.BMP
INSPECT	BMP	502	11-02-91	5:42p	INSPECT.BMP
LOADREC	BMP	502	11-03-91	12:36p	LOADREC.BMP
MIX	BMP	502	11-03-91	12:11p	MIX.BMP
OPENFIL2	BMP	502	11-19-91	10:55p	OPENFIL2.BMP
OPENFILE	BMP	502	11-03-91	12:46p	OPENFILE.BMP
PAGE	BMP	502	01-07-92	11:00p	PAGE.BMP
PRESENT	BMP	502	01-07-92	10:54p	PRESENT.BMP
PRINT	BMP	502	11-12-91	11:09p	PRINT.BMP
PROGRAM	BMP	502	11-22-91	7:53p	PROGRAM.BMP
REPORT	BMP	502	01-07-92	10:58p	REPORT.BMP
RR	BMP	1,870	02-17-97	2:55p	RR.BMP
SAMPLE	BMP	526	07-20-93	11:56a	SAMPLE.BMP
SAMPLER	BMP	502	01-07-92	11:18p	SAMPLER.BMP
SAVEREC	BMP	502	11-03-91	12:36p	SAVEREC.BMP
SEQUENC	BMP	502	11-03-91	11:40a	SEQUENC.BMP
SETUP	BMP	502	11-26-91	9:03p	SETUP.BMP
SPACE	BMP	502	05-06-93	11:22a	SPACE.BMP
STOP	BMP	502	08-03-93	5:25p	STOP.BMP
TEMPLATE	BMP	502	11-02-91	2:14p	TEMPLATE.BMP
VDATA	BMP	502	11-17-91	11:01a	VDATA.BMP
VRESULT	BMP	502	11-17-91	11:03a	VRESULT.BMP
WASH	BMP	502	11-03-91	12:06p	WASH.BMP
ZMLEFT	BMP	502	11-02-91	4:46p	ZMLEFT.BMP
ZMRIGHT	BMP	502	11-02-91	4:47p	ZMRIGHT.BMP
MDATA	PAG	6,138	01-02-96	10:18a	MDATA.PAG
MDIFF	PAG	6,791	12-15-95	9:43a	MDIFF.PAG
MFIT	PAG	7,095	12-15-95	9:41a	MFIT.PAG
MPARAM	PAG	10,997	12-15-95	9:43a	MPARAM.PAG
MRESHIST	PAG	12,112	12-15-95	9:34a	MRESHIST.PAG
MRESMEAS	PAG	11,710	12-15-95	9:30a	MRESMEAS.PAG
MRESSIVA	PAG	12,026	12-15-95	9:39a	MRESSIVA.PAG
MRESSIVB	PAG	12,035	12-15-95	9:39a	MRESSIVB.PAG
MRESSIVI	PAG	11,951	12-15-95	9:44a	MRESSIVI.PAG
MRESSTAT	PAG	12,841	12-15-95	9:38a	MRESSTAT.PAG
MSHAPE	PAG	3,388	12-14-95	5:30p	MSHAPE.PAG
MSTATS	PAG	4,359	12-15-95	9:42a	MSTATS.PAG
MTROMP	PAG	11,039	12-15-95	9:47a	MTROMP.PAG
COLOURS	PAG	510	07-21-91	10:17p	COLOURS.PAG
MDATA	REP	6,815	01-02-96	10:19a	MDATA.REP
MDIFF	REP	8,142	08-07-95	1:28p	MDIFF.REP
MFIT	REP	8,352	08-09-95	4:20p	MFIT.REP
MPARAM	REP	21,440	11-01-95	11:37a	MPARAM.REP
MRESHIST	REP	13,362	08-07-95	1:35p	MRESHIST.REP
MRESMEAS	REP	12,693	08-07-95	1:39p	MRESMEAS.REP
MRESSIVA	REP	13,239	08-07-95	2:10p	MRESSIVA.REP
MRESSIVB	REP	13,308	11-01-95	10:53a	MRESSIVB.REP
MRESSIVI	REP	13,249	08-07-95	2:13p	MRESSIVI.REP
MRESSTAT	REP	14,148	08-07-95	2:19p	MRESSTAT.REP
MSHAPE	REP	5,499	12-14-95	5:22p	MSHAPE.REP

Appendix C to QSpec Software Update Notification for version 2.18 software

```
MSTATS   REP           4,675  02-13-95  2:59p  MSTATS.REP
MTROMP   REP          14,183  08-07-95  2:22p  MTROMP.REP
COLOURS  REP             510   07-21-91 10:17p  COLOURS.REP
HEADER   WMF          11,126  07-05-94  1:21p  HEADER.WMF
SIZER    WMF           7,766  08-04-91  7:39p  SIZER.WMF
IHSD204  IH          24,078  12-07-95  4:17p  IHSD204.IH
MICRO105 IH           5,228  05-23-95  9:17a  MICRO105.IH
        106 file(s)      1,393,429 bytes
Total files listed:
        108 file(s)      1,424,969 bytes
         3 dir(s)        13,312 bytes free
```

Disk 2

```
Directory of A:\
SIZER    <DIR>          02-25-98  3:51p  sizer
        0 file(s)          0 bytes
```

```
Directory of A:\sizer
.        <DIR>          02-25-98  3:51p  .
..       <DIR>          02-25-98  3:51p  ..
SIZER    HLP          1,229,802 09-04-97  4:08p  sizer.hlp
        1 file(s)          1,229,802 bytes
Total files listed:
        1 file(s)          1,229,802 bytes
         3 dir(s)          227,328 bytes free
```

Disk 3

```
Directory of A:\
SIZER    <DIR>          02-25-98  3:51p  sizer
        0 file(s)          0 bytes
Directory of A:\sizer
.        <DIR>          02-25-98  3:51p  .
..       <DIR>          02-25-98  3:51p  ..
4$$D     PM           37,245 11-15-93 11:18a 4$$D.PM
4BHF     PM           37,245 04-24-95  1:12a 4BHF.PM
4CDF     PM           37,245 04-24-95  3:13a 4CDF.PM
4CFF     PM           37,245 04-05-95 12:56p 4CFF.PM
4CHE     PM           37,245 04-05-95  2:12p 4CHE.PM
4CHF     PM           37,245 04-05-95  3:33p 4CHF.PM
4GAD     PM           37,245 04-05-95  4:47p 4GAD.PM
4GAE     PM           37,245 04-05-95  6:05p 4GAE.PM
4GDD     PM           37,245 04-06-95  8:57a 4GDD.PM
4GFD     PM           37,245 04-06-95 10:10a 4GFD.PM
4GHD     PM           37,245 04-06-95 11:23a 4GHD.PM
4IAD     PM           37,245 04-06-95 12:36p 4IAD.PM
4IAE     PM           37,245 04-06-95  1:58p 4IAE.PM
4IDD     PM           37,245 04-06-95  3:19p 4IDD.PM
4IDF     PM           37,245 04-06-95  5:13p 4IDF.PM
4IFD     PM           37,245 04-06-95  6:26p 4IFD.PM
4IFF     PM           37,245 04-07-95  8:44a 4IFF.PM
4IHD     PM           37,245 04-07-95  9:56a 4IHD.PM
4IHE     PM           37,245 04-07-95 11:11a 4IHE.PM
4IHF     PM           37,245 04-07-95 12:30p 4IHF.PM
4JAD     PM           37,245 11-12-93  2:38a 4JAD.PM
4JDD     PM           37,245 04-07-95  1:43p 4JDD.PM
```

Appendix C to QSpec Software Update Notification for version 2.18 software

```

4JFD      PM          37,245  04-07-95  2:55p  4JFD.PM
4JHD      PM          37,245  04-07-95  4:38p  4JHD.PM
4KAF      PM          37,245  04-10-95  8:05a  4KAF.PM
4KHF      PM          37,245  04-10-95  9:26a  4KHF.PM
4LAD      PM          37,245  11-12-93  12:06a 4LAD.PM
4LAF      PM          37,245  04-10-95  10:47a 4LAF.PM
4LDD      PM          37,245  04-10-95  12:00p 4LDD.PM
4LDF      PM          37,245  04-10-95  1:20p  4LDF.PM
4LFD      PM          37,245  04-24-95  12:31p 4LFD.PM
4LFF      PM          37,245  04-10-95  3:19p  4LFF.PM
4LHD      PM          37,245  04-11-95  8:12a  4LHD.PM
4LHF      PM          37,245  04-11-95  9:31a  4LHF.PM
4MAC      PM          37,245  04-11-95  11:53a 4MAC.PM
4MAD      PM          37,245  04-11-95  1:06p  4MAD.PM
4MAE      PM          37,245  04-11-95  2:22p  4MAE.PM
4MAF      PM          37,245  04-11-95  3:43p  4MAF.PM
      38 file(s)          1,415,310 bytes
Total files listed:
      38 file(s)          1,415,310 bytes
      3 dir(s)            35,840 bytes free

```

Disk 4

```

Directory of A:\
SIZER      <DIR>          02-25-98  3:54p  sizer
      0 file(s)          0 bytes

```

```

Directory of A:\sizer
.          <DIR>          02-25-98  3:54p  .
..         <DIR>          02-25-98  3:54p  ..
4MDD      PM          37,245  04-12-95  8:01a  4MDD.PM
4MDF      PM          37,245  04-24-95  8:19a  4MDF.PM
4MFC      PM          37,245  04-24-95  8:19a  4MFC.PM
4MFD      PM          37,245  04-24-95  8:19a  4MFD.PM
4MFE      PM          37,245  04-24-95  8:19a  4MFE.PM
4MFF      PM          37,245  04-24-95  8:19a  4MFF.PM
4MHC      PM          37,245  04-24-95  8:19a  4MHC.PM
4MHD      PM          37,245  04-24-95  8:19a  4MHD.PM
4MHE      PM          37,245  04-24-95  8:19a  4MHE.PM
4MHF      PM          37,245  04-24-95  8:19a  4MHF.PM
4NAD      PM          37,245  11-12-93  1:22a  4NAD.PM
4NAE      PM          37,245  04-13-95  12:10p 4NAE.PM
4NAF      PM          37,245  04-13-95  1:31p  4NAF.PM
4NBD      PM          37,245  04-13-95  2:43p  4NBD.PM
4NDD      PM          37,245  04-13-95  3:56p  4NDD.PM
4NDE      PM          37,245  04-13-95  5:12p  4NDE.PM
4NDF      PM          37,245  04-13-95  6:32p  4NDF.PM
4NED      PM          37,245  04-18-95  6:30a  4NED.PM
4NFD      PM          37,245  04-18-95  7:43a  4NFD.PM
4NFE      PM          37,245  04-18-95  9:13a  4NFE.PM
4NFF      PM          37,245  04-18-95  10:36a 4NFF.PM
4NGD      PM          37,245  04-18-95  11:49a 4NGD.PM
4NGE      PM          37,245  04-18-95  1:05p  4NGE.PM
4NHD      PM          37,245  04-18-95  2:20p  4NHD.PM
4NHE      PM          37,245  04-18-95  4:41p  4NHE.PM
4NHF      PM          37,245  04-18-95  6:01p  4NHF.PM
4NJD      PM          37,245  04-19-95  7:15a  4NJD.PM
4NJE      PM          37,245  04-19-95  10:16a 4NJE.PM
4NJF      PM          37,245  04-19-95  1:32p  4NJF.PM

```

Appendix C to QSpec Software Update Notification for version 2.18 software

4OAD	PM	37,245	11-11-93	8:10p	4OAD.PM
4OAE	PM	37,245	04-19-95	4:27p	4OAE.PM
4ODD	PM	37,245	04-19-95	5:39p	4ODD.PM
4OED	PM	37,245	10-21-94	1:23p	4OED.PM
4OFD	PM	37,245	09-05-94	6:50p	4OFD.PM
4OGD	PM	37,245	04-20-95	6:38a	4OGD.PM
4OHD	PM	37,245	11-11-93	9:25p	4OHD.PM
4OHE	PM	37,245	04-20-95	7:53a	4OHE.PM

Appendix C to QSpec Software Update Notification for version 2.18 software

```

40ID      PM      37,245  04-20-95 10:04a 40ID.PM
          38 file(s)      1,415,310 bytes
Total files listed:
          38 file(s)      1,415,310 bytes
          3 dir(s)        35,840 bytes free
    
```

Disk 5

```

Directory of A:\
SIZER      <DIR>          02-25-98  3:57p  sizer
          0 file(s)          0 bytes
    
```

```

Directory of A:\sizer
.          <DIR>          02-25-98  3:57p  .
..         <DIR>          02-25-98  3:57p  ..
40JD      PM      37,245  04-20-95 12:39p 40JD.PM
40JE      PM      37,245  04-20-95  3:48p 40JE.PM
4OKD      PM      37,245  04-20-95  5:17p 4OKD.PM
4PAD      PM      37,245  11-11-93  6:54p 4PAD.PM
4PAE      PM      37,245  04-21-95  8:05a 4PAE.PM
4PDD      PM      37,245  04-21-95 12:11p 4PDD.PM
4PDE      PM      37,245  04-21-95  2:27p 4PDE.PM
4PED      PM      37,245  04-21-95  3:42p 4PED.PM
4PFD      PM      37,245  12-05-94  4:46p 4PFD.PM
4PFE      PM      37,245  04-21-95  8:05p 4PFE.PM
4PFF      PM      37,245  04-21-95 10:08p 4PFF.PM
4PHD      PM      37,245  05-19-94 10:52a 4PHD.PM
4PHE      PM      37,245  04-22-95 12:03a 4PHE.PM
4PHF      PM      37,245  04-22-95  2:05a 4PHF.PM
4PID      PM      37,245  04-22-95  3:56a 4PID.PM
4PJD      PM      37,245  04-22-95  5:51a 4PJD.PM
4PJE      PM      37,245  04-22-95  7:49a 4PJE.PM
4PKD      PM      37,245  04-22-95 10:05a 4PKD.PM
4QAD      PM      37,245  04-22-95 11:58a 4QAD.PM
4QAE      PM      37,245  04-22-95  1:57p 4QAE.PM
4QDD      PM      37,245  04-22-95  3:49p 4QDD.PM
4QFD      PM      37,245  04-22-95  5:41p 4QFD.PM
4QHD      PM      37,245  04-22-95  7:34p 4QHD.PM
4QHE      PM      37,245  04-22-95  9:31p 4QHE.PM
4QJD      PM      37,245  04-22-95 11:27p 4QJD.PM
4RAD      PM      37,245  04-23-95  1:23a 4RAD.PM
4RAE      PM      37,245  04-23-95  3:24a 4RAE.PM
4RFD      PM      37,245  04-23-95  5:19a 4RFD.PM
4RHD      PM      37,245  01-10-94  3:50p 4RHD.PM
4RKD      PM      37,245  04-23-95  7:36a 4RKD.PM
4SHD      PM      37,245  04-23-95  9:35a 4SHD.PM
4SKD      PM      37,245  04-23-95 11:54a 4SKD.PM
4THD      PM      37,245  04-23-95  1:58p 4THD.PM
4TJD      PM      37,245  11-11-93 10:51p 4TJD.PM
4TKD      PM      37,245  04-23-95  4:22p 4TKD.PM
4UHD      PM      37,245  04-23-95  6:33p 4UHD.PM
4UKD      PM      37,245  04-23-95  9:01p 4UKD.PM
4_METAL  PM      37,245  04-23-95 11:11p 4_METAL.PM
          38 file(s)      1,415,310 bytes
Total files listed:
          38 file(s)      1,415,310 bytes
          3 dir(s)        35,840 bytes free
    
```

10. Appendix D

Complete listing of disks and files for version 2.18 Software
(Part Number PSW0005/2.18) for the Mastersizer Microplus

Disk 1

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,884	02-19-98	11:17a	SIZER.INF
SIZER	<DIR>		02-20-98	8:59a	sizer
		2 file(s)			31,572 bytes

Directory of A:\sizer

.	<DIR>		02-20-98	8:59a	.
..	<DIR>		02-20-98	8:59a	..
SIZER	EXE	651,728	01-26-98	12:06p	sizer.exe
SMG	EXE	109,008	11-06-95	5:12a	smg.exe
PRESENT	DLL	48,000	11-06-95	4:52a	present.dll
MSYAN	DLL	53,920	01-26-98	11:44a	msyan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	7,982	02-19-98	11:39a	SIZER.INI
PAINT	EXE	86,528	03-11-92	8:50a	PAINT.EXE
USR_RI	INI	532	04-26-95	10:50a	USR_RI.INI
STD_RI	INI	413	11-08-95	11:10a	STD_RI.INI
DIMSZS	DD	1,040	07-14-93	12:37p	DIMSZS.DD
R0300RFP	SD	254	10-26-94	10:43a	R0300RFP.SD
LOG100	SIZ	1,420	03-07-95	4:50p	LOG100.SIZ
CTL3D	DLL	14,448	02-08-93	4:04p	CTL3D.DLL
DEFAULT	SAM	1,536	11-11-97	9:57a	DEFAULT.SAM
STANDARD	SHA	2,104	02-13-95	3:08p	STANDARD.SHA
DDELINK	BSC	268	10-28-92	2:19p	DDELINK.BSC
LOGFILEH	BSC	947	10-28-92	2:25p	LOGFILEH.BSC
LOGFILEL	BSC	772	10-28-92	2:30p	LOGFILEL.BSC
LOGREMH	BSC	651	10-28-92	2:32p	LOGREMH.BSC
LOGREML	BSC	652	10-28-92	2:35p	LOGREML.BSC
SETUPALL	BSC	35	12-09-91	6:54a	SETUPALL.BSC
MTROMP	BSC	2,104	11-11-94	11:13a	MTROMP.BSC
8255	BSC	1,182	10-28-92	1:54p	8255.BSC
PROMSAM	BSC	63	05-23-95	10:15a	PROMSAM.BSC
1	BMP	502	01-20-92	9:25p	1.BMP
2	BMP	502	01-20-92	9:29p	2.BMP
3	BMP	502	01-20-92	9:31p	3.BMP
4	BMP	502	01-20-92	9:35p	4.BMP
5	BMP	502	01-20-92	9:36p	5.BMP
6	BMP	502	01-20-92	9:37p	6.BMP
7	BMP	502	01-20-92	9:38p	7.BMP
8	BMP	502	01-20-92	9:38p	8.BMP
ALIGN	BMP	502	11-02-91	5:14p	ALIGN.BMP
BACKGND	BMP	502	01-07-92	11:04p	BACKGND.BMP
CALCULA2	BMP	502	11-19-91	11:00p	CALCULA2.BMP
CALCULAT	BMP	502	11-02-91	5:23p	CALCULAT.BMP
CHECK	BMP	502	01-17-92	10:54p	CHECK.BMP
CLEARGRF	BMP	502	11-02-91	5:01p	CLEARGRF.BMP
DISKLOAD	BMP	502	11-12-91	11:00p	DISKLOAD.BMP
DISKSAVE	BMP	502	11-12-91	11:00p	DISKSAVE.BMP
DOCUMENT	BMP	502	11-03-91	11:27a	DOCUMENT.BMP

Appendix D to QSpec Software Update Notification for version 2.18 software

EXIT	BMP	502	04-07-93	2:48p	EXIT.BMP
GO	BMP	502	11-03-91	11:46a	GO.BMP
GRFDOWN	BMP	502	11-02-91	4:56p	GRFDOWN.BMP
GRFLEFT	BMP	502	11-02-91	4:39p	GRFLEFT.BMP
GRFRIGHT	BMP	502	11-02-91	4:37p	GRFRIGHT.BMP
GRFUP	BMP	502	11-02-91	4:55p	GRFUP.BMP
HARDWARE	BMP	502	01-08-92	8:32a	HARDWARE.BMP
HELP	BMP	502	01-17-92	10:37p	HELP.BMP
INSPECT	BMP	502	11-02-91	5:42p	INSPECT.BMP
LOADREC	BMP	502	11-03-91	12:36p	LOADREC.BMP
MIX	BMP	502	11-03-91	12:11p	MIX.BMP
OPENFIL2	BMP	502	11-19-91	10:55p	OPENFIL2.BMP
OPENFILE	BMP	502	11-03-91	12:46p	OPENFILE.BMP
PAGE	BMP	502	01-07-92	11:00p	PAGE.BMP
PRESENT	BMP	502	01-07-92	10:54p	PRESENT.BMP
PRINT	BMP	502	11-12-91	11:09p	PRINT.BMP
PROGRAM	BMP	502	11-22-91	7:53p	PROGRAM.BMP
REPORT	BMP	502	01-07-92	10:58p	REPORT.BMP
RR	BMP	1,870	02-17-97	2:55p	RR.BMP
SAMPLE	BMP	526	07-20-93	11:56a	SAMPLE.BMP
SAMPLER	BMP	502	01-07-92	11:18p	SAMPLER.BMP
SAVEREC	BMP	502	11-03-91	12:36p	SAVEREC.BMP
SEQUENC	BMP	502	11-03-91	11:40a	SEQUENC.BMP
SETUP	BMP	502	11-26-91	9:03p	SETUP.BMP
SPACE	BMP	502	05-06-93	11:22a	SPACE.BMP
STOP	BMP	502	08-03-93	5:25p	STOP.BMP
TEMPLATE	BMP	502	11-02-91	2:14p	TEMPLATE.BMP
VDATA	BMP	502	11-17-91	11:01a	VDATA.BMP
VRESULT	BMP	502	11-17-91	11:03a	VRESULT.BMP
WASH	BMP	502	11-03-91	12:06p	WASH.BMP
ZMLEFT	BMP	502	11-02-91	4:46p	ZMLEFT.BMP
ZMRIGHT	BMP	502	11-02-91	4:47p	ZMRIGHT.BMP
MDATA	PAG	6,138	01-02-96	10:18a	MDATA.PAG
MDIFF	PAG	6,791	12-15-95	9:43a	MDIFF.PAG
MFIT	PAG	7,095	12-15-95	9:41a	MFIT.PAG
MPARAM	PAG	10,997	12-15-95	9:43a	MPARAM.PAG
MRESHIST	PAG	12,112	12-15-95	9:34a	MRESHIST.PAG
MRESMEAS	PAG	11,710	12-15-95	9:30a	MRESMEAS.PAG
MRESSIVA	PAG	12,026	12-15-95	9:39a	MRESSIVA.PAG
MRESSIVB	PAG	12,035	12-15-95	9:39a	MRESSIVB.PAG
MRESSIVI	PAG	11,951	12-15-95	9:44a	MRESSIVI.PAG
MRESSTAT	PAG	12,841	12-15-95	9:38a	MRESSTAT.PAG
MSHAPE	PAG	3,388	12-14-95	5:30p	MSHAPE.PAG
MSTATS	PAG	4,359	12-15-95	9:42a	MSTATS.PAG
MTROMP	PAG	11,039	12-15-95	9:47a	MTROMP.PAG
COLOURS	PAG	510	07-21-91	10:17p	COLOURS.PAG
MDATA	REP	6,815	01-02-96	10:19a	MDATA.REP
MDIFF	REP	8,142	08-07-95	1:28p	MDIFF.REP
MFIT	REP	8,352	08-09-95	4:20p	MFIT.REP
MPARAM	REP	21,440	11-01-95	11:37a	MPARAM.REP
MRESHIST	REP	13,362	08-07-95	1:35p	MRESHIST.REP
MRESMEAS	REP	12,693	08-07-95	1:39p	MRESMEAS.REP
MRESSIVA	REP	13,239	08-07-95	2:10p	MRESSIVA.REP
MRESSIVB	REP	13,308	11-01-95	10:53a	MRESSIVB.REP
MRESSIVI	REP	13,249	08-07-95	2:13p	MRESSIVI.REP
MRESSTAT	REP	14,148	08-07-95	2:19p	MRESSTAT.REP
MSHAPE	REP	5,499	12-14-95	5:22p	MSHAPE.REP
MSTATS	REP	4,675	02-13-95	2:59p	MSTATS.REP

Appendix D to QSpec Software Update Notification for version 2.18 software

MTROMP	REP	14,183	08-07-95	2:22p	MTROMP.REP
COLOURS	REP	510	07-21-91	10:17p	COLOURS.REP
HEADER	WMF	11,126	07-05-94	1:21p	HEADER.WMF
SIZER	WMF	7,766	08-04-91	7:39p	SIZER.WMF
IHSD204	IH	24,078	12-07-95	4:17p	IHSD204.IH
MICRO105	IH	5,228	05-23-95	9:17a	MICRO105.IH
		105 file(s)			1,377,856 bytes
Total files listed:					
		107 file(s)			1,409,428 bytes
		3 dir(s)			29,696 bytes free

Disk 2

```
Directory of A:\
SIZER          <DIR>          02-20-98  9:25a  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          02-20-98  9:25a  .
..            <DIR>          02-20-98  9:25a  ..
SIZER      HLP          1,229,802  09-04-97  4:08p  sizer.hlp
SHAPEMP    BSC           19,321    02-17-97  2:54p  SHAPEMP.BSC
                2 file(s)          1,249,123 bytes
```

```
Total files listed:
                2 file(s)          1,249,123 bytes
                3 dir(s)          207,872 bytes free
```

Disk 3

```
Directory of A:\
SIZER          <DIR>          02-20-98  9:08a  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          02-20-98  9:08a  .
..            <DIR>          02-20-98  9:08a  ..
5$$$D        PM          37,245    10-26-94  2:26p  5$$$D.PM
5BHF         PM          37,245    03-02-95  8:23p  5BHF.PM
5CDF         PM          37,245    03-02-95  9:23p  5CDF.PM
5CFF         PM          37,245    03-02-95  10:23p 5CFF.PM
5CHE         PM          37,245    03-02-95  11:19p 5CHE.PM
5CHF         PM          37,245    03-03-95  12:19a 5CHF.PM
5GAD         PM          37,245    02-28-95  3:46p  5GAD.PM
5GAE         PM          37,245    03-03-95  3:21a  5GAE.PM
5GDD         PM          37,245    02-28-95  5:31p  5GDD.PM
5GFD         PM          37,245    02-28-95  6:23p  5GFD.PM
5GHD         PM          37,245    02-28-95  7:16p  5GHD.PM
5IAD         PM          37,245    02-28-95  1:08p  5IAD.PM
5IAE         PM          37,245    03-03-95  4:19a  5IAE.PM
5IDD         PM          37,245    03-03-95  5:14a  5IDD.PM
5IDF         PM          37,245    03-03-95  6:14a  5IDF.PM
5IFD         PM          37,245    02-28-95  2:01p  5IFD.PM
5IFF         PM          37,245    03-03-95  7:13a  5IFF.PM
5IHD         PM          37,245    03-03-95  8:08a  5IHD.PM
5IHE         PM          37,245    03-03-95  9:05a  5IHE.PM
5IHF         PM          37,245    03-03-95  10:04a 5IHF.PM
5JAD         PM          37,245    02-27-95  4:01p  5JAD.PM
5JDD         PM          37,245    03-03-95  12:57p 5JDD.PM
5JFD         PM          37,245    02-28-95  2:53p  5JFD.PM
5JHD         PM          37,245    02-27-95  4:53p  5JHD.PM
5KAF         PM          37,245    02-27-95  5:51p  5KAF.PM
5KHF         PM          37,245    02-27-95  6:48p  5KHF.PM
5LAD         PM          37,245    02-27-95  7:41p  5LAD.PM
5LAF         PM          37,245    03-03-95  1:56p  5LAF.PM
5LDD         PM          37,245    03-03-95  2:49p  5LDD.PM
5LDF         PM          37,245    03-03-95  3:47p  5LDF.PM
5LFD         PM          37,245    03-03-95  4:40p  5LFD.PM
```

Appendix D to QSpec Software Update Notification for version 2.18 software

```

5LFF      PM          37,245  03-03-95  5:38p  5LFF.PM
5LHD      PM          37,245  03-03-95  6:31p  5LHD.PM
5LHF      PM          37,245  02-28-95  12:15p 5LHF.PM
5MAC      PM          37,245  03-03-95  8:27p  5MAC.PM
5MAD      PM          37,245  02-27-95  8:35p  5MAD.PM
5MAE      PM          37,245  03-03-95  9:23p  5MAE.PM
5MAF      PM          37,245  03-03-95  10:22p 5MAF.PM
      38 file(s)      1,415,310 bytes
Total files listed:
      38 file(s)      1,415,310 bytes
      3 dir(s)        35,840 bytes free
    
```

Disk 4

```

Directory of A:\
SIZER      <DIR>          02-20-98  9:11a  sizer
      0 file(s)          0 bytes
    
```

```

Directory of A:\sizer
.          <DIR>          02-20-98  9:11a  .
..         <DIR>          02-20-98  9:11a  ..
5MDD      PM          37,245  03-03-95  11:15p  5MDD.PM
5MDF      PM          37,245  03-04-95  12:12a  5MDF.PM
5MFC      PM          37,245  03-04-95  1:04a   5MFC.PM
5MFD      PM          37,245  03-04-95  1:57a   5MFD.PM
5MFE      PM          37,245  03-04-95  2:52a   5MFE.PM
5MFF      PM          37,245  03-04-95  3:50a   5MFF.PM
5MHC      PM          37,245  03-04-95  4:42a   5MHC.PM
5MHD      PM          37,245  03-04-95  5:34a   5MHD.PM
5MHE      PM          37,245  03-04-95  6:29a   5MHE.PM
5MHF      PM          37,245  03-04-95  7:27a   5MHF.PM
5NAD      PM          37,245  10-27-94  2:11a   5NAD.PM
5NAE      PM          37,245  02-27-95  9:30p   5NAE.PM
5NAF      PM          37,245  03-04-95  9:21a   5NAF.PM
5NBD      PM          37,245  03-04-95  10:14a  5NBD.PM
5NDD      PM          37,245  02-27-95  10:22p  5NDD.PM
5NDE      PM          37,245  03-04-95  11:09a  5NDE.PM
5NDF      PM          37,245  03-04-95  12:07p  5NDF.PM
5NED      PM          37,245  03-04-95  1:00p   5NED.PM
5NFD      PM          37,245  03-04-95  1:53p   5NFD.PM
5NFE      PM          37,245  03-04-95  2:48p   5NFE.PM
5NFF      PM          37,245  03-04-95  3:45p   5NFF.PM
5NGD      PM          37,245  03-04-95  4:38p   5NGD.PM
5NGE      PM          37,245  03-04-95  5:33p   5NGE.PM
5NHD      PM          37,245  01-20-95  2:00p   5NHD.PM
5NHE      PM          37,245  02-27-95  11:16p  5NHE.PM
5NHF      PM          37,245  03-04-95  6:31p   5NHF.PM
5NJD      PM          37,245  03-04-95  7:25p   5NJD.PM
5NJE      PM          37,245  02-28-95  12:11a  5NJE.PM
5NJF      PM          37,245  03-04-95  8:23p   5NJF.PM
5OAD      PM          37,245  10-26-94  8:29p   5OAD.PM
5OAE      PM          37,245  03-04-95  10:27p  5OAE.PM
5ODD      PM          37,245  02-28-95  1:03a   5ODD.PM
5OED      PM          37,245  03-04-95  11:20p  5OED.PM
5OFD      PM          37,245  10-27-94  12:48a  5OFD.PM
5OGD      PM          37,245  03-05-95  12:13a  5OGD.PM
5OHD      PM          37,245  10-26-94  7:05p   5OHD.PM
5OHE      PM          37,245  02-28-95  1:57a   5OHE.PM
    
```

Appendix D to QSpec Software Update Notification for version 2.18 software

```
5OID      PM      37,245  03-05-95  1:06a  5OID.PM
          38 file(s)      1,415,310 bytes
Total files listed:
          38 file(s)      1,415,310 bytes
          3 dir(s)        35,840 bytes free
```

Disk 5

```
Directory of A:\
SIZER          <DIR>          02-20-98  9:14a  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          02-20-98  9:14a  .
..            <DIR>          02-20-98  9:14a  ..
5OJD         PM           37,245  02-28-95  2:51a  5OJD.PM
5OJE         PM           37,245  02-28-95  3:47a  5OJE.PM
5OKD         PM           37,245  03-05-95  2:10a  5OKD.PM
5PAD         PM           37,245  10-26-94  9:53p  5PAD.PM
5PAE         PM           37,245  03-05-95  3:06a  5PAE.PM
5PDD         PM           37,245  02-28-95  4:39a  5PDD.PM
5PDE         PM           37,245  03-05-95  4:02a  5PDE.PM
5PED         PM           37,245  03-05-95  4:55a  5PED.PM
5PFD         PM           37,245  02-28-95  5:32a  5PFD.PM
5PFE         PM           37,245  03-05-95  5:50a  5PFE.PM
5PFF         PM           37,245  03-05-95  6:49a  5PFF.PM
5PHD         PM           37,245  02-28-95  6:24a  5PHD.PM
5PHE         PM           37,245  03-05-95  7:44a  5PHE.PM
5PHF         PM           37,245  03-05-95  8:42a  5PHF.PM
5PID         PM           37,245  03-05-95  9:35a  5PID.PM
5PJD         PM           37,245  02-28-95  8:10p  5PJD.PM
5PJE         PM           37,245  03-05-95  10:32a 5PJE.PM
5PKD         PM           37,245  03-05-95  11:36a 5PKD.PM
5QAD         PM           37,245  03-05-95  12:31p 5QAD.PM
5QAE         PM           37,245  03-05-95  1:28p  5QAE.PM
5QDD         PM           37,245  03-05-95  2:22p  5QDD.PM
5QFD         PM           37,245  03-05-95  3:16p  5QFD.PM
5QHD         PM           37,245  02-28-95  7:17a  5QHD.PM
5QHE         PM           37,245  03-05-95  5:08p  5QHE.PM
5QJD         PM           37,245  03-05-95  7:05p  5QJD.PM
5RAD         PM           37,245  03-05-95  8:01p  5RAD.PM
5RAE         PM           37,245  03-05-95  8:59p  5RAE.PM
5RFD         PM           37,245  03-05-95  10:51p 5RFD.PM
5RHD         PM           37,245  02-28-95  8:11a  5RHD.PM
5RKD         PM           37,245  03-06-95  5:41a  5RKD.PM
5SHD         PM           37,245  03-06-95  10:44a 5SHD.PM
5SKD         PM           37,245  03-06-95  4:52p  5SKD.PM
5THD         PM           37,245  02-28-95  9:10a  5THD.PM
5TJD         PM           37,245  10-26-94  11:26p 5TJD.PM
5TKD         PM           37,245  02-28-95  10:17a 5TKD.PM
5UHD         PM           37,245  03-07-95  1:15a  5UHD.PM
5UKD         PM           37,245  02-28-95  9:19p  5UKD.PM
5_METAL     PM           37,245  02-28-95  11:18a 5_METAL.PM
                38 file(s)          1,415,310 bytes
Total files listed:
                38 file(s)          1,415,310 bytes
                3 dir(s)           35,840 bytes free
```