
Software Update Notification



Filename: QSpec Software Update Notification for version 2.19 Doct. Version No: 1
software.doc
Date: 14th June 1999 Author: R . M. Jones

Mastersizer (S, X, Micro, Microplus) version 2.19 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 PSW0002/2.19)

Mastersizer X

(Part Number PSW0003/2.19)

Mastersizer Micro

(Part Number PSW0004/2.19)

Mastersizer Microplus

(Part Number PSW0005/2.19)

Contents

Mastersizer S	1
(Part Number PSW0002/2.19)	1
Mastersizer X.....	1
(Part Number PSW0003/2.19)	1
Mastersizer Micro.....	1
(Part Number PSW0004/2.19)	1
Mastersizer Microplus.....	1
(Part Number PSW0005/2.19)	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.....	5
3.6 Distribution Media	7
4. Installing Software.....	7
5. Directory Structure.....	7
7. Appendix A.....	8
8. Appendix B.....	13
9. Appendix C.....	19
10. Appendix D.....	25

1. Introduction

- 1.1** This update notification describes the version 2.19 release of Mastersizer software for Microsoft Windows and describes the software bugs addressed by this release. This release addresses some important on-line process control applications. When ordering this version of software you will be supplied with a copy of this 'QSpec Update Notification' document and the software on CD-ROM. (The default. Floppy disks are available if specified.) No user manuals are supplied (as the originals remain valid and unchanged) when updating from any prior version (2.xx) of software.

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 CD-ROM or 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 bug fixes were carried out on this release.

Software Change Request Number	Description
267	<p>The phone number has changed for Malvern Instruments SA</p> <p>REPLACE THE LINE Subs-Malvern Instruments SA (Orsay, France) Tel. + (33) 1 60 19 02 00 Fax + (33) 1 60 19 13 26 by Subs-Malvern Instruments SA (Orsay, France) Tel. + (33) 1 69 35 18 00 Fax + (33) 1 60 19 13 26</p>
548	<p>Change required in Help for Date\$ function</p> <p>Basic Date\$ and Year 2000 It has been pointed out by the Malvern Helpdesk that the description of the Malvern Basic Date\$ function, as described in the Mastersizer v2.18 Help, is not correct - it states that the year number 00 refers to 1900 when it could be 2000. The description should be:</p> <p>Application General function.</p> <p>Syntax Date\$</p> <p>Notes This function returns the current date as MM:DD:YY where MM is the month number, DD is the day number and YY the year number. The example below assumes that any year number less than 80 specifies a year in the 21st century</p> <pre>Months\$ = "JanFebMarAprMayJunJulAugSepOctNovDec" D\$ = Mid\$ (Date\$, 4, 2) M = Val (Left\$ (Date\$, 2)) M\$ = Mid\$ (Months\$, 3 * M - 2, 3) Y\$= Right\$ (Date\$, 2) Y = Val(Y\$) ' if year number is greater than 80 then assume year is 19xx If Y > 80 Then Fmt\$ = "Today is \$\$ \$\$\$ 19\$\$" Else Fmt\$ = "Today is \$\$ \$\$\$ 20\$\$" End If</pre>

	Print Status Using, Fmt\$, D\$, M\$, Y\$;
--	---

627	Failure of RunMode function In a process control application in a minerals process plant, the software hangs, owing to the way in which the RunMode error trapping functions.
713	Changes needed to implement revised error reporting for minerals process plant operation problem reported in SCR 713

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 PSW0002/2.19)	Appendix A
For Mastersizer X (Part Number PSW0003/2.19)	Appendix B
For Mastersizer Micro (Part Number PSW0004/2.19)	Appendix C
For Mastersizer Microplus (Part Number PSW0005/2.19)	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 2.19 Software
(Part Number PSW0002/2.19) for the Mastersizer S

Disk 1

Volume in drive A has no label
Volume Serial Number is 160A-3D34

Directory of A:\

INSTALL	EXE	26,688	28/09/95	16:04	INSTALL.EXE
SIZER	INF	4,176	04/05/99	12:28	SIZER.INF
SIZER	<DIR>		04/05/99	15:09	sizer
		2 file(s)			30,864 bytes

Directory of A:\sizer

.	<DIR>		04/05/99	15:09	.
..	<DIR>		04/05/99	15:09	..
SIZER	EXE	651,808	23/03/99	18:04	sizer.exe
SMG	EXE	109,008	06/11/95	6:12	smg.exe
PRESENT	DLL	48,000	06/11/95	5:52	present.dll
MSYAN	DLL	53,920	26/01/98	12:44	msyan.dll
SMG	HLP	45,474	19/12/94	5:11	smg.hlp
SIZER	INI	8,358	23/03/99	17:25	SIZER.INI
PAINT	EXE	86,528	11/03/92	8:50	PAINT.EXE
USR_RI	INI	532	26/04/95	10:50	USR_RI.INI
STD_RI	INI	438	22/09/95	17:20	STD_RI.INI
DIMSZS	DD	1,040	14/07/93	12:37	DIMSZS.DD
R0300F	SD	206	10/03/93	8:31	R0300F.SD
R0300RF	SD	266	10/03/93	8:31	R0300RF.SD
R1000F	SD	186	10/03/93	8:33	R1000F.SD
LOG100	SIZ	1,420	14/11/94	8:58	LOG100.SIZ
STND300R	SHA	2,206	03/11/95	10:25	STND300R.SHA
STND300	SHA	1,695	03/11/95	10:25	STND300.SHA
STND1000	SHA	1,526	03/11/95	10:25	STND1000.SHA
CTL3D	DLL	14,448	08/02/93	16:04	CTL3D.DLL
DEFAULT	SAM	1,536	11/11/97	9:52	DEFAULT.SAM
CEMENTM	BSC	3,012	28/10/92	14:09	CEMENTM.BSC
DDELINK	BSC	268	28/10/92	14:19	DDELINK.BSC
LOGFILEH	BSC	947	28/10/92	14:25	LOGFILEH.BSC
LOGFILEL	BSC	772	28/10/92	14:30	LOGFILEL.BSC
LOGREMH	BSC	651	28/10/92	14:32	LOGREMH.BSC
LOGREML	BSC	652	28/10/92	14:35	LOGREML.BSC
PCDEMOC	BSC	1,717	28/10/92	14:44	PCDEMOC.BSC
SETUPALL	BSC	35	09/12/91	6:54	SETUPALL.BSC
SPRAY	BSC	1,218	23/11/92	9:28	SPRAY.BSC
MTROMP	BSC	2,104	11/11/94	11:13	MTROMP.BSC
8255	BSC	1,182	28/10/92	13:54	8255.BSC
SHAPES	BSC	19,578	17/02/97	16:16	SHAPES.BSC
1	BMP	502	20/01/92	21:25	1.BMP
2	BMP	502	20/01/92	21:29	2.BMP
3	BMP	502	20/01/92	21:31	3.BMP
4	BMP	502	20/01/92	21:35	4.BMP
5	BMP	502	20/01/92	21:36	5.BMP

Appendix A to QSpec Software Update Notification for version 2.19 software

6	BMP	502	20/01/92	21:37	6.BMP
7	BMP	502	20/01/92	21:38	7.BMP
8	BMP	502	20/01/92	21:38	8.BMP
ALIGN	BMP	502	02/11/91	17:14	ALIGN.BMP
BACKGND	BMP	502	07/01/92	23:04	BACKGND.BMP
CALCULA2	BMP	502	19/11/91	23:00	CALCULA2.BMP
CALCULAT	BMP	502	02/11/91	17:23	CALCULAT.BMP
CHECK	BMP	502	17/01/92	22:54	CHECK.BMP
CLEARGRF	BMP	502	02/11/91	17:01	CLEARGRF.BMP
DISKLOAD	BMP	502	12/11/91	23:00	DISKLOAD.BMP
DISKSAVE	BMP	502	12/11/91	23:00	DISKSAVE.BMP
DOCUMENT	BMP	502	03/11/91	11:27	DOCUMENT.BMP
EXIT	BMP	502	07/04/93	14:48	EXIT.BMP
GO	BMP	502	03/11/91	11:46	GO.BMP
GRFDOWN	BMP	502	02/11/91	16:56	GRFDOWN.BMP
GRFLEFT	BMP	502	02/11/91	16:39	GRFLEFT.BMP
GRFRIGHT	BMP	502	02/11/91	16:37	GRFRIGHT.BMP
GRFUP	BMP	502	02/11/91	16:55	GRFUP.BMP
HARDWARE	BMP	502	08/01/92	8:32	HARDWARE.BMP
HELP	BMP	502	17/01/92	22:37	HELP.BMP
INSPECT	BMP	502	02/11/91	17:42	INSPECT.BMP
LOADREC	BMP	502	03/11/91	12:36	LOADREC.BMP
MIX	BMP	502	03/11/91	12:11	MIX.BMP
OPENFIL2	BMP	502	19/11/91	22:55	OPENFIL2.BMP
OPENFILE	BMP	502	03/11/91	12:46	OPENFILE.BMP
PAGE	BMP	502	07/01/92	23:00	PAGE.BMP
PRESENT	BMP	502	07/01/92	22:54	PRESENT.BMP
PRINT	BMP	502	12/11/91	23:09	PRINT.BMP
PROGRAM	BMP	502	22/11/91	19:53	PROGRAM.BMP
REPORT	BMP	502	07/01/92	22:58	REPORT.BMP
RR	BMP	1,870	17/02/97	14:55	RR.BMP
SAMPLE	BMP	526	20/07/93	11:56	SAMPLE.BMP
SAMPLER	BMP	502	07/01/92	23:18	SAMPLER.BMP
SAVEREC	BMP	502	03/11/91	12:36	SAVEREC.BMP
SEQUENC	BMP	502	03/11/91	11:40	SEQUENC.BMP
SETUP	BMP	502	26/11/91	21:03	SETUP.BMP
SPACE	BMP	502	06/05/93	11:22	SPACE.BMP
STOP	BMP	502	03/08/93	17:25	STOP.BMP
TEMPLATE	BMP	502	02/11/91	14:14	TEMPLATE.BMP
VDATA	BMP	502	17/11/91	11:01	VDATA.BMP
VRESULT	BMP	502	17/11/91	11:03	VRESULT.BMP
WASH	BMP	502	03/11/91	12:06	WASH.BMP
ZMLEFT	BMP	502	02/11/91	16:46	ZMLEFT.BMP
ZMRIGHT	BMP	502	02/11/91	16:47	ZMRIGHT.BMP
MDATA	PAG	6,138	02/01/96	10:18	MDATA.PAG
MDIFF	PAG	6,791	15/12/95	9:43	MDIFF.PAG
MFIT	PAG	7,095	15/12/95	9:41	MFIT.PAG
MPARAM	PAG	10,997	15/12/95	9:43	MPARAM.PAG
MRESHIST	PAG	12,112	15/12/95	9:34	MRESHIST.PAG
MRESMEAS	PAG	11,710	15/12/95	9:30	MRESMEAS.PAG
MRESSIVA	PAG	12,026	15/12/95	9:39	MRESSIVA.PAG
MRESSIVB	PAG	12,035	15/12/95	9:39	MRESSIVB.PAG
MRESSIVI	PAG	11,951	15/12/95	9:44	MRESSIVI.PAG
MRESSTAT	PAG	12,841	15/12/95	9:38	MRESSTAT.PAG
MSHAPE	PAG	3,388	14/12/95	17:30	MSHAPE.PAG
MSTATS	PAG	4,359	15/12/95	9:42	MSTATS.PAG
MTROMP	PAG	11,039	15/12/95	9:47	MTROMP.PAG
COLOURS	PAG	510	21/07/91	22:17	COLOURS.PAG

Appendix A to QSpec Software Update Notification for version 2.19 software

```

MDATA      REP          6,815  02/01/96  10:19  MDATA.REP
MDIFF      REP          8,142  07/08/95  13:28  MDIFF.REP
MFIT       REP          8,352  09/08/95  16:20  MFIT.REP
MPARAM     REP         21,440  01/11/95  11:37  MPARAM.REP
MRESHIST  REP         13,362  07/08/95  13:35  MRESHIST.REP
MRESMEAS  REP         12,693  07/08/95  13:39  MRESMEAS.REP
MRESSIVA  REP         13,239  07/08/95  14:10  MRESSIVA.REP
MRESSIVB  REP         13,308  01/11/95  10:53  MRESSIVB.REP
MRESSIVI  REP         13,249  07/08/95  14:13  MRESSIVI.REP
MRESSTAT  REP         14,148  07/08/95  14:19  MRESSTAT.REP
MSHAPE    REP          5,499  14/12/95  17:22  MSHAPE.REP
MSTATS    REP          4,675  13/02/95  14:59  MSTATS.REP
MTROMP    REP         14,183  07/08/95  14:22  MTROMP.REP
COLOURS   REP           510   21/07/91  22:17  COLOURS.REP
HEADER    WMF         11,126  05/07/94  13:21  HEADER.WMF
SIZER     WMF          7,766  04/08/91  19:39  SIZER.WMF
IHSD204   IH         24,078  07/12/95  16:17  IHSD204.IH
          111 file(s)      1,402,298 bytes
    
```

Total files listed:

```

          113 file(s)      1,433,162 bytes
           3 dir(s)       3,584 bytes free
    
```

Disk 2

Volume in drive A has no label
Volume Serial Number is 16CF-1710

Directory of A:\

```

SIZER      <DIR>          04/05/99  15:14  sizer
          0 file(s)          0 bytes
    
```

Directory of A:\sizer

```

.          <DIR>          04/05/99  15:14  .
..         <DIR>          04/05/99  15:14  ..
SIZER     HLP         1,228,717  23/03/99  17:25  sizer.hlp
DE1000    SHA          1,526   17/02/97  14:54  DE1000.SHA
DE300     SHA          1,695   17/02/97  14:54  DE300.SHA
DE300RF   SHA          2,206   17/02/97  14:54  DE300RF.SHA
ASPECT    PAG         11,748   17/02/97  14:55  ASPECT.PAG
ASPECT    REP         12,791   17/02/97  14:55  ASPECT.REP
RRT       PAG          5,758   17/02/97  14:55  RRT.PAG
RRT       REP          5,758   17/02/97  14:55  RRT.REP
ASPECT2   BSC          1,456   17/02/97  14:55  ASPECT2.BSC
RORA2     BSC          2,818   17/02/97  14:55  RORA2.BSC
ALI1000   BSC          5,864   29/04/99   9:28  ALI1000.BSC
          11 file(s)      1,280,337 bytes
    
```

Total files listed:

```

          11 file(s)      1,280,337 bytes
           3 dir(s)      174,592 bytes free
    
```

Disk 3

Volume in drive A has no label
Volume Serial Number is 16CF-2C50

Directory of A:\

```
SIZER          <DIR>          04/05/99  15:15  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          04/05/99  15:15  .
..            <DIR>          04/05/99  15:15  ..
3$$1         PM          111,485  06/07/94  13:06  3$$1.PM
3$$A         PM          111,485  04/07/94  17:12  3$$A.PM
3$$D         PM          111,485  04/07/94  17:02  3$$D.PM
3_METAL     PM          111,485  15/11/94  12:15  3_METAL.PM
3JAD        PM          111,485  24/06/94  23:25  3JAD.PM
3JHD        PM          111,485  14/11/94  15:23  3JHD.PM
3KAF        PM          111,485  25/06/94   2:48  3KAF.PM
3KHF        PM          111,485  25/06/94   6:07  3KHF.PM
3LAD        PM          111,485  14/11/94  16:13  3LAD.PM
3MAD        PM          111,485  25/06/94   9:12  3MAD.PM
3NAD        PM          111,485  25/06/94  12:16  3NAD.PM
3NAE        PM          111,485  25/06/94  15:28  3NAE.PM
3NDD        PM          111,485  14/11/94  17:04  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

Volume in drive A has no label
Volume Serial Number is 16D0-0611

Directory of A:\

```
SIZER          <DIR>          04/05/99  15:17  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          04/05/99  15:17  .
..            <DIR>          04/05/99  15:17  ..
3NHD        PM          111,485  25/06/94  18:29  3NHD.PM
3NHE        PM          111,485  25/06/94  21:38  3NHE.PM
3NJE        PM          111,485  14/11/94  17:57  3NJE.PM
3OAD        PM          111,485  24/06/94  20:20  3OAD.PM
3ODD        PM          111,485  23/09/94  10:12  3ODD.PM
3OFD        PM          111,485  26/06/94   0:40  3OFD.PM
3OHD        PM          111,485  24/06/94  15:50  3OHD.PM
3OHE        PM          111,485  14/11/94  18:50  3OHE.PM
3OJD        PM          111,485  14/11/94  19:42  3OJD.PM
3OJE        PM          111,485  15/11/94   9:20  3OJE.PM
```

Appendix A to QSpec Software Update Notification for version 2.19 software

```
3PAD      PM      111,485  26/06/94   3:46  3PAD.PM
3PDD      PM      111,485  15/11/94  10:10  3PDD.PM
3PFD      PM      111,485  26/06/94   6:49  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

Volume in drive A has no label
Volume Serial Number is 16D0-1B04

Directory of A:\

```
SIZER      <DIR>      04/05/99  15:19  sizer
      0 file(s)      0 bytes
```

Directory of A:\sizer

```
.          <DIR>      04/05/99  15:19  .
..         <DIR>      04/05/99  15:19  ..
3PHD      PM      111,485  26/06/94   9:52  3PHD.PM
3QAA      PM      111,485  30/09/94  11:55  3QAA.PM
3QHD      PM      111,485  14/10/94   0:57  3QHD.PM
3RAA      PM      111,485  26/06/94  12:26  3RAA.PM
3RHA      PM      111,485  26/06/94  14:57  3RHA.PM
3RHD      PM      111,485  14/10/94   7:32  3RHD.PM
3RJA      PM      111,485  26/06/94  17:35  3RJA.PM
3SAA      PM      111,485  26/06/94  20:13  3SAA.PM
3SHA      PM      111,485  26/06/94  22:48  3SHA.PM
3SJA      PM      111,485  27/06/94   1:30  3SJA.PM
3THD      PM      111,485  27/06/94   4:55  3THD.PM
3TJD      PM      111,485  27/06/94   8:23  3TJD.PM
3TKD      PM      111,485  15/11/94  11:16  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 2.19 Software
(Part Number PSW0003/2.19) for the Mastersizer X

Disk 1

Volume in drive A has no label
Volume Serial Number is 15DF-371E

Directory of A:\

INSTALL	EXE	26,688	28/09/95	16:04	INSTALL.EXE
SIZER	INF	4,636	04/05/99	12:34	SIZER.INF
SIZER	<DIR>		04/05/99	14:18	sizer
		2 file(s)			31,324 bytes

Directory of A:\sizer

.	<DIR>		04/05/99	14:18	.
..	<DIR>		04/05/99	14:18	..
SIZER	EXE	651,808	23/03/99	18:04	sizer.exe
SMG	EXE	109,008	06/11/95	6:12	smg.exe
PRESENT	DLL	48,000	06/11/95	5:52	present.dll
MSXAN	DLL	39,504	07/11/97	9:41	msxan.dll
SMG	HLP	45,474	19/12/94	5:11	smg.hlp
SIZER	INI	8,316	24/03/99	9:11	SIZER.INI
PAINT	EXE	86,528	11/03/92	8:50	PAINT.EXE
USR_RI	INI	533	30/05/95	14:33	USR_RI.INI
STD_RI	INI	438	22/09/95	17:20	STD_RI.INI
DIMSZX	DD	626	12/12/91	9:43	DIMSZX.DD
R0045	SD	138	07/11/91	13:19	R0045.SD
R0300	SD	138	07/11/91	13:19	R0300.SD
R1000	SD	138	07/11/91	13:20	R1000.SD
R0100	SD	138	07/11/91	13:19	R0100.SD
LOG100	SIZ	1,420	09/06/95	11:23	LOG100.SIZ
CTL3D	DLL	14,448	08/02/93	16:04	CTL3D.DLL
DEFAULT	SAM	1,536	11/01/93	12:29	DEFAULT.SAM
STND45	SHA	1,116	03/11/95	10:19	STND45.SHA
STND100	SHA	1,117	03/11/95	10:19	STND100.SHA
STND300	SHA	1,117	03/11/95	10:20	STND300.SHA
STND1000	SHA	1,118	03/11/95	10:20	STND1000.SHA
CEMENTM	BSC	3,012	28/10/92	14:09	CEMENTM.BSC
DDELINK	BSC	268	28/10/92	14:19	DDELINK.BSC
LOGFILEH	BSC	947	28/10/92	14:25	LOGFILEH.BSC
LOGFILEL	BSC	772	28/10/92	14:30	LOGFILEL.BSC
LOGREMH	BSC	651	28/10/92	14:32	LOGREMH.BSC
LOGREML	BSC	652	28/10/92	14:35	LOGREML.BSC
PCDEMOC	BSC	1,717	28/10/92	14:44	PCDEMOC.BSC
SETUPALL	BSC	35	09/12/91	6:54	SETUPALL.BSC
MTROMP	BSC	2,104	11/11/94	11:13	MTROMP.BSC
SPRAY	BSC	1,218	23/11/92	9:28	SPRAY.BSC
8255	BSC	1,182	28/10/92	13:54	8255.BSC
SHAPEX	BSC	19,527	17/02/97	14:54	SHAPEX.BSC
1	BMP	502	20/01/92	21:25	1.BMP
2	BMP	502	20/01/92	21:29	2.BMP
3	BMP	502	20/01/92	21:31	3.BMP

Appendix B to QSpec Software Update Notification for version 2.19 software

4	BMP	502	20/01/92	21:35	4.BMP
5	BMP	502	20/01/92	21:36	5.BMP
6	BMP	502	20/01/92	21:37	6.BMP
7	BMP	502	20/01/92	21:38	7.BMP
8	BMP	502	20/01/92	21:38	8.BMP
ALIGN	BMP	502	02/11/91	17:14	ALIGN.BMP
BACKGND	BMP	502	07/01/92	23:04	BACKGND.BMP
CALCULA2	BMP	502	19/11/91	23:00	CALCULA2.BMP
CALCULAT	BMP	502	02/11/91	17:23	CALCULAT.BMP
CHECK	BMP	502	17/01/92	22:54	CHECK.BMP
CLEARGRF	BMP	502	02/11/91	17:01	CLEARGRF.BMP
DISKLOAD	BMP	502	12/11/91	23:00	DISKLOAD.BMP
DISKSAVE	BMP	502	12/11/91	23:00	DISKSAVE.BMP
DOCUMENT	BMP	502	03/11/91	11:27	DOCUMENT.BMP
EXIT	BMP	502	07/04/93	14:48	EXIT.BMP
GO	BMP	502	03/11/91	11:46	GO.BMP
GRFDOWN	BMP	502	02/11/91	16:56	GRFDOWN.BMP
GRFLEFT	BMP	502	02/11/91	16:39	GRFLEFT.BMP
GRFRIGHT	BMP	502	02/11/91	16:37	GRFRIGHT.BMP
GRFUP	BMP	502	02/11/91	16:55	GRFUP.BMP
HARDWARE	BMP	502	08/01/92	8:32	HARDWARE.BMP
HELP	BMP	502	17/01/92	22:37	HELP.BMP
INSPECT	BMP	502	02/11/91	17:42	INSPECT.BMP
LOADREC	BMP	502	03/11/91	12:36	LOADREC.BMP
MIX	BMP	502	03/11/91	12:11	MIX.BMP
OPENFIL2	BMP	502	19/11/91	22:55	OPENFIL2.BMP
OPENFILE	BMP	502	03/11/91	12:46	OPENFILE.BMP
PAGE	BMP	502	07/01/92	23:00	PAGE.BMP
PRESENT	BMP	502	07/01/92	22:54	PRESENT.BMP
PRINT	BMP	502	12/11/91	23:09	PRINT.BMP
PROGRAM	BMP	502	22/11/91	19:53	PROGRAM.BMP
REPORT	BMP	502	07/01/92	22:58	REPORT.BMP
RR	BMP	1,870	17/02/97	14:55	RR.BMP
SAMPLE	BMP	526	20/07/93	11:56	SAMPLE.BMP
SAMPLER	BMP	502	07/01/92	23:18	SAMPLER.BMP
SAVEREC	BMP	502	03/11/91	12:36	SAVEREC.BMP
SEQUENC	BMP	502	03/11/91	11:40	SEQUENC.BMP
SETUP	BMP	502	26/11/91	21:03	SETUP.BMP
SPACE	BMP	502	06/05/93	11:22	SPACE.BMP
STOP	BMP	502	03/08/93	17:25	STOP.BMP
TEMPLATE	BMP	502	02/11/91	14:14	TEMPLATE.BMP
VDATA	BMP	502	17/11/91	11:01	VDATA.BMP
VRESULT	BMP	502	17/11/91	11:03	VRESULT.BMP
WASH	BMP	502	03/11/91	12:06	WASH.BMP
ZMLEFT	BMP	502	02/11/91	16:46	ZMLEFT.BMP
ZMRIGHT	BMP	502	02/11/91	16:47	ZMRIGHT.BMP
MDATA	PAG	6,138	02/01/96	10:18	MDATA.PAG
MDIFF	PAG	6,791	15/12/95	9:43	MDIFF.PAG
MFIT	PAG	7,095	15/12/95	9:41	MFIT.PAG
MPARAM	PAG	10,997	15/12/95	9:43	MPARAM.PAG
MRESHIST	PAG	12,112	15/12/95	9:34	MRESHIST.PAG
MRESMEAS	PAG	11,710	15/12/95	9:30	MRESMEAS.PAG
MRESSIVA	PAG	12,026	15/12/95	9:39	MRESSIVA.PAG
MRESSIVB	PAG	12,035	15/12/95	9:39	MRESSIVB.PAG
MRESSIVI	PAG	11,951	15/12/95	9:44	MRESSIVI.PAG
MRESSTAT	PAG	12,841	15/12/95	9:38	MRESSTAT.PAG
MSHAPE	PAG	3,388	14/12/95	17:30	MSHAPE.PAG
MSTATS	PAG	4,359	15/12/95	9:42	MSTATS.PAG

Appendix B to QSpec Software Update Notification for version 2.19 software

MTROMP	PAG	11,039	15/12/95	9:47	MTROMP.PAG
COLOURS	PAG	510	21/07/91	22:17	COLOURS.PAG
MDATA	REP	6,815	02/01/96	10:19	MDATA.REP
MDIFF	REP	8,142	07/08/95	13:28	MDIFF.REP
MFIT	REP	8,352	09/08/95	16:20	MFIT.REP
MPARAM	REP	21,440	01/11/95	11:37	MPARAM.REP
MRESHIST	REP	13,362	07/08/95	13:35	MRESHIST.REP
MRESMEAS	REP	12,693	07/08/95	13:39	MRESMEAS.REP
MRESSIVA	REP	13,239	07/08/95	14:10	MRESSIVA.REP
MRESSIVB	REP	13,308	01/11/95	10:53	MRESSIVB.REP
MRESSIVI	REP	13,249	07/08/95	14:13	MRESSIVI.REP
MRESSTAT	REP	14,148	07/08/95	14:19	MRESSTAT.REP
MSHAPE	REP	5,499	14/12/95	17:22	MSHAPE.REP
MSTATS	REP	4,675	13/02/95	14:59	MSTATS.REP
MTROMP	REP	14,183	07/08/95	14:22	MTROMP.REP
COLOURS	REP	510	21/07/91	22:17	COLOURS.REP
HEADER	WMF	11,126	05/07/94	13:21	HEADER.WMF
SIZER	WMF	7,766	04/08/91	19:39	SIZER.WMF
IHSD204	IH	24,078	07/12/95	16:17	IHSD204.IH
		113 file(s)			1,386,311 bytes

Total files listed:

115 file(s)	1,417,635 bytes
3 dir(s)	17,408 bytes free

Disk 2

Volume in drive A has no label
Volume Serial Number is 15E0-231C

Directory of A:\

SIZER	<DIR>	04/05/99	14:23	sizer
	0 file(s)			0 bytes

Directory of A:\sizer

.	<DIR>	04/05/99	14:23	.
..	<DIR>	04/05/99	14:23	..
SIZER	HLP	1,228,717	23/03/99	17:25 sizer.hlp
ALI1000	BSC	5,864	29/04/99	9:28 ALI1000.BSC
		2 file(s)		1,234,581 bytes

Total files listed:

2 file(s)	1,234,581 bytes
3 dir(s)	222,208 bytes free

Disk 3

Volume in drive A has no label
 Volume Serial Number is 1505-3513

Directory of A:\

```
SIZER          <DIR>          04/05/99  14:24  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          04/05/99  14:24  .
..            <DIR>          04/05/99  14:24  ..
2$$1         PM           50,045  29/03/93   2:00  2$$1.PM
2$$A         PM           50,045  21/02/92   9:52  2$$A.PM
2$$D         PM           50,045  20/02/92  17:27  2$$D.PM
2AKD         PM           50,045  18/02/92  14:05  2AKD.PM
2GAD         PM           50,045  29/01/92   4:44  2GAD.PM
2GDD         PM           50,045  03/03/92   3:52  2GDD.PM
2GFD         PM           50,045  28/01/92   5:54  2GFD.PM
2GHD         PM           50,045  28/01/92  17:13  2GHD.PM
2IAD         PM           50,045  11/03/92   4:12  2IAD.PM
2IDD         PM           50,045  10/03/92  17:42  2IDD.PM
2IFD         PM           50,045  10/03/92   7:25  2IFD.PM
2IHD         PM           50,045  09/03/92  21:08  2IHD.PM
2IHF         PM           50,045  29/09/93  21:01  2IHF.PM
2JAD         PM           50,045  25/01/92  16:47  2JAD.PM
2JDD         PM           50,045  02/03/92  17:34  2JDD.PM
2JFD         PM           50,045  02/02/92  14:26  2JFD.PM
2JHD         PM           50,045  01/02/92   4:35  2JHD.PM
2LAD         PM           50,045  15/12/91  14:26  2LAD.PM
2LAF         PM           50,045  23/12/93  21:59  2LAF.PM
2LDD         PM           50,045  26/02/92   5:24  2LDD.PM
2LDF         PM           50,045  24/12/93   1:59  2LDF.PM
2LFD         PM           50,045  02/02/92   3:09  2LFD.PM
2LFF         PM           50,045  23/12/93  23:59  2LFF.PM
2LHD         PM           50,045  01/02/92  15:52  2LHD.PM
2LHF         PM           50,045  23/12/93  19:58  2LHF.PM
2MAD         PM           50,045  04/02/92   5:33  2MAD.PM
2MDD         PM           50,045  25/02/92  19:06  2MDD.PM
2MFD         PM           50,045  04/02/92  16:51  2MFD.PM
2MHD         PM           50,045  05/02/92   4:08  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

Volume in drive A has no label
 Volume Serial Number is 1506-0F15

Directory of A:\

```
SIZER          <DIR>          04/05/99  14:26  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          04/05/99  14:26  .
..            <DIR>          04/05/99  14:26  ..
2NAD          PM           50,045  22/01/92   5:39  2NAD.PM
2NDD          PM           50,045  04/03/92  18:18  2NDD.PM
2NFD          PM           50,045  25/01/92   5:15  2NFD.PM
2NFE          PM           50,045  04/07/92  13:37  2NFE.PM
2NGE          PM           50,045  08/11/93  21:13  2NGE.PM
2NHD          PM           50,045  14/12/91  15:37  2NHD.PM
2NHE          PM           50,045  13/10/93   0:37  2NHE.PM
2NJF          PM           50,045  30/11/93  21:46  2NJF.PM
2OAD          PM           50,045  17/12/91   1:09  2OAD.PM
2ODD          PM           50,045  04/03/92   5:21  2ODD.PM
2OFD          PM           50,045  22/01/92  16:59  2OFD.PM
2OHD          PM           50,045  13/12/91   4:02  2OHD.PM
2OHE          PM           50,045  12/10/93  21:19  2OHE.PM
2OID          PM           50,045  24/07/93   5:33  2OID.PM
2OKD          PM           50,045  24/07/93   9:25  2OKD.PM
2PAD          PM           50,045  13/12/91  15:36  2PAD.PM
2PDD          PM           50,045  03/03/92  18:59  2PDD.PM
2PFD          PM           50,045  06/02/92  20:09  2PFD.PM
2PFF          PM           50,045  30/06/93   0:31  2PFF.PM
2PHD          PM           50,045  15/12/91   2:58  2PHD.PM
2PHF          PM           50,045  29/06/93  21:06  2PHF.PM
2PJD          PM           50,045  24/07/93   2:24  2PJD.PM
2PKD          PM           50,045  04/07/94  20:25  2PKD.PM
2QAA          PM           50,045  19/02/92  14:10  2QAA.PM
2QAD          PM           50,045  07/02/92   7:55  2QAD.PM
2QDD          PM           50,045  06/03/92  11:07  2QDD.PM
2QFD          PM           50,045  08/02/92  14:05  2QFD.PM
2QHD          PM           50,045  17/12/91  22:51  2QHD.PM
2QJD          PM           50,045  23/07/93  23:07  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

Volume in drive A has no label
Volume Serial Number is 1506-244A

Directory of A:\

```
SIZER          <DIR>          04/05/99  14:27  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          04/05/99  14:27  .
..            <DIR>          04/05/99  14:27  ..
2RAA        PM           50,045  24/02/92  17:19  2RAA.PM
2RAD        PM           50,045  09/02/92   2:08  2RAD.PM
2RAE        PM           50,045  02/12/93   0:33  2RAE.PM
2RDD        PM           50,045  06/03/92   0:33  2RDD.PM
2RFD        PM           50,045  16/02/92  23:38  2RFD.PM
2RHA        PM           50,045  15/12/91  23:46  2RHA.PM
2RHD        PM           50,045  03/02/92   2:14  2RHD.PM
2RID        PM           50,045  02/09/93  20:44  2RID.PM
2RJA        PM           50,045  15/11/93  23:42  2RJA.PM
2SAA        PM           50,045  21/01/92   0:13  2SAA.PM
2SFA        PM           50,045  08/02/92   2:32  2SFA.PM
2SFD        PM           50,045  28/09/94  19:04  2SFD.PM
2SHA        PM           50,045  16/12/91   9:23  2SHA.PM
2SHF        PM           50,045  24/06/93  22:03  2SHF.PM
2SKE        PM           50,045  19/05/93  12:53  2SKE.PM
2SKF        PM           50,045  29/06/93   1:41  2SKF.PM
2TDD        PM           50,045  07/07/93   0:47  2TDD.PM
2TFD        PM           50,045  06/07/93  21:12  2TFD.PM
2THA        PM           50,045  20/02/92   3:36  2THA.PM
2THD        PM           50,045  14/12/91   4:21  2THD.PM
2TJD        PM           50,045  17/05/93  15:18  2TJD.PM
2TKD        PM           50,045  19/02/92   4:51  2TKD.PM
2UHD        PM           50,045  16/08/93  21:41  2UHD.PM
2UJA        PM           50,045  15/11/93  20:56  2UJA.PM
2UJD        PM           50,045  17/08/93   1:51  2UJD.PM
2UKA        PM           50,045  29/11/93  21:13  2UKA.PM
2VHF        PM           50,045  07/12/93  20:32  2VHF.PM
2VJF        PM           50,045  07/12/93  23:11  2VJF.PM
2VKF        PM           50,045  08/12/93   2:02  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 2.19 Software
(Part Number PSW0004/2.19) for the Mastersizer Micro

Disk 1

Volume in drive A has no label

Directory of A:\

INSTALL	EXE	26,688	28/09/95	16:04	INSTALL.EXE
SIZER	INF	4,852	24/03/99	15:21	SIZER.INF
SIZER	<DIR>		24/03/99	15:51	sizer
	2 file(s)		31,540 bytes		

Directory of A:\sizer

.	<DIR>		24/03/99	15:51	.
..	<DIR>		24/03/99	15:51	..
SIZER	EXE	651,808	23/03/99	17:04	sizer.exe
SMG	EXE	109,008	06/11/95	5:12	smg.exe
PRESENT	DLL	48,000	06/11/95	4:52	present.dll
MSYAN	DLL	47,728	07/11/97	8:49	msyan.dll
SMG	HLP	45,474	19/12/94	4:11	smg.hlp
SIZER	INI	8,001	24/03/99	8:56	SIZER.INI
PAINT	EXE	86,528	11/03/92	8:50	PAINT.EXE
USR_RI	INI	532	26/04/95	10:50	USR_RI.INI
STD_RI	INI	413	08/11/95	11:09	STD_RI.INI
DIMSZS	DD	1,040	14/07/93	12:37	DIMSZS.DD
R0100RF	SD	190	11/11/93	15:47	R0100RF.SD
LOG100	SIZ	1,420	25/04/95	12:28	LOG100.SIZ
CTL3D	DLL	14,448	08/02/93	16:04	CTL3D.DLL
DEFAULT	SAM	4,608	19/11/93	11:10	DEFAULT.SAM
STANDARD	SHA	1,559	30/11/93	14:23	STANDARD.SHA
DDELINK	BSC	268	28/10/92	14:19	DDELINK.BSC
LOGFILEH	BSC	947	28/10/92	14:25	LOGFILEH.BSC
LOGFILEL	BSC	772	28/10/92	14:30	LOGFILEL.BSC
LOGREMH	BSC	651	28/10/92	14:32	LOGREMH.BSC
LOGREML	BSC	652	28/10/92	14:35	LOGREML.BSC
SETUPALL	BSC	35	09/12/91	6:54	SETUPALL.BSC
MTROMP	BSC	2,104	11/11/94	11:13	MTROMP.BSC
8255	BSC	1,182	28/10/92	13:54	8255.BSC
PROMSAM	BSC	62	19/02/97	17:27	PROMSAM.BSC
SHAPEM	BSC	19,284	17/02/97	14:54	SHAPEM.BSC
1	BMP	502	20/01/92	21:25	1.BMP
2	BMP	502	20/01/92	21:29	2.BMP
3	BMP	502	20/01/92	21:31	3.BMP
4	BMP	502	20/01/92	21:35	4.BMP
5	BMP	502	20/01/92	21:36	5.BMP
6	BMP	502	20/01/92	21:37	6.BMP
7	BMP	502	20/01/92	21:38	7.BMP
8	BMP	502	20/01/92	21:38	8.BMP
ALIGN	BMP	502	02/11/91	17:14	ALIGN.BMP
BACKGND	BMP	502	07/01/92	23:04	BACKGND.BMP
CALCULA2	BMP	502	19/11/91	23:00	CALCULA2.BMP
CALCULAT	BMP	502	02/11/91	17:23	CALCULAT.BMP

Appendix C to QSpec Software Update Notification for version 2.19 software

CHECK	BMP	502	17/01/92	22:54	CHECK.BMP
CLEARGRF	BMP	502	02/11/91	17:01	CLEARGRF.BMP
DISKLOAD	BMP	502	12/11/91	23:00	DISKLOAD.BMP
DISKSAVE	BMP	502	12/11/91	23:00	DISKSAVE.BMP
DOCUMENT	BMP	502	03/11/91	11:27	DOCUMENT.BMP
EXIT	BMP	502	07/04/93	14:48	EXIT.BMP
GO	BMP	502	03/11/91	11:46	GO.BMP
GRFDOWN	BMP	502	02/11/91	16:56	GRFDOWN.BMP
GRFLEFT	BMP	502	02/11/91	16:39	GRFLEFT.BMP
GRFRIGHT	BMP	502	02/11/91	16:37	GRFRIGHT.BMP
GRFUP	BMP	502	02/11/91	16:55	GRFUP.BMP
HARDWARE	BMP	502	08/01/92	8:32	HARDWARE.BMP
HELP	BMP	502	17/01/92	22:37	HELP.BMP
INSPECT	BMP	502	02/11/91	17:42	INSPECT.BMP
LOADREC	BMP	502	03/11/91	12:36	LOADREC.BMP
MIX	BMP	502	03/11/91	12:11	MIX.BMP
OPENFIL2	BMP	502	19/11/91	22:55	OPENFIL2.BMP
OPENFILE	BMP	502	03/11/91	12:46	OPENFILE.BMP
PAGE	BMP	502	07/01/92	23:00	PAGE.BMP
PRESENT	BMP	502	07/01/92	22:54	PRESENT.BMP
PRINT	BMP	502	12/11/91	23:09	PRINT.BMP
PROGRAM	BMP	502	22/11/91	19:53	PROGRAM.BMP
REPORT	BMP	502	07/01/92	22:58	REPORT.BMP
RR	BMP	1,870	17/02/97	14:55	RR.BMP
SAMPLE	BMP	526	20/07/93	11:56	SAMPLE.BMP
SAMPLER	BMP	502	07/01/92	23:18	SAMPLER.BMP
SAVEREC	BMP	502	03/11/91	12:36	SAVEREC.BMP
SEQUENC	BMP	502	03/11/91	11:40	SEQUENC.BMP
SETUP	BMP	502	26/11/91	21:03	SETUP.BMP
SPACE	BMP	502	06/05/93	11:22	SPACE.BMP
STOP	BMP	502	03/08/93	17:25	STOP.BMP
TEMPLATE	BMP	502	02/11/91	14:14	TEMPLATE.BMP
VDATA	BMP	502	17/11/91	11:01	VDATA.BMP
VRESULT	BMP	502	17/11/91	11:03	VRESULT.BMP
WASH	BMP	502	03/11/91	12:06	WASH.BMP
ZMLEFT	BMP	502	02/11/91	16:46	ZMLEFT.BMP
ZMRIGHT	BMP	502	02/11/91	16:47	ZMRIGHT.BMP
MDATA	PAG	6,138	02/01/96	10:18	MDATA.PAG
MDIFF	PAG	6,791	15/12/95	9:43	MDIFF.PAG
MFIT	PAG	7,095	15/12/95	9:41	MFIT.PAG
MPARAM	PAG	10,997	15/12/95	9:43	MPARAM.PAG
MRESHIST	PAG	12,112	15/12/95	9:34	MRESHIST.PAG
MRESMEAS	PAG	11,710	15/12/95	9:30	MRESMEAS.PAG
MRESSIVA	PAG	12,026	15/12/95	9:39	MRESSIVA.PAG
MRESSIVB	PAG	12,035	15/12/95	9:39	MRESSIVB.PAG
MRESSIVI	PAG	11,951	15/12/95	9:44	MRESSIVI.PAG
MRESSTAT	PAG	12,841	15/12/95	9:38	MRESSTAT.PAG
MSHAPE	PAG	3,388	14/12/95	17:30	MSHAPE.PAG
MSTATS	PAG	4,359	15/12/95	9:42	MSTATS.PAG
MTROMP	PAG	11,039	15/12/95	9:47	MTROMP.PAG
COLOURS	PAG	510	21/07/91	22:17	COLOURS.PAG
MDATA	REP	6,815	02/01/96	10:19	MDATA.REP
MDIFF	REP	8,142	07/08/95	13:28	MDIFF.REP
MFIT	REP	8,352	09/08/95	16:20	MFIT.REP
MPARAM	REP	21,440	01/11/95	11:37	MPARAM.REP
MRESHIST	REP	13,362	07/08/95	13:35	MRESHIST.REP
MRESMEAS	REP	12,693	07/08/95	13:39	MRESMEAS.REP
MRESSIVA	REP	13,239	07/08/95	14:10	MRESSIVA.REP

Appendix C to QSpec Software Update Notification for version 2.19 software

MRESSIVB	REP	13,308	01/11/95	10:53	MRESSIVB.REP
MRESSIVI	REP	13,249	07/08/95	14:13	MRESSIVI.REP
MRESSTAT	REP	14,148	07/08/95	14:19	MRESSTAT.REP
MSHAPE	REP	5,499	14/12/95	17:22	MSHAPE.REP
MSTATS	REP	4,675	13/02/95	14:59	MSTATS.REP
MTROMP	REP	14,183	07/08/95	14:22	MTROMP.REP
COLOURS	REP	510	21/07/91	22:17	COLOURS.REP
HEADER	WMF	11,126	05/07/94	13:21	HEADER.WMF
SIZER	WMF	7,766	04/08/91	19:39	SIZER.WMF
IHSD204	IH	24,078	07/12/95	16:17	IHSD204.IH
MICRO105	IH	5,228	23/05/95	9:17	MICRO105.IH
		106 file(s)	1,393,509 bytes		

Total files listed:

108 file(s)	1,425,049 bytes
3 dir(s)	12,800 bytes free

Disk 2

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>	24/03/99	15:56	sizer
		0 file(s)	0 bytes	

Directory of A:\sizer

.	<DIR>	24/03/99	15:56	.
..	<DIR>	24/03/99	15:56	..
SIZER	HLP	1,228,717	23/03/99	16:25 sizer.hlp
		1 file(s)	1,228,717 bytes	

Total files listed:

1 file(s)	1,228,717 bytes
3 dir(s)	228,352 bytes free

Disk 3

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>	24/03/99	15:57	sizer
		0 file(s)	0 bytes	

Directory of A:\sizer

.	<DIR>	24/03/99	15:57	.
..	<DIR>	24/03/99	15:57	..
4\$\$D	PM	37,245	15/11/93	11:18 4\$\$D.PM
4BHF	PM	37,245	24/04/95	1:12 4BHF.PM
4CDF	PM	37,245	24/04/95	3:13 4CDF.PM
4CFF	PM	37,245	05/04/95	12:56 4CFF.PM
4CHE	PM	37,245	05/04/95	14:12 4CHE.PM
4CHF	PM	37,245	05/04/95	15:33 4CHF.PM
4GAD	PM	37,245	05/04/95	16:47 4GAD.PM

Appendix C to QSpec Software Update Notification for version 2.19 software

```

4GAE      PM      37,245  05/04/95  18:05 4GAE.PM
4GDD      PM      37,245  06/04/95   8:57 4GDD.PM
4GFD      PM      37,245  06/04/95  10:10 4GFD.PM
4GHD      PM      37,245  06/04/95  11:23 4GHD.PM
4IAD      PM      37,245  06/04/95  12:36 4IAD.PM
4IAE      PM      37,245  06/04/95  13:58 4IAE.PM
4IDD      PM      37,245  06/04/95  15:19 4IDD.PM
4IDF      PM      37,245  06/04/95  17:13 4IDF.PM
4IFD      PM      37,245  06/04/95  18:26 4IFD.PM
4IFF      PM      37,245  07/04/95   8:44 4IFF.PM
4IHD      PM      37,245  07/04/95   9:56 4IHD.PM
4IHE      PM      37,245  07/04/95  11:11 4IHE.PM
4IHF      PM      37,245  07/04/95  12:30 4IHF.PM
4JAD      PM      37,245  12/11/93   2:38 4JAD.PM
4JDD      PM      37,245  07/04/95  13:43 4JDD.PM
4JFD      PM      37,245  07/04/95  14:55 4JFD.PM
4JHD      PM      37,245  07/04/95  16:38 4JHD.PM
4KAF      PM      37,245  10/04/95   8:05 4KAF.PM
4KHF      PM      37,245  10/04/95   9:26 4KHF.PM
4LAD      PM      37,245  12/11/93   0:06 4LAD.PM
4LAF      PM      37,245  10/04/95  10:47 4LAF.PM
4LDD      PM      37,245  10/04/95  12:00 4LDD.PM
4LDF      PM      37,245  10/04/95  13:20 4LDF.PM
4LFD      PM      37,245  24/04/95  12:31 4LFD.PM
4LFF      PM      37,245  10/04/95  15:19 4LFF.PM
4LHD      PM      37,245  11/04/95   8:12 4LHD.PM
4LHF      PM      37,245  11/04/95   9:31 4LHF.PM
4MAC      PM      37,245  11/04/95  11:53 4MAC.PM
4MAD      PM      37,245  11/04/95  13:06 4MAD.PM
4MAE      PM      37,245  11/04/95  14:22 4MAE.PM
4MAF      PM      37,245  11/04/95  15:43 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

Volume in drive A has no label

Directory of A:\

```

SIZER      <DIR>      24/03/99  16:01 sizer
0 file(s)      0 bytes
    
```

Directory of A:\sizer

```

.          <DIR>      24/03/99  16:01 .
..         <DIR>      24/03/99  16:01 ..
4MDD      PM      37,245  12/04/95   8:01 4MDD.PM
4MDF      PM      37,245  24/04/95   8:19 4MDF.PM
4MFC      PM      37,245  24/04/95   8:19 4MFC.PM
4MFD      PM      37,245  24/04/95   8:19 4MFD.PM
4MFE      PM      37,245  24/04/95   8:19 4MFE.PM
4MFF      PM      37,245  24/04/95   8:19 4MFF.PM
4MHC      PM      37,245  24/04/95   8:19 4MHC.PM
4MHD      PM      37,245  24/04/95   8:19 4MHD.PM
    
```

Appendix C to QSpec Software Update Notification for version 2.19 software

4MHE	PM	37,245	24/04/95	8:19	4MHE.PM
4MHF	PM	37,245	24/04/95	8:19	4MHF.PM
4NAD	PM	37,245	12/11/93	1:22	4NAD.PM
4NAE	PM	37,245	13/04/95	12:10	4NAE.PM
4NAF	PM	37,245	13/04/95	13:31	4NAF.PM
4NBD	PM	37,245	13/04/95	14:43	4NBD.PM
4NDD	PM	37,245	13/04/95	15:56	4NDD.PM
4NDE	PM	37,245	13/04/95	17:12	4NDE.PM
4NDF	PM	37,245	13/04/95	18:32	4NDF.PM
4NED	PM	37,245	18/04/95	6:30	4NED.PM
4NFD	PM	37,245	18/04/95	7:43	4NFD.PM
4NFE	PM	37,245	18/04/95	9:13	4NFE.PM
4NFF	PM	37,245	18/04/95	10:36	4NFF.PM
4NGD	PM	37,245	18/04/95	11:49	4NGD.PM
4NGE	PM	37,245	18/04/95	13:05	4NGE.PM
4NHD	PM	37,245	18/04/95	14:20	4NHD.PM
4NHE	PM	37,245	18/04/95	16:41	4NHE.PM
4NHF	PM	37,245	18/04/95	18:01	4NHF.PM
4NJD	PM	37,245	19/04/95	7:15	4NJD.PM
4NJE	PM	37,245	19/04/95	10:16	4NJE.PM
4NJF	PM	37,245	19/04/95	13:32	4NJF.PM
4OAD	PM	37,245	11/11/93	20:10	4OAD.PM
4OAE	PM	37,245	19/04/95	16:27	4OAE.PM
4ODD	PM	37,245	19/04/95	17:39	4ODD.PM
4OED	PM	37,245	21/10/94	13:23	4OED.PM
4OFD	PM	37,245	05/09/94	18:50	4OFD.PM
4OGD	PM	37,245	20/04/95	6:38	4OGD.PM
4OHD	PM	37,245	11/11/93	21:25	4OHD.PM
4OHE	PM	37,245	20/04/95	7:53	4OHE.PM
4OID	PM	37,245	20/04/95	10:04	4OID.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

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>	24/03/99	16:03	sizer
		0 file(s)	0 bytes	

Directory of A:\sizer

.	<DIR>	24/03/99	16:03	.
..	<DIR>	24/03/99	16:03	..
4OJD	PM	37,245	20/04/95	12:39 4OJD.PM
4OJE	PM	37,245	20/04/95	15:48 4OJE.PM
4OKD	PM	37,245	20/04/95	17:17 4OKD.PM
4PAD	PM	37,245	11/11/93	18:54 4PAD.PM
4PAE	PM	37,245	21/04/95	8:05 4PAE.PM
4PDD	PM	37,245	21/04/95	12:11 4PDD.PM
4PDE	PM	37,245	21/04/95	14:27 4PDE.PM
4PED	PM	37,245	21/04/95	15:42 4PED.PM
4PFD	PM	37,245	05/12/94	16:46 4PFD.PM

Appendix C to QSpec Software Update Notification for version 2.19 software

4PFE	PM	37,245	21/04/95	20:05	4PFE.PM
4PFF	PM	37,245	21/04/95	22:08	4PFF.PM
4PHD	PM	37,245	19/05/94	10:52	4PHD.PM
4PHE	PM	37,245	22/04/95	0:03	4PHE.PM
4PHF	PM	37,245	22/04/95	2:05	4PHF.PM
4PID	PM	37,245	22/04/95	3:56	4PID.PM
4PJD	PM	37,245	22/04/95	5:51	4PJD.PM
4PJE	PM	37,245	22/04/95	7:49	4PJE.PM
4PKD	PM	37,245	22/04/95	10:05	4PKD.PM
4QAD	PM	37,245	22/04/95	11:58	4QAD.PM
4QAE	PM	37,245	22/04/95	13:57	4QAE.PM
4QDD	PM	37,245	22/04/95	15:49	4QDD.PM
4QFD	PM	37,245	22/04/95	17:41	4QFD.PM
4QHD	PM	37,245	22/04/95	19:34	4QHD.PM
4QHE	PM	37,245	22/04/95	21:31	4QHE.PM
4QJD	PM	37,245	22/04/95	23:27	4QJD.PM
4RAD	PM	37,245	23/04/95	1:23	4RAD.PM
4RAE	PM	37,245	23/04/95	3:24	4RAE.PM
4RFD	PM	37,245	23/04/95	5:19	4RFD.PM
4RHD	PM	37,245	10/01/94	15:50	4RHD.PM
4RKD	PM	37,245	23/04/95	7:36	4RKD.PM
4SHD	PM	37,245	23/04/95	9:35	4SHD.PM
4SKD	PM	37,245	23/04/95	11:54	4SKD.PM
4THD	PM	37,245	23/04/95	13:58	4THD.PM
4TJD	PM	37,245	11/11/93	22:51	4TJD.PM
4TKD	PM	37,245	23/04/95	16:22	4TKD.PM
4UHD	PM	37,245	23/04/95	18:33	4UHD.PM
4UKD	PM	37,245	23/04/95	21:01	4UKD.PM
4_METAL	PM	37,245	23/04/95	23:11	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 2.19 Software
(Part Number PSW0005/2.19) for the Mastersizer Microplus

Disk 1

Volume in drive A has no label

Directory of A:\

```
INSTALL  EXE           26,688  28/09/95  16:04  INSTALL.EXE
SIZER    INF           4,884  24/03/99  15:15  SIZER.INF
SIZER    <DIR>          24/03/99  15:39  sizer
          2 file(s)          31,572 bytes
```

Directory of A:\sizer

```
.          <DIR>          24/03/99  15:39  .
..         <DIR>          24/03/99  15:39  ..
SIZER     EXE       651,808  23/03/99  17:04  sizer.exe
SMG       EXE       109,008  06/11/95   5:12  smg.exe
PRESENT   DLL        48,000  06/11/95   4:52  present.dll
MSYAN     DLL        53,920  26/01/98  11:44  msyan.dll
SMG       HLP        45,474  19/12/94   4:11  smg.hlp
SIZER     INI         7,982  24/03/99   9:02  SIZER.INI
PAINT     EXE       86,528  11/03/92   8:50  PAINT.EXE
USR_RI    INI         532    26/04/95  10:50  USR_RI.INI
STD_RI    INI         413    08/11/95  11:10  STD_RI.INI
DIMSZS    DD          1,040  14/07/93  12:37  DIMSZS.DD
R0300RFP  SD           254    26/10/94  10:43  R0300RFP.SD
LOG100    SIZ          1,420  07/03/95  16:50  LOG100.SIZ
CTL3D     DLL       14,448  08/02/93  16:04  CTL3D.DLL
DEFAULT   SAM          1,536  11/11/97   9:57  DEFAULT.SAM
STANDARD  SHA          2,104  13/02/95  15:08  STANDARD.SHA
DDELINK   BSC           268    28/10/92  14:19  DDELINK.BSC
LOGFILEH  BSC           947    28/10/92  14:25  LOGFILEH.BSC
LOGFILEL  BSC           772    28/10/92  14:30  LOGFILEL.BSC
LOGREMH   BSC           651    28/10/92  14:32  LOGREMH.BSC
LOGREML   BSC           652    28/10/92  14:35  LOGREML.BSC
SETUPALL  BSC            35    09/12/91   6:54  SETUPALL.BSC
MTROMP    BSC          2,104  11/11/94  11:13  MTROMP.BSC
8255      BSC          1,182  28/10/92  13:54  8255.BSC
PROMSAM   BSC            63    23/05/95  10:15  PROMSAM.BSC
1         BMP            502    20/01/92  21:25  1.BMP
2         BMP            502    20/01/92  21:29  2.BMP
3         BMP            502    20/01/92  21:31  3.BMP
4         BMP            502    20/01/92  21:35  4.BMP
5         BMP            502    20/01/92  21:36  5.BMP
6         BMP            502    20/01/92  21:37  6.BMP
7         BMP            502    20/01/92  21:38  7.BMP
8         BMP            502    20/01/92  21:38  8.BMP
ALIGN     BMP            502    02/11/91  17:14  ALIGN.BMP
BACKGND   BMP            502    07/01/92  23:04  BACKGND.BMP
CALCULA2  BMP            502    19/11/91  23:00  CALCULA2.BMP
CALCULAT  BMP            502    02/11/91  17:23  CALCULAT.BMP
CHECK     BMP            502    17/01/92  22:54  CHECK.BMP
```

Appendix D to QSpec Software Update Notification for version 2.19 software

CLEARGRF	BMP	502	02/11/91	17:01	CLEARGRF.BMP
DISKLOAD	BMP	502	12/11/91	23:00	DISKLOAD.BMP
DISKSAVE	BMP	502	12/11/91	23:00	DISKSAVE.BMP
DOCUMENT	BMP	502	03/11/91	11:27	DOCUMENT.BMP
EXIT	BMP	502	07/04/93	14:48	EXIT.BMP
GO	BMP	502	03/11/91	11:46	GO.BMP
GRFDOWN	BMP	502	02/11/91	16:56	GRFDOWN.BMP
GRFLEFT	BMP	502	02/11/91	16:39	GRFLEFT.BMP
GRFRIGHT	BMP	502	02/11/91	16:37	GRFRIGHT.BMP
GRFUP	BMP	502	02/11/91	16:55	GRFUP.BMP
HARDWARE	BMP	502	08/01/92	8:32	HARDWARE.BMP
HELP	BMP	502	17/01/92	22:37	HELP.BMP
INSPECT	BMP	502	02/11/91	17:42	INSPECT.BMP
LOADREC	BMP	502	03/11/91	12:36	LOADREC.BMP
MIX	BMP	502	03/11/91	12:11	MIX.BMP
OPENFIL2	BMP	502	19/11/91	22:55	OPENFIL2.BMP
OPENFILE	BMP	502	03/11/91	12:46	OPENFILE.BMP
PAGE	BMP	502	07/01/92	23:00	PAGE.BMP
PRESENT	BMP	502	07/01/92	22:54	PRESENT.BMP
PRINT	BMP	502	12/11/91	23:09	PRINT.BMP
PROGRAM	BMP	502	22/11/91	19:53	PROGRAM.BMP
REPORT	BMP	502	07/01/92	22:58	REPORT.BMP
RR	BMP	1,870	17/02/97	14:55	RR.BMP
SAMPLE	BMP	526	20/07/93	11:56	SAMPLE.BMP
SAMPLER	BMP	502	07/01/92	23:18	SAMPLER.BMP
SAVEREC	BMP	502	03/11/91	12:36	SAVEREC.BMP
SEQUENC	BMP	502	03/11/91	11:40	SEQUENC.BMP
SETUP	BMP	502	26/11/91	21:03	SETUP.BMP
SPACE	BMP	502	06/05/93	11:22	SPACE.BMP
STOP	BMP	502	03/08/93	17:25	STOP.BMP
TEMPLATE	BMP	502	02/11/91	14:14	TEMPLATE.BMP
VDATA	BMP	502	17/11/91	11:01	VDATA.BMP
VRESULT	BMP	502	17/11/91	11:03	VRESULT.BMP
WASH	BMP	502	03/11/91	12:06	WASH.BMP
ZMLEFT	BMP	502	02/11/91	16:46	ZMLEFT.BMP
ZMRIGHT	BMP	502	02/11/91	16:47	ZMRIGHT.BMP
MDATA	PAG	6,138	02/01/96	10:18	MDATA.PAG
MDIFF	PAG	6,791	15/12/95	9:43	MDIFF.PAG
MFIT	PAG	7,095	15/12/95	9:41	MFIT.PAG
MPARAM	PAG	10,997	15/12/95	9:43	MPARAM.PAG
MRESHIST	PAG	12,112	15/12/95	9:34	MRESHIST.PAG
MRESMEAS	PAG	11,710	15/12/95	9:30	MRESMEAS.PAG
MRESSIVA	PAG	12,026	15/12/95	9:39	MRESSIVA.PAG
MRESSIVB	PAG	12,035	15/12/95	9:39	MRESSIVB.PAG
MRESSIVI	PAG	11,951	15/12/95	9:44	MRESSIVI.PAG
MRESSTAT	PAG	12,841	15/12/95	9:38	MRESSTAT.PAG
MSHAPE	PAG	3,388	14/12/95	17:30	MSHAPE.PAG
MSTATS	PAG	4,359	15/12/95	9:42	MSTATS.PAG
MTROMP	PAG	11,039	15/12/95	9:47	MTROMP.PAG
COLOURS	PAG	510	21/07/91	22:17	COLOURS.PAG
MDATA	REP	6,815	02/01/96	10:19	MDATA.REP
MDIFF	REP	8,142	07/08/95	13:28	MDIFF.REP
MFIT	REP	8,352	09/08/95	16:20	MFIT.REP
MPARAM	REP	21,440	01/11/95	11:37	MPARAM.REP
MRESHIST	REP	13,362	07/08/95	13:35	MRESHIST.REP
MRESMEAS	REP	12,693	07/08/95	13:39	MRESMEAS.REP
MRESSIVA	REP	13,239	07/08/95	14:10	MRESSIVA.REP
MRESSIVB	REP	13,308	01/11/95	10:53	MRESSIVB.REP

Appendix D to QSpec Software Update Notification for version 2.19 software

MRESSIVI	REP	13,249	07/08/95	14:13	MRESSIVI.REP
MRESSTAT	REP	14,148	07/08/95	14:19	MRESSTAT.REP
MSHAPE	REP	5,499	14/12/95	17:22	MSHAPE.REP
MSTATS	REP	4,675	13/02/95	14:59	MSTATS.REP
MTROMP	REP	14,183	07/08/95	14:22	MTROMP.REP
COLOURS	REP	510	21/07/91	22:17	COLOURS.REP
HEADER	WMF	11,126	05/07/94	13:21	HEADER.WMF
SIZER	WMF	7,766	04/08/91	19:39	SIZER.WMF
IHSD204	IH	24,078	07/12/95	16:17	IHSD204.IH
MICRO105	IH	5,228	23/05/95	9:17	MICRO105.IH
		105 file(s)			1,377,936 bytes

Total files listed:

107 file(s)	1,409,508 bytes
3 dir(s)	29,184 bytes free

Disk 2

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>		24/03/99	15:43	sizer
		0 file(s)			0 bytes

Directory of A:\sizer

.	<DIR>		24/03/99	15:43	.
..	<DIR>		24/03/99	15:43	..
SIZER	HLP	1,228,717	23/03/99	16:25	sizer.hlp
SHAPEMP	BSC	19,321	17/02/97	14:54	SHAPEMP.BSC
		2 file(s)			1,248,038 bytes

Total files listed:

2 file(s)	1,248,038 bytes
3 dir(s)	208,896 bytes free

Disk 3

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>		24/03/99	15:44	sizer
		0 file(s)			0 bytes

Directory of A:\sizer

.	<DIR>		24/03/99	15:44	.
..	<DIR>		24/03/99	15:44	..
5\$\$D	PM	37,245	26/10/94	14:26	5\$\$D.PM
5BHF	PM	37,245	02/03/95	20:23	5BHF.PM
5CDF	PM	37,245	02/03/95	21:23	5CDF.PM
5CFF	PM	37,245	02/03/95	22:23	5CFF.PM
5CHE	PM	37,245	02/03/95	23:19	5CHE.PM
5CHF	PM	37,245	03/03/95	0:19	5CHF.PM
5GAD	PM	37,245	28/02/95	15:46	5GAD.PM

Appendix D to QSpec Software Update Notification for version 2.19 software

```

5GAE      PM      37,245  03/03/95   3:21  5GAE.PM
5GDD      PM      37,245  28/02/95  17:31  5GDD.PM
5GFD      PM      37,245  28/02/95  18:23  5GFD.PM
5GHD      PM      37,245  28/02/95  19:16  5GHD.PM
5IAD      PM      37,245  28/02/95  13:08  5IAD.PM
5IAE      PM      37,245  03/03/95   4:19  5IAE.PM
5IDD      PM      37,245  03/03/95   5:14  5IDD.PM
5IDF      PM      37,245  03/03/95   6:14  5IDF.PM
5IFD      PM      37,245  28/02/95  14:01  5IFD.PM
5IFF      PM      37,245  03/03/95   7:13  5IFF.PM
5IHD      PM      37,245  03/03/95   8:08  5IHD.PM
5IHE      PM      37,245  03/03/95   9:05  5IHE.PM
5IHF      PM      37,245  03/03/95  10:04  5IHF.PM
5JAD      PM      37,245  27/02/95  16:01  5JAD.PM
5JDD      PM      37,245  03/03/95  12:57  5JDD.PM
5JFD      PM      37,245  28/02/95  14:53  5JFD.PM
5JHD      PM      37,245  27/02/95  16:53  5JHD.PM
5KAF      PM      37,245  27/02/95  17:51  5KAF.PM
5KHF      PM      37,245  27/02/95  18:48  5KHF.PM
5LAD      PM      37,245  27/02/95  19:41  5LAD.PM
5LAF      PM      37,245  03/03/95  13:56  5LAF.PM
5LDD      PM      37,245  03/03/95  14:49  5LDD.PM
5LDF      PM      37,245  03/03/95  15:47  5LDF.PM
5LFD      PM      37,245  03/03/95  16:40  5LFD.PM
5LFF      PM      37,245  03/03/95  17:38  5LFF.PM
5LHD      PM      37,245  03/03/95  18:31  5LHD.PM
5LHF      PM      37,245  28/02/95  12:15  5LHF.PM
5MAC      PM      37,245  03/03/95  20:27  5MAC.PM
5MAD      PM      37,245  27/02/95  20:35  5MAD.PM
5MAE      PM      37,245  03/03/95  21:23  5MAE.PM
5MAF      PM      37,245  03/03/95  22:22  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

Volume in drive A has no label

Directory of A:\

```

SIZER      <DIR>      24/03/99  15:47  sizer
0 file(s)      0 bytes
    
```

Directory of A:\sizer

```

.          <DIR>      24/03/99  15:47  .
..         <DIR>      24/03/99  15:47  ..
5MDD      PM      37,245  03/03/95  23:15  5MDD.PM
5MDF      PM      37,245  04/03/95   0:12  5MDF.PM
5MFC      PM      37,245  04/03/95   1:04  5MFC.PM
5MFD      PM      37,245  04/03/95   1:57  5MFD.PM
5MFE      PM      37,245  04/03/95   2:52  5MFE.PM
5MFF      PM      37,245  04/03/95   3:50  5MFF.PM
5MHC      PM      37,245  04/03/95   4:42  5MHC.PM
5MHD      PM      37,245  04/03/95   5:34  5MHD.PM
    
```

Appendix D to QSpec Software Update Notification for version 2.19 software

5MHE	PM	37,245	04/03/95	6:29	5MHE.PM
5MHF	PM	37,245	04/03/95	7:27	5MHF.PM
5NAD	PM	37,245	27/10/94	2:11	5NAD.PM
5NAE	PM	37,245	27/02/95	21:30	5NAE.PM
5NAF	PM	37,245	04/03/95	9:21	5NAF.PM
5NBD	PM	37,245	04/03/95	10:14	5NBD.PM
5NDD	PM	37,245	27/02/95	22:22	5NDD.PM
5NDE	PM	37,245	04/03/95	11:09	5NDE.PM
5NDF	PM	37,245	04/03/95	12:07	5NDF.PM
5NED	PM	37,245	04/03/95	13:00	5NED.PM
5NFD	PM	37,245	04/03/95	13:53	5NFD.PM
5NFE	PM	37,245	04/03/95	14:48	5NFE.PM
5NFF	PM	37,245	04/03/95	15:45	5NFF.PM
5NGD	PM	37,245	04/03/95	16:38	5NGD.PM
5NGE	PM	37,245	04/03/95	17:33	5NGE.PM
5NHD	PM	37,245	20/01/95	14:00	5NHD.PM
5NHE	PM	37,245	27/02/95	23:16	5NHE.PM
5NHF	PM	37,245	04/03/95	18:31	5NHF.PM
5NJD	PM	37,245	04/03/95	19:25	5NJD.PM
5NJE	PM	37,245	28/02/95	0:11	5NJE.PM
5NJF	PM	37,245	04/03/95	20:23	5NJF.PM
5OAD	PM	37,245	26/10/94	20:29	5OAD.PM
5OAE	PM	37,245	04/03/95	22:27	5OAE.PM
5ODD	PM	37,245	28/02/95	1:03	5ODD.PM
5OED	PM	37,245	04/03/95	23:20	5OED.PM
5OFD	PM	37,245	27/10/94	0:48	5OFD.PM
5OGD	PM	37,245	05/03/95	0:13	5OGD.PM
5OHD	PM	37,245	26/10/94	19:05	5OHD.PM
5OHE	PM	37,245	28/02/95	1:57	5OHE.PM
5OID	PM	37,245	05/03/95	1:06	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

Volume in drive A has no label

Directory of A:\

SIZER	<DIR>	24/03/99	15:49	sizer
0 file(s)		0 bytes		

Directory of A:\sizer

.	<DIR>	24/03/99	15:49	.
..	<DIR>	24/03/99	15:49	..
5OJD	PM	37,245	28/02/95	2:51 5OJD.PM
5OJE	PM	37,245	28/02/95	3:47 5OJE.PM
5OKD	PM	37,245	05/03/95	2:10 5OKD.PM
5PAD	PM	37,245	26/10/94	21:53 5PAD.PM
5PAE	PM	37,245	05/03/95	3:06 5PAE.PM
5PDD	PM	37,245	28/02/95	4:39 5PDD.PM
5PDE	PM	37,245	05/03/95	4:02 5PDE.PM
5PED	PM	37,245	05/03/95	4:55 5PED.PM
5PFD	PM	37,245	28/02/95	5:32 5PFD.PM

Appendix D to QSpec Software Update Notification for version 2.19 software

5PFE	PM	37,245	05/03/95	5:50	5PFE.PM
5PFF	PM	37,245	05/03/95	6:49	5PFF.PM
5PHD	PM	37,245	28/02/95	6:24	5PHD.PM
5PHE	PM	37,245	05/03/95	7:44	5PHE.PM
5PHF	PM	37,245	05/03/95	8:42	5PHF.PM
5PID	PM	37,245	05/03/95	9:35	5PID.PM
5PJD	PM	37,245	28/02/95	20:10	5PJD.PM
5PJE	PM	37,245	05/03/95	10:32	5PJE.PM
5PKD	PM	37,245	05/03/95	11:36	5PKD.PM
5QAD	PM	37,245	05/03/95	12:31	5QAD.PM
5QAE	PM	37,245	05/03/95	13:28	5QAE.PM
5QDD	PM	37,245	05/03/95	14:22	5QDD.PM
5QFD	PM	37,245	05/03/95	15:16	5QFD.PM
5QHD	PM	37,245	28/02/95	7:17	5QHD.PM
5QHE	PM	37,245	05/03/95	17:08	5QHE.PM
5QJD	PM	37,245	05/03/95	19:05	5QJD.PM
5RAD	PM	37,245	05/03/95	20:01	5RAD.PM
5RAE	PM	37,245	05/03/95	20:59	5RAE.PM
5RFD	PM	37,245	05/03/95	22:51	5RFD.PM
5RHD	PM	37,245	28/02/95	8:11	5RHD.PM
5RKD	PM	37,245	06/03/95	5:41	5RKD.PM
5SHD	PM	37,245	06/03/95	10:44	5SHD.PM
5SKD	PM	37,245	06/03/95	16:52	5SKD.PM
5THD	PM	37,245	28/02/95	9:10	5THD.PM
5TJD	PM	37,245	26/10/94	23:26	5TJD.PM
5TKD	PM	37,245	28/02/95	10:17	5TKD.PM
5UHD	PM	37,245	07/03/95	1:15	5UHD.PM
5UKD	PM	37,245	28/02/95	21:19	5UKD.PM
5_METAL	PM	37,245	28/02/95	11:18	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