
Software Update Notification



Filename: QSpec Update Notification for v2.17 software.doc Doct. Version No: 1
Date: 17th December 1997 Author: R . M. Jones

Mastersizer (S, X, Micro, Microplus) version 2.17 Software
for use with (English) Microsoft Windows version 3.1,
Windows for Workgroups version 3.11 and Windows95

Mastersizer S

(Part Number PSW 0002/2.17)

Mastersizer X

(Part Number PSW 0003/2.17)

Mastersizer Micro

(Part Number PSW 0004/2.17)

Mastersizer Microplus

(Part Number PSW 0005/2.17)

Contents

Mastersizer S	1
(Part Number PSW 0002/2.17)	1
Mastersizer X	1
(Part No .PSW 0003/2.17)	1
Mastersizer Micro	1
(Part Number PSW 0004/2.17)	1
Mastersizer Microplus	1
(Part Number PSW 0005/2.17)	1
Contents	2
1. Introduction	2
2. Instrument compatibility	2
3. Functional Features	2
4. Installing Software	4
5. Directory Structure	4

1. Introduction

- 1.1** This note describes the version 2.17 release of Mastersizer software for Windows and describes the software bugs addressed by this release.

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

3.3.1 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.

Version 2.16 was only issued under concession and therefore, for completeness, all of the bug fixes in that release have been included in this list.

Malvern Software Change Request Number	Version in which Bug was spotted	Description
11	2.16	The DDE Request appears not to work; when it did work in V2.15
18	2.15	If system runs with software version 2.15 under remote control, the time-out on obscuration triggering doesn't work.
46	2.16	1) Align display causes crash in Japanese version and the icon disappears in "help about" dialog the icon in the "help about dialog" disappears sometimes.
48	2.15	When the date is set to 2000 or above the software shows the date as "100" for 2000 e.g. 10/2/100 for 10th February 2000. This only occurs for the short version of the date. The long version does not show any problems.
49	2.15	DDE poke problem
52	1.2b	In Malvern Basic, when the command "experiment delay type" is written or "experiment pulselength n,type" is used: The available options can be "first" or "all" pulses but whatever is typed, "first" is always shown no matter what has been typed.
53	2.00 Beta	Memory is not recovered if the sizer software crashes
55	1.26	Values in the data and fit tables sometimes show a small difference.
56	2.15	Results are sometimes different if the calculation is performed with the sizer window minimised.
58	2.15	Memory leak when using "print string" command
59	2.16a1	Crash after printing
60	2.15	There is a problem with the shutdown of the Sizer with Windows 3.1 This manifests itself as follows: 1. Generate a new result by checking measure – calculate 2. Minimise sizer program 3. Edit Windows by closing Program Manager 4. When message from sizer appears asking to save configuration - click "yes" Sizer icon flashes but does not respond to double click. Other parts of Windows do not respond and Windows has to be restarted.
61	2.15	The μ sign needs to be replaced by the word "micro".
62	2.15	The test for more than 50 variables in the Basic language is incorrect.
63	2.15	The Rosin-Rammler distribution generated by the Edit-Generate Result command should produce linear undersize curves when plotted on the Rosin-Rammler axes. If the distribution has a component below 0.1 micron the curve is no longer linear.
65	2.15	The diameter alarm can only be set between the default 5 and 20microns
85	2.16	When using a file with a large number of records, reopening with the basic editor visible causes a stack overflow.

3.5 Distribution Media

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

For Mastersizer S (Part Number PSW 0002/2.17)	Appendix A
For Mastersizer X (Part Number PSW 0003/2.17)	Appendix B
For Mastersizer Micro (Part Number PSW 0004/2.17)	Appendix C
For Mastersizer Microplus (Part Number PSW 0005/2.17)	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\PRESNT	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\PRESNT	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\PRESNT	Malvern Presentation Matrices
\SIZERMU\PROGRAMS	Malvern and User measurement and analysis procedures

For Mastersizer Microplus

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

Complete listing of disks and files for version 2.17 Software
(Part Number PSW 0002/2.17) for the Mastersizer S**Disk 1**

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,113	11-07-97	12:24p	SIZER.INF
SIZER	<DIR>		11-07-97	4:03p	sizer
		2 file(s)			30,801 bytes

Directory of A:\sizer

.	<DIR>		11-07-97	4:03p	.
..	<DIR>		11-07-97	4:03p	..
SIZER	EXE	651,728	11-19-97	11:45a	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	55,456	11-06-97	5:32p	msyan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	8,358	09-04-97	3:20p	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	11-19-97	12:35p	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
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

Appendix A to QSpec Release Note for version 2.17 software

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
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) 1,403,754 bytes

Total files listed:

113 file(s) 1,434,555 bytes

Disk 2

```
Directory of A:\
SIZER          <DIR>          11-07-97  4:13p  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-07-97  4:13p  .
..            <DIR>          11-07-97  4:13p  ..
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
```

Disk 3

```
Directory of A:\
SIZER          <DIR>          11-07-97  4:19p  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-07-97  4:19p  .
..            <DIR>          11-07-97  4:19p  ..
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
```

Disk 4

```
Directory of A:\
SIZER          <DIR>          11-07-97  4:26p  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-07-97  4:26p  .
..            <DIR>          11-07-97  4:26p  ..
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
```

Appendix A to QSpec Release Note for version 2.17 software

```
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
```

Disk 5

```
Directory of A:\
SIZER      <DIR>      11-07-97  4:31p sizer
          0 file(s)      0 bytes
```

```
Directory of A:\sizer
.          <DIR>      11-07-97  4:31p .
..         <DIR>      11-07-97  4:31p ..
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
```


Complete listing of disks and files for version 2.17 Software
(Part Number PSW 0003/2.17) for the Mastersizer X

Disk 1

Directory of A:\

```

INSTALL  EXE          26,688  09-28-95  4:04p  INSTALL.EXE
SIZER    INF          4,575  11-07-97 12:30p  SIZER.INF
SIZER    <DIR>                11-10-97 12:12p  sizer
                2 file(s)                31,263 bytes
    
```

Directory of A:\sizer

```

.          <DIR>                11-10-97 12:12p  .
..         <DIR>                11-10-97 12:12p  ..
SIZER     EXE          651,728  11-19-97 11:45a  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  09-04-97  3:23p  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
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 B to QSpec Release Note for version 2.17 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
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

Appendix B to QSpec Release Note for version 2.17 software

Total files listed:
115 file(s) 1,417,494 bytes

Disk 2

Directory of A:\
SIZER <DIR> 11-10-97 12:20p sizer
0 file(s) 0 bytes

Directory of A:\sizer
. <DIR> 11-10-97 12:20p .
.. <DIR> 11-10-97 12:20p ..
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

Disk 3

Directory of A:\
SIZER <DIR> 11-10-97 12:25p sizer
0 file(s) 0 bytes

Directory of A:\sizer
. <DIR> 11-10-97 12:25p .
.. <DIR> 11-10-97 12:25p ..
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
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

Disk 4

Directory of A:\
SIZER <DIR> 11-10-97 12:30p sizer
QSpec Software Update 17th December 1997
Notification for version 2.17
software.doc

Appendix B to QSpec Release Note for version 2.17 software

0 file(s) 0 bytes

Directory of A:\sizer

```

.          <DIR>          11-10-97 12:30p .
..         <DIR>          11-10-97 12:30p ..
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
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

```

Disk 5

```

Directory of A:\
SIZER      <DIR>          11-10-97 12:40p sizer
0 file(s) 0 bytes

```

Directory of A:\sizer

```

.          <DIR>          11-10-97 12:40p .
..         <DIR>          11-10-97 12:40p ..
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

```

Appendix B to QSpec Release Note for version 2.17 software

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	

Complete listing of disks and files for version 2.17 Software
(Part Number PSW 0004/2.17) for the Mastersizer Micro**Disk 1**

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,852	11-07-97	12:28p	SIZER.INF
SIZER	<DIR>		11-10-97	12:46p	sizer
	2 file(s)				31,540 bytes

Directory of A:\sizer

.	<DIR>		11-10-97	12:46p	.
..	<DIR>		11-10-97	12:46p	..
SIZER	EXE	651,728	11-19-97	11:45a	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	09-04-97	3:28p	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
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

Appendix C to QSpec Release Note for version 2.17 software

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
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	

Disk 2

Directory of A:\

Appendix C to QSpec Release Note for version 2.17 software

```
SIZER          <DIR>          11-10-97 12:52p sizer
                0 file(s)          0 bytes
```

Directory of A:\sizer

```
.              <DIR>          11-10-97 12:52p .
..            <DIR>          11-10-97 12:52p ..
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
```

Disk 3

Directory of A:\

```
SIZER          <DIR>          08-08-97  9:50a sizer
                0 file(s)          0 bytes
```

Directory of A:\sizer

```
.              <DIR>          08-08-97  9:50a .
..            <DIR>          08-08-97  9:50a ..
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
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
```


Disk 4

```
Directory of A:\
SIZER          <DIR>          11-10-97 12:55p sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-10-97 12:55p .
..            <DIR>          11-10-97 12:55p ..
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
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
4OID         PM           37,245 04-20-95 10:04a 4OID.PM
                38 file(s)          1,415,310 bytes
Total files listed:
                38 file(s)          1,415,310 bytes
```

Disk 5

```
Directory of A:\
SIZER          <DIR>          11-10-97 12:58p sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-10-97 12:58p .
..            <DIR>          11-10-97 12:58p ..
4OJD         PM           37,245 04-20-95 12:39p 4OJD.PM
4OJE         PM           37,245 04-20-95  3:48p 4OJE.PM
4OKD         PM           37,245 04-20-95  5:17p 4OKD.PM
4PAD         PM           37,245 11-11-93  6:54p 4PAD.PM
```

Appendix C to QSpec Release Note for version 2.17 software

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	

Complete listing of disks and files for version 2.17 Software
(Part Number PSW 0005/2.17) for the Mastersizer Microplus**Disk 1**

Directory of A:\

INSTALL	EXE	26,688	09-28-95	4:04p	INSTALL.EXE
SIZER	INF	4,884	11-07-97	12:26p	SIZER.INF
SIZER	<DIR>		11-10-97	2:08p	sizer
	2 file(s)				31,572 bytes

Directory of A:\sizer

.	<DIR>		11-10-97	2:08p	.
..	<DIR>		11-10-97	2:08p	..
SIZER	EXE	651,728	11-19-97	11:45a	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	55,456	11-06-97	5:32p	msyan.dll
SMG	HLP	45,474	12-19-94	4:11a	smg.hlp
SIZER	INI	7,982	09-04-97	3:26p	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	11-19-97	12:39p	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
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

Appendix D to QSpec Release Note for version 2.17 software

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
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)		1,379,392 bytes	
Total files listed:					
		107 file(s)		1,410,964 bytes	

Disk 2

Directory of A:\

Appendix D to QSpec Release Note for version 2.17 software

```
SIZER          <DIR>          11-10-97  2:14p  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          11-10-97  2:14p  .
..            <DIR>          11-10-97  2:14p  ..
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
```

Disk 3

Directory of A:\

```
SIZER          <DIR>          11-10-97  2:15p  sizer
                0 file(s)                0 bytes
```

Directory of A:\sizer

```
.              <DIR>          11-10-97  2:15p  .
..            <DIR>          11-10-97  2:15p  ..
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
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
```

Disk 4

```
Directory of A:\
SIZER          <DIR>          08-08-97  9:16a  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          08-08-97  9:16a  .
..            <DIR>          08-08-97  9:16a  ..
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
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
```

Disk 5

```
Directory of A:\
SIZER          <DIR>          11-10-97  2:47p  sizer
                0 file(s)          0 bytes
```

```
Directory of A:\sizer
.              <DIR>          11-10-97  2:47p  .
..            <DIR>          11-10-97  2:47p  ..
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
```

Appendix D to QSpec Release Note for version 2.17 software

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	