VisionMate SR User Manual



WEEE Compliance:

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on Thermo Fisher Scientific's compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at <u>www.thermo.com/WEEERoHS</u>

Conformité WEEE:

Ce produit doit être conforme à la directive européenne (2002/96/EC) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est marqué par le symbole suivant:



Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur <u>www.thermo.com/</u>

WEEERoHS

WEEE Konformität:

Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96/ EC erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet:



Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs- / Entsorgungs-anlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Scientific Produkte zu identifizieren, die unter diese RoHS Anweisung fallen, finden Sie unter <u>www.thermo.com/WEEERoHS</u>

Product specifications are subject to change without prior notice. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

See the latest version at www.thermo.com

Les spécifications des produits peuvent être modifiées sans préavis. Voir la dernière version sur **www.thermo.com**

Änderung von Produktbeschreibungen ist vorbehalten ohne vorherige Benachrichtigungt. Sehen Sie die neueste Version an www.thermo.com

1 VisionMate SR Overview

1.1 Specifications				
Dimensions (W x D x	55 x 280 x 160 mm			
Н)	2.17" x 11" x 6.3"			
Weight	Approx 0.8 kg (1.76 lbs)			
Power Supply	110-240 V, 50-60 Hz			
Speed	Approx 5-12 seconds per rack depending on PC performance.			
Decoding Capacity	1 up to 96 tubes.			
Supported Barcode	Any Thermo Scientific 2D Barcoded Storage Tube as well as virtually			
Туре	any other tube with high contrast 2D codes			
Computer	1 GB RAM PC with Microsoft® Windows® 2000, Windows XP or higher			
Requirements	rements (including Windows Vista®) with 2 free USB2 ports.			
Connection Method	Via supplied USB cable to USB2 interface.			
Power Consumption	12 W (operation)			
Technology	Colour CCD image sensor			

- 1.2 Software
 - 1.2.1 So after I've installed the software... what do I do next?
 - Place your rack on the scanner
 - Enable the racks you are using in the software by clicking the enable buttons, then hit scan.
 - 1.2.2 What does the Software do?
 - The VisionMate SR software allows you to scan a rack of Thermo Scientific 2D data matrix-coded tubes and save the results to output files in one simple operation.
 - 1.2.3 How can I store my data?
 - You can save data to Text file, CSV file, MS Excel file, COM Port, XML file, ODBC database and via a TCP/IP connection.
 - The storage of data is designed to be very flexible and easy to configure.
 - Data is exported as soon as a scan is complete without the need for user information.
 - 1.2.4 What configuration changes do I need to make?
 - None for decoding racks.
 - The only configuration required is for export function which is typically a case of selecting what type of file to create and where you want output files to be stored.
 - 1.2.5 Can I use my old software with a new VisionMate SR?
 - No. The hardware is not compatible with older versions of Thermo software.
 - 1.2.6 Can I run the new VisionMate SR on Windows Vista, or Windows 7?
 - Yes. This software is compatible with both Windows Vista and Windows 7.

- 1.3 Hardware
 - 1.3.1 Can I run a SmartScan Solo or VisionMate High Speed reader on the same computer as the VisionMate SR reader?
 - Yes. But you only need to use the security key already included in the VisionMate reader (so the separate key for the Solo is not necessary).
 - 1.3.2 Can I run two VisionMate SRs on the same computer?
 - No
 - 1.3.3 Can I run the VisionMate SR on a Mac?
 - In some cases.....with the latest Mac models and utilizing the latest version of the Mac OS running Windows XP in a virtual machine this is possible.
 - Older models and software or running straight off of the Mac OS is not possible.
 - 1.3.4 Can I use a USB 1.X port to connect the VisionMate SR Scanner?
 - No. The scanner requires a USB 2 port. Using a USB 1.X port will cause very slow scan speeds and could cause erratic performance and application errors.
 - 1.3.5 Do I need to perform maintenance on my VisionMate SR scanner?
 - No, apart from keeping the glass clean and dust-free, including ensuring the corners do not accumulate debris, there is no maintenance necessary.
 - 1.3.6 Can I use an external USB hub to connect the scanner and dongle?
 - Yes, this is normally fine if you use a good quality hub, but if you do begin to experience problems with image quality then the scanner should be connected directly to the PC USB port.
- 1.4 Speed
 - 1.4.1 What affects the speed of my VisionMate SR scanner?
 - The speed of the scanner is generally influenced by the computer hardware.
 - Increasing memory and processor speed will often decrease scanning times.
 - Additionally, using the intensive scan mode will decrease the scan cycle time, due to the increased amount of data being processed by the software.
 - 1.4.2 How long does it take to scan a rack?
 - The scanner will typically read a single rack in 5 to 12 seconds.
 - 1.4.3 What affects decoding speed?
 - The main factor which affects decoding speed is the processing performance of the PC.
 - An unloaded PC which meets the recommended specification should scan a good rack of 96 tubes in seven to eight seconds.

- 1.5 Scanning Different Tube Types/Conditions:
 - 1.5.1 How do I scan different rack sizes?
 - Each rack has a separate rack size configuration box.
 - Either use the drop-down menu to select a common rack size e.g. 96 or 48 tubes, or alternatively use the rows and columns options to set the correct rack configuration.
 - Supported rack sizes are from 1x1 tubes up to 24x16 tubes.
 - 1.5.2 How do I scan racks from different manufacturers?
 - The scanner is configured to read all rack types on the same settings.
 - 1.5.3 Which manufacturers' tubes are supported?
 - All tubes produced by Thermo Scientific are supported as well as most other manufacturers tubes.
 - 1.5.4 Does this reader work with REMP tubes?
 - No. We do not support or promote this product as being compatible with REMP's type of 2D code.
 - 1.5.5 Can the new VisionMate read Matrix 12 mL tubes? 2.0 mL or 5 mL Abgene tubes? Matrix Glass tubes? 14 x 14 codes? Nunc screw cap tubes? 14x14 and 12x12 codes in the same rack?
 - Yes, yes, yes, yes, yes, yes!
 - 1.5.6 Can the VisionMate SR read Matrix 384 tubes?
 - No, Other Thermo Scientific readers are available if you need to read Matrix 384 tubes.
 - 1.5.7 Can I scan very cold racks (from VPLN) on the VisionMate SR scanner?
 - Yes. The scanner contains protection against internal condensation, though if the base of the tubes is highly frosted over we would recommend wiping the base of the tubes with a pre-saturated isopropyl wipe.

2 Parts

- 2.1 Included in this packaging you should find the following:
 - 1) VisionMate SR Scanner
 - 2) Blue License Dongle
 - 3) Installation CD
 - 4) Printed User Manual
 - 5) Power leads (1 each) for US, UK and European use
 - 6) Gray plastic rack locator/template
- 2.2 If you are missing any of the above parts, please contact your Thermo Fisher Scientific representative or their authorized distributor.

3 Installation

3.1 Insert CD ROM and double click on the installer.



3.2 Click "Next" when prompted.

🕏 Thermo VisionMate Flatbed Setup					
	Welcome to the Thermo VisionMate Flatbed Setup Wizard This wizard will guide you through the installation of Thermo VisionMate Flatbed . It is recommended that you close all other applications before starting Setup. This will make it possible to update relevant system files without having to reboot your computer. Click Next to continue.				
	Next > Cancel				

3.3 Read license agreement and click "I Agree".

🐨 Thermo VisionMate SR Setup	
License Agreement Please review the license terms before installing Thermo VisionMate SR.	
Press Page Down to see the rest of the agreement.	
Thermo	^
END USER LICENCE AGREEMENT	
NOTICE TO USER: PLEASE READ THIS AGREEMENT ("AGREEMENT") CAREFULLY. BY USING ALL OR ANY PORTION OF THE SOFTWARE YOU ACCEPT ALL OF THE TERMS AND CONDITIONS OF THIS AGREEMENT. YOU AGREE THAT THIS AGREEMENT IS ENFORCEABLE LIKE ANY WRITTEN NEGOTIATED CONTRACT SIGNED BY YOU. IF YOU DO NOT AGREE, DO NOT USE THIS SOFTWARE. IF YOU ACQUIRED THIS SOFTWARE ON TANGIABLE MATERIAL (EG. CD) WITHOUT AN OPPORTUNITY TO REVIEW THIS LICENCE AND YOU DO NOT ACCEPT THE TERM OF THIS AGREEMENT, YOU MAY	▶
If you accept the terms of the agreement, click I Agree to continue. You must accept the agreement to install Thermo VisionMate SR.	
Thermo — < Back I Agree Cance	el

3.4. Choose the directory and click "Next"

😚 Thermo VisionMate SR Setup	
Choose Install Location Choose the folder in which to install Thermo VisionMate SR.	(and a second se
Setup will install Thermo VisionMate SR in the following folder. To install in a differen click Browse and select another folder. Click Install to start the installation.	t folder,
Destination Folder D:\Program Files\Thermo\Thermo VisionMate SR Browse	
Space required: 6.0KB Space available: 329.7GB	
Thermo	Cancel

3.5 Once the installation is finished, plug in the Dongle and scanner hardware.

🗟 Thermo VisionMate SR Setup					
	Welcome to the Thermo VisionMate SR Setup Wizard				
	This wizard will guide you through the installation of Thermo VisionMate SR.				
	It is recommended that you close all other applications before starting Setup. This will make it possible to update relevant system files without having to reboot your computer.				
57	Click Next to continue.				
	Next > Cancel				

3.6 Follow the Windows hardware wizard, clicking "Next" and "Continue Anyway" when prompted,

Hardwa	Hardware Installation					
1	The software you are installing for this hardware: AVA6 Plus has not passed Windows Logo testing to verify its compatibility with Windows XP. (<u>Tell me why this testing is important</u> .) Continuing your installation of this software may impair or destabilize the correct operation of your system either immediately or in the future. Microsoft strongly recommends that you stop this installation now and contact the hardware vendor for software that has passed Windows Logo testing.					
	Continue Anyway					

3.7 Click "Finish" and now you are now ready to run the software.

Found New Hardware Wizard				
Completing the Found New Hardware Wizard The wizard has finished installing the software for: AVA6 Plus				
	Click Finish to close the wizard.			
	Keinish Cancel			

4 Configuration

- 4.1 Setting the ROI (region of interest) Alignment
 - 4.1.1 The ROI alignment should normally be accurate without the need to modify the default settings.
 - 4.1.2 However, if the ROI alignment looks incorrect or if you experience read failures due to codes not being completely inside their grid squares then it may be necessary to adjust the ROI alignment.
 - 4.1.3 To set the ROI alignment place a rack on the scanner so that you can align the grid to the tubes in the rack.
 - 4.1.4 Open the 'Config' menu, then click 'Set Rack ROI' to show the 'Scan Area Configuration' screen.



4.1.5 Open the 'System' menu, then click 'Scan'. The scanner will scan the full glass area of the scanner.

- 4.1.6 When the scan has completed, the scanned image will be displayed, and the ROI alignment for the Rack will be displayed as a red grid.
- 4.1.7 Change the location and size of the grid, using the mouse, as necessary until all of the codes are centrally located in their grid squares as shown in the image below.



- 4.1.8 When you are satisfied with the alignment of the grid, open the 'System' menu and click 'Save and close' which will save the new settings and close the 'Scan Area Configuration' screen. Alternatively click 'Discard and close' to discard any changes you have made.
- 4.1.9 The changes that you have made will be implemented with immediate effect.
- 4.1.10 The ROI can be manipulated and saved for each rack configuration type (24, 48, 96, etc).

- 4.2 Setting the Image Control Properties
 - 4.2.3 The image control properties allow you to manipulate the scanned image and also manage the way the decoder decodes the scanned image.
 - 4.2.4 It is not normally necessary to change these properties, but if you experience difficulties in scanning racks it is normally possible to improve readability by modifying these parameters.
 - 4.2.5 To show the image control properties tab open the 'Config' menu.
 - 4.2.6 Click the 'Image Control' option. This will display a new tab at the bottom of the screen called 'Image Control'.
 - 4.2.7 The image control tab has two sections: One each for specifying the decoder and scanner properties.

Rack 1 Rack 2 Rack3 Results List Image Control			
8 2↓ □			
Decoder			
Enable Image Filtering:	True		
Enable Varying Threshold:	True		
Varying Threshold Lower Limit:	70		
Varying Threshold Upper Limit:	130		
E Scanner			
Brightness:	-50		
Contrast:	-100		
Resolution Intensive:	400		
Resolution Normal:	200		

Resolution Normal: Scanner Resolution (dpi) for use on normal setting: Max Value is 9600

4.2.8	The decoder properties affect the way the image is decoded once it
	has already been acquired by the scanner.

Property	Default Value	Range	Description	
Enable Image Filtering	False	True or False	This function cyclically sharpens and blurs the image. In some cases, this improves the reading of difficult racks, but in some cases it can prolong the read-time of otherwise quick reading racks. It may be worth enabling and disabling this function to see the effect that it has on your particular racks.	
Enable Varying Threshold	True	True or False	 the effect that it has on your particular racks. Varying threshold is a decoding mode that applies filters to the image to brighten and darken the image. This effect can increase the readability of racks by increasing contrast and reducing the effects of unintentional brightness variations. There is an upper and lower threshold for this mode which can be set as wide as 1 to 255 or can be narrowed depending on the image being decoded. If you dramatically change the brightness or the contrast of the scanned image, then the thresholds may need to also be changed. If varying threshold is not enabled, the scan times for 'easy' to read racks will be quicker, but it is more likely that 	
Varying Threshold Lower Limit	1	1 to 127	The lower threshold for varying threshold mode. The values relate to the darkness of the image with 1 being the darkest.	
Varying Threshold Upper Limit	250	128 to 255	T he upper threshold for varying threshold mode. The values relate to the darkness of the image 255 being the brightest.	

4.2.9 The scanner properties affect the way that the image is acquired by the scanner. These are properties that are sent to the scanner each time a scan is started.

Property	Default Value	Range	Description	
Brightness	200	-999 to 999	This is the brightness of the scanned image.	
Contrast	600	-999 to 999	This is the contrast of the scanned image.	
Resolution (Intensive)	400	100 to 600	This is the resolution (dpi) of the scanner which will be used when intensive scanning mode is selected. Setting the resolution higher will result in more certain decoding at the cost of higher scan and decode times.	
Resolution (Normal)	300	100 to 600	 decoding at the cost of higher scan and decode times. This is the resolution (dpi) of the scanner which will be used when normal scanning mode is selected. Setting the resolution higher will result in more certain decoding at the cost of higher scan and decode times. If you are experiencing no-reads, then try increasing this value in increments of 50 until reliable reading is achieved. 	

- 4.3 Resetting the Software to Default Configuration
 - 4.3.3 It is possible to undo any changes that have been made to the software so that it is configured with the default settings just as straight after installation
 - 4.3.4 To do this shutdown the VisionMate SR software.
 - 4.3.5 Open 'Windows Explorer' or 'My Computer' and navigate to the installation directory which will be something like "C:\Program Files\Thermo\Thermo VisionMate SR"
 - 4.3.6 Within the installation directory, locate the "Config" folder and delete all of the files contained in the config directory (but not the config directory itself)
 - 4.3.7 Restart the VisionMate SR software which will automatically recreate all of the default settings.

5 **Operation**

- 5.1 Rack positions
 - 5.1.1 The scanner has one rack position. Place any rack type on to the scanner insuring the A1 position is in the top left.



- 5.2 Specifying the Rack Type.
 - 5.2.1 Racks are specified by the number of tubes which the rack is designed to hold.

Rows:	Cols:	Quio	k Select
8 😂	12 拿	96	*
Expl	t'd Tube	s: 9	6 🛟

- 5.3 Quick Select
 - 5.3.1 The easiest way of specifying a rack is use the 'Quick Select' selector.
 - 5.3.2 This allows a single-click selection of 24, 48, 96 or 384 tube rack types. Selecting one of these options automatically sets the correct number of rows and columns and sets the expected number of tubes to the full rack size.
- 5.4 Rows and Cols
 - 5.4.1 The Rows and Cols selector allows the specification of any rack size from a 1x1 rack (single tube) up to a 24x16 rack (384 tubes).
- 5.5 Expected Number of Tubes (Expt'd Tubes)
 - 5.5.1 The expected number of tubes function is optional and it is not necessary to set this option.
 - 5.5.2 However, if you are scanning racks which are not full, setting the expected number of tubes to the actual number of tubes in the rack will result in quicker scan times as the software will not waste time looking for tubes which are not present.
- 5.6 Setting the Rack Name
 - 5.6.1 The Rack Name is used to identify each rack both on screen and also optionally in any exported data.
 - 5.6.2 There are five available methods for setting the rack Name. You choose which method you wish to use by selecting it from the 'RackName' menu.
 - 5.6.3 Option 1: Request Before Scan
 - This option displays a rack Name entry box when a scan is started so that the user may manually enter a rack name for all of the enabled racks.
 - The rack Name may optionally be typed in using the keyboard or could alternatively be entered using a keyboard-wedge style barcode scanner.
 - Click the green check button, which allows the scan to proceed using the entered rack Name.
 - Clicking the red X button will cancel the scan operation

📁 Rack Na	ne Entry	X
Rack 1	Thermo Rack	
		:

5.6.4 Option 2: Request After Scan

- This option displays a rack Name entry box when a scan is finished so that the user may manually enter a rack name for all of the enabled racks.
- The rack Name may optionally be typed in using the keyboard or could alternatively be entered using a keyboard-wedge style barcode scanner.
- Click the green check button, which allows the scan to proceed using the entered rack Name.
- Clicking the red X button will cancel the scan operation
- 5.6.5 Option 3: Use Date and Time.
 - This option requires no user intervention during scanning and ensures that each rack is given a unique rack name.
 - The format of the rack name will be similar to '20091228-121242'
 - Where the first part is the date (YYYYMMDD) and the second part is the time (hhmmss).
- 5.6.6 Option 4: Set Via TCPIP.
 - This option should be selected if you are setting the rack name by using the TCPIP server connection.
- 5.6.7 Option 5 Use Fixed Rack Name
 - This option allows you to set a fixed rack name which will be used for every scan.
- 5.7 Starting a Scan
 - 5.7.1 Once the rack is placed on the scanner you can initiate a scan by pressing the start scan button.



- 5.7.2 The scanner will scan the rack, and then the software will load the image and decode the rack.
- 5.7.3 The live decoding of the rack will be displayed in the rack window.
- 5.7.4 Once decoding is complete, the tube-codes will be shown in the results grid at the bottom of the screen and the decoding statistics are shown to the right of the rack window. The decoding time is also displayed in the bottom right of the screen.

- 5.8 Displaying the decoding in full-size window mode.
 - 5.8.1 It is possible to show the decoding window for the rack in a full-size window.
 - 5.8.2 This can be done by clicking the Zoom button above the rack window. This option may also be selected by clicking 'Enable full-screen decoding' on the 'System' menu.
 - 5.8.3 The zoom function has no effect on the decoding performance, but is available only as a tool for visual inspection and informational purposes.
- 5.9 Changing the Scan Resolution
 - 5.9.1 The scanner resolution affects the read performance and the scan speed of the scanner.
 - 5.9.2 The scanner has two default resolutions:

Name	Default Resolution	Description
Normal Resolution	300 dpi	Normal resolution results in reliable decoding of most racks coupled with a fast scan time. If you are experiencing incorrect no-reads, then Intensive resolution should be selected.
Intensive Resolution	400 dpi	Intensive resolution results in very reliable decoding, but with a slight increase in decoding time. If read time is not important, but reliable decoding is essential then this is a good option.

- 5.9.3 The scan resolution is selected using the 'Decode Type' button located on the menu bar at the top of the screen.
- 5.9.4 Pressing the button toggles the resolution between:
 - normal mode (button is displayed in blue)



- intensive mode (button is displayed in orange)
- 5.9.5 It is also possible to change the decode resolution mode by using the menu shortcuts on the 'System' menu.
- 5.9.6 Note: It is possible to change the values for the two resolution modes by changing the values in the image control properties. See Section xxx

- 5.10 Scan Statistics
 - 5.10.1 When a scan is complete then the statistics for the number of tubes read, number of tubes not read and the scan and decode time are shown on screen.

Reads:	96
No Reads:	0
No Tubes:	0

- 5.10.2 The statistics for each rack are shown to the right of the rack window. This shows the number of reads (tubes successfully decoded), the number of 'No Reads' (tubes not decoded), and the number of 'No Tubes' (empty rack location).
- 5.10.3 If the number of reads equals the expected number of tubes then the results are shown in green, otherwise they are displayed in red.

Decode Type: Normal Scan time: 8.391 sec. Decode time: 2.015 sec.

5.10.4 The decode type of Normal or Intensive along with the time taken for the scan (image acquisition) and decoding are displayed at the bottom right of the screen.

	0		0	-	٥	0		4	0	Ø	٥
L.,	Q	8	9	-	3		9	M	-	CI.	0
0	•	0	0	0		-	0	0	0	Q	Ø
Q	۲		B	Q)	8	٢	87	0	0	D	Ø
	3	0		M	0		120		Ø	6	
Ð	\diamond	0	0	-		-	1	8	192		8
8	0		0	٩	8	0	8	0	0		0
TA	-	8	8	Ø	•		0		Ø	Ø	0

5.11 Scan Results: Upon completion of a scan, the scan results are shown in the results tab at the bottom of the screen.

- 5.12 Grid Results.
 - 5.12.1 Rack locations that have been read successfully are shown in green.
 - 5.12.2 Rack locations which resulted in a 'No Read' are shown in orange.
 - 5.12.3 Rack locations which resulted in a 'No Tube' are shown in red.

ſ.	Rack	1 (2009	13603-043627-1	Image Conti	rol									
			1	2	3	4	5	6	7	8	9	10	11	12
	۲.	A	0082862310	0082862309	0082862308	0082860753	0082860747	0082860748	0082862304	0082860745	0082860744	0082862301	0082862300	0082862299
		В	0082862287	0082862288	0082862289	0082862290	0082862291	0082862292	0082862293	0082862294	0082862295	0082862296	0082862297	0082860746
		С	0082862286	0082862285	0082862284	0082862283	0082862282	0082862281	0082862280	0082862279	0082862278	0082862277	0082862276	0082862275
		D	0082862263	0082862264	0082862265	0082862266	0082862267	0082862268	0082862269	0082862270	0082862271	0082862272	0082862273	0082862274
		E	0082860752	0082862261	0082862260	0082862259	0082862258	0082862257	0082862256	0082862255	0082862254	0082862253	0082862252	0082862251
		F	0082862239	0082862240	0082862241	0082862242	0082862243	0082862244	0082862245	0082862246	0082862247	0082862248	0082862249	0082862250
		G	0082862238	0082862237	0082862236	0082862235	0082862234	0082862233	0082862232	0082862231	0082862230	0082862229	0082862228	0082862227
		н	0082860734	0082860733	0082862217	0082862218	0082862219	0082862220	0082862221	0082862222	0082862223	0082862224	0082862225	0082859038

5.13 List Results: After a successful scan the 'Results List' is shown on the far right of the software window, the results are displayed in a list format.

Bow	Column	Code	<u> </u>
A	01	0082862310	
A	02	0082862309	
A	03	0082862308	
A	04	0082860753	
A	05	0082860747	
A	06	0082860748	
A	07	0082862304	
A	08	0082860745	
A	09	0082860744	
A	10	0082862301	
A	11	0082862300	
A	12	0082862299	
В	01	0082862287	
В	02	0082862288	
В	03	0082862289	
В	04	0082862290	
В	05	0082862291	
В	06	0082862292	
В	07	0082862293	
В	08	0082862294	
В	10	0082862295	
Б	10	0002052235	
D	10	0002062237	
D C	12	0002050745	
L	01	0002002200	~

5.14 Printing Scan Results

- 5.14.1 It is possible to print the results for any decoded rack by selecting 'Print', from the 'System Menu.
- 5.14.2 The resulting print-out will be arranged in rack layout similar to the sample output displayed below.

Rack ID 07/09/2009-2

	1	2	3	4	5	6	7	8	9	10	11	12
A	100032664	100032664	100032663	100032640	100032640 7	100032640	100032640	100032640	100032640	100032640	1000326613	100032649
В	100032664	100032641 3	100032664 0	100032649	100032664 5	100032641 6	100032648 3	100032641 4	100032662 0	100032663 5	1000326647	100032664 8
С	100032642	100032641 8	100032643 3	100032643	100032643	100032643 0	100032642 9	100032642 8	100032642 7	100032642 6	1000326470	100032644 9
D	100032644	100032663 8	100032643 8	100032643 9	100032644 0	100032644 1	100032644 2	100032644 3	100032644 4	100032644 5	1000326446	100032645 0
E	100032645	100032645 8	100032648 4	100032645 6	100032645 5	100032645 4	100032645 3	100032645 2	100032647 6	100032645 1	1000326469	100032647 4
F	100032646	100032646 1	100032646 2	100032646 3	100032646 4	100032646 5	100032646 6	100032646 7	100032647 5	100032647 1	1000326491	100032644 7
G	100032649 0	100032648 2	100032667 6	100032648 0	100032647 9	100032647 8	100032647 7	100032646 8	100032643 4	100032645 7	1000326637	100032664 3
н	100032647	100032641	100032664	100032648 7	100032648	100032648	100032641 7	100032649 4	100032649	100032641	25SLHE1A B00168571	100032647

Print Date: 07 September 2009 17:10

Created by Thermo VisionMate Flatbed Scanner

5.14.3 If you wish to print the results in a list layout, simply copy and paste the results from the results list tab into a text editor such as Notepad and print it from there.

5.15 Tube Locator

- 5.15.1 The tube locator function allows you to search for a single tube within the decoded results of the previously scanned rack of tubes.
- 5.15.2 To start the tube locator, select 'Tube Locator' from the 'System' menu.

📁 Tube Locator					
Find Tube:	1000326454				
Result:	Rack: 2 Location: E:6				

- 5.15.3 To locate a tube, enter the full tube code into the 'Find Tube' text box, then click the search button.
- 5.15.4 If the tube is found, its rack number and cell location will be displayed. The tube will also be highlighted in the results grid.

6 Data Export

- 6.1 Background
 - 6.1.1 To enable an Export method, simply click on it in the list of export options on the 'Export To' menu.
 - 6.1.2 When an export method is enabled, its configuration tab will be displayed at the bottom of the screen after the results tabs allowing you to specify exactly how you want the data to be presented.
 - 6.1.3 The available data export methods are:

Export Method	Description
Text File	This function exports to various types of text file including standard text files and comma-delimited text files (*.csv)
Microsoft Excel File	This function exports to native MS Excel Worksheet files (*.xls) which may then be opened directly in Excel. It is not necessary for MS Excel to be installed for this function to work,
	although of course MS Excel is necessary to view the files.
COM Port	This function exports the data to a Serial port.
XML File	This function exports to a XML standard file.
TCP/IP Connection	This function exports data via a TCP/IP connection, allowing communication with another application on the same PC or across the intranet or internet.
ODBC Database	This function exports data to an existing ODBC database. Possible database technologies include MS Access, SQL Server, Oracle and DSN.
VisionMate Compatible CSV File	This function exports data to a CSV file formatted to be compatible with older styles of VisionMate reader output files.

- 6.1.4 It is possible to use a single export method, or any number of different export methods simultaneously.
- 6.1.5 The data is automatically exported at the end of each scan cycle without the need for user intervention, so you should configure the export option for your selected export method before starting a scan.
- 6.1.6 Please see the relevant section of this help manual for a complete description of each export function.
- 6.2 Text File Export
 - 6.2.4 This export function allows the data to be decoded to any of the following text file types:
 - Comma-delimited file (*.csv for Microsoft Excel compatibility)
 - Text file (*.txt for Notepad compatibility)
 - Log file (*.log)
 - 6.2.5 The main difference between the files is the application they normally open with in Windows.
 - 6.2.6 To enable the CSV / Text File Export Function, open the "Export" menu.
 - 6.2.7 Open the "Export To" sub-menu.

6.2.8 Select the "Text File (*.csv)" option. The text file configuration tab will be displayed at the bottom of the screen.

False
False
True
False
True
False
yyyyMM dd
, (Comma)
Row
hh:mm:ss
False
C:\
Use Date and Time
.CSY
Thermo Rack

6.2.9 Choose the appropriate export parameters according to the details below.

6.2.10 The data output is performed automatically at the end of each scan.

DATA EXPORT GROUP

Item	Description
Export Date to each line	Exports the date to each line in the output file based on the date
	format.
Export Rack Name to each Line	Exports the Name to each line in the output file.
Export Rack Name to top of file	Exports the rack name to the top of the file on a single line.
Export Time to each line	Exports the time to each line in the output file based on the time
	format.
Export Tube rack location to each	Exports the tube locations to each line in the output file.
line	
Export User Name to each line	Exports the current user name to each line in the output file.

DATA FORMAT GROUP

Item	Description
Date Format	A drop down box allows the selection of a date format to use when
	exporting the date to the output file. (y = year; M = month; d = day;
	MMM = month abbreviation; MMMM = month full name)
Delimiter Character	A drop down box allows the selection of the delimiter character.
Order Results By	Option to order the results by Column or Row in the output file.
Time Format	A drop down box allows the selection of a time format to use when
	exporting the date to the output file. (h = hour; m = minute; s =
	second; t = AM/PM)
Use Delimiter Character at End of	This option allows the user to include or not include the delimiter
Line	character at the end of each line in the output file.

FILE FORMAT GROUP

Item	Description
Directory	Use the browse button to select a folder in which the output files are to be created.
File Naming Method	 There are three options for naming the csv file: Use Date and Time: This option uses the date and time at the point of file creation to name the file. The date format is as specified in the "Date Format" property. Use Rack Name: This option uses the rack name which was given to the rack during the scan and is probably the most useful option. Use Fixed Filename: This option always names the file with the filename specified in the "Fixed Filename" property. If the file exists, it is overwritten.
Filename Extension	 This property allows you to select the filename extension for the output file. The available options are *.csv (comma-delimited file), *.txt (text file) and *.log (log text file) The *.csv file will normally open automatically in MS Excel where available. The *.txt file will normally open automatically in MS Notepad. The *.log file is not normally associated with any program.
Filename Prefix	It is possible to add a text prefix for the filename here. E.g., if the file naming method is "Use Date and Time" and a "Filename Prefix" of "Test_Rack" is used, then the output file would be named e.g. "Test_Rack01012008.csv" instead of just "01012008.csv"
Fixed File Name	If you select the "File Naming Method" of "Fixed Filename" then you should enter the fixed filename here. E.g. "Test_Rack".

6.3 MS Excel File Export

- 6.3.4 This export function allows the data to be exported to a Microsoft Excel file (*.xls).
- 6.3.5 The export function will create files even if MS Excel is not installed on the system.
- 6.3.6 To enable the MS Excel File Export Function, open the "Export" menu.
- 6.3.7 Open the "Export To" sub-menu.
- 6.3.8 Select the "MS Excel File" option. The configuration tab will be displayed at the bottom of the screen.

2+ E		
Export Date to each line	False	^
Export Rack Name to each line	False	
Export Rack Name to specified cell (below)	False	
Export Rack Name to specified cell: Column	N	
Export Rack Name to specified cell: Row	1	
Export Time to each line	False	
Export Tube Rack Location to each line	True	
🗆 Data Format		
A1 Tube Column	A	
A1 Tube Row	1	
Data Layout Style	Row Layout	
Date format	ddMMyyyy	
Time Format	HH:mm:ss	
🗄 File Format		
Add Sheet if workbook exists	False	
Directory	C:\WisionMate Results	
File Naming Method	Use Rack Name	
Filename Prefix		
Fixed File Name	Thermo Rack	
Overwrite workbook if it exists	False	×
		×
Averwrite workbook if it eviete		

- 6.3.9 Choose the appropriate export parameters according to the details below:
- 6.3.10 The data output is performed automatically at the end of each scan.

DATA EXPORT GROUP

	Item	Description
ſ	Export Date to each line	Exports the date to each line in the output file based on the
		date format.
	Export Rack Name to each Line	Exports the rack Name to each line in the output file.
	Export Rack Name to specified cell	This option enables / disables the option to put the rack name in a cell in the Excel sheet. If enabled, the rack name will be
ļ		put in the cell specified by the column and row positions below.
	Export Rack Name to specified cell:	This is the Excel sheet column in which the rack name should
	Column	be placed.
	Export Rack Name to specified cell: Row	This is the Excel sheet row in which the rack name should be
		placed.
	Export Time to each line	Exports the time to each line in the output file based on the time format. (Only compatible when exporting in either row or column format)
	Export Tube rack location to each line	Exports the tube locations to each line in the output file.
	Export User Name to each line	Exports the current user name to each line in the output file.

DATA FORMAT GROUP

Description	
A1 This option allows you to shift the layout of the data in the Excel sheet. E.g if the	value is "1" 🔹 🗧 – – 🗍 Formatted Table
Tube then the data will be written starting in the first column of the sheet. If the value	s "3" then the
Column data will be written starting in the third column of the sheet, giving you two empty	columns to the
left of the worksheet which could be used for other external data.	
A1 This option allows you to shift the layout of the data in the Excel sheet. E.g if the	value is "A"
Tube then the data will be written starting in the first row of the sheet. If the value is "C	" then the data
Row will be written starting in the third row of the sheet, giving you two empty rows at	he top of the
worksheet which could be used for other external data.	
Data This drop-down option allows the choice of either "Rack Layout" or "Column Lay	out" as a method
Layout of displaying the data in the Excel Worksheet.	
• Rack Layout: The data is displayed in same layout as if they were in the rac	k. Therefore
the tube in position E4 in the rack will be displayed in cell E4 in the workshe	et. (subject to
the above two properties being set to a start cell of A1)	
 Column Layout: The data is displayed in vertical columns ordered by the co 	umn letter of
the tube in ascending order. Each column containing separate information f	elds, e.g., code,
date, tube location.	
Row Layout: The data is displayed in vertical columns ordered by the row n	umber of the
tube in ascending order. Each worksheet column contains separate information	tion fields, e.g.
code, date, tube location.	
Formatted Rack Layout: The data is displayed in the same orientation as in	the physical
rack. The data includes column and row headers	
Date A drop down box allows the selection of a date format to use when exporting the	date to the
Format output file. (y = year; M = month; d = day; MMM = month abbreviation; MMMM =	month full
name)	
Time A drop down box allows the selection of a time format to use when exporting the	date to the
Format output file. (h = hour; m = minute; s = second; t = PM/AM)	

FILE FORMAT GROUP

Item	Description
Add sheet	If this option is set to true then each time a rack is exported to an existing MS Excel workbook
if	a new worksheet page will be added to store the data,
workbook	
exists	
Directory	Use the browse button to select a folder in which the output files are to be created.
File	There are three options for naming the Excel file:
Naming	Use Date and Time: This option uses the date and time at the point of file creation to
Method	name the file. The date format is as specified in the "Date Format" property.
	Use Rack Name: This option uses the rack name which was given to the rack during the
	scan and is probably the most useful option.
	• Use Fixed Filename: This option always names the file with the filename specified in the
	"Fixed Filename" property.
Filename	It is possible to add a text prefix for the filename here. E.g., if the file naming method is "Use
Prefix	Date and Time" and a "Filename Prefix" of "Test_Rack" is used, then the output file would be
	named e.g. "Test_Rack01012008.xls" instead of just "01012008.xls"
Fixed File	If you select the "File Naming Method" of" Fixed Filename" then you should enter the fixed
Name	filename here. E.g., "Test_Rack".
Overwrite	If this option is true, then if a rack is exported to an existing MS Excel workbook, the existing
workbook	MS Excel file will be deleted and recreated with the new rack data.
if it exists	

6.4 COM Port Export

- 6.4.4 This export function allows the data to be outputted to an onboard COM Port
- 6.4.5 To enable the COM Port Function, open the "Export" menu.
- 6.4.6 Open the "Export To" sub-menu.6.4.7 Select the "COM Port" option to bring up the COM port configuration tab.

Rack 1 Rack 2 Rack 3 Results List COM Port	
81 2↓ 🖾	
🗆 Data Export	
Export Date to each line	True
Export Rack Name to each line	False
Export Rack Name to top of file	True
Export Time to each line	True
Export Tube Rack Location to each line	True
Export User Name to each line	False
🗆 Data Format	
Date format	ууууMMdd
Delimiter Character	, (Comma)
Order results by	Column
Prefix Character	[13]
Suffix Character	[10]
Time Format	hh:mm:ss
Use Delimiter Character at line end	True
🗆 File Format	
File Naming Method	Use Date and Time
Rackname Prefix	
Port Config	
Baud Rate	9600
COM Port	COM1
Data Bits	8
Handshaking	None
Parity	None
Stop Bits	1
Baud Rate Baud rate setting for the CDM Port	

- 6.4.8 Choose the appropriate export parameters according to the details below.
- 6.4.9 The data output is performed automatically at the end of each scan.

DATA EXPORT GROUP

Item	Description
Export Date to each line	Exports the date to each line in the output file based on the date format.
Export Rack Name to	Exports the rack name to each line in the output file.
each Line	
Export Rack Name to	Exports the rack name to the top of the file on a single line.
top of file	
Export Time to each line	Exports the time to each line in the output file based on the time format.
Export Tube rack	Exports the tube locations to each line in the output file.
location to each line	
Export User Name to	Exports the current user name to each line in the output file.
each line	

DATA FORMAT GROUP

Item	Description
Date Format	A drop down box allows the selection of a date format to use when exporting the
	date to the output file. (y = year; M = month; d = day; MMM = month
	abbreviation; MMMM = month full name)
Delimiter Character	A drop down box allows the selection of the delimiter character.
Order Results By	Option to order the results by Column or Row in the output file.
Prefix Character	If you wish to prefix the output with a character then it should be entered here.
	The prefix can be plain text e.g. "ABC", or it can be an ASCII character code
	(see www.asciitable.com). To specify an ASCII character, enclose the ASCII
	value with square brackets. E.g., [13] would add a carriage return as the prefix.
Suffix Character	If you wish to suffix the output with a character then it should be entered here.
	The suffix can be plain text e.g. "ABC", or it can be an ASCII character code (see
	www.asciitable.com). To specify an ASCII character, enclose the ASCII value
	with square brackets. E.g., [13] would add a carriage return as the suffix.
Time Format	A drop down box allows the selection of a time format to use when exporting the
	date to the output file. (h = hour; m = minute; s = second; t = PM/AM)
Use Delimiter Character	This option allows the user to include or remove a delimiter character at the end
at End of Line	of each line in the output file.

FILE FORMAT GROUP

Item	Description
File Naming Method	There are three options for naming the csv file:
	Use Date and Time: This option uses the date and time at the point of file
	creation to name the file. The date format is as specified in the "Date Format"
	property.
	Use Rack Name: This option uses the rack name which was given to the rack
	during the scan and is probably the most useful option.
	Use Fixed Filename: This option always names the file with the filename
	specified in the "Fixed Filename" property. If the file exists, it is overwritten.
Rackname Prefix	It is possible to add a text prefix for the rack name here. E.g., if the file naming method is "Use Date and Time" and a "Filename Prefix" of "Test Rack" is used
	then the output file would be named e.g. "Test_Rack01012008.csv" instead of
	just "01012008.csv".

PORT CONFIG GROUP

Item	Description
Baud Rate	This is the transmission speed used for the COM Port communication. This should match the baud rate of the receiving COM port.
COM Port	This is the physical COM port to use for data export. If you change the COM port that you are using, you must then disable and re-enable the Export to COM port option to make the change.
Data Bits	This is the number of data bits used in the data communication structure. This should match the number of data bits used by the receiving COM port.
Handshaking	This is the type of handshaking, or flow control used by the COM port. The handshaking method should match the type used by the receiving COM port.
Parity	This the error detection method used by the COM port. The parity method should match the type used by the receiving COM port.
Stop Bits	This is the number of stop bits used in the data communication structure. This should match the number of stop bits used by the receiving COM port.

- 6.5 XML File Export
 - 6.5.4 This export function allows the data to be exported to an "Extensible Markup Language file" (*.xml).
 - 6.5.5 XML files are text based files conforming to a standard and are increasingly used to share information between platforms in a common format.
 - 6.5.6 To enable the XML file Function, Open the "Export" menu.
 - 6.5.7 Open the "Export To" sub-menu.
 - 6.5.8 Select the "XML" option to bring up the XML configuration tab.

utput Files
and Time
sk
0

- 6.5.9 Choose the appropriate export parameters according to the details below.
- 6.5.10 The data output is performed automatically at the end of each scan.

DATA EXPORT GROUP

Item	Description
Export Date to each line	Exports the date to each line in the output file based on the date format.
Export Rack Name to each	Exports the rack name to each line in the output file.
Line	
Export Rack Name to top of	Exports the rack name to the top of the file on a single line.
file	
Export Time to each line	Exports the time to each line in the output file based on the time format.
Export Tube rack location to	Exports the tube locations to each line in the output file.
each line	
Export User Name to each	Exports the current user name to each line in the output file.
line	

DATA FORMAT GROUP

Item	Description
Order Results By	Option to order the results by Column or Row in the output file.
Date Format	A drop down box allows the selection of a date format to use when exporting the date to the output file. ($y = year$; $M = month$; $d = day$; MMM = month abbreviation; MMMM = month full name)
Time Format	A drop down box allows the selection of a time format to use when exporting the date to the output file. (h = hour; m = minute; s = second; t = PM/AM)

FILE FORMAT GROUP

Item	Description
Directory	Use the browse button to select a folder in which the output files are to
	be created.
File Naming Method	There are three options for naming the Excel file:
	Use Date and Time: This option uses the date and time at the point of
	file creation to name the file. The date format is as specified in the
	"Date Format" property.
	Use Rack Name: This option uses the rack name which was given to
	the rack during the scan and is probably the most useful option.
	Use Fixed Filename: This option always names the file with the
	filename specified in the "Fixed Filename" property.
Filename Prefix	It is possible to add a text prefix for the filename here. E.g., if the file
	naming method is "Use Date and Time" and a "Filename Prefix" of
	"Test_Rack" is used, then the output file would be named
	"Test_Rack01012008.xls" instead of just "01012008.xls".
Fixed File Name	If you select the "File Naming Method" of "Fixed Filename" then you
	should enter the fixed filename here. E.g., "Test_Rack".

Remote (TCP/IP) Operation 7

- 7.1
- Viewing the TCP/IP Configuration Tab
 7.1.1 Open the "Server" menu.
 7.1.2 Select the "Show Config Tab" option to bring up both the TCP/IP configuration tab and the TCP/IP Log tab.

Rack 1 TCPIP Config TCPIPLog	
🗆 Command Set	
Acknowledge Text	OK
Change Product	P
Get / Set Expected No of Tubes	A
Get Current Product	C
Get Data	D
Get Decode Time	ĸ
Get Number of No Reads	6
Get Number of No Tubes	н
Get Number of Reads	F
Get Rack Name	B
Get Scanner Status	L
Get String Status	J
Set Rack Name	I contraction of the second seco
Start Scan	S
Data Export	
Export Date to each line	False
Export Rack Name to each line	False
Export Time to each line	False
Export Tube Rack Location to each line	True
🗆 Data Format	
Date format	ууууMMdd
Delimiter Character	, (Comma)
Echo received commands	True
Order results by	Column
Prefix Character	
Send Data Length	False
Suffix Character	
Time Format	hh:mm:ss
Use Delimiter Character at line end	True
Port Configuration	
IP Address mode	Automatic IP Address
Port Number	8000
User Defined IP Address	
Acknowledge Text This text is returned by the scanner to acknowledge commands	

COMMAND SET GROUP

Item	Description
Acknowledge	This is the response which will be returned by the VisionMate Server to
Text	acknowledge a command from the client. The response format is [Acknowledge
	Text][Command]. E.g., if you send an "S" the response will be "OKS"
Change Product	This is a command to change and load a new product. The format of the
	command is [Change Product Command] [Number of Rows][Number of Columns].
	E.g., To change to a 48 tube rack, the command is "P0806".
Get / Set	This command sets the expected number of tubes. The format of the command is
Expected No of	[Command][Number of tubes].
Tubes	E.g. A96 will set the expected number of tubes to 96.
Get Current	The currently loaded product is returned
Product	The response structure is [Number of Rows][Number of Columns]
	E.g., The response for a 96 tube rack would be "0806".
Get Data	The last scanned rack data is returned.
	The format of the data is dependant on the settings made in the data format group.
	Command Structure: DE.g.'D' returns the data for rack.
Get Decode Time	This returns the decode time of the last scan in seconds.
	Command Structure: K
	E.g The response may be OKK2.431
Get Number of No	This returns the number of No Reads for the last scanned rack.
Reads	Command Structure: G
	E.g. The reponse for 2 no reads might ne OKG2
Get Number of No	This returns the number of No Tubes for the last scanned rack.
Tubes	Command Structure: H
	E.g. The reponse for 2 no tubes might ne OKH2
Get Number of	This returns the number of tubes read for the last scanned rack.
Reads	Command Structure: F
	E.g. The reponse for 96 reads might ne OKF96
Get Rack Name	The Rack name is returned
	Command Structure: G
	E.g. G might return 'OKGThermo_Rack' if the rack name is 'Thermo_Rack'
Get Scanner	A byte value is returned indicating the scanner status based on status bits as
Status	described below
	0 = Initialised
	1 = Scanning
	2 = Finished scan
	3 = Data ready
	4 = Data sent
	5 = Rack96
	6 = Empty
	7 = Error
	E.g., if the scanner is initialized and is scanning a 96 rack, but the data is not ready or sent and there is no error, then the value would be 100011. The returned value is then 35 (which is the decimal value of 10011). Then the scan is finished and the data is ready, the bits would be 101101 and the return value is 45.
	 6 = Empty 7 = Error E.g., if the scanner is initialized and is scanning a 96 rack, but the data is not ready or sent and there is no error, then the value would be 100011. The returned value is then 35 (which is the decimal value of 10011). Then the scan is finished and the data is ready, the bits would be 101101 and the return value is 45.

Get String Status	 This returns a string representation of the status of the scanner. The possible return values are: Error : The scanner is in an error state. Ready: The scanner is ready to scan after initialization. Scanning: The scanner is currently scanning and decoding a rack. DataReady: The scan is complete and data is ready. The scanner is ready to perform the next scan Command Structure: J E.j. Whilst scanning the command might return OKJScanning
Set Rack Name	This command sets the Rack name. Command Structure: I[Rack Name] E.g.' IThermo_Rack' sets the name of the rack to 'Thermo_Rack'
Start Scan	The command to start scanning a rack.

DATA EXPORT GROUP

Item	Description
Export Date to each line	Exports the date to each line based on the date format.
Export Rack Name to each	Exports the rack name to each line in the output file.
Line	
Export Rack Name to top of file	Exports the rack name to the top of the file on a single line.
Export Time to each line	Exports the time to each line in the output file based on the time format.
Export Tube rack location	Exports the tube locations to each line in the output file.
to each line	

DATA FORMAT GROUP

Item	Description
Date Format	A drop down box allows the selection of a date format to use when
	exporting the date to the output. ($y = year$; $M = month$; $d = day$; MMM
	= month abbreviation; MMMM = month full name)
Delimiter Character	A drop down box allows the selection of the delimiter character.
Order Results By	Option to order the results by Column or Row in the output.
Prefix Character	If you wish to prefix the output with a character then it should be
	entered here.
	The prefix can be plain text e.g. "ABC", or it can be an ASCII
	character code (see www.asciitable.com). To specify an ASCII
	character, enclose the ASCII value with square brackets. E.g., [13]
	would add a carriage return as the prefix.
Suffix Character	If you wish to suffix the output with a character then it should be
	entered here. The suffix can be plain text e.g. "ABC", or it can be an
	ASCII character code (see www.asciitable.com). To specify an ASCII
	character, enclose the ASCII value with square brackets. E.g., [13]
	would add a carriage return as the suffix.
Time Format	A drop down box allows the selection of a time format to use when
	exporting the date. (h = hour; m = minute; s = second; t = PM/AM)
Use Delimiter Character at End	This option allows the user to include or remove a delimiter character
of Line	at the end of each line in the output.

PORT CONFIGURATION GROUP

Item	Description
IP Address Mode	 This is the IP address which the on which the server will listen for incoming connections. There are two options: Loopback Address: This uses the local loopback address 127.0.0.1. This can only be used, and is the recommended method, if the client is on the same machine. Automatic IP Address: This option automatically retrieves and uses the primary IP address of the server PC.
Port Number	This is port on which the server will listen for incoming connections. The valid range is 1024-65536.

- 7.2 Enabling the TCP/IP Server
 - 7.2.1 Open the "Export" menu.
 - 7.2.2 Select the "Enable TCP/IP Server" option.
 - The TCP/IP server will start and wait for a client connection.
 - The IP address and port number in use will be displayed in the TCP/IP Log screen along with a message confirming that the server has started.

Rack 1 [08/09/2009-1] Rack 2 Rack3 Results List TCPIP Config TCPIPLog	
I.P. Address: 192.168.1.12 Port: 8000	
198/09/2009 01:07:52 Client Connected: 192.168.1.12 08/09/2009 01:07:43 Server Started	
	~

- 7.3 Testing the TCP/IP Server with HyperTerminal
 - 7.3.1 Within the VisionMate software, display the TCP/IP configuration and log tabs as detailed in section 10.1 above.
 - 7.3.2 Enable the TCP/IP server as shown in section 10.2 above.
 - 7.3.3 Note the IP Address and port number which are being used. This information is shown on the TCP/IP Log tab.
 - 7.3.4 Start HyperTerminal.
 - go to Windows Start Menu All Programs Accessories Communications- HyperTerminal
 - 7.3.5 Enter a meaningful name for the connection (E.g., "VisionMate Test"), and select "OK".

7.3.6 Choose to connect using TCP/IP (Winsock), enter the IP Address and port number previously noted in step 3 above and select "OK".

Connect To	? 🛛
🧞 VisionMa	ate Test
Enter details for	the host that you want to call:
<u>H</u> ost address:	192.168.1.7
Port nu <u>m</u> ber:	8000
Co <u>n</u> nect using:	TCP/IP (Winsock)
	OK Cancel

7.3.7 HyperTerminal will connect to the server and the VisionMate TCP/IP log tab will display a "Client Connected" message.



- 7.3.8 Within the HyperTerminal window, type an uppercase "S" and press return. The server will acknowledge the command by responding with "OKS" and the scan will start.
- 7.3.9 To check that the scan has finished, send the Get Status command by typing "L" and pressing enter. When the scan has finished, the response will be "OK45".
- 7.3.10 Retrieve the data from the rack by typing "D" and pressing enter. The results will be returned.
 - Note: If an unknown command is received, the response will be "OK?[Unknown Command Text]".

8 Maintenance

- 8.1 The only maintenance required is to remove any debris from reader surface
 - 8.1.1 Remove the Rack locator template
 - 8.1.2 Using a wet, non-abrasive cloth, remove any debris wiping gently
 - 8.1.3 Alternately canned air can be used to remove any debris
- 8.2 Service/Maintenance
 - 8.2.1 For technical assistance or to reorder contact your local Thermo Fisher Scientific representative or authorized distributor.

9 <u>Disposal</u>

9.1 The disposal of this equipment is responsibility of the customer, and must be in an environmentally responsible manner.

10 Warranty

10.1 A 12 month warranty is included within the price of the unit.

11 Troubleshooting

- 11.1 I am getting no-reads or unreliable reading
 - <u>Is the scanner glass clean</u>? If not, remove the plastic rack locator and clean the glass scan area.
 - <u>Is the plastic rack locator placed correctly on the scanner glass and are the</u> <u>racks correctly within the rack locator holes</u>?
 - <u>Are the bottoms of the tubes dirty, damaged or obscured by frost</u>? If the tubes are frosted it is advised to wipe them with an alcohol wipe. If the tubes are damaged then they may have to be replaced.
 - <u>Are the codes aligned well within the grid</u>? If not, see the section on ROI alignment.
 - If you are using normal decode mode, do the no-reads disappear if you use <u>intensive decoding mode</u>? If so, try increasing the resolution for the normal decoding mode. See the section on setting the Image Control Properties. If not try using intensive decoding again after increasing the intensive resolution in the image control properties.
 - <u>Is the PC struggling to cope and running very slowly</u>? If so, new hardware or a hardware upgrade may be required. Try closing other running applications if they are using a lot of processing power.
 - <u>Does the image look very bright or very dark</u>? If so, try changing the brightness and / or contrast levels. See the section on setting the Image Control Properties.
- 11.2 I cannot get the TCP/IP Server to work
 - <u>Is the TCP/IP Server enabled</u>? Check in the TCP/IP Log tab to make sure that an I.P address and port are shown and that the TCP/IP Server has started.
 - <u>Are you using the correct IP address and Port number</u>? Check in the TCP/IP Log tab to make sure that the I.P address and port being used match the ones that you are connecting with.
 - <u>Can you 'ping' the IP address</u>? Open a dos command prompt and ping the ip address by entering ping [IP address] and pressing enter. E.g. if the TCP/IP server is using IP address 192.168.1.1 then type 'ping 192.168.1.1'

into the command prompt and press enter. You should receive a number of replies. If you don't receive replies, then there is a problem either with the TCP/IP network or firewalls which must be solved before the TCP/IP Server will be able to function.

- <u>Does the scanner respond if you use HyperTerminal</u>? Try following the example which explains how to use HyperTerminal to control the scanner.
- <u>Are you sending a carriage-return character after the messages</u>? The TCP/IP server looks for an ASCII character 13 [CR] as a message terminator.

12 Contact Details

North America Contact (Main Office): Thermo Fisher Scientific 22 Friars Drive Hudson, NH 03051 PH: (603) 595-0505 FAX: (603) 595-0106

Europe Contact: Thermo Fisher Scientific Unit 2 Lower Meadow Road Brooke Park Handforth Wilmslow SK9 3LP England PH: +44 (0) 161 486 2110

Asia Pacific Contact: matrix.ap.info@thermofisher.com