

# **Fluoromark™ Microplate Fluorometer and Fluoromark Software**

## **Instruction Manual**

**Catalog Numbers  
170-6941, 170-6966,  
170-6970, 170-6942,  
170-6967, 170-6971,  
170-6943, 170-6968,  
170-6972, 170-6944,  
170-6969 and 170-6973**



# Warranty and Regulatory Notices

## Warranty Statement

This warranty may vary outside of the continental United States. Contact your local Bio-Rad office for the exact terms of your warranty.

Bio-Rad Laboratories warrants that the Fluoromark system will be free from defects in material and workmanship, and will meet all performance specifications for the period of 1 year from the date of shipment. This warranty covers all parts and labor.

In the event that the instrument must be returned to the factory for repair under warranty, the instrument must be packed for return in the original packaging.

Bio-Rad shall not be liable for any incidental, special, or consequential loss, damage, or expense directly or indirectly arising from the use of the Fluoromark system. Bio-Rad makes no warranty whatsoever in regard to products or parts furnished by third parties, such being subject to the warranty of their respective manufacturers. Service under this warranty shall be requested by contacting your nearest Bio-Rad office.

This warranty does not extend to any instruments or parts thereof that have been subject to misuse, neglect, or accident, or that have been modified by anyone other than Bio-Rad or that have been used in violation of Bio-Rad instructions.

The foregoing obligations are in lieu of all other obligations and liabilities including negligence and all warranties, of merchantability, fitness for a particular purpose or otherwise, expressed or implied in fact or by law, and state Bio-Rad's entire and exclusive liability and buyer's exclusive remedy for any claims or damages in connection with the furnishing of goods or parts, their design, suitability for use, installation, or operation. Bio-Rad will in no event be liable for any special, incidental, or consequential damages whatsoever, and Bio-Rad's liability under no circumstances will exceed the contract price for the goods for which liability is claimed.

## Regulatory Notices

**Important:** This Bio-Rad instrument is designed and certified to meet EN55011, EN50082-1, and IEC 1010-1 requirements, which are internationally accepted electrical safety standards. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will:

Void the manufacturer's warranty.

Void the regulatory certifications.

Create a potential safety hazard.

**Note:** This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

# Table of Contents

<b>Section 1</b>	<b>Introduction</b> .....	<b>1</b>
1.1	General Information .....	1
1.2	Specifications .....	1
<b>Section 2</b>	<b>Equipment Set-up and Installation</b> .....	<b>2</b>
2.1	Unpacking the Fluoromark Microplate Fluorometer .....	2
2.2	RS232 Connection to PC .....	2
2.3	Power Cord Connection .....	2
2.4	Location of Instrument .....	2
<b>Section 3</b>	<b>Maintenance</b> .....	<b>3</b>
3.1	Disinfection Procedure .....	3
<b>Section 4</b>	<b>Software Installation</b> .....	<b>4</b>
4.1	Installing Excel 5.0 Software .....	4
4.2	Installing the Fluoromark Software .....	5
<b>Section 5</b>	<b>Using the Fluoromark Windows PC Software</b> .....	<b>8</b>
5.1	Hints for Operation .....	8
5.2	General View of the Menus and the Corresponding Icons .....	8
5.3	Performing a Measurement .....	11
5.4	Incubating Plates (if installed) .....	14
5.5	Defining a Layout .....	15
5.6	Defining a Test .....	20
<b>Section 6</b>	<b>Configuration of the Fluoromark Microplate Fluorometer and Definition of Microplates</b> .....	<b>22</b>
6.1	Serial Port for Communication PC to Fluoromark .....	22
6.2	Adjusting the Mechanical Offset .....	23
6.3	Definition of the Mechanical Dimensions of Microplates .....	24
6.4	Definition of the Filter Configuration .....	26
<b>Section 7</b>	<b>Microsoft Excel Data Reduction Routines</b> .....	<b>27</b>
7.1	General Hints for Operation .....	27
7.2	Data Reduction in EXCEL .....	27
<b>Section 8</b>	<b>Error Codes</b> .....	<b>32</b>
<b>Section 9</b>	<b>Product Information</b> .....	<b>33</b>
9.1	Product Information .....	33
9.2	Filter Table .....	34
9.3	Disinfection Certificate .....	34

# Section 1

## General Information

### 1.1 Introduction

The Fluoromark microplate fluorometer provides all the flexibility you need for virtually any fluorescent assay. You can use it to perform top and bottom epifluorescence for detection in solution assays or cell monolayers bound to well bottoms, and use 6, 12, 24, 48, 96, and 384 well plate formats. With the widest working wavelength range of any microplate fluorometer (250–900 nm), the Fluoromark fluorometer lets you employ fluorophores such as dansylamide for protein quantitation or Cy® dyes for DNA quantitation. You can run both time-resolved studies using lanthanide chelates and standard fluorescent determinations. For assays requiring temperature control, the optional incubator features a range from ambient to 45 °C. Features of the Fluoromark system include

- Perform both standard and time-resolved fluorescent measurements
- Read either from the top or the bottom of the plate
- Use any microplate format including 6, 12, 24, 96, and 384 well plates
- Optional incubation capabilities
- Extended ultraviolet read capability down to 250 nm
- Cleaner, economical, more sensitive alternative to radioactive assays

### 1.2 Specifications

Light source	High energy Xenon flash lamp
Detector	Side window, current type photomultiplier tube
Filters	Excitation filter wheel for 8 interference filters Emission filter wheel for 8 interference filters
Excitation wavelength range	290–900 nm (250 nm optional)
Emission wavelength range	300–900 nm
Raw result range	0–9999
Sensitivity	5 pg/well sodium fluorescein
Incubation system	Ambient to 42 °C (optional)
Interface	RS232 bi-directional
Power	150 VA 100/120 V and 220/250 V Fuses: T2.5 A 250 V for main power 230 V T1.6 A 250 V for main power 115 V
Physical	20 x 44 x 48 cm (h x w x d)
Weight	23 Kg
Operating Conditions	Operating temperature 15–30 °C Storage temperature -10–50 °C Humidity 20%–80%

This instrument conforms to Overvoltage Category II and Contamination Class II

## Section 2

### Equipment Set-up and Installation

#### 2.1 Unpacking the Fluoromark Microplate Fluorometer

1. Before opening the box, check it for any outer damage. Make a note of any damages, if present.
2. Place the box upright with the TOP label facing up, and open the container. Remove the instrument from the box in this position only.

**Warning:** Grasp the instrument firmly under the right and left sides before attempting to lift or carry it.

3. Check the instrument for damaged or broken parts. Note specifically any damages, if they are present.
4. Each Fluoromark instrument includes the following accessories. Verify that they are present.
  - 1 power cord
  - 1 RS232 computer cable
  - 1 instruction manual
  - 3 Software diskettes
  - 1 box containing 2 spare fuses
  - 3 Allen wrenches
  - Excel Software
5. Verify that the serial number on the back of the instrument is the same as that on the outside of the box. Check also that the main voltage shown on the back of the instrument corresponds to the local voltage in your country.

#### 2.2 RS232 Cable Connection

The RS232 port is on the back of the instrument. The included interface cable will connect the Fluoromark microplate fluorometer to COM1 of your PC.

**Note:** Communication parameters of the PC interface are automatically configured by the software.

#### 2.3 Power Connection

1. Plug the included power cord into the main power connector port. Do not switch power on yet. Make sure that the instrument is connected to a grounded conductor of the main power outlet.
2. Before powering the instrument up for the first time, allow it to warm to room temperature for 6 hours, to prevent any condensation problems.

#### 2.4 Location of the Instrument

The Fluoromark microplate fluorometer is a precise optical measuring instrument. The following conditions must be met to insure maximum sensitivity.

1. Place the instrument on a flat and vibration free surface.
2. Do not place the instrument in direct sunlight.
3. Insure that the location is dust free.

## **Section 3 Maintenance**

### **3.1 Disinfection Procedure**

All parts of this instrument that can come in contact with patient sera or positive samples must be handled as hazardous items. Bio-Rad recommends using gloves when performing maintenance or working with the instrument.

**Note:** It is very important that the instrument undergo thorough disinfection before performing maintenance or before physical removal of the instrument from the laboratory. The instrument must be disinfected before you send it back to Bio-Rad for service. For safety reasons, you must fill out a Disinfection Certificate and include it with the instrument. Without a Disinfection Certificate, Bio-Rad will not accept any returned instruments.

Bio-Rad recommends using a solution of 10% formaldehyde, 70% alcohol, and 20% deionized water. Make sure to observe all national regulations for safe handling of formaldehyde. The disinfection procedure must be performed by authorized persons wearing one way gloves and protective clothes. The location must possess adequate ventilation.

To disinfect Fluoromark, use the following procedure.

**Note:** Be sure that you wear disposable gloves.

1. Disconnect the instrument from the main power supply.
2. Remove the RS232 cable from its port.
3. Clean all the outside surfaces of the instrument carefully with a wad of cotton which has been soaked in the formaldehyde solution.
4. Place the instrument in a sealed plastic bag.
5. Repeat the procedure for disinfection on any accessories which are also being returned with the instrument.
6. Complete the Certificate of Disinfection and send it in with the instrument

## Section 4 Software Installation

The Fluoromark Excel software runs on IBM compatible PC processors under Microsoft Windows® Version 3.1, 3.11 and '95. The CPU must be a 486 or higher with minimum 4 Mb RAM. The data reduction is done in the Microsoft spread sheet program Excel version 5.0c. Before installing Fluoromark software, verify proper installation of Excel 5.0c.

### 4.1 Installing Excel 5.0 Software

The data base engine, Microsoft Query, must be installed during Excel installation.

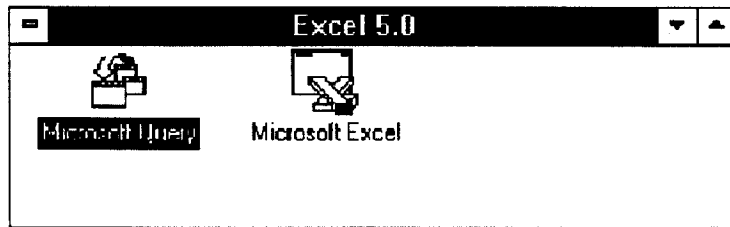


Fig. 4.1. Required components of Excel software.

If Microsoft Query is not installed on your computer, you must reinstall Excel specifying Query, or Data Access during custom installation.

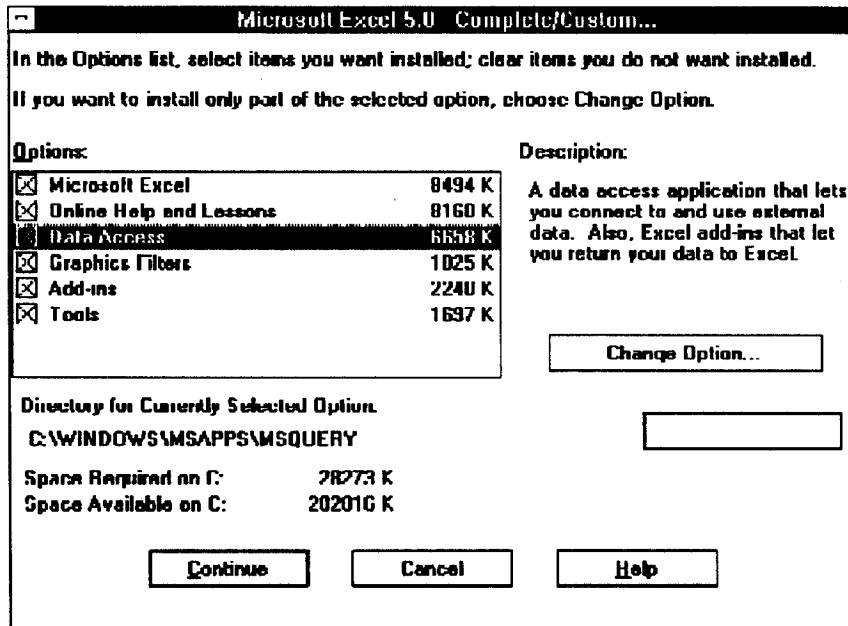
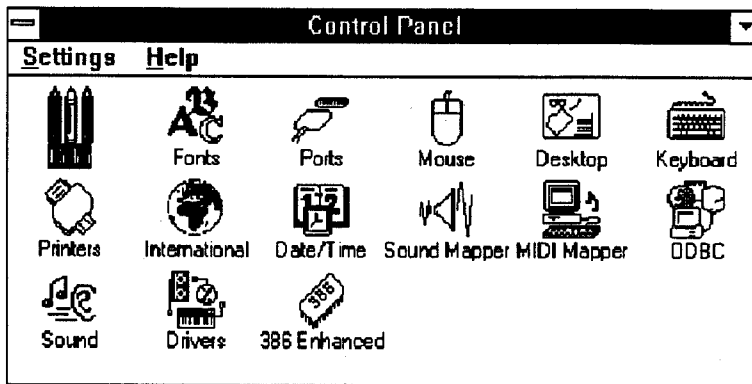


Fig. 4.2. Reinstallation of the data base handler Microsoft Query.

After proper installation of Excel; the Microsoft Query icon should appear in the Excel program group, and the data base driver ODBC should appear in the System Control program group.



Changes the Windows screen colors

Fig. 4.3. ODBC driver in system control.

## 4.2 Installing the Fluoromark Software

After verifying installation of Excel 5.0c, install the Fluoromark software by placing the installation diskette in the drive, and running the Setup program from the Program Manager.

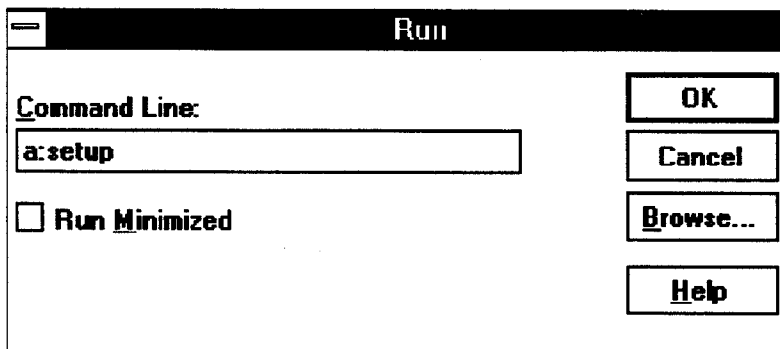


Fig. 4.4. Starting the Fluoromark installation.

The installation program will prompt for the directory in which to install the Borland database tools.

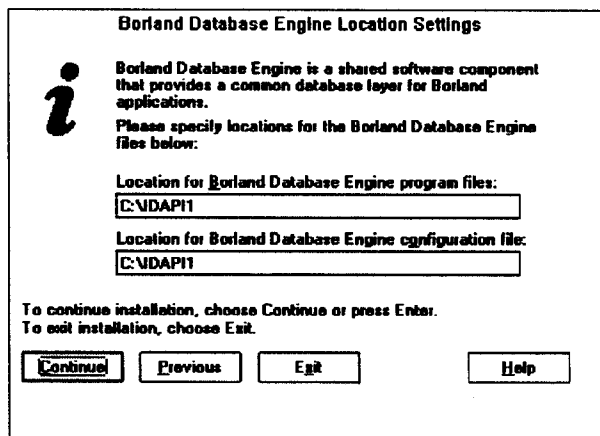


Fig. 4.5. Directory for the database.

The database of the Fluoromark software will be installed in the selected directory, from diskettes 1 and 2. Next install the Fluoromark reader software by running the Setup program on diskette 3.

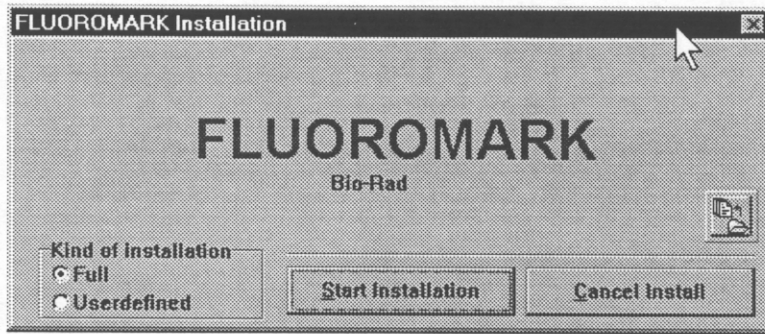


Fig. 4.6. Fluoromark installation screen.

After selecting the Start Installation option, the setup program prompts for the directory in which to install the Fluoromark program files.

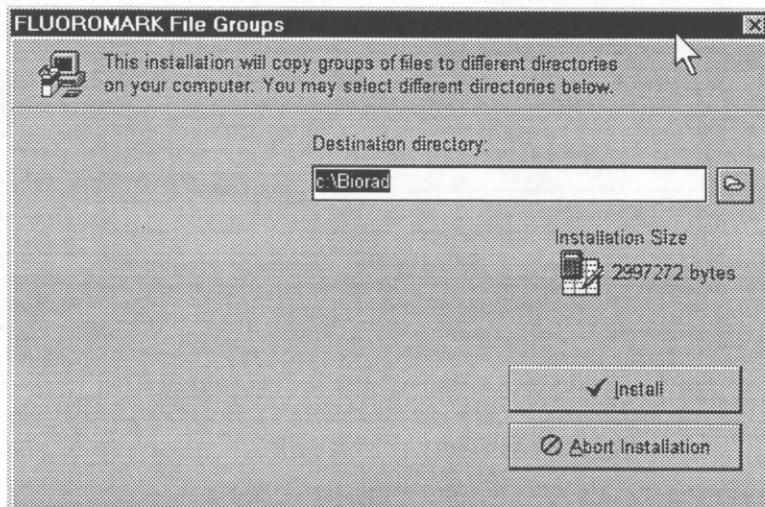


Fig. 4.7. Directory for the Fluoromark software.

After installation of the Fluoromark software, a Fluoromark icon will be created in a Program Group named Fluoromark.

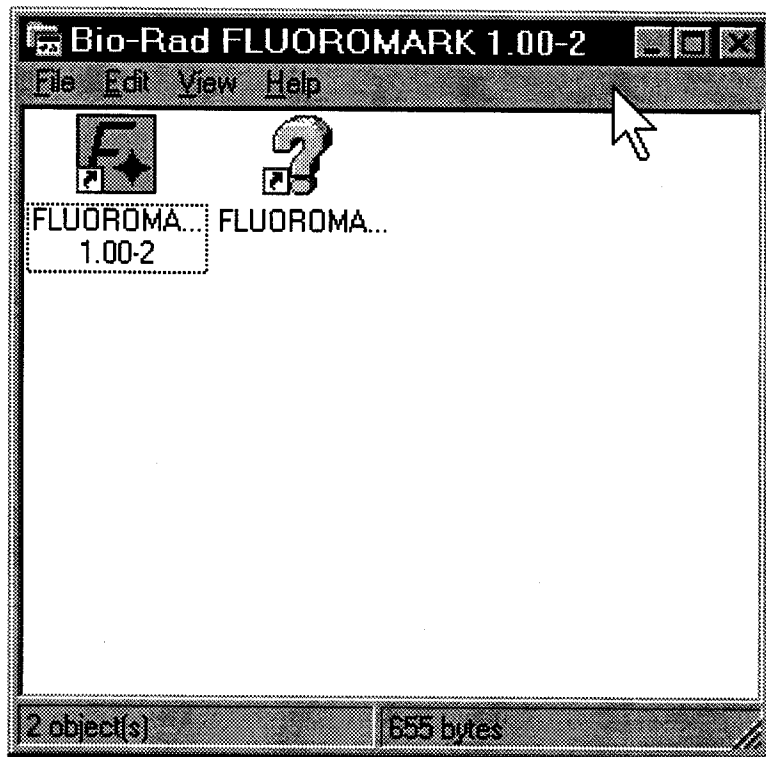


Fig. 4.8. Screen to start Fluoromark software.

You are now ready to begin using the Fluoromark package.

## Section 5

### Using the Fluoromark Windows PC Software

#### 5.1 Hints for Operation

The Fluoromark software runs under Microsoft Windows version 3.1, 3.11 (Workgroups), or '95. While any computer capable of running Windows will run the Fluoromark software, a Pentium processor with a minimum of 16 Mbytes of RAM is recommended.

The red transport locking pin must be removed to free up the plate carrier after the instrument is installed.

The Offset values from the back of the instrument must be entered on the Instrument Setup, Offset, screen before a measurement can be made.

All screen savers must be disabled while using the Fluoromark PC program, as the system can lose data across the serial line from Fluoromark to the PC when screen savers are in effect.

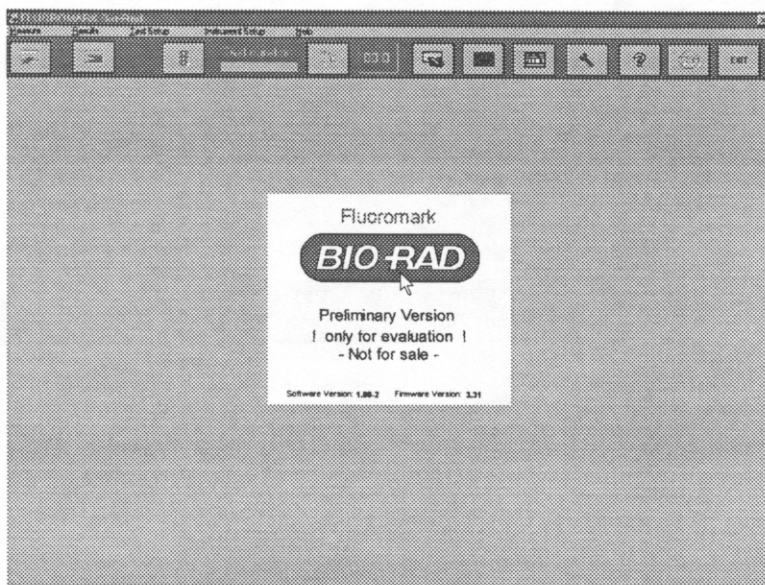


Fig. 5.1. Startup screen with menus and symbols.

#### 5.2. General View of the Menus and the Corresponding Icons

Frequently used menu selections can be accessed directly from the Tool Bar. Tool Bar icons perform the same functions as the menu items listed below.

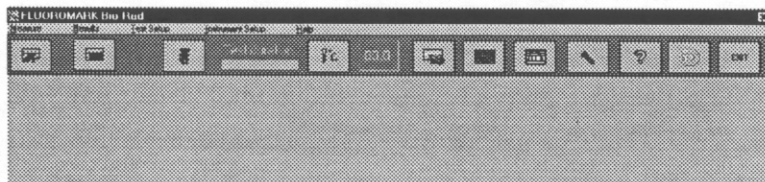
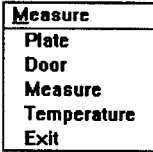









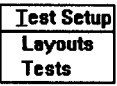


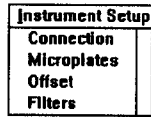

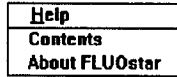



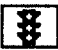


Fig. 5.2. Menus and symbols.

Icon	Description	Action
	Measure	Menu for Fluoromark hardware control and for starting test runs. The Fluoromark microplate fluorometer has to be switched on and connected to the PC for this menu to be active.
	Plate out	Opens the Fluoromark microplate fluorometer door and ejects plate carrier. If the plate carrier is already out, no action is taken.
	Plate in	Brings the plate carrier in. If the plate carrier is already in, no action is taken.
	Door open	Open the door
	Door close	Close the door
	Measure	Performs a measurement on a microplate, using a pre-defined test procedure
	Real-time graphic	During a test run the measure icon will change to the real-time graphics icon. This opens the real-time graphic screen.
	Temperature	Sets the target temperature of the incubator 25 °C to 45 °C in 0.1 °C steps, and begins incubating
	Results	Transfers control to Excel data reduction template for measurement results
	Results	Transfers control to Excel data reduction template for measurement results
	Test setup	Menu to define microplate sample positions and test protocols
	Layouts	Menu to define the positions of Standards, Blanks, Unknowns, Empty wells, and concentrations
	Tests	Menu to define measurement timing, shaking, and filters for excitation and emission

Icon	Description	Action
	Instrument setup	Defines hardware configurations and plate types
	Connection	Menu to define com port (1 to 4) for connecting Fluoromark microplate fluorometer to PC
	Microplates	Menu to define the dimensions of microplate
	Offset	Menu to set the "home position" for each Fluoromark instrument
	Filters	Menu to define excitation and emission filters
	Help	Menu for the integrated help function of the program with information about the version number of the software
	About Fluoromark	Software version, version number software of the PC software. Firmware version, version number of the Fluoromark software.
	Stop button	Stops an active test run, saving any data collected
	End of program	Ends the program, sends init command to the Fluoromark fluorometer

## 5.3 Performing a Measurement

Test procedures are performed by selecting the  symbol or by selecting Measure in the measurement menu. After initiating the test run, a message is displayed prompting for insertion of a microplate.

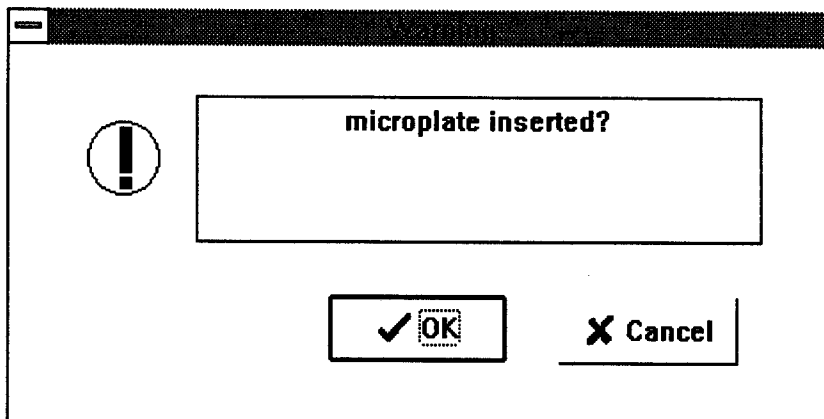


Fig. 5.3. Warning before starting the test run.

The warning screen (Figure 5.3) prompts for insertion of a microplate before a test run.

### Select buttons

**OK** proceeds to the next menu for selecting a pre-defined test

**Cancel** aborts measurement, returns to main menu

After pressing OK button, all pre-defined tests will be listed.

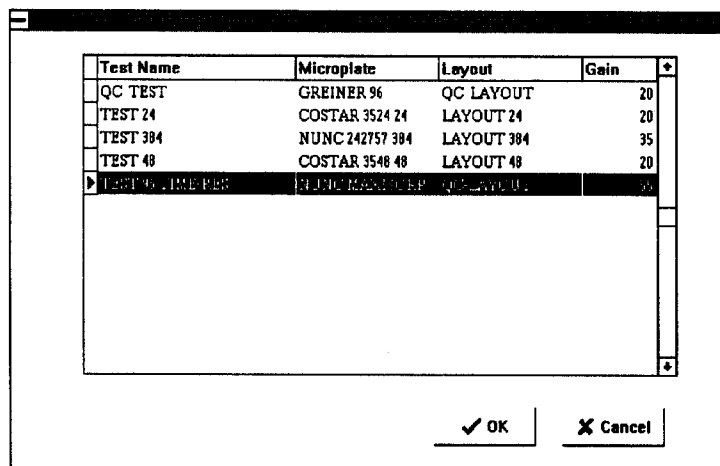


Fig. 5.4. Listing of all pre-defined tests for selection.

The test screen (Figure 5.4) lists all pre-defined tests, showing the following information.

**Test Name** previously defined test procedures

**Microplate** microplate type for the corresponding test

**Layout** layout name for the corresponding test

**Gain** defined sensitivity (gain value) for the corresponding test

Select a test procedure with a mouse click or the cursor keys, and start the selected test by pressing the OK button.

**Select buttons**

**OK** proceeds to the next screen, for selection of the plate ID

**Cancel** aborts measurement, returns to main menu

After specifying a test procedure, an alpha numeric identifier for the experiment may be entered. This name will be saved with the results of the measurement, and may be used to recall, calculate, and export the data.

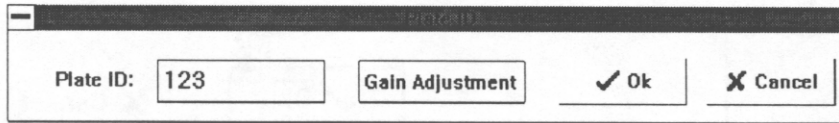


Fig. 5.5. Enter an identification number before test start.

The enter an identifier for the test run screen (Figure 5.5) lets you enter an identification number. The name or number will not be used in calculations.

**Select buttons**

**Gain Adjustment** proceeds to the next screen, see Figure 5.8, for running Gain adjustment

**OK** begin measuring the selected test

**Cancel** aborts measurement, returns to main menu

After starting a test, a time gauge will be displayed in the tool bar, indicating the elapsed time and time left for completion of the measurement. This gauge is updated dynamically during the test run. If temperature control has been selected, the actual temperature will also be displayed.

The measure icon will change to the real-time graphics icon in the Tool Bar.



Fig. 5.6. Menu and symbols during test run.

After selecting the real-time graphics icon, the following screen will appear.

96	1	2	3	4	5	6	7	8	9	10	11	12
A												
B												
C												
D												
E												
F												
G												
H												

Max:     
 Min:

Cycle:     
 Intervals:

Auto Scale  
 Manual Scale

Fig. 5.7. Screen for real-time graphics.

The real-time graphics screen (Figure 5.7) displays real time graphic in defined microplate format (384, 96, 48, 24). During the test run all measured values will be automatically displayed as points. If you select Auto Scale minimum and maximum limits will be updated automatically. If you select Manual Scale minimum and maximum limits can be specified manually. You can specify the number of points to be shown in a well by manipulating the number of intervals.

**Select buttons**

**Max** maximum limit for displayed values

**Min** minimum limit for displayed values

**Intervals** number of points to be shown. Default value is number of cycles defined in test definition.

**OK** closes real-time graphics screen and saves measured values if test run is finished

Selected Test: QC TEST      Layout: QC LAYOUT

96	1	2	3	4	5	6	7	8	9	10	11	12
A												
B	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1
C	S2											S2
D	S3											S3
E	S4											S4
F	S5											S5
G	S6											S6
H	B											B

Required Value:

Gain:

Fig. 5.8. Screen for Gain Adjustment.

The gain adjustment screen (Figure 5.8) performs gain adjustment for the plate. Specify the required value up to 9999. (Bio-Rad recommends using 9500). The gain adjustment button will determine the optimum gain for the selected well (for example the highest known standard).

**Select buttons**

**Required value** required value of the result in relative fluorescent units

**Gain optimum** gain for required result


**Gain Adjustment** performs gain adjustment for the selected well in the layout above

**Adjustment**

**OK** closes gain adjustment screen and saves gain value in corresponding test definition

**Cancel** aborts gain adjustment

### 5.4 Incubating Plates (optional equipment)

Incubation (if installed) is implemented with either the  icon, or by selecting the temperature option in the measurement menu.

After selecting this function the following screen will appear.

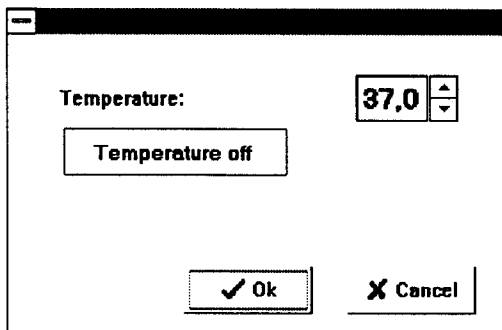


Fig. 5.9. Screen temperature selection.

Specify the target temperature of the incubator (25 °C to 45 °C), in steps of 0.1 °C using either the spinner or by typing directly into the selection box. The Temperature On button toggles to Temperature Off during incubation.

**Select buttons**

**Temperature** defines the target incubation temperature


**Temperature off** de-activates incubation

**OK** closes temperature screen

**Cancel** aborts last action, returns to main menu

If you have selected incubation, the measured temperature will be displayed in the tool bar. If the temperature is below the target temperature, the field will be red. If the target temperature has been reached, the field will be green.

## 5.5 Defining a Layout

Microplate layouts are defined using the  icon, or by selecting Layouts in the test setup menu. After selecting this function, all previously defined layouts will be listed.

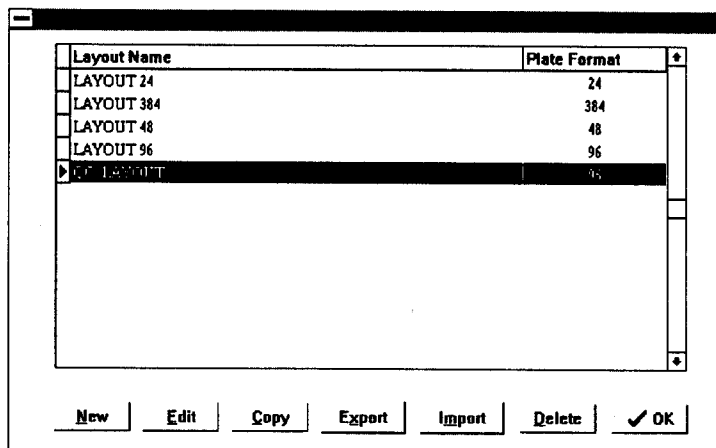


Fig. 5.10. List of all defined layouts.

The layouts screen (Figure 5.10) lists all previously defined and stored layouts.

### Select buttons

- New** creates a new microplate layout
- Edit** modifies an existing microplate layout
- Copy** copies an existing microplate layout
- Export** exports an existing microplate layout
- Import** imports an existing microplate layout
- Delete** deletes an existing microplate layout
- OK** closes the layout screen, and returns to main menu

When a new layout is defined, the plate format (384, 96, 48, or 24 wells) must be defined first.

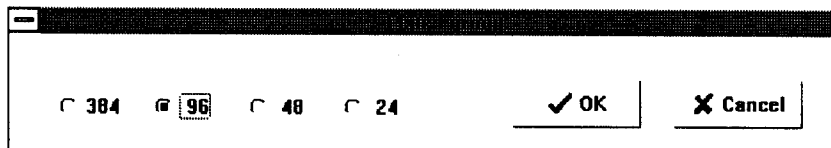


Fig. 5.11. Defining the plate format of a new layout.

The plate format screen (Figure 5.11) defines the plate format for the new microplate layout.

### Select buttons

- 384** selects plate format to 384 wells
- 96** selects plate format to 96 wells
- 48** selects plate format to 48 wells
- 24** selects plate format to 24 wells
- OK** begins definition of a new layout of the selected plate format
- Cancel** aborts definition, returns to layouts screen

After selecting the plate format for the new microplate layout, the software proceeds to the corresponding layout definition screen.

Layout Name:

Content:

24	1	2	3	4	5	6
A	B	X1		B	X5	X6
B	S1	S1	S2	S2		
C	S3	S3	S4	S4	X1	X2
D	B			B	X3	X4

Index:

Constant  
 Increase

Replicates:

Horizontal  
 Vertical

Fig. 5.12. Screen for modifying or defining a 24 well layout.

Layout Name:

Content:

48	1	2	3	4	5	6	7	8
A	B	B	B	B				
B	S1	S1	S2	S2				
C	S3	S3	S4	S4				
D	B	B	B	B				
E	X1	X2	X3	X4				
F	X5	X6	X7	X8				

Index:

Constant  
 Increase

Replicates:

Horizontal  
 Vertical

Fig. 5.13. Screen for modifying or defining a 48 well layout.

Layout Name:

Content:

Index:

Constant  
 Increase

Replicates:

Horizontal  
 Vertical

96	1	2	3	4	5	6	7	8	9	10	11	12
A												
B	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1	S1
C	S2											S2
D	S3											S3
E	S4											S4
F	S5											S5
G	S6											S6
H	B											B

Fig. 5.14. Screen for modifying or defining a 96 layout.

Layout Name:

Content:

Index:

Constant  
 Increase

Replicates:

Horizontal  
 Vertical

384	1	2	3	4	5	6	7	8	9	10	11	12
A	B	B	B	B	B	B	B	B	B	B	B	B
B	B	B	B	B	B	B	B	B	B	B	B	B
C	S1	S1	S1	S1	S2	S2	S2	S2	S3	S3	S3	S3
D	S7	S7	S7	S7	S8	S8	S8	S8	S9	S9	S9	S9
E	S13	S13	S13	S13	S14	S14	S14	S14	S15	S15	S15	S15
F	S19	S19	S19	S19	S20	S20	S20	S20	S21	S21	S21	S21
G	X1	X2	X3	X4	X5	X6	X7	X8	X9	X10	X11	X12
H	B	B	B	B	B	B	B	B	B	B	B	B

Fig. 5.15. Screen for modifying or defining a 384 layout.

The layout screens (Figure 5.13 through 5.15) are used to modify or to define a layout.

**Select buttons**

- Content** Allows selection of well type (Empty, Samples, Blanks, Standards)
- Index** Selects the number of the next well defined. (Define Auto Increment to automatically increase the index, or Constant Index for continuous replicates.)
- Replicates** Selects the default number of replicates. Horizontal or Vertical determines the order of the replicates.
- Concentrations** Switches to definition of standard concentrations (Figure 5.16)
- OK** Saves all modifications and returns to Figure 5.10 (layouts)
- Cancel** Returns to Figure 5.10 without saving

## How to Modify A Layout

Define the appropriate type of sample for a given group of wells in the Content drop-down box. After defining the well type, select the well locations using the left mouse button, and dragging across all wells of that type. Wells can also be defined with the keyboard, by holding the shift key and highlighting the appropriate wells with the cursor keys. Wells can be defined by row or column by clicking the row or column label (A to H and 1 to 12). All wells of the microplate can be selected by clicking the field in the upper left border.

The concentrations screen (Figure 5.16) shows concentrations for the defined layout. Concentrations can be defined for individual wells by typing directly or by using the auto function. To use the auto function, define start value (concentration) and select factor, increment, or decrement and type in the offset or factor. Then select concentration by using the right mouse button and dragging across the fields you want to change. Fields can also be changed with the keyboard, by holding the shift key and highlighting the appropriate fields with the cursor keys.

### Select buttons

**Concentration** defines the concentration of the available standards with the auto function

**Layout** saves all modifications and returns back to Figure 5.10

### Definition

**Cancel** returns to Figure 5.10 (layout definition) without saving

Layout Name: QC LAYOUT

Concentration

Factor

Increment

Decrement

Content	Concentration
S1	3000
S2	2000
S3	1000
S4	500
S5	250
S6	125
B	

Fig. 5.16. Screen for defining concentrations.

The copy screen (Figure 5.17) is used to copy layouts, tests, or microplates

### Select buttons

**OK** begin copying the selected layout, test or microplate

**Cancel** aborts copy function

Copy

To

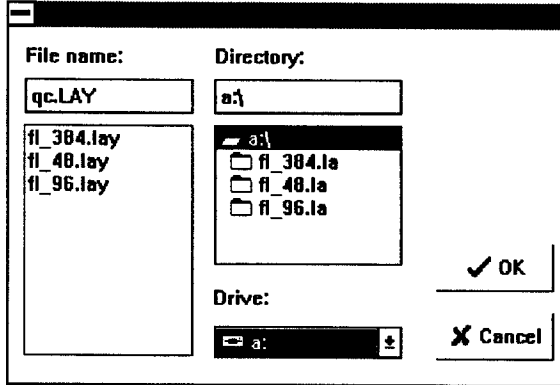
Fig. 5.17. Screen for copying layouts, tests or microplates.

The export screen (Figure 5.18) is to export layouts, tests or microplates.

**Select buttons**

**OK** begin exporting the selected layout, test or microplate

**Cancel** aborts export function



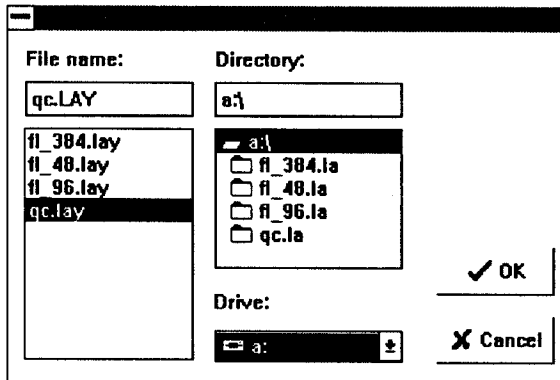
**Fig. 5.18. Screen for exporting layouts, tests, or microplates.**

The import screen (Figure 5.19) is to import layouts, tests, or microplates

**Select buttons**

**OK** begin importing the selected layout, test or microplate

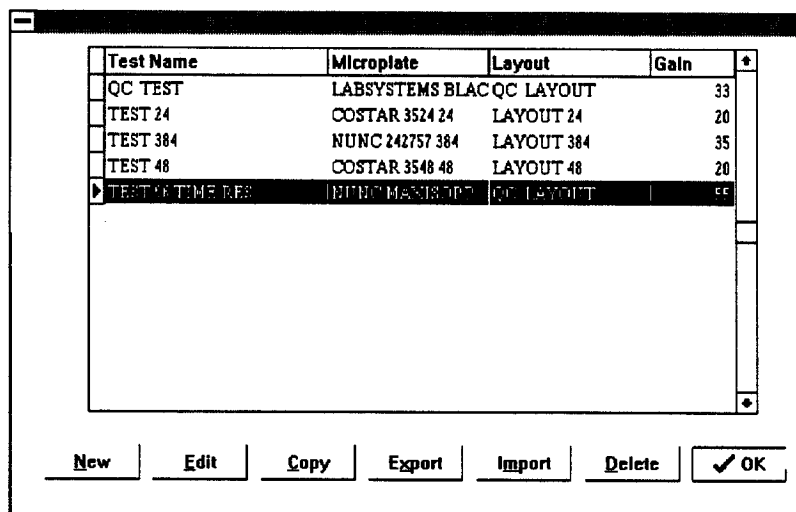
**Cancel** aborts import function



**Fig. 5.19. Screen for importing layouts, tests, or microplates.**

## 5.6 Defining a Test

Test procedures are defined using the  icon on the tool bar, or by selecting tests from the test setup menu. After selecting this function all defined tests will be listed.



Test Name	Microplate	Layout	Gain	+
QC TEST	LABSYSTEMS BLAC	QC LAYOUT	33	
TEST 24	COSTAR 3524 24	LAYOUT 24	20	
TEST 384	NUNC 242757 384	LAYOUT 384	35	
TEST 48	COSTAR 3548 48	LAYOUT 48	20	
TEST 96 TIME RES	NUNC MANIP 96	QC LAYOUT	66	

Buttons: New, Edit, Copy, Export, Import, Delete, OK

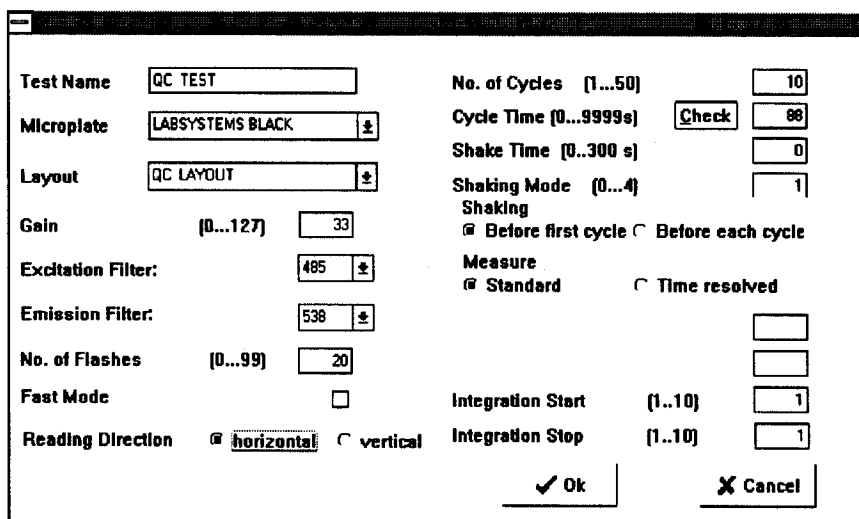
Fig. 5.20. List of all defined tests.

The tests screen (Figure 5.20) gives a list of all pre-defined tests.

### Select buttons

- New** defines a new test (Figure 5.21, test definition)
- Edit** modifies a previously defined test (Figure 5.21, test definition)
- Copy** copies an existing test (Figure 5.17, layout)
- Export** exports an existing test (Figure 5.18, layout)
- Import** imports an existing test (Figure 5.19, layout)
- Delete** deletes an existing test
- OK** returns to main menu, saving all modifications

When a new test procedure is defined, the software proceeds to the test definition screen to define the measurement procedure.



Test Name: QC TEST

Microplate: LABSYSTEMS BLACK

Layout: QC LAYOUT

Gain: [0...127] 33

Excitation Filter: 485

Emission Filter: 538

No. of Flashes: [0...99] 20

Fast Mode:

Reading Direction:  horizontal  vertical

No. of Cycles [1...50]: 10

Cycle Time [0...9999s]:  88

Shake Time [0..300 s]: 0

Shaking Mode [0...4]: 1

Shaking:  Before first cycle  Before each cycle

Measure:  Standard  Time resolved

Integration Start [1..10]: 1

Integration Stop [1..10]: 1

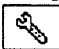
Buttons: Ok, Cancel

Fig. 5.21. Screen to define test.

<b>Test screen (Figure 5.20)</b>	to define a test
<b>Test name</b>	alpha numeric identifier for the test procedure
<b>Microplate</b>	selects the appropriate microplate type from a pre-defined list (default plates are included, and custom plates may also be defined in setup)
<b>Layout</b>	selects a pre-defined template, which designates the positions of Samples, Standards, etc. The Layout is used both for data reduction and to determine which wells should be measured.
<b>Gain</b>	adjusts voltage to the PMT, to provide optimum sensitivity for each assay
<b>Excitation filter</b>	selects the appropriate excitation filter type from a pre-defined list (defined in setup)
<b>Emission filter</b>	selects the appropriate emission filter type from a pre-defined list (defined in setup)
<b>Fast mode</b>	for small kinetic test runs you can select fast mode operation. Normally after inserting your microplate the transport system will be calibrated before measurement will begin. In fast mode no calibration will be done.
<b>Reading direction</b>	selects reading in horizontal or vertical direction
<b>No. of cycles</b>	defines the number of times the entire plate is measured, from 1 to 50 cycles; for every cycle, each well defined in the layout will be measured
<b>Cycle time</b>	defines the time interval between the start of one plate measurement to the start of the next cycle; this time includes the time to position the plate carrier, shake the plate (if used), the measurement time for each well, and a waiting time if necessary. If a Fluoromark fluorometer is connected, the check cycle time button is activated to validate the Cycle Time directly from the Fluoromark fluorometer.
<b>Shake time</b>	defines length of time to shake the plate before a cycle, from 0 to 300 seconds
<b>Shaking Mode</b>	defines mode for shaking. Mode 0 means high speed and low diameter (96 well plate) and mode 4 low speed and high diameter (24 well plate)
<b>Shaking before first cycle / Before each cycle</b>	selects if shaking is done only before the first or before each cycle
<b>Measure standard/ Time resolved</b>	selects mode for measure. Standard is done with fixed values for integration delay and integration time. For Time Resolved the values for integration delay and integration time can be defined in a range from 10 $\mu$ s to 1,534 $\mu$ s.
<b>Integration start/ stop</b>	in addition to using a measurement point from a single cycle for data analysis, the Excel template also has the ability to handle data from measurement points from several cycles; Integration Start/Stop selects the measurement points that will be used for data reduction (these integration limits can also be changed in Excel)
<b>Select buttons</b>	
<b>OK</b>	returns to test screen (Figure 5.20) saving all changes to the test definition
<b>Cancel</b>	returns to test screen (Figure 5.20) without saving any changes to the test definition

## Section 6

### Configuration of the Fluoromark and Definition of Microplates

Configuration of the communication port, Offset, definition of microplates, and configuration of the filters is initiated with the  symbol, or by selecting the Instrument Setup menu.

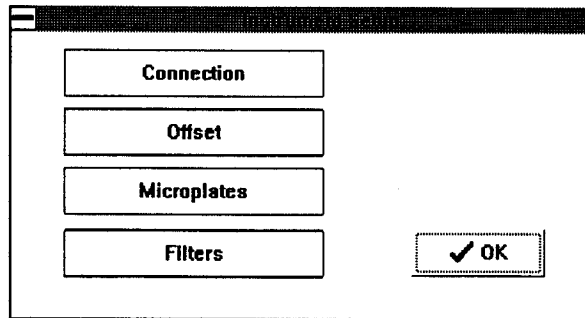


Fig. 6.1. Screen for instrument settings.

#### 6.1 Serial Port For Communication PC to Fluoromark Microplate Fluorometer

To change the PC serial communication port, select Connection in the instrument setup menu.

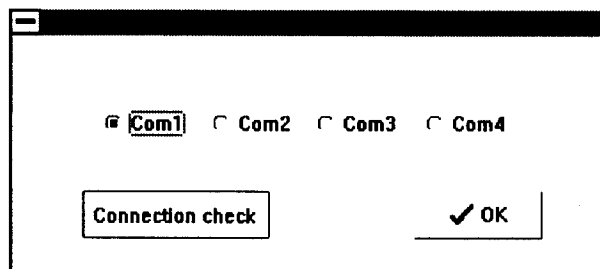


Fig. 6.2. Screen for selecting the com port.

Screen to select com port for the communication between PC and Fluoromark.

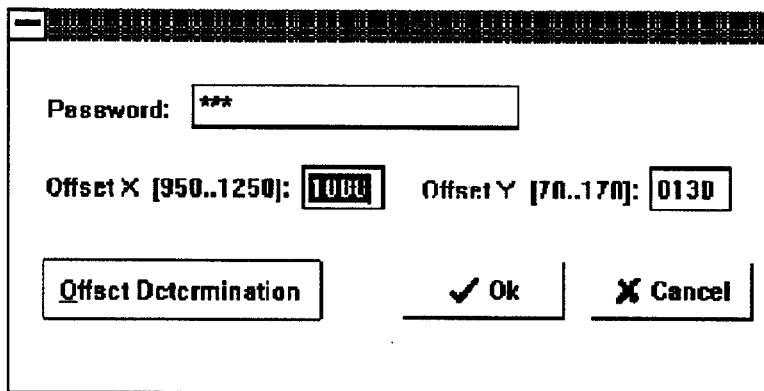
#### Select buttons

<b>Com1</b>	to select com port 1
<b>Com2</b>	to select com port 2
<b>Com3</b>	to select com port 3
<b>Com4</b>	to select com port 4
<b>OK</b>	return to main menu, saving changes
<b>Connection Check</b>	check the selected com port to verify the presence of a Fluoromark instrument

## 6.2 Definition of the Mechanical Offset

Every Fluoromark instrument is calibrated with a different offset position. This value is normally determined at the factory and is not changed. If service is performed to the plate transport or to the measuring head, the new offset position must be determined by a service technician. The calibrated offsets are posted on the back of each instrument.

Select Offset in the instrument setup menu to enter the offsets.



The screenshot shows a graphical user interface for setting mechanical offsets. At the top, there is a password field labeled 'Password:' containing three asterisks. Below this are two input fields: 'Offset X [950..1250]:' with the value '1000' and 'Offset Y [70..170]:' with the value '0130'. At the bottom, there are three buttons: 'Offset Determination', 'Ok', and 'Cancel'.

Fig. 6.3. Screen to define the offset.

Valid Offset calibration values are vital to optimum performance. Optical crosstalk and variability of results, can occur if incorrect offset values are used. Type Bio-Rad as the password, and enter the X and Y offset values from the back of the instrument.

If service has been performed, or if the offset values are not available, instruments can be field calibrated. To determine new offset values, fix the plate with the red locking pin. Then remove the pin without changing the position of the plate transport, and press Offset Determination button. The plate transport will move to the home position, and the new calibrated Offset values will be placed in the appropriate fields.

Password requires you to enter the password BMG. When the password is entered, the offset values can be changed, or automatically determined

**Offset X** field to enter the offset in x direction

**Offset Y** field to enter the offset in y direction

### Select buttons

**Offset Determination** starts the automatic offset determination. Before you can start this, you have to type in the correct password. Then you can position the plate carrier under the measure head and check the right position with the stick. After this take out the stick and start the offset determination function.

**OK** saves the offset values, back to main menu

### 6.3 Definition of the Mechanical Dimensions of the Microplates

The Fluoromark software is supplied with the dimensions of standard microplates from all major microplate manufacturers. While most plates have the standard spacing and footprint, some plates have slightly different dimensions, and must be positioned accordingly for optimum results. In addition, new plates can easily be defined by selecting Microplates in the menu instrument setup.

Microplate	Length	Width	X (1)	Y (1)	X (n)	Y (n)	Plate Format
CORNING 25860 96	1276	855	144	112	1132	742	96
CORNING 25880 96	1276	855	144	112	1132	742	96
COSTAR 3524 24	1278	853	175	140	1140	715	24
COSTAR 3548 48	1277	854	180	99	1097	755	48
COSTAR 96	1280	856	145	113	1138	744	96
DYNATECH 96	1280	858	144	112	1136	741	96
FALCON 24	1275	854	140	138	1104	717	24
FALCON 3072 96	1276	858	143	114	1134	744	96
FALCON 3078 48	1278	856	187	105	1092	752	48
GREINER 24	1280	858	165	144	1115	714	24
GREINER 96	1280	855	144	112	1136	743	96
LINBRO 24	1275	860	109	110	1176	750	24
NEW 124256 96	1276	855	144	112	1132	742	96

Fig. 6.4. List of defined microplates.

**Microplate screen (Figure 6.4)** lists all defined microplates with their specific mechanical dimensions.

#### Select buttons

- New** proceeds to microplate screen (Figure 6.5) to create a new microplate name with no pre-defined dimensions
- Edit** proceeds microplate screen (Figure 6.5) to change a defined microplate
- Copy** copies an existing microplate layout screen (Figure 5.17)
- Export** exports an existing microplate layout screen (Figure 5.18)
- Import** imports an existing microplate layout screen (Figure 5.19)
- Delete** deletes the currently selected microplate definition
- OK** returns to the main menu, saving changes

Select New or Edit to proceed to microplate screen (Figure 6.3) for definition of the mechanical dimensions of a microplate, in steps of 0.1 mm.

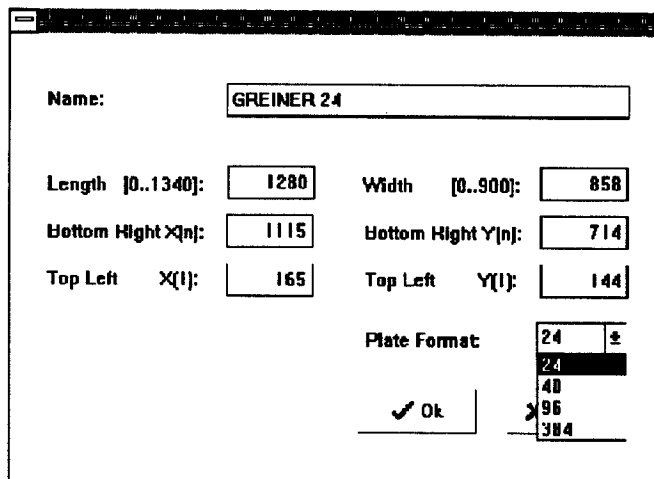


Fig. 6.5. Screen to define the mechanical dimensions of microplates.

Microplate screen (Figure 6.5) defines the plate format (384, 96, 48, or 24 wells) and the mechanical dimensions of microplates, in steps of 0.1 mm.

<b>Name</b>	name for the definition of the microplate
<b>Length</b>	outer length of the entire microplate, in the 1–12 dimension, in steps of 0.1 mm
<b>Width</b>	outer width of the entire microplate, in the A–H dimension, in steps of 0.1 mm
<b>X(n)</b>	distance from the center of the lower right well (normally H12) to the left outer edge of the microplate, in steps of 0.1 mm
<b>Y(n)</b>	distance from the center of the lower right well (normally H12) to the top outer edge of the microplate, in steps of 0.1 mm
<b>X(1)</b>	distance from the center of the upper left well (normally A1), to the left outer edge of the microplate, in steps of 0.1 mm
<b>Y(1)</b>	distance from the center of the upper left well (normally A1) to the top outer edge of the microplate, in steps of 0.1 mm

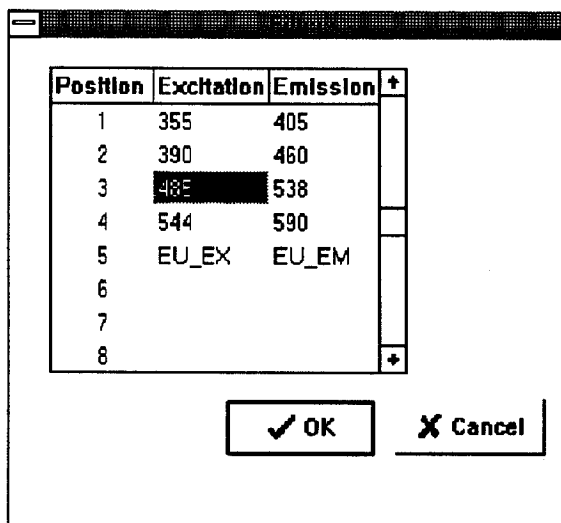
**Plate format** selects the plate format of the Microplate (384, 96, 48 or 24 wells)

**Select buttons**

<b>OK</b>	accepts current microplate name and dimensions, and returns to microplates screen.
<b>Cancel</b>	returns to microplates screen without saving changes

## 6.4 Definition of the Filter Configuration

The Fluoromark can be configured with 8 excitation and 8 emission filters. To define the wavelength and position of the filters in your Fluoromark select Filters in the instrument setup menu.



Position	Excitation	Emission	↑
1	355	405	
2	390	460	
3	485	538	
4	544	590	
5	EU_EX	EU_EM	
6			
7			
8			+

✓ OK      ✗ Cancel

Fig. 6.6. Screen to define the filter configuration.

### Select buttons

**OK**            saves all positions and wavelength of the filters, and returns to main menu

**Cancel**        returns to main menu without saving changes


## Section 7

### Microsoft Excel Data Reduction Routines

**Note:** Data Access must be installed with Excel, in order for the database functions to be enabled.

#### 7.1 General Hints for Operation

Although no software can incorporate every conceivable feature, modern spreadsheet packages provide powerful data reduction with virtually unlimited flexibility. Pre-written data reduction templates in Microsoft's Excel display raw data, plot flash curves for single or multiple wells, calculate the average of replicates, perform blanking, calculate the concentrations of unknown samples, and graph standard curves.

Open Excel by pressing the  icon from the tool bar. Starting Excel this way, rather than by using Windows Program Manager, will automatically load the Fluoromark.XLT template. When the template is loaded, the name is changed to Fluoromark1.XLS to prevent accidentally overwriting the original template. (**Note:** the XLT extension denotes a template, and the XLS extension indicates a spreadsheet.)

Fluoromark1.XLS is actually several worksheets bound together in a workbook. The data from each experiment is divided into separate groups, and each sheet displays a different set of data. The name of each sheet appears on the tab at the bottom of the sheet, and users activate individual sheets by selecting the appropriate tab with the mouse.

The first sheet, called testruns, is a database of all previously measured plates, which lists the name, plate type, layout, and other parameters of each experiment. Data is recalled from the computer's hard drive by double-clicking anywhere on the line corresponding to the measurement of choice. Sheets for displaying raw data, graphs, and final data can be selected in any order.

An additional pull down menu Fluoromark in the Excel menu line allows tests to be imported into the database, or exported from the database to a separate file, add comment information to the test run, copy or delete a test run and a menu option switch back to Fluoromark software.

#### 7.2 Performing Data Reduction in Excel

When the Testruns sheet is selected, a database of all previously measured plates is displayed.

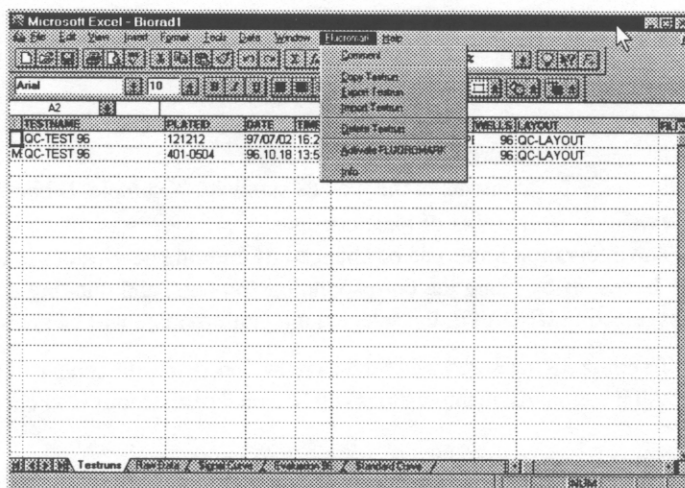


Fig. 7.1. Sheet 1 (testruns) in Excel.

Testruns screen (Figure 7.1) gives a list of all test runs with following information

<b>Test Name</b>	name of the corresponding test setup
<b>Plate ID</b>	measurement identification name of number
<b>Date</b>	date of the test run
<b>Time</b>	time of the test run
<b>Microplate</b>	name of the corresponding microplate definition
<b>Wells</b>	plate format of the corresponding microplate
<b>Layout</b>	name of the corresponding layout definition
<b>Filename</b>	file name of the test results

Select a test run by double clicking on the line corresponding to the measurement of interest.

Test runs can be copied, exported, imported and deleted with the menu item Fluoromark, and selecting the appropriate menu option. Additionally it is possible to add a comment to every test run. With the menu option Activate Fluoromark, you can switch back to the Fluoromark program. After selecting a test run, the program will automatically switch to sheet 2 (raw data).

Well	Cont.	Cycles	1	2	3	4	5	6	7	8	9	10	11	12
B01	S1	2170	2	2	2	2	2	2	2	2	2	2	2	2
B02	S1	218	3	3	3	3	3	3	3	3	3	3	3	3
B03	S1	220	3	3	3	3	4	8	7	9	8			
B04	S1	222	2	2	2	2	3	7	8	10	7			
B05	S1	215	2	2	2	2	2	29	50	61	66			
B06	S1	215	2	2	2	2	8	24	52	52	60			
B07	S1	215	2	2	2	2	7	222	448	468	422			
B08	S1	217	2	2	2	2	51	279	481	491	460			
B09	S1	215	2	2	2	2	38	1894	2821	4916	4221			
B10	S1	218	2	2	2	2	637	1782	4060	4737	4605			
B11	S1	218	2	2	2	2	412	22970	43270	44060	41060			
B12	S1	218	2	2	2	2	3	19100	46430	11900	42150			
C01	S2	218	2	2	2	2	2	2	2	2	2			
C12	S2	215	3	3	3	3	3	3	3	3	3			
D01	S3	218	2	2	2	2	2	2	2	2	2			
D12	S3	215	3	3	3	3	3	3	3	3	3			
E01	S4	215												
E12	S4	217												
F01	S6	216												
F12	S6	218												
G01	S8	220												
G12	S8	214												
H01	S	215												
H12	S	215												

Fig. 7.2. List of all raw data for the selected test run AEQUORIN TEST.

Raw data screen (Figure 7.2) for the Fluoromark data analysis worksheet, displays the fluorescence signal data for each measured well. Relative fluorescence units for as many as 50 time intervals are shown after the well descriptions.

Each line gives the following information for each measured well.

<b>Well</b>	microplate well location
<b>Cont.</b>	contents of the well (S standard, X sample, B blank)
<b>Interval</b>	number of plate cycle

The integration limits can be changed by entering new values in the integration Start and Stop boxes, or by using the corresponding spinners, and clicking on the UPDATE button. After selecting UPDATE, the selected integration range will be shown in red.

To remove a measured well for data reduction, delete the content of the well and click on the UPDATE button. All manipulations will be stored permanently after clicking on the Save button.

**Select button**

**UPDATE** updates and saves your changes on integration start and stop and the selected wells for the following sheets

**Save** saves all manipulations in the test run file

To view a plot of the fluorescence curve of a single well or group of wells, highlight the well(s) of interest, select UPDATE, and switch to the Signal Curve sheet by clicking on the corresponding tab at the bottom of the sheet. A range of wells can be selected by clicking and dragging, and individual wells can be chosen by clicking on individual wells with the CTRL key depressed.

Signal curve screen (Figure 7.3) selects the signal curve of the selected wells from sheet 2. All formatting and mathematical functions of Excel are enabled, and can be used without restrictions.

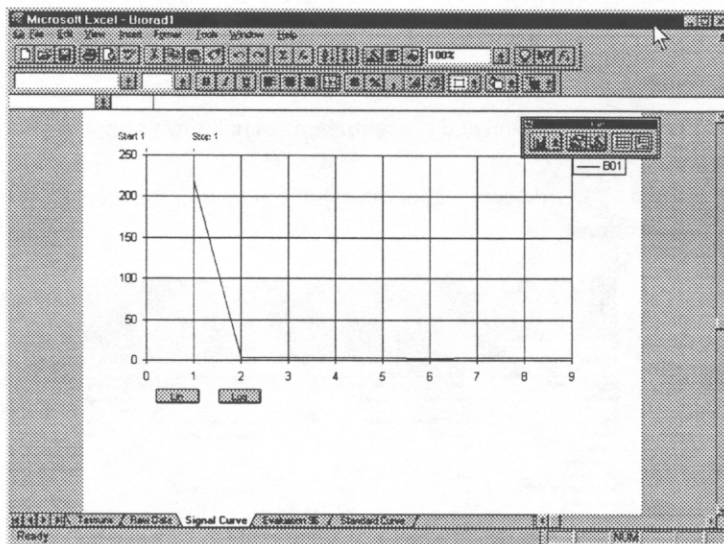


Fig. 7.3. Linear display of the signal curve.

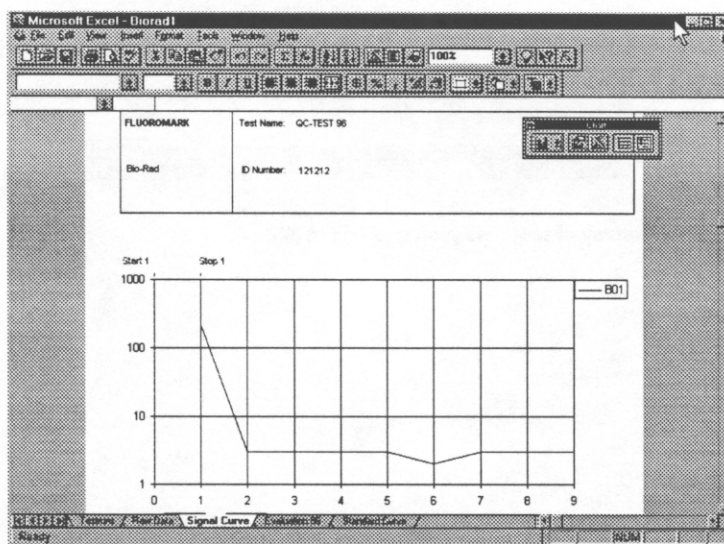


Fig. 7.4. Logarithmic display of the signal curve.

### Switch buttons

- **Lin** changes Y axis (RFU) of signal curve to linear
- **Log** changes Y axis (RFU) of signal curve to logarithmic

Sheet 4, evaluation384, evaluation96, evaluation48, evaluation24, displays test definitions and calculated data. For displayed data, three tables in the microplate format are shown. The type of data (Raw, Blanked, Concentrations, etc.), can be selected using the drop-down menu on each table. The format of this sheet depends on the microplate used for the selected test run.

### Select button

- Layout** displays the microplate layout of the selected plate
- Raw results** integral of the raw data from start to stop interval
- Averages** displays the average integrals of replicate wells
- Blank** displays averages the Blank values subtracted
- Correction (RLU)**
- Unknown concentrations** calculated concentrations of unknown samples
- Standard concentrations** theoretical concentrations of known standards

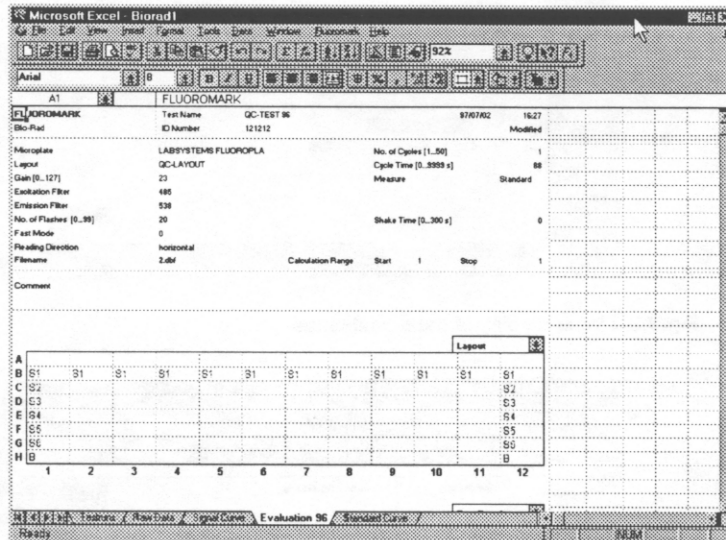


Fig. 7.5. Display of test definition and layout (96).

Sheet 5, standard curve, plots the measured RFU values for the standard concentrations, defined in the Test Setup, as a Standard Curve, or Calibration Curve. The Standard Curve can be displayed by selecting the Standard Curve tab at the bottom of the sheet. The Curve can be shown in linear or logarithmic scale.

**Select buttons**

**Lin** changes Y axis (RFU) and X axis (Conc.) to linear

**Log** correlation curve in logarithmic display

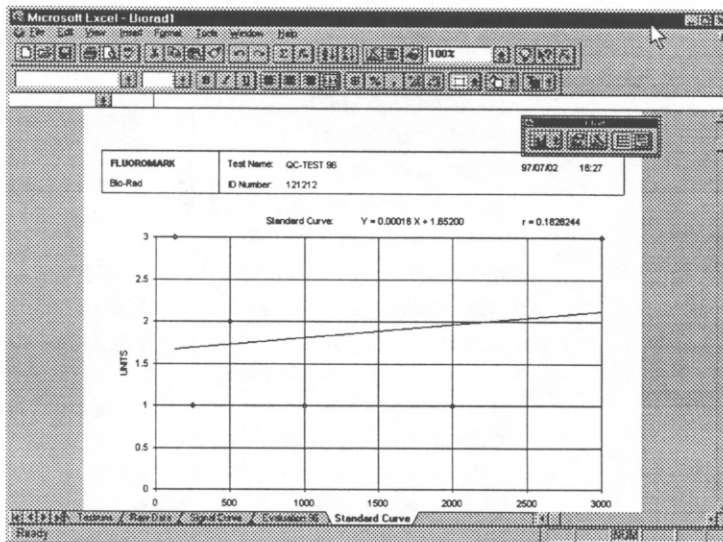


Fig. 7.6. Correlation curve in linear display.

## Section 8 Error Codes

<b>Error Code</b>	<b>Error Type</b>	<b>Error Message</b>
0100	wrong command format	Interface error
0200	command character not OK	Invalid command
0300	command parameter not OK	Invalid parameter
0501	Error during writing in Fluo.init	
0502	no temperature defined	
0503	offset X not defined	
0504	offset Y not defined	
0505	no com port selected	
0506	no microplate def. for this test	
0507	no layout definition for this test	
0811	calibration error X-motor	initialization error transp. system
0812	calibration error Y-motor	initialization error transp. system
0813	calibration error X- and Y-motor; home sensor defect	initialization error transp. system
0814	calibration error ex. filter wheel	Initialization error ex. filter wheel
0815	calibration error em. filter wheel	Initialization error em filter wheel
0821	right XY area limit error	transport error
0822	left XY area limit error	transport error
0823	forward XY area limit error	transport error
0824	back XY area limit error	transport error
0851	error temperature sensor 1	signal temperature sensor 1 missing
0852	error temperature sensor 2	signal temperature sensor 2 missing
0853	temperature 1 higher than maximum	temperature 1 higher than maximum
0854	temperature 2 higher than maximum	temperature 2 higher than maximum

## Section 9 Product Information

### 9.1 Product Information

Catalog Number	Product Description
170-6941	<b>Fluoromark Microplate Fluorometer, 110 V</b> , includes software, power cord, instruction manual, 5 excitation filters (355,390, 485, 544 nm, and time-resolved filter*), 5 emission filters (405, 460, 538, 590 nm, and time-resolved filter*)
170-6966	<b>Fluoromark Fluorometer, 100 V</b> (see description of 170-6941)
170-6970	<b>Fluoromark Fluorometer, 220 V EC version</b> (see description of 170-6941)
170-6942	<b>Fluoromark Microplate Fluorometer with built-in incubator</b> , includes the same accessory items as above.
170-6967	<b>Fluoromark Fluorometer with built-in incubator, 100 V</b> (see descriptions above)
170-6971	<b>Fluoromark Fluorometer with built-in incubator 220 V EC version</b> (see descriptions above)
170-6943	<b>Fluoromark Microplate Fluorometer with low UV read option down to 250 nm</b> , includes the same accessory items as above.
170-6968	<b>Fluoromark Fluorometer 220 V with low UV</b> , Japanese version (see descriptions above)
170-6972	<b>Fluoromark Fluorometer 220 V with low UV</b> , EC version (see descriptions above)
170-6944	<b>Fluoromark Microplate Fluorometer with built-in incubator and low UV read option down to 250 nm</b> , includes the same accessory items as above.
170-6969	<b>Fluoromark Fluorometer with built-in incubator and low UV</b> , 220 V Japanese version (see descriptions above)
170-6973	<b>Fluoromark Fluorometer with built-in incubator and low UV</b> , 220 V EC version (see descriptions above)
170-6946	<b>Microplate Manager Fluoromark Software with manual</b> , includes EXCEL version 5.0c
170-6947	<b>355 nm Excitation Filter</b>
170-6948	<b>390 nm Excitation Filter</b>
170-6949	<b>485 nm Excitation Filter</b>
170-6950	<b>544 nm Excitation Filter</b>
170-6951	<b>Time-Resolved Excitation Filter</b>
170-6958	<b>320 nm Excitation Filter</b>
170-6959	<b>515 nm Excitation Filter</b>
170-6960	<b>584 nm Excitation Filter</b>
170-6952	<b>405 nm Emission Filter</b>
170-6953	<b>460 nm Emission Filter</b>
170-6954	<b>538 nm Emission Filter</b>
170-6955	<b>590 nm Emission Filter</b>
170-6956	<b>Time-Resolved Emission Filter Eu</b>
170-6957	<b>Time-Resolved Emission Filter Sm</b>
170-6961	<b>555 nm Emission Filter</b>
170-6962	<b>612 nm Emission Filter</b>
170-6963	<b>96 Well Fluorescence Microplate</b> , solid black, 25
170-6964	<b>96 Well Fluorescence Microplate</b> , black with clear bottom, 25
170-6965	<b>384 Well Fluorescence Microplate</b> , solid black, 25

\* Standard time-resolved excitation and emission filters support use of Europium and Samarium fluorophores.

## 9.2 Filter Table

<b>Excitation</b>	<b>Emission</b>
355 nm (standard)	405 nm (standard)
390 nm (standard)	460 nm (standard)
485 nm (standard)	538 nm (standard)
544 nm (standard)	590 nm (standard)
Time-resolved filter (standard)	555 nm (optional)
320 nm (optional)	612 nm (optional)
515 nm (optional)	Europium time-resolved filter (optional)
	Samarium time-resolved filter (optional)

Bio-Rad also offers filters for other custom applications. Contact your sales representative or Bio-Rad Laboratories directly at 1 (800) 4BIORAD.

## 9.3 Disinfection Certificate

This instrument and its inventory has 1) Never been in contact with any dangerous biological material, or 2) been disinfected according to the instructions of the operating manual of this instrument.

Name: \_\_\_\_\_  
(Print) \_\_\_\_\_

Firm: \_\_\_\_\_  
(Print) \_\_\_\_\_  
\_\_\_\_\_  
\_\_\_\_\_

Signature: \_\_\_\_\_ Date: \_\_\_\_\_

Windows and Excel are trademarks of Microsoft.  
Cy is a trademark of Amersham.

**BIO-RAD**

## Declaration of Conformity

Bio-Rad Laboratories, Inc., 1000 Alfred Nobel Drive, Hercules, California, 94547, U.S.A., declares that the product  
Fluoromark Microplate Reader (Catalog nos. 170-6941, 170-6942, 170-6943, 170-6944, 170-6970, 170-6971, 170-6972, 170-6973)

---

to which this declaration relates, is in conformity to the following standards or normative documents

EN61010 Part 1/03.94, IEC 1010-1/09.90 & Amendment 1/09.92, EU Low Voltage (73/23EWG), EN 60555, EN 55011, EN 50082-1, VDE Cert. No. 90337 G, checked for conformity with DIN EN 61010 Part 1/03.94, Classification VDE 0411 Part 1/03.94, UL Listing No. 171095-E168510, and with UL 3101-1, CAN/CSA C22.2 No. 1010-1

---

following the provisions of the Low Voltage, EMC Directive.

The product is imported into the EU by Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead (London area), Hertfordshire HP2 7TD England.

July 30, 1997

date of issue

---

Norman Schwartz  
Vice President

DC1706973 Rev A

**BIO-RAD****Bio-Rad  
Laboratories****Life Science  
Group**

**Website** [www.bio-rad.com](http://www.bio-rad.com) **Bio-Rad Laboratories Main Office** 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800  
**Also in:** **Australia** Ph. 02-9914-2800, Fx. 02-9914-2889 **Austria** Ph. (1)-877 89 01, Fx. (1) 876 56 29 **Belgium** Ph. 09-385 55 11, Fx. 09-385 65 54  
**Canada** Ph. (905) 712-2771, Fx. (905) 712-2990 **China** Ph. (86-10) 2046622, Fx. (86-10) 2051876 **Denmark** Ph. 39 17 9947, Fx. 39 27 1698  
**Finland** Ph. 90 804 2200, Fx. 90 804 1100 **France** Ph. (1) 43 90 46 90, Fx. (1) 46 71 24 67 **Germany** Ph. 089 31884-0, Fx. 089 31884-100  
**Hong Kong** Ph. 7893300, Fx. 7891257 **India** Ph. 91-11-461-0103, Fx. 91-11-461-0765 **Israel** Ph. 03 951 4127, Fx. 03 951 4129  
**Italy** Ph. 02-21609.1, Fx. 02-21609.399 **Japan** Ph. 03-5811-6270, Fx. 03-5811-6272 **The Netherlands** Ph. 0313 18-540666, Fx. 0313 18-542216  
**New Zealand** Ph. 09-443 3099, Fx. 09-443 3097 **Singapore** Ph. (65) 272-9877, Fx. (65) 273-4838 **Spain** Ph. (91) 661 70 85, Fx. (91) 661 96 98  
**Sweden** Ph. 46 (0) 8 627 50 00, Fx. 46 (0) 8 627 54 00 **Switzerland** Ph. 01-809 55 55, Fx. 01-809 55 00  
**United Kingdom** Ph. 0800 181134, Fx. 01442 259118