



# Configuring SpectraMax<sup>®</sup> and FlexStation<sup>®</sup> Multi-Mode Microplate Readers for IMAP<sup>®</sup> Assays

## Set Up Guide

## Instrument settings for IMAP Fluorescence Polarization (FP) with FAM-labeled substrates

Parameter	SpectraMax i3x*	SpectraMax Paradigm*	SpectraMax iD5*	SpectraMax M5/M5e FlexStation 3
Optical Configuration	FP-FLUO	FP-FLUO	N/A	N/A
Detection Cartridge (if required)	Fluorescence Polarization Detection Cartridge (Fluorescein)	Fluorescence Polarization Detection Cartridge (Fluorescein)	N/A	N/A
Read Mode	FP	FP	FP	FP
Read Type	Endpoint	Endpoint	Endpoint	Endpoint
Wavelengths	Excitation: 485 nm Emission: 535 nm	Excitation: 485 nm Emission: 535 nm	Excitation: <input checked="" type="checkbox"/> Use Filter 485 nm Emission: <input checked="" type="checkbox"/> Use Filter 535 nm	Excitation: 485 nm Emission: 525 nm Cutoff: 515 nm
Plate Type	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm
Read Area	[user-defined]	[user-defined]	[user-defined]	[user-defined]
PMT and Optics	On the Fly Detection: Off – Stop and Go Integration Time: 100–500 ms Read Height: [user-optimized]**	On the Fly Detection: Off – Stop and Go Integration Time: 100–500 ms Read Height: [user-optimized]**	PMT Gain: Automatic or Medium Integration Time: 100–500 ms Read Height: [user-optimized]**	PMT Gain: High or Medium Flashes per read: 100
More Settings	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	<input checked="" type="checkbox"/> Calibrate <input checked="" type="checkbox"/> Settling Time: 100 ms
G factor (set in Data Reduction)	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]

\*See also the section "Accessories required for IMAP detection with SpectraMax readers" on page 6.

\*\*Optimize microplate and read height by checking the box next to "Show Pre-Read Optimization Options" in More Settings; after clicking Read, follow the instructions as prompted.

## Instrument settings for IMAP FP with TAMRA-labeled substrates

Parameter	SpectraMax i3x*	SpectraMax Paradigm*	SpectraMax iD5*	SpectraMax M5/M5e FlexStation 3
Optical Configuration	FP-RHOD	FP-RHOD	N/A	N/A
Detection Cartridge (if required)	Fluorescence Polarization Detection Cartridge (Rhodamine)	Fluorescence Polarization Detection Cartridge (Rhodamine)	N/A	N/A
Read Mode	FP	FP	FP	FP
Read Type	Endpoint	Endpoint	Endpoint	Endpoint
Wavelengths	Excitation: 535 nm Emission: 595 nm	Excitation: 535 nm Emission: 595 nm	Excitation: <input checked="" type="checkbox"/> Use Filter 535 nm Emission: <input checked="" type="checkbox"/> Use Filter 595 nm	Excitation: 530 nm Emission: 590 nm Cutoff: 570 nm
Plate Type	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm
Read Area	[user-defined]	[user-defined]	[user-defined]	[user-defined]
PMT and Optics	On the Fly Detection: Off – Stop and Go Integration Time: 100–500 ms Read Height: [user-optimized]**	On the Fly Detection: Off – Stop and Go Integration Time: 100–500 ms Read Height: [user-optimized]**	PMT Gain: Automatic or Medium Integration Time: 100–500 ms Read Height: [user-optimized]**	PMT Gain: High or Medium Flashes per read: 100
More Settings	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	<input checked="" type="checkbox"/> Calibrate <input checked="" type="checkbox"/> Settling Time: 100 ms
G factor (set in Data Reduction)	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]

\*See also the section “Accessories required for IMAP detection with SpectraMax readers” on page 6.

\*\*Optimize microplate and read height by checking the box next to “Show Pre-Read Optimization Options” in More Settings; after clicking Read, follow the instructions as prompted.

## Instrument settings for IMAP TR-FRET with FAM-labeled substrates

Parameter	SpectraMax i3x*	SpectraMax Paradigm*	SpectraMax iD5*	SpectraMax M5/M5e FlexStation 3
Optical Configuration (Cartridge)	[Custom Order]	[Custom Order]	N/A	N/A
Detection Cartridge (if required)	[Custom Order]	[Custom Order]	N/A	N/A
Read Mode	TR-FRET	TR-FRET	TR-FRET	TRF
Read Type	Endpoint	Endpoint	Endpoint	Endpoint
Wavelengths	Excitation: 340 nm Emission 1: 490 nm Emission 2: 520 nm	Excitation: 340 nm Emission 1: 490 nm Emission 2: 520 nm	Excitation: <input checked="" type="checkbox"/> Use Filter 340 nm Emission: <input checked="" type="checkbox"/> Use Filter 490 nm <input checked="" type="checkbox"/> Use Filter 520 nm	Excitation: Lm1: 330 nm Lm2: 330 nm Emission: Lm1: 490 nm Lm2: 520 nm Cutoff: Lm1: 475 nm Lm2: 515 nm
Plate Type	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm
Read Area	[user-defined]	[user-defined]	[user-defined]	[user-defined]
TRF Settings				Integration Delay: 200 $\mu$ s Integration Time: 1000 $\mu$ s
PMT and Optics	On the Fly Detection: Off – Stop and Go Number of Pulses: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1 ms Read Height: [user-optimized]**	On the Fly Detection: Off – Stop and Go Number of Pulses: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1 ms Read Height: [user-optimized]**	Flashes per read: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1.0 ms Read Height: [user-optimized]**	Flashes per read: 100
More Settings	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	<input checked="" type="checkbox"/> Calibrate
G factor (set in Data Reduction)	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]	1.0 [or user-defined]

\*See also the section “Accessories required for IMAP detection with SpectraMax readers” on page 6.

\*\*Optimize microplate and read height by checking the box next to “Show Pre-Read Optimization Options” in More Settings; after clicking Read, follow the instructions as prompted.

## IMAP TR-FRET with TAMRA-labeled substrates

Parameter	SpectraMax i3x*	SpectraMax Paradigm*	SpectraMax iD5*	SpectraMax M5/M5e FlexStation 3
Optical Configuration (Cartridge)	[Custom Order]	[Custom Order]	N/A	N/A
Detection Cartridge (if required)	[Custom Order]	[Custom Order]	N/A	N/A
Read Mode	TR-FRET	TR-FRET	TR-FRET	TRF
Read Type	Endpoint	Endpoint	Endpoint	Endpoint
Wavelengths	Excitation: 340 nm Emission 1: 545 nm Emission 2: 570 nm	Excitation: 340 nm Emission 1: 545 nm Emission 2: 570 nm	Excitation: <input checked="" type="checkbox"/> Use Filter 340 nm Emission: <input checked="" type="checkbox"/> Use Filter 535 nm <input checked="" type="checkbox"/> Use Filter 595 nm	Excitation: Lm1: 330 nm Lm2: 330 nm Emission: Lm1: 545 nm Lm2: 570 nm Cutoff: Lm1: 530 nm Lm2: 550 nm
Plate Type	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm [user-optimized]**	384-well: 384 Well Corning flatbtm
Read Area	[user-defined]	[user-defined]	[user-defined]	[user-defined]
TRF Settings				Integration Delay: 200 $\mu$ s Integration Time: 1000 $\mu$ s
PMT and Optics	On the Fly Detection: Off – Stop and Go Number of Pulses: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1 ms Read Height: [user-optimized]**	On the Fly Detection: Off – Stop and Go Number of Pulses: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1 ms Read Height: [user-optimized]**	Flashes per read: 100 Excitation Time: 0.05 ms Measurement Delay: 0.2 ms Integration Time: 1.0 ms Read Height: [user-optimized]**	Flashes per read: 100
More Settings	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	Read Order: Row or Column Show Pre-Read Optimization Options <input checked="" type="checkbox"/>	<input checked="" type="checkbox"/> Calibrate

\*See also the section "Accessories required for IMAP detection with SpectraMax readers" on page 6.

\*\*Optimize microplate and read height by checking the box next to "Show Pre-Read Optimization Options" in More Settings; after clicking Read, follow the instructions as prompted.

## Accessories required for IMAP detection with SpectraMax readers

### SpectraMax i3x or SpectraMax Paradigm reader

#### IMAP FP

- FAM detection requires a Fluorescence Polarization (FP) (Fluorescein Detection Cartridge (Molecular Devices P/N 0200-7009)
- TAMRA detection requires a Fluorescence Polarization (FP) (Rhodamine) Detection Cartridge (Molecular Devices P/N 0200-7010)

#### IMAP TR-FRET

- FAM or TAMRA detection requires a custom order. A custom detection cartridge request form is available through sales representatives.

### SpectraMax iD5 reader

#### IMAP FP

- Filters for FAM detection:
  - Set of 2 Fluorescence Polarization Filters 485 nm BW 25 nm Polarized Vertical & Horizontal (Molecular Devices P/N 6590-0136)
  - Set of 2 Fluorescence Polarization Filters 535 nm BW 25 nm Polarized Vertical & Horizontal (Molecular Devices P/N 6590-0137)
- Filters for TAMRA detection:
  - Set of 2 Fluorescence Polarization Filters 535 nm BW 25 nm Polarized Vertical & Horizontal (Molecular Devices P/N 6590-0137)
  - Set of 2 Fluorescence Polarization Filters 595 nm BW 25 nm Polarized Vertical & Horizontal (Molecular Devices P/N 6590-0139)

#### IMAP TR-FRET

- Enhanced TRF Module (Molecular Devices P/N 0200-7030)
- Filters for FAM detection:
  - Time Resolved Filter with holder 340nm BW 70nm (Molecular Devices P/N 6590-0080)
  - Time Resolved Filter with holder 490nm BW 10nm (Molecular Devices P/N 6590-0095)
  - Time Resolved Filter with holder 520nm BW 15nm (Molecular Devices P/N 6590-0098)
- Filters for TAMRA detection:
  - Time Resolved Filter with holder 340nm BW 70nm (Molecular Devices P/N 6590-0080)
  - Fluorescence Filter with holder 535nm BW 25nm (Molecular Devices P/N 6590-0099)
  - Fluorescence Filter with holder 595nm BW 35nm (Molecular Devices P/N 6590-0105)

## IMAP® microplate reader compatibility

Product Name	Part Number	SpectraMax® M5/M5e	FlexStation®	SpectraMax iD5	SpectraMax i3x	SpectraMax Paradigm	SpectraMax M4
<b>Demo Kit</b>							
IMAP® Evaluation Demo Kit	R8166	✓	✓	✓	✓	✓	(TR-FRET only)
<b>Evaluation Kits</b>							
IMAP FP Evaluation Kit	R8155	✓	✓	✓	✓	✓	
IMAP TR-FRET Evaluation Kit	R8161	✓	✓	✓	✓	✓	✓
IMAP PDE FP Evaluation Kit	R8175	✓	✓	✓	✓	✓	
IMAP PDE TR-FRET Evaluation Kit	R8176	✓	✓	✓	✓	✓	✓
<b>Screening Express Kits</b>							
IMAP FP Screening Express Kit	R8127	✓	✓	✓	✓	✓	
IMAP TR-FRET Screening Express Kit	R8160	✓	✓	✓	✓	✓	✓

## Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)  
Check our website for a current  
listing of worldwide distributors.

## Regional Offices

USA and Canada	<a href="tel:+18006355577">+1.800.635.5577</a>	Taiwan/Hong Kong	<a href="tel:+886226567585">+886.2.2656.7585</a>
United Kingdom	<a href="tel:+441189448000">+44.118.944.8000</a>	Japan	<a href="tel:+81363629109">+81.3.6362.9109</a>
Europe*	<a href="tel:0080066532860">00800.665.32860</a>	South Korea	<a href="tel:+82234719531">+82.2.3471.9531</a>
China	<a href="tel:+864008203586">+86.4008203586</a>	India	<a href="tel:+917386611198">+91.73.8661.1198</a>

*\*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom*