# FocalCheck<sup>™</sup> Fluorescence Microscope Test Slides

## F36909 FocalCheck™ fluorescence microscope test slide #1 F36913 FocalCheck™ fluorescence microscope test slide #2 F36914 FocalCheck™ fluorescence microscope test slide #3

Table 1. Contents and storage information for FocalCheck™ fluorescence microscope test slide #1

	1	2	3	4	5	Storage and stability	
Row A	15 μm	6 µm	4 µm	1 μm	0.5 μm	<ul> <li>Room temperature</li> <li>Avoid prolonged exposure to light</li> <li>When stored as directed, the slides should be stable for at least 12 months.</li> </ul>	
	Green/orange/far red thin rings	Blue throughout/ green ring	Blue/green/red/ far red	Blue/green/red/ far red	Blue/green/red/ far red		
Row B (Intensity Series)	6 µm	6 µm	6 µm	6 µm	6 µm		
	Green	Green	Green	Green	Green		
	100% intensity	33% intensity	10% intensity	3% intensity	0.667% intensity		

Table 2. Contents and storage information for FocalCheck™ fluorescence microscope test slide #2

	1	2	3	4	5	Storage and stability	
	6 μm	6 µm	6 µm	6 µm		• Room	
	Green	Orange	Red	Far red	Mixture of all eight reference	temperature • Avoid	
	Bead 1 503/511 nm Bead 2 511/524 nm	Bead 1 541/555 nm Bead 2 545/565 nm	Bead 1 578/605 nm Bead 2 589/613 nm	Bead 1 657/676 nm Bead 2 671/692 nm	beads	prolonged exposure to light	
Row B	6 μm	6 µm	6 µm	6 µm			
	Green	Orange	Red	Far red	Mixture of all	When stored as directed, the slides should be stable for at least 12 months.	
	Ring 503/511 nm Core 511/524 nm	Ring 541/555 nm Core 545/565 nm	Ring 589/613 nm Core 578/605 nm	Ring 671/692 nm Core 657/676 nm	four double- stained beads		

Table 3. Contents and storage information for FocalCheck™ fluorescence microscope test slide #3

	1	2	3	4	5	Storage and stability	
Row A	2.5 μm						
	Blue	Green	Orange	Red	Deep red		
	350/440 nm	505/515 nm	540/560 nm	580/605 nm	633/660 nm		
	Bright	Bright	Bright	Bright	Bright	<ul><li> Room temperature</li><li> Avoid prolonged</li></ul>	
	100% intensity	exposure to light					
Row B	2.5 μm	When stored as directed,					
	Blue	Green	Orange	Red	Deep red	the slides should be stable for at least 12 months.	
	350/440 nm	505/515 nm	540/560 nm	580/605 nm	633/660 nm		
	Dim	Dim	Dim	Dim	Dim		
	1% intensity						

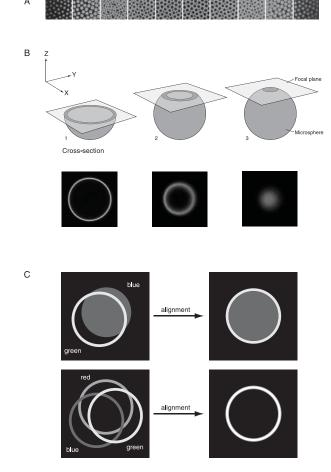
FocalCheck<sup>™</sup> fluorescence microscope test slides #1, #2, and #3 are specifically designed for calibrating fluorescence microscope systems and evaluating system and filter performance. The slides each contain 10 sample areas coated with proprietary fluorescent microspheres designed specifically for microscopy applications. The microspheres are mounted in optical cement (refractive index of ~1.52) for maximum stability and are intended for use on all fluorescence microscope systems.

- **FocalCheck<sup>™</sup> fluorescence microscope test slide #1** is ideal for routine checking and calibration of confocal and widefield fluorescence microscopes.
- FocalCheck<sup>™</sup> fluorescence microscope test slide #2 provides a robust, reproducible method of evaluating the performance of spectral imaging systems, as well as the ability to discriminate closely overlapping spectra. The slide consists of 6-µm-diameter microspheres labeled with a series of spectrally overlapping dyes.
- FocalCheck<sup>™</sup> fluorescence microscope test slide #3 is useful for basic evaluation of filter performance and as a general practice slide for fluorescence microscopy and digital imaging.

## **Experimental Protocols**

The basic procedure for using the FocalCheck<sup>™</sup> fluorescence microscope test slides is to image the beads using standard techniques. Below are several general applications.

**Color registration** For multicolor experiments, particularly on widefield microscopes, switching between different filter configurations can result in misregistration between the various color channels. Using the multicolor beads on FocalCheck<sup>™</sup> fluorescence microscope test slide #1 (sample positions A1–5), the extent of the misregistration can be digitally corrected independently of the sample (Figure 1).



**Figure 1.** Confocal laser-scanning microscope optical cross-sectioning and alignment with FocalCheck<sup>™</sup> microspheres. A) Serial optical sectioning from top-to-bottom along the Z-axis of ring-stained microspheres reveals a continuous pattern of disc-to-ring-to-disc images. B) The diameter of the fluorescent ring (or disc) is dependent on the depth of the optical focal plane. C) In the confocal laser-scanning microscope, separate light paths exist for UV and visible wavelengths. Also, emitted fluorescence is detected by different photomultipliers. Proper optical alignment may be obtained with either of two types of FocalCheck<sup>™</sup> microspheres. For example, the microspheres with a green ring stain that are blue-fluorescent throughout the bead allow UV/visible wavelength alignment in three dimensions upon aligning the green ring with the blue disc. Focal alignment is also possible simultaneously in three colors by aligning the red, blue, and green rings of the FocalCheck<sup>™</sup> microspheres containing fluorescent red, blue, and green ring stains.

**Chromatic distortions** Modern microscopes use highly color-corrected (achromatic) optical components. However, it is possible that certain optical components may introduce wavelength-dependent image shifts that can affect the correct registration of multicolor labels. Such aberrations are critical to avoid in colocalization experiments. FocalCheck<sup>™</sup> fluorescence microscope slide #1 is ideal for ensuring that the imaging system is free from chromatic distortions.

**X- and Y-axis calibration** For distance and other morphometric measurements, microscope imaging should be calibrated for each objective and magnification setting using a stage micrometer. The larger 15-μm beads on FocalCheck<sup>™</sup> fluorescence microscope test slide #1, sample positions A1 and A2, provide ideal targets for assessing the calibration of the microscope in three dimensions. When measured at the equator, the beads are nominally 15 μm in diameter; this should correspond with the measured X and Y dimensions by imaging.

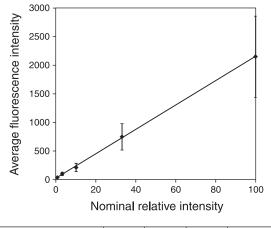
**Z-axis calibration** Positions A1 and A2 on FocalCheck<sup>™</sup> fluorescence microscope test slide #1 can be used to assess the performance of the motorized focus control of an imaging system. To do this, acquire Z-series stacks from bottom-to-top and top-to-bottom. The refractive index of the slide mounting media is close to 1.515, which minimizes spherical aberration when using

an oil immersion lens. When viewed in orthogonal projections, the bead images should appear relatively circular. Any distortions could indicate problems in the Z-stepping of the microscope's focus system or a refractive index mismatch.

Lamp stability Most standard mercury lamps are rated for 200–300 hours of use. While many can last longer, there is often little warning that the lamp is failing. Bead samples on any of the FocalCheck<sup>™</sup> fluorescence microscope test slides can be used to monitor the stability of the excitation source on the microscope. Periodic imaging of the same sample, using the same acquisition parameters, can indicate changes in lamp output by plotting the intensity of the beads over time.

Intensity calibration FocalCheck<sup>™</sup> fluorescence microscope test slide #1 contains a series of intensity beads. By acquiring a series of images, you can create a relative intensity curve against which samples can be evaluated (Figure 2). The intensity of the beads can be used to generate curves for different instruments, thus providing a basis for assessing the relative fluorescence intensity of a sample in a system-independent manner.

FocalCheck<sup>™</sup> fluorescence microscope test slide #1 can also be used to examine the linearity and dynamic range of the detection system on the microscope under various settings. The intensity range of the beads covers the approximate dynamic range of a typical 12-bit CCD camera (4096 gray levels) when imaged at a single exposure setting.



Nominal relative intensity	1.00%	5.00%	10.00%	40.00%	100%
Actual intensity	18.81	88.70	201.91	738.01	2138.48
Determined relative intensity	0.88%	4.15%	9.44%	34.51%	100%

**Figure 2.** Average fluorescence intensity of the intensity beads series on FocalCheck<sup>™</sup> microscope test slide #1. The intensity bead series was imaged on a Nikon Eclipse 800 microscope using a 100-watt Hg-arc lamp and a 60×/1.4NA objective. The images were acquired using a Hamamatsu Orca-ER 12-bit CCD camera. The average fluorescence intensity and standard deviation of the bead targets were determined and plotted against nominal relative intensity as determined by flow cytometry. In this particular configuration, all five intensities were imaged using a single exposure setting. The measured average intensities shown in the table give relative intensities that correspond closely to the nominal values measured by flow cytometry.

### Spectral scanning

FocalCheck<sup>™</sup> fluorescence microscope test slide #2 can be used on spectral imaging instruments to help ensure that the wavelength detection system is working correctly. This slide can also be used to perform many of the general tests described above.

### Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat no.	Product name	Unit size
F36909	FocalCheck™ fluorescence microscope test slide #1 *for alignment, intensity, and calibration*	each
F36913	FocalCheck™ fluorescence microscope test slide #2 *for spectral imaging systems*	each
F36914	FocalCheck™ fluorescence microscope test slide #3 *5 colors, high and low intensities*	each
P7220	PS-Speck™ Microscope Point Source Kit *blue, green, orange, and deep red fluorescent beads*	1 kit

# **Contact Information**

#### **Corporate Headquarters**

5791 Van Allen Way Carlsbad, CA 92008 USA Phone: +1 760 603 7200 Fax: +1 760 602 6500 Email: techsupport@lifetech.com

### **European Headquarters**

Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF UK Phone: +44 141 814 6100 Toll-Free Phone: 0800 269 210 Toll-Free Tech: 0800 838 380 Fax: +44 141 814 6260 Tech Fax: +44 141 814 6117 Email: euroinfo@invitrogen.com Email Tech: eurotech@invitrogen.com

#### Japanese Headquarters

LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Japan Phone: +81 3 5730 6509 Fax: +81 3 5730 6519 Email: jpinfo@invitrogen.com

Additional international offices are listed at www.lifetechnologies.com.

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

#### **Obtaining Support**

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- · Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- · Search through frequently asked questions (FAQs)
- · Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations. and other product support documents
- Obtain information about customer training
- Download software updates and patches

#### SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

#### Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to **www.lifetechnologies.com/support** and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

#### **Limited Warranty**

Invitrogen (a part of Life Technologies Corporation) is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Support Representatives.

All Invitrogen products are warranted to perform according to specifications stated on the certificate of analysis. The Company will replace, free of charge, any product that does not meet those specifications. This warranty limits the Company's liability to only the price of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. The Company reserves the right to select the method(s) used to analyze a product unless the Company agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore the Company makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Support Representatives.

Life Technologies Corporation shall have no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

#### Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

©2011 Life Technologies Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.