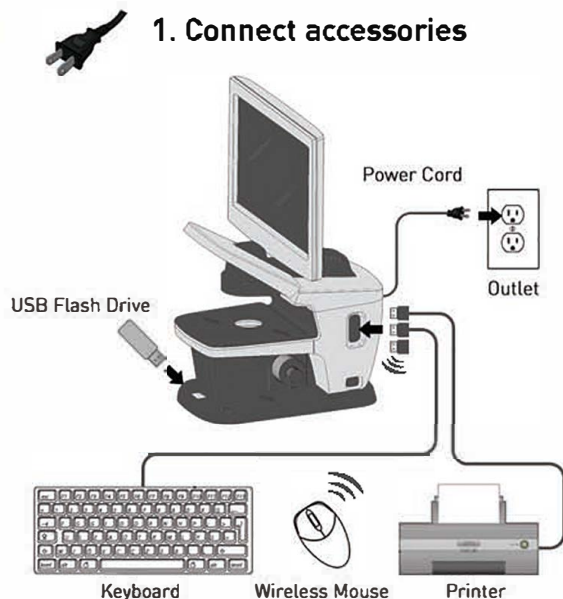


# FLoid® Cell Imaging Station quick start guide

Online Specials

## What's in the box?

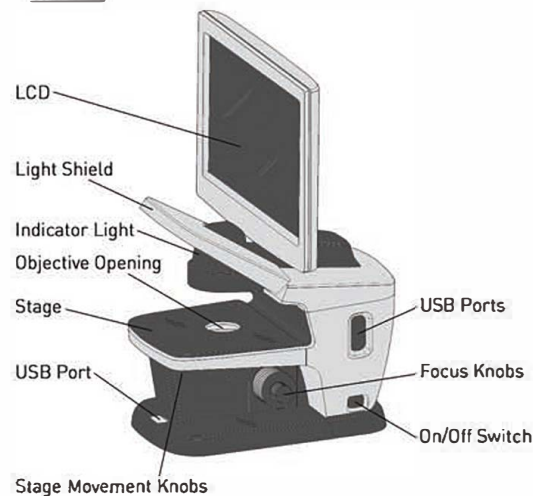
- FLoid® Cell Imaging Station  
[Cat. no. 4471136]
- Power cord
- Keyboard
- Mouse
- USB flash drive  
(with user manual.)
- Quick start guide
- Printer [Cat. no. 4472351]



## 1. Connect accessories



## 2. Power on



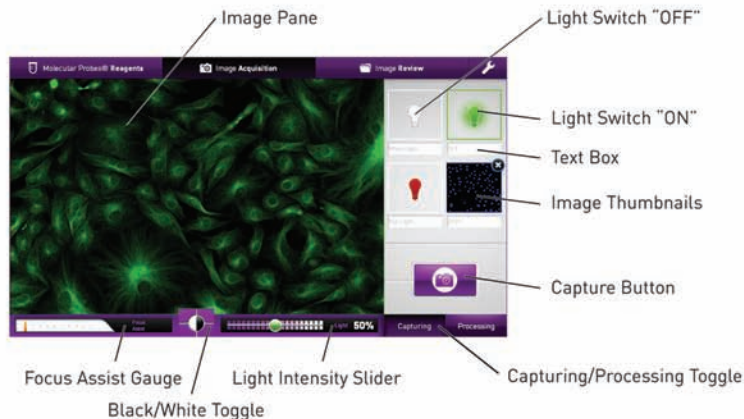
# FLoid® Cell Imaging Station

## quick start guide



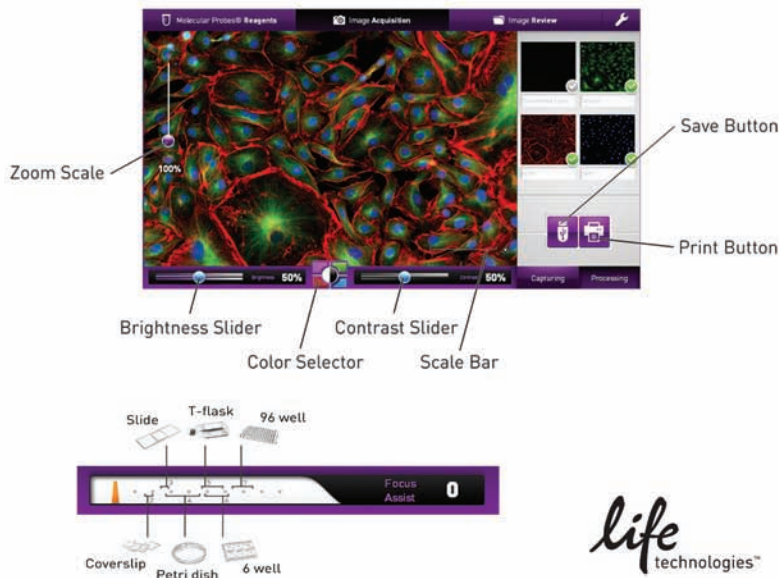
### 3. Capture Image

- Place sample on stage over objective opening
- Turn on light by selecting light bulb icon
- Focus image using focus knobs
- Capture image(s) then press "Processing" button



### 4. Process Image

- Adjust brightness and contrast as needed
- Save image to USB flash drive and print image



#### Limited Product Warranty

Life Technologies Corporation and/or its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](mailto:www.lifetechnologies.com/support).

**DISCLAIMER** - LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATES DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT, TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATES BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners.

3 December 2012

life  
technologies™