

# Cellometer<sup>®</sup>

## Calcein-AM/PI Vitality and Viability Kit

Product Number: CSK-0118

Sample Kit: CSK-0118-S (Not available for purchase)

Cell Type Markers  
Apoptosis

Cell Cycle  
Proliferation  
Metabolism  
**Viability**



This product is for RESEARCH USE ONLY and is not approved for diagnostic or therapeutic use

Page left blank intentionally

# Table of Contents

## 1. Introduction

1.1 Assay Description.....	Page 2
1.2 Materials and Reagents.....	Page 2
1.3 Instrument and Software.....	Page 3

## 2. Assay Protocol

2.1 Preparation of Adherent Cells.....	Page 3
2.2 Preparation of Reagent.....	Page 3
2.3 Staining Procedure.....	Page 3
2.4 Staining Procedure for Clinical Samples.....	Page 3
2.5 Data Acquisition.....	Page 3

## 3. Technical Support..... Page 7

## 4. Software Settings

4.1 Calcein-AM Counting Options Screen.....	Page 7
4.2 Default Calcein-AM / PI Assay and Cell Type Settings.....	Page 8
4.3 Software Settings for Clinical Samples.....	Page 12

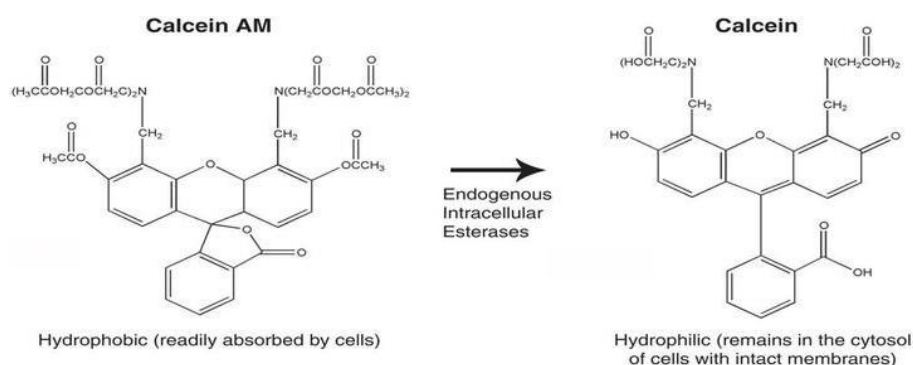
## 5. Importing a New Assay Type, Cell Type, and Templates..... Page 14

5.1 Import Result and Print Templates.....	Page 15
5.2 Import Cell Type.....	Page 15
5.3 Import Assay Type.....	Page 16

## 1.0 Introduction

### 1.1 Assay Description

Calcein AM (Calcein acetoxymethyl ester) is a cell permeable, non-fluorescent compound. Upon crossing the cell membrane, Calcein AM is rapidly hydrolyzed by cellular esterases inside live cells. The hydrolysis cleaves the AM group, converting the non-fluorescent Calcein AM to a strongly green fluorescing Calcein. The more hydrophilic Calcein is trapped inside the cell (1). Cells that do not possess active cytoplasmic esterases are unable to convert Calcein AM to Calcein, and therefore do not fluoresce green. This allows for a quick and easy detection of metabolically-active cells in a sample.



Stains such as propidium iodide (PI), 7-AAD, and ethidium bromide (EB), are membrane exclusion dyes that are frequently used to stain non-viable nucleated cells with compromised membranes. Acridine orange freely diffuses across the cell membrane and stains DNA in all nucleated cells. When AO and PI are combined it is possible to determine % viability for nucleated cells. When Calcein AM is used in conjunction with PI, it is possible to determine % vitality / viability based on the number of metabolically-active (green fluorescent) and non-viable (red fluorescent) cells in a sample.

Since Calcein AM does not require DNA binding, it stains all metabolically-active cells and can be used to measure metabolic activity in non-nucleated cells, such as platelets (4). Calcein-AM is also a good alternative for analysis of adipocytes, as the AO dye has shown some non-specific binding of lipid droplets that does not occur with Calcein AM (5). Because Calcein AM is photostable, shows low cytotoxicity, does not affect cellular functions, and requires cellular esterases for conversion to green-fluorescing Calcein, it is a popular stain for the examination of cell vitality and viability. (1,2,3).

1. Braut-Boucher, F. et al. *Journal of Immunological Methods*. Vol. 178, Issue 41 (1995).
2. Luc S. De Clerck. et al. *Journal of Immunological Methods*. Vol. 172, Issue 1, (1994).
3. Parish, CR. *Immunology and Cell Biology*. Vol. 77 (1999)
4. Verheul, HW. et al. *Blood*. Vol. 96 No. 13 (2000)
5. Kilroy, G. et al. *PLoS One*. Vol.4, Issue 9 (2009)

## 1.2 Materials Included

- Calcein-AM/PI Vitality and Viability Kit  
**Cat. # CSK-0118 (100 tests)**
  - (Component A) Cellometer Propidium Iodide Staining Solution
  - (Component B) Cellometer Calcein-AM Staining Solution

## 1.3 Additional Materials required, but not included

- dH<sub>2</sub>O
- Trypsin EDTA (if working with adherent cells)

## 1.4 Instrument and Software Requirement

- Cellometer Vision or Vision CBA Image Cytometry System
- Cellometer Vision or Vision CBA Software
- Fluorescence Optics Module VB-535-402 and VB-660-502

## 2.0 Assay Protocol for CSK-0118

### Preparation of Adherent Cells for Staining

1. Using 1 x Trypsin-Versene (EDTA), trypsinize cells until they have lifted off the plate (approximately 15 minutes).
2. Use the Cellometer Sample Adjustment Calculator to determine the sample volume required to obtain a concentration of 2-3 x 10<sup>6</sup> cells/ mL.
3. Spin down cells at 1,000 to 2,000 rpm for five minutes.
4. Decant the supernatant and re-suspend cells in 1ml of 1 x PBS or culture media in which the cells were grown.

### 2.2 Preparation of Calcein AM Reagent

1. Pipette 2 µl Calcein-AM (Nexcelom Part# CSK-0118 Component A) into 18 µl of dH<sub>2</sub>O. This is now Calcein-AM Solution A. Mix by pipetting up and down at least 15 times or vortex.

### 2.3 Staining Procedure for Cultured Cells

1. Add 5 µl of Calcein-AM Solution A and 5 µl of PI Staining Solution (Nexcelom Part# CSK-0118 Component B) to 40 µl of cell sample.
2. Gently pipette the sample up and down ten times, then incubate for 20 min at 37°C in the dark.
3. After the 20 minute incubation, the sample is ready for analysis. Proceed to step 2.5.

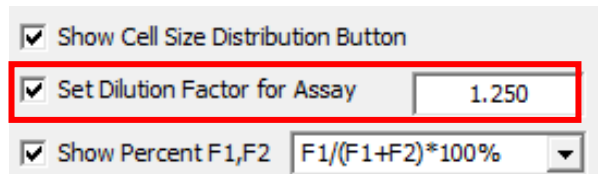
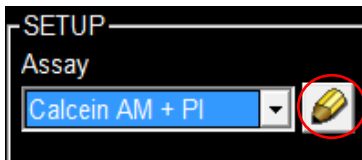
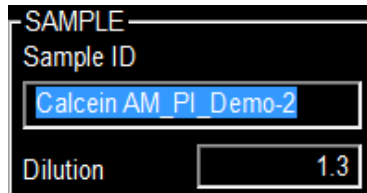
### 2.4 Staining Procedure for Whole Blood, Cord Blood, and Bone Marrow Clinical Samples.

1. Pipette 10 µl of fresh blood sample into 70 µl of 1 x PBS
2. Pipette 40 µl of the diluted blood sample into a new eppendorf tube.
3. Add 5 µl of Calcein-AM Solution A and 5 µl of PI Staining Solution(Nexcelom Part# CSK-0118 Component B) to 40 µl of cell sample.
4. Gently pipette the sample up and down ten times, then incubate for 20 min at 37°C in the dark. After the 20 minute incubation, the sample is ready for analysis. Proceed to step 2.5.

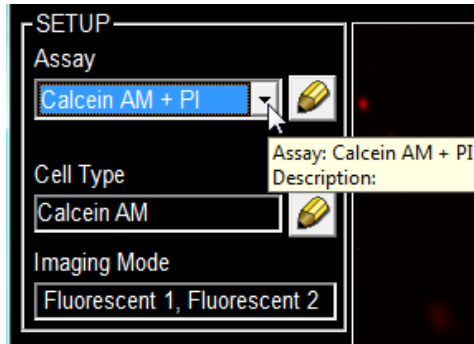
## 2.5 Data Acquisition

**NOTE: Please review the auto save set-up to make sure the acquired data is properly saved (See step 16).**

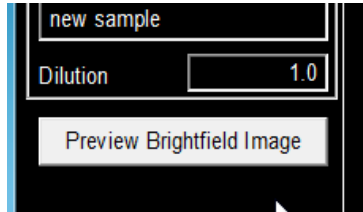
1. Gently mix the cell sample by pipetting up and down at least ten times, then load 20 µL into the Cellometer counting chamber and insert into the Cellometer instrument.
2. Wait 60 seconds for the cells to settle in the chamber
3. Type a name for your sample into the Sample ID text box
4. Verify that the default dilution factor for the Calcein AM + PI (For cultured cells) assay is 1.25 by clicking on the pencil icon and locating the “Set Dilution Factor for Assay” in the dialog pop-up screen. The default dilution factor for the Calcein AM + PI CS (Clinical Samples) assay is 10. If a higher sample dilution was performed in step 2.4, or the cell culture sample was pre-diluted in step 2.3, adjust the dilution factor accordingly. Please note that dilution factor on the main screen is rounded up for display purposes and will read as 1.3.



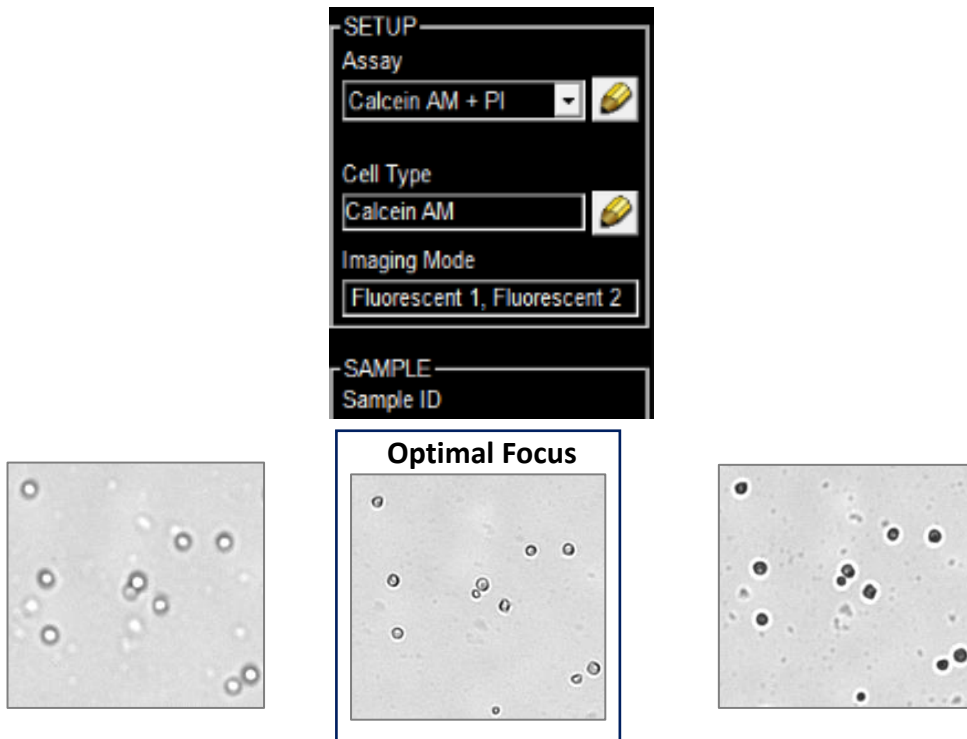
5. Select the **Calcein AM + PI or Calcein AM + CS Assay** from the Assay drop-down menu in the upper left corner of the main Vision CBA software screen. If this assay is not present in your drop-down menu, import the Calcein AM + PI assay files using the instructions in section 5.0.



6. Unless you are testing clinical samples, the Calcein AM + PI assay should be run with the default software settings. If you suspect that the settings may have been changed, review the default software settings in section 4.0. To update the Calcein AM + PI assay to the Calcein AM + PI CS assay for whole blood, cord blood, and bone marrow clinical samples see section 4.3.
7. Click Preview Brightfield Image at the bottom left of the main Vision screen.



8. Turn the focus knob and adjust focus for the bright field image. Cells in focus for the Calcein-AM/PI assay will have a bright center and dark outline. There should be a crisp contrast between background and the cell membrane. See Focus Guide below.

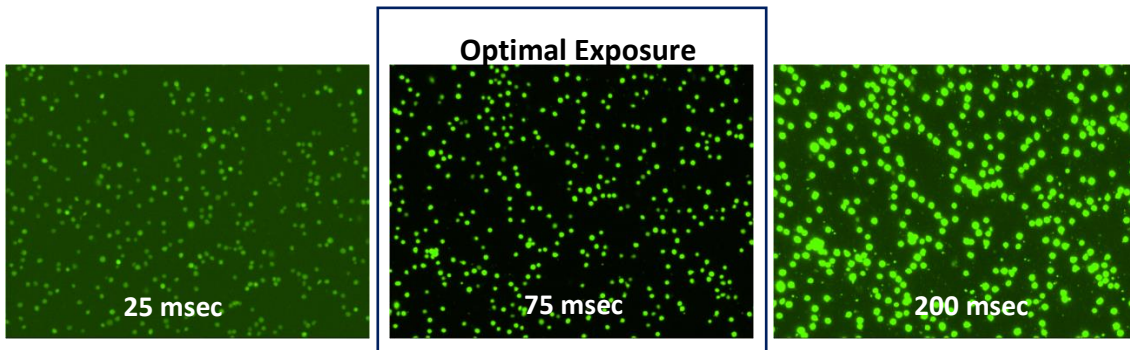
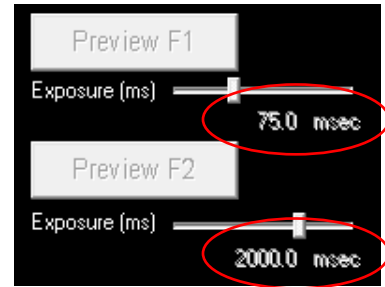




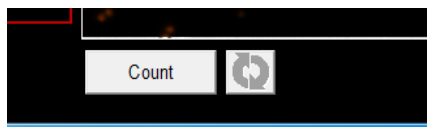
9. Click Stop Preview
10. Click the Preview F1 Image button (bottom left of screen) and verify that the fluorescence signal displays as 100% of range.

For the Calcein AM + PI Assay the default exposure time is 75 milliseconds (msec). For the Calcein AM + PI CS Assay the optimal exposure is 200 milliseconds (msec). Optimal exposure time will generate a bright image with well-defined fluorescent spots.

Under-exposure will yield dark images with weak spots, like the 25 msec image below, with insufficient fluorescent signal. Over-exposure will yield images that are too bright with fluorescent spots that are large and sometimes overlapping, as shown in the 200 ms image below. Spots are also less distinct from background.



11. Click Preview F2. The F2 (PI) channel should be set to 2000 msec exposure for the Calcein AM + PI assay and 4000 msec for the Calcein AM + PI SC assay.
12. Click the Count button at the bottom of the screen



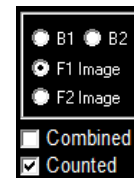
13. When counting is complete, an initial Results Table will appear on the screen. For optimal counting results, the total cell concentration should be between  $3 \times 10^5$  cells/mL and  $1 \times 10^7$  cells/mL (see circled value in the report below). Recount a concentrated or diluted cell sample if necessary.

<b>Assay: Calcein AM + PI</b>		<b>Date: 10/25/2012 14:02:18</b>	Show Size Distribution
F1 Channel: Calcein AM - Green F2 Channel: Propidium Iodide - Red			Intensity Distribution
Sample ID: 10uM_75msec_3-2 Dilution: 1.30		Instrument Serial #: Vision-301-0088 Instrument Optics: X050-Vx-535-402-Vx-660-502	
<b>Results:</b>			
Count	Concentration	Mean Diameter	
=====	=====	=====	
Calcein positive: 1724	$3.17 \times 10^6$ cells/mL	10.7 micron	
PI positive cells: 1560	$2.86 \times 10^6$ cells/mL	9.3 micron	
Total cells: 3284	$6.03 \times 10^6$ cells/mL		
Vitality/Viability: 52.6%			

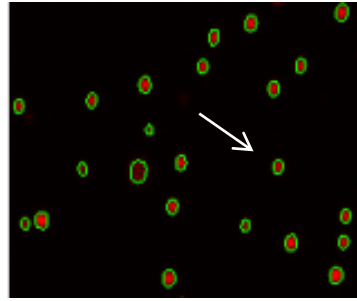
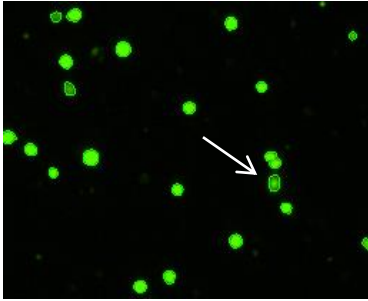
The default report displays:

- **Calcein-positive (metabolically-active) cells** counted, cell concentration, and mean cell diameter
- **PI-positive (non-viable) cells** counted, cell concentration, and mean cell diameter
- **Total cells** counted and total cell concentration
- **% Vitality/Viability** for the cell sample: Calcein-positive cells / (Calcein-positive cells + PI-positive cells)

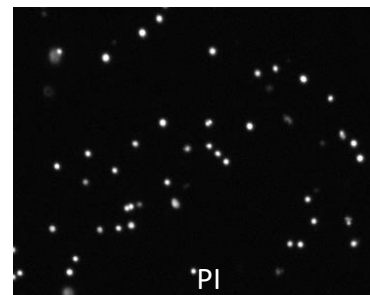
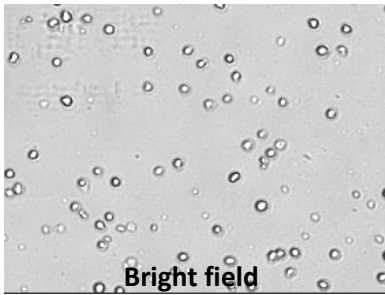
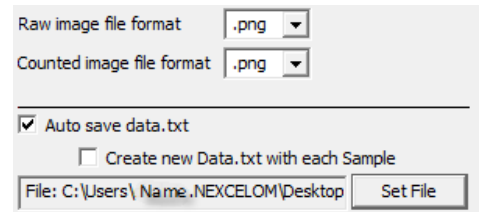
14. Click the **Close** button at the bottom right corner of the Counting Results table. Select the F1 Image and check Counted at the right-hand side of the screen.



15. Review the counted image to confirm that the Calcein-positive cells are being counted correctly. Individual cells within clumps should be circled in green, indicating that they are being counted individually. Click F2 Image to confirm that the PI stained cells are being counted correctly as well. If cells are not being counted correctly, please contact Nexcelom Technical Support for assistance with optimization of counting parameters (see Section 3.0).



16. To view bright field and fluorescent cell images that have been saved, open the image folder where the data has been saved. Bright field and fluorescent images are captured for the Calcein AM / PI assay and may be saved automatically. Saved **Raw** images may be opened in the Cellometer software for re-analysis. Images are only saved if “Auto save data.txt” is selected in the save options menu. The options menu is found by selecting options, then “save options” at the top of the main screen.



A screen-capture software may be used to save both uncounted and counted colored fluorescent images for presentation and publication.

### 3.0 Technical Support

4.0 Nexcelom Technical Support is available from 9am to 5pm EST.

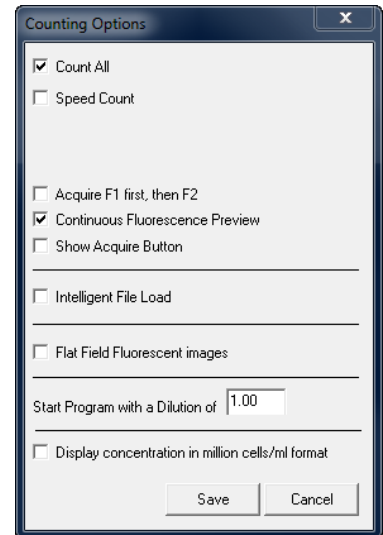
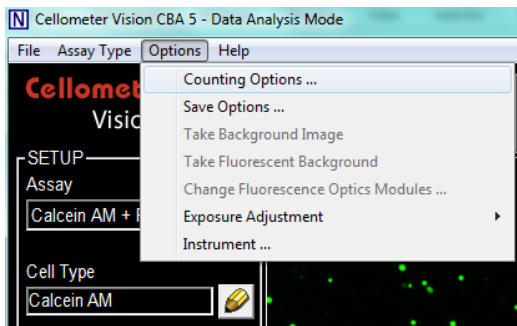
5.0 E-mail: [support@nexcelom.com](mailto:support@nexcelom.com)

6.0 Phone: 978-327-5340

### 4.0 Software Settings

#### 4.1 Review Calcein AM/PI Counting Options Screen

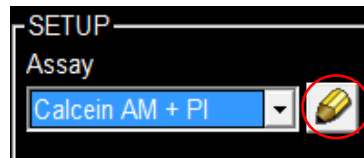
Click on the Options Page and select Counting Options. If you suspect that default settings may have been changed, verify that all selections on the instrument screen match the default settings below.



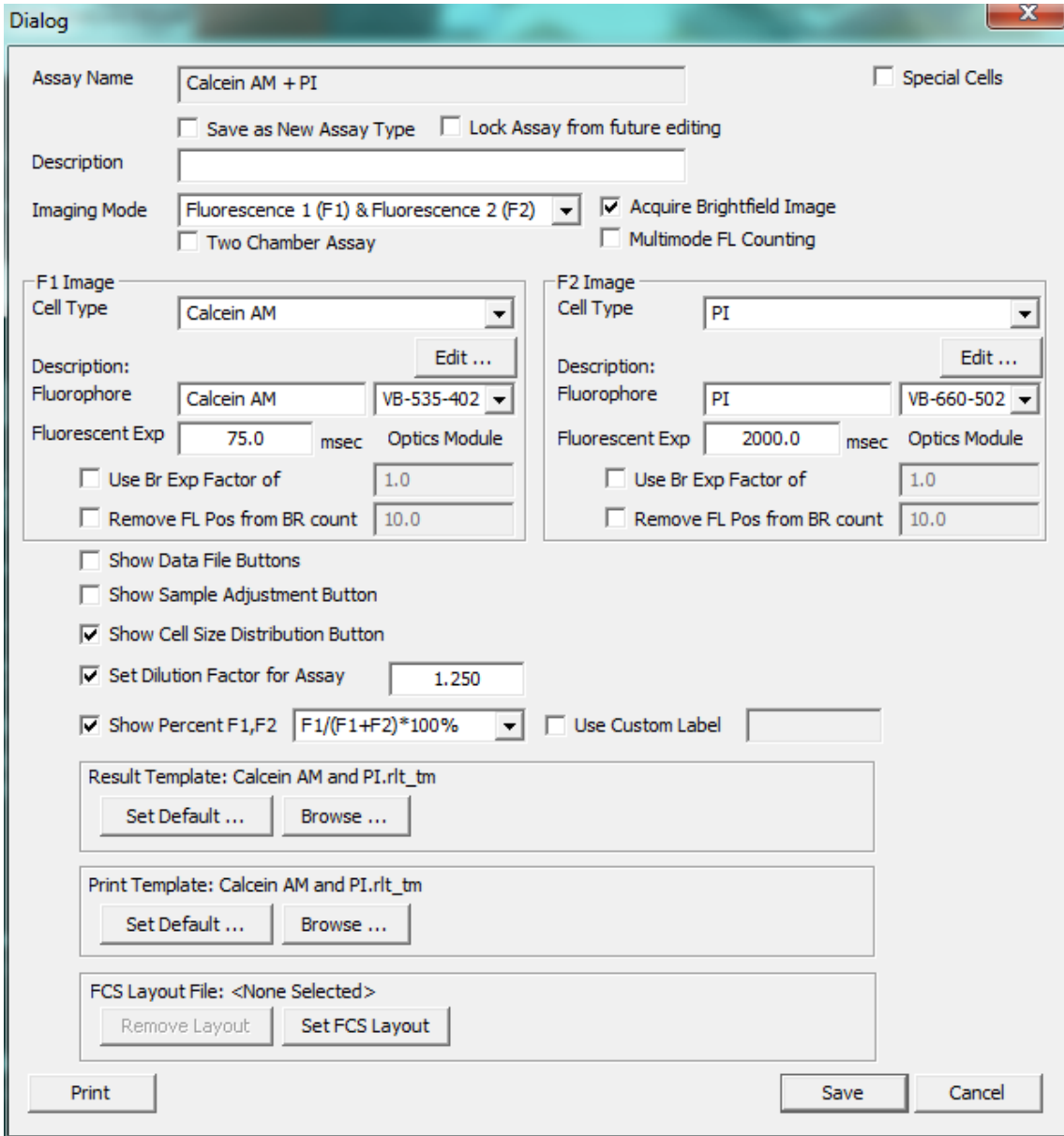
#### 4.2 Default (Calcein AM / PI) Software Settings

##### 4.2.1 Check Dialog Screen Settings

1. Click on the pencil icon under Assay on the main Vision CBA screen.

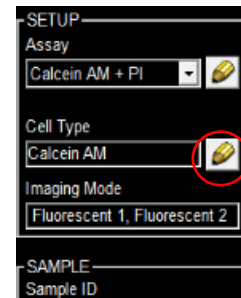


- Verify that all selections on the instrument screen match the default settings below.



#### 4.2.2 Check Cell Type Settings

- Click on the pencil icon under Cell Type on the main Vision CBA software screen to check the Calcein AM cell type settings.



- Verify that all selections for the bright field (BR) tab on the instrument screen match the default settings below.

Cell Type

Cell Type Name   Save as New Cell Type  
 Lock from future editing

Detailed Description

Brightfield (BR) | Fluorescence (FL)

	Minimum		Maximum	
Cell Diameter	<input type="text" value="6.0"/>	micron	<input type="text" value="30.0"/>	micron
Roundness	<input type="text" value="0.10"/>	default: 0.10; range: 0 - 1.0; 1.0 for perfect circle		
Contrast Enhancement	<input type="text" value="0.40"/>	default: 0.40; range: 0 - 0.8; high value for light cells		

Deduster Parameters

Do not decluster dumps

Deduster Edge Factor	<input type="text" value="0.5"/>	default: 0.5; range 0 - 1.0; higher value for more edge enhancement
Deduster Th Factor	<input type="text" value="1.0"/>	default: 1.0; range 0 - 1.0; higher value for more sensitivity
Background Adjustment	<input type="text" value="1.0"/>	default: 1.0; range 0 - 1.0; lower value to pick up dim cells

Trypan Blue Viability Parameters

	Minimum		Maximum	
Dead Cell Diameter	<input type="text" value="3.0"/>	micron	<input type="text" value="50.0"/>	micron
Sensitivity	<input type="text" value="1.0"/>	default: 1.0; range 0 - 6.0; higher value to pick up more dead cells		
Uniformity	<input type="text" value="150"/>	default: 150; range 100 - 255 higher value for non-uniform dead cells		

Very Dim Dead Cells

Print Save Cancel

- Click on the Fluorescence (FL) tab.

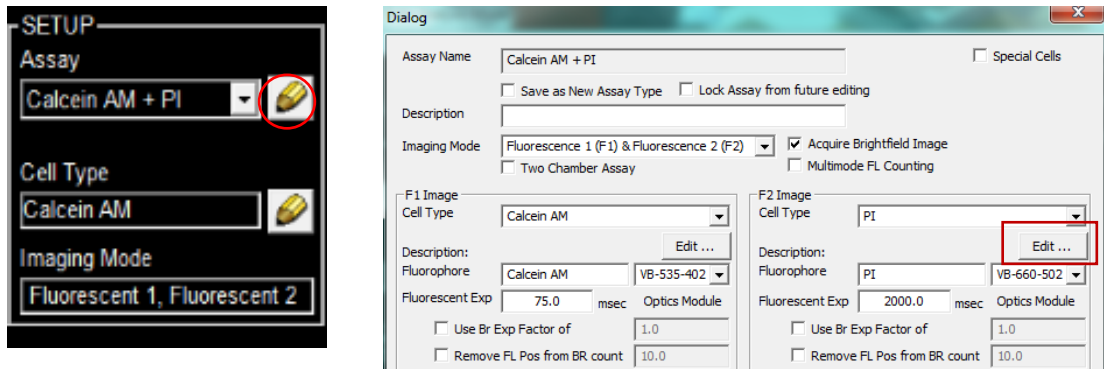
The screenshot shows the 'Cell Type' dialog box with the 'Fluorescence (FL)' tab selected. The 'Cell Type Name' is 'Calcein AM' and the 'Detailed Description' field is empty. The 'Brightfield (BR)' and 'Fluorescence (FL)' tabs are visible at the bottom.

- Verify that all selections for the Fluorescence (FL) tab on the instrument screen match the default settings below.

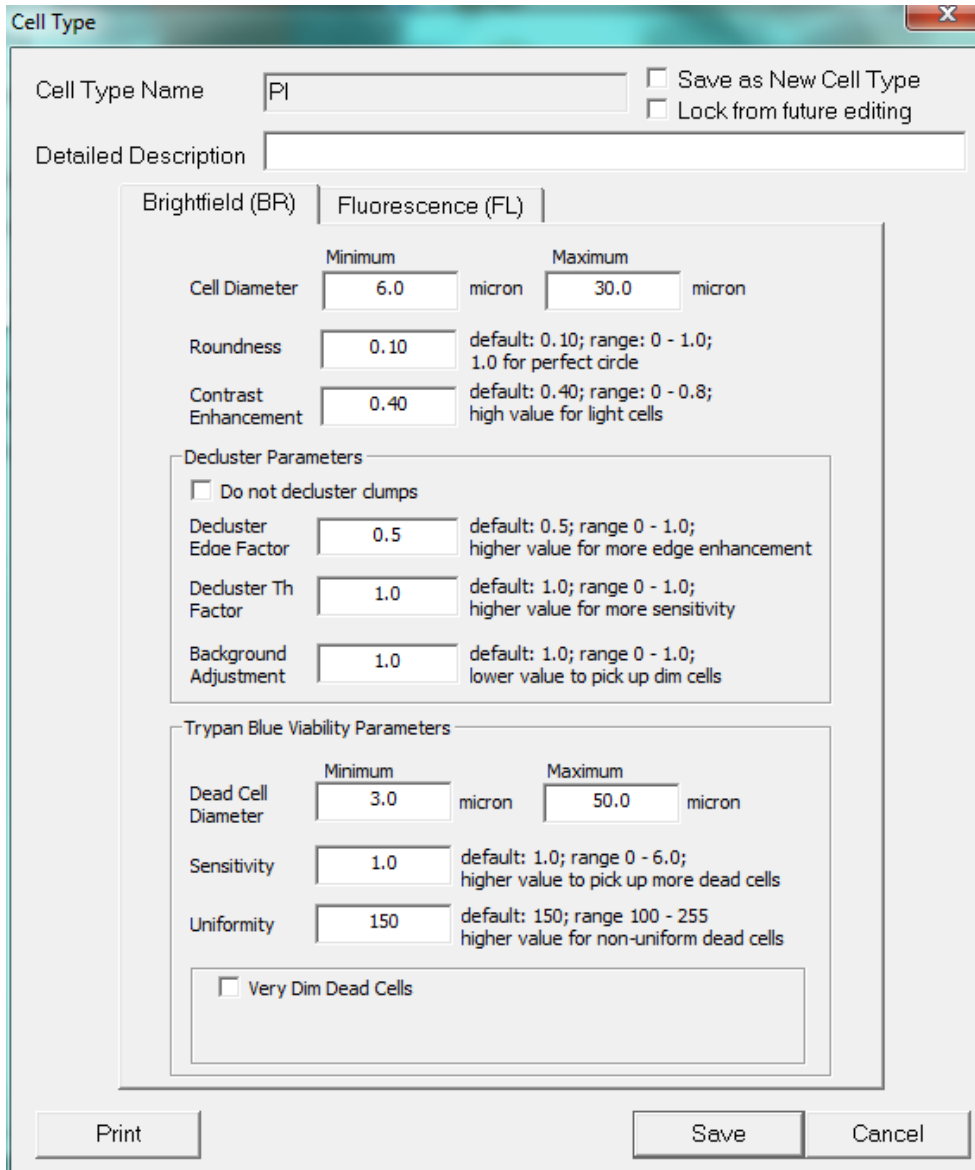
The screenshot shows the 'Cell Type' dialog box with the 'Fluorescence (FL)' tab selected. The 'Cell Type Name' is 'Calcein AM' and the 'Detailed Description' field is empty. The 'Brightfield (BR)' and 'Fluorescence (FL)' tabs are visible at the bottom. The 'Fluorescence (FL)' tab is active, showing the following settings:

- Cell Type Name:** Calcein AM
- Detailed Description:** (empty)
- Save as New Cell Type:**
- Lock from future editing:**
- Brightfield (BR) / Fluorescence (FL) tabs:** Fluorescence (FL) is selected.
- Description:** Calcein AM
- Cell Diameter:**
  - Minimum: 6.0 micron
  - Maximum: 30.0 micron
- Normalize intensity for cell size:**
- Non-uniform cell fluorescence:**
- Roundness:** 0.10 (default: 0.10 range: 0 - 1.0; 1.0 for perfect circle)
- Do not count free nuclei:**
- Advanced BR/F mode:**
- Fluorescence Threshold Parameters:**
  - Auto Threshold Fluorescent:**  (value: 10.0) \* Count range 0 - 100% of brightest cell. Lower values count dimmer cells.
  - Manual Threshold Fluorescent:**  (value: 20.0) \* Count range 0 - 100% of image max. Lower values count dimmer cells.
  - Deduster Th Factor:** 1.00 (default: 0.9; range 0 - 1.0; lower value for better deduster)
- Buttons:** Print, Save, Cancel

- Click on the pencil icon under Assay on the main Vision CBA screen, then click the Edit button on the right-hand side of the dialog box.



- Verify the Brightfield (BR) settings for the F2 image (PI).





7. Verify the fluorescent settings for the F2 image, by selecting the Fluorescence (FL) tab.

The screenshot shows the 'Cell Type' dialog box with the 'Fluorescence (FL)' tab selected. The 'Cell Type Name' is 'PI'. The 'Detailed Description' field is empty. The 'Fluorescence (FL)' tab is active, showing a 'Description' of 'Propidium Iodide (PI)'. The 'Cell Diameter' is set with a minimum of 4.0 and a maximum of 30.0 microns. There are checkboxes for 'Normalize intensity for cell size' and 'Non-uniform cell fluorescence', both of which are unchecked. The 'Roundness' is set to 0.10, with a default of 0.10 and a range of 0 to 1.0. There are checkboxes for 'Do not count free nuclei' and 'Advanced BR/F mode', both of which are unchecked. The 'Fluorescence Threshold Parameters' section has two options: 'Auto Threshold Fluorescent' (unchecked) and 'Manual Threshold Fluorescent' (checked). Both options have a threshold value of 10.0. The 'Decuster Th Factor' is set to 1.00, with a default of 0.9 and a range of 0 to 1.0. At the bottom of the dialog are 'Print', 'Save', and 'Cancel' buttons.

Cell Type

Cell Type Name   Save as New Cell Type  
 Lock from future editing

Detailed Description

Brightfield (BR) | Fluorescence (FL) |

Description

Cell Diameter Minimum  micron Maximum  micron

Normalize intensity for cell size  
 Non-uniform cell fluorescence

Roundness  default: 0.10 range: 0 - 1.0;  
1.0 for perfect circle

Do not count free nuclei  Advanced BR/F mode

Fluorescence Threshold Parameters

Auto Threshold Fluorescent  
 \* Count range 0 - 100% of brightest cell  
Lower values count dimmer cells

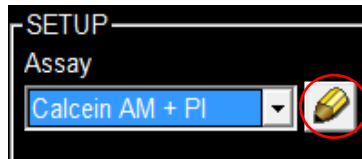
Manual Threshold Fluorescent  
 \* Count range 0 - 100% of image max  
Lower values count dimmer cells

Decuster Th Factor  default: 0.9; range 0 - 1.0;  
lower value for better decuster

Print Save Cancel

### 4.3 Software Settings for Whole Blood, Cord Blood, and Bone Marrow Clinical Samples (Calcein AM + PI CS)

1. Click on the pencil icon under Assay on the main Vision CBA screen.



2. Adjust the current settings.
  - 2.1 Check the Save as New Assay Type box
  - 2.2 Rename Assay to Calcein AM + PI CS
  - 2.3 Change the Fluorescent Exp for Calcein AM from 75.0 to 200.0
  - 2.4 Change the Fluorescent Exp for PI from 2000.0 to 4000.0
  - 2.5 Change the Dilution Factor for Assay from 1.25 to 10

## 2.6 Click Save

Dialog

Assay Name:   Special Cells

Save as New Assay Type  Lock Assay from future editing

Description:

Imaging Mode:   Acquire Brightfield Image  
 Two Chamber Assay  Multimode FL Counting

**F1 Image**

Cell Type:

Description:

Fluorophore:

Fluorescent Exp:  msec

Use Br Exp Factor of

Remove FL Pos from BR count

Show Data File Buttons

Show Sample Adjustment Button

Show Cell Size Distribution Button

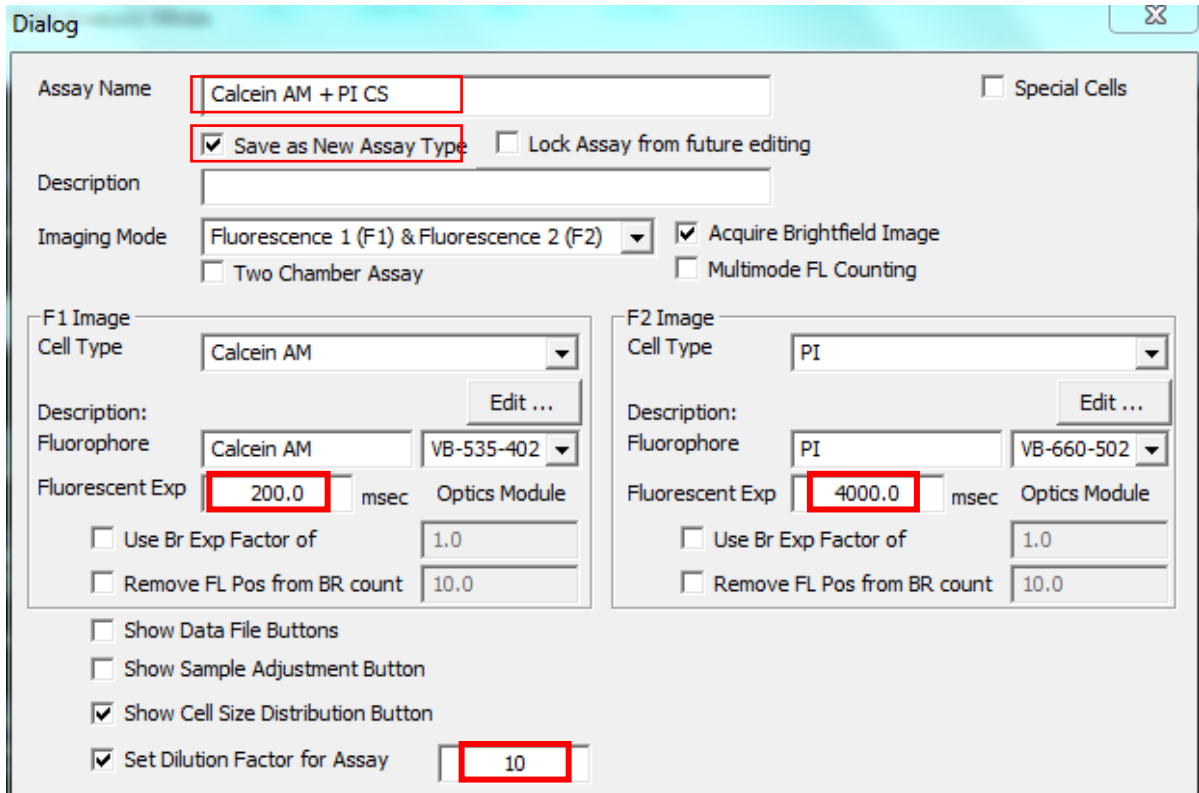
Set Dilution Factor for Assay

Show Percent F1,F2   Use Custom Label

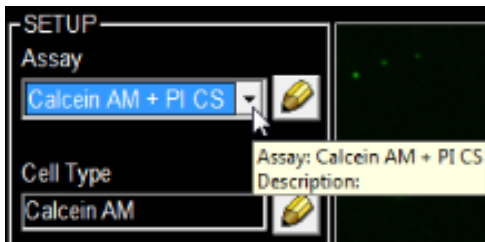
Result Template: Calcein AM and PI.rlt\_tm

Print Template: Calcein AM and PI.rlt\_tm

FCS Layout File: <None Selected>

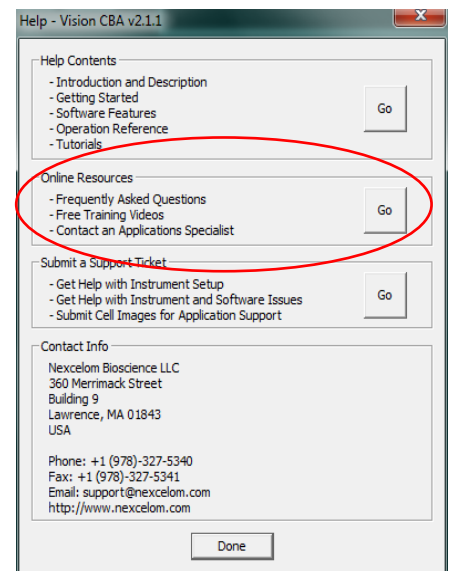


3. The updated setting should match those shown below.
4. You may now select the Calcein AM + PI CS assay from the drop down menu and proceed with data acquisition in section 2.5.



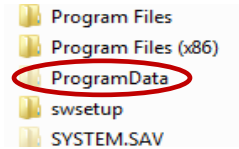
## 5.0 Importing a New Cell Type, Assay Type, and Templates

1. Click the question mark in the right-hand corner to access the help menu.
2. Click on the "Go" button in the "Online Resources" section. This will automatically load the Cellometer Vision CBA Online Resources webpage.
3. Under the "Assay Files" tab, locate and download the Calcein AM + PI files onto the desktop.

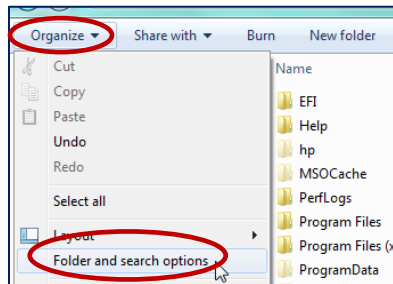


## 5.1 Import Result and Print templates

1. Navigate to the START menu and select Computer.
2. Double click on the C: Drive and locate the ProgramData folder (shown at right), then proceed to Step 3. IF the ProgramData folder is not present, it may be hidden. Follow the instructions below to show hidden folders.



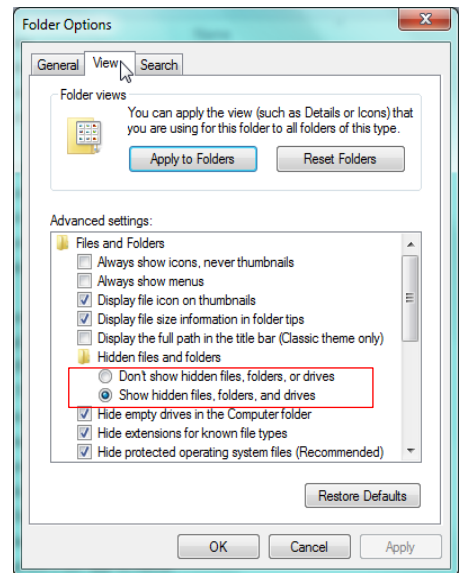
- 2.1 Click on “Organize” (top left of screen) and select “Folder and search options”



- 2.2 A Folder Options menu will pop up. Select View.

- 2.3 Under the “Hidden files and folders” file, select “Show hidden files, folders, and drives”

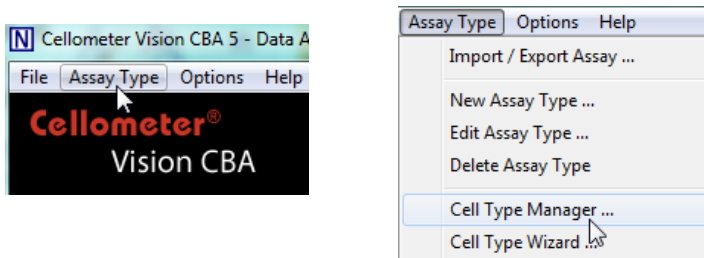
- 2.4 Click OK.



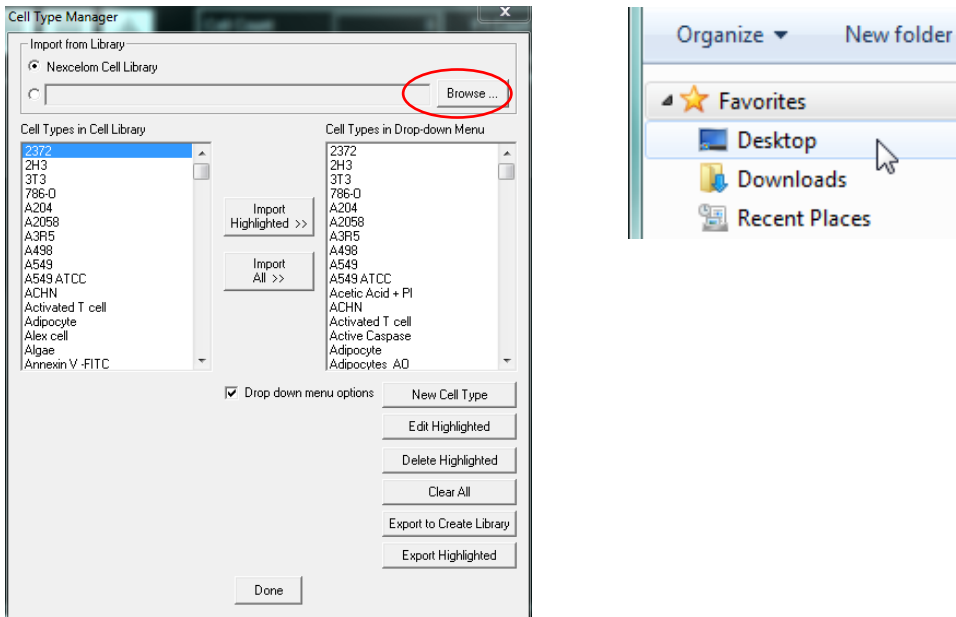
3. Open ProgramData folder, then open the Nexcelom\_VisionCBA folder.
4. Open the Template folder. Copy the result template (Calcein AM and PI.rlt\_tm) and print template (Calcein AM and PI .prn\_tm) files from the new folder on the desktop and paste them into the Template folder.

## 5.2 Import Cell Type

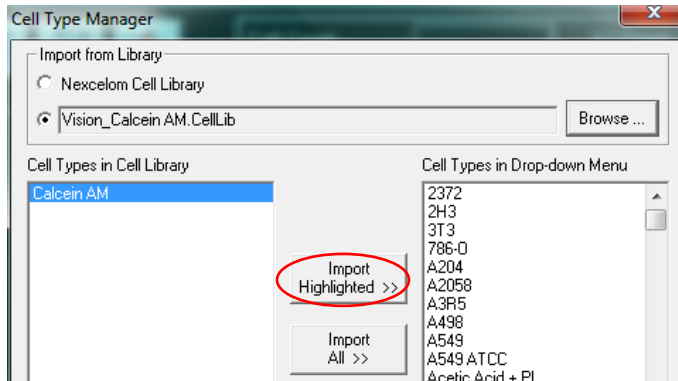
1. Locate and click on the “Assay Type” at the top of the screen. Followed by Cell Type Manager.



- Once the Cell Type Manager appears, click on Browse, navigate to the desktop and locate the “Vision\_Calcein AM.CellLib” file that has been downloaded onto your desktop.



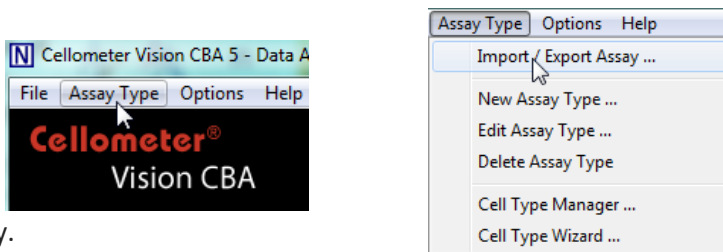
- Select Vision\_Calcein AM.CellLib on the desktop, and click Import Highlighted.



- Repeat step 1-4 to download “Vision\_PI.CellLib”

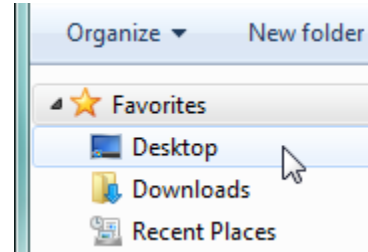
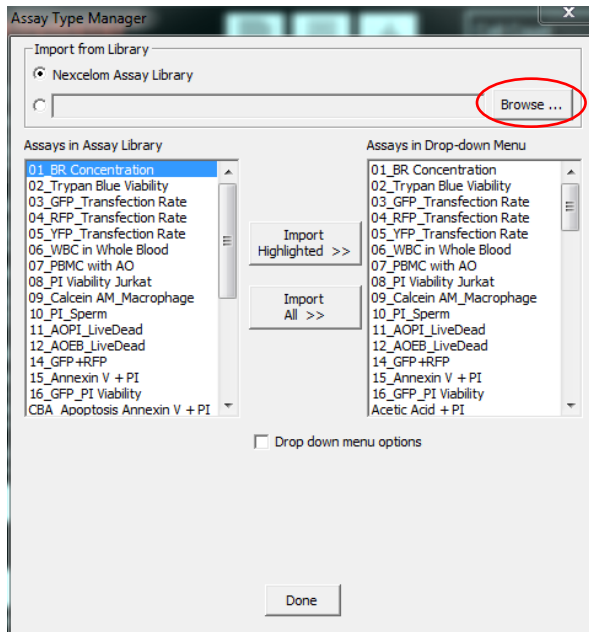
### 5.3 Import Assay Type

- Locate and click on the “Assay Type” at the top of the screen. Followed by Import / Export

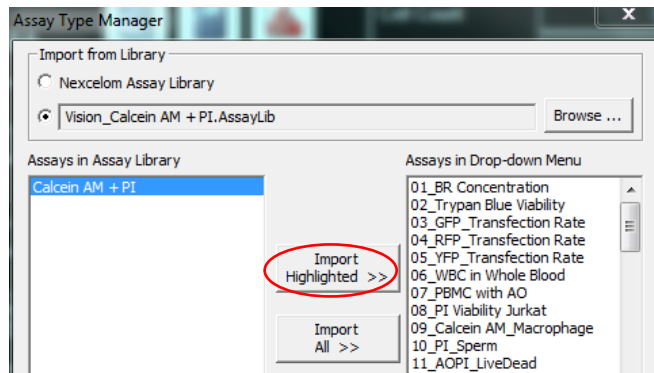


Assay.

- Once the Assay Type Manager appears, click on Browse, navigate to the desktop and locate the “Vision\_Calcein AM.CellLib” file that has been downloaded onto your desktop.



- Select “Vision\_Calcein AM + PI.AssayLib” on the desktop, and click Import Highlighted.



- This assay is now available for use. See section 2.3.4 to select and use assay.

## 4.0 Additional Resources

### 4.1 Technical Support

- Nexcelom Technical Support is available from 9am to 5:00pm EST.  
E-mail: [support@nexcelom.com](mailto:support@nexcelom.com)  
Phone: 978-327-5340

### 4.3 Storage and Handling

- For long term storage, store CSK-0118 Component A between 2°C and 8°C .
- For long term storage, store CSK-0118 Component B between -24°C and -16°C.

### 4.4 Warranty

This product is for RESEARCH USE ONLY and is not approved for diagnostic or therapeutic use. Product is warranted to meet the specifications outlined in the Certificate of Analysis when stored and used according to the manufacturer's instructions. No other warranty, expressed or implied (such as merchantability, fitness for a particular purpose, or non-infringement) is granted. Warranty is valid until the expiration date stated on the product label. If no expiration is listed, the warranty is valid for 6 months from the date of product receipt. Warranty will be void if product is stored incorrectly, the recommended protocol is not followed, or the product is used for a different application.

## 5.0 Ordering Information

### 5.1 How to Reorder

#### **For orders shipping to destinations in the United States:**

- When ordering with a Purchase order
  - Fax a copy of your order to 978-327-5341
  - Email a copy of your order to [sales@nexcelom.com](mailto:sales@nexcelom.com)
- When ordering with a Credit Card
  - Visit [www.shop.nexcelom.com](http://www.shop.nexcelom.com) and place your order

#### **For orders shipping to destinations outside the United States:**

Contact your local distributor or Nexcelom Representative to place your order