

MetaMatters

NEW in MetaMorph® Software version 7.7.1...

Support for:

- The new Hamamatsu ORCA®-Flash 2.8 camera with the Scientific CMOS Image Sensor FL-280!
- The Olympus Soft Imaging Solutions IX2-FCB controller
- The Hamamatsu C9300-221 high speed VGA interline CCD camera
- The Zeiss Definite Focus z-drift compensator
- The Sutter Lambda XL extended life illumination source

Improved support for:

- Hamamatsu C9100-13 high dynamic range EMCCD camera

To download this update, you must already be at version 7.x and have a current Software Maintenance Plan. Go [HERE](#) to initiate download.

For MetaMorph Software support, contact our Technical Support at 800-635-5577 x1820 or email us at support.dtn@moldev.com

For more information, contact info@moldev.com

MetaMorph Software Webinar Series: "Performing FRET with MetaMorph Software"

August 10, 2010

10:30 AM, EDT

Recording will be posted to our website for those who cannot attend live.

Look for your invitation in your email in-box!

Inside this issue:

MetaMorph Software Incentives	2
MetaMorph Software News	1, 4
FOCUS: Quantifying Cell Motility Using Tracking Applications	2, 3
Upcoming Events	4

MetaTool Tips:

Update vs. Upgrade

- An update is available to accounts with current software maintenance and is only good within a single version (i.e. 7.1 » 7.7).
- An upgrade is available to all customers who wish to go from an older version to a newer version (i.e. 6.x » 7.x).

Amazing Software Incentives!

Submit a 500 - 800 word
MetaMorph® Software method and
receive a FREE software upgrade or
application module.

Submit a 100 - 200 word
MetaMorph® Software tip
and receive a FREE 12 month software
maintenance agreement.

The method description and tip paragraph will be
published in *MetaMatters*.

*The free MetaMorph®
Software Basics
Training Course is
September 21 & 22,
2010!*

*Email:
Mary.David@moldev.com
for more information on
incentive programs.
For training courses click here:
[MetaMorph Software Training](#)*

*The MetaMorph®
Software Advanced
Training Course is
September 23 & 24,
2010!*

FOCUS: Quantifying Cell Motility Using Tracking Applications

*George McNamara, Ph.D., Image Core Manager
Analytical Imaging Core Facility, University of Miami*

In part one, I described making a panoramic movie of the classic timelapse film by David Rogers of a neutrophil chasing bacterium. The movie was acquired using phase contrast microscopy – a great method for visualizing transparent cells but a nightmare for fully automated tracking. I added the “human touch” by using the Trace Region tool and Display – Graphics - Paint Region, intensity level 255, to paint the neutrophils white in every plane. Prior to painting, I first used Process – Arithmetic; subtract constant 10, to make sure only the pixels I painted would be intensity level 255. Here, I will compare different tracking applications in MetaMorph Software:

1. Track Points (manual)
2. Track Objects (using Threshold algorithm)
3. MultiDimensional Motion Analysis (using either of two algorithms)
4. Integrated Morphometry Analysis (IMA) / Morphometry - Measure Objects (MO)

I used the original panoramic stack with Track Points, after Log – Open Data Log to enable saving to Microsoft Excel 2007 by Dynamic Data Exchange. I attempted to track the “center of mass” (centroid in image analysis terms). *Figure 1* shows the three planes from the original data, overlaid with Track points performed three times. The overlay was made by using Process – Arithmetic, maximum, with 24-bit result between bitmap image saved from Track Points – Graph Data and an image made by copying/pasting the three planes with regions of interest. If I did perfectly reproducible tracking, only one color cross would be visible. The positions are not identical and in some areas, wanders more than others.

(continued on page 3)

FOCUS: MetaMorph® Software Method - Making a Panoramic Movie from a 16 mm film

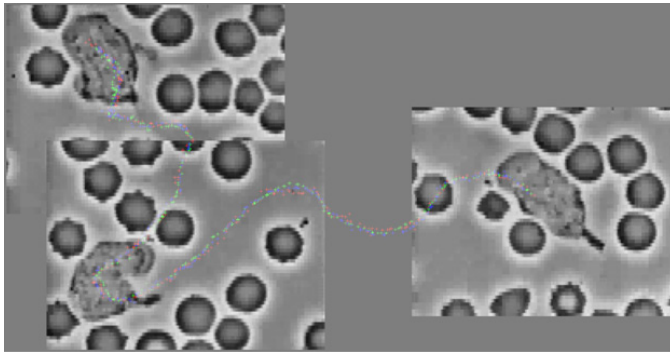


Figure 1. First (#1), middle (#220) and last (#439) planes merged into composite image, and overlaid with Track Points positions manually tracked three times (red, green, blue).

MultiDimensional Motion Analysis has several algorithm options. The Threshold Result option and Template Match option use the Track Objects commands. The Auto-Find Objects algorithm, with simple threshold segmentation method, works fine with threshold (intensity 255) painted objects; though it is somewhat slower than Threshold Result. Auto-Find using the current IMA (Integrated Morphometry Analysis) morphometry setting is faster than simple threshold; though fine tuning settings is hardly needed if you have high contrast objects such as region painted cells as here.

The three Apps commands discussed above provide identical data display and graphing tools. If you do not need these tools, or do not have the trackobj drop-in, you can track using the morphometry commands. The Measure - Integrated Morphometry Analysis (IMA) and the Morphometry - Measure Objects (MO) are two interfaces to the same engine. When tracking objects with IMA/MO, I recommend having only one object per tracking cycle. If you are starting with more than one object present, you can often use Process - Binary Operations - Binarize, to make a binary stack, then Process - Stack Arithmetic, Maximum, to make a binary path image. If you need to set up to have only one object/track at one time, you can use a region of interest with Paint Region, intensity level 0, to black out all but one path or part of one path (Figure 3). You can then use the Process - Arithmetic, Logical And (stack all planes, binary path mask), to blackout everything but the object of immediate interest.

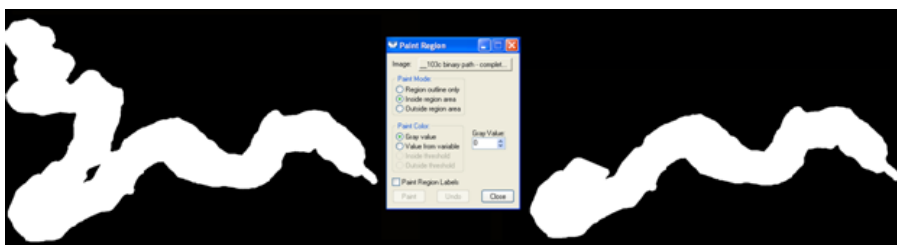


Figure 3. Complete binary path (left), Paint Region dialog box, and path truncated by blacking out part.

(continued from page 2)

For Track Objects, I used Search Options-Algorithm = Threshold, with the tracking regions (40x40, 60x60 pixels) entirely inside the thresholded object in plane 1. Figure 2 shows the path generated by Track Objects. The path is smoother than the manual Track Points method. Only one track is shown since multiple runs will generate the same result.

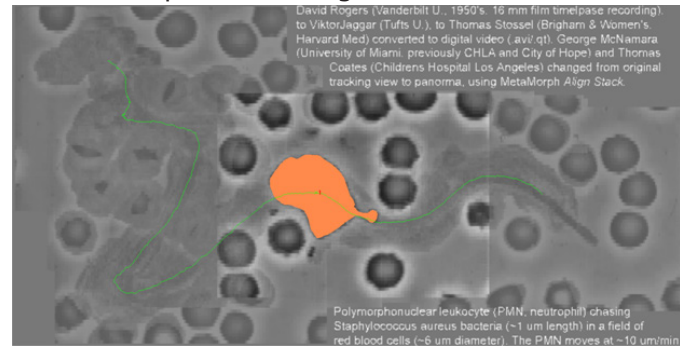


Figure 2. Panoramic movie plane #110 with threshold (orange) on for intensity 255 of the neutrophil; green trace is the Track Objects path, green X is the current position, and red "1" is Track Objects designating this as track 1.

Once you have your one object stack, look closely at the IMA Measurements tab to make sure your parameter limits are usable. You may also want to enable more (or fewer!) parameters, by using the Select Measurements button.

In part 3, I will discuss IMA/MO in more detail. Here I just summarize the key points:

1. Pick your parameters.
2. Pay attention to Preferences.
3. Test the parameters make sense.
4. You will need to record IMA or MO into a journal in order to "loop for all planes".
5. Open log files to separate Excel worksheet tabs (or text files).



Improved throughput is just a click away!

Is your MetaMorph® acquisition system overwhelmed with users?

Are users waiting in line to analyze their data?

Adding a dedicated MetaMorph offline analysis system provides improved productivity to your laboratory or core facility.

Accelerate throughput and save time by moving image analysis off the acquisition system and onto the separate dedicated workstation. This is especially helpful for analysis that requires high computational power such as deconvolution or performing segmentation and measurements on large data sets.

We offer three offline products:

- MetaMorph Offline Basic Software: Includes standard measurement and analysis capabilities such as thresholding, segmentation, filtering and morphometry.
- MetaMorph Offline Premier Software: Enhanced version of Basic Software plus advanced application modules and the 4D viewer application.
- MetaMorph Offline Premier Software Network Licensing: Equivalent to offline Premier version but in a cost effective bundle with attractive per seat license pricing.

These MetaMorph packages can be purchased through any of our valued distribution partners or directly from Molecular Devices, Inc. Click [HERE](#) to find a distributor near you.

402 Boot Road
Downingtown, PA 19335

Phone: 800-635-5577
Fax: 610-873-5499
meta.admin@moldev.com
support.dtn@moldev.com
training.dtn@moldev.com

**MetaMorph® Software...
the gold standard in
research imaging**

**We're on the web!
MetaMorph.com**

Upcoming Training, Courses and Conferences

AUGUST 10, 2010

Performing FRET within
MetaMorph Software
WEBINAR @ 10:30 AM, EDT

SEPTEMBER 21 – 22, 2010

Fundamentals of
MetaMorph Software
Downingtown, PA

SEPTEMBER 23 – 24, 2010

Advanced Topics in
MetaMorph Software
Downingtown, PA

OCTOBER 13 – 26, 2010

Immunocytochemistry, In Situ
& Live Cell Imaging
Cold Spring Harbor

NOVEMBER 13 – 17, 2010

Society for Neuroscience
San Diego, CA

DECEMBER 11 – 15, 2010

American Society for Cell
Biology
Philadelphia, PA



MetaMorph® products @ QFM,
2008