



November 16th, 2022

Subject: Discontinuation of MetaMorph Microscopy Automation and Image Analysis Software Suite

Dear Valued MetaMorph Customer,

For 30 years, MetaMorph Microscopy Automation and Image Analysis Software, including MetaFluor and MetaVue, has enabled Scientists like you to make landmark discoveries in biology. Over 25,000 scholarly papers cite MetaMorph Software, exemplifying the learning that it has helped unlock. We thank you for being our customer and for your contributions to science and advancing discovery!

After careful consideration, we have decided to cease offering the MetaMorph Software Suite, including MetaFluor and MetaVue. This decision aligns with our strategy of providing new innovative products with improved security, design, and functionality in support of scientific research. The intent of this letter is to serve as an Advanced Notice of Discontinuation to help you plan accordingly.

## Important dates for purchase and support:

- <u>Last day to purchase MetaMorph Software is June 30, 2023</u>. After this date, it will no longer be available for purchase.
- Last day to purchase a maintenance plan for MetaMorph Software is June 30, 2023. All maintenance plan coverage will end by December 29, 2023. After this date, all service and support will be on a reasonable effort basis. The software maintenance plan provides access to product support and access to software updates. To check your expiration, please see <a href="http://www.meta.moleculardevices.com/maintenance/">http://www.meta.moleculardevices.com/maintenance/</a>

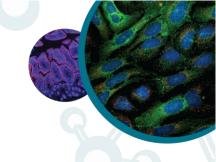
The strength of MetaMorph lives on in our leading microscopy instrumentation and we encourage you to investigate what we have to offer. Our systems for high-content imaging (HCI) and high-content analysis (HCA) provide flexible scalability making it easy to evolve your system alongside your research. If you know MetaMorph you will easily be able to grow into the high-content image acquisition and analysis software MetaXpress designed for a range of applications optimized for ImageXpress instruments.

Additionally, IN Carta image analysis software solves the challenges of complex image analysis by providing easy-to-use workflows and machine learning to generate insights quickly from ImageXpress imaging systems. And if you are looking for an easy to learn and use system for walk-up microscopy, please consider our ImageXpress® Pico Automated Cell Imaging System.

ImageXpress Confocal HT.ai High-Content Imaging System: <a href="https://www.moleculardevices.com/products/cellular-imaging-systems/high-content-imaging/imagexpress-confocal-ht-ai">https://www.moleculardevices.com/products/cellular-imaging-systems/high-content-imaging/imagexpress-confocal-ht-ai</a>

MetaXpress High-Content Image Acquisition and Analysis Software: <a href="https://www.moleculardevices.com/products/cellular-imaging-systems/acquisition-and-analysis-software/metaxpress#Overview">https://www.moleculardevices.com/products/cellular-imaging-systems/acquisition-and-analysis-software/metaxpress#Overview</a>

IN Carta Image Analysis Software: <a href="https://www.moleculardevices.com/products/cellular-imaging-systems/acquisition-and-analysis-software/in-carta-image-analysis-software">https://www.moleculardevices.com/products/cellular-imaging-systems/acquisition-and-analysis-software/in-carta-image-analysis-software</a>





ImageXpress® Pico Automated Cell Imaging System: <a href="https://www.moleculardevices.com/products/cellular-imaging-systems/high-content-imaging/imagexpress-pico">https://www.moleculardevices.com/products/cellular-imaging-systems/high-content-imaging/imagexpress-pico</a>

We encourage you to contact our team, using the information provided below. This is a good time to ensure that all your MetaMorph systems are fully up to date with the latest version of the software, which is 7.10.5 (released in 2022) – please provide your MetaMorph License ID(s) so our team can help you with any update requests.

For more information, please contact our team: <a href="mailto:support.dtn@moldev.com">support.dtn@moldev.com</a>

As always, we remain committed to providing you with the highest quality of products and services. Thank you for the opportunity to continue to support and help advance your research and discovery.

Please let us know if we can answer any questions.

Sincerely,

KD Modi

Director of Product Management