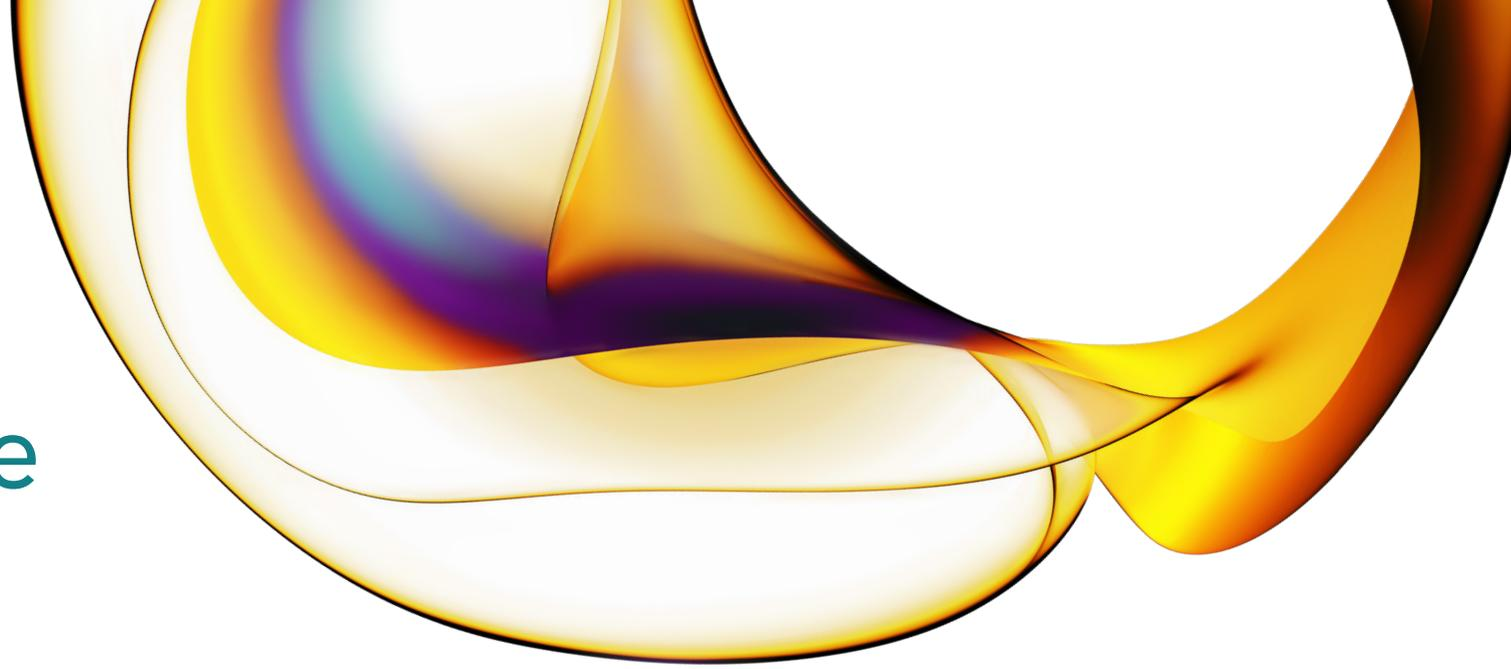


revvity

Next-generation image data management.



Signals Image Artist: Giving you more options to explore your data and gain new insight

Storing, managing, and sharing large volumes of image data can be challenging, but with Signals Image Artist™ image analysis and management platform, it's easier than ever.

All your image data from the Celigo® image cytometer and other major cellular imaging systems is stored, accessed, and managed in one place. With its powerful image processing capabilities and ready-made image analysis building blocks, the Signals Image Artist platform can be used alongside the Celigo system's acquisition, visualization, and analysis software.

Benefit from machine learning and AI capabilities to train the software to develop analysis algorithms – giving you answers faster. You can even access data and perform analysis remotely via a web browser.





Benefits of Signals Image Artist:

- **Compatible with Celigo** image cytometer and major cell imaging systems
- **A central location** to store and access image data from the lab, office, or remotely
- **Advanced image analysis** capabilities, including machine learning and AI
- **Multiuser solution** that can support your entire lab without compromising performance
- **Scalable data storage** to expand with your labs needs over time
- **Easy IT management** with single system and installation across image data types
- **Cloud and on-premise** options to choose from
- **Web browser application** to access data and perform analysis remotely

The Signals Image Artist platform and the Celigo image cytometer are sold separately.

revvity