

# IncuCyte® ZOOM Software Version Feature Comparison

	2013B	2014A	2015A	2016A	2016B	2018A
<b>Data Storage</b>						
Data Archiving	●	●	●	●	●	●
IncuStore™ Expansion Data Storage Compatibility	●	●	●	●	●	●
<b>Whole Well Imaging</b>						
Points of Interest (POIs) - Digital Monitoring Tools	●	●	●	●	●	●
Automated vessel marking - IncuCyte® ZOOM Marking Tool	●	●	●	●	●	●
35mm Dishes, 6-, 12-, 24-, 96-Well Enabled	●	●	●	●	●	●
384-Well Enabled			●	●	●	●
<b>General Features</b>						
New Movie Formats (MPEG-4, AVI uncompressed)	●	●	●	●	●	●
Improved Scheduling Flexibility			●	●	●	●
Fluorescence Analysis Tools for Co-culture Scratch Wound Assays			●	●	●	●
Enhanced NeuroTrack™ Fluorescence Tools			●	●	●	●
Enhanced autofocus & additional plate support for single spheroid applications in round bottom ULA plates					●	●
<b>Data Processing</b>						
Background Fluorescence Subtraction		●	●	●	●	●
Fluorescent NeuroTrack™ Analysis		●	●	●	●	●
Import/Export Image Processing Definitions		●	●	●	●	●
Coincident Object Analysis				●	●	●
Archive Analysis						●
<b>IncuCyte® ZOOM Rack System</b>						
Space Saving Rack - Holds 2 instruments and 2 IncuStore™ Units		●	●	●	●	●
Uninterruptable Power Supply		●	●	●	●	●
'In The Lab' Access to User Interface		●	●	●	●	●
<b>Enabled Applications</b>						
Angiogenesis (additional software module required)	●	●	●	●	●	●
Apoptosis	●	●	●	●	●	●
Cell Culture QC	●	●	●	●	●	●
Chemotaxis (additional software module required)			●	●	●	●
Clonal Dilution	●	●	●	●	●	●
Cytotoxicity	●	●	●	●	●	●
Immune Cell Killing	●	●	●	●	●	●
Immune Cell Proliferation & Clustering	●	●	●	●	●	●
Neurite Dynamics - Label-free (additional software module required)	●	●	●	●	●	●
Neuronal Co-Culture - Fluorescence (additional software module required)		●	●	●	●	●
Proliferation - Cell Count	●	●	●	●	●	●
Proliferation - Confluence	●	●	●	●	●	●
Reporter Gene	●	●	●	●	●	●
Scratch Wound Migration (additional software module required)	●	●	●	●	●	●
Spheroids	●	●	●	●	●	●
Stem Cell Monitoring & Reprogramming	●	●	●	●	●	●
Transfection Efficiency	●	●	●	●	●	●

For additional information on this and other products, please contact us at [sales@essenbio.com](mailto:sales@essenbio.com)