Dilution cloning using IncuCyte[®] ZOOM System **Real-time verification of clonality**

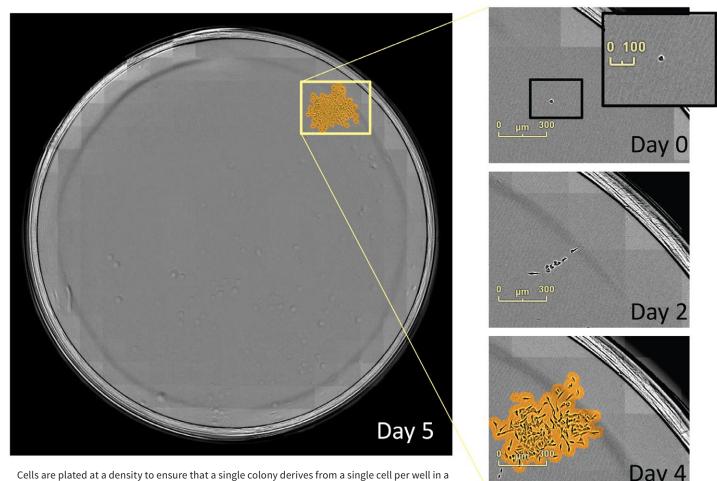


Spend more time inspecting and analyzing at your desk while your cells spend more time in the incubator.

Inspect colony morphology, monitor growth characteristics and validate clonality in real time while cells remain undisturbed in the incubator.

Pinpoint rare events to track and monitor morphology and growth rate over time.

Analyze data on-the-fly simple and easily share them to facilitate collaboration with multiple user access.



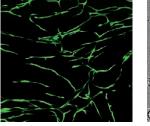
Cells are plated at a density to ensure that a single colony derives from a single cell per well in a 96-well plate. HD-phase contrast images of the whole well are acquired daily.



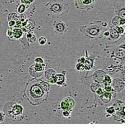
Key applications using the IncuCyte® ZOOM Live-Cell Analysis System



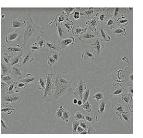
3D-Spheroids



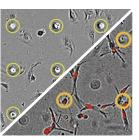
Angiogenesis



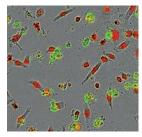
Apoptosis



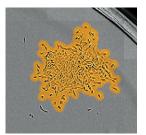
Cell Culture QC



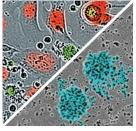
Chemotaxis Migration & Invasion



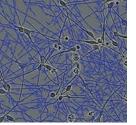
Cytotoxicity



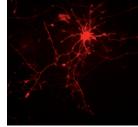
Dilution Cloning



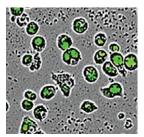
Immune Cell Killing, Clustering & Proliferation



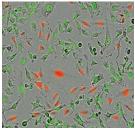
Neurite Dynamics -Label-Free



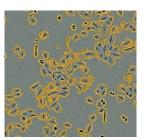
Neuronal Co-Culture -Fluorescence



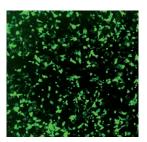
Phagocytosis



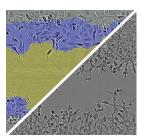
Proliferation -Cell Count



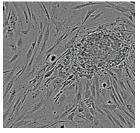
Proliferation -Confluence



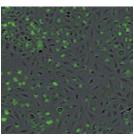
Reporter Gene



Scratch Wound Migration & Invasion



Stem Cell Monitoring & Reprogramming



Transfection Efficiency

Learn more at **essenbioscience.com/applications**. Or contact us at **sales@essenbio.com**.

Essen BioScience, 300 West Morgan Road, Ann Arbor, Michigan, USA 48108

© 2016 Essen BioScience. All rights reserved. All trademarks are the property of Essen BioScience unless otherwise specified.

