

Cell Monitoring and QC using IncuCyte® ZOOM System

Real-time monitoring of cell morphology and growth



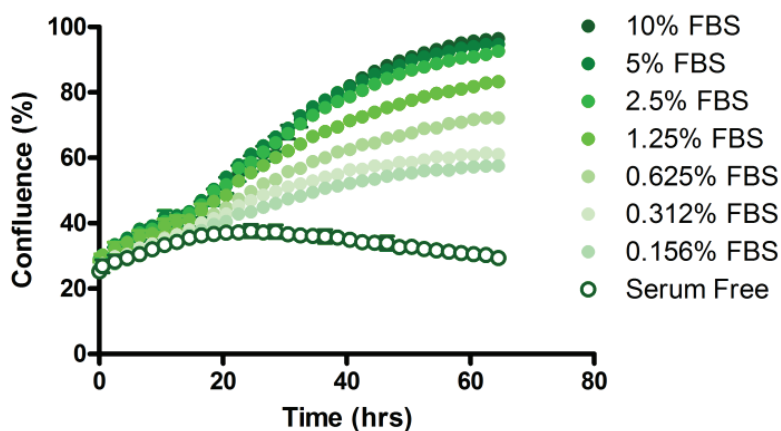
Minimize variation, improve consistency and predictability via real-time cell monitoring over time.

Assess health and phenotype of cells by monitoring cell morphology, growth properties and differentiation state in their native environment.

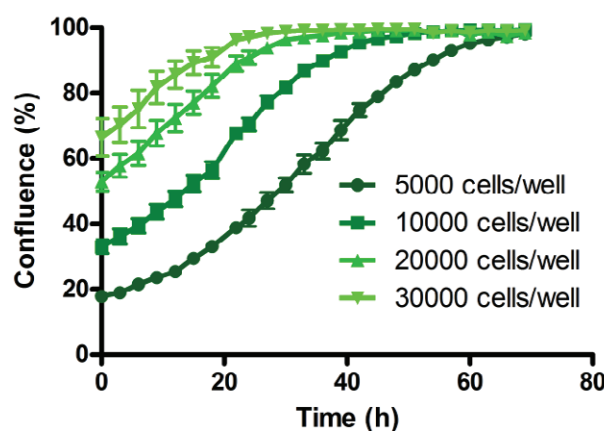
Assess yield and consistency by monitoring cell numbers and plating density via non-invasive and non-perturbing approach.

Enable data-driven decisions while the experiment is still in progress.

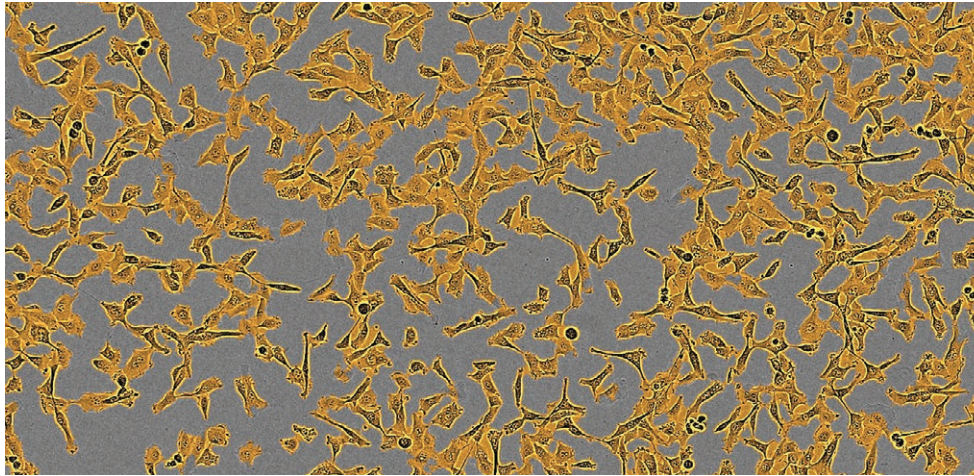
Access data remotely at your convenience.



HUVECs were cultured in basal media containing several growth factors and supplemented with decreasing amounts of FBS. The optimal concentration of FBS for normal propagation of HUVEC was determined to be >2.5%.

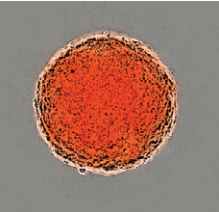
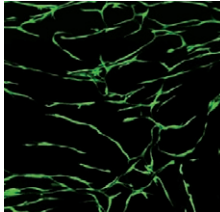
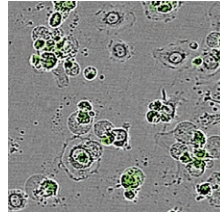
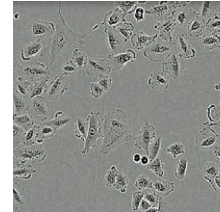
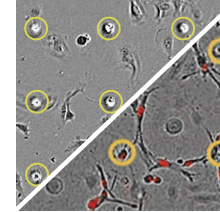
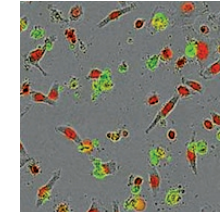
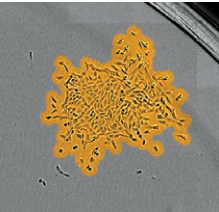
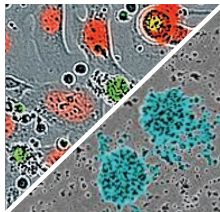
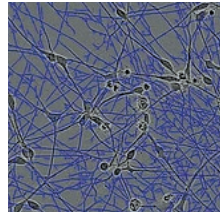
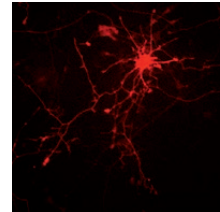
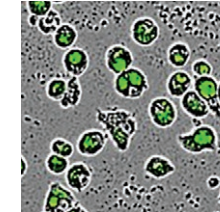
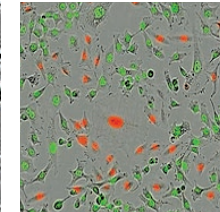
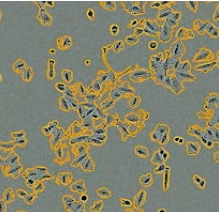
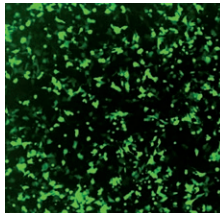
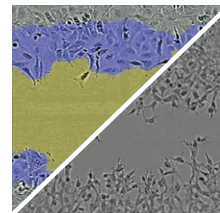
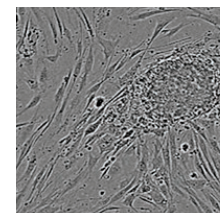
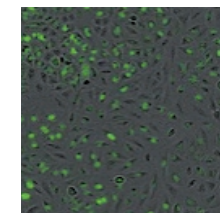


HT-1080 cells, plated at the indicated cell densities, were imaged every 3 hours for a total of 3 days. The IncuCyte® ZOOM Basic Analyzer software was used to obtain real-time updates on cell confluence over time.



The IncuCyte® ZOOM Basic Analyzer provides real-time updates on cell confluence, based on segmentation of HD-phase contrast images (segmentation mask illustrated in yellow).

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.