

Abstract:

Being fundamental and ubiquitous features of proper cellular function, cytoskeletal rearrangement and cell cycle progression are sensitive indicators of cell growth and cell stress. Previously, we developed a number of multiplexed, quantitative, cell-based, high-content screening (HCS) assays to monitor these key cellular functions using a Thermo Scientific Cellomics ArrayScan® HCS Reader. For example, the Thermo Scientific Cellomics Cytoskeletal Rearrangement HCS Kit simultaneously detects changes in cytoskeletal structures (F-actin and microtubule fibers), nuclear DNA content, and nuclear and cellular morphologies using bright fluorescent probes. The Thermo Scientific Cellomics Cell Cycle I HCS Kit characterizes cell cycle progression by simultaneously quantifying DNA content, BrdU incorporation and Histone H3 phosphorylation. Here we demonstrate the value of using these two multiparameter kits together. By assessing the cell cycle using BrdU and phospho-Histone H3 antibodies while simultaneously monitoring cytoskeletal rearrangements and cell morphological changes, we established a multiparameter profile of different drugs affecting cell cycle, cell growth, cell differentiation and cell death. A variety of cells were treated with different actin or tubulin affecting-drugs or cell cycle inhibitors, and then monitored the drug effect on cellular and nuclear shape, cytoskeletal structure, and cell cycle phases. Evaluation of drugs using these multiplexed cytoskeletal and cell cycle assays is a simple and easy-to-implement method for compound toxicity profiling in high content screening.

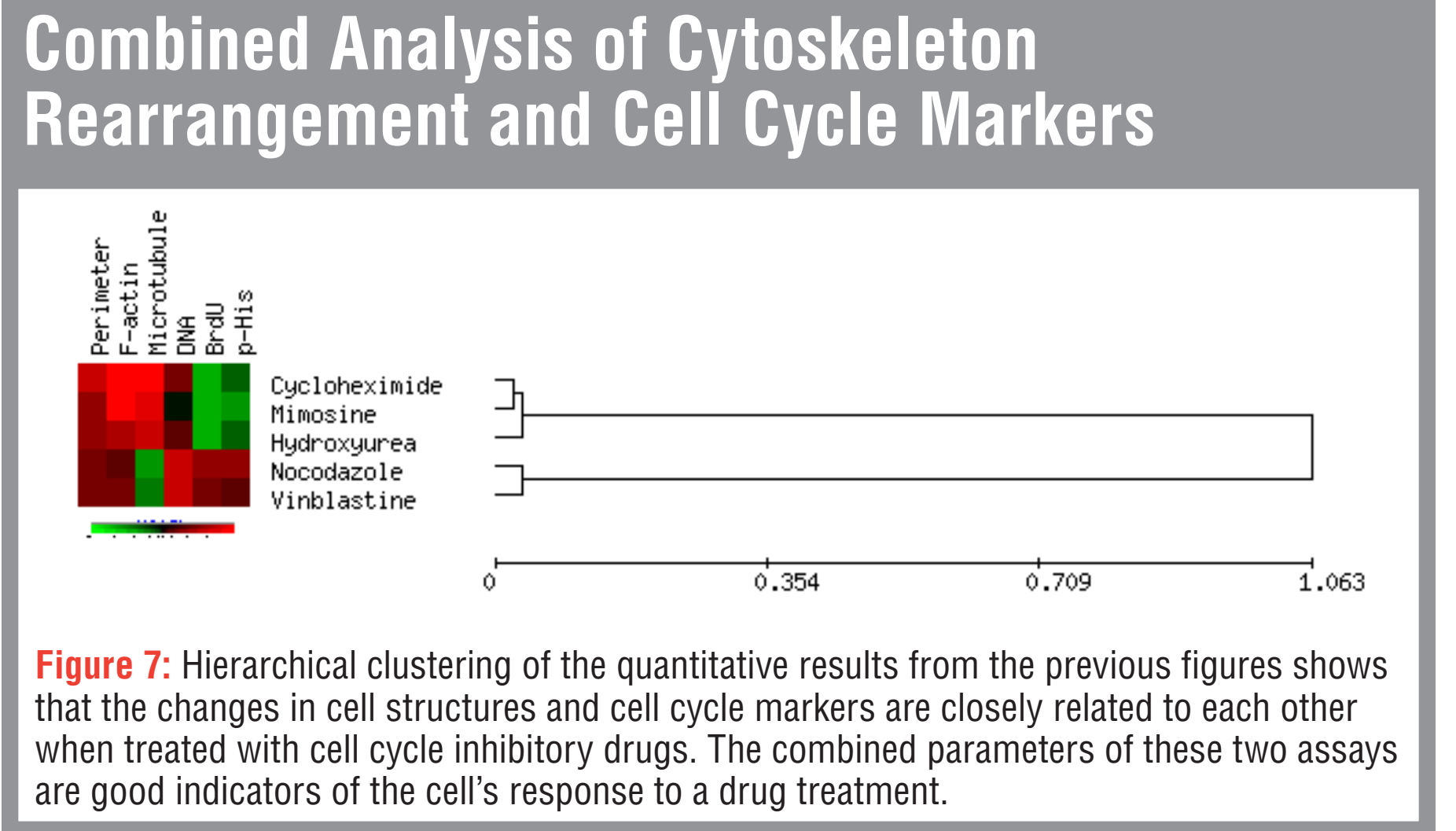
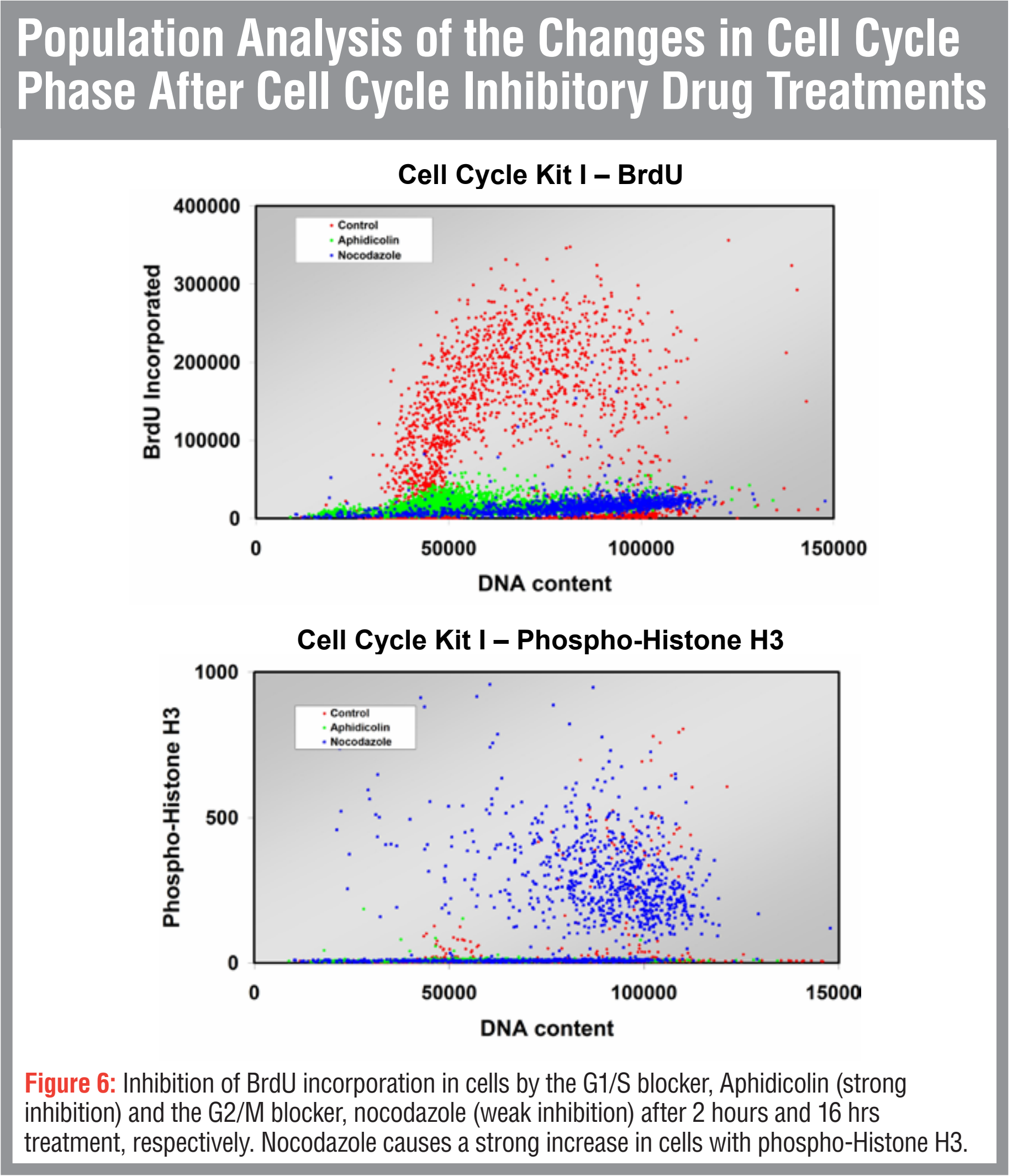
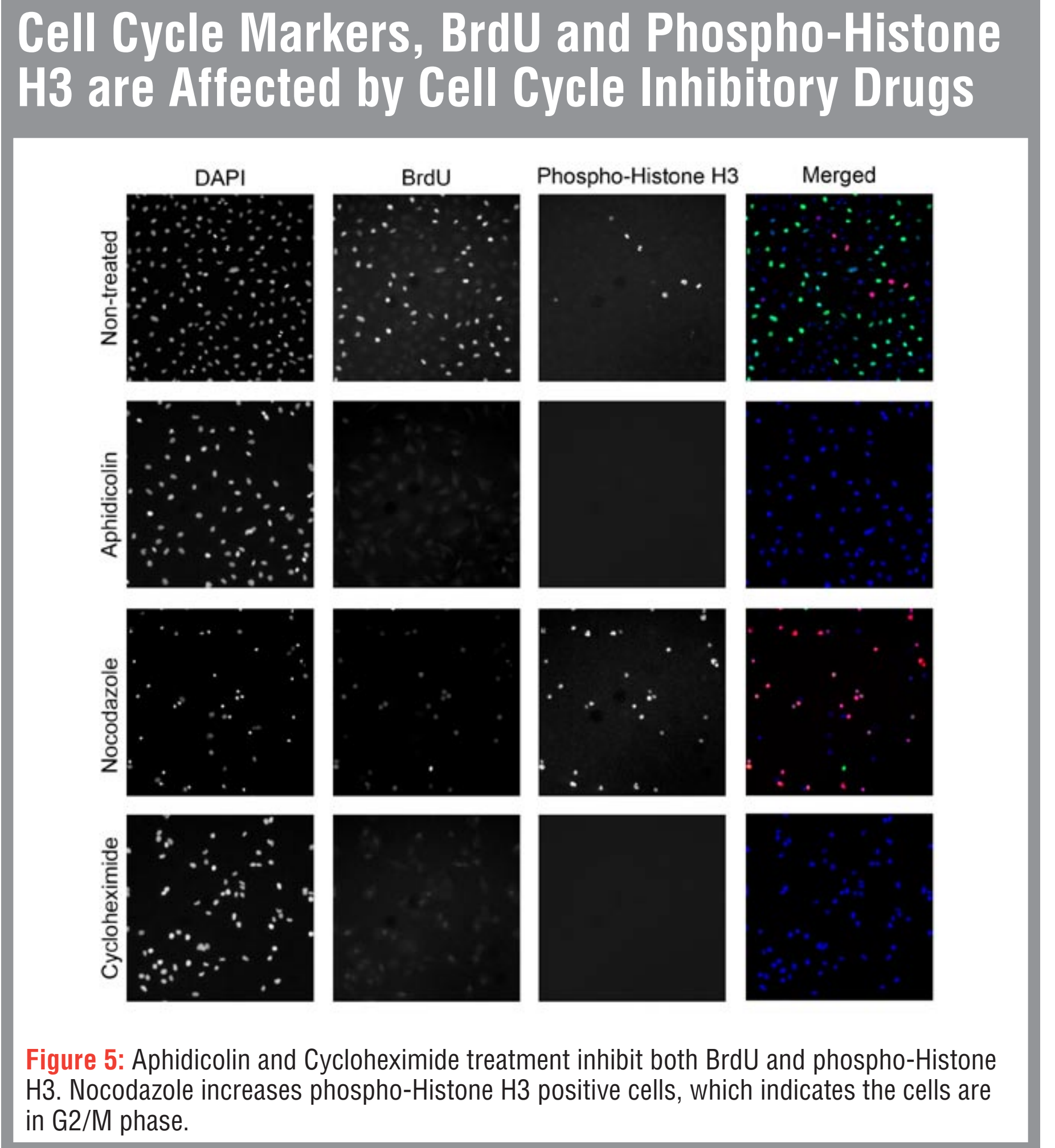
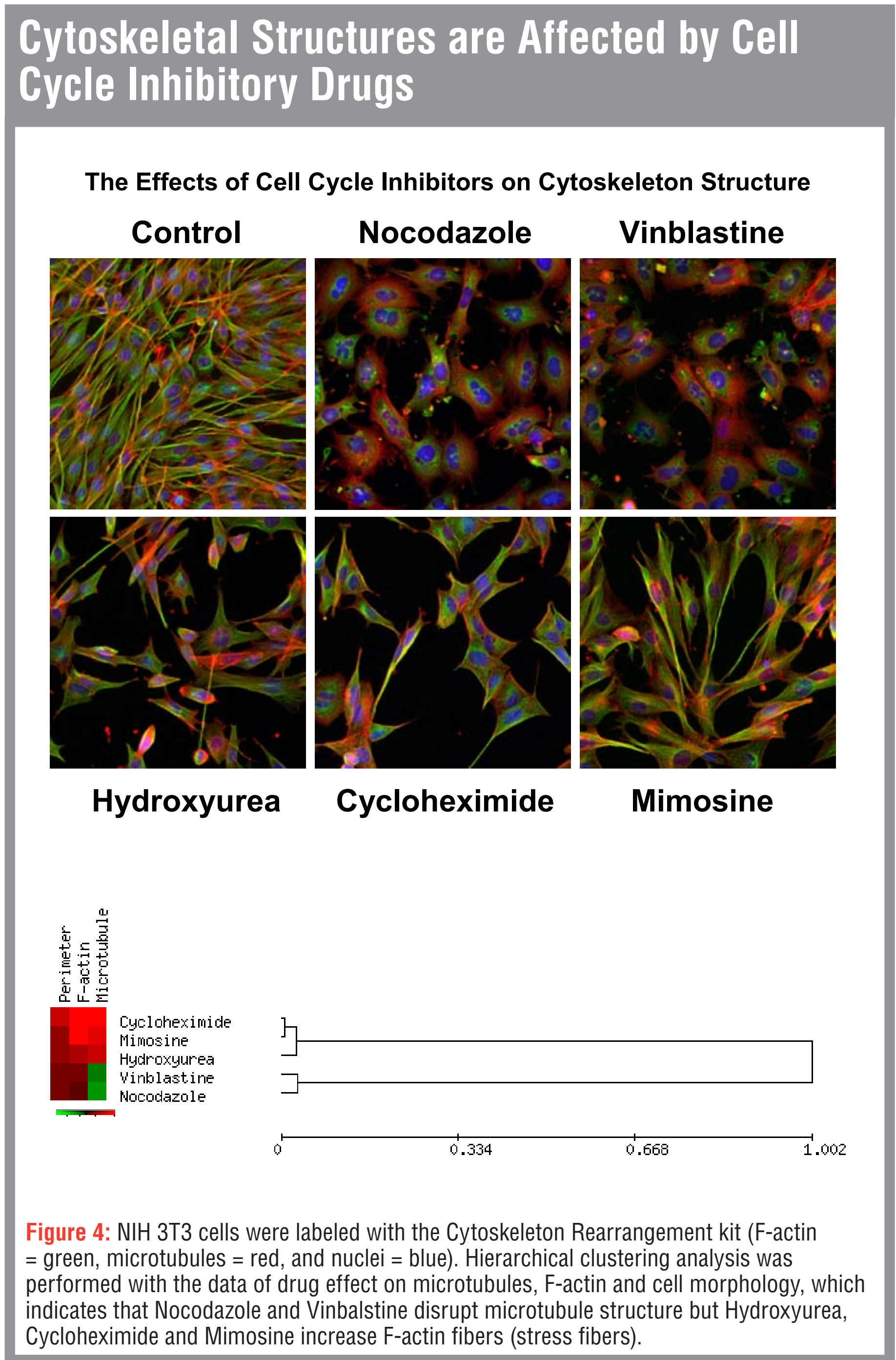
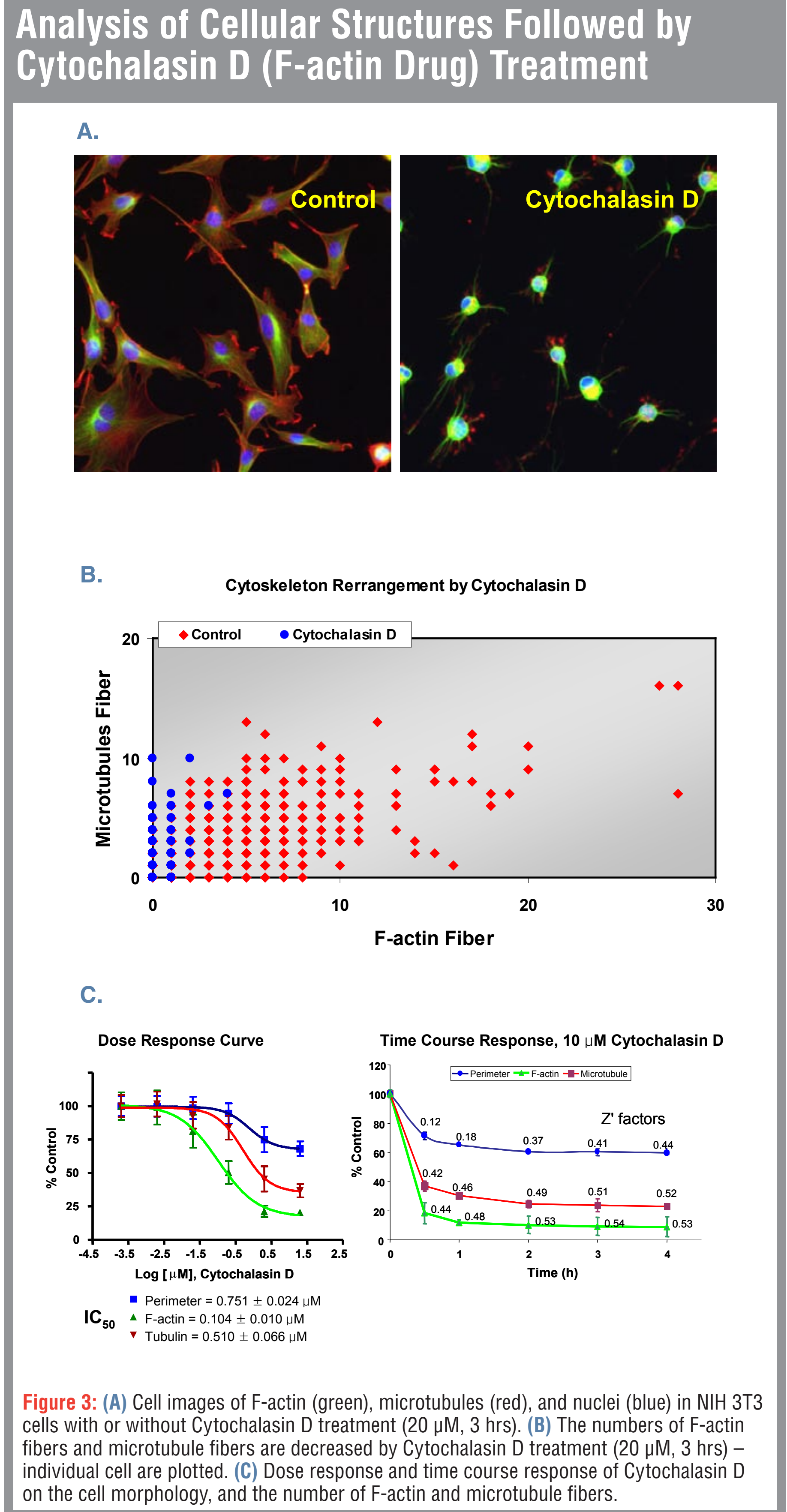
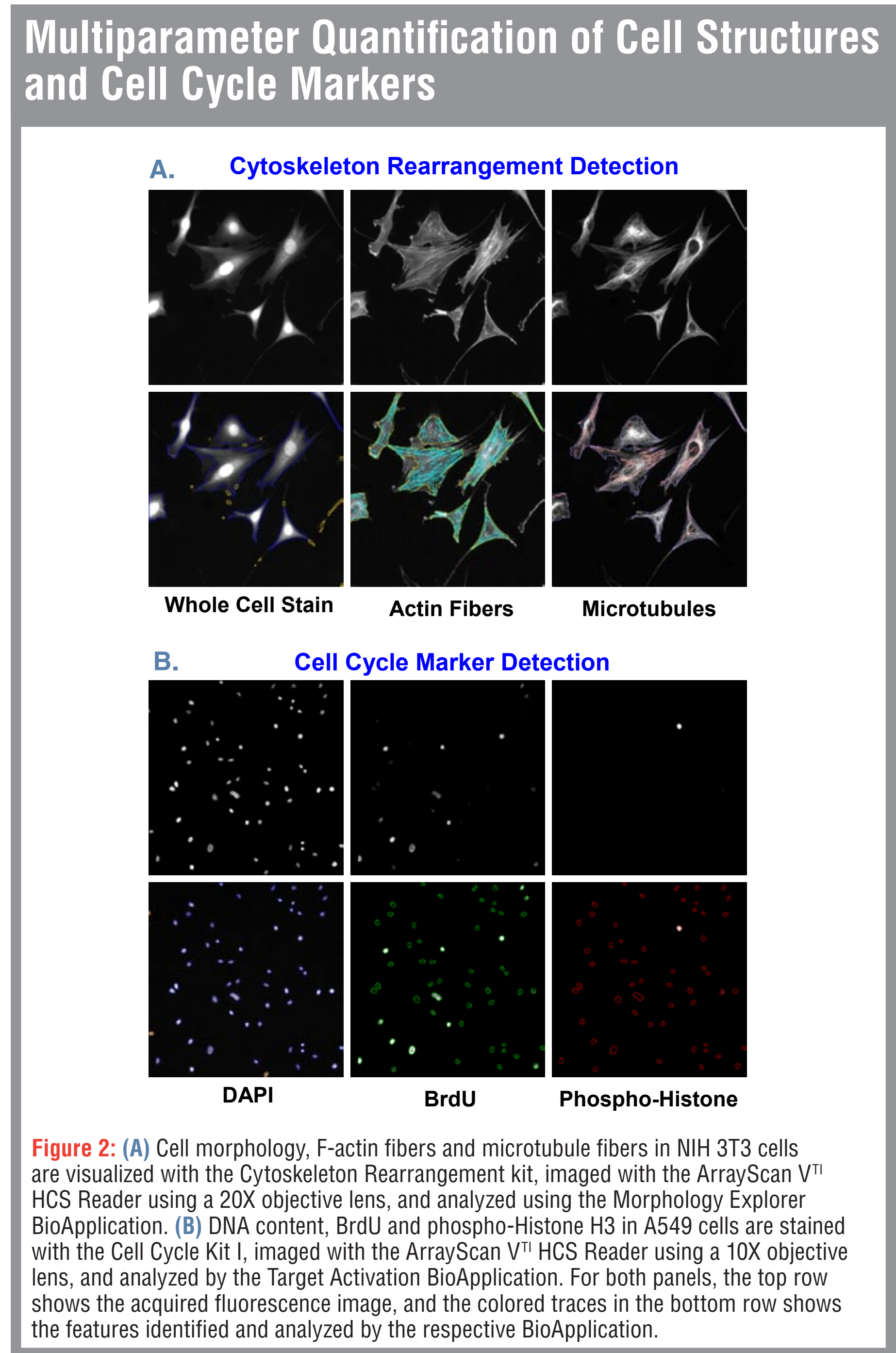
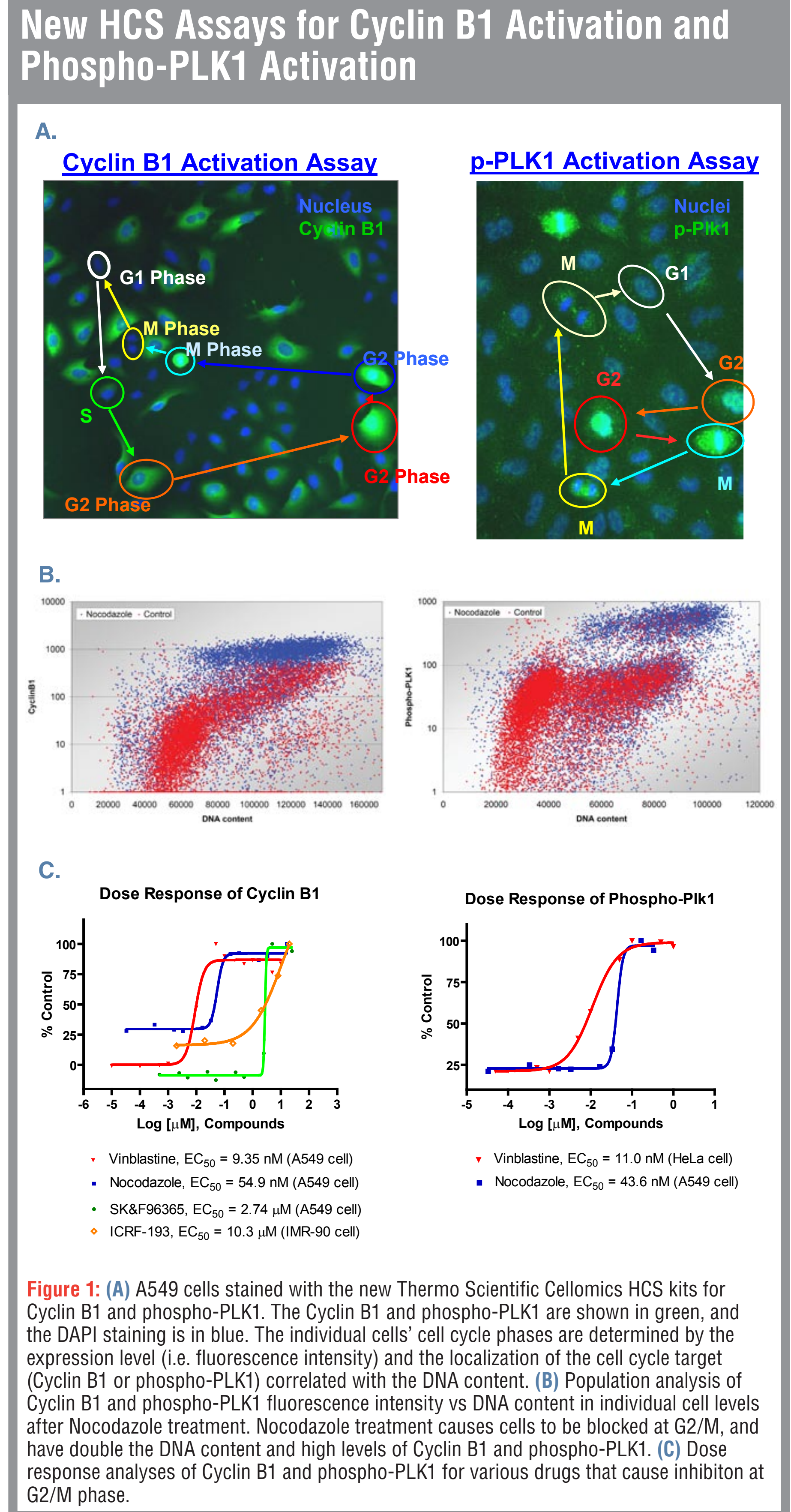
Introduction:

Cell stress, cell proliferation and the characterization of agents that either promote or inhibit cell proliferation are particularly of interest in cell biology and drug-discovery research. High content analysis (HCA) involves a fluorescence cell-based assay where the cells are automatically imaged and analyzed using quantitative fluorescence microscopy. HCA can be used to quantify the cell proliferation activity by measuring their DNA content, the state of cell cycle-associated proteins, and morphological and structural changes in individual cells and in cell populations.

We have developed a portfolio of Cellomics HCS kits to identify individual cells' cell cycle phase and proliferation state, using a fixed end-point HCS assay based on immunofluorescence detection in cells grown on standard high-density microplates. These kits all contain the DNA binding dye DAPI which is used to determine the cell cycle phase in individual cells by DNA content, and this is verified by correlating the expression level or phosphorylation state of other cell cycle associated indicators. Cells which have incorporated BrdU into DNA can be quickly detected by using monoclonal antibody directed against BrdU and fluorophore-conjugated secondary antibody. BrdU staining facilitates the identification of cells that have progressed through the S-phase of the cell cycle during the BrdU-labeling period. The Thermo Scientific Cellomics BrdU Cell Proliferation HCS kits have been specifically designed to enable simple BrdU detection in nuclear DNA without extreme treatments to the cell. This enables multiplexed detection of BrdU with other targets in the same cells. We also developed several other cell cycle assay targets, which include Cyclin B1, phosphorylated polo-like kinase 1 (phospho-PLK1), phosphorylated retinoblastoma protein (phospho-Rb) and phosphorylated Histone H3. These kits enable different options to directly determine individual cells' cell cycle phase, such as the G1, S phase (DNA content & BrdU) and G2/M phase (Cyclin B1, Phospho-PLK1, Phospho-Rb and mitosis specific Histone H3 phosphorylation).

The intracellular cytoskeletal meshwork is responsible for maintaining cell shape, cell movement, cytokinesis, and the organization of organelles within the cell. The dynamic network of the cytoskeleton also facilitates proper function of other cellular proteins through direct binding, transportation, repositioning and sequestration of these proteins. The structure of the cytoskeleton is controlled by bundling, aligning and repositioning of the filaments through cytoskeleton-associated proteins in response to the external signaling. Alterations of cytoskeletal structure are often associated with pathologies and cell death. Signaling defects in conjunction with cytoskeletal rearrangement can also contribute to increased cell proliferation rate and tumor cell motility which result in metastasis. Thus, the key players controlling cytoskeletal rearrangement in the cell are potential therapeutic targets.

Thermo Scientific, Inc. provides an integrated set of products that work together to deliver a “total solution” platform for HCS. Integral components of this set are the Thermo Scientific Cellomics HCS Reagent Kits. These kits provide easy-to-use methods and reagents for preparing high-quality samples for automated cell-based imaging assays. Using Thermo Scientific Cellomics Cytoskeleton Rearrangement kit and other Cell Cycle kits, we have developed quantitative HCS assays for assessing the cytoskeletal rearrangement and the cell cycle phases in cells treated with different compounds. The Cellomics HCS kits makes these assays easy to implement, and provides a powerful set of multiplexed parameters for monitoring the response of the cell due to drug treatment in these key functional areas. All images and data shown were acquired and quantitatively analyzed using Thermo Scientific's Cellomics ArrayScan® HCS Reader with associated Cellomics BioApplication image processing software modules.



Summary:

- **New Cell Cycle Marker analysis provide comprehensive information on cell cycle phases (Kits: Cyclin B1:#8404401&2, p-PLK1:#8404801&2, p-Rb:#8404501&2 and BrdU/p-Histone H3:#8404601&2).**
- **Cytoskeletal rearrangement (#8402401&2) and cell cycle marker detection allow to evaluate and profile the multiparameter cellular effect of cell cycle inhibitory drugs.**
- **Thermo Scientific Cellomics HCS Reagent Kits and ArrayScan HCS Reader offer the powerful tools for High Content Cell-Based Screening and Analysis**