

Development of a 3-color assay for mitosis and apoptosis using the laser-scanning IsoCytTM and comparison with the microscope-based INCell Analyzer 1000

Nadya Smotrov^{1*}, Grace K.Y.Chan^{1*}, Alexander A. Szwczak¹, Nathan Bays¹, Steven C. Miller², Evan F.Cromwell², Ilona Kariv¹

*Both authors contributed equally

¹Automated Lead Optimization, Merck Research Laboratories, Boston, MA, USA, ²Blueshift Biotechnologies, Inc., Sunnyvale, CA



Abstract

High Content Screening (HCS) has become increasingly popular for compound profiling in support of lead optimization because it allows quantification of multiple markers at the cellular level, providing information-rich analysis of complex biological systems. However, this information often comes at the price of slower assay speed. Here we describe the evaluation of the IsoCytTM from Blueshift Biotechnologies, a new instrument which has the capability to combine the best of both High Throughput Screening (HTS) and HCS.

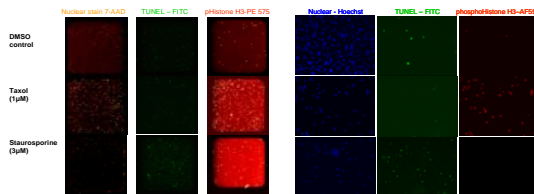
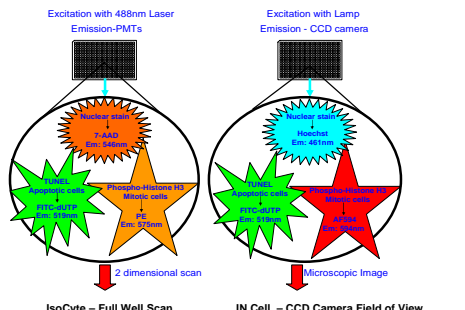
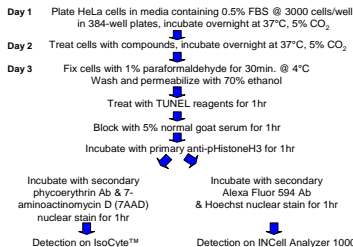
The IsoCytTM is a fast laser scanning fluorimeter capable of imaging all cells in a single well at a rate closer to HTS requirements, while still providing enhanced cell-by-cell analysis. Using the IsoCytTM equipped with one 488 nm laser we developed a three-color fluorescence cellular assay with detection of apoptosis by terminal deoxynucleotide transferase dUTP nick end labeling (TUNEL) and detection of mitosis via quantification of phospho-histone H3 levels. Data from the IsoCyt was compared with data obtained using the INCell Analyzer 1000 (GE Healthcare), a traditional automated microscope-based imager. Despite the different methods and modes of detection, the results from both instruments were very comparable. Our results demonstrate that the IsoCytTM system provides multiplexed intensity measurements with similar accuracy, yet requires significantly less imaging time compared to the INCell Analyzer.

Introduction

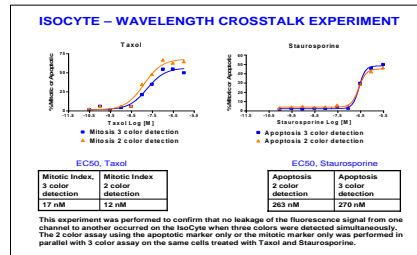
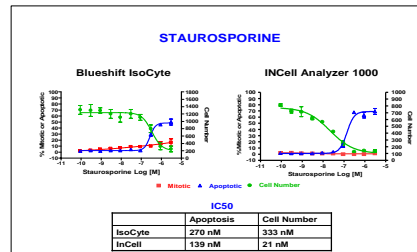
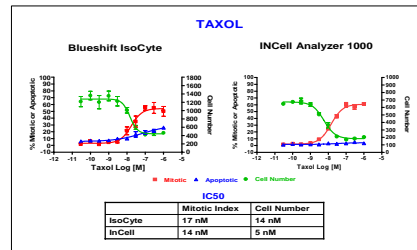
Modulation of the cell cycle and processes leading to apoptosis are common targets for pharmacological interventions in drug discovery. Applying high content screening technology, we developed a rapid assay to simultaneously measure these important cellular events using the Blueshift IsoCytTM and validated it against the INCell Analyzer 1000, a more traditional high content imaging platform.

We measured the mitotic index by monitoring levels of the mitotic marker phospho-histone H3 (phospho-Ser28), while apoptotic cells were identified using the TUNEL method to detect single-strand DNA breaks. At the same time, a nuclear stain allowed cell counting as an indication of proliferation.

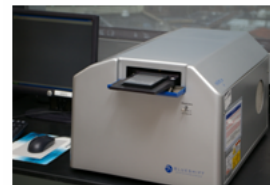
Method and Detection



Results



BlueShift IsoCyt



InCell Analyzer 1000



INSTRUMENT COMPARISON

	IsoCyt TM	INCell Analyzer 1000
Acquisition Speed	~ 2 min / 384 well plate	~ 105 min / 384 plate (3 image fields/well)
Acquisition Area	Whole well in 1 scan	Fraction of a well per image field
Analysis	Simultaneous with acquisition	Separate analysis required
Applications	Best suited for intensity-based measurements	Best suited for measuring morphology changes and subcellular events

Conclusion

We have developed a rapid multi-parametric high content assay for simultaneous detection of apoptosis, mitotic arrest and proliferation in a single microplate well using the IsoCyt laser-based fluorimeter. Our data demonstrates that the IsoCyt is comparable in accuracy to traditional microscope-based imaging system for cell fluorescence intensity measurements, but has the added advantages of fast acquisition with concurrent image analysis. It also images an entire well, collecting data on a much larger population of cells. Based on our evaluation, this type of high content assay on the IsoCyt can be utilized not only for lead optimization but meets the needs of high throughput primary screening. Together the IsoCyt and INCell 1000 are two complementary instruments that allow an HTS group to perform a broad range of high content experiments -- from detecting subcellular features to rapidly analyzing large populations of cells.

C2 The Merck logo within the grey bar is part of a new branding policy and should not be altered

Campbelj, 7/28/2004