

# Multiparametric analysis of cell health status using flow cytometry

Quentin Low, Jolene Bradford, Veronica Calderon and Leticia Montoya  
Thermo Fisher Scientific, 29851 Willow Creek Road, Eugene, OR 97402

## ABSTRACT

Cellular responses to drug treatment can be varied and heterogeneous. Often, a multiparametric approach is necessary to identify cellular pathways that are responding or affected. To maximize the number of parameters analysed per sample, we aimed to combine several functional sensors into one assay. In addition, we set out to streamline the workflow focusing on live cell responses. To this end, we designed a flow cytometry-based detection system centered on a single incubation condition for sensors that have minimal spectral overlap. Analysing Jurkat cells treated with camptothecin, a topoisomerase inhibitor that triggers cellular apoptosis, we were able to simultaneously detect the loss of cells in G2/M phase using the stoichiometric DNA dye, Hoechst 33342; an increase in apoptotic cells with a sensor for activated caspase 3/7 (CellEvent Caspase-3/7 Green); the loss of mitochondrial membrane potential with tetramethylrhodamine, methyl ester (TMRM); and an increase in dead cells as well as cells in late apoptotic stage, using a SYTOX viability dye. We also tested another four-sensor combination that focused on detecting cellular stress. Upon prolonged camptothecin induction, we were able to detect an increase in mitochondrial stress with MitoSOX, a sensor that detects mitochondrial oxidative stress; an increase in cellular stress with a CellROX reagent specific for reactive oxygen species. There was also a loss in mitochondrial membrane potential as well as a greater loss of cellular viability. In summary, using a model of cellular response to drug treatment, we performed multiparametric cell status analysis with a combination of cellular detection reagents on a flow cytometer.

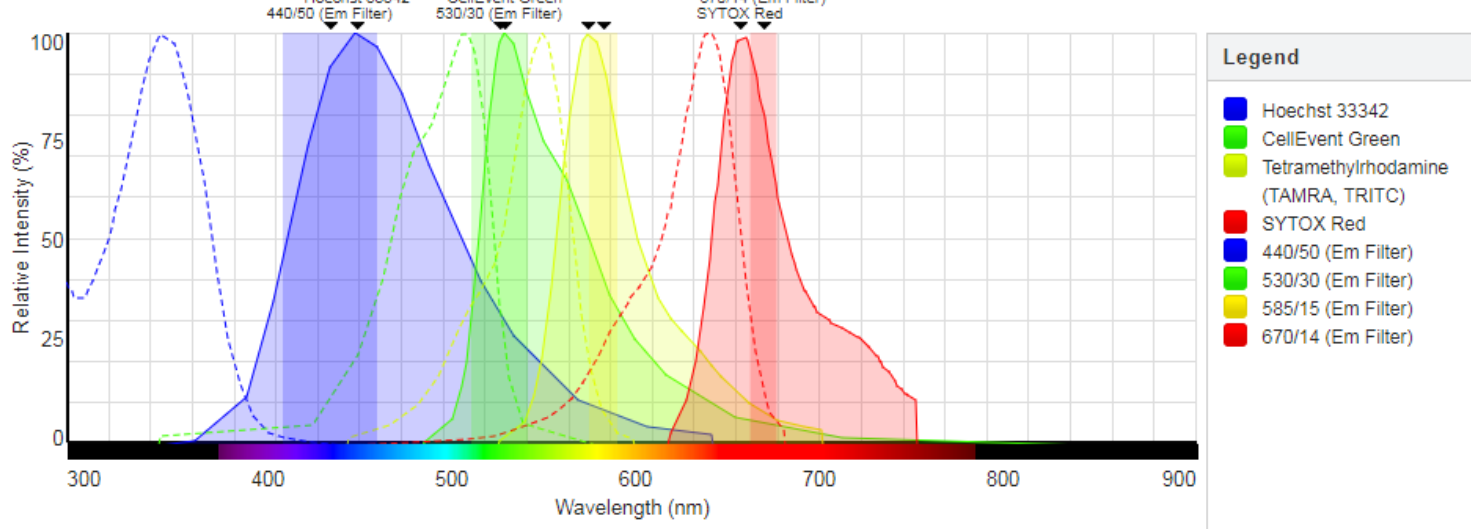
## INTRODUCTION

Determination of cell status has often been limited to single parameter analysis. This approach limits the number of parameters examined, usually resulting in sequential testing using different reagents that target distinct biological readouts. Compounding the limited throughput of a single readout assay, single parameter analysis have the potential to miss out on other confounding biological effects. Flow cytometry leads itself to multiparametric analysis. In this study, we highlight the use of flow cytometric analysis in detecting four different cell status parameters in a single analytical run. One set of sensor will detect cell cycle state, induction of apoptosis, mitochondrial health and cell viability. The second set will determine the stress state of the cell by measuring mitochondrial stress, cellular stress, mitochondrial health and cellular viability. These sensors are marketed as single parameter reagents, but in this study, we show that four different sensors are compatible in live cell staining to demonstrate cellular health. Furthermore, we are able to identify hits from a compound library using a combination of these sensors on a flow cytometer.

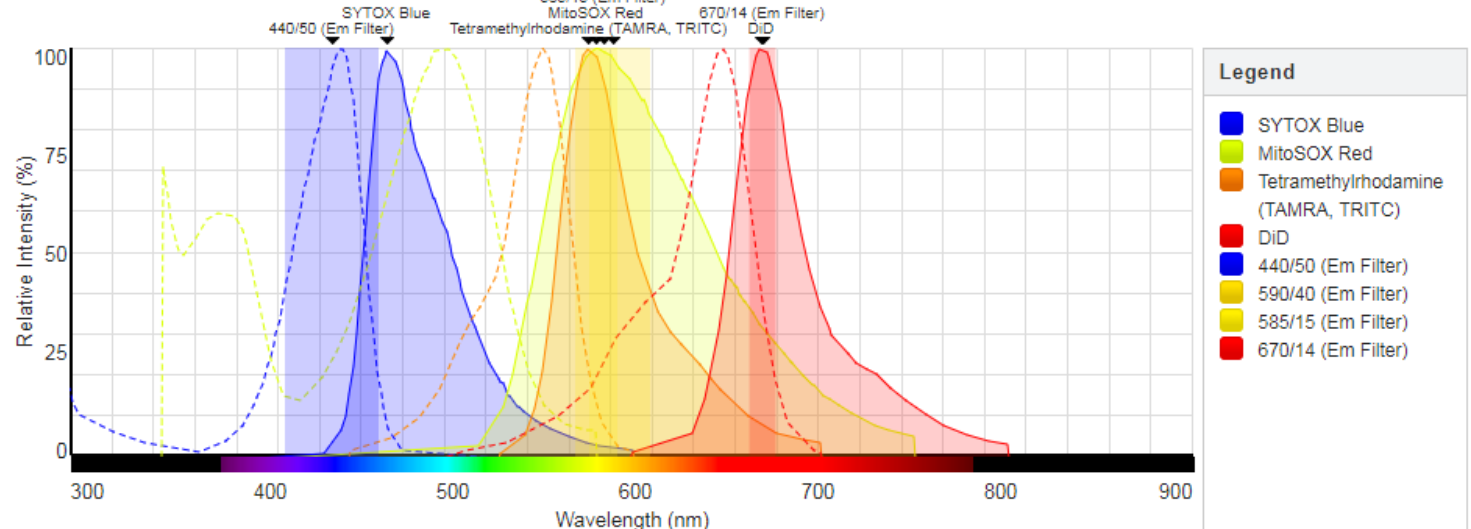
## MATERIALS AND METHODS

**Verification testing:** Jurkat cells were resuspended at  $1 \times 10^6$  cells/ml, and 1 ml of cells were stained with sensors in FACS tubes. A concentration range from 0.1  $\mu$ M to 10  $\mu$ M of camptothecin was used to stimulate the cells for 18 hours, in 37°C with 5% CO<sub>2</sub>. Individual sensors outlined in table 1 was added either individually or in combination. Sensors were incubated with the cells for 30 minutes at 37°C with 5% CO<sub>2</sub>. After that period, cellular responses were analyzed using an Attune NxT Acoustic Cytometer.

**Treatment for compound library analysis:** For this test, 10,000 cells in 80  $\mu$ l of RPMI 1640 + 10% fetal bovine serum were plated into each well of a 96 well plate. Compounds from one plate of the Tocriscreen Mini (Tocris, Cat No.2890) were used. The initial 10  $\mu$ M stock was diluted to 100  $\mu$ M. This allowed for a 1:10 addition of the compounds to 90  $\mu$ l of cell to result in the final compound concentration of 10  $\mu$ M. Incubation of the compounds with the cells were carried out in 37°C with 5% CO<sub>2</sub> for 18 hours. After which, components of MultiKit 1 and MultiKit 2 were added to the final assay concentrations. Cells were returned to a cell culture incubator for 30 mins prior to flow cytometric analysis, using an Attune NxT Acoustic Cytometer with the attached Autosampler.



MultiKit 1	SKU	Assay Concentration
Hoechst 33342	H21492	10 $\mu$ g/ml
CellEvent Caspase-3/7 Green Flow Cytometry Assay Kit	C10427	500 nM
MitoProbe TMRM Assay Kit for Flow Cytometry	M20036	20 nM
SYTOX Red Dead Cell Stain, for 633 or 653 nm excitation	S34859	5 nM



MultiKit 2	SKU	Assay Concentration
SYTOX Blue Dead Cell Stain, for flow cytometry	S35857	1 $\mu$ M
MitoSOX Red Mitochondrial Superoxide Indicator, for live-cell imaging	M36008	1 $\mu$ M
MitoProbe TMRM Assay Kit for Flow Cytometry	M20036	20 nM
CellROX Deep Red Flow Cytometry Assay Kit	C10491	500 nM

## RESULTS

Figure 1. Individual cell status sensors can be used in combination

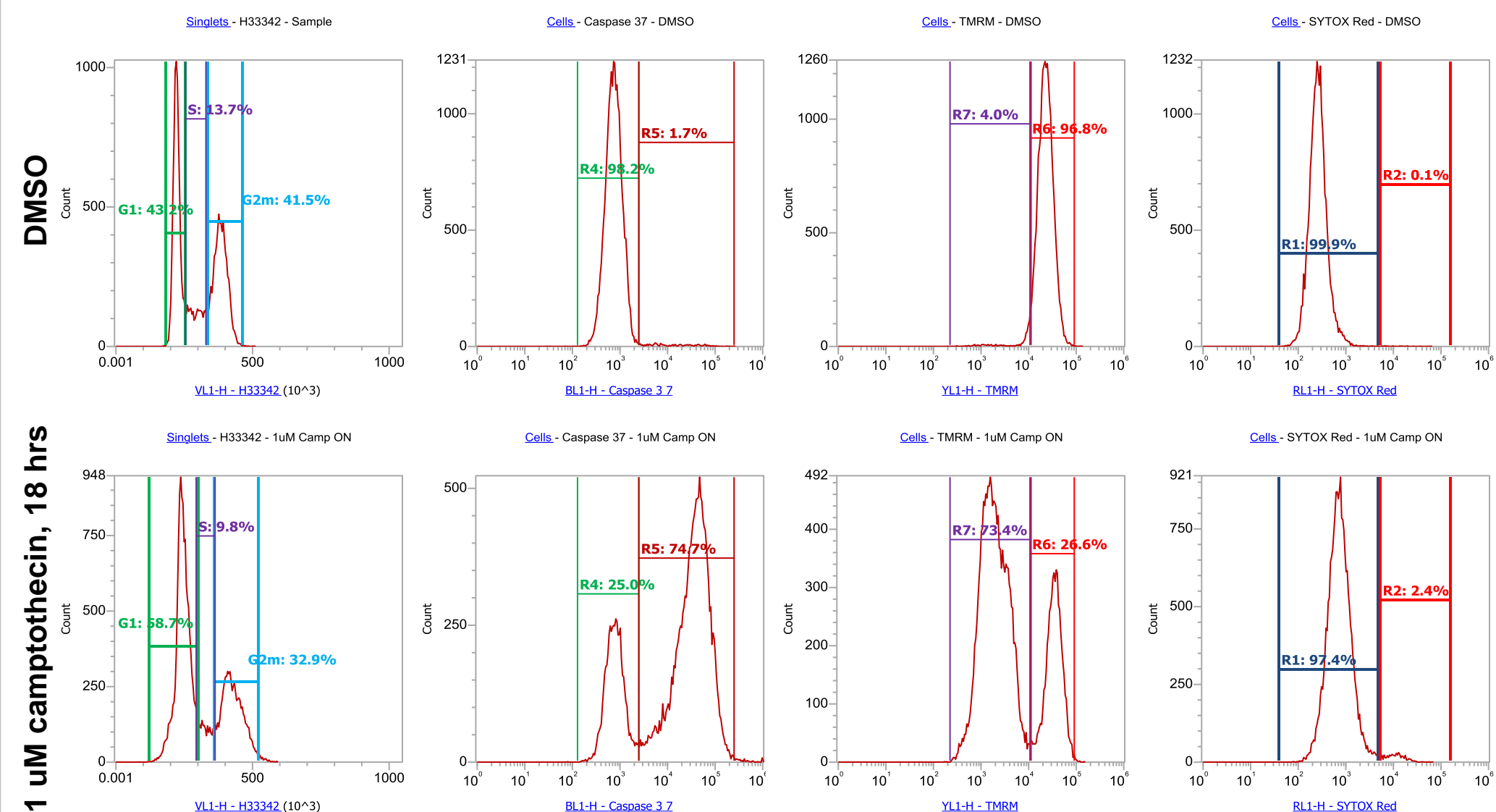


Figure 1a. Individually added sensors detect the concomitant changes in apoptotic state and mitochondrial health, along with an increase of cells in the G1 phase and a small increase in cell death resulting from camptothecin treatment.

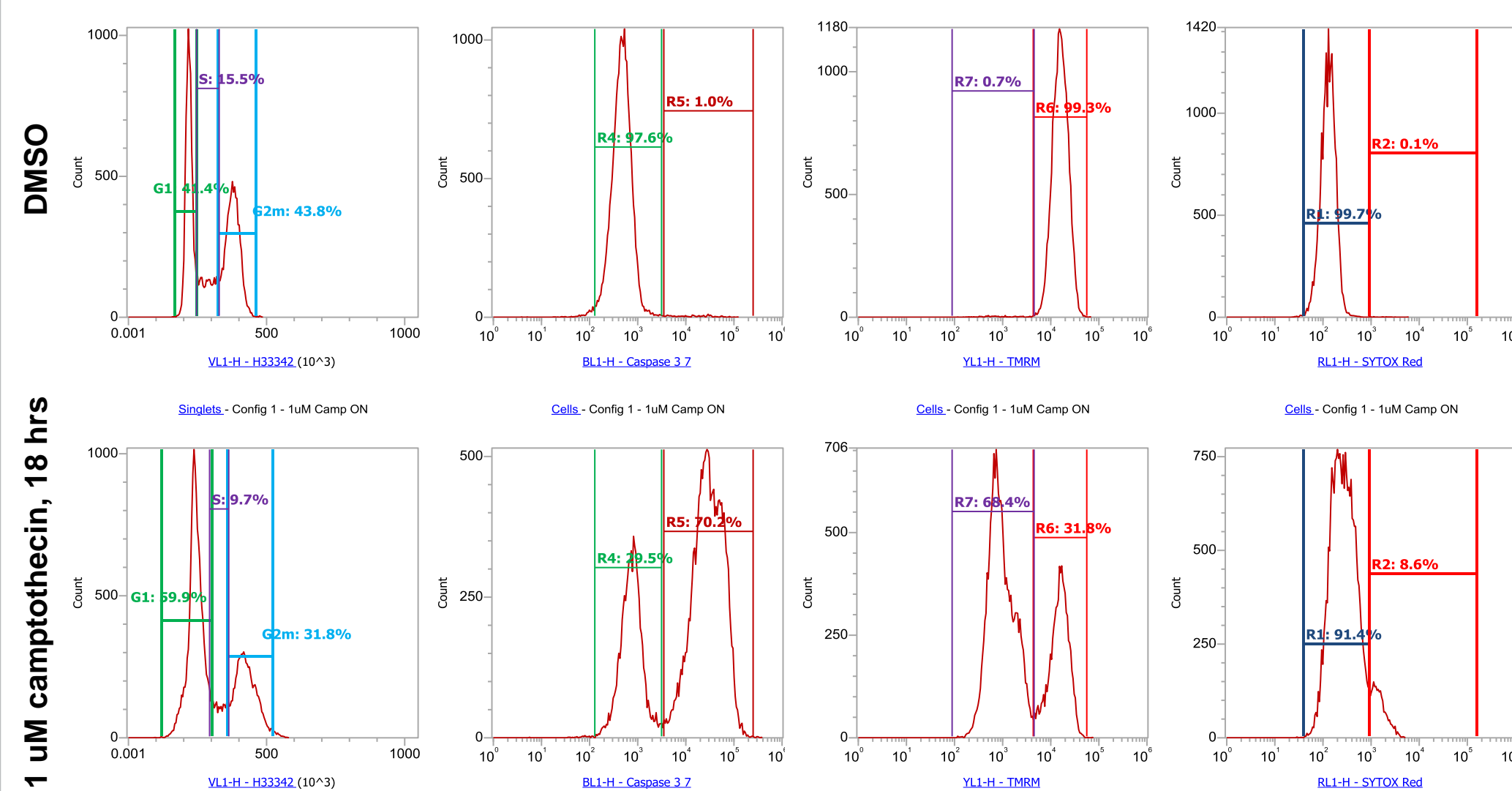


Figure 1b. Sensors added in combination detected camptothecin-related effects similar to Figure 1a. Apoptotic state (70% vs 75%), mitochondrial health (68% vs 73%), increase of cells in the G1 phase (60% vs 59%) and the increase in cell death (8.6% vs 2.4%). Detection was performed without fluorescence compensation.

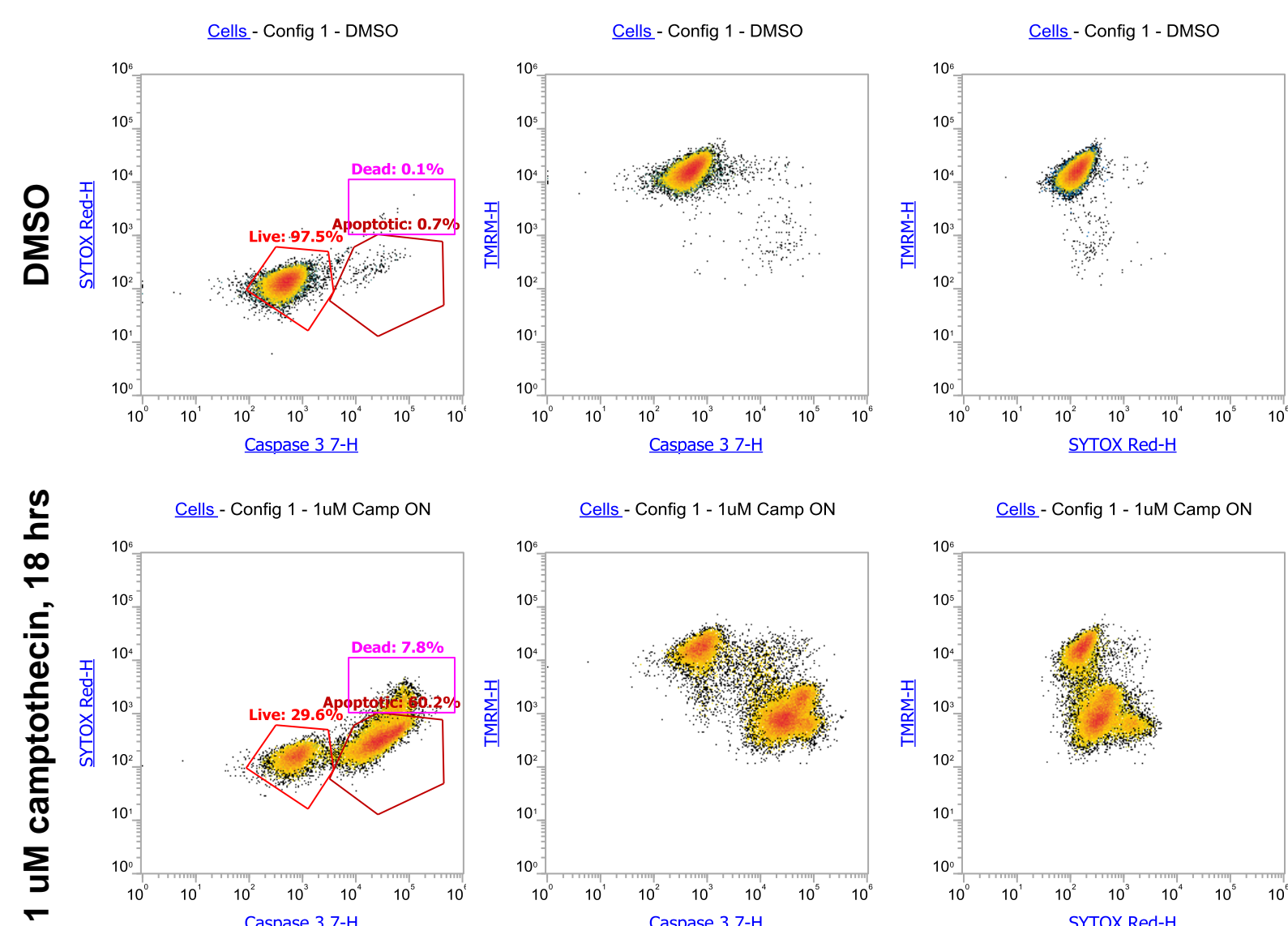


Figure 1c. Multiparameter analysis of Jurkat cells treated with 1  $\mu$ M camptothecin for 18 hours show cells undergoing apoptosis with some that are dead/late apoptotic (Left Plots). Cell that undergo apoptotic changes have a reduce mitochondrial function as is evident in the loss of TMRM staining (Middle Plots). Loss of TMRM staining and gain of SYTOX staining is consistent with what is expected for dead/late apoptotic cells (Right Plots).

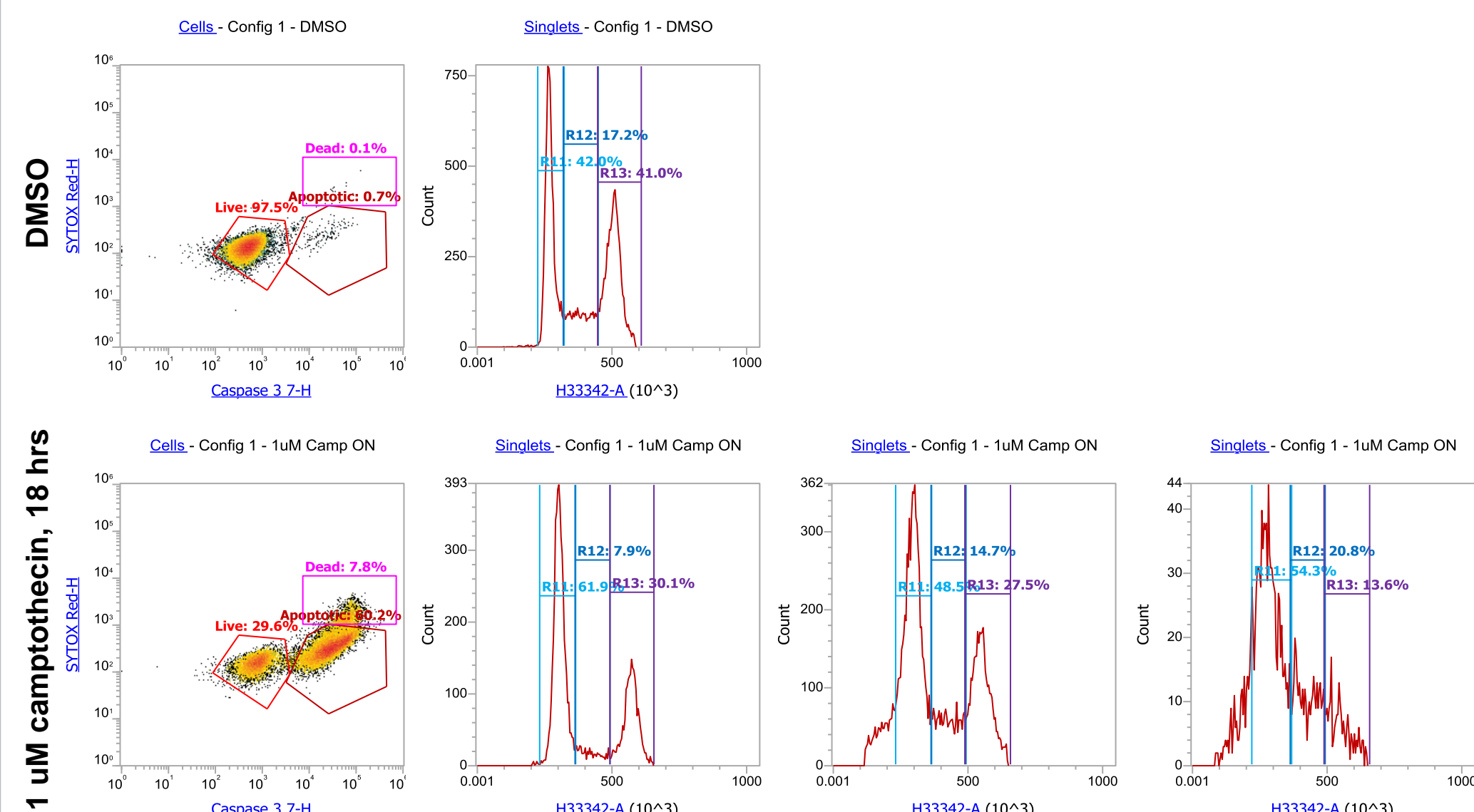


Figure 1d. Multiparameter cell cycle analysis of Jurkat cells treated with 1  $\mu$ M camptothecin for 18 hours. Relative to DMSO treated cells, camptothecin treatment resulted in loss of cells in G2/M phase (Bottom Histograms). The loss corresponded to the cellular states; treated live cells show a loss of S phase cells, with an increasing loss of G2/M phase cells as the cells undergo apoptosis.

Figure 2. Individual cell status sensors to detect cellular stress

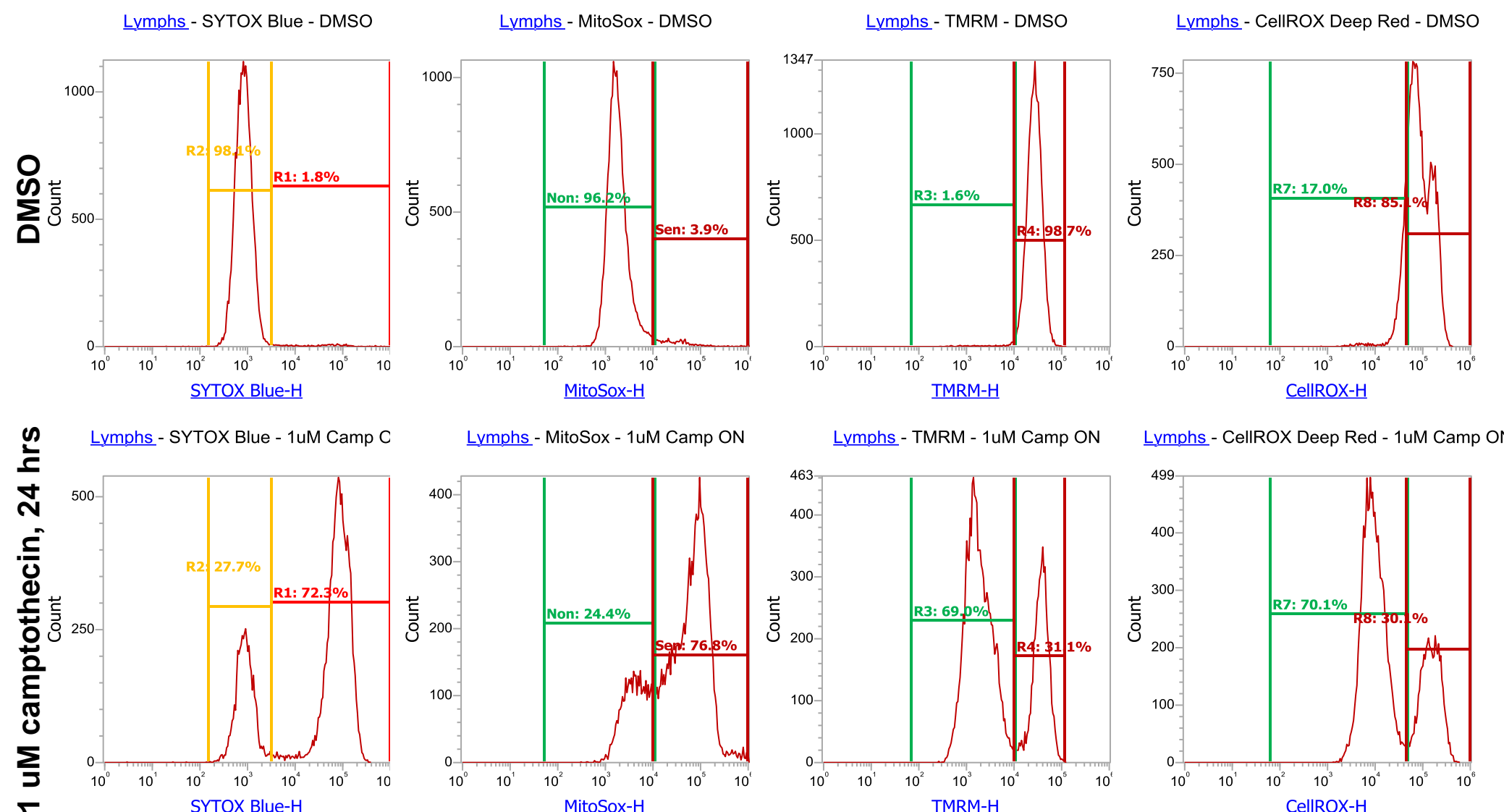


Figure 2a. Camptothecin treatment of Jurkat cells for 24 hours resulted in significant cell death, a loss of mitochondrial membrane potential and increases in mitochondrial oxidative stress.

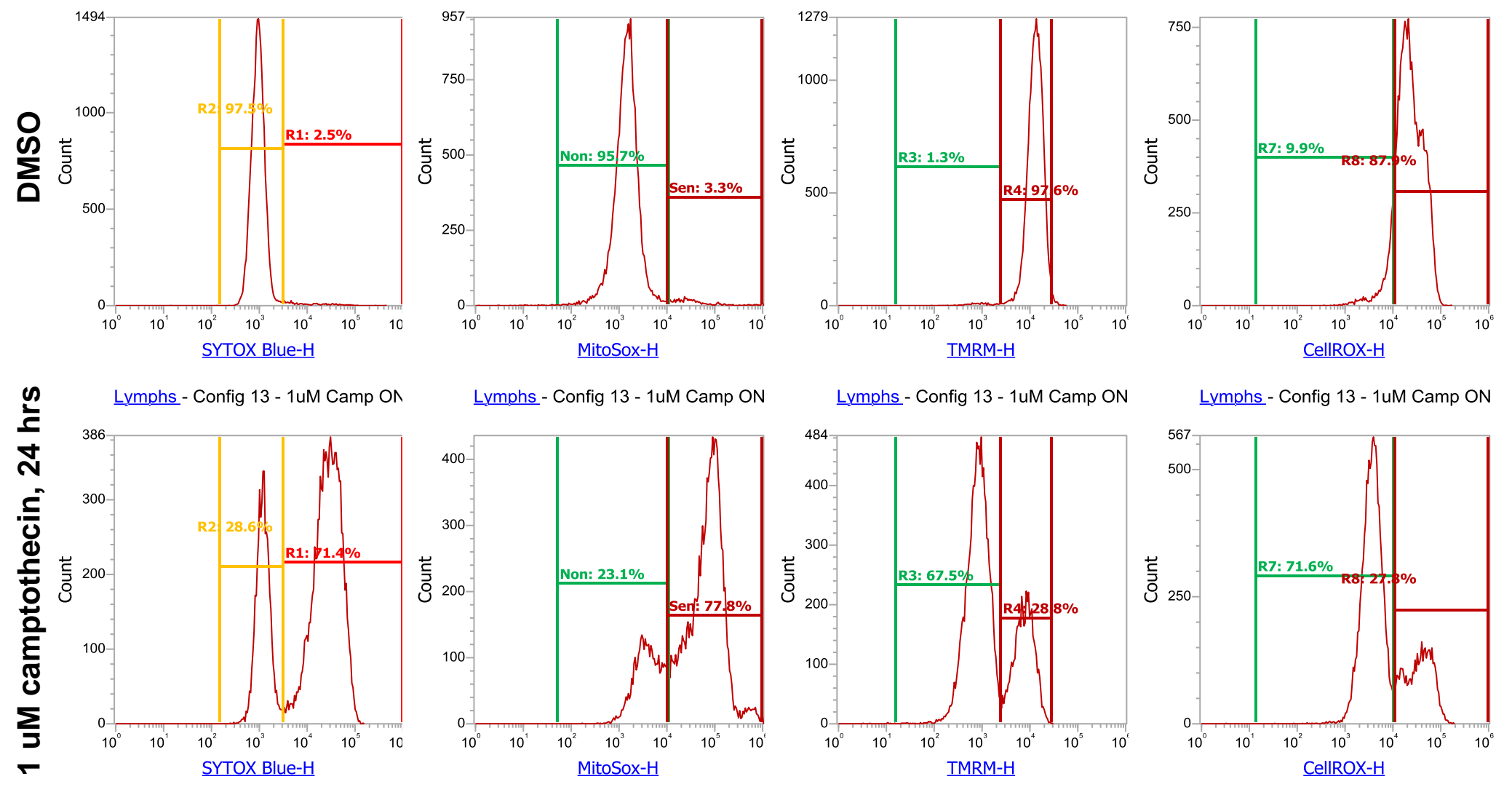


Figure 2b. Using MultiKit 2, the effects of camptothecin on Jurkat cells were detected in a multiparametric analysis. Consistent with the results in Figure 2a, treated Jurkat cells showed an increase in cell death, a loss of mitochondrial membrane potential and increases in mitochondrial oxidative stress. Unlike MultiKit 1, using MultiKit 2 required fluorescence compensation. Uncompensated data shown in lower, boxed row.

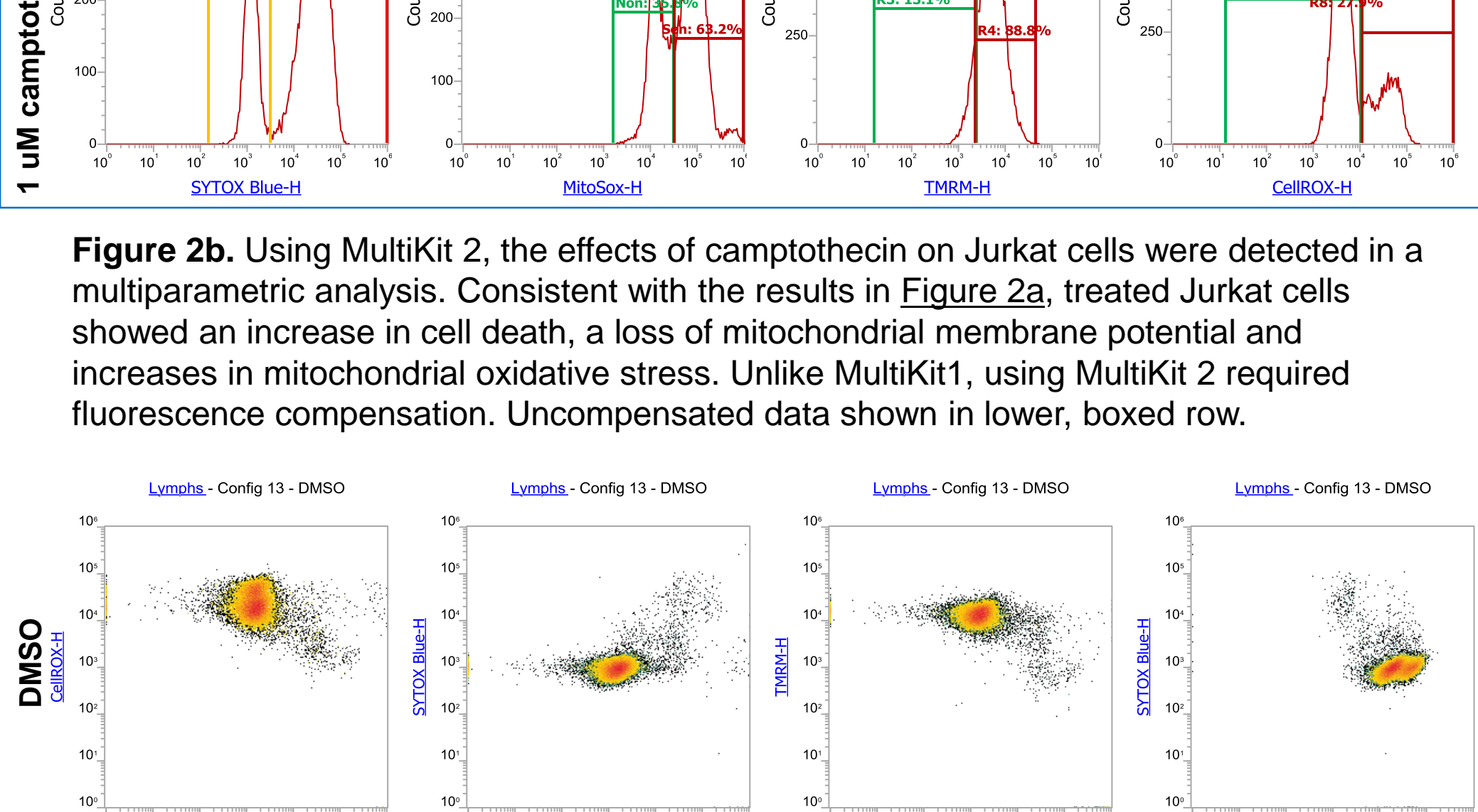


Figure 2c. Multiparameter analysis of Jurkat cells treated with 1  $\mu$ M camptothecin for 24 hours. There was no induction of cellular stress by camptothecin treatment, however, mitochondrial stress was detected along with a loss in mitochondrial membrane potential.

Figure 3. Screening with Tocriscreen Mini compound library

Position	Batch Specific Product Name	Reagent Description	Position	Batch Specific Product Name	Reagent Description
A01	AMPA	Selective NMDA agonist	E01	AMPA	Selective NMDA agonist
A02	AMPA	Selective NMDA agonist	E02	AMPA	Selective NMDA agonist
A03	AMPA	Selective NMDA agonist	E03	AMPA	Selective NMDA agonist
A04	AMPA	Selective NMDA agonist	E04	AMPA	Selective NMDA agonist
A05	AMPA	Selective NMDA agonist	E05	AMPA	Selective NMDA agonist
A06	AMPA	Selective NMDA agonist	E06	AMPA	Selective NMDA agonist
A07	AMPA	Selective NMDA agonist	E07	AMPA	Selective NMDA agonist
A08	AMPA	Selective NMDA agonist	E08	AMPA	Selective NMDA agonist
A09	AMPA	Selective NMDA agonist	E09	AMPA	Selective NMDA agonist
A10	AMPA	Selective NMDA agonist	E10	AMPA	Selective NMDA agonist
A11	AMPA	Selective NMDA agonist	E11	AMPA	Selective NMDA agonist
A12	AMPA	Selective NMDA agonist	E12	AMPA	Selective NMDA agonist
A13	AMPA	Selective NMDA agonist	E13	AMPA	Selective NMDA agonist
A14	AMPA	Selective NMDA agonist	E14	AMPA	Selective NMDA agonist
A15	AMPA	Selective NMDA agonist	E15	AMPA	Selective NMDA agonist
A16	AMPA	Selective NMDA agonist	E16	AMPA	Selective NMDA agonist
A17	AMPA	Selective NMDA agonist	E17	AMPA	Selective NMDA agonist
A18	AMPA	Selective NMDA agonist	E18	AMPA	Selective NMDA agonist
A19	AMPA	Selective NMDA agonist	E19	AMPA	Selective NMDA agonist
A20	AMPA	Selective NMDA agonist	E20	AMPA	Selective NMDA agonist
A21	AMPA	Selective NMDA agonist	E21	AMPA	Selective NMDA agonist
A22	AMPA	Selective NMDA agonist	E22	AMPA	Selective NMDA agonist
A23	AMPA	Selective NMDA agonist	E23	AMPA	Selective NMDA agonist
A24	AMPA	Selective NMDA agonist	E24	AMPA	Selective NMDA agonist
A25	AMPA	Selective NMDA agonist	E25	AMPA	Selective NMDA agonist
A26	AMPA	Selective NMDA agonist	E26	AMPA	Selective NMDA agonist
A27	AMPA	Selective NMDA agonist	E27	AMPA	Selective NMDA agonist
A28	AMPA	Selective NMDA agonist	E28	AMPA	Selective NMDA agonist
A29	AMPA	Selective NMDA agonist	E29	AMPA	Selective NMDA agonist
A30	AMPA	Selective NMDA agonist	E30	AMPA	Selective NMDA agonist
A31	AMPA	Selective NMDA agonist	E31	AMPA	Selective NMDA agonist
A32	AMPA	Selective NMDA agonist	E32	AMPA	Selective NMDA agonist
A33	AMPA	Selective NMDA agonist	E33	AMPA	Selective NMDA agonist
A34	AMPA	Selective NMDA agonist	E34	AMPA	Selective NMDA agonist
A35	AMPA	Selective NMDA agonist	E35	AMPA	Selective NMDA agonist
A36	AMPA	Selective NMDA agonist	E36	AMPA	Selective NMDA agonist
A37	AMPA	Selective NMDA agonist	E37	AMPA	Selective NMDA agonist
A38	AMPA	Selective NMDA agonist	E38	AMPA	Selective NMDA agonist
A39	AMPA	Selective NMDA agonist	E39	AMPA	Selective NMDA agonist
A40	AMPA	Selective NMDA agonist	E40	AMPA	Selective NMDA agonist
A41	AMPA	Selective NMDA agonist	E41	AMPA	Selective NMDA agonist
A42	AMPA	Selective NMDA agonist	E42	AMPA	Selective NMDA agonist
A43	AMPA	Selective NMDA agonist	E43	AMPA	Selective NMDA agonist
A44	AMPA	Selective NMDA agonist	E44	AMPA	Selective NMDA agonist
A45	AMPA	Selective NMDA agonist	E45	AMPA	Selective NMDA agonist
A46	AMPA	Selective NMDA agonist	E46	AMPA	Selective NMDA agonist
A47	AMPA	Selective NMDA agonist	E47	AMPA	Selective NMDA agonist
A48	AMPA	Selective NMDA agonist	E48	AMPA	Selective NMDA agonist
A49	AMPA	Selective NMDA agonist	E49	AMPA	Selective NMDA agonist
A50	AMPA	Selective NMDA agonist	E50	AMPA	Selective NMDA agonist
A51	AMPA	Selective NMDA agonist	E51	AMPA	Selective NMDA agonist
A52	AMPA	Selective NMDA agonist	E52	AMPA	Selective NMDA agonist
A53	AMPA	Selective NMDA agonist	E53	AMPA	Selective NMDA agonist
A54	AMPA	Selective NMDA agonist	E54	AMPA	Selective NMDA agonist
A55	AMPA	Selective NMDA agonist	E55	AMPA	Selective NMDA agonist
A56	AMPA	Selective NMDA agonist	E56	AMPA	Selective NMDA agonist
A57	AMPA	Selective NMDA agonist	E57	AMPA	Selective NMDA agonist
A58	AMPA	Selective NMDA agonist	E58	AMPA	Selective NMDA agonist
A59	AMPA	Selective NMDA agonist	E59	AMPA	Selective NMDA agonist
A60	AMPA	Selective NMDA agonist	E60	AMPA	Selective NMDA agonist
A61	AMPA	Selective NMDA agonist	E61	AMPA	Selective NMDA agonist
A62	AMPA	Selective NMDA agonist	E62	AMPA	Selective NMDA agonist
A63	AMPA	Selective NMDA agonist	E63	AMPA	Selective NMDA agonist
A64	AMPA	Selective NMDA agonist	E64	AMPA	Selective NMDA agonist
A65	AMPA	Selective NMDA agonist	E65	AMPA	Selective NMDA agonist
A66	AMPA	Selective NMDA agonist	E66	AMPA	Selective NMDA agonist
A67	AMPA	Selective NMDA agonist	E67	AMPA	Selective NMDA agonist
A68	AMPA	Selective NMDA agonist	E68	AMPA	Selective NMDA agonist
A69	AMPA	Selective NMDA agonist	E69	AMPA	Selective NMDA agonist
A70	AMPA	Selective NMDA agonist	E70	AMPA	Selective NMDA agonist
A71	AMPA	Selective NMDA agonist	E71	AMPA	Selective NMDA agonist
A72	AMPA	Selective NMDA agonist	E72	AMPA	Selective NMDA agonist
A73	AMPA	Selective NMDA agonist	E73	AMPA	Selective NMDA agonist
A74	AMPA	Selective NMDA agonist	E74	AMPA	Selective NMDA agonist
A75	AMPA	Selective NMDA agonist	E75	AMPA	Selective NMDA agonist
A76	AMPA	Selective NMDA agonist	E76	AMPA	Selective NMDA agonist
A77	AMPA	Selective NMDA agonist	E77	AMPA	Selective NMDA agonist
A78	AMPA	Selective NMDA agonist	E78	AMPA	Selective NMDA agonist
A79	AMPA	Selective NMDA agonist	E79	AMPA	Selective NMDA agonist
A80	AMPA	Selective NMDA agonist	E80	AMPA	Selective NMDA agonist
A81	AMPA	Selective NMDA agonist	E81	AMPA	Selective NMDA agonist
A82	AMPA	Selective NMDA agonist	E82	AMPA	Selective NMDA agonist
A83	AMPA	Selective NMDA agonist	E83	AMPA	Selective NMDA agonist
A84	AMPA	Selective NMDA agonist	E84	AMPA	Selective NMDA agonist
A85	AMPA	Selective NMDA agonist	E85	AMPA	Selective NMDA agonist
A86	AMPA	Selective NMDA agonist	E86	AMPA	Selective NMDA agonist
A87	AMPA	Selective NMDA agonist	E87	AMPA	Selective NMDA agonist
A88	AMPA	Selective NMDA agonist	E88	AMPA	Selective NMDA agonist
A89	AMPA	Selective NMDA agonist	E89	AMPA	Selective NMDA agonist
A90	AMPA	Selective NMDA agonist	E90	AMPA	Selective NMDA agonist
A91	AMPA	Selective NMDA agonist	E91	AMPA	Selective NMDA agonist
A92	AMPA	Selective NMDA agonist	E92	AMPA	Selective NMDA agonist
A93	AMPA	Selective NMDA agonist	E93	AMPA	Selective NMDA agonist
A94	AMPA	Selective NMDA agonist	E94	AMPA	Selective NMDA agonist
A95	AMPA	Selective NMDA agonist	E95	AMPA	Selective NMDA agonist
A96	AMPA	Selective NMDA agonist	E96	AMPA	Selective NMDA agonist
A97	AMPA	Selective NMDA agonist	E97	AMPA	Selective NMDA agonist
A98	AMPA	Selective NMDA agonist	E98	AMPA	Selective NMDA agonist
A99	AMPA	Selective NMDA agonist	E99	AMPA	Selective NMDA agonist
A100	AMPA	Selective NMDA agonist	E100	AMPA	Selective NMDA agonist

Figure 4. Responses from compound library analysis

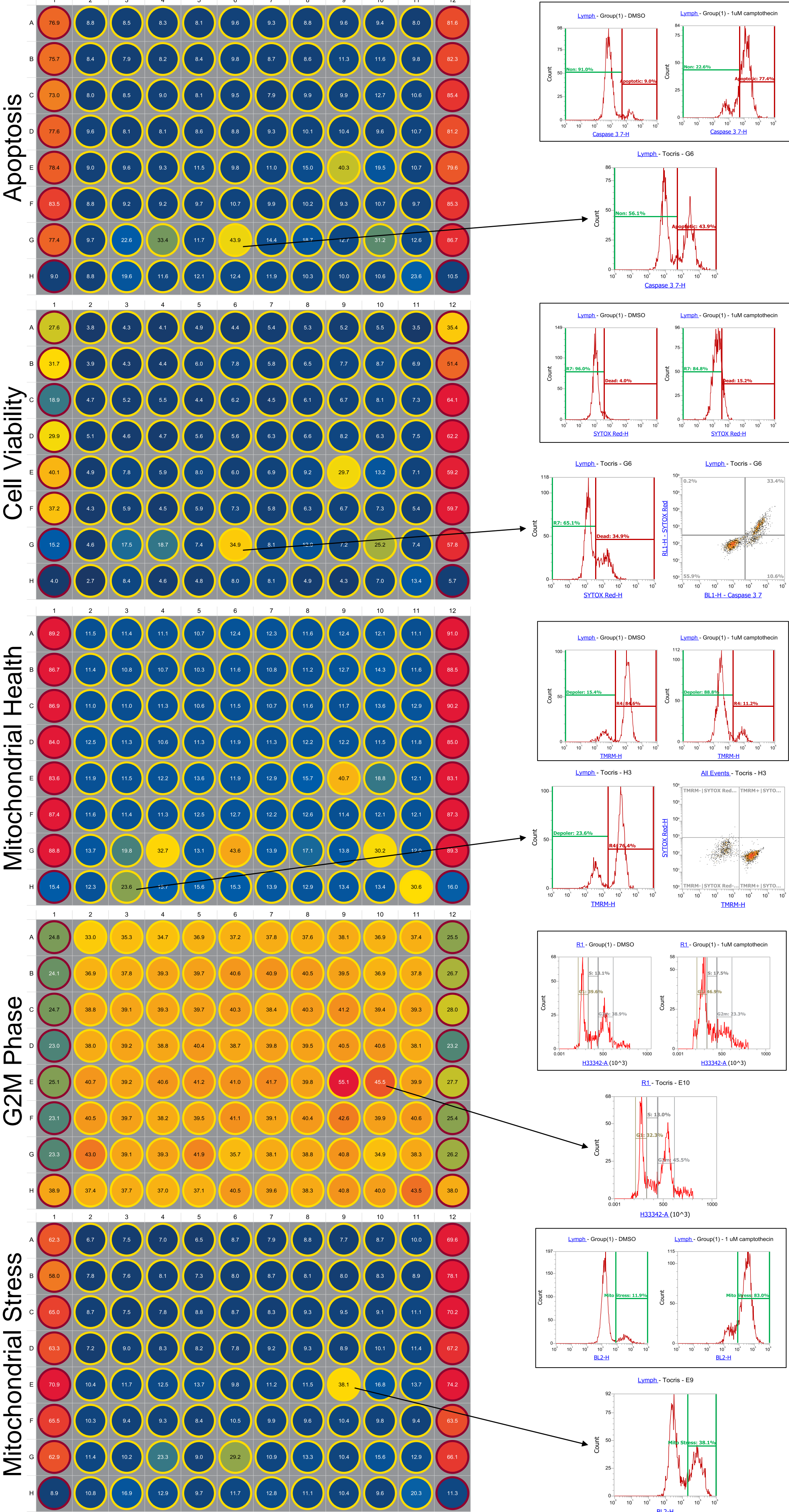


Figure 4. Multiparametric readouts from Jurkat cells in response to incubation with compounds from the Tocriscreen library. Control readouts are shown in boxed histograms. Heat map shows potential hits, and specific responses from those wells are highlighted (arrows).

## CONCLUSIONS

- Single parameter detection reagents can be used in combination to create multiparameter kits
- Spectrally distinct reagents allows for experimental setup without employing fluorescence compensation
- Multiparameter kits can be tailored to determining cellular responses in drug compound library analyses.

## REFERENCES

Sklar LA, Carter MB, Edwards BS. Curr Opin Pharmacol. 2007 Oct;7(5):527-34. Epub 2007 Jul 24.  
Feng Y, Mitchison TJ, Bender A, Young DW, Tallarico JA. Nat Rev Drug Discov. 2009 Jul;8(7):567-78. doi: 10.1038/nrd2876.

## LEGAL INFORMATION

For Research Use Only. Not for use in diagnostic procedures.

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others

**Thermo Fisher**  
SCIENTIFIC