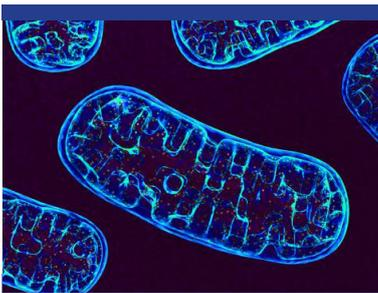


# Mitochondrial Oxidative Stress

## Background Information



'ROS trafficking between mitochondria could constitute a positive-feedback mechanism resulting in an elevated production of ROS that could be propagated throughout the cell and may cause perceptible mitochondrial and cellular injury.'

<sup>1</sup>Zorov DB *et al.* (2014) *Physiol Rev* **94**(3); 909-950

- Mitochondria consume the majority of cellular oxygen and regulate redox-signalling<sup>1</sup>.
- Toxic drugs can induce mitochondrial reactive oxygen species (mROS) causing mitochondrial/cellular damage which has been linked to the pathology of many diseases<sup>2</sup>.
- The Cyprotex mitochondrial oxidative stress assay detects selective mROS production caused by the toxicity of novel compounds.

### Protocol

#### Cell Line

HepG2 cell line; other cell types on request

#### Analysis Platform

Cellomics ArrayScan® CX7, XTI and VTI (Thermo Scientific)

#### Test Compound Concentrations

8 point dose response curve with top concentration based on 100x C<sub>max</sub> or solubility limit\*

#### Compound Requirements

50 µL of a solution to achieve 100x C<sub>max</sub> (200x top concentration to maintain 0.5% DMSO) or equivalent amount in solid compound

#### Number of Replicates

3 replicates per concentration\*

#### Time Points

24 hour\*

#### Quality Controls

Negative control: 0.5% DMSO (vehicle)

Positive controls: 2 known mROS-inducing compounds

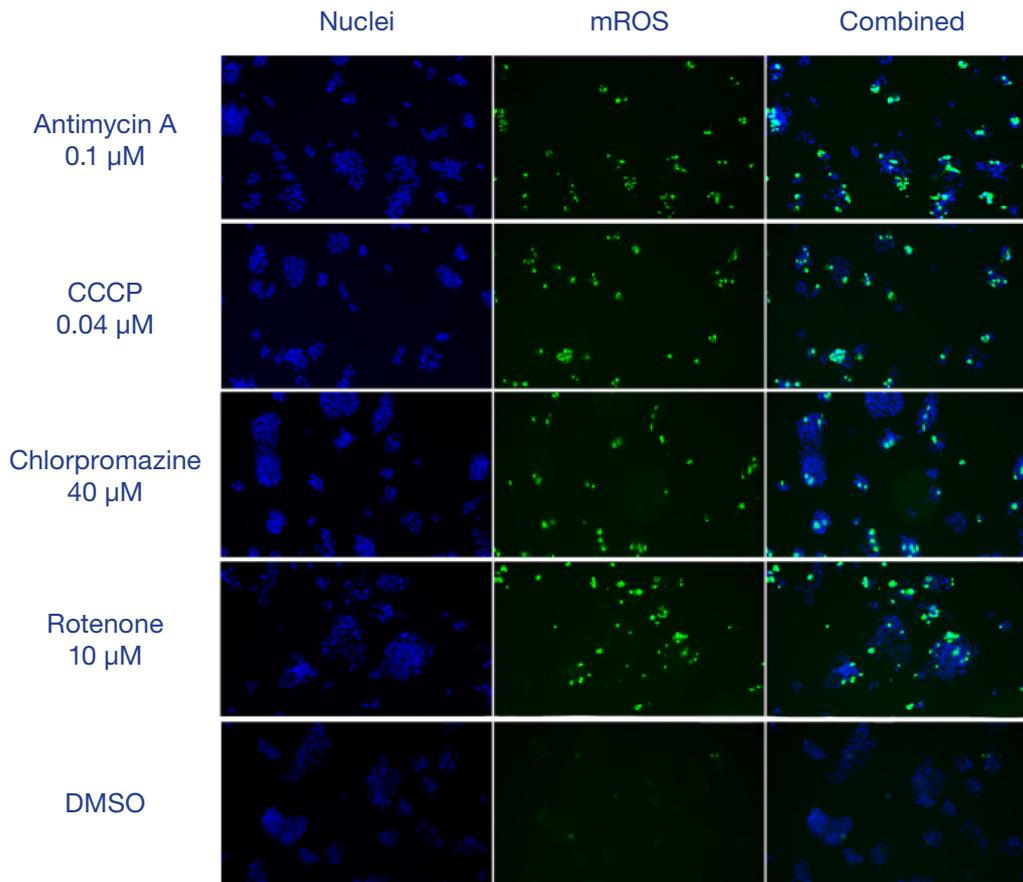
#### Data Delivery

Minimum effective concentration (MEC) and AC<sub>50</sub> value for each measured parameter, mitochondrial oxidative stress (mROS), and cell health (cell count, nuclear size, DNA structure and cellular ATP)

\*Other options available on request

**Figure 1**

Representative HCS images of four mROS inducing test compounds alongside vehicle control; nuclei (blue), mROS (green).

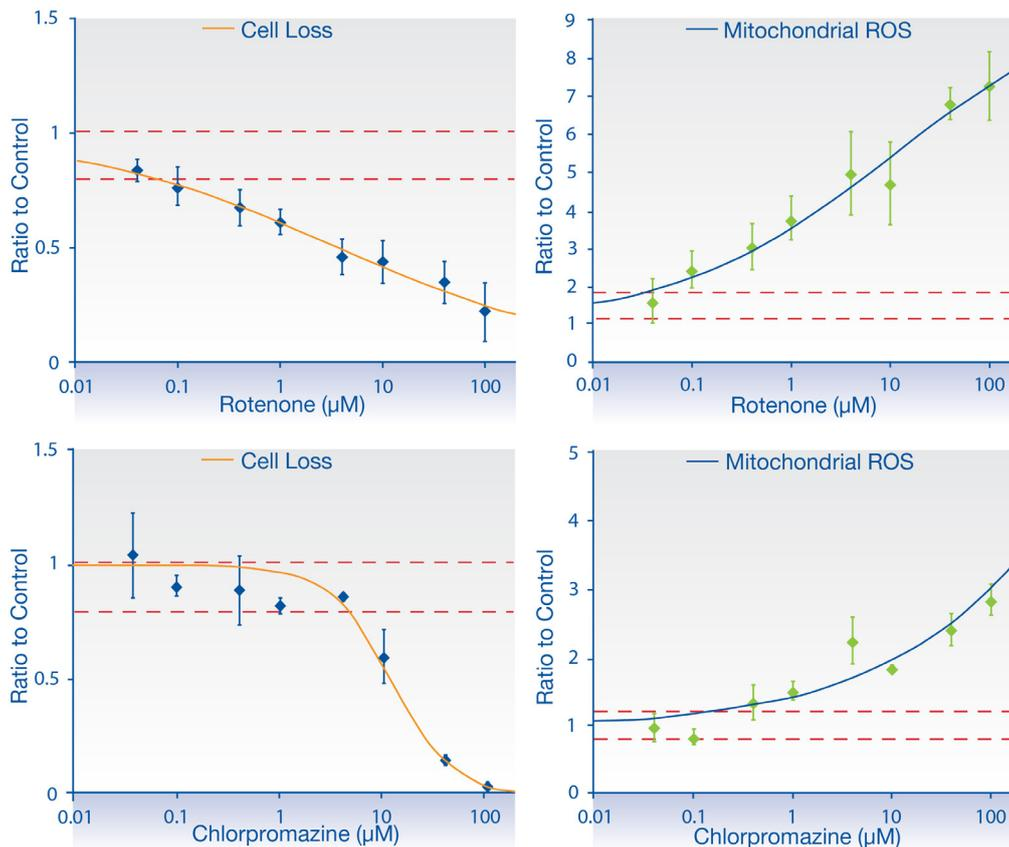


HepG2 cells were plated on tissue culture treated black walled clear bottomed polystyrene plates. The cells were dosed with test compound at a range of concentrations.

At the end of the incubation period (24 hours), the cells were labelled with Hoechst (nuclei) and mitoSox® (mROS) then imaged using an automated fluorescent cellular imager, CellInsight CX7 High-Content Screening (HCS) Platform (Thermo Scientific Celloomics).

**Figure 2**

Representative dose-response curves for rotenone and chlorpromazine displaying increased mROS formation.



#### References

- Zorov DB *et al.*, (2014) Mitochondrial Reactive Oxygen Species (ROS) and ROS-Induced ROS Release. *Physiol Rev* **94**(3); 909-950
- Ježek P *et al.*, (2020) Redox Signaling from Mitochondria: Signal Propagation and its Targets. *Biomolecules* **10**(1); 93

Read online:

