

## Introduction

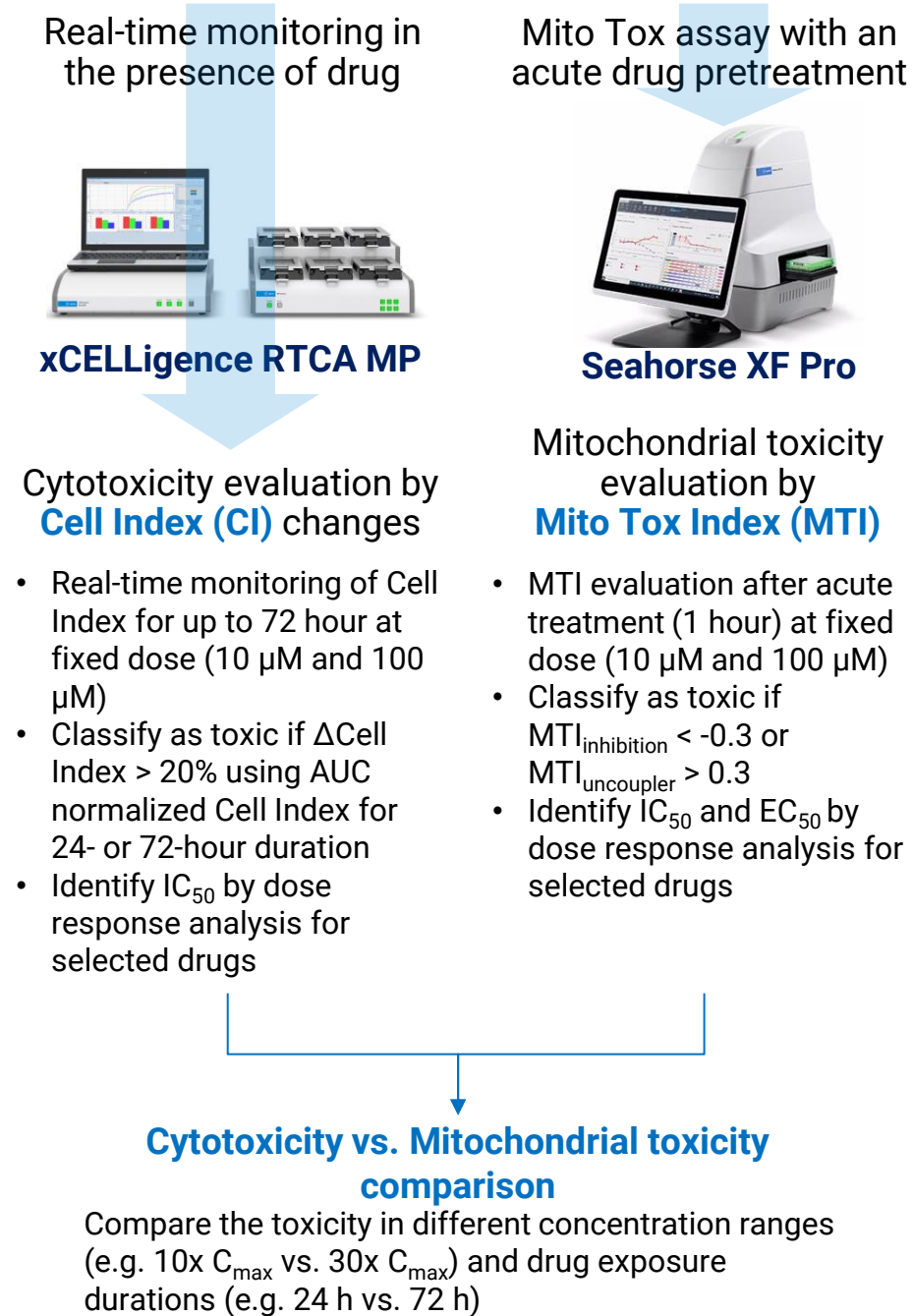
Label-free measurement of cellular impedance in Cell Index using Agilent xCELLigence RTCA provides insight into the cytotoxic effects of drugs or compounds via real-time monitoring of cell status, including cell number, viability, and morphology.

The Mito Tox Index (MTI) measured by Agilent Seahorse XF Pro analyzer quantifies the extent to which substances disrupt mitochondrial function by distinguishing between ETC inhibition and uncoupling.

This study aims to optimize and validate a drug safety testing workflow that integrates mitochondrial toxicity assessment using MTI with cell viability assessment via the Cell Index. Fourteen drugs, known to cause varying degrees of hepatotoxic adverse effects, with or without impacting mitochondrial function, were tested.

The cytotoxicity of the drugs was quantitatively evaluated using the Cell Index, while mitochondrial toxicity was assessed using an MTI threshold of 0.3, both in the HepG2 cell line. Drugs were then categorized based on the type of toxicity determined by these two sets of data.

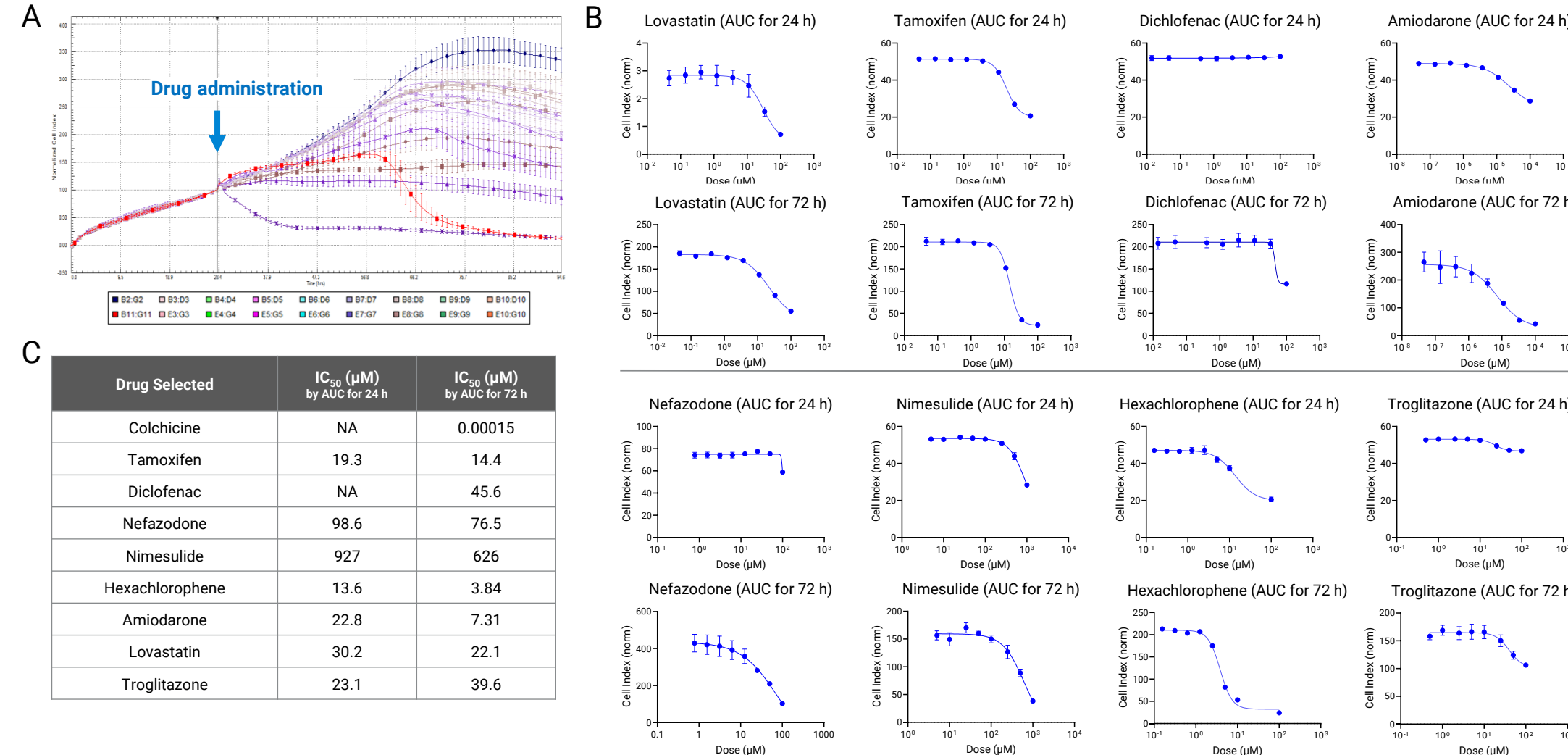
## Assay Design



**Figure 1. Toxicity evaluation workflow using dual cell analysis platforms, Agilent xCELLigence RTCA MP and Seahorse XF Pro**

## Results and Discussion

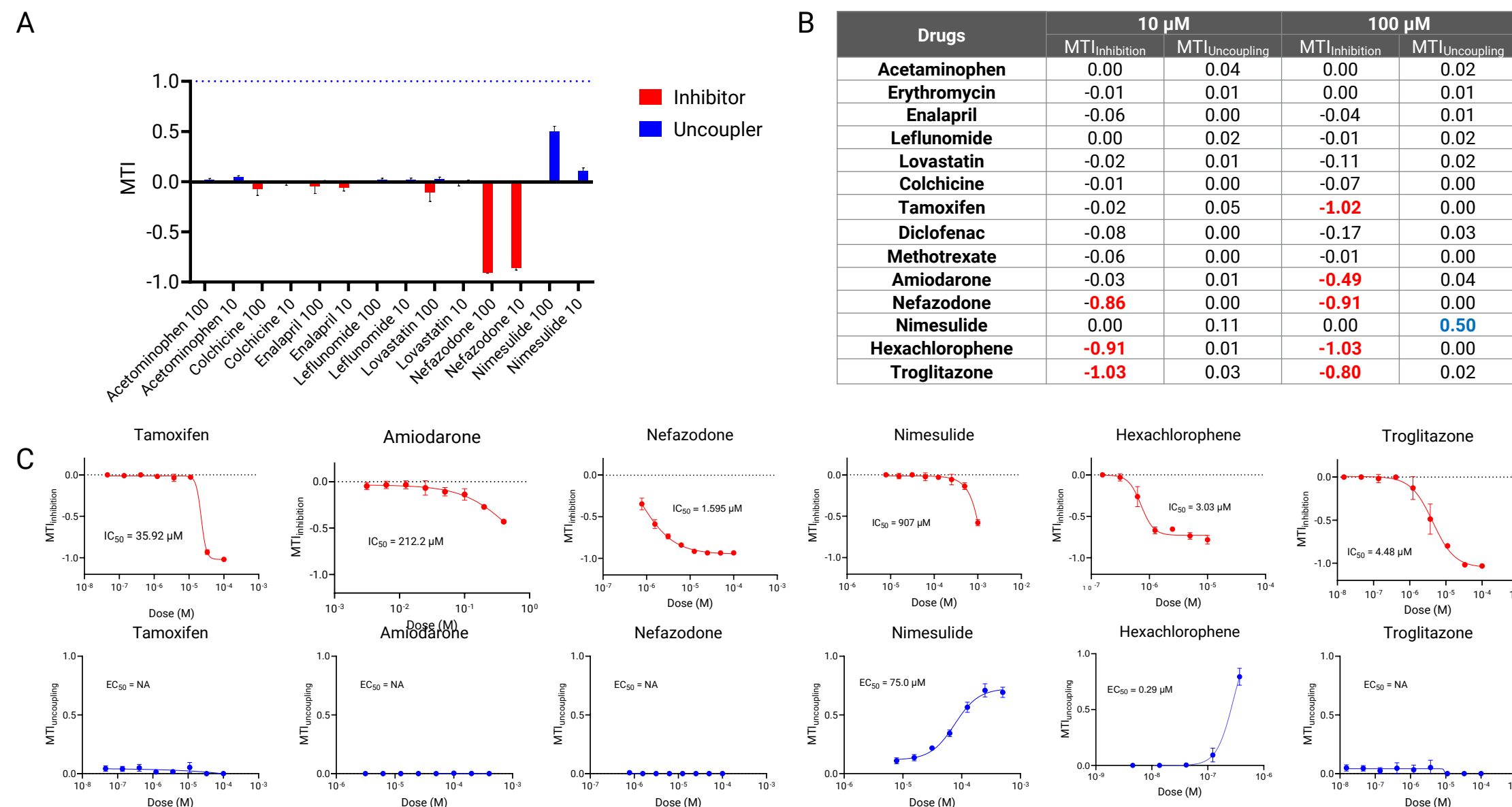
### Cytotoxicity Evaluation



**Figure 2. Cytotoxicity evaluation by Cell Index comparison**

A. A representative changes in Cell Index of HepG2 cells measured for up to 72 hours in the presence of drugs at 10 or 100  $\mu$ M.  
B. Dose-response of normalized Cell Index in AUC for the indicated time to drugs selected  
C.  $IC_{50}$  summary of the drugs selected.

### Mitochondrial Toxicity Evaluation



**Figure 3. Mitochondrial toxicity evaluation by MTI comparison**

A. A representative MTI data from HepG2 cells acutely treated for 1 hour with drugs at 10 or 100  $\mu$ M.  
B. Summary of  $MTI_{inhibition}$  and  $MTI_{uncoupling}$  for drugs tested.  
C. Dose-response of MTI to drugs show  $|0.3|$  or higher at 100  $\mu$ M.

## Summary

### Toxicity Evaluation Results

Toxicity at 30x $C_{max}$				Toxicity at 10x $C_{max}$			
Drugs	MitoTox <sup>1)</sup>	CytoTox <sup>2)</sup>		Drugs	MitoTox <sup>1)</sup>	CytoTox <sup>2)</sup>	
Name	MTI  > 0.3	$\Delta AUC_{24h}$ > 20%	$\Delta AUC_{72h}$ > 20%	Name	MTI  > 0.3	$\Delta AUC_{24h}$ > 20%	$\Delta AUC_{72h}$ > 20%
Acetaminophen	NA	NA	NA	Acetaminophen	No	No	No
Erythromycin	NA	NA	NA	Erythromycin	No	No	No
Enalapril	No	No	No	Enalapril	No	No	No
Leflunomide	No	No	No	Leflunomide	No	No	No
Lovastatin	No	No	Yes	Lovastatin	No	No	No
Colchicine	No	No	Yes	Colchicine	No	No	No
Tamoxifen	No	No	Yes	Tamoxifen	No	No	No
Diclofenac	No	No	Yes	Diclofenac	No	No	Yes
Methotrexate	No	No	Yes	Methotrexate	No	No	Yes
Amiodarone	Yes	Yes	Yes	Amiodarone	No	Yes	Yes
Nefazodone	Yes	Yes	Yes	Nefazodone	Yes	No	Yes
Nimesulide <sup>3)</sup>	Yes	Yes	Yes	Nimesulide <sup>3)</sup>	Yes	No	Yes
Hexachlorophene <sup>3)</sup>	Yes	Yes	Yes	Hexachlorophene <sup>3)</sup>	Yes	Yes	Yes
Troglitazone	Yes	Yes	Yes	Troglitazone	Yes	Yes	Yes

1) Classified as toxic if  $\Delta$ Cell Index > 20% using AUC normalized Cell Index for 24 or 72-hour duration  
2) Classified as toxic if  $MTI_{inhibition} < -0.3$  or  $MTI_{uncoupler} > 0.3$   
3) Uncoupler

Nontoxic  
No evidence of mitochondrial toxicity  
Attention required for mitochondrial toxicity  
Toxic

### Impact of mitochondrial toxicity

Drug Selected	Cytotoxicity		Mitochondrial Toxicity	
	$IC_{50}$ ( $\mu$ M, 24 h)	$IC_{50}$ ( $\mu$ M, 72 h)	MTI $IC_{50}$ ( $\mu$ M)	MTI $EC_{50}$ ( $\mu$ M)
Tamoxifen	19.3	14.4	39.9	NA
Amiodarone	22.8	7.31	212	NA
Nefazodone	98.6	76.5	1.60	NA
Nimesulide	927	626	907	75.0
Hexachlorophene	13.6	3.84	3.03	0.29
Troglitazone	23.1	39.6	4.48	NA

## Conclusion

By collecting data from both Agilent xCELLigence and Seahorse XF technologies, a rich dataset is generated that goes beyond simple cytotoxicity measures.

Cellular impedance measurements provide insights into cell viability, proliferation, and morphological changes in real-time. The ability to track these changes over time results in sensitive and comprehensive data sets.

Mitochondrial toxicity assessment offers quantitative mechanistic information, which may be informative for decisions and predictions on whether drug candidates should or shouldn't be removed from a drug pipeline.

This study highlights the complexity of drug-induced cytotoxicity, particularly in identifying those associated with mitochondrial dysfunction, and the importance of considering both dose and time-dependent effects.

In vitro models using cell lines like HepG2 have limitations in accurately predicting liver toxicity and are often insufficient to be physiologically relevant. The enriched data from dual platforms may offer more comprehensive insights into this complexity. Further research with optimized workflows and improved data integration methods is anticipated to enhance these predictions.

## References

- Huang *et al.* (2021) Assessing Water Cytotoxicity with an Impedance-Based, Real-Time, and Label-Free Cellular Assay. Agilent Technologies, application note, publication number 5994-3712EN
- Kam *et al.* (2022) A Customized XF Workflow for Detection and Characterization of Mitochondrial Toxicity. Agilent Technologies, application note, publication number 5994-4778EN
- Schulz, *et al.* (2012) Therapeutic and toxic blood concentrations of nearly 1,000 drugs and other xenobiotics. Critical Care 16:R136