

- ✓ Impedance
- ✓ High-throughput (96 well plate)
- ✓ Long-term measurements
- ✓ Extracellular field potential (EFP)
- ✓ Electrical and optical stimulation
- ✓ Force-frequency relationship
- ✓ Label-free
- ✓ Incubation chamber
- ✓ 2D and 3D culture

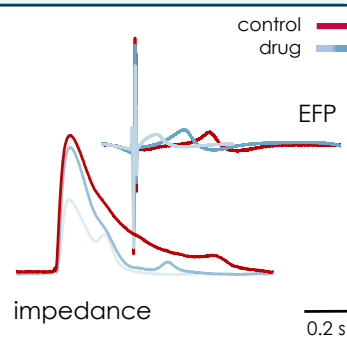


## Cardiac applications:

- Drug screening: contractility and electrophysiology readout
- Test acute and chronic drug effect
- Induce cardiomyocyte maturation through long-term pacing
- Investigate compound-induced cardiotoxicity



*hiPSC-CMs, hESC-CMs, neonatal mice and rat CMs*

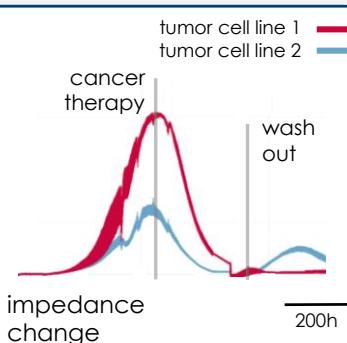


## Oncology applications:

- Monitor cancerous cell proliferation and viability
- Monitor regrowth of tumor cells after treatment
- Investigate cytotoxicity effect of cancer therapies



*Mammary carcinoma cell line (H8N8)*

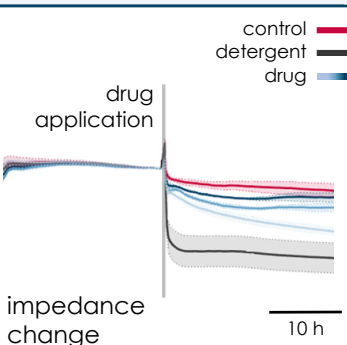


## Hepatology applications:

- Investigate compound-induced hepatotoxicity
- Monitor membrane integrity and tight junctions



*Monocyte-derived hepatocyte-like (MH): MetaHeps®  
Liver sinusoidal endothelial cells and hepatocytes: Upcyte®*



## Other applications:

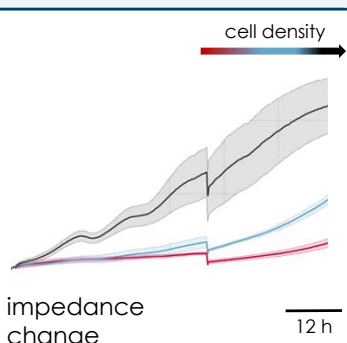
- Monitor proliferation, shape change and death of several cell types, induced not only by test reagents but for example by photoswitchable compounds or hypoxia



*HeLa cells*



*Chinese hamster ovary (CHO) cells*



...and find out more