

>> Standards for hiPSC-derived cardiomyocyte electrophysiology using the MEA assay

Mike Clements^{1,✉}, Cristina Altrocchi², Devon Guerrelli³, Adam P. Hill⁴, Nikki Posnack³, Chris Strock⁵, & Takashi Yoshinaga⁶

¹Axion BioSystems, GA, USA; ²Johnson & Johnson, Belgium; ³Children's National Hospital, DC, USA; ⁴Victor Chang Cardiac Research Institute, Australia; ⁵Cyprotex, MA, USA; ⁶Eisai Co., Ltd, Japan. ✉ Email: mclements@axionbio.com

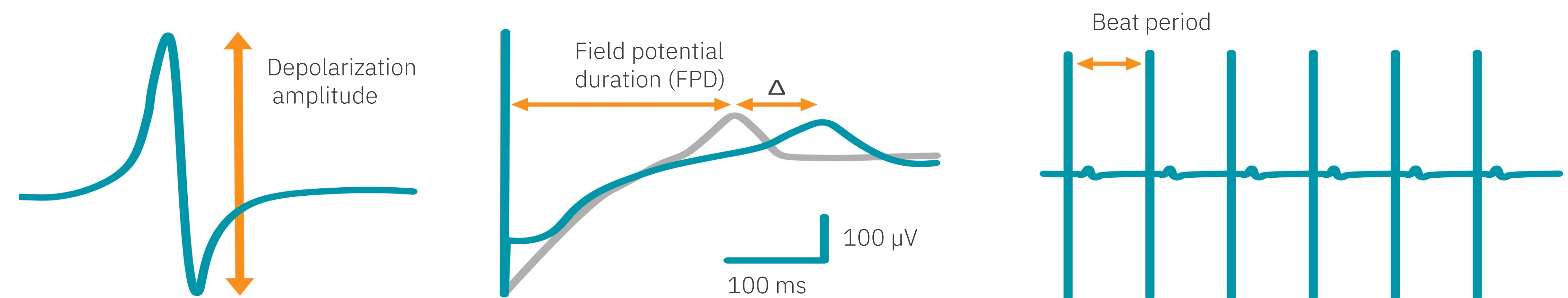
Advancing hiPSC-CM assay adoption will require standardization for data comparison and validation

Background

Cardiac stem cell models are transforming research and discovery—but a **lack of standardized criteria** to assess functional activity can lead to inconsistent results. Over the last decade, the multielectrode array (MEA) assay has become a popular tool for characterizing hiPSC-cardiomyocyte (CM) batches, studying disease models, screening therapeutics, and evaluating drug-induced cardiotoxicity.

The goal of this project is to **set the minimum acceptance criteria** for a spontaneous beating wild type hiPSC-ventricular cardiomyocyte field potential assay **for compound testing and/or disease modeling**.

Identifying key cardiac metrics



Leveraging in-house experience in academia and industry, published data, and international consortia (CiPA and JiCSA) validating hiPSC-cardiomyocyte assays, we have developed the **Axion iPSC Model Standards (AIMS)** framework. This proposed standard focuses on the **spontaneous beat rate**, features of the cardiac waveform (**depolarization spike amplitude** and **field potential duration**), and the synchronization of activity in the syncytia.

AIMS #CM01: Defining minimal acceptance criteria for wild-type hiPSC-ventricular cardiomyocyte

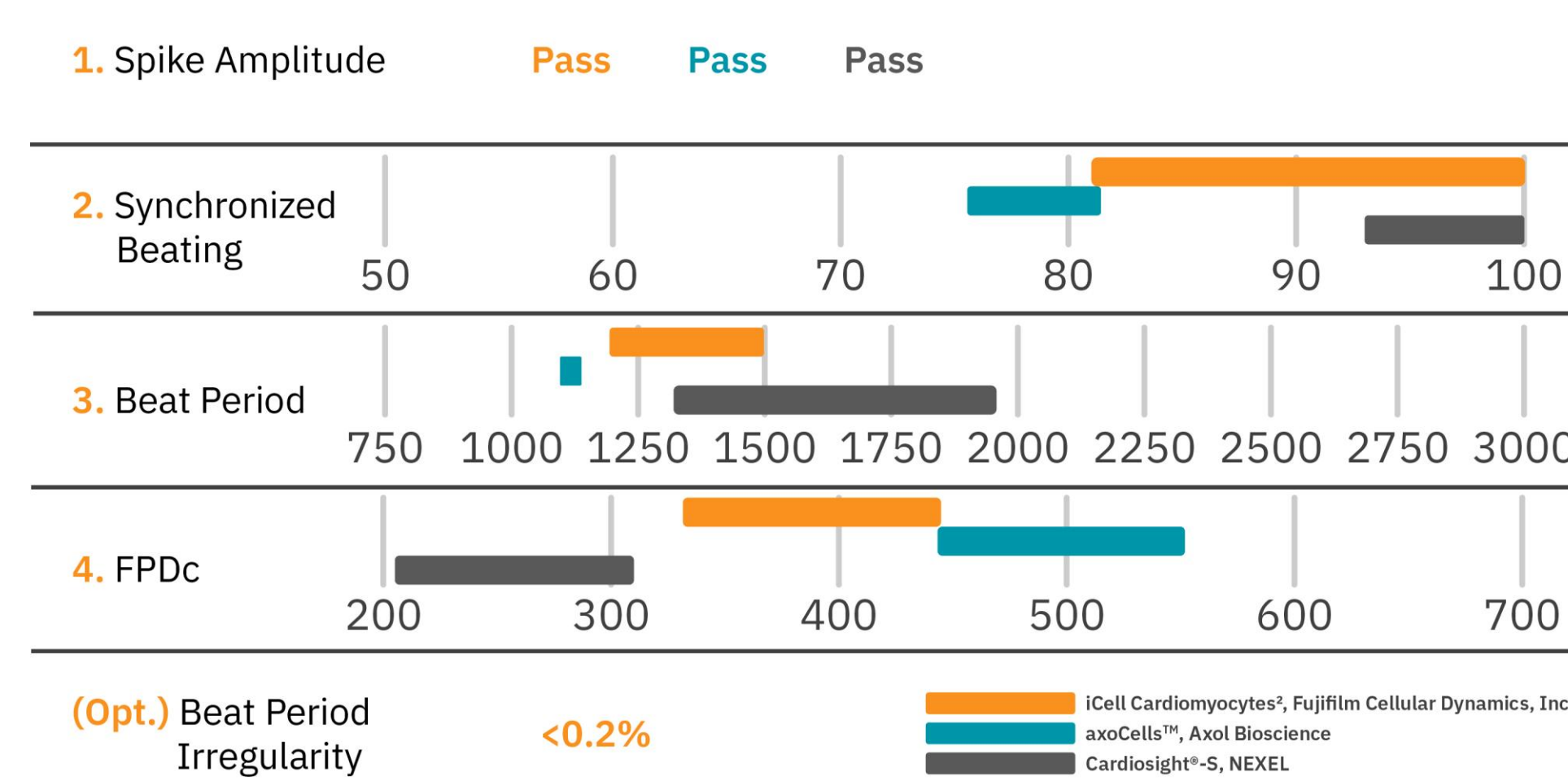
AIMS #CM01: Proposed standards

An hiPSC-CM cell source is considered to meet standards when **≥80% of MEA wells** meet all four of the following well acceptance criteria:

1. Spike amplitude of **≥0.5 mV** on ≥50% electrodes .
2. Synchronized beating across **≥50% electrodes**.
3. Spontaneous beat period of **750-3000 ms** (i.e. 20-80 beats per minute, BPM).
4. Corrected field potential duration (FPDc) of **200-700 ms**. Fridericia correction: $FPDc = FPD / (\text{Beat Period})^{1/3}$

(Optional) Beat period irregularity **<0.2%**

Standards accommodate a range of cell sources



This figure shows three vendor sources where the specification was met. Specification range was inferred from Millard et. al. 2018¹. Although basal activity of vendor cell sources differed significantly, **consistent concentration-dependent effects were observed**¹.

Important considerations

- >> Reproducibility and quality is dependent on culture protocols. Participating vendor protocols must specify key details (see AIMS website).
- >> The synchronized beating should initiate from a single point of origin. Competing pacemakers can confound analysis
- >> For compound-induced effects on spike amplitude, ≥1.0 mV would provide a larger assay window.
- >> Primary ventricular CMs do not spontaneously beat, and this behavior in iPSC-CMs is often attributed to “immaturity.” As models develop, a paced hiPSC-CM assay could be the subject of a future AIMS but is beyond the scope of this AIMS.

Select References

1. Millard D et al. Cross-Site Reliability of Human Induced Pluripotent stem cell-derived Cardiomyocyte Based Safety Assays Using Microelectrode Arrays: Results from a Blinded CiPA Pilot Study. *Toxicol Sci.* 2018 Aug 1;164(2):550-562.
2. Blinova K et al. International Multisite Study of Human-Induced Pluripotent Stem Cell-Derived Cardiomyocytes for Drug Proarrhythmic Potential Assessment. *Cell Rep.* 2018 Sep 25
3. Cooper B et al. Comparative cardiotoxicity assessment of bisphenol chemicals and estradiol using human induced pluripotent stem cell-derived cardiomyocytes. *Toxicol Sci.* 2024 Apr 2;198:273-287
4. Thorpe, J et al. Development of a robust induced pluripotent stem cell atrial cardiomyocyte differentiation protocol to model atrial arrhythmia. *Stem Cell Res Ther* 14, 183 (2023).
5. Clements M et al. High-Throughput Multi-Parameter Profiling of Electrophysiological Drug Effects in Human Embryonic Stem Cell Derived Cardiomyocytes using Multi-Electrode Arrays. *Toxicol Sci.* 2014 Aug 1;140(2):445-61.

More references available at www.axionbiosystems.com/axion-ipsc-model-standards-aims/aims-cm01

Acknowledgements

Thanks to FCDI, Axol Biosciences, and NEXEL for sharing their hiPSC-cardiomyocyte electrophysiological characterization data.

Additional thanks to Daniel Millard, Stacie Chvatal, Anthony Nicolini, Parker Ellingson, and Rika Yamazaki for AIMS review.



**Learn more
about AIMS**



Scan the QR code to visit our website for more information about AIMS, iPSC model standardization, and provide feedback.