

### Introduction

the symptomatology of almost all disease areas.

discovery giving rise to better lead molecules.

simultaneous recordings for interpretation of the cellular response.



Fig. 1: Qube 384. With 384 parallel measurement sites and 10–15 minutes per plate run, the Qube enables testing of more than 1,500 compounds on different cells in one hour. With full automation features for unattended operation – a plate stacker and a cell preparation unit –, the system provides walk-away functionality for more than 6,000 compounds before user interventions is needed.

### Conclusion

Here we show experiments in voltage and current clamp modes on Qube 384 using HL-1 mouse atrial cardiomyocytes and Axiogenesis matured COR.4U iPS cell-derived cardiomyocytes. In both cell types it was possible to evoke action potentials by injecting current pulses whereas HL-1 cells also showed spontaneous action potentials. In both HL-1 and iPS CMs it was possible to evoke inward sodium currents and in a fraction of the cells also calcium currents (not shown). A distinguished feature of the Qube is the possibility to combine both current and voltage clamp recordings in the same sweep making it possible to carry out very fine-tuned experiments.

# Voltage- and current clamp on induced pluripotent cardiomyocytes with automated patch clamp

Daniel Sauter, Anders Lindqvist, Rasmus Bjørn Jacobsen, Jens Henneke, Mads P.G. Korsgaard



![](_page_0_Figure_15.jpeg)

### References

Claycomb WC, Lanson NA Jr, Stallworth BS, Egeland DB, Delcarpio JB, Bahinski A, Izzo NJ Jr. "HL-1 cells: a cardiac muscle cell line that contracts and retains phenotypic characteristics of the adult cardiomyocyte". Proc Natl Acad Sci U S A. 95:2979-84, 1998.

## Materials and methods

HL-1 cells were cultured as described in Claycomb et al. (1998). iPS-derived cardiomyocytes (CMs) were kindly provided by Axiogenesis AG (Cologne, Germany) and cultured according to instructions provided by Axiogenesis. Experiments with iPSderived CMs were carried out at 20 degrees and experiments with HL-1 at ambient temperature.

Sophion Bioscience A/S Baltorpvej 154, 2750 Ballerup, Denmark Phone: +45 4460 8800 info@sophion.com www.sophion.com

![](_page_0_Picture_23.jpeg)