

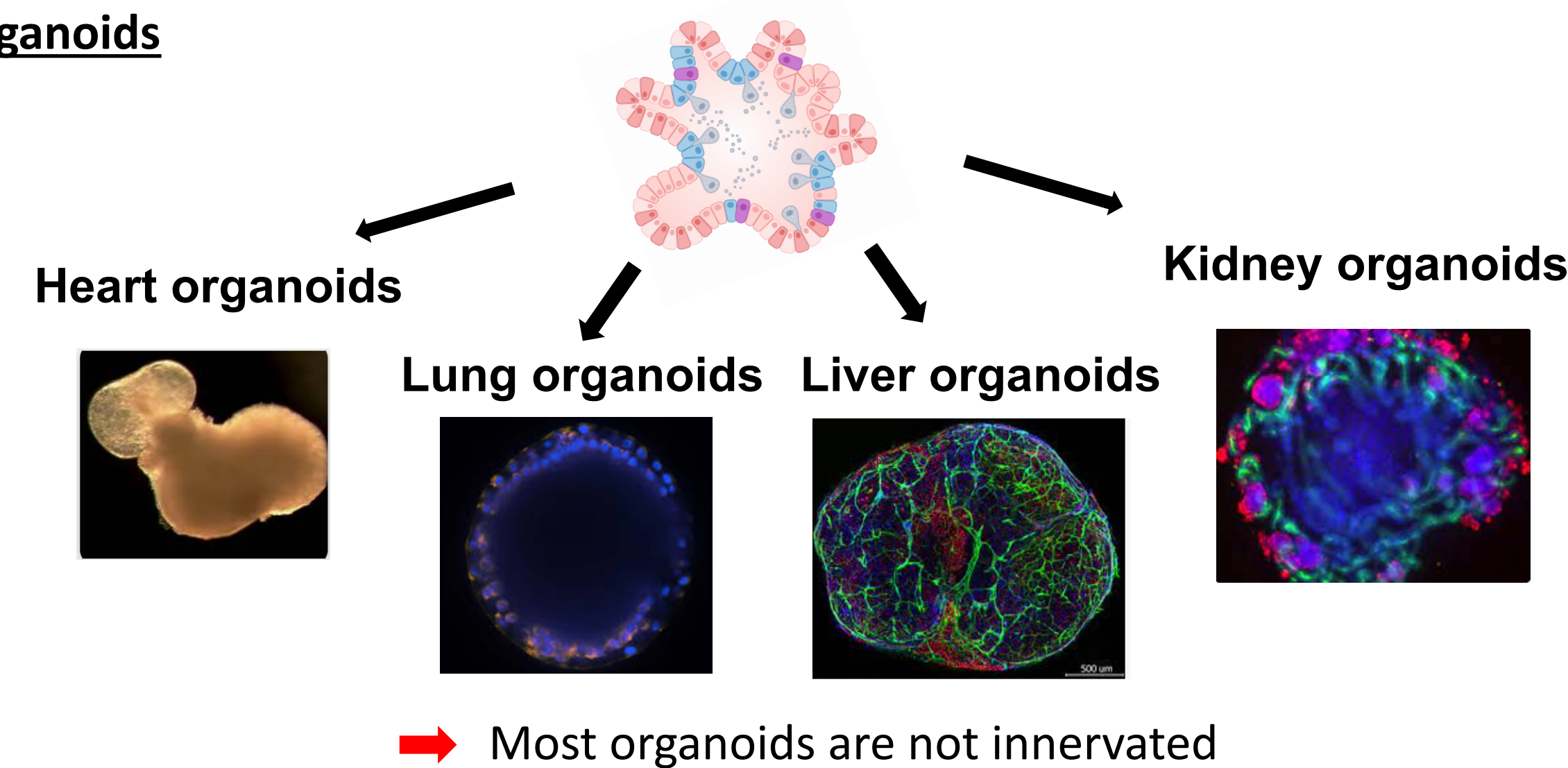


Abstract

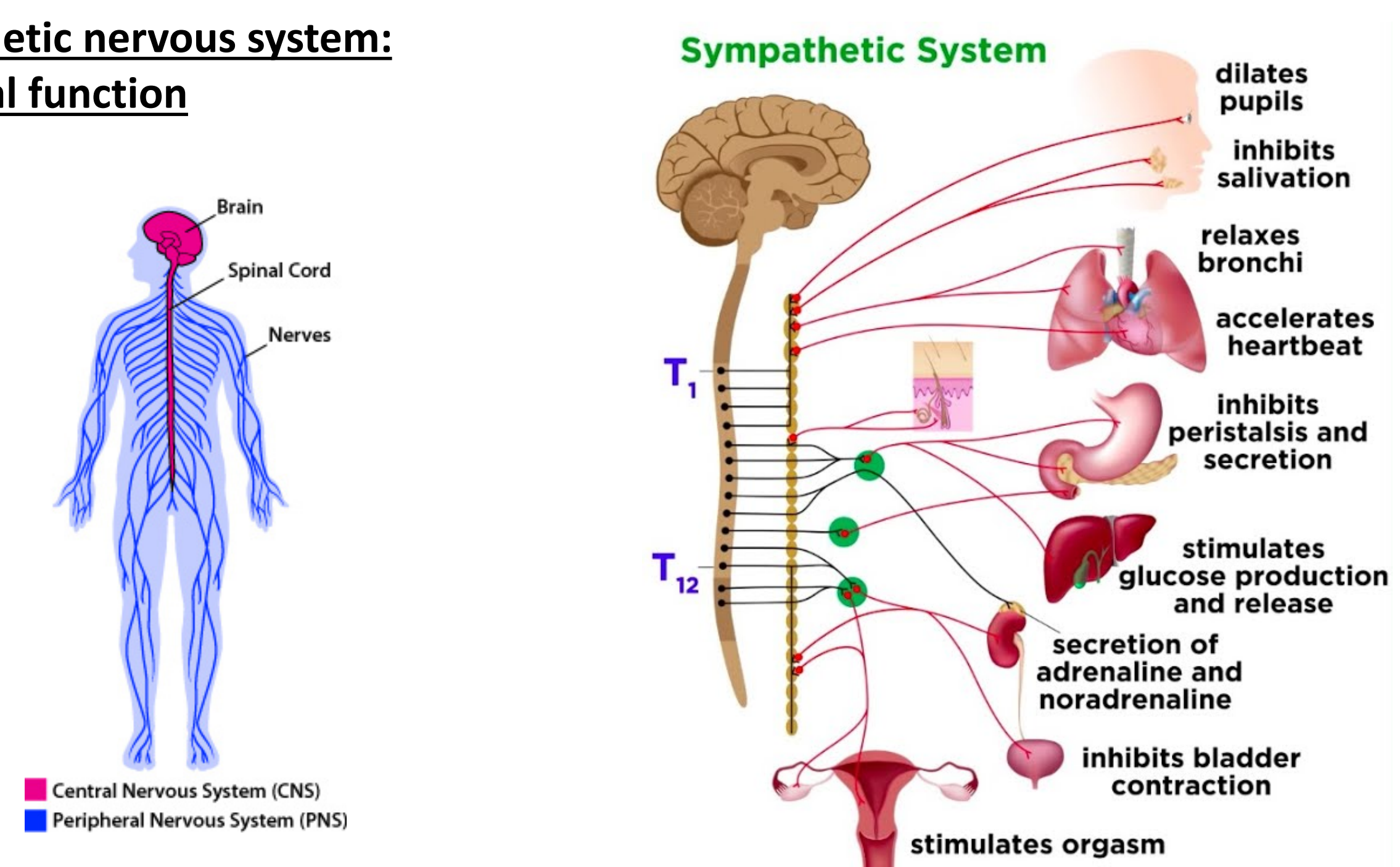
The peripheral sympathetic nervous system (SNS) innervates and regulates the maturation and functionality of almost all organs in the body. One of the most vital organs that the SNS regulates is the heart. Sympathetic neuron (symN) signaling facilitates cardiac development, maturation, and increases heartbeat. SNS dysregulation has been linked to cardiac dysfunction, such as arrhythmia and myocardial infarction. Human organoids derived from human pluripotent stem cells (hPSCs) are valuable tools in studying organ development and functionality in healthy and diseased states. However, despite a wide availability of cardiac organoid protocols, none of these organoids are symN-innervated, thus lacking the neurocardiac interaction. We have previously reported a well-established symN protocol using hPSCs, that has been applied to modeling multiple SNS diseases. Here, we developed an efficient strategy to make sympathetic neuron (symN)-innervated cardiac assembloids (hSCAs) are self-organized, and showed cardiomyocyte maturation, cardiac cavity formation, atrial to ventricular patterning, and spontaneous beating. In hSCAs, we also observed symN innervation with neurotransmitter release, and regulation of the cardiomyocyte beating rate, which could be manipulated pharmacologically or optogenetically. Using this platform, we modeled symN-mediated early heart development and myocardial infarction. This easy-to-access and highly versatile modular platform will facilitate the study of neuron-organ interaction *in vitro* and may be applicable to make more assembloid models of variable organs (such as kidney and lung) with diverse peripheral neurons (such as parasympathetic and sensory neurons).

Introduction

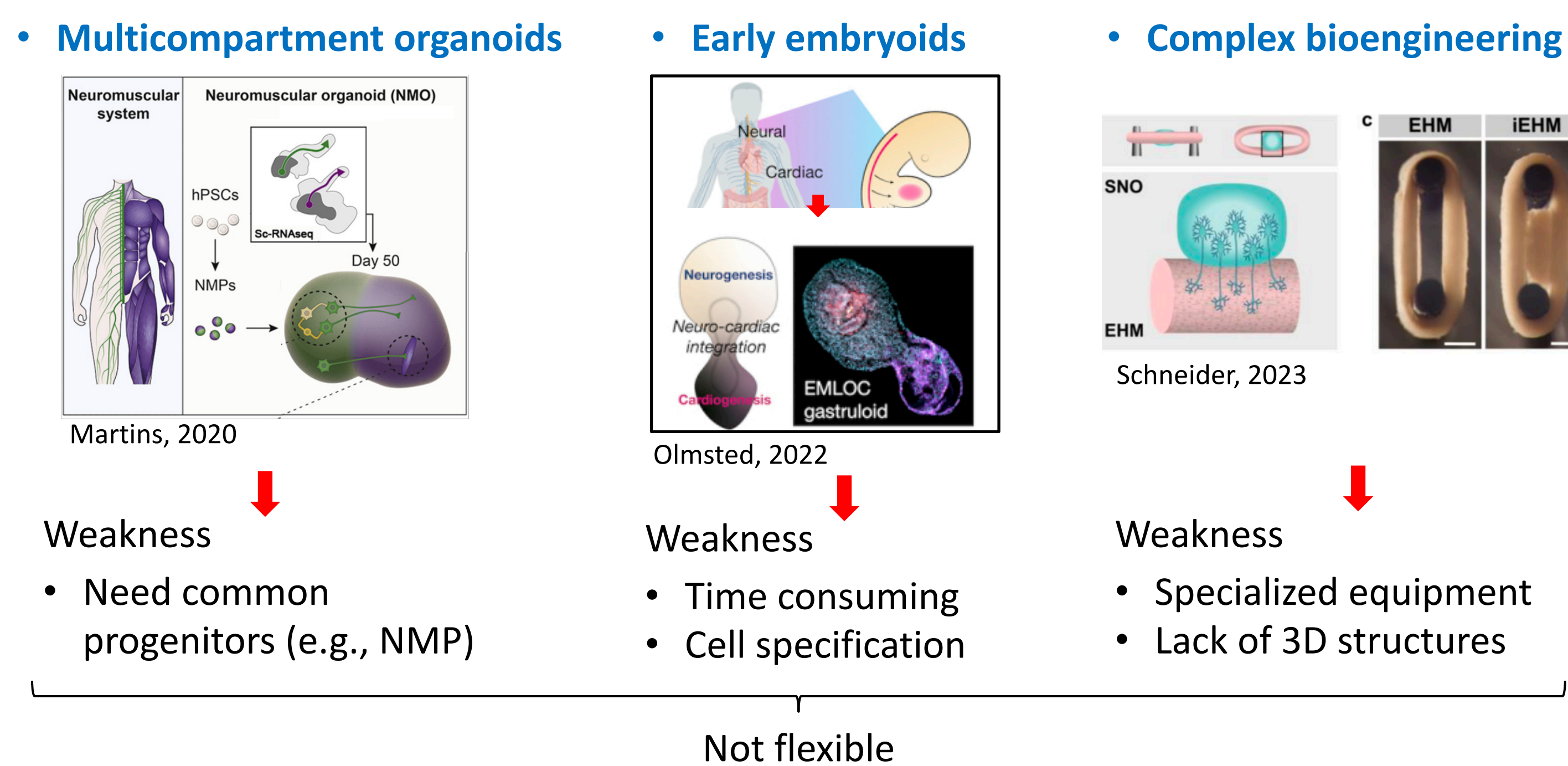
3D organoids



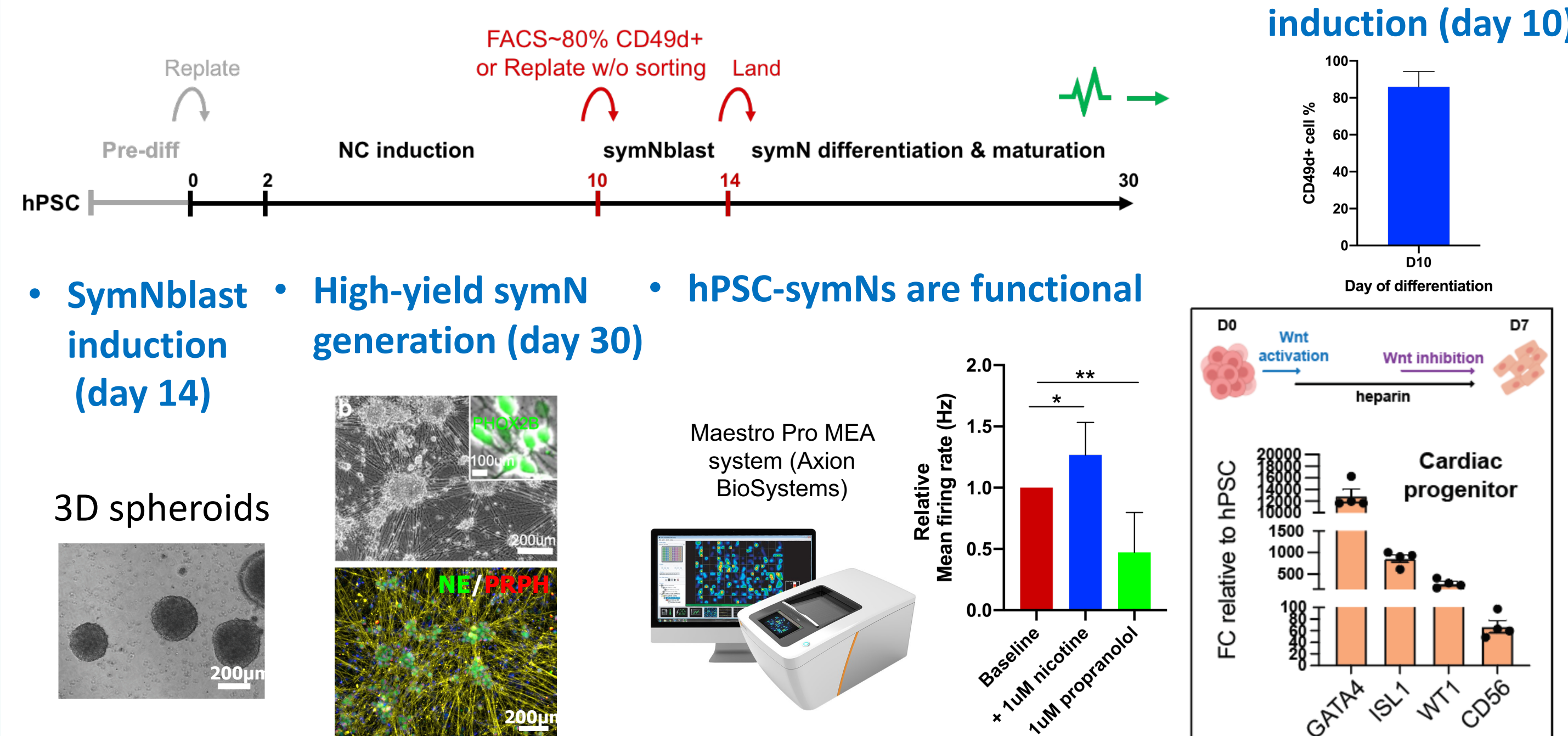
Sympathetic nervous system: biological function



Current methods to make innervated organoids

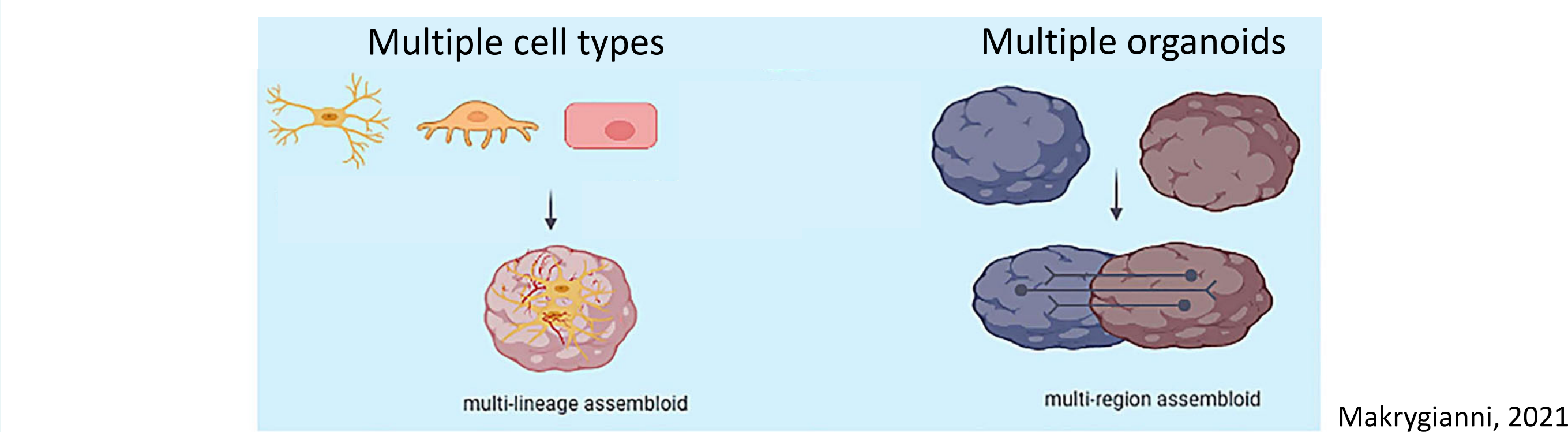


Human pluripotent stem cell-derived symNs and cardiac progenitors



Specific aims

- To develop ANS-innervated organ models
 - A simple, easy to reproduce, efficient, modular platform
- => Assembloid technology: self-organizing admixtures of ≥ 2 cell types/organoids (Kanton, 2022)



Results

Figure 1. Building the organoid using the assembloid technology

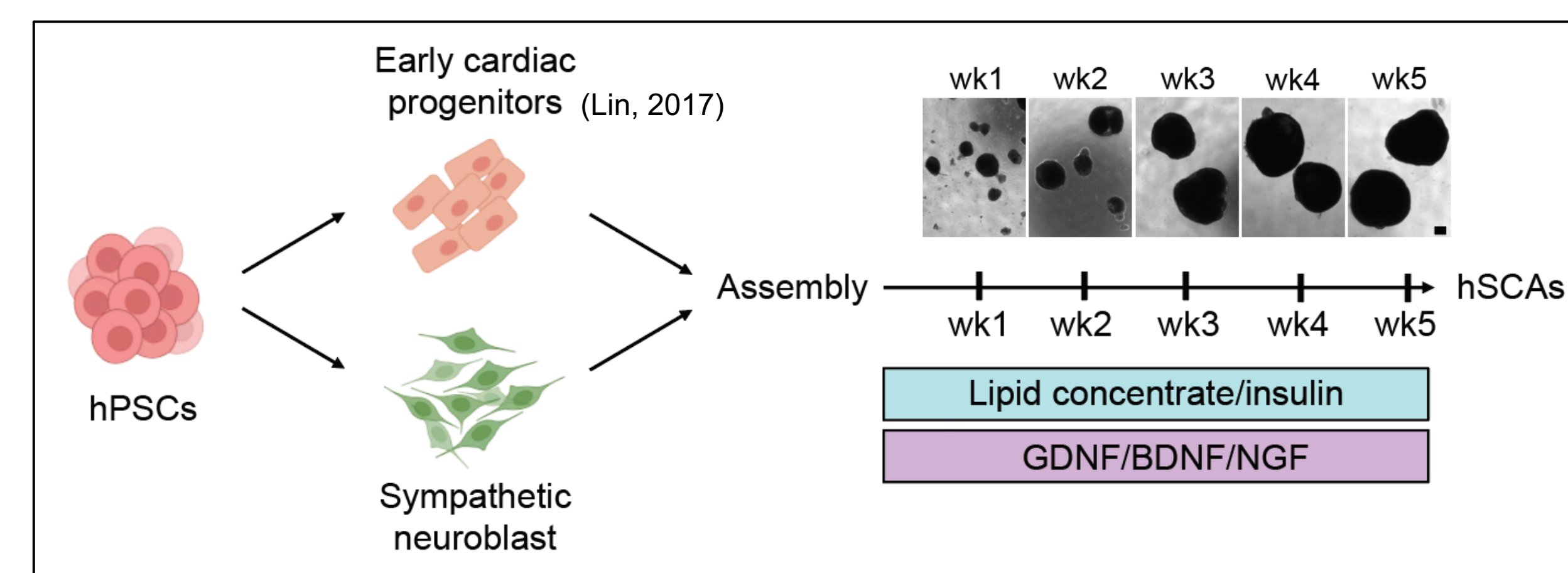
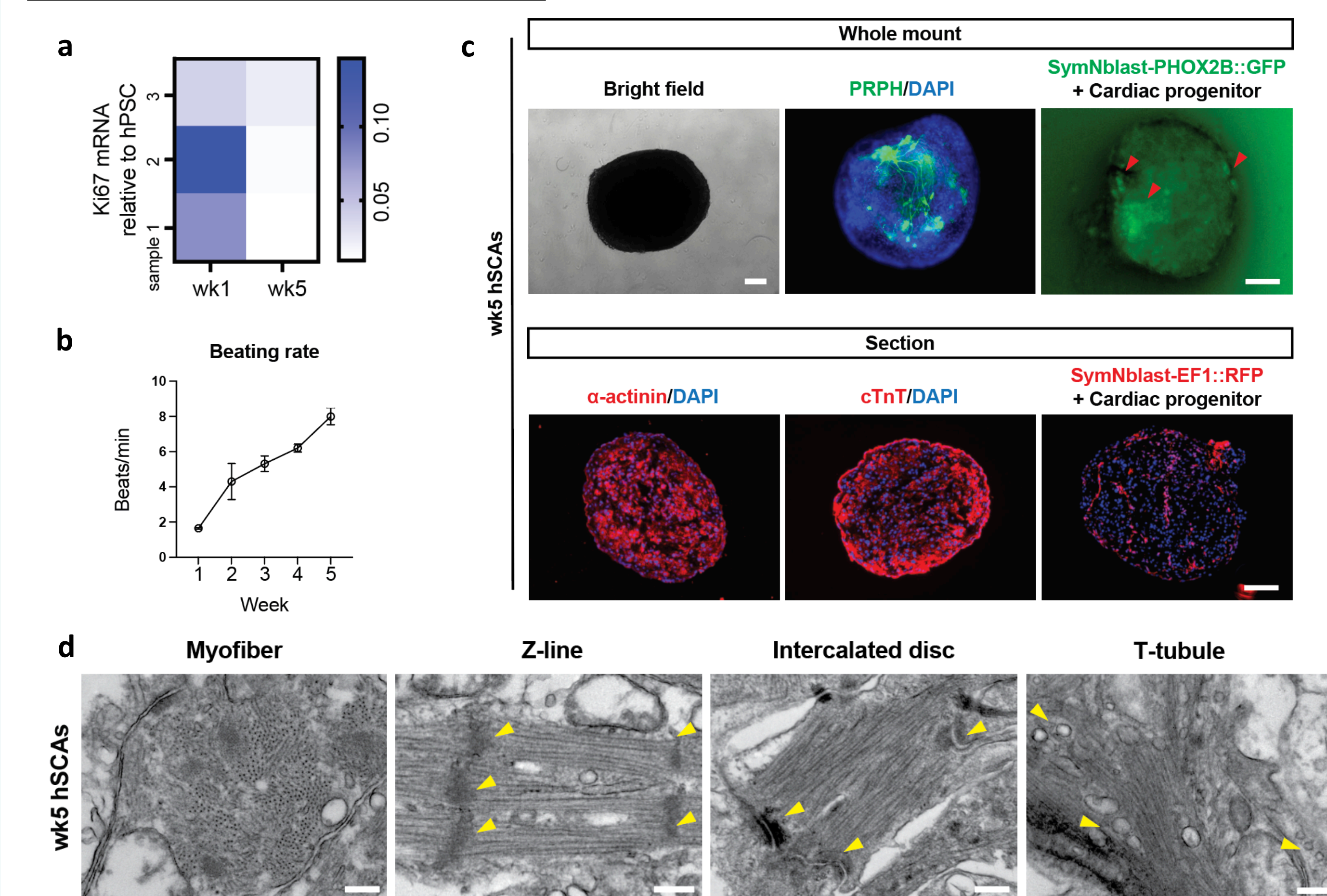


Figure 2. Characterization of hSCAs



(a) Ki67 mRNA expression. (b) hSCAs beating over time. (c) IF staining. (d) Mature features by TEM

Efficient NC induction (day 10)

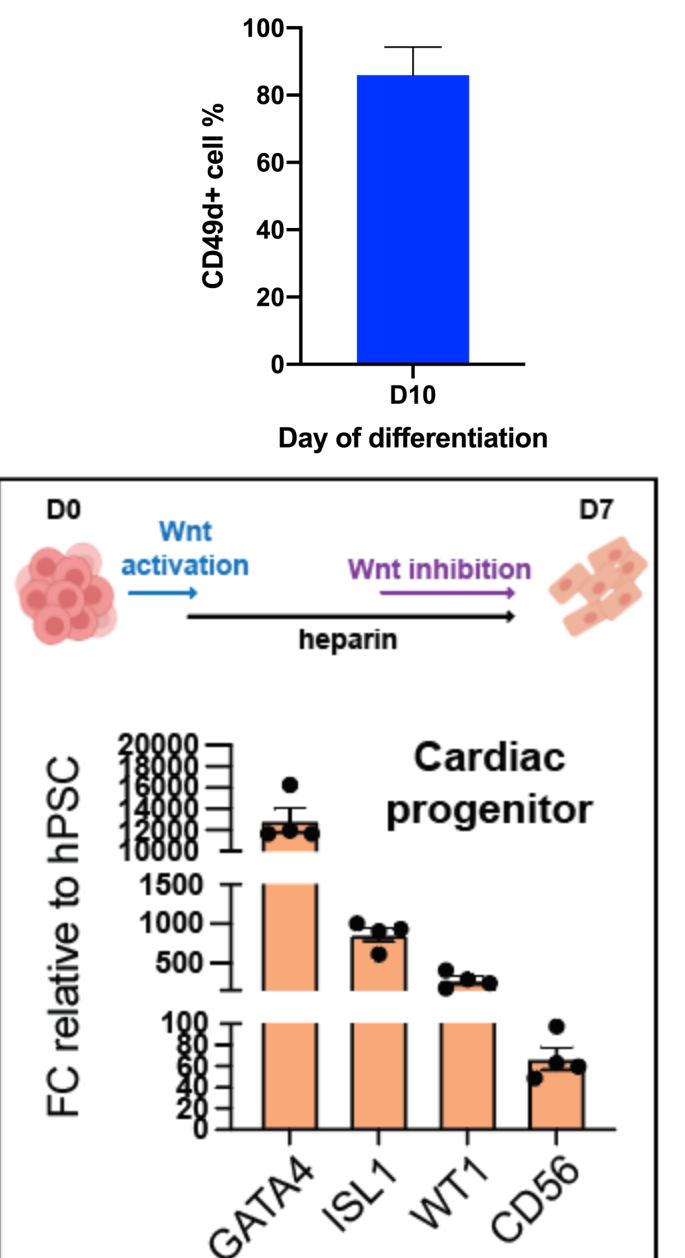


Figure 3. hSCAs are self-organized

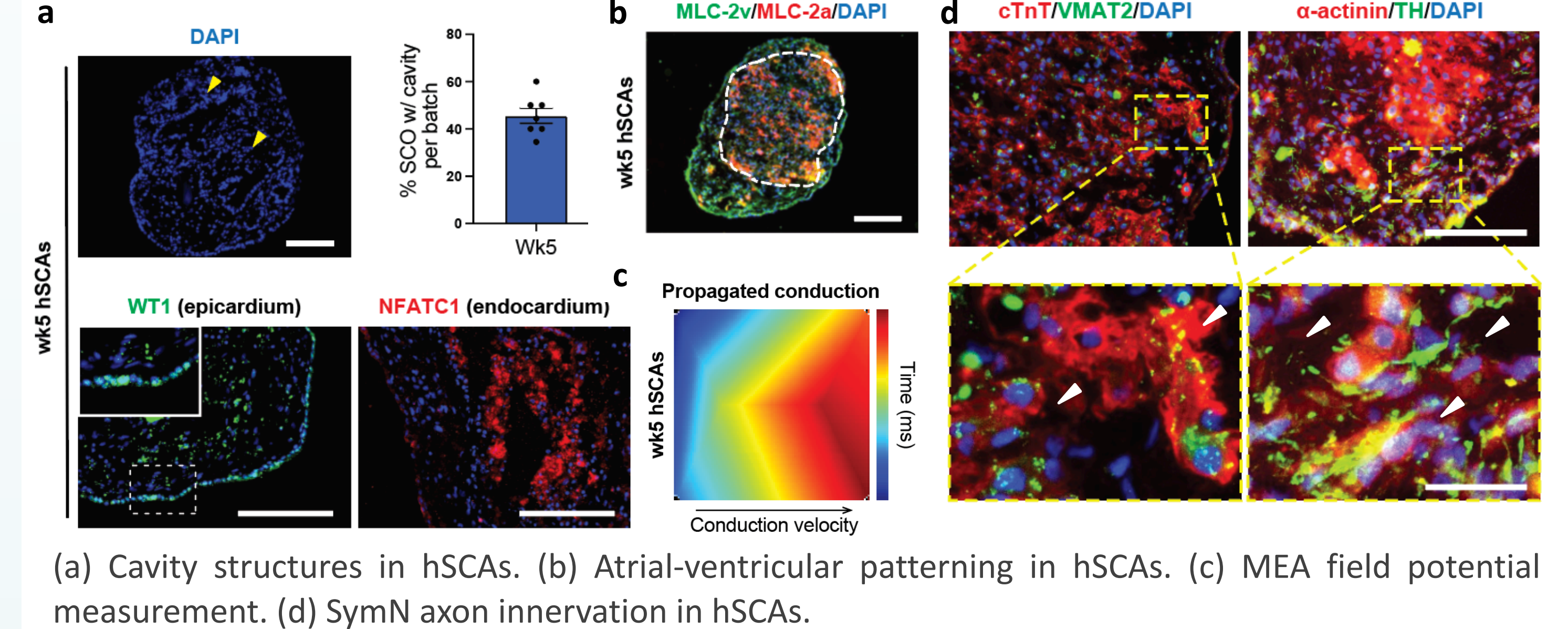
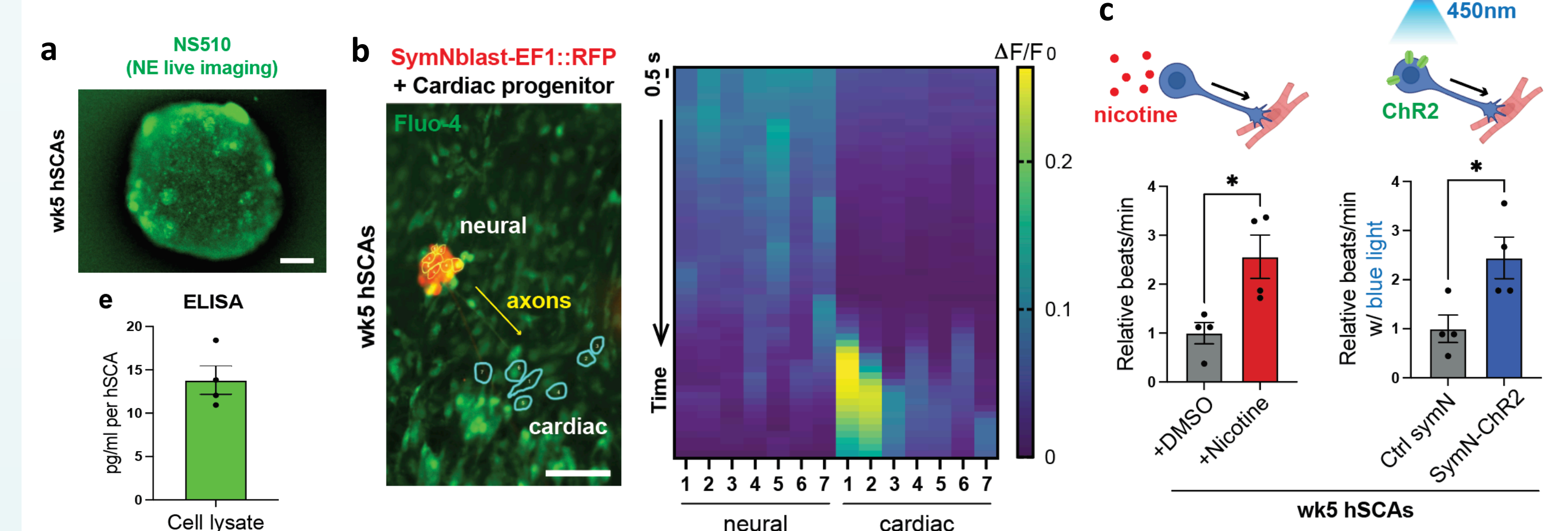
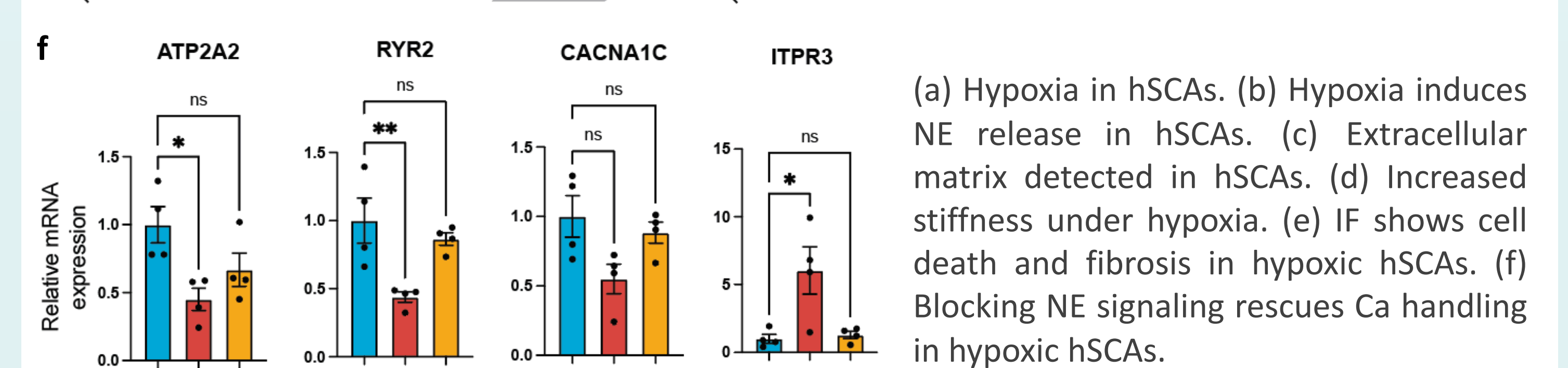
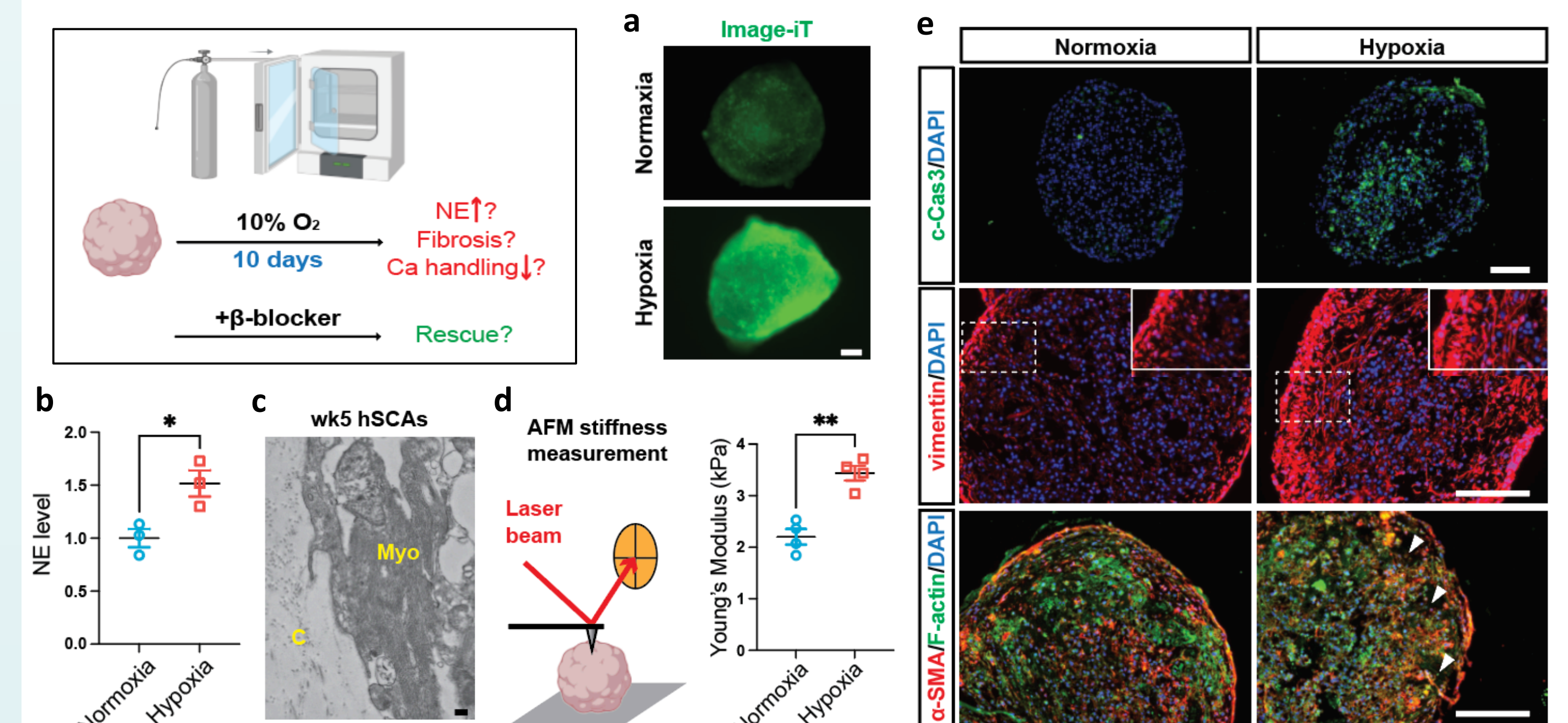


Figure 4. SymN regulation in hSCAs



(a) NE release in hSCAs. (b) SymN signaling regulates hSCA beating. (c) SymN activation increases hSCA beating.

Figure 5. hSCAs model myocardial infarction



(a) Hypoxia in hSCAs. (b) Hypoxia induces NE release in hSCAs. (c) Extracellular matrix detected in hSCAs. (d) Increased stiffness under hypoxia. (e) IF shows cell death and fibrosis in hypoxic hSCAs. (f) Blocking NE signaling rescues Ca handling in hypoxic hSCAs.

Conclusion

Here, we developed a simple and flexible strategy to generate symN-innervated cardiac assembloids from hPSCs. The hSCAs are functional, and may be applied to generate more ANS-innervated organoids in the future. Such human-based *in vitro* systems may significantly facilitate the understanding of peripheral network regulation in diseases.

References

1. Wu et al. Efficient Differentiation of Postganglionic Sympathetic Neurons using Human Pluripotent Stem Cells under Feeder-free and Chemically Defined Culture Conditions. *J. Vis. Exp.* (159), e60843, doi:10.3791/60843 (2020).
2. Wu et al. Generation of postganglionic sympathetic neurons using human pluripotent stem cells for electrophysiological and functional assessments. STAR protocol. (2024)