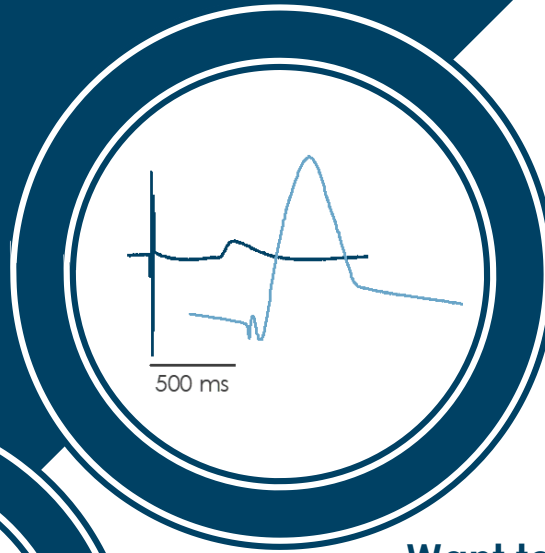


CardioExcyte 96 FLEXcyte 96

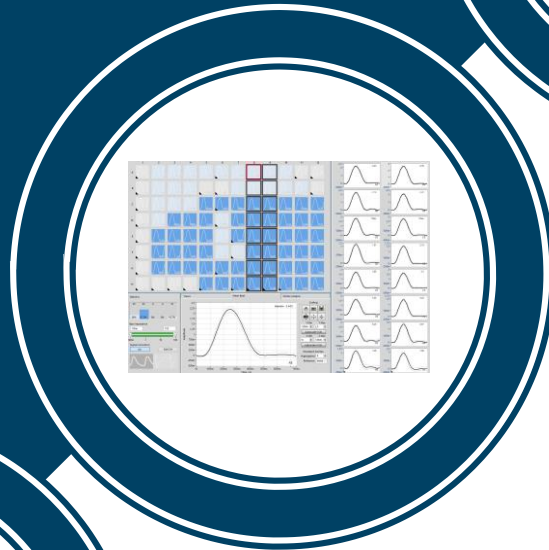
Electrophysiology
Contractility
Viability

nanjion



Want to monitor activity of your cells?

We offer sensitive, multiparametric readouts on any timescale.



Want to optimize your lab time?

We provide the software to speed up your analysis.



What to increase insight in your cells' behavior?

We provide overall solutions for your research.

ONLINE DEMOS



Book yours now!
info@nanion.de