

## Introducing the PatchXpress<sup>®</sup> Plus Automated Parallel Patch Clamp System

### Current-clamp upgrade kits now available for PatchXpress Systems

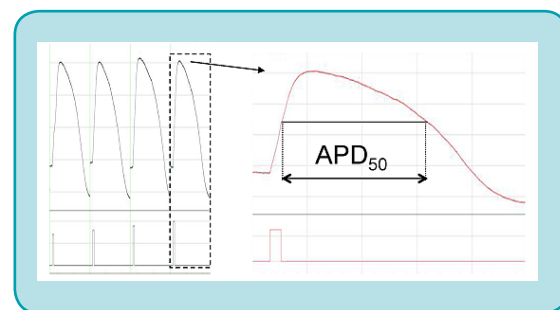
- Monitor resting membrane potentials in the absence and presence of test compounds
- Inject current to modify membrane potentials or to elicit action potentials

### Measure action potentials in excitable cells including stem-cell cardiomyocytes and neurons

- Detect action potentials during data acquisition

### Measure Action Potential Durations (APDs) in DataXpress<sup>®</sup> Analysis Software

- APDs can be measured based on user selectable parameters (APD<sub>30</sub>, APD<sub>50</sub>, etc.)



Action potential recording on the PatchXpress Plus System from nodal stem-cell cardiomyocyte.