

SyncroPatch 384

IV analysis tool

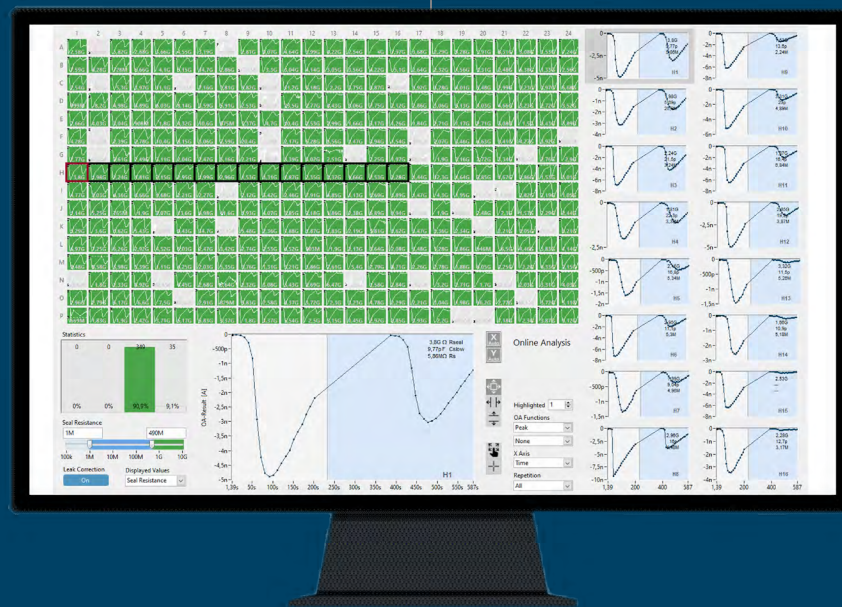
nanion

Analysis of compound effects on current-voltage relationships with DataControl 384

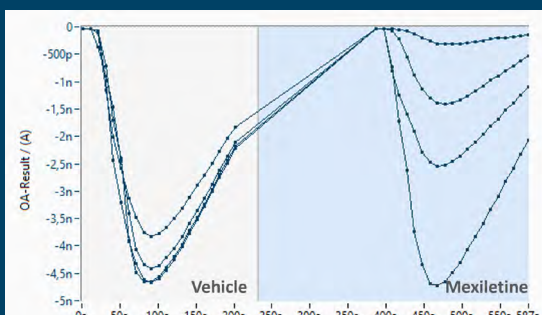
- Seamless generation of activation and inactivation curves pre- and post-compound.
- Estimate concentration response relationships of compound effects on e.g. V_{half} or G_{max} .
- Generate IV curves from averaged data from multiple wells.

Get the most out of your data

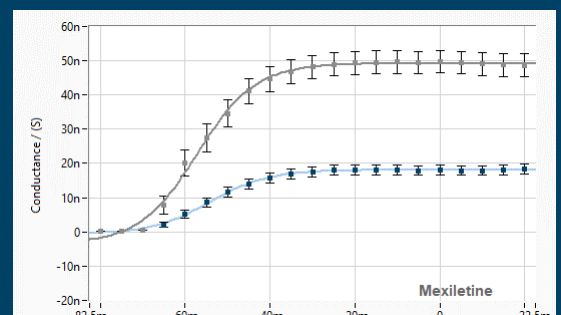
- Normalize and overlay IV curves to easily spot effects or differences.
- Use heatmaps for quick overview of IV derived parameters.
- Use IV data as QC parameter for automated filtering.



www.nanion.de



Average IV plots recorded in different conditions can be displayed – here 3 different concentrations of mexiletine (blue) and the corresponding vehicle (grey) are displayed.

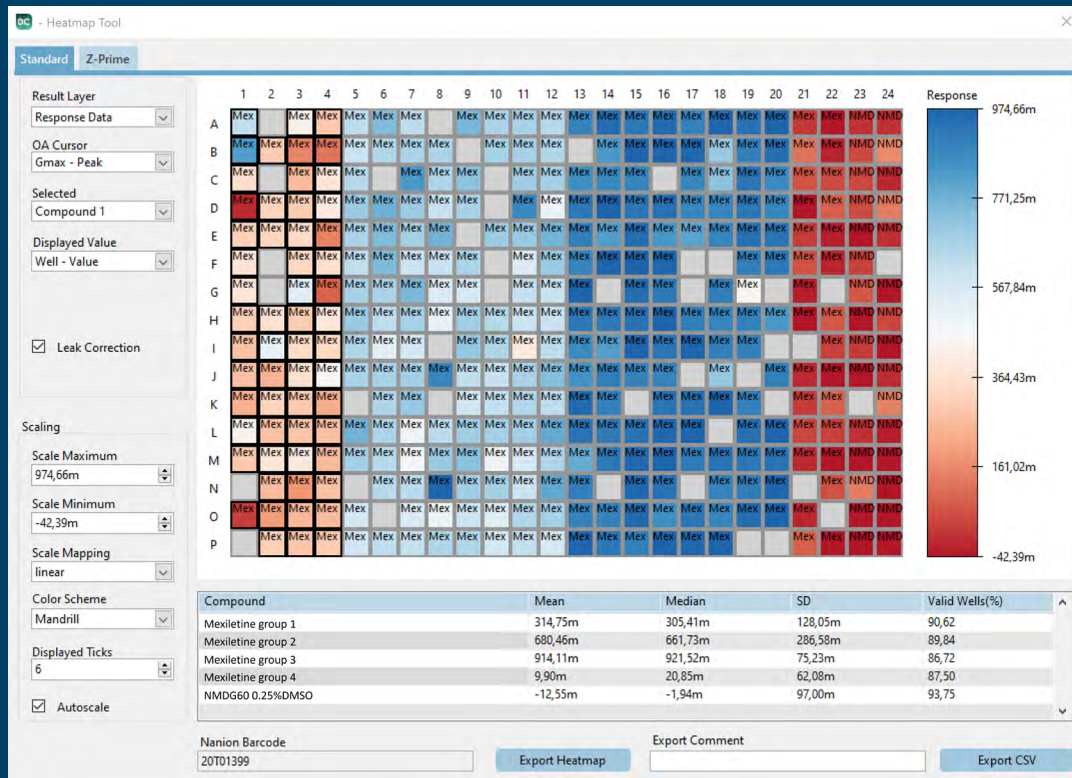


Automatically generate conductance-voltage (GV) curves for better clarity. Here the curves from reference and one concentration of mexiletine is shown.

SyncroPatch 384

IV analysis tool

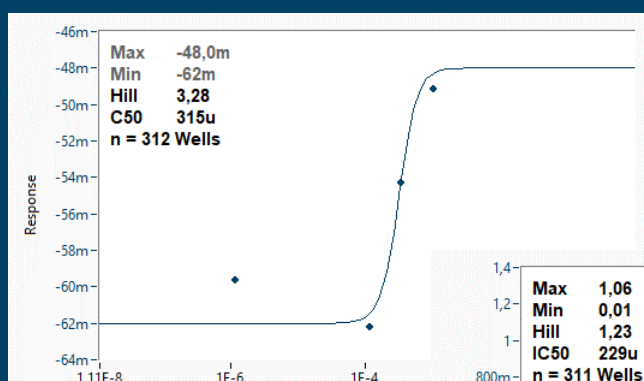
nanjion



The data derived from the IV curves (Gmax, Gmin, Vhalf and slope) can be viewed as a heatmap for a quick overview. Here the effect of mexiletine on Gmax is shown across the plate.

Full Flexibility

Use different parameters to generate CRC



The IC_{50} of compounds can be calculated using a variety of parameters e.g. Vhalf (left) or Gmax (below).

