



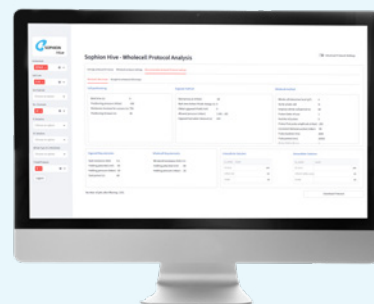
Sophion Hive assay development software

Simplify setup, accelerate your research.

Sophion Hive is a data-driven web-based platform built on decades of experimental data. It helps new and experienced users design and optimize whole-cell protocols to maximize experiment throughput. Transfer settings directly to QPatch or QPatch Compact for faster, reproducible, high-quality results.

With Sophion Hive, you simplify setup, optimize your whole-cell protocols, and configure your experiments for success.

- Accelerate your research with over 60 pre-configured protocol settings, and access a proprietary database of more than 23,000 ready-to-use whole-cell protocols for rapid setup and onboarding
- Harness over 25 years of Sophion experiment data to optimize whole-cell protocols for consistent, reproducible, high-quality results
- Smart filtering by experimental conditions, ion channel and cell type – even the most challenging cell types like hiPSC-cardiomyocytes, neurons, etc.
- Transfer optimized settings directly to QPatch and QPatch Compact to minimize platform setup time and reduce manual errors



Get instant recommendations for whole-cell protocol settings based on your selected filters.

GD46287-1

Get started today.

Discover more: sophion.com

 **SOPHION** | Discover more
BIOSCIENCE