

Genedata Ready-to-Run

Scale Up Your Data Analysis to Match Your Instrument

Genedata and Sophion Bioscience® have partnered to enable the direct analysis of raw data from Sophion instruments in Genedata Screener®, allowing a smooth and easy execution of automated patch clamp (APC) workflows.

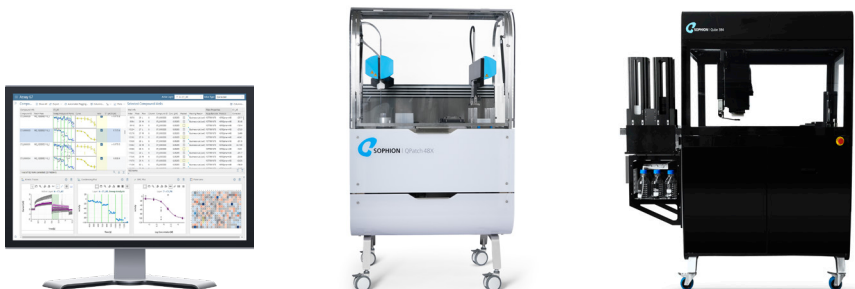
The integration provides automated and standardized data analysis and ensures that data can easily be imported and analyzed in Genedata Screener, using current and future Genedata Screener software versions.

The joint solution offers:

- ▶ Direct import of current and voltage traces data from Sophion instruments into Genedata Screener
- ▶ Automated, standardized, and scalable analysis of APC data in a screening context
- ▶ Easy comparison of ion channel and toxicity results with results from other assays in a data warehouse

The Ready-to-Run integration covers:

- ▶ Sophion Qube 384
- ▶ Sophion QPatch II



Activate your Genedata-Sophion Ready-to-Run integration:

Contact your Genedata Screener scientific account manager or your Sophion account manager.

If you have one solution and would like to add the other, contact:

info@sophion.com
screener@genedata.com

Genedata Ready-to-Run

Genedata Ready-to-Run program aims to create seamless connections by providing and maintaining instrument- and software-specific integrations for key workflows, saving our common customers time and effort when setting up or expanding their operations.

Go to www.genedata.com/ready-to-run for the latest information on instruments integrated with Genedata Screener.



Genedata transforms data into intelligence with innovative software solutions and domain-specific consulting services that automate complex, large-scale experimental processes and enable organizations to maximize the ROI in their R&D, spanning early discovery all the way to the clinic. Founded in 1997, Genedata is headquartered in Basel, Switzerland with additional offices in Boston, London, Munich, San Francisco, Singapore, and Tokyo.

www.genedata.com