

# Genedata Ready-to-Run

## Scale Up Your Data Analysis to Match Your Instrument

Genedata and Nanion® have partnered to offer a Ready-to-Run integration, allowing you to directly analyze your APC data in a screening context.

The joint solution offers:

- ▶ Direct import of current and voltage traces, directly from the binary file
- ▶ Complete data analysis from current traces to IC50s in one environment
- ▶ Integration of your Nanion instrument into your screening workflow
- ▶ Analysis and combination of multiple events

The integration ensures that data from Nanion instruments can easily be analyzed in Genedata Screener, using both current and future software versions.

The Ready-to-Run integration covers:

- ▶ SyncroPatch 384PE

Go to [www.genedata.com/ready-to-run](http://www.genedata.com/ready-to-run) for the latest information on instruments integrated with Genedata Screener.



## Activate your Genedata-Nanion Ready-to-Run integration:

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

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

[info@nanion.de](mailto:info@nanion.de)

[screener@genedata.com](mailto:screener@genedata.com)

## Genedata Ready-to-Run

Genedata Ready-to-Run provides out-of-the-box integrations for complex screening technologies. The integrations require little to no configuration, are instrument specific, come with defined functionality, and are maintained by Genedata as part of a software license agreement.



Genedata Screener® is part of the Genedata portfolio of advanced software solutions that serve the evolving needs of drug discovery, industrial biotechnology, and other life sciences.

Basel | Boston | Munich | San Francisco | Tokyo

[www.genedata.com](http://www.genedata.com) | [screener@genedata.com](mailto:screener@genedata.com)

© 2016 Genedata AG. All rights reserved. Genedata Screener is a registered trademark of Genedata AG. All other product and service names mentioned are the trademarks of their respective companies. 01S16

