

# PerkinElmer Signals Notebook



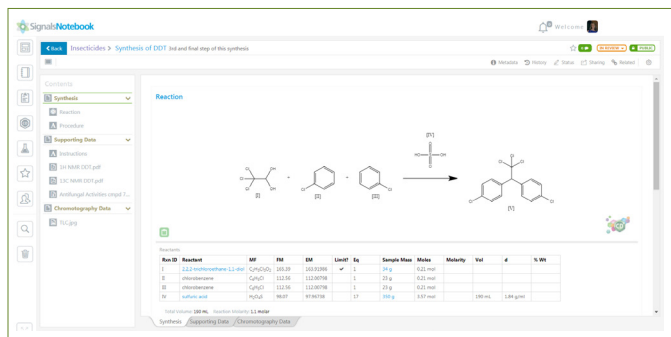
## Signals Notebook - Which Version Is Right For Me?

PerkinElmer Signals™ Notebook is the cloud-based electronic lab notebook (ELN) and scientific collaboration platform for today's scientists. With embedded ChemDraw to support synthetic chemistry workflows, drag and drop capture of a wide variety of electronic data files, lightning fast data entry and search, and a modern, highly responsive design, Signals Notebook is quickly becoming the preferred cloud-based ELN for discerning chemists and researchers alike.

**PerkinElmer Signals Notebook Standard** is designed to provide scientists with the tools they need to design, record and share their experiments. Signals Notebook provides organizations with peace of mind, delivering a business friendly framework to support the security, configurability and administrative capabilities demanded by today's leading research and industrial organizations. Organizations have control over sharing of experiments both throughout their enterprise and external to their organization, such as allowing controlled access to external partners. Administrators can define metadata to be captured along with experiments, and the access rules based upon them such as setting up project or auto-numbering rules for experiments. Organizations also retain ownership of the data when researchers move on.

**PerkinElmer Signals Notebook Private Cloud** provides all of the functionality offered in the Standard version, yet provides control over mandatory updates. Software updates are not executed automatically, enabling organizations to achieve greater control over when and how changes are made, as typically required for a validated environment. Signals Notebook Private Cloud seamlessly supports the validation of processes used in controlled workflows. Additional test environments can also be provided.

**PerkinElmer Signals Notebook Individual Edition** provides access to the same data capture and scientific workflow capabilities as PerkinElmer Signals Notebook Standard, but is designed to facilitate a more flexible sharing and collaborative model where individual researchers maintain ownership and control of their own data. Ideally suited for researchers who want to share with a wide variety of users inside and outside of their organizations (such as researchers at universities), Signals Notebook Individual Edition is only available to licensees of ChemOffice® Professional, our premiere chemical sketching and scientific productivity suite. Signals Notebook Individual Edition provides a lower cost entry into the modern scientific collaboration platform of PerkinElmer Signals Notebook Standard for users who do not require organizational control.



Designing Reactions with Signals Notebook Individual Edition

Top Above: Reaction interpretation in PerkinElmer Signals Notebook Individual Edition

Capability	Signals Notebook Private Cloud	Signals Notebook Standard (Preferred by most)	Signals Notebook Individual Edition
Cloud based application, no server required (SaaS solution)	✓	✓	✓
Embedded sketcher, with optional stoichiometric table for chemical reactions	✓	✓	✓
Comment on experiments	✓	✓	✓
Configure email and app notifications	✓	✓	✓
ChemDraw round-trip editing	✓	✓	✓
Access to ChemACX Explorer	✓	✓	✓
Define timeout period	✓	✓	✓
Public Tenant (Share content with any other Individual Edition user)			✓
Included with ChemOffice Professional			✓
Per user licensing model	✓	✓	
Private Tenant (Dedicated URL, logo)	✓	✓	
Administrators create and remove users, manage usergroups	✓	✓	
Transfer user data to another user	✓	✓	
Create system rules for sharing	✓	✓	
Create rules for notebook and experiment auto-numbering	✓	✓	
Create new metadata for experiments (projects, protocols etc.)	✓	✓	
Restrict access via IP range	✓	✓	
Configurable SAML authentication (shared desktop credentials)	✓	✓	
Alternate geographical location for data storage	✓	✓	
Configure sign, close and witness workflows	✓	✓	
Define user alias for auto-naming	✓	✓	
Create rules for trashing notebooks and experiments	✓	✓	
Define autotext (lists, text snippets)	✓	✓	
Private Environment (Dedicated architecture)	✓		
Optionally validated system	✓		
Automatic updates		✓	✓

## Common Questions

***I am a user on Signals Notebook Individual Edition. Can I move my data to my organization's Signals Notebook Standard?***

Yes, we can execute a bulk data migration from the public tenant to a private Signals Notebook Standard tenant. Please reach out to your PerkinElmer account representative for more information.

***I am a user on Signals Notebook Individual Edition. Can I share my data with someone else on the system?***

Yes, you can share your experiments and notebooks with anyone else, assuming that person is also a user of Signals Notebook Individual Edition.

***I am a user of Signals Notebook Standard. Can I share my data with someone else on the system?***

You can only share your experiments and notebooks with other users of your organization's system. Your organization can choose to set up CROs and other outside users with whom you can share, but you cannot share with users of another organization's Signals Notebook or Signals Notebook Individual Edition. Your organization may also choose to limit sharing to align with their internal business rules.

***Can I purchase a single license to Signals Notebook Individual Edition?***

No. Currently, Signals Notebook Individual is only provided via a license to ChemOffice Professional.

***I bought ChemOffice Professional. How do I get access to Signals Notebook Individual Edition?***

You should have received an email with your activation details for ChemOffice Professional. There is an activation link in that email for Signals Notebook Individual. If you no longer have the email please contact support at [informatics.support@perkinelmer.com](mailto:informatics.support@perkinelmer.com)

***I bought ChemOffice Professional and my company bought Signals Notebook Standard. Which Notebook should I use?***

There are cases where you may have eligibility to both an organizational Signals Notebook Standard and Signals Notebook Individual Edition. Access to each is defined by the URL you use to access them. You should use each appropriately for your work. Data entered into one is not available in the other.

***I had access to PerkinElmer Signals Notebook for ChemDraw. What has changed?***

Nothing except the name. Signals Notebook Individual Edition is identical to Signals Notebook for ChemDraw and still has all your data.

To register for a free trial of Signals Notebook Standard Edition for you and your colleagues, please visit <https://signalsnotebook.perkinelmer.cloud/trial/>

To download your free trial of Signals Notebook Individual Edition via ChemOffice Professional please visit <http://www.perkinelmer.com/product/chemoffice-professional-chemofficepro>

Contact us at 1-800-762-4000 to speak with a PerkinElmer Representative

For more information on PerkinElmer Signals Notebook Standard and PerkinElmer Signals Notebook Individual Edition, visit <http://www.perkinelmer.com/signalsnotebook>

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2017-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.