

# ChemOffice® 17.1 Suite of Products



PerkinElmer is collaborating with Chemical Abstracts Service to combine the power of two leading chemistry solutions, ChemDraw® and SciFinder®.

New Features	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
Hotkey Enhancements	Win/Mac	X	X	X
High-DPI Monitor Support	Win/Mac	X	X	X
Facilitated Copy/Pasting	Win/Mac	X	X	X
Stereochemistry Handling Improvements	Win/Mac	X	X	X
Support for HELM Notation	Win/Mac		X	X
CAS RN to Structure from ChemACX.com	Win/Mac		X	X
IUPAC name-based Atom Numbering	Win/Mac		X	X
PerkinElmer Signals™ Notebook Individual Edition	Win/Mac*			X
Mnova ChemDraw Edition	Win/Mac			X
ChemDraw Add-ins	Win/Mac			X

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
ChemDraw	Win/Mac	X	X	X
Multiple ChemDraw Items Folders	Win/Mac	X	X	X
Save and Read Graphic Files	Win/Mac	X	X	X
Save and Read Chemical Files	Win/Mac	X	X	X
Printing Options	Win/Mac	X	X	X
Chemical Templates	Win/Mac	X	X	X
Equipment Templates	Win/Mac	X	X	X
Analyze/Check Structures	Win/Mac	X	X	X
Insert OLE Object in ChemDraw	Win	X	X	X
In-place OLE Editing of ChemDraw Objects	Win	X	X	X
Show Stereochemistry	Win/Mac	X	X	X
Relative Stereochemistry (ISIS compatibility)	Win/Mac	X	X	X
Reaction Interpretation	Win/Mac	X	X	X

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
Reaction Mapping	Win/Mac	X	X	X
Calculate Properties	Win/Mac	X	X	X
Document Tagging	Win/Mac	X	X	X
Manual spectrum/structure assignments	Win/Mac	X	X	X
Chemical Polymer Tools	Win/Mac	X	X	X
Structure Clean Up	Win/Mac	X	X	X
Hotkeys	Win/Mac	X	X	X
Expand/Contract Labels	Win/Mac	X	X	X
Create/Use Nicknames	Win/Mac	X	X	X
Expand Generic Structure	Win/Mac	X	X	X
Multicenter Attachments	Win/Mac	X	X	X
TLC/GEP Tools	Win/Mac	X	X	X
Fragmentation Tools	Win/Mac	X	X	X
ChemDraw Active X Plugin	Win	X	X	X
Copy/Paste as SMILES	Win/Mac	X	X	X
Copy/Paste as SYBYL (SLN)	Win/Mac	X	X	X
Copy/Paste as InChI	Win/Mac	X	X	X
Copy/Paste as Molfile/Mol3000	Win/Mac	X	X	X
Copy/Paste as CDXML	Win/Mac	X	X	X
pKa LogP LogS	Win/Mac	X	X	X
tPSA	Win/Mac	X	X	X
Advanced Retrosynthesis Tool	Win/Mac		X	X
Auto-numbering of multiple structures	Win/Mac		X	X
Search SciFinder	Win/Mac		X	X
Name = Structure/Structure = Name	Win/Mac		X	X
cLogP	Win/Mac		X	X
Biopolymer Toolbar	Win/Mac		X	X
BioDraw	Win/Mac		X	X
Reaction Stoichiometry Grid	Win/Mac		X	X
Calculate <sup>1</sup> H <sup>13</sup> C NMR Spectra	Win/Mac		X	X
Query Features	Win/Mac		X	X
Query Tools	Win/Mac		X	X
Advanced Stereochemistry	Win/Mac		X	X
Paste as HELM	Win/Mac		X	X
Create Sequence	Win/Mac		X	X
Create New Monomer	Win/Mac		X	X
Copy as HELM	Win/Mac		X	X
ChemDraw Cloud	Win/Mac*		X	X
ChemFinder (Std in CD Prof; Ultra in CO)	Win		X	X
ChemDraw for Excel	Win		X	X
Name=Struct for ChemDraw for Excel	Win		X	X
ChemScript + Python	Win		X	X

\*Access to ChemDraw Cloud and Signals Notebook is provided for one year and can be renewed if current with maintenance.

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
CombiChem for Excel	Win		X	X
3D Search	Win		X	X
Chem3D (Pro in CDPro; Ultra in CO)	Win		X	X
Chem3D Hotlink	Win		X	X
Chem3D Active X Plugin	Win		X	X
GAMESS	Win			X
Interface to MOPAC 2016	Win			X
Interface to Gaussian	Win			X
Interface to Conflex	Win			X
Interface to Autodock	Win			X
ChemFinder / Oracle	Win			X
ChemFinder for Office	Win			X
BioViz in ChemFinder Ultra	Win			X
Compound Profiles in ChemFinder Ultra	Win			X
Clustering in ChemFinder Ultra	Win			X
Combine ChemFinder Query Hit Lists	Win			X
ChemFinder exports to MS Word / Excel	Win			X

Visit [www.cambridgesoft.com/software/overview.aspx](http://www.cambridgesoft.com/software/overview.aspx) for more information.

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 © 2015-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.