

# ChemOffice+ Cloud 20.0 Suite of Products



New features	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice+ Cloud
Atom Hotkey Enhancements	Win/Mac	X	X	X
New "Enter" Molecule Hotkey	Win/Mac	X	X	X
3D Clean-up	Win/Mac		X	X
Enhanced 3D Display	Win/Mac		X	X
Atom/Bonds Color Highlights	Win/Mac		X	X
Google Patents Add-in	Win/Mac			X
ChemOffice+ Individual (requires Signals Notebook Login)	Win/Mac			X

ChemOffice+ Features	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice+ Cloud
Browse & Drill-down Into ChemDraw Files	Win/Mac			X
Browse .mol/.sdf Files	Win/Mac			X
Browse & Drill-down ChemDraw Files Inside MS Word	Win			X
Browse & Drill-down ChemDraw Files Inside MS Powerpoint	Win			X
Copy to Structure(s) to Clipboard	Win/Mac			X
Create Collection	Win/Mac			X
Adding Properties to Collection	Win/Mac			X
Edit Properties	Win/Mac			X
Edit Multiple Chemical Structures	Win/Mac			X
Combine Collections	Win/Mac			X
Export to SD file (v2000, v3000)	Win/Mac			X
Create Powerpoint Report Slide	Win/Mac			X

Recent Additions	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice+ Cloud
Aromatic cycle display toggle	Win/Mac	X	X	X
Hotkey Enhancements	Win/Mac	X	X	X
Smart Copy/Paste	Win/Mac	X	X	X
Stereochemistry Handling Improvements	Win/Mac	X	X	X
Rotation Shortcut	Win/Mac	X	X	X
4K Display Support	Win/Mac	X	X	X
Ring Fill Coloring	Win/Mac		X	X
HELM 2.0 Support	Win/Mac		X	X
Search into Reaxys	Win/Mac		X	X
Search into SciFinder-n	Win/Mac		X	X
IUPAC-based Atom Numbering	Win/Mac		X	X
PubChem GHS Safety Add-in	Win/Mac			X
Add-ins Dynamic Download	Win/Mac			X
ChemACX Explorer	Win/Mac			X
Shared HELM Libraries	Win/Mac			X

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice+ Cloud
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
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
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
ChemDraw JS (with Site Subscription Only)	Win/Mac			X

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice+ Cloud
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
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
Interface to Conflex	Win			X
Interface to Autodock	Win			X
ChemFinder / Oracle	Win			X
Interface to GAMESS 18	Win			X
Interface to Gaussian 16W	Win			X
Interface to MOPAC 2016	Win			X
ChemFinder for Office	Win			X
Explorer Window View in ChemFinder Ultra	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

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

Visit [www.perkinelmer.com/category/chemdraw](http://www.perkinelmer.com/category/chemdraw) 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 © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

125790 (77247) PKI