

ChemOffice® 18.0 Suite of Products



New Features	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
Search into Reaxys	Win/Mac		X	X
ChemACX Explorer	Win/Mac			X
Shared HELM Libraries	Win/Mac			X

Recent Additions	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
ChemDraw JS (with Site Subscription Only)	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

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
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
Reaction Stoichiometry Grid	Win/Mac		X	X
Calculate ¹ H ¹³ 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

Includes	Platform	ChemDraw Prime	ChemDraw Professional	ChemOffice Professional
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 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

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

Visit 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



For a complete listing of our global offices, visit 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.

012204Q_01

PKI