

Q Do end-users need to install any software?

A No. All the end-user needs to run is a compliant web browser with network access to the ChemDraw JS server.

Q What is “extended copy-paste”?

A Modern browsers place restrictions on what data can be read from the clipboard. Most chemical drawing programs place data on the clipboard in a format that cannot be read by any web browser. Extended copy-paste enables ChemDraw JS to read chemical structures from the clipboard, but require installing an additional browser applet or extension.

Q Can I copy/paste from ChemDraw JS into other applications?

A Yes. With an additional, browser-specific clipboard extension, it is possible to copy (Ctrl+C) ChemDraw JS objects and paste them into ChemDraw or a MS Office document. After modifying them, you can paste them from ChemDraw or MS Word/Powerpoint back into ChemDraw JS without loss of information. (Windows only, does not work with MS Office Online documents).

Q What features require access to ChemDraw JS services?

A End user tools/actions:

- Name-to-Structure
- Structure-to-Name

JavaScript API calls:

- findReactions
- getProperties
- getSVG
- getMolecularWeight
- nameToStructure
- structureToName

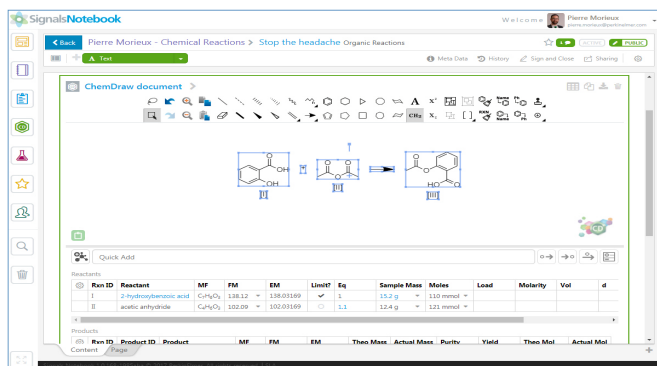


Figure 3. ChemDraw JS Sketcher in PerkinElmer Signals™ Notebook.