



## CHEMDRAW JS SHARING

ChemDraw JS generates and reads the same file formats used by ChemDraw Desktop so that files can be easily shared. Additional formats are also available for sharing with other applications:

- CDXML
- molfile and rxnfile
- SMILES
- InChI and InChIKey
- Chemical drawings copied from ChemDraw Desktop, or from other third-party editors, can be directly pasted into ChemDraw JS for immediate exchange of structures and reactions between desktop and web browsers.
- Chemical drawings embedded in MS Office applications (Windows) can be copied/pasted into ChemDraw JS without loss of information (requires browser extension).

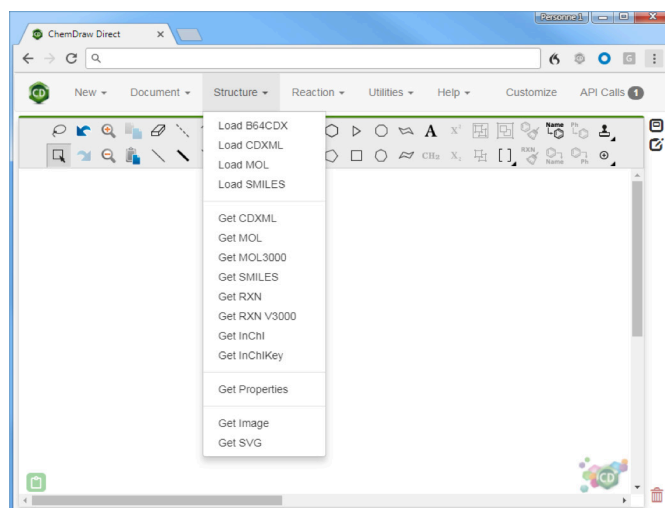
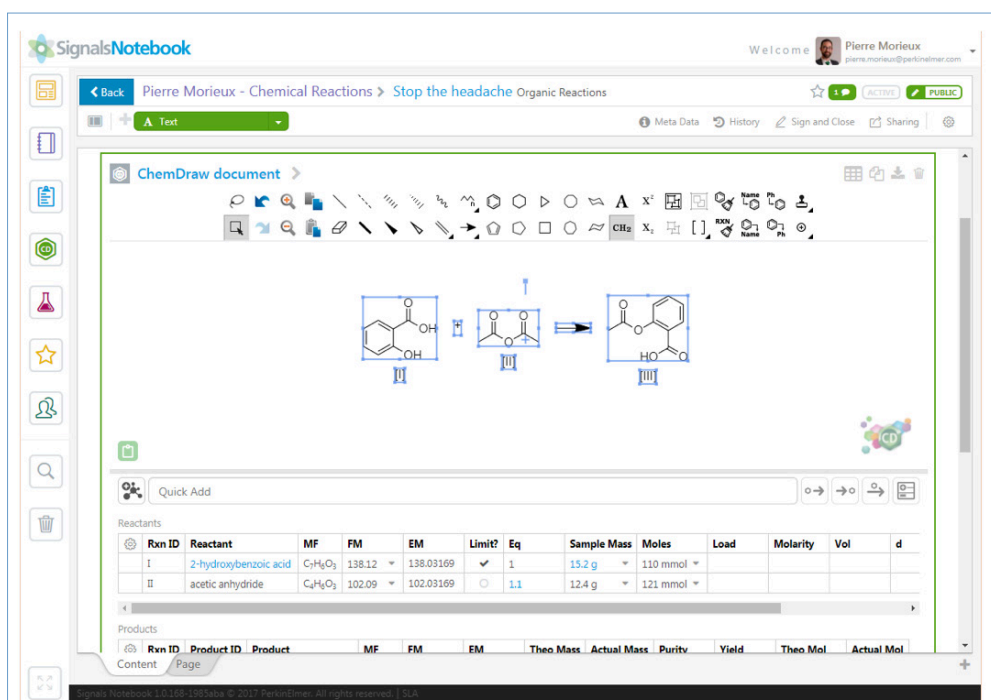


Figure 2. Available API calls for ChemDraw JS.

## CHEMDRAW JS DEPLOYMENT

ChemDraw JS is supplied with a fully documented User's Guide, Installation Guide and Developers' Guide. The latter includes instructions on how to load and to embed ChemDraw JS into an application, and details of how to make API calls to generate the various file formats that are available.

A screenshot of the PerkinElmer Signals Notebook interface. The main window displays a chemical reaction scheme involving 2-hydroxybenzoic acid and acetic anhydride. Below the reaction is a table of reactants with columns for Rxn ID, Reactant, MF, FM, EM, Limit?, Eq, Sample Mass, Moles, Load, Molarity, Vol, and d. The table contains two rows of data for reactants I and II.

Rxn ID	Reactant	MF	FM	EM	Limit?	Eq	Sample Mass	Moles	Load	Molarity	Vol	d
I	2-hydroxybenzoic acid	C <sub>7</sub> H <sub>6</sub> O <sub>3</sub>	138.12	138.03169	✓	1	15.2 g	110 mmol				
II	acetic anhydride	C <sub>4</sub> H <sub>6</sub> O <sub>3</sub>	102.09	102.03169	○	1.1	12.4 g	121 mmol				

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

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.

012318G\_01

PKI

DRAFT