

ChemDraw Add-in Architecture

An Add-in architecture has been developed that allows you to develop custom add-ins and extend the capabilities of ChemDraw, by building integrations with other research applications using the latest web technologies. The Add-in platform gives access to a ChemDraw JS API that can be called by external applications to access the various functions of ChemDraw. Add-ins can now be dynamically downloaded from a local folder or a remote URL for a more controlled access of additional capabilities.

Available with ChemOffice Professional

ChemDraw JS

The JavaScript version of ChemDraw allows your custom applications to become "ChemDraw Smart." Built on JavaScript and HTML5, ChemDraw JS can be plugged into web browsers and internal applications that need to be chemicalized. Inside its lightweight packages, ChemDraw JS carries the core functionality of the ChemDraw family including: your favorite Hotkeys/Shortcuts, Name to Structure, Structure to Name, Templates and Structure Query Tools.

Available in ChemOffice Professional Site License, Internal usage only.

For more information please visit www.perkinelmer.com/product/chemdraw-and-chemoffice-chemdraw

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 ©2018-2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014423B_01

PKI

DRAFT