

HUMAN HEALTH

ENVIRONMENTAL HEALTH

ChemOffice Version 19.0

Hardware & Software Guide

Last Updated: September 20, 2019



Table of Contents

1	Introduction.....	3
2	Hardware Requirements	3
3	Software Requirements.....	3
4	Technical Support.....	4

1 Introduction

This document is intended for general use with PerkinElmer Desktop 19.0 customers.

2 Hardware Requirements

Hardware Component	Minimum Specification	Recommended Specification
Memory (RAM)	For 32-bit OS 1 GB For 64-bit OS 2 GB	For 32-bit OS 1 GB For 64-bit OS 2 GB
Processor	1.6 GHz or higher	3.0 GHz or higher dual core processor
Free Space Required on Hard Drive	300 MB - 1 GB (1 GB is for installing the full ChemOffice Professional). This is for installation only and does not include any creation/expansion of any databases.	10 GB
Screen resolution	1024 x 768 or higher	1024 x 768 or higher

3 Software Requirements

Software	Windows	Macintosh (ChemDraw Only)
Operating System	<ul style="list-style-type: none"> Windows 7 Professional and Ultimate (32-bit) Windows 7 Professional and Ultimate (64-bit) Windows 8.1 (64-bit) Windows 10 (64-bit)* 	<ul style="list-style-type: none"> Mac OS X 10.15*, 10.14 and 10.13 <p>* Qualified against the beta candidate.</p>
Microsoft Office (Standard, Professional, or Enterprise)	<ul style="list-style-type: none"> Microsoft Office 2016 (32-bit and 64-bit)* Microsoft Office 2019 (32-bit and 64-bit)* Office 365* <p>*The latest update used for testing is 1908 (11929.20300).</p> <p>Note: We test with the latest versions of MS Office available at the time, but that is no guarantee of full compatibility with future updates.</p>	<ul style="list-style-type: none"> Microsoft Office 2016* Microsoft Office 2019* Office 365* <p>* The latest update used for testing is 16.29.1 (19091700).</p>
.Net Framework	<ul style="list-style-type: none"> .Net Framework 4.x 	<ul style="list-style-type: none"> N/A
Browser ActiveX controls:	<ul style="list-style-type: none"> Microsoft Internet Explorer 10.x (32-bit and 64-bit) Microsoft Internet Explorer 11.x (32-bit and 64-bit) 	<ul style="list-style-type: none"> N/A <p>Note: The ChemDraw & Chem3D Plugins are no longer part of any ChemOffice/ChemDraw package. The ChemDraw & Chem3D ActiveX controls are still included with windows packages.</p>

4 Technical Support

To access FAQs, KB articles, support forums and more, please visit:

<https://informatics.perkinelmer.com/Support/>

For support contact options, please visit:

<https://informatics.perkinelmer.com/Support/Contact/>

PerkinElmer Inc. | For the Better
HUMAN HEALTH | ENVIRONMENTAL HEALTH
940 Winter Street | Waltham, MA 02451