

HUMAN HEALTH

ENVIRONMENTAL HEALTH



Signals Notebook

Hardware and Software Guide

Last Updated: July 25, 2018

1 Introduction

This document describes the supported and qualified software components to seamlessly use Signals Notebook application. It also describes the minimum and recommended hardware components required to access Signals Notebook application.

2 Hardware Requirements

Any mid-range client machine with minimum 8GB of RAM should be sufficient.

Note: Signals Notebook is not qualified currently on mobile platforms (tablets, phones).

3 Software Requirements

We specifically qualify Internet Explorer 11 and support most modern browsers such as Chrome, Firefox and Safari 11.

For the ChemDraw Web Clipboard extension, Java Runtime Environment 1.7+ may be required according to end user's Operating System and browser.

Editing in ChemDraw Desktop requires ChemDraw 17.1 or above for PC and ChemDraw 17.1.1 for Mac. ChemDraw Edit in Desktop 17.1 is supported in Sierra and High Sierra operating system on Mac and Windows 7 and 10.

3.1 Supported Environments

Following environments are supported:

Browser	Operating System	JAVA
Google Chrome	Windows or macOS	Not required
Mozilla Firefox 50 or later	Windows or macOS	Not required
Apple Safari 10 or later	macOS 10.12 or later	Not required
Microsoft Internet Explorer 11	Windows	Required, Applet enabled
Mozilla Firefox 49 or earlier	Windows or macOS	Required, Applet enabled
Apple Safari 9 or earlier	macOS 10.11 or earlier	Required, Applet enabled

3.2 MS Office

Signals Notebook is qualified to work with MSOffice 2013 Service Pack 1 and MSOffice 2016.

To work with Online MSOffice option, you need to have a valid Office 365 account with Office Online subscription.

The following table covers the file size limits for viewing and editing documents in Office Online:

Application	Size Limit in View Mode	Size Limit in Edit Mode
Excel Online	5 MB	5 MB
PowerPoint Online	None	300 MB
Word Online	None	None