

HUMAN HEALTH

ENVIRONMENTAL HEALTH



Signals Notebook

Hardware and Software Guide

Last Updated: Sep 30, 2020

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 support most modern browsers such as Chrome, Firefox, Safari 11, and Chromium-based Microsoft Edge.
- 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.
- ChemDraw Edit in Desktop 18.2 is supported in MacOS High Sierra and Mojave and Windows 7 (32-bit and 64-bit), 8.1 (64-bit) and 10 (64-bit).
- Editing in Tibco Spotfire requires Tibco Spotfire Analyst LTS 10.3 or 10.10. The add-in **SignalsNotebookSpotfireOpener-v1.x.spk** must also be installed on the Spotfire server and distributed to the user. Editing in Spotfire is supported in Windows 10 system only. The file size limit that is supported for opening and saving dxp files in Spotfire is 50 MB. Users can modify larger files via Windows Registry Editor, for more details please refer to: <https://support.microsoft.com/en-sg/help/2668751/you-cannot-download-more-than-50-mb-or-upload-large-files-when-the-upl>.

3.1 Supported Environments

Following environments are supported:

Browser	Operating System
Google Chrome	Windows or macOS
Mozilla Firefox 50 or later	Windows or macOS
Apple Safari 10 or later	macOS 10.12 or later
Chromium-based Microsoft Edge	Windows or macOS

3.2 MS Office

Signals Notebook is qualified to work with MSOffice 2016 and MSOffice 2019.

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	100 MB

Note: The overall Signals Notebook limit is set to 2 GB.

3.3 SAML

Signals Notebook supports SAML 2.0 IDPs.

PKI requires from the customer	PKI provides to the customer
<ul style="list-style-type: none">• Login URL for the IDP• SAML Key/IDP certificate	<ul style="list-style-type: none">• Asset URL• Application ID