

HUMAN HEALTH

ENVIRONMENTAL HEALTH

Lead Discovery™ 8.1

System Requirements

Powered by TIBCO Spotfire®

Last Updated: Oct 26, 2018

PerkinElmer Lead Discovery™ 8.1 System Requirements

Software Requirements	
PerkinElmer Lead Discovery™ 8.1 for TIBCO Spotfire 7.7, 7.8, 7.9, 7.10 and 7.11	
Category	Requirement
TIBCO Spotfire	<p>TIBCO Spotfire Client:</p> <p>TIBCO Spotfire Analyst 7.7.x TIBCO Spotfire Business Author/Consumer 7.7.x TIBCO Spotfire Analyst 7.8.x TIBCO Spotfire Business Author/Consumer 7.8.x TIBCO Spotfire Analyst 7.9.x TIBCO Spotfire Business Author/Consumer 7.9.x TIBCO Spotfire Analyst 7.10.x TIBCO Spotfire Business Author/Consumer 7.10.x TIBCO Spotfire Analyst 7.11.x LTS TIBCO Spotfire Business Author/Consumer 7.11.x LTS</p> <p>TIBCO Spotfire Server:</p> <p>TIBCO Spotfire Server 7.7.x TIBCO Spotfire Server 7.8.x TIBCO Spotfire Server 7.9.x TIBCO Spotfire Server 7.10.x TIBCO Spotfire Server 7.11.x LTS</p>
Data Sources	<p>PerkinElmer CSCartridge 13.1 and 15.1 PerkinElmer ChemSearch Module for Attivio 1.0 Accelrys Accord Chemistry Cartridge 7.1 ChemAxon JChem Cartridge 5.2 and 6.2 IDBS ChemXtra 7.2 Accelrys Direct 5.1, 7.0, 8.0 and 9.0 GGA Bingo 1.6</p>

	<p>NOTE: Lead Discovery is designed to work with any Oracle-based cartridge, but this release of Lead Discovery is only tested and qualified with some listed above.</p>
Data Source Database	Oracle 10g or higher
Renderer	<p>ChemDraw ActiveX Enterprise 17.1 and earlier versions, 32 bit</p> <p>NOTE: ChemDraw ActiveX 12 has limitations for rendering and other functionalities in Lead Discovery, thus a higher version of ChemDraw ActiveX is recommended to be used.</p> <p>Accelrys Accord Chemistry Control 6.x and 7.0 ChemAxon Marvin View 17.x, 18.x, both 32 bit and 64 bit IDBS ChemIQ 7.2 BIOVIA Draw 2017, 2018, both 32 bit and 64 bit</p> <p>NOTE: Using Lead Discovery with the TIBCO Spotfire Web Player requires the use of a renderer that supports the Operating System the Web Player runs on (note that you also need to use the same renderer that is used in TIBCO Spotfire Analyst to generate the analyses that are then to be opened in the Web Player).</p> <p>Make sure to check the system requirements for the renderer before installing. For renderer system requirements. Please refer to the respective renderer provider.</p>
Editor	<p>For TIBCO Spotfire Analyst:</p> <p>ChemDraw ActiveX Enterprise 17.1 and earlier versions, 32 bit ChemAxon Marvin Sketch 17.x, 18.x, both 32 bit and 64 bit PerkinElmer ChemDraw JS 17.1 BIOVIA Draw 2017, 2018, both 32 bit and 64 bit</p> <p>For TIBCO Spotfire Web Client:</p> <p>PerkinElmer ChemDraw JS 17.1</p> <p>Note: By default, the ChemDraw JS in AWS hosted by PerkinElmer is used.</p> <p>ChemAxon Marvin JS 17.24 LTS</p>

Chemical Structure Formats	Lead Discovery supports embedded SMILES, CDX (CDXML), CHIME and MOLFILES strings, depending on the support of the renderer/editor.
Microsoft .NET Framework	Microsoft .NET Framework 4.5 or higher Note: Spotfire also has minimum .NET framework requirement. Refer to the system requirements of Spotfire Client for more information.
Browser	Please refer to the Spotfire Web Client System Requirements as reference for Lead Discovery in Spotfire Web Client. MS Internet Explorer 9 or higher is expected to support Trellis Card Visualization in Spotfire Analyst.

Hardware Requirements

As an extension of Spotfire, Lead Discovery uses the same hardware requirements as Spotfire.

Please refer to the following website for the Spotfire hardware requirements:

<http://support.spotfire.com/sr.asp>