



# **Spectral Database**

## Installation Instructions





# Installation Instructions

Review the system recommendations at [www.knowitall.com/minimum\\_system](http://www.knowitall.com/minimum_system).

1. Plug the Installation USB into your computer. From the USB, open the **KnowItAll Installation** folder, double-click **KnowItAllInstall.exe** and follow the instructions. If you are unable to install from the USB, download the software from [www.knowitall.com/download](http://www.knowitall.com/download).
2. Start the KnowItAll ID Expert software and follow the instructions to activate your license. The **Product Registration Code** can be found on the back of the KnowItAll packaging.
3. From the Installation USB, open the **Individually Licensed Databases** folder and copy the files for the spectral database(s) you purchased to your computer.

## What's Included in Your Spectral Database Purchase

Your spectral database purchase also includes	
<b>KnowItAll® ID Expert™</b>	Spectral identification software with single-component, multi-component, and functional group analysis. KnowItAll ID Expert is easy to use: simply open a spectrum and view the results. See demo at <a href="http://www.knowitall.com/idexpertvideo">www.knowitall.com/idexpertvideo</a> .
<b>ATR-IR Solvents - Bio-Rad Sadtler Database</b>	629 high-quality ATR-IR spectra of organic solvents
<b>KnowItAll IR Spectral Library (Trial)</b>	30-day access to over 235,000 IR spectra
<b>KnowItAll Raman Spectral Library (Trial)</b>	30-day access to over 15,000 Raman spectra

**English**   日本語   **Deutsch**   **Français**   中文

PDF versions of this document are available in English, Japanese, German, French, & Chinese on the KnowItAll Software & Database Installation USB.

このドキュメントは英語、日本語、ドイツ語、フランス語、中国語それぞれPDF形式でインストールUSBに格納されています。

PDF-Versionen dieses Dokuments finden sie auf Englisch, Japanisch, Deutsch, Französisch und Chinesisch auf dem Installations-USB-Stick für die KnowItAll-Software und Datenbanken.

Des versions PDF de ce document sont disponibles en anglais, japonais, allemand, français & chinois sur la clé USB d'installation du logiciel et des bases de données KnowItAll.

本文档有英、德、法、日语及中文PDF版。

**BIO-RAD**

**Bio-Rad**  
**Laboratories, Inc.**

*Informatics Division*

**Website:** [www.knowitall.com](http://www.knowitall.com)   **Contact Us:** [www.knowitall.com/contactus](http://www.knowitall.com/contactus)



210407-REV20160825