



Spectral Database

Installation Instructions





Installation Instructions

Review the system recommendations at 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.
2. Start the KnowItAll ID Expert software and follow the instructions to activate your license. The **License Verification Code** can be found on the back of the KnowItAll packaging or email if downloaded.
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	
KnowItAll® ID Expert™	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 www.knowitall.com/idexpertvideo .
ATR-IR Solvents - Bio-Rad Sadtler Database	629 high-quality ATR-IR spectra of organic solvents
KnowItAll IR Spectral Library (Trial)	30-day access to over 250,000 IR spectra
KnowItAll Raman Spectral Library (Trial)	30-day access to over 22,000 Raman spectra

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

PDF versions of this document are available in English, Japanese, German, French, & Chinese at www.knowitall.com/training

本マニュアルのPDFファイル（英語、日本語、ドイツ語、フランス語、中国語）は www.knowitall.com/training からダウンロードできます。

PDF-Versionen dieses Dokuments finden sie auf Englisch, Japanisch, Deutsch, Französisch und Chinesisch auf www.knowitall.com/training

Des versions PDF de ce document sont disponibles en anglais, japonais, allemand, français & chinois sur page www.knowitall.com/training

本文档有英、德、法、日语及中文 PDF 版，请看 www.knowitall.com/training

BIO-RAD

Bio-Rad
Laboratories, Inc.

Informatics Division

Website: www.knowitall.com **Contact Us:** www.knowitall.com/contactus



210407-REV20180202