



Databases Available with KnowItAll® U

Below is a listing of all of the databases that are available in Bio-Rad's KnowItAll U Spectroscopy Library (subscription for colleges / universities) and their codes.* For more information, visit www.knowitall.com

Database	Code
IR - Pesticides & Agricultural Chemicals - Bio-Rad Sadtler	AX
IR - Commonly Abused Drugs (Acid) - Bio-Rad Sadtler	AAX
IR - Commonly Abused Drugs (Base) - Bio-Rad Sadtler	ABX
MS - Bio-Rad Sadtler AAFS Toxicology Section Mass Spectra of Drugs	AF
IR - Adhesives & Sealants (Subset) - Bio-Rad Sadtler	ALX
IR - Automobile Paint Chips	APX
IR - Adhesives & Sealants - Bio-Rad Sadtler	ASX
IR - Polyols - Bio-Rad Sadtler	BX
IR - Polymers & Monomers (Basic) 2 - Bio-Rad Sadtler	BMX
IR - Polymers & Monomers (Basic) 3 - Bio-Rad Sadtler	BM2X
IR - Polymers & Monomers (Basic) 1 - Bio-Rad Sadtler	BPX
ATR-IR - Polymers & Monomers (Basic) 1 - Bio-Rad Sadtler	BWX
ATR-IR - Polymers & Monomers (Basic) 2 - Bio-Rad Sadtler	BW2X
ATR-IR - Polymers & Monomers (Basic) 3 - Bio-Rad Sadtler	BW3X
ATR-IR - Polymers & Monomers (Basic) 4 - Bio-Rad Sadtler	BW4X
IR - Surfactants (Comprehensive) - Bio-Rad Sadtler	CAX
IR - Surfactants (Basic) - Bio-Rad Sadtler	CBX
IR - Coating Chemicals - Bio-Rad Sadtler	CCX
IR - Surfactants (Subset) 1 - Bio-Rad Sadtler	CLX
IR - Surfactants (Subset) 2 - Bio-Rad Sadtler	CMX
IR - Epoxy Resins, Curing Agents & Additives - Bio-Rad Sadtler	CRX
IR - Polymers & Monomers (Comprehensive) - Bio-Rad Sadtler	DAX
IR - Polymers, Controlled Pyrolyzates - Bio-Rad Sadtler	DDX
Multi-Technique Demo Database - Bio-Rad Sadtler	DEMO
IR - Controlled and Prescription Drugs 1 - Bio-Rad Sadtler	DIX
IR - Georgia State Crime Lab	DLX
IR - Polymer Additives - Bio-Rad Sadtler	DPX
ATR-IR - Controlled & Prescription Drugs 1 - Bio-Rad Sadtler	DWX
ATR-IR - Controlled & Prescription Drugs 2 - Bio-Rad Sadtler	DW2X
ATR-IR - Controlled & Prescription Drugs 3 - Bio-Rad Sadtler	DW3X

IR - Plasticizers - Bio-Rad Sadtler	EX
IR - EPA Vapor Phase - Bio-Rad Sadtler	ELX
ATR-IR - Plasticizers - Bio-Rad Sadtler	EWX
IR - Food Additives - Bio-Rad Sadtler	FX
IR - Food Additives (Revised) - Bio-Rad Sadtler	FAX
IR - Flavors, Fragrances & Oils - Bio-Rad Sadtler	FFX
Raman - Flavors & Fragrances - Bio-Rad Sadtler	FFRX
ATR-IR - Flavors & Fragrances - Bio-Rad Sadtler	FFWX
IR - Fibers by Microscope - Bio-Rad Sadtler	FMX
IR - Flavors & Fragrances - Bio-Rad Sadtler	FIRX
IR - Flame Retardants - Bio-Rad Sadtler	FRX
IR - Fats, Waxes & Derivatives - Bio-Rad Sadtler	GX
ATR-IR - Nutraceuticals - Bio-Rad Sadtler	GNX
ATR-IR - Nutraceuticals 2 - Bio-Rad Sadtler	GN2X
IR - Gases & Vapors - Bio-Rad Sadtler	GSX
IR - Lubricants 1 - Bio-Rad Sadtler	HX
IR - Lubricants 2 - Bio-Rad Sadtler	H2X
IR - Surfactants, Hummel - Wiley	HAX
¹ H NMR - Bio-Rad Sadtler	HB
IR - Polymers, Hummel Industrial Polymers - Wiley	HCX
IR - Polymers, Hummel Industrial Monomers - Wiley	HEX
IR - Polymer Additives, Hummel Industrial - Wiley	HFX
IR - Polymers, Hummel Defined Basic - Wiley	HDX
IR - Industrial Chemicals, Pure Organic Compounds - Wiley	HIX
IR - Industrial Chemicals, Basic Organic Compounds - Wiley	HLX
IR - Polymers, Hummel Defined - Wiley	HMX
IR - Polymers, Hummel Industrial - Wiley	HPX
IR - Polymer Processing Chemicals - Bio-Rad Sadtler Scholl	HSX
IR - Polymers, Hummel - Bio-Rad Sadtler	HUX
IR - Rubber Chemicals - Bio-Rad Sadtler	JX
IR - Coating Chemicals (Revised) - Bio-Rad Sadtler	JLX
IR - Fibers & Textile Chemicals - Bio-Rad Sadtler	KLX
IR - Solvents - Bio-Rad Sadtler	LX
IR - Lubricant Additives 1 - Bio-Rad Sadtler	LA1X
IR - Lubricant Additives 2 - Bio-Rad Sadtler	LA2X
IR - Solvents (Basic) - Bio-Rad Sadtler	LLX
IR - Solvents (Vapor Phase) - Bio-Rad Sadtler	LUX
IR - Intermediates - Bio-Rad Sadtler	MX
IR - Merck - Bio-Rad Sadtler	MBX
MS - Wiley Drugs	MD

NMR - Metabolites - Bio-Rad Sadtler	MET
MS - Wiley Geochemicals, Petrochemicals & Biomarkers	MG
IR - Intermediates (Basic) - Bio-Rad Sadtler	MLX
IR - Minerals & Clays - Bio-Rad Sadtler	MNX
MS - Wiley Industrial Chemicals	MT
MS - Wiley Androstanes, Estrogens & other Steroids	MU
MS - Wiley Volatile Compounds in Food	MV
IR - Organometallics - Bio-Rad Sadtler	NX
ATR-IR - NIOSH Pocket Guide to Chemical Hazards Compounds - Bio-Rad Sadtler	NAX
¹ H NMR - Wiley HNMR Spectra of Organic Compounds	NB
¹³ C NMR - Bio-Rad Sadtler	NC
IR - Electric Power Plant Materials - Bio-Rad Sadtler	NEX
¹³ C NMR - Flavors & Fragrances - Wiley	NF
¹ H NMR - NIOSH Pocket Guide to Chemical Hazards Compounds - Bio-Rad Sadtler	NH
IR - NIOSH Pocket Guide to Chemical Hazards Compounds - Bio-Rad Sadtler	NIX
¹³ C NMR - Polymers & Monomers - Bio-Rad Sadtler	NM
¹³ C NMR - NIOSH Pocket Guide to Chemical Hazards Compounds - Bio-Rad Sadtler	NN
¹³ C NMR - Organic Compounds -Wiley	NO
¹³ C NMR - Natural Products - Wiley	NP
¹ H NMR - Chemical Shifts - Bio-Rad Sadtler	NQ
IR - NIOSH Pocket Guide to Chemical Hazards Compounds - (Vapor Phase) - Bio-Rad Sadtler	NVX
ATR-IR - Organometallics 1 - Bio-Rad Sadtler	NWX
ATR-IR - Organometallics 2 - Bio-Rad Sadtler	NW2X
¹ H NMR - Organic Compounds (Comprehensive) - Wiley	NY2X
IR - Polymers & Monomers (Subset) 1 - Bio-Rad Sadtler	OLX
IR - Polymers & Monomers (Subset) 2 - Bio-Rad Sadtler	OMX
IR - Organosilicons	OSX
IR - Petroleum Chemicals - Bio-Rad Sadtler	PX
IR - Prepared & Prescription Drugs (Acid) - Bio-Rad Sadtler	PAX
IR - Prepared & Prescription Drugs (Base) - Bio-Rad Sadtler	PBX
IR - Polymer Additives (Revised) - Bio-Rad Sadtler	PLX
IR - Standards (Subset) 1 - Bio-Rad Sadtler	QX
IR - Standards (Subset) 2 - Bio-Rad Sadtler	QFX
IR - Standards (Organic & Polymeric Compounds Subset) - Bio-Rad Sadtler	QPX
Raman - Polymers & Monomers (Basic) - Bio-Rad Sadtler	QRX
Raman - Polymers & Monomers (Basic) 2- Bio-Rad Sadtler	QR2X
IR - Pharmaceuticals - Bio-Rad Sadtler	RX
Raman - Polymers & Processing Chemicals - Bio-Rad Sadtler	RAX

¹¹ B NMR - Wolfgang Robien	RBX
IR - Rubber Chemicals - Bio-Rad Sadtler	RCX
¹⁹ F NMR - Wolfgang Robien	RFX
Raman - Nutraceuticals - Bio-Rad Sadtler	RGNX
Raman - Forensic - HORIBA	RHX
Raman - Inorganics - Bio-Rad Sadtler	RIX
Raman - JASCO	RJX
Raman - Biomaterials - HORIBA	RLX
Raman - Minerals - HORIBA	RMX
¹⁵ N NMR - Wolfgang Robien	RNX
¹⁷ O NMR - Wolfgang Robien	ROX
Raman - Organometallics - Bio-Rad Sadtler	RORX
³¹ P NMR - Wolfgang Robien	RPX
Raman - Standards 1 - Bio-Rad Sadtler	RST1X
Raman - Standards 2 - Bio-Rad Sadtler	RST2X
Raman - Standards 3 - Bio-Rad Sadtler	RST3X
Raman - Standards 4 - Bio-Rad Sadtler	RST4X
Raman - Standards 5 - Bio-Rad Sadtler	RST5X
Raman - Standards 6 - Bio-Rad Sadtler	RST6X
²⁹ Si NMR - Wolfgang Robien	RSX
Raman - Minerals (FT) - HORIBA	RTX
Raman - Semiconductor Materials - HORIBA	RVX
Raman - Controlled & Prescription Drugs - Bio-Rad Sadtler	RZX
Raman - Controlled & Prescription Drugs 2 - Bio-Rad Sadtler	RZ2X
Raman - Controlled & Prescription Drugs 3 - Bio-Rad Sadtler	RZ3X
IR - Steroids 2 - Bio-Rad Sadtler	SX
IR - Standards (Comprehensive) - Bio-Rad Sadtler	SAX
IR - Standards (Selected Subset) - Bio-Rad Sadtler	SLX
IR - Demo Database - Bio-Rad Sadtler	SRX
IR - Standards 1 - Bio-Rad Sadtler	STIR1X
IR - Standards 2 - Bio-Rad Sadtler	STIR2X
IR - Standards 3 - Bio-Rad Sadtler	STIR3X
IR - Standards 4 - Bio-Rad Sadtler	STIR4X
IR - Standards 5 - Bio-Rad Sadtler	STIR5X
ATR-IR - Standards 1 - Bio-Rad Sadtler	STW1X
ATR-IR - Standards 2 - Bio-Rad Sadtler	STW2X
ATR-IR - Standards 3 - Bio-Rad Sadtler	STW3X
ATR-IR - Standards 4 - Bio-Rad Sadtler	STW4X

ATR-IR - Standards 5 - Bio-Rad Sadtler	STW5X
ATR-IR - Standards 6 - Bio-Rad Sadtler	STW6X
SWGDRUG Library – Bio-Rad Sadtler	SWGX
IR - Steroids, Androgens, Progestins, & Estrogens - Bio-Rad Sadtler	SYX
R - Tutorial Database - Bio-Rad Sadtler	TUX
IR - Priority Pollutants (Vapor Phase) - Bio-Rad Sadtler	UGX
IR - Priority Pollutants - Bio-Rad Sadtler	ULX
IR - University Standards - Bio-Rad Sadtler	USX
UV-Vis - 200 to 350 nm - Bio-Rad Sadtler	UW
UV-Vis - 200 to 500 nm - Bio-Rad Sadtler	UX
UV-Vis - 200 to 800 nm - Bio-Rad Sadtler	UY
IR - Standards (Vapor Phase Comprehensive) - Bio-Rad Sadtler	VAX
IR - Standards (Vapor Phase Selected Subset) - Bio-Rad Sadtler	VLX
ATR-IR - Polymers - Bio-Rad Sadtler	WX
ATR-IR - Solvents - Bio-Rad Sadtler	WLX
MS - Wiley Mass Spectral Library of Lipids	WMLX
MS - Wiley Mass Spectra of Pesticides with Retention Indices, 2 nd Edition	WMP3X
MS - Wiley Registry of Mass Spectral Data 11 th Edition	WMS3X
MS - Wiley Registry of Mass Spectra Data 11 th Edition Deleted Removed Spectra	WMSD3X
MS - Wiley Registry of Mass Spectra Data 11 th Edition Replicate Spectra	WMSR3X
¹³ C NMR - Wolfgang Robien	WRX
IR - Water Treatment Chemicals - Bio-Rad Sadtler	WTX
IR - Dyes, Pigments & Stains - Bio-Rad Sadtler	XX
¹⁹ F NMR - Wiley	XF
IR - Dyes - Bio-Rad Sadtler	XLX
¹⁵ N NMR - Wiley	XN
¹⁷ O NMR - Wiley	XO
³¹ P NMR - Wiley	XP
²⁹ Si NMR - Wiley	XS
IR - Inorganics - Bio-Rad Sadtler	YX
IR - Inorganics (Subset) - Bio-Rad Sadtler	YLX
ATR-IR - Inorganics 1 - Bio-Rad Sadtler	YWX
ATR-IR - Inorganics 2 - Bio-Rad Sadtler	YW2X

* Databases accessible at each school depends upon the license the school has chosen.

