

ChemACX Vendor Program

ChemACX Database and Vendor Program

ChemACX is a database of available chemicals that provides a unified, up-to-date, structure-searchable source of chemicals from over 550 Vendors worldwide. The ChemACX Database can be used as a standalone product as well as integrated with CambridgeSoft's Inventory and E-Notebook products. Participation in the ChemACX Vendor Program puts your products in front of qualified chemical buyers and researchers in a range of industries, government laboratories, and academic institutions.

Add your catalog to the ChemACX Database

The Vendor Program allows Vendors to have their catalog appear in the ChemACX database as well as receive communications from CambridgeSoft regarding update schedule, data needs, etc. The program is free. To participate in the ChemACX Vendor Program, we would ask that you send CambridgeSoft your updated catalog information on your regular update schedule (ChemACX is updated quarterly) and include the information outlined in the data required section. We would encourage you to send all of the data types outlined, although we understand that this is not always possible.

We prefer to receive the data in an SDfile containing structures and text data. If you have difficulty creating SDfiles, we can provide products and services to help produce these files. We also accept data in Excel spreadsheets, ChemFinder databases, ISIS/Base databases or tables of Access databases.

Data Required

Required Data	
	Company information
	Vendor Name
	E-mail address for sales inquiries
	Website URL
	Mailing address
	DUNS Number
	Telephone and FAX numbers
	Contact person for this catalog information
	Catalog or collection name
	Catalog or Collection version/update/release number
	Chemical Product Information (for each product)
	Chemical name (or product name, if not a single compound)
	CAS Registry Number, if known
	Catalog number
	UN/CEFACT unit of measure codes for packages (EA = Each, PK = Package, CS = Case, etc).

All specifications subject to change without notice.

US 1 800 315-7300 **INT'L** 1 617 588-9300 **FAX** 1 617 588-9390 **EMAIL** info@cambridgesoft.com

EU 00 800 875 20000 **UK** +44 1223 464900 **JP** 0120 146 700 **WWW** www.cambridgesoft.com

MAIL CambridgeSoft Corporation 100 CambridgePark Drive Cambridge, Massachusetts 02140 USA

¹ ChemBioOffice, ChemBioDraw, ChemBio3D & ChemBioFinder are trademarks of CambridgeSoft Corporation ©2008

CambridgeSoft[®]
www.cambridgesoft.com

ChemACX Vendor Program

Preferred Data			
	Data Type	Preferred Format	Units
	Structure	ChemDraw CDX file	
	MSDS	PDF	
	Package sizes available		
	Prices in all currencies quoted	Dollar and Yen	
	Density (when applicable)	numeric values only	g/ml
	Wt% (i.e. purity)	numeric values only	
	Molarity (when applicable) (Dissolution solvent)	numeric values only	M
	Risk Values	R1, R2– R68	
	Safety Values	S1, S2– S64	
	Melting Point		Degree Celsius
	Boiling Point		Degree Celsius
	Synonyms for chemical name		
	Product description (e.g., purity, concentration, branding information)		

Optional Data			
	Data Type	Preferred Format	Units
	Molecular weight		
	Molecular formula		
	Environmental health and safety data		
	Optical Rotation (when applicable)	positive and negative numeric values only	Degrees

Contact Us

To contact us with questions and comments regarding the ChemACX Database and the Vendor Program or to submit your catalog, please email to ACXcatalogs@cambridgesoft.com. We look forward to hearing from you!

All specifications subject to change without notice.

US 1 800 315-7300 **INT'L** 1 617 588-9300 **FAX** 1 617 588-9390 **EMAIL** info@cambridgesoft.com

EU 00 800 875 20000 **UK** +44 1223 464900 **JP** 0120 146 700 **WWW** www.cambridgesoft.com

MAIL CambridgeSoft Corporation 100 CambridgePark Drive Cambridge, Massachusetts 02140 USA

ChemBioOffice, ChemBioDraw, ChemBio3D & ChemBioFinder are trademarks of CambridgeSoft Corporation ©2008

CambridgeSoft®
www.cambridgesoft.com