

ChemACX.cambridgesoft.com Monthly Update

February 2019

Updated Catalogs

AK SCIENTIFIC, INC.

ALDRICH

ARZA BIOSCIENCE LTD.

BIO FARMA LTD.

CHEMCIA SCIENTIFIC, LLC

MATRIX SCIENTIFIC

PRINCETON BIOMOLECULAR RESEARCH, INC.
(SCREENING COMPOUNDS)

SIGMA ALDRICH LIBRARY OF RARE CHEMICALS

SUPELCO (A SIGMA-ALDRICH BRAND)

SYNCHEM LABORGEMEINSCHAFT OHG

UNITED STATES PHARMACOPEIA (A SIGMA-ALDRICH
BRAND)

New Vendors

AVANTI (A SIGMA -ALDRICH BRAND)

MM (A SIGMA-ALDRICH BRAND)

SIGALD (A SIGMA-ALDRICH BRAND)

All specifications subject to change without notice.

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

EU 00 800 875 20000 UK +44 1223 464900 JP 0120 73 800 WWW www.cambridgesoft.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2018 PerkinElmer Inc. All Rights Reserved.

