

ChemACX Release Statistics

19.19.1 Update Statistics

18 Updated Vendors

FOR COMPLETE LIST PLEASE SEE LIST OF VENDORS
POSTED ON PERKINELMER.COM

4 New Vendors

AMBEED, INC.

CHEMSCENE LLC (BUILDING BLOCKS)

CHEMSCENE LLC (REFERENCE COMPOUNDS)

MEDCHEMEXPRESS (SCREENING LIBRARIES)

19.19.1 Data Information

TOTAL NUMBER OF VENDORS: 748

NUMBER OF PRODUCTS: 23,284,594

NUMBER OF SUBSTANCES: 11,072,574

NUMBER OF SUBSTANCES WITH STRUCTURES:
10,999,002

NUMBER OF PACKAGES: 17,528,896

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.co

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.

