

ChemACX Release Statistics

21.21.3 Update Statistics

23 Updated Vendors

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

1 New Vendor

ASINEX (SCREENING COMPOUNDS)

21.21.3 Data Information

TOTAL NUMBER OF VENDORS: 779

NUMBER OF PRODUCTS: 38,969,931

NUMBER OF SUBSTANCES: 21,262,276

NUMBER OF SUBSTANCES WITH STRUCTURES:
21,178,153

NUMBER OF PACKAGES: 31,864,267

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 1998-2021 PerkinElmer Inc. All Rights Reserved.

