

# ChemACX Release Statistics

## 21.21.2 Update Statistics

### 17 Updated Vendors

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

### 1 New Vendor

AARON CHEMICALS LLC

### 21.21.2 Data Information

TOTAL NUMBER OF VENDORS: 780

NUMBER OF PRODUCTS: 37,417,690

NUMBER OF SUBSTANCES: 20,283,333

NUMBER OF SUBSTANCES WITH STRUCTURES:  
20,199,113

NUMBER OF PACKAGES: 30,154,688

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](mailto:informatics.info@perkinelmer.com)

EU 00 800 875 20000 UK +44 1223 464900 JP 0120 73 800 WWW [www.cambridgesoft.com](http://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.

