



## GET YOUR PRODUCTS NOTICED

### Join the ChemACX Vendor Program

The ChemACX Vendor Program puts your products in front of thousands of qualified chemical buyers and researchers in a range of industries, government laboratories, and academic institutions.

ChemACX is a database of available chemicals that provides a unified, up-to-date, structure-searchable source of chemicals from over 780 vendors worldwide. The database can be used as a standalone product, or can be integrated with our ChemDraw<sup>®</sup>, PerkinElmer Signals<sup>™</sup> Notebook, E-Notebook<sup>™</sup>, and Inventory software.

#### Additional Benefits of the ChemACX Vendor Program

- Your catalog appears in the ChemACX database
- You'll receive communications regarding scheduling, data needs, and more
- Your membership is FREE

To participate, send us your current catalog information on your regular update schedule, and include the following:

#### Required Company Information

- |                   |   |  |
|-------------------|---|--|
| ✓ Vendor name     | ✓ Data Universal Numbering System (DUNS) number | ✓ Catalog or collection name (e.g. building blocks, screening compounds) |
| ✓ Email address   | ✓ Telephone and FAX numbers                     | ✓ Catalog or collection version/ update/release number                   |
| ✓ Website URL     | ✓ Contact person for catalog information        |  |
| ✓ Mailing address |   |  |

#### Required Chemical Product Information

- |   |   |   |
|---|---|---|
| ✓ Chemical name (or product name, if not a single compound) | ✓ Product description   | ✓ Optional: chemical, physical, or biological properties or safety data, if any |
| ✓ Synonyms  | ✓ Catalog number  |   |
| ✓ CAS registry number                                       | ✓ United Nations Standard Products and Services Code (UNSPSC) |   |

#### Required Package Information

- |  |   |
|--|---|
| ✓ Prices for given package sizes in all available currencies | ✓ Available currencies include USD, EUR, GBP, JPY, CHF, CNY |
|--|---|

**Acceptable data formats:** An SDfile containing structures and text data, (we can provide products and services to produce these); Excel spreadsheets; ChemFinder databases; ISIS/Base databases; Access databases

It may also be possible, by special arrangement, to have live links from your product listings on our website to your website. Contact us for more details.

For more information, please email [Informatics.ACXCatalogs@perkinelmer.com](mailto:Informatics.ACXCatalogs@perkinelmer.com).

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)

**PerkinElmer**  
For the Better

For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2019 PerkinElmer, Inc. All rights reserved. PerkinElmer<sup>®</sup> is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.