

# ChemACX Vendor List

**A B C D E F G H I J K L M N O P Q R S T U V W X Y Z**

**-#-**

**1K SCIENTIFIC**

**3A PHARMATECH**

**3B SCIENTIFIC CORP.**

**3RCHEM**

**8S LABORATORIES LLC**

**-A-**

**A&A LIFE SCIENCE INC. DBA  
AVA CHEM SCIENTIFICS**

**A&C PHARMTECH, INC**

**A&J PHARMATECH**

**A.G. SCIENTIFIC**

**A1 BIOCHEM LABS**

Updated Jun. 2021

**A2B CHEM LLC**

Updated May 2021

**AA BLOCKS LLC**

**AB CHEM TECHNOLOGIES LLC**

**ABCR GMBH & Co. KG**

**ABSCO LTD**

**ABX ADVANCED BIOCHEMICAL  
COMPOUNDS**

**AECHEM**

**AK SCIENTIFIC, INC.**

Updated Jun. 2021

**AKL RESEARCH LLP**

**AOB CHEM USA**

**AOKBIO Co., LTD**

**AOKCHEM Co., LTD.**

**APAC PHARMACEUTICAL**

**AQCHEM LLC**

**ASW MEDCHEM, INC.**

**AURUM PHARMATECH LLC**

**AARON CHEMICALS LLC**

Updated Jun. 2021

**AARON CHEMISTRY**

**ABAMACHEM LTD**

**ABBAY COLOR**

**ABBOTT LABORATORIES CHEMICAL  
DIVISION**

**ABBY PHARMATECH LLC**

**ABERJONA LABORATORIES**

**ABOV CHEM LLC**

**ABSOLUTE CHIRAL**

Updated Jul. 2021

**ACADECHEM COMPANY LIMITED**

**ACADEMIQUE PHYTOCANADA**

**ACCELA CHEMBIO Co., LTD.**

Updated Aug. 2020

**ACCELEDEV CHEMICAL LLC**

**ACCUSTANDARD, INC.**

**ACE SYNTHESIS LLC**

**ACENTEX SCIENTIFIC, INC.**

**ACESCHEM INC.**

**ACESYS PHARMATECH**

**ACETO CORPORATION**

**ACHEMICA**

**ACINOPEPTIDE Co. LTD**

**ACME BIOSCIENCE, INC.**

**ACROS ORGANICS**

**ACROS ORGANICS – USA**

**ACTIVATE SCIENTIFIC GMBH**

**ADESIS INC.**

**ADVAMACS**

**ADVANCE RESEARCH CHEMICALS, INC.**

**ADVANCED AROMATICS, L.P.**

**ADVANCED ASYMMETRICS, INC.**

**ADVANCED CHEMBLOCKS INC.**

Updated Jun. 2021

**ADVANCED CHEMTECH**

**ADVANCED MOLECULAR  
TECHNOLOGIES PTY LTD**

**ADVANCED SCIENTIFIC  
INTERNATIONAL, LLC**

**ADVANCED SYNTHESIS  
TECHNOLOGIES, S.A.**

**ADVANCED TECHNOLOGY &  
INDUSTRIAL Co., LTD.**

**AIMS FINE CHEMICALS INC.**

**AIR PRODUCTS & CHEMICALS, INC.**

**AJINOMOTO - AMINO ACID  
DEPARTMENT**

**ALBANY MOLECULAR RESEARCH, INC.**

**ALBEMARLE CORPORATION**

**ALBRIGHT & WILSON AMERICAS**

**ALCHEM INTERNATIONAL LIMITED**

**ALCHEM PHARMTECH, INC.**

**ALDLAB CHEMICALS, LLC**

**ALDRICH**

**ALFA AESAR**

**ALFA CHEMISTRY**

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 800

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

## ALFA CHEMISTRY (MATERIALS)

New Aug. 2020

## ALFA PYRIDINES

## ALICHEM INC.

Updated May. 2021

## ALLIEDSIGNAL INC. SPECIALTY CHEMICALS

## AMBEED, INC.

Updated Jun. 2021

## AMERICAN ADVANCED ORGANICS (DIV OF ALBANY MOLECULAR)

## AMERICAN ADVANCED SCIENTIFIC INC.

## AMERICAN BIOANALYTICAL

## AMERICAN BIOCHEMICALS

## AMERICAN CUSTOM CHEMICAL CORP

## AMERICAN PEPTIDE COMPANY

## AMERICAN RADIOLABELED CHEMICALS, INC.

## AMERICHEM, INC.

## AMRESCO INC.

## ANGEL PHARMATECH, LTD.

## ANGELINI FINE CHEMICALS

## ANGENE INTERNATIONAL LIMITED

## ANGUS CHEMICAL COMPANY

## ANNKER ORGANICS

## ANTIBIOTICOS SPA

## ANWARD CHEMICALS

## APEIRON SYNTHESIS

## APEX SCIENTIFIC INC.

## APEXMOL TECHNOLOGY

## APITOXIN

## APOLLO SCIENTIFIC LTD. (BIOCHEMICALS)

## APOLLO SCIENTIFIC LTD. (FLUORO CHEMICALS)

## APOLLO SCIENTIFIC LTD. (INTERMEDIATES)

## APOLLO SCIENTIFIC LTD. (SUPPORT RESEARCH LIBRARY)

## APPLICHEM

## ARBOŃOVA OY AB

## ARCTOM CHEMICALS

## ARK PHARM, INC.

## ARKÉ ORGANICS SRL

## AROMA CIRCLE

## ARONIS

## ARRAN CHEMICAL CO.

## ARYLCHEM CO.

## ARZA BIOSCIENCE LIMITED

Updated Jul. 2021

## ASAHI GLASS COMPANY - FLUORO CHEMICALS DIVISION

## ASCENT SCIENTIFIC

## ASHWOOD CHEMICALS LTD.

## ASINEX LTD. (SCREENING COMPOUNDS)

New Jul. 2021

## ASISCHEM INC.

## ASTATECH, INC.

## ASTAR PHARMACEUTICALS, LLC

## ATOMOLE SCIENTIFIC CO., LTD.

## AUGUSTUS OILS LTD

## AURORA FINE CHEMICALS LTD

Updated Feb. 2021

## AVANTI (A SIGMA-ALDRICH BRAND)

## AVIABOR FINE BORON COMPOUNDS

## AVONCHEM/CHROMOS EXPRESS LTD

## AXISPHARM

New Aug. 2020

## AXITEC CO., LTD

## AZEPINE LTD

## -B-

## BACHEM AG

## BASF CHEMICAL INTERMEDIATES

## BE-ASSOCIATES, LLC

## BLD PHARMATECH LTD

Updated Jan. 2021

## BOC SCIENCES

## BSC CHEMICALS

## BAIHUA BIO-PHARMACEUTICAL CO. LTD.

## BALAJI AMINES LTD

## BAYER CORPORATION

## BAYER INORGANICS

## BEDOUKIAN RESEARCH, INC.

## BEIJING CENTURY RICHAP CHEMISTRY CO. LTD

## BEIJING GOLDEN OLIVE COMPANY

## BEPHARM LTD.

## BESTIPHARMA

## BETA PHARMA, INC.

## BEWE LEEPHARMA INC.

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

BIDDLE SAWYER CORPORATION

BIDE PHARMATECH

BIO FARMA LTD.

BIO-RESEARCH PRODUCTS, INC.

BIOBLOCKS, INC.

BIOQUADRANT INC.

BIOVISION, INC.

BIOCHEMIE GMBH

BIOCHEMPEG

BIOCORE PHARMTECH LIMITED

BIOFINE INTERNATIONAL INC.

Updated Feb. 2021

BIOMOL

BIOSYNTH AG

BIOTIUM, INC.

BIOTREND CHEMIKALIEN GMBH  
(NEUROCHEMICALS)

BIOTREND CHEMIKALIEN GMBH  
(RADIOCHEMICALS)

BOERCHEM PHARMTECH CO., LTD

BOROCHEM

BORON MOLECULAR

BOSCHE SCIENTIFIC LLC

BOULDER SCIENTIFIC COMPANY

BRIDGE ORGANICS CO.

BRIGHT EVERGREEN

BROADPHARM

BURDICK & JACKSON, INC.

BUSH BOAKE ALLEN INC.

BUTT PARK LTD

**-C-**

C/D/N ISOTOPES INC.

CF PLUS CHEMICALS S.R.O.

CGENETECH, INC.

CMS CHEMICALS LTD --  
CARBOHYDRATES

CMS CHEMICALS LTD -- ORGANICS

CNH TECHNOLOGIES, INC.

CU CHEMIE UETIKON

CACHESYN INC.

CALAIRE CHIMIE S.A.

CALLERY CHEMICAL COMPANY

CAMBREX CORPORATION

CAMBRIDGE ISOTOPE LABS

CAMBRIDGE MAJOR LABORATORIES,  
INC.

CANCERPHARM GROUP

CAPOT CHEMICAL COMPANY LIMITED

CARBOMER INC.

CARBOSYNTH LTD.

CAYMAN CHEMICAL COMPANY

CELANESE, LTD.

CENTEC INTERNATIONAL

CENTRAL GLASS CO. LTD.- FINE  
CHEMICAL DEPARTMENT

CERILLIANT CORP.

CHANGZHOU SHUOCHENG  
BIOLOGICAL TECHNOLOGY CO. LTD

CHANGZHOU WELTON CHEMICAL  
CO.,LTD.

CHARKIT CHEMICAL CORPORATION  
CHATTEM CHEMICALS, INC.

CHEM-IMPEX INTERNATIONAL, INC.

CHEMALONG LABS LLC

CHEMBASICS PTY LTD

CHEMBRIDGE CORPORATION  
(BUILDING BLOCKS)

Updated Feb. 2021

CHEMBRIDGE CORPORATION  
(SCREENING COMPOUNDS)

Updated May. 2021

CHEMCONTRACT RESEARCH INC.

CHEMDESIGN CORP. (A BAYER  
COMPANY)

CHEMDIV, INC.

Updated Jun. 2021

CHEMDIV (DISCOVERY CHEMICALS)

Updated Jul. 2021

CHEMFOCUS, LLC

CHEMPACIFIC CORPORATION

CHEMSAMP Co, LLC

CHEMSERVICE, INC.

CHEMADA

CHEMATEK S.P.A.

CHEMBO PHARMA Co LTD

CHEMCIA SCIENTIFIC, LLC

CHEMFUTURE PHARMATECH

CHEMGENX LLC

CHEMI S.P.A.

CHEMICAL TECHNOLOGIES &  
INVESTIGATIONS, LTD.

CHEMICREA INC.

CHEMIELIVA PHARMACEUTICAL CO.,  
LTD

Updated Sep. 2020

CHEMINISCHE LABORTORIEN  
DR. SOENKE PETERSEN

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

CHEMISPHERE LIMITED

CHEMIVATE LIMITED

CHEMIZON INC.

CHEMODEX LTD (BUILDING BLOCKS)

CHEMODEX LTD (SCREENING COMPOUNDS)

CHEMSCENE LLC (BUILDING BLOCKS)

CHEMSCENE LLC (REFERENCE COMPOUNDS)

CHEMTEC LEUNA - SPECIALTY CHEMICALS AND FINE CHEMICALS

CHEMTOS, LLC

CHEMTURA CORPORATION

CHEMZON SCIENTIFIC

CHENGDU GT BIOCHEM Co.,LTD.

CHESS GMBH

CHEVRON PHILLIPS CHEMICAL COMPANY LLC

CHIREX LTD.

CHINA NATIONAL CHEMICAL CONSTRUCTION CORPORATION

CHIRACON GMBH

CHIRAGENE CORPORATION (FORMERLY CELGENE)

CHIRALSTAR INC

CHIROBLOCK GMBH

CHIROTECH TECHNOLOGY LIMITED

CHIRON AS

Updated Oct. 2020

CHONGQING JIULONGPO YUHUA CHEMICALS Co., LTD.

CHUGAI KASEI Co., LTD.

CILAG AG

CITY CHEMICAL LLC

CIVENTI CHEM

CLARIANT CORPORATION

CLASSIC FLAVORS & FRAGRANCES

CLEARSYNTH INC.

COALITE CHEMICALS

COMBI-BLOCKS, INC.

COMBIPHOS CATALYSTS, INC.

COOL PHARM, LTD.

CORVINUS CHEMICALS

CREANOVA, INC. (FORMERLY HÜLS AMERICA, INC.)

CUSCHEM INC.

CYMTEK LABORATORIES, LIMITED

CYTH CHEMICAL TRADE Co., LTD.

**-D-**

D&O CHEMICALS, INC.

D.F. GOLDSMITH CHEMICALS & METALS Co.

DC CHEMICALS

New Aug. 2020

DSM FINE CHEMICALS INC.

DAICEL CHEMICAL INDUSTRIES, LTD.

DAIKIN INDUSTRIES LTD

DAISO Co., LTD.

DALTON PHARMA SERVICES INC.

DANNIER CHEMICAL, INC.

DANYANG HENGAN CHEMICAL Co., LTD.

DAVOS CHEMICAL CORP.

DAYBIOCHEM TECHNOLOGY Co., LTD  
DE VOIT ASSOCIATES

DEBYE SCIENTIFIC

Updated Jun. 2021

DEEPAK CHEMICALS GROUP

DEEPWATER CHEMICALS

DEGUSSA CORP.

DHRUVTARA CHEMICALS PVT LTD

DIAGNOSTIC CHEMICALS LIMITED

DIGITAL SPECIALTY CHEMICALS, INC.

DINAMITE DIPHARMA SPA

DIOSYNTH B.V.

DISCOVERY MOLECULES

DISHMAN USA

DIVERCHIM

DIXIE CHEMICAL COMPANY, INC.

DOTTIKON EXCLUSIVE SYNTHESIS AG

DYSTAR LP

DYNAMIT NOBEL GMBH

**-E-**

ECA INTERNATIONAL CORPORATION

EMD CHEMICALS INC.

EMKA-CHEMIE

EASTAR CHEMICAL CORPORATION

EASTMAN CHEMICAL COMPANY

ECHELON BIOSCIENCES INC.

ELAN CHEMICAL COMPANY, INC.

ELECTRON MICROSCOPY SERVICES

ELEMENT STORE

ELKIMIA INC.

ELPAR INTERNATIONAL INC.

ENAMINE LTD.

Updated Jul. 2021

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

ENDOTHERM GMBH

ENNOVA MEDCHEM GROUP, INC.

E NOVATION CHEMICALS LLC

ERNESTO VENTOS S.A.

ERREGIERRE INDUSTRIA CHIMICA  
S.P.A

ESPRIT CHEMICAL COMPANY

ESSTECH, INC.

EUROLABS LIMITED

EUROPEAN TECHNOLOGIES INC.

EVOBLOCKS

EXCENEN PHARMATECH (BIOCHEMICALS)  
Updated Jan 2021

EXCHEM ORGANICS

EXIMED LABORATORY

EXCLUSIVE CHEMISTRY LTD

EXPANSIA S.A.

**-F-**

F2 CHEMICALS LTD.

FCH GROUP

FTECH, INC.

FABBRICA ITALIANA SINTETICI

FAR RESEARCH

FARCHEMIA DIVISION OF  
TESSENDERLO CHEMIE

FARMAK, A.S.

FINE CHEMICAL Co., LTD. LANZHOU  
BELFAST

FINE CHEMICALS CORPORATION  
(PTY) LTD.

FINE ORGANICS LTD., SEAL SANDS

FINETECH INDUSTRY LTD

FINORNIC CHEMICALS (I) PVT. LTD.

FIRMENICH SA

FIRST CHEMICAL CORP.

FISHER SCIENTIFIC (U.S.)

FISHER SCIENTIFIC UK

FLEXCHEM BV

FLUKA CHEMICAL CORP.

FLUOROCEM GROUP -- SCALE UP  
SOLUTIONS

FLUOROCEM UK

FLUOROUS TECHNOLOGIES, INC.

FOCUS SYNTHESIS LLC

FRAMOCEM (CHAYSECHEM, NA  
DISTRIBUTOR)

FRINTON LABORATORIES

FRONTIER SCIENTIFIC SERVICES INC.  
Updated Apr. 2021

FRONTIER SCIENTIFIC, INC.

FRUTAROM (LABORATORY CHEMICALS)

FRUTAROM LTD (AROMA CHEMICALS)

FUJI HUNT SPECIALTY PRODUCTS

**-G-**

G&J RESEARCH CHEMICALS LTD.

G.J. CHEMICAL COMPANY, INC.

GFS CHEMICALS, INC.

GNI CHEMICALS CORPORATION

GANES CHEMICALS INC.

GAYLORD CHEMICAL CORPORATION

GELEST INC.

GENERAL INTERMEDIATES OF

CANADA

GENERICHEM CORPORATION

GENOSYS (A SIGMA ALDRICH BRAND)

GENZYME PHARMACEUTICALS

GEORGE UHE COMPANY, INC.

GREATFIELD MINERAL AND  
CHEMICAL INC.

GREENCHEM USA

GUANGZHOU TOPWORK  
INTERNATIONAL CO. LTD.

GWALIOR CHEMICAL INDUSTRIES  
LTD.

**-H-**

H&S CHEMICAL CO. INC.

HAISO PHARMCHEM

HALOCARBON PRODUCTS CORP.

HANDE SCIENCES US

HANGZHOU APICHEM TECHNOLOGY  
Co., LTD.  
Updated Aug. 2020

HANGZHOU PHARMA & CHEM. CO. LTD

HANGZHOU TRYLEAD CHEMICAL  
TECHNOLOGY Co., LTD

HANGZHOU YUHAO CHEMICAL  
TECHNOLOGY Co., LTD

HANSHAN AGILEBIOCHEM

HETCAT

HETEROCYCLICS INC.

HEXAGON LABS INC.

HIGH PERFORMANCE PRODUCT  
ENGINEERING

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

**HOKKO CHEMICAL INDUSTRY CO., LTD.**  
**- FINE CHEMICALS DIVISION**

**HOVIONE GROUP**

**HUHUGENE INC.**

## **-I-**

**IAN HETEROCYCLE CO., LTD.**

**ISP FINE CHEMICALS (FORMERLY GAF)**

**IVD (A SIGMA-ALDRICH BRAND)**

**ICON SERVICES**

**IHARA CHEMICAL INDUSTRY CO., LTD.**

**IHARA NIKKEI CHEMICAL INDUSTRY CO., LTD.**

**INFARMATIK, INC.**

**INABATA AMERICA CORP.**

**INDOFINE CHEMICAL COMPANY, INC.**

**INDUSTRIA CHIMICA MILANESE S.P.A.**

**INNOCEHEM CHEMICAL CO., LTD.**

**INSIGHT CHEMICAL SOLUTIONS**

**INTERBIOSCREEN LTD.**  
**(BIOLOGICALS)**

**INTERBIOSCREEN LTD. (BUILDING BLOCKS)**

**INTERBIOSCREEN LTD (SCREENING COMPOUNDS)**

**INTERCHEM CORP.**

**INTERNATIONAL FLAVORS & FRAGRANCES INC.**

**IOLiTEC GMBH & Co KG**

**IRIS BIOTECH GMBH**

**ISOCHEM, (CHAYSECHEM, NA DISTRIBUTOR)**

**ISOTEC INC.**

## **-J-**

**J&K SCIENTIFIC LTD**

**J&W PHARMLAB, LLC**  
Updated May 2021

**J.P. PHARMA-CHEM INDUSTRIES**

**J.H. CALO COMPANY**

**J.T. BAKER**

**JPM2 PHARMACEUTICALS, INC.**

**JRD FLUOROCEMICALS LTD**  
Updated Sep. 2020

**JIANGSU INSTITUTE OF ECOMONES**

**JIANGSU ZHONGDAN CHEMICAL IMP. & EXP. CO., LTD.**

**JOST CHEMICAL**

**JUNSEI CHEMICAL CO., LTD**

**JUPITER BIOSCIENCES LTD**

**JUPITER CHEMICALS, INC.**

## **-K-**

**K-I CHEMICAL INDUSTRY CO., LTD.**

**KADEN BIOCHEMICALS GMBH**

**KANEKA CORPORATION - FINE CHEMICALS DIVISION**

**KANTO CHEMICAL CO., INC.**

**KATWIJK CHEMIE BV**

**KAIRONKEM**

**KEMIRA FINE CHEMICALS**

**KEY MOLECULAR CORPORATION**

**KEY ORGANICS BIONET (BIOCHEMICALS)**

**KEY ORGANICS BIONET (BUILDING BLOCKS)**  
Updated Jun. 2021

**KEY ORGANICS BIONET (COMBINED-SCREENING-FRAGMENTS)**  
Updated Jul. 2021

**KEY ORGANICS BIONET (FRAGMENT LIBRARY)**  
Updated May 2021

**KEYINTERMEDIATE**

**KINENTIA BIOSCIENCES**

**KINGCHEM INC.**

**KINGSCIENTIFIC**

**KINGSH CHEMICALS INC.**

**KISHIDA BUILDING BLOCKS**

**KISHIDA CHEMICAL CO., LTD**

**KLAUS F. MEYER GMBH**

**KOEI CHEMICAL COMPANY, LTD**

**KYOWA HAKKO KOGYO CO., LTD.**

## **-L-**

**LC LABORATORIES**

**LGM PHARMA**

**LKT LABORATORIES, INC.**

**LABNETWORK**

**LABORATORI MAG SPA**

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588- FAX 1 617 588- EMAIL  
EU 00 800 875 UK +44 1223 JP 0120 73 800 informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

**LAN PHARMATECH LTD.**

Updated Apr. 2021

**LANCASTER SYNTHESIS INC.**

**LANSDOWNE CHEMICALS PLC**

**LANXESS AG**

**LANZHOU TECH BIOCHEMICAL CO. LTD.**

**LAPORTE PERFORMANCE CHEMICALS (FORMERLY INSPEC GROUP)**

**LATOXAN**

**LEADGEN LABS, LLC**

**LIFE CHEMICALS INC. (BUILDING BLOCKS)**

**LIFE CHEMICALS INC. (SCREENING COMPOUNDS)**

**LIVCHEM LOGISTICS GMBH**

**LOBA FEINCHEMIE AG**

**LONZA LTD.**

**LUKNOVA INC**

**LUMAC\*LSC B.V.**

**LYONDELL CHEMICAL COMPANY**

**-M-**

**MM (A SIGMA-ALDRICH BRAND)**

**MP BIOMEDICALS**

**MV LABORATORIES, INC.**

**MACFARLAN SMITH**

**MALLINCKRODT LABORATORY CHEMICALS**

**MANCHESTER ORGANICS LTD**

Updated Mar. 2021

**MARKER GENE TECHNOLOGIES, INC**

**MARSHALLTON RESEARCH LABORATORIES, INC.**

**MATERIA INC.**

**MATRIX SCIENTIFIC**

**MAYBRIDGE CHEMICAL COMPANY, LTD.**

**MCULE**

Updated Nov. 2020

**MCULE ULTIMATE**

Updated Jun. 2021

**MEDCHEMEXPRESS LLC (SCREENING LIBRARIES)**

**MEDKOO BIOSCIENCES**

**MEDICAL ISOTOPES, INC.**

**MEGAZYME INTERNATIONAL**

**MELFORD LABORATORIES LTD.**

**MENAI ORGANICS**

**MERCK SCHUCHARDT OHG**

**METALL-CHEMIE GOERRIG GMBH & CO. KG**

**MICROCOMBICHEM E.K.**

**MILESTONE PHARMTECH LLC**

**MILLIONLINK (TIANJIN) INTERNATIONAL TRADE CO., LTD.**

**MINOL ACIDS & CHEMICALS PVT. LTD.**

**MITCHELL COTTS CHEMICALS**

**MITENI SPA**

**MITSUBISHI RAYON CO., LTD.**

**MITSUI CHEMICALS, INC. MOLBRIDGE**

**MOLE CHEMICAL CORPORATION**

**MOLECULAR BIOSCIENCES**

**MOLECULAR PROBES, INC.**

**MOLEKULA**

**MOLEPEDIA Co., LTD.**

**MORAVEK BIOCHEMICALS**

**MORFLEX, INC.**

**MORRE-TEC INDUSTRIES, INC.**

**MORTON INTERNATIONAL, INC.**

**MUSECHEM**

Updated Jan. 2021

**-N-**

**NSC TECHNOLOGIES**

**NACALAI TESQUE INC**

**NAMIKI-SHOJI CO., LTD.**

Updated Sep. 2020

**NANJING DAXIAN CHEMICAL INSTITUTE LTD.**

**NANJING PHARMACEUTICAL FACTORY CO., LTD.**

**NANJING RAYMON BIOTECH Co., LTD**

**NANJING REDLINE LTD.**

**NEOMPS, INC.**

**NEOSYN LABORATORIES INC.**

**NETQEM**

**NEW ZEALAND PHARMACEUTICALS LIMITED**

**NICHIMEN AMERICA INC.**

**NIGU CHEMIE GMBH**

**NIHON NOHYAKU Co., LTD.**

**NIPA HARDWICKE INC.**

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 800

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights



# ChemACX Vendor List

NIPPOH CHEMICALS CO., LTD.

NIPPON SHOKUBAI CO. LTD.

NORCHIM S.A.

NORQUAY TECHNOLOGY INC. USA

NORTHEAST PHARMACEUTICAL  
GROUP CO. LTD

NOVABIOCHEM (EMD BIOSCIENCES,  
INC.)

NOVASYN ORGANICS PVT. LTD.

NOVELKORE

NOVOCHEMY LTD.

**-O-**

OAKWOOD PRODUCTS, INC.

OCCIDENTAL PETROLEUM  
CORPORATION

OMEGACHEM, INC.

ONTARIO PHARMA RESEARCH  
CHEMICALS

ONYX SCIENTIFIC LIMITED

OPTIMA CHEMICAL GROUP LLC

ORGANICA FEINCHEMIE GMBH  
WOLFEN

ORGANIX, INC.

ORGANOCEM LTD

ORGAPHARM-AXYNTIS GROUP

ORGASYNTH

ORIBASE PHARMA

ORION CORPORATION FERMION

OSAKA ORGANIC CHEMICAL INDUSTRY,  
LTD.

OTAVA

OXCHEM CORPORATION

**-P-**

PFL FINE CHEMICALS GMBH

PK CHEM

PNK UNITED, INC.

PPG-SIPSY

PALINDENT ORGANIC  
INTERMEDIATES

PANCHIM

PARADIGM SCIENTIFIC

PARISH CHEMICAL COMPANY

PARKWAY SCIENTIFIC LLC

PEAKDALE FINE CHEMICALS LIMITED

PENG TENG SHANGHAI FINE CHEMICAL  
CO.

PENN SPECIALTY CHEMICALS, INC.

PENTAGON CHEMICALS

PEPTeCH CORPORATION

PERKINELMER CHEMICALS

PFALTZ & BAUER, INC.

PFANSTIEHL LABORATORIES, INC.

PHARMABLOCK R&D Co. LTD.

PHARMACORE, INC.

PHARMABRIDGE

PHARMACEUTICAL WORKS POLPHARMA  
S.A.

PHARMACIA & UPJOHN COMPANY

PHARMCO PRODUCTS INC.

PHARMTEN CHEMICAL Co.,  
LIMITED

PHOENIX CHEMICALS

PHYTOLAB GMBH & Co. KG

PLATTE VALLEY SCIENTIFIC, LLC

POLI INDUSTRIA CHIMICA S.P.A.

POLIUM TECHNOLOGIES, INC.

POLYCARBON INDS., INC.

POLYMER LABORATORIES LTD

POLYORGANIX, INC.

POLYSCIENCES, INC.

PORSE FINE CHEMICAL Co., LTD.

PRATAP ORGANICS

PRECISION BIOCHEMICALS, INC.

PRECURSOR CHEMICALS, INC.

PRINCETON BIOMOLECULAR  
RESEARCH, INC. (BUILDING BLOCKS)

PRINCETON BIOMOLECULAR  
RESEARCH, INC. (SCREENING  
COMPOUNDS)

Updated Oct. 2020

PROCHEM, INC.

PRODOTTI CHIMICI E ALIMENTARI  
SPA

PURE CHEMISTRY SCIENTIFIC INC.

**-Q-**

QUANTA BIODESIGN, LTD.

QUEEN'S UNIVERSITY OF BELFAST

**-R-**

R&S CHEMICALS, INC.

R.I. CHEMICAL, INC.

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights



# ChemACX Vendor List

RCA KG (REUTER CHEMISCHE  
APPARATEBAU KG)

RSP AMINO ACID ANALOGUES, INC.

RARE CHEMICALS GMBH  
Updated Oct. 2020

RARE CHEMICALS GMBH (SCREENING  
LIBRARY)

RASCHIG AG

REAXON GBR

REILLY INDUSTRIES, INC.

RELAX LTD.

RELIABLE BIOPHARMACEUTICAL  
CORPORATION

RESEARCH ORGANICS

RHODIA FINE ORGANICS

RICHMAN CHEMICAL INC.

RIEDEL-DE HAËN

RIEKE METALS, INC.  
Updated Jun. 2021

RIEKE METALS, INC. (NOVEL  
CHEMICAL SOLUTIONS)

ROBINSON BROTHERS LIMITED

ROCHE

ROCKWOOD LITHIUM GMBH

ROHM & HAAS PROCESS CHEMICALS

ROHNER LTD.

ROMIL LTD

RUETGERS-NEASE CORPORATION

RYAN SCIENTIFIC, INC.

RYSCOR SCIENCE, INC.

**-S-**

S.I.M.S. S.R.L.

S.K. CHEMICAL INDUSTRIES

SEAC

SK ENERGY AND CHEMICAL, INC.

SKW CHEMICALS

SNPE CHEMICALS, GROUPE SNPE

SALFORD ULTRAFINE CHEMICALS &  
RESEARCH LTD.

SAMCHULLY PHARMACY COMPANY,  
LTD.

SANFUR USA, INC.

SANOCHEMIA PHARMAZEUTIKA AG

SANTAI LABS

SANTIAGO

SAUDI BASIC INDUSTRIES  
CORPORATION

SCHERING AG

SELENA CHEMICALS

SELLECKCHEM

SENN CHEMICALS

SENSACHEM, INC.

SHAANXI BONO BIOTECH Co., LTD

SEQUOIA RESEARCH PRODUCTS LTD

SHANDONG YOUBANG BIOCHEMICAL  
TECHNOLOGY Co., LTD.

SHANGHAI ACT CHEMICAL LIMITED

SHANGHAI AOBO BIO-PHARMACEUTICAL  
TECHNOLOGY Co., LTD.

SHANGHAI CHEMHERE Co., LTD

SHANGHAI ELITTES ORGANICS Co., LTD.

SHANGHAI EXCELLENT CHEMICAL Co., LTD.

SHANGHAI FWD CHEMICALS Co., LTD.

SHANGHAI HAUYUAN CHEMEXPRESS  
Co., LTD

SHANGHAI INTERNATIONAL  
PHARMACEUTICAL Co., LTD

SHANGHAI PI CHEMICALS LTD.

SHANGHAI PHARMA TECH Co., LTD.  
(PARLING)

SHANGHAI PU-JIE FRAGRANCE Co.,  
LTD.

SHANGHAI RACE CHEMICAL Co., LTD

SHANGHAI REITO CHEMICAL  
TECHNOLOGY Co., LTD.

SHANGHAI RONGKANG FINE  
CHEMICAL Co., LTD

SHANGHAI SINOFLUORO SCIENTIFIC  
CORPORATION LTD.

SHANGHAI SPECBIOCHEM CHEMISTRY  
GROUP LIMITED

SHELL CHEMICAL LP

SICHUAN TONGSHENG AMINO ACID  
Co., LTD

SIGALD (A SIGMA-ALDRICH BRAND)

SIGMA ALDRICH DISCOVERY  
CHEMICALS

SIGMA CHEMICAL COMPANY

SIGMA-ALDRICH BRAND CHEMICALS

SIGMA-ALDRICH JAPAN K.K.

SILAR LABORATORIES - A DIVISION OF  
WRIGHT CORPORATION

SIMAFEX

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 80C

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

SINOCHEM HEBEI TIANJIN IMPORT - EXPORT CORPORATION	SUZHOU SINOERA CHEM CO., LTD	SÄUREFABRIK SCHWEIZERHALL (SF-CHEM)
SINON CORPORATION	SUZHOU YIELD PHARMACEUTICAL R&D Co., LTD	-T-
SMALL MOLECULES	SYNAFFIX BV	
SOHVE CORPORATION	SYNCHEM, INC.	TC CHEMICALS Co., LTD.
SOLTEC VENTURES, INC.	SYNFINE RESEARCH	TOSLAB
SPECS (BUILDING BLOCKS)	SYNHET INC. Updated Apr. 2021	TAIZHOU KANGDUO SEA BIOCHEMICAL & PHARMACEUTICAL Co., LTD.
SPECS (NATURAL PRODUCTS)	SYNNNOVA	TAKASAGO INTERNATIONAL CORP.
SPECS (SCREENING COMPOUNDS)	SYNQUEST LABORATORIES, INC. Updated Jun. 2021	TAOKA CHEMICAL Co., LTD.
SPECTRUM CHEMICAL & LABORATORY PRODUCTS	SYNCHEM LABORGEMEINSCHAFT UG & Co. KG	TECOLAND CORPORATION
SPECTRUM INFO LTD.	SYNCOZYMES (SHANGHAI) Co., LTD.	TENOVA PHARMACEUTICALS INC
SPHEROTECH, INC.	SYNERGY-SCIENTIFIC LTD.	TEVA GROUP API DIVISION
SPIROCHEM AG	SYNGANIX SCIENCE INC.	TEXAS BIOCHEMICALS INC.
STANFORD MATERIALS COMPANY	SYNOVEL LABORATORY Co., LTD.	THE NEW CENTURY CHEMICAL & PHARMACEUTICAL ENTERPRISE
STREM CHEMICALS, INC.	SYNOVO GMBH	THE SHEPHERD CHEMICAL COMPANY
SUMIKIN CHEMICAL Co. LTD.	SYNTAI CHEMICALS & PHARMACEUTICALS, LTD.	TOCRIS BIOSCIENCE
SUMITOMO SEIKA CHEMICALS Co., LTD.	SYNTASTIC AB	TOKUYAMA CORPORATION
SUN BIOCHEMICALS, INC.	SYNTECH SOLUTION	TOKYO CHEMICAL INDUSTRY Co., LTD.
SUNOCO CHEMICALS	SYNTESIA CHEMIE GMBH	TOMAH PRODUCTS, INC.
SUNSHINE CHEMICALS COMPANY	SYNTHE163	TORAY INDUSTRIES, INC.
SUNTTO CHEMICAL	SYNTHETECH	TORONTO RESEARCH CHEMICALS, INC.
SUPELCO (A SIGMA ALDRICH BRAND)	SYNTHON CHEMICALS GMBH & Co. KG	TOSOH CORPORATION
SUPRIYA CHEMICALS	SYNTHON CHIRAGENICS CORPORATION	TRACTUS COMPANY LIMITED
SUZHOU CHONTECH PHARMACHEM TECHNOLOGY Co., LTD.	SYNTHONIX CORPORATION Updated Jun. 2021	TRANS WORLD CHEMICALS INC.
SUZHOU ROVATHIN PHARMA TECH Co., LTD.	SYNTIDE MEDI-CHEM TECHNOLOGY Co., LTD.	TRATAMIENTOS GUADALQUIVIR S.A. (TRAGUSA)
SUZHOU RYAN PHARMACHEM TECHNOLOGY Co., LTD.		

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588- FAX 1 617 588- EMAIL  
EU 00 800 875 UK +44 1223 JP 0120 73 800 informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights

# ChemACX Vendor List

TRIFARMA SPA

TRINAPCO, INC.

TRIOMPHE FINE CHEMISTRY CO.

TRIPLEBOND CORP.

TYGER SCIENTIFIC INC.

## -U-

USDAC COMPANY

UBE INDUSTRIES, LTD. - FINE  
CHEMICALS DIVISION

UBICHEM PLC

UNAVERA CHEMLAB GMBH

UNITED STATES PHARMA COPEIA (A  
SIGMA ALDRICH BRAND)

## -V-

VWR INTERNATIONAL GMBH

VWR INTERNATIONAL, LLC

VANDEMARK GROUP

VARSAL INSTRUMENTS, INC.

VESINO INDUSTRIAL CO.

VETEC QUIMICA FINA LTDA

VIJAYA PHARMACEUTICALS

VIS FARMACEUTICI SPA

VITAS-M LABORATORY, LTD.  
Updated Aug. 2020

VITAS-M LABORATORY, LTD.  
(SCREENING COMPOUNDS)  
Updated Apr. 2021

VIVA CORPORATION

VIWIT PHARMATECH CO., LTD.

## -W-

WACKER CHEMICALS (USA)

WAKO PURE CHEMICAL INDUSTRIES,  
LTD.

WATANABE CHEMICAL INDUSTRIES,  
LTD.  
Updated Jun. 2021

WATEC LABORATORIES, INC.

WATERSTONE TECHNOLOGY, LLC

WHYTE CHEMICALS LIMITED

WILLIAM BLYTHE LTD.

WILSHIRE CHEMICAL CO., INC.

WINNER GROUP INTERNATIONAL LTD.

WONDA SCIENCE INC.

WONDERCHEM Co., LIMITED

WUXI APPTec Co., LTD

WUHAN ZHONGFU BIO-TECH Co.,  
LTD.

WUJIN LINCHUAN CHEMICAL  
FACTORY

WUXI D-STONE INTERNATIONAL Co.,  
LTD.

WYCHEM LIMITED

## -X-

XPR INDUSTRIAL (CHINA) LIMITED

XiMO HUNGARY, LTD.

XINCHEM COMPANY

## -Y-

YAMAKAWA CHEMICAL

YUANHONG SINO CHEMICAL

YUHAN CORPORATION

## -Z-

ZAMBON GROUP

ZENECA SPECIALTIES

ZEON CORP

ZERENEX BIOACTIVES

ZERENEX MOLECULAR LIMITED  
Updated Mar. 2021

ZHEJIANG KANGYU PHARMACEUTICAL  
CO.,LTD

ZHEJIANG OUHUA CHEMICAL  
IMP.&EXP.CO.,LTD

ZHENGTAI PHARMA & CHEM Co., LTD.

ZYLEXA PHARMA LTD

All specifications subject to change

US 1 800 315-7300 INT'L 1 617 588-

EU 00 800 875 UK +44 1223

FAX 1 617 588-

JP 0120 73 800

EMAIL

informatics.info@perkinelmer.com

MAIL PerkinElmer 940 Winter St. Waltham, MA 02451 USA

ChemBioOffice, ChemBioDraw, Chem3D, ChemBioFinder & ChemInfo Copyright 1998-2021 PerkinElmer Inc. All Rights